

Maurice, Maurice C., Maurice S. and MauriceFlex

User Guide for Maurice, Maurice C., Maurice S. and MauriceFlex

Copyright © 2023. All Rights Reserved.

PN 046-295

Revision D, March 2023

For research use only. Not for use in diagnostic procedures

Patents and Trademarks

ProteinSimple's Maurice and iCE technology is covered by issued and pending patents in the U.S. and other countries. For more information please see <https://www.bio-techne.com/support/instrument-support/intellectual-property>.

ProteinSimple and the ProteinSimple logo are trademarks or registered trademarks of ProteinSimple. Other marks appearing in these materials are marks of their respective owners.

Table of Contents

Chapter 1: Let's Get Started 16

Welcome	17
Maurice Systems	17

Chapter 2: Getting Your Lab Ready 18

Introduction	19
Space Requirements.....	19
Physical Specifications	19
Electrical Requirements	20
Environmental Requirements.....	20
Software and Computer Requirements	20
Using Maurice with Compass for iCE.....	20
Using Maurice with Empower®.....	21
General Guidelines and Information	21
Intended Use	21
Lifting and Moving the System: Lift Maurice Correctly	21

Chapter 3: Maurice 22

Maurice Systems	23
External Components.....	24
System Door	24
Status Light	24
Internal Components.....	25
Cartridge Slot.....	25
Sample and Reagent Platform	26
Fractionation Adapter.....	29

Filter.....	31
Using the Plate Lid.....	31
Rear Panel.....	32
System Labels.....	33
Computer Workstation.....	34

Chapter 4: Compass for iCE Overview..... 35

Launching Compass for iCE	36
Compass for iCE Overview	36
Changing the Screen View	37
Batch Screen	37
Run Summary Screen.....	39
Analysis Screen	41
Screen Panes	42
Title Bar.....	43
Main Menu	43
Instrument Status Bar.....	43
Screen Tab.....	44
View Bar.....	44
Compass for iCE Status Bar	44
Software Menus	44
File Menu	45
Edit Menu	45
View Menu	45
Instrument Menu	46
Window Menu	46
Help Menu.....	47

Changing the Compass for iCE	
Main Window Layout.....	47
Resizing the Main Compass for iCE Window	47
Resizing the Screen Tab.....	47
Resizing Screen Panes	48
Changing the Location of Screen Panes	49
Restoring the Main Window to the Default Layout.....	50
Viewing the User Guide.....	50
Visiting the Bio-Techne Academy.....	50
Checking for and Installing New Versions of Compass for iCE	50
Viewing Release Notes.....	51
Sending Run Files to Technical Support	51
Compass for iCE Version Information.....	52
Directory and File Information.....	53
File Types.....	54

Chapter 5:	
cIEF Batches.....	55
Overview	56
Batch Screen Overview	56
Batch Screen Panes.....	56
Software Menus Active in the Batch Screen	57
Opening a Batch.....	58
Creating a New Batch	59
Step 1 - Open a Template Batch	59
Step 2 - Assign Your Samples.....	60
Step 3 - Assign Your Method Parameters.....	63
Step 4 - Set Up Your Injections	67
Step 5 - Setup On-Board Mixing (Optional)	70
Step 6 - Add Programmed Pauses and Stops (Optional).....	73

Step 7 - Add Batch Notes (Optional)	74
Step 8 - Modify Default Analysis Parameters (Optional)	75
Step 9 - Modify Default Analysis View Parameters (Optional).....	76
Step 10 - Save Your Batch	76
Viewing Replicate Injections	77
On-board Mixing (Maurice and Maurice C.).....	78
Batch History	80
Making Changes to a Batch	81
Viewing and Editing Batches in Completed Runs.....	82
Copying and Pasting Injection Names or Sample IDs from other Documents	82
Importing and Exporting Injections	84
Exporting Injections.....	85
Creating an Injections File	85
Importing Injections	86
Importing and Exporting Methods.....	86
Importing a Method	86
Exporting a Method.....	87
Batch Reports.....	87

Chapter 6:	
Running cIEF Applications	90
Overview	91
Before You Throw the Switch.....	91
Power Up	91
Running cIEF Applications	91
What You'll Need	91
Step 1: Prep Your Markers, Samples and Reagents.....	92
Step 2: Prep the cIEF Cartridge	96
Step 3: Install the Cartridge.....	97

Step 4: Load Samples and Reagents	99
Step 5: Check for Cartridge Alerts.....	101
Step 6: Create a Batch.....	104
Step 7: Start the Batch	116
Editing a Running Batch	119
Post-batch Procedures	120
Checking Your Data	123
Step 1: Select Your Detection Mode.....	123
Step 2: Check Your pI Markers.....	124
Step 3: Checking Sample Peaks.....	126
Step 4: Assigning Peak Names.....	128

Chapter 7: MauriceFlex cIEF Batches..... 129

Overview	130
Batch Screen Overview	130
Batch Screen Panes.....	130
Software Menus Active in the Batch Screen	131
Opening a Batch.....	133
Creating a New Batch	133
Step 1 - Open a Template Batch	133
Step 2 - Assign Your Samples.....	134
Step 3 - Assign Your Method Parameters.....	137
Step 4 - Set Up Your Injections.....	140
Step 5 - Add Programmed Pauses and Stops (Optional).....	142
Step 6 - Add Batch Notes (Optional)	144
Step 7 - Modify Default Analysis Parameters (Optional)	144
Step 8 - Modify Default Analysis View Parameters (Optional).....	145
Step 9 - Save Your Batch	145
Viewing Replicate Injections.....	146

Batch History	147
Making Changes to a Batch	148
Viewing and Editing Batches in Completed Runs.....	148
Copying and Pasting Injection Names or Sample IDs from other Documents	149
Importing and Exporting Injections	150
Exporting Injections.....	151
Creating an Injections File	151
Importing Injections	152
Importing and Exporting Methods.....	152
Importing a Method	152
Exporting a Method.....	153
Batch Reports.....	153

Chapter 8: Running MauriceFlex cIEF Applications..... 156

Overview	157
Before You Throw the Switch.....	157
Power Up	157
Running MauriceFlex cIEF Applications	157
What You'll Need	157
Step 1: Prep Your Markers, Samples and Reagents.....	159
Step 2: Load Samples and Reagents	161
Step 3: Prep the cIEF Fractionation Cartridge	163
Step 4: Install the Cartridge.....	165
Step 5: Check for Cartridge Alerts.....	166
Step 6: Create a Batch.....	168
Step 7: Start the Batch	178
Editing a Running Batch	180
Post-batch Procedures	182
Checking Your Data	184
Step 1: Check Your pI Markers.....	185

Step 2: Checking Sample Peaks.....	187
Step 3: Assigning Peak Names.....	189

Chapter 9: MauriceFlex Fractionation Batches 190

Overview	191
Batch Screen Overview	191
Batch Screen Panes.....	191
Software Menus Active in the Batch Screen	192
Opening a Batch.....	194
Creating a New Batch	194
Step 1 - Open a Template Batch	195
Step 2 - Assign Your Samples.....	196
Step 3 - Assign Your Method Parameters.....	200
Step 4 - Set Up Your Injection.....	204
Step 5 - Add Programmed Pauses and Stops (Optional).....	206
Step 6 - Add Batch Notes (Optional)	207
Step 7 - Modify Default Analysis Parameters (Optional)	208
Step 8 - Modify Default Analysis View Parameters (Optional).....	209
Step 9 - Save Your Batch	209
Viewing Fractions To Be Collected.....	210
Batch History	211
Making Changes to a Batch	212
Viewing and Editing Batches in Completed Runs.....	212
Copying and Pasting an Injection Name or Sample ID from other Documents	213
Importing and Exporting an Injection	214
Exporting an Injection.....	214
Creating an Injection File.....	215
Importing an Injection.....	215

Importing and Exporting Methods.....	216
Importing a Method	216
Exporting a Method.....	216
Batch Reports.....	217

Chapter 10: Running MauriceFlex Fractionation Applications... 220

Overview	221
Before You Throw the Switch.....	221
Power Up	221
MauriceFlex Fractionation Workflow	222
Running MauriceFlex Fractionation Applications	223
What You'll Need	223
Step 1: Prep Your Markers, Samples and Reagents.....	224
Step 2: Load Samples and Reagents	226
Step 3: Prep the cIEF Fractionation Cartridge	229
Step 4: Install the Cartridge.....	231
Step 5: Check for Cartridge Alerts.....	232
Step 6: Create a Batch.....	234
Step 7: Start the Batch	244
Editing a Running Batch	246
Post-batch Procedures	248
Checking Your Data	250
Step 1: Check the pI Markers for the Injected Sample.....	251
Step 2: Checking Sample Peaks.....	254
Step 3: Assigning Sample Peak Names.....	256
Step 4: Checking Peaks in the Mobilization Electropherograms.....	256
Step 5: Assigning Peak Names in the Mobilization Electropherograms.....	258

Chapter 11: CE-SDS PLUS Batches 259

Overview	260
Batch Screen Overview	260
Batch Screen Panes.....	260
Software Menus Active in the Batch Screen	261
Opening a Batch	262
Creating a New Batch	263
Step 1 - Open a Template Batch	263
Step 2 - Assign Your Samples	264
Step 3 - Assign Your Method Parameters.....	267
Step 4 - Set Up Your Injections	269
Step 5 - Add Programmed Pauses and Stops (Optional).....	272
Step 6 - Add Batch Notes (Optional)	274
Step 7 - Modify Default Analysis Parameters (Optional)	274
Step 8 - Modify Default Analysis View Parameters (Optional).....	275
Step 9 - Save Your Batch	275
Viewing Replicate Injections	276
Batch History	277
Making Changes to a Batch	278
Viewing and Editing Batches in Completed Runs.....	278
Copying and Pasting Injection Names or Sample IDs from other Documents	279
Importing and Exporting Injections	281
Exporting Injections.....	281
Creating an Injections File	282
Importing Injections	282
Importing and Exporting Methods.....	283
Importing a Method	283
Exporting a Method.....	283

Batch Reports.....	284
--------------------	-----

Chapter 12: Running CE-SDS PLUS Applications..... 287

Overview	288
Before You Throw the Switch.....	288
Power Up	288
Running CE-SDS PLUS Applications	288
What You'll Need	288
Step 1: Prep Your Internal Standard, Samples and Reagents.....	289
Step 2: Prep the Cartridge.....	295
Step 3: Install the Cartridge.....	297
Step 4: Load Samples and Reagents	299
Step 5: Check for Cartridge Alerts.....	300
Step 6: Create a Batch.....	303
Step 7: Start the Batch	312
Editing a Running Batch	315
Post-batch Procedures	316
Checking Your Data	319
Step 1: Check Your Internal Standard	319
Step 2: Set Your Molecular Weight (MW) Markers	322
Step 3: Checking Sample Peaks.....	325
Step 4: Assigning Peak Names.....	327

Chapter 13: Turbo CE-SDS Batches 328

Overview	329
Batch Screen Overview	329
Batch Screen Panes.....	329
Software Menus Active in the Batch Screen	330
Opening a Batch	331
Creating a New Batch	332

Step 1 - Open a Template Batch	332
Step 2 - Assign Your Samples	333
Step 3 - Assign Your Method Parameters.....	336
Step 4 - Set Up Your Injections	338
Step 5 - Add Programmed Pauses and Stops (Optional).....	341
Step 6 - Add Batch Notes (Optional)	343
Step 7 - Modify Default Analysis Parameters (Optional)	343
Step 8 - Modify Default Analysis View Parameters (Optional).....	344
Step 9 - Save Your Batch	344
Viewing Replicate Injections	345
Batch History	346
Making Changes to a Batch	346
Viewing and Editing Batches in Completed Runs.....	347
Copying and Pasting Injection Names or Sample IDs from other Documents	348
Importing and Exporting Injections	350
Exporting Injections.....	350
Creating an Injections File	350
Importing Injections	351
Importing and Exporting Methods.....	351
Importing a Method	351
Exporting a Method.....	352
Batch Reports.....	353
Chapter 14:	
Running Turbo CE-SDS Applications	356
Overview	357
Before You Throw the Switch.....	357
Power Up	357
Running Turbo CE-SDS Applications.....	357

What You'll Need	357
Step 1: Prep Your Internal Standard, Samples and Reagents	358
Step 2: Prep the Cartridge.....	364
Step 3: Install the Cartridge.....	366
Step 4: Load Samples and Reagents	368
Step 5: Check for Cartridge Alerts.....	369
Step 6: Create a Batch.....	371
Step 7: Start the Batch	381
Editing a Running Batch	384
Post-batch Procedures	385
Checking Your Data	388
Step 1: Check Your Internal Standard.....	388
Step 2: Set Your Molecular Weight (MW) Markers	391
Step 3: Checking Sample Peaks.....	394
Step 4: Assigning Peak Names.....	396

Chapter 15:

Cartridge and Adapter Alerts 397

Overview	398
Cartridge Alerts	398
Cartridge Maintenance	398
Cartridge Prep	402
Cartridge Type.....	402
Adapter and Insert Alerts	403
Alerts when Starting a Batch.....	403
Alerts when Pausing a Batch	404
Compass Alerts	404

Chapter 16:

Run Status 406

Run Summary Screen Overview	407
Run Summary Screen Panes	407

Software Menus Active in the Run Summary Screen.....	409
Opening Run Files	411
Opening One Run File.....	411
Opening Multiple Run Files	411
Batch Injection Information.....	413
Injection Flags	414
Run Status Information.....	415
Viewing the Injection Focus Series (cIEF, MauriceFlex cIEF, MauriceFlex Fractionation Only)	417
Viewing the Fractionation Focus Series (MauriceFlex Fractionation Only)	418
Viewing the Separation Plot (CE-SDS PLUS and Turbo CE-SDS Only)	420
Current and Voltage Plots.....	420
Run History	422
Viewing Multiple Events.....	423
Copying History Info	424
Viewing Run Errors or Warnings.....	424
Injection Reports	425
Example Analysis and Injection Report: CE-SDS.....	427
Example Analysis and Injection Report: cIEF	431
Example Analysis and Injection Report: MauriceFlex Fractionation	434
Switching Between Open Run Files.....	437
Closing Run Files	437

Chapter 17: Controlling Maurice, Maurice C., Maurice S. and MauriceFlex 438

Instrument Control.....	439
Starting a Run	439
Stopping a Run.....	442

Emergency Stop.....	443
Status Modes.....	444
Shutdown	444
Instrument Software (Embedded) Updates.....	445
Instrument Self Test.....	446
Viewing and Changing System Properties.....	447
Checking Cartridge Status	448
Cartridge Self-Test.....	449
Instrument Reports.....	452
Example Instrument Report	453
Viewing Log Files	467
Runs Log.....	467
Sending a Log File to Technical Support.....	468
Instrument Log	469
Self Test Logs.....	470
Command Log	471

Chapter 18: cIEF Data Analysis 473

Analysis Screen Overview	474
Analysis Screen Panes	474
Software Menus Active in the Analysis Screen.....	475
Opening Run Files	477
Opening One Run File.....	477
Opening Multiple Run Files	477
How Run Data is Displayed	478
Experiment Pane: Batch Injection Information	478
Graph Pane: Electropherogram Data	479
Peaks Pane: Calculated Results.....	480
Injections Pane: User-Specified Peak Names.....	482
Analysis Options Pane	483
Viewing Run Data.....	484

Switching Between Samples and Markers Data Views	484	Changing the Electropherogram View	507
Selecting and Displaying Injection Data	486	Autoscaling the Electropherogram	508
Switching Between Selected and All Views of Injections	488	Customizing the Data Display	509
Hiding Injection Data	490	Stacking Multiple Electropherograms	514
Switching Between Absorbance and Fluorescence Exposures (For Standard cIEF Runs only)	490	Overlaying Multiple Electropherograms	515
Data Notifications and Warnings	492	Overlaying Fluorescence and Absorbance Electropherograms (standard cIEF runs only)	515
Checking Your Results	493	Zooming	517
Naming Peaks	493	Selecting Data Viewing Options	518
Adding New Peak Names in the Analysis Options Pane	493	Adjusting the Baseline	522
Adding Peak Names from the Graph or Peaks Table	496	Selecting the Graph X-axis Range	526
Changing Peak Name Information	498	Closing Run Files	528
Group Statistics	498	Analysis Settings Overview	528
Using Groups	498	Detection Settings	529
Viewing Sample Injection Groups	499	Changing the Detection Method (Standard Maurice cIEF runs only)	530
Viewing Statistics	500	Changing the Detection Exposure	531
Hiding or Removing Injections in Group Analysis	501	Peak Names Settings	532
Copying Results Tables and Graphs	501	Peak Names Analysis Settings Groups	533
Copying Results Tables	502	Creating a Peak Names Group	533
Copying the Graph	502	Modifying a Peak Names Group	536
Saving the Graph as an Image File	502	Deleting a Peak Names Group	536
Exporting Run Files	503	Applying Peak Names Groups to Run Data	536
Exporting Results Tables	503	Peak Fit Analysis Settings	539
Exporting Raw Sample Electropherogram Data	503	Range Settings	540
Changing Sample Protein Identification	504	Baseline Settings	541
Adding or Removing Sample Data	504	Peak Find Settings	541
Hiding Sample Data	505	Peak Fit Analysis Settings Groups	542
Changing Peak Names for Sample Data	506	Creating a New Peak Fit Group	543
		Changing the Default Peak Fit Group	544
		Modifying a Peak Fit Group	545
		Deleting a Peak Fit Group	545

Applying Peak Fit Groups to Specific Run Data	546
Manual Peak Integration	547
Advanced Analysis Settings	551
pI Markers Settings	552
Advanced Analysis Settings Groups	552
Creating a New Analysis Group	553
Changing the Default Analysis Group	554
Modifying an Analysis Group	555
Deleting an Analysis Group	555
Applying Analysis Groups to Specific Run Data	556
pI Markers Analysis Settings	558
pI Markers Analysis Settings Groups	558
Creating a New Markers Group	559
Changing the Default Markers Group	561
Modifying a Markers Group	561
Deleting a Markers Group	562
Applying Markers Groups to Specific Run Data	562
Injection Reports	564
Example Analysis and Injection Report	566
Importing and Exporting Analysis Settings	567
Importing Analysis Settings	567
Exporting Analysis Settings	568
Chapter 19:	
MauriceFlex Fractionation Data Analysis	569
Analysis Screen Overview	570
Analysis Screen Panes	570
Software Menus Active in the Analysis Screen	571
Opening Run Files	573
Opening One Run File	573
Opening Multiple Run Files	573
How Run Data is Displayed	574
Experiment Pane: Batch Injection Information	575
Graph Pane: Electropherogram Data	575
Peaks Pane: Calculated Results	576
Injections Pane: User-Specified Peak Names	578
Fractions Pane: Fraction Peak Predictor	579
Analysis Options Pane	580
Viewing Run Data	580
Switching Between Samples and Fractions Data Views	580
Selecting and Displaying Injection Data	583
Switching Between Selected and	
All Electropherogram Views (Fractions View)	585
Data Notifications and Warnings	587
Checking Your Results	589
Naming Peaks	589
Adding New Peak Names in the	
Analysis Options Pane	589
Adding Peak Names from the Graph or	
Peaks Table	592
Changing Peak Name Information	593
Copying Results Tables and Graphs	594
Copying Results Tables	594
Copying the Graph	594
Saving the Graph as an Image File	595
Exporting Run Files	595
Exporting Results Tables	595
Exporting Raw Sample Electropherogram Data	596
Changing Sample Protein Identification	596
Adding or Removing Sample Data	596
Hiding Sample Data	598
Changing Peak Names for Sample Data	599
Adding or Removing Fractions Data	599

Changing the Electropherogram View.....	601	Advanced Analysis Settings Groups	642
Autoscaling the Electropherogram	602	Creating a New Analysis Group.....	643
Customizing the Data Display	603	Changing the Default Analysis Group	644
Stacking Multiple Electropherograms	608	Modifying an Analysis Group	645
Overlaying Multiple Electropherograms	609	Deleting an Analysis Group	645
Zooming	609	Applying Analysis Groups to Specific Run Data	645
Selecting Data Viewing Options	610	pI Markers Analysis Settings	647
Adjusting the Baseline.....	615	pI Markers Analysis Settings Groups	648
Selecting the Graph X-axis Range.....	618	Creating a New Markers Group	649
Closing Run Files	620	Changing the Default Markers Group.....	651
Analysis Settings Overview	620	Modifying a Markers Group	651
Detection Settings.....	621	Deleting a Markers Group	651
Peak Names Settings	622	Applying Markers Groups to Specific Run Data.....	652
Peak Names Analysis Settings Groups.....	622	Injection Reports	653
Creating a Peak Names Group.....	623	Example Analysis and Injection Report	655
Modifying a Peak Names Group	626	Importing and Exporting Analysis Settings	657
Deleting a Peak Names Group	626	Importing Analysis Settings	657
Applying Peak Names Groups to Run Data	626	Exporting Analysis Settings.....	657
Peak Fit Analysis Settings	629	Chapter 20:	
Range Settings	630	CE-SDS Data Analysis	659
Baseline Settings	631	Analysis Screen Overview	660
Peak Find Settings.....	631	Analysis Screen Panes	660
Peak Fit Analysis Settings Groups.....	632	Software Menus Active in the Analysis Screen.....	661
Creating a New Peak Fit Group	633	Opening Run Files	663
Changing the Default Peak Fit Group.....	634	Opening One Run File.....	663
Modifying a Peak Fit Group	634	Opening Multiple Run Files	663
Deleting a Peak Fit Group	635	How Run Data is Displayed	664
Applying Peak Fit Groups to Specific Run Data	635	Experiment Pane: Batch Injection Information	665
Manual Peak Integration	637	Graph Pane: Electropherogram Data	665
Advanced Analysis Settings	642	Lane Pane: Virtual Gel-Like Image Data	666
pI Markers Settings.....	642	Peaks Pane: Calculated Results	667

Injections Pane: User-Specified Peak Names.....	669	Changing Peak Names for Sample Data.....	693
Analysis Options Pane: Peak Names	671	Changing the Electropherogram View.....	694
Viewing Run Data.....	671	Autoscaling the Electropherogram	694
Switching Between Samples and		Customizing the Data Display	695
Standards Data Views.....	671	Stacking Multiple Electropherograms	699
Selecting and Displaying Injection Data	673	Overlaying Multiple Electropherograms	700
Switching Between Single and		Zooming	701
Multiple Views of Injections	675	Selecting Data Viewing Options	702
Hiding Injection Data	676	Adjusting the Baseline.....	706
Data Notifications and Warnings.....	678	Selecting the Graph X-axis Range.....	709
Checking Your Results	679	Changing the Virtual Gel View	711
Naming Peaks.....	679	Adjusting the Contrast	712
Adding New Peak Names in the		Inverting the Virtual Gel.....	712
Analysis Options Pane	679	Selecting Lane Labels	713
Adding Peak Names from the Graph or		Closing Run Files	713
Peaks Table	683	Analysis Settings Overview.....	713
Changing Peak Name Information	684	Markers Analysis Settings.....	715
Group Statistics	684	Markers Settings.....	716
Using Groups	684	Changing the Injection Used for the	
Viewing Sample Injection Groups	685	CE-SDS MW Markers	716
Viewing Statistics	686	Standards Analysis Settings Groups.....	716
Hiding or Removing Injections in		Creating a New Standards Group	717
Group Analysis.....	688	Changing the Default Standards Group.....	720
Copying Data Views and Results Tables	688	Modifying a Standards Group	720
Copying Results Tables	688	Deleting a Standards Group	720
Copying Data Views	688	Applying Standards Groups to	
Saving the Data Views as an Image File	688	Specific Run Data	721
Exporting Run Files	689	Peak Names Settings	723
Exporting Results Tables	689	Peak Names Analysis Settings Groups.....	723
Exporting Raw Sample Electropherogram Data	689	Creating a Peak Names Group.....	724
Changing Sample Protein Identification.....	690	Modifying a Peak Names Group	727
Adding or Removing Sample Data	690	Deleting a Peak Names Group	727
Hiding Sample Data.....	691		

Applying Peak Names Groups to Run Data	727
Peak Fit Analysis Settings	731
Range Settings	731
Baseline Settings	732
Peak Find Settings.....	732
Peak Fit Analysis Settings Groups.....	734
Creating a New Peak Fit Group	734
Changing the Default Peak Fit Group.....	736
Modifying a Peak Fit Group	736
Deleting a Peak Fit Group	737
Applying Peak Fit Groups to Specific Run Data	737
Manual Peak Integration	739
Advanced Analysis Settings	743
Internal Standard Settings.....	743
Advanced Analysis Settings Groups	743
Creating a New Analysis Group.....	744
Changing the Default Analysis Group	745
Modifying an Analysis Group	746
Deleting an Analysis Group	746
Applying Analysis Groups to Specific Run Data	747
Signal Processing Settings.....	749
Reference Channel Correction	749
Baseline Smoothing	750
Injection Reports	752
Example Analysis and Injection Report	754
Importing and Exporting Analysis Settings	756
Importing Analysis Settings	756
Exporting Analysis Settings.....	757

Chapter 21: **Setting Your Preferences 758**

Customize Your Preferences	759
----------------------------------	-----

Enabling Access Control	760
Setting Data Export Options	760
Setting up Automatic Injection Reports	761
Selecting Custom Plot Colors for Graph Overlay.....	762
Grouping Options.....	764
Setting Up Maurice Systems to Send Tweets.....	764

Chapter 22: **Compass Access Control and** **21 CFR Part 11 Compliance..... 768**

Overview	769
Compass Authorization Server for iCE	769
Server Administration.....	769
Adding Non-admin Users	771
Adding Admin Users	775
Resetting User Passwords.....	776
Audit Trail	777
Password Policy Settings	780
LDAP Settings.....	780
Encryption Details.....	781
Enabling Access Control	781
Changing the Software Inactivity Auto Lock	782
Logging In to Compass for iCE.....	782
Locking and Unlocking the Application.....	783
Resolving Log In Issues.....	784
Saving Changes.....	785
Signing Files.....	785
Exporting Uncontrolled Files.....	786
Instrument Command Log.....	787
Run File History	788
Troubleshooting Problems and	
Suggested Solutions.....	788

Chapter 23:
Maintenance and Troubleshooting..... 790

Cartridge Handling and Care.....	791
cIEF Cartridge.....	792
cIEF Fractionation Cartridge	793
CE-SDS PLUS Cartridges	793
Turbo CE-SDS Cartridges	794
Installing or Replacing the System Filter	795
Maintenance	798
Daily	798
Yearly.....	798
Changing the Fuse	798
Spare Parts	799
Software Updates	800
Instrument Software (Embedded) Updates.....	800
Frequently Asked Questions: General.....	800
Frequently Asked Questions: cIEF Applications	801
Frequently Asked Questions: cIEF Fractionation Applications	805
Frequently Asked Questions: CE-SDS Applications	807
Best Practices	808
Maurice cIEF	808
MauriceFlex cIEF and MauriceFlex Fractionation	808
Maurice CE-SDS	808
General Tips	809
Troubleshooting	809
Maurice System Troubleshooting	810
cIEF Application Troubleshooting.....	810
cIEF Fractionation Application Troubleshooting	812
CE-SDS Application Troubleshooting.....	813

Chapter 24:
General Information..... 815

Compliance	816
Safety Guidelines	817
Chemical Hazards.....	817
Chemical Waste Hazards	818
Waste Production and Disposal.....	818
Safety Data Sheets	819
Instrument Safety Labels	820
Consumables and Reagents.....	820
Maurice CE-SDS Consumables, Kits and Reagents	820
Maurice cIEF Consumables, Kits and Reagents	822
MauriceFlex cIEF Fractionation Consumables, Kits and Reagents	825
Customer Service and Technical Support	826
Legal Notices	827
Maurice Disclaimer of Warranty	827
Compass Software and Compass Authorization Server for iCE License Agreement.....	827

Chapter 1:

Let's Get Started

Chapter Overview

- Welcome
- Maurice Systems

Welcome

Congratulations on bringing Maurice into your lab! We welcome you as a new user and are excited to be a part of your work. This user guide will provide you with details on system hardware, operating the system, how to use Compass for iCE software, maintenance procedures and other useful information.

To help you get the most from your new lab addition, we've added some attention phrases to guide you through the user guide:

NOTE	Points out useful information.
IMPORTANT	Indicates information necessary for proper operation of Maurice systems.
CAUTION	Cautions you about potentially hazardous situations that could result in injury to you or damage to the system.
!WARNING!	Warns you that serious physical injury can result if the listed precautions aren't followed.

Maurice Systems

Maurice, Maurice C., Maurice S. and MauriceFlex systems give you identity, purity and heterogeneity data on your biologics, and get you to results faster with short development times and simple workflows!

- **They're fluent in cIEF and CE-SDS.** They take cIEF up a notch, and CE-SDS is a breeze. You'll get pI and charge heterogeneity data in less than 10 minutes flat — with the added bonus of same-time absorbance and native fluorescence for sensitivity down to 0.7 µg/mL. Their size applications have the high res and wide molecular weight range you need and they're done in as little as 5.5 minutes.
- **They collect fractions of the charge variants:** The feature of icIEF fractionation enables you to isolate the protein charge variants for in-depth characterization by mass spectrometry, and the offline fraction collection gives you the freedom to carry out the intact mass, reduced and subunit mass, and peptide mapping analyses on any vendor's mass spectrometers.
- **They make it easy.** Just pop in a ready-to-go cartridge, drop in your sample vials or a 96-well plate in the appropriate adaptor, and hit start — they'll do the rest!
- **They're time-savers.** Develop methods fast so you get to results even faster. Your cIEF and CE-SDS methods are done in a day. The icing? You can develop platform methods and use them for multiple molecules. No maintenance and clean-up needed between the two applications.
- **They're dependable.** Get reproducible results with tight CVs day in and day out. Your data is reliable no matter what — across samples, users, instruments or labs.

Chapter 2:

Getting Your Lab Ready

Chapter Overview

- Introduction
- Space Requirements
- Physical Specifications
- Electrical Requirements
- Environmental Requirements
- Software and Computer Requirements
- General Guidelines and Information

Introduction

This chapter will help you prepare the lab for Maurice. Please have the space, electrical and environmental requirements ready prior to scheduling your installation.

NOTE: Please wait for an authorized ProteinSimple Field Service Engineer to unpack and install Maurice for you. Don't try doing this yourself. Handling Maurice incorrectly could cause injury to yourself or damage to the system.

Space Requirements

You need a lab bench or table that can support 100 lb (46 kg) and has enough space for both Maurice and his computer. There should be sufficient clearance for both heat ventilation and to provide access if Maurice needs service.

IMPORTANT

Maurice needs a stable surface and must remain level to work properly. The lab bench or table can't shift or wobble under heavy weight. Don't use anti-vibration tables either, since Maurice may not stay level while he's working.

Dimension	Meters	Feet
Width	1.5	5.0
Depth	0.8	2.5
Height	0.5	1.5

Recommended space requirements for Maurice.

Physical Specifications

Description	Specification
Maurice's Dimensions (Door Closed)	0.44 m x 0.42 m x 0.61 m (H x W x D) 1.46' x 1.38' x 2.0' (H x W x D)
Maurice's Dimensions (Door Open)	0.44 m x 0.57 m x 0.61 m (H x W x D) 1.46' x 2.43' x 2.0' (H x W x D)
Maurice's Weight	46 kg (100 lb)
Computer Workstation Dimensions	Computer Workstation Dimensions

For indoor use only. Use up to altitudes of 1524 meters (5000 feet).

Electrical Requirements

Maurice requires a dedicated, grounded circuit capable of delivering the appropriate current and voltage for your country. The power requirements for all three Maurice systems are 100 V–240 V (AC), 50/60 Hz, 500 W.

In addition to these requirements, Maurice needs the grounded circuits terminate at the receptacles, and receptacles must be located within 10 ft (3 m) of the instrument.

Environmental Requirements

Maurice likes a consistent temperature in the lab (not too hot – not too cold). He works best when conditions stay within these ranges:

Requirement	Specification
Operating temperature range	8–25 °C (64–77 °F)
Operating humidity range	20–80% relative, non-condensing

Software and Computer Requirements

Maurice brings his own computer to the lab with Compass for iCE software pre-installed.

Using Maurice with Compass for iCE

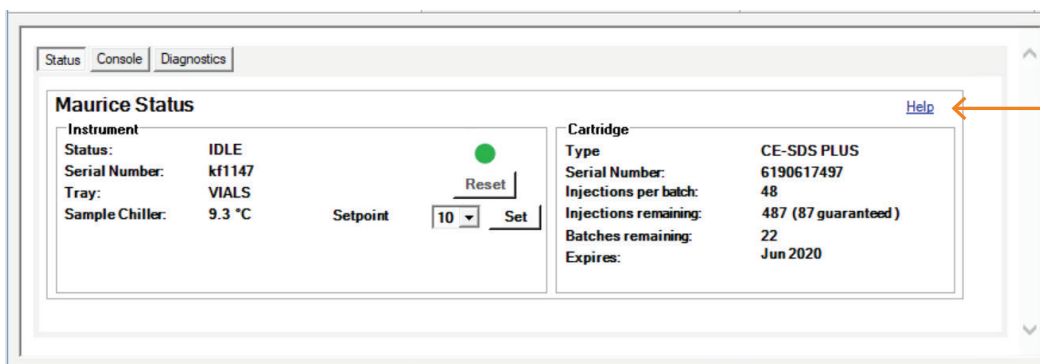
Provided with all Maurice systems, Compass for iCE is used to run cIEF and CE-SDS applications on Maurice, cIEF applications on Maurice C., CE-SDS applications on Maurice S. and cIEF, CE-SDS, MauriceFlex cIEF and MauriceFlex Fractionation applications on MauriceFlex, and analyze resulting data. Just in case you need it, a CD containing Compass for iCE software also comes in the box. If you don't want to analyze your data at Maurice's workstation in the lab, Compass for iCE software can also be installed on a separate workstation, such as your desktop computer. Your computer must meet the recommended requirements listed below to run Compass for iCE software and process data.

Component	Minimum Requirements
Operating System	Windows 10
Processor	Core i5
Memory	8 GB
Free Disk Space	250 GB
Ethernet Ports	2 - One is required to connect to Maurice or MauriceFlex, the other is used for network access
USB Ports	2 - To connect the keyboard and mouse

Using Maurice with Empower®

The optional Maurice Empower® Control Kit lets you collect, manage and report Maurice results with the Waters™ Empower® 3 Chromatography Data Software. You can use Empower to control a Maurice, Maurice C. or Maurice S. when performing a Maurice cIEF or CE-SDS PLUS batch. To use Empower®, contact ProteinSimple to install the instrument software if not previously done. Then install the Maurice ICS driver software following the instructions in the Maurice with Empower® Installation Guide on the CD that came in your Maurice Empower® Control kit.

To learn how to create Maurice method and sample sets, run samples and view data in Empower®, open the Run Samples window and click **Help** in the top right corner of the Maurice Control Panel.



General Guidelines and Information

Intended Use

NOTE: Maurice is for research use only. Not for use in diagnostic procedures.

Lifting and Moving the System: Lift Maurice Correctly

IMPORTANT

Take all the standard precautions when lifting or moving Maurice. Since Maurice systems weigh 46 kg (100 lb), you should not lift him by yourself. Two people should lift him onto the lab bench.

Chapter 3:

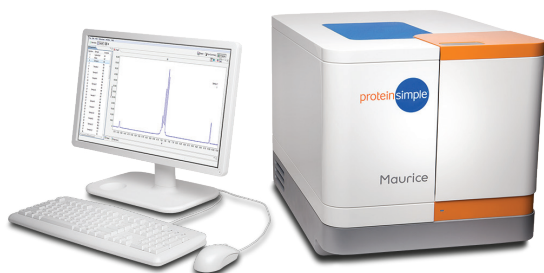
Maurice

Chapter Overview

- Maurice Systems
- External Components
- Internal Components
- Computer Workstation

Maurice Systems

Maurice, Maurice C., Maurice S. and MauriceFlex systems include the instrument, computer workstation, Compass for iCE software and cIEF, CE-SDS PLUS, Turbo CE-SDS or cIEF Fractionation Cartridges.



Maurice with Computer Workstation



CE-SDS, cIEF and cIEF Fractionation Cartridges

All systems have the same hardware components, computer and software, the only difference between them are the applications you can run:

Application	Cartridge	Maurice	Maurice C.	Maurice S.	MauriceFlex
cIEF	cIEF	●	●		●
cIEF with on-board mixing	cIEF	●	●		
CE-SDS	CE-SDS PLUS or Turbo CE-SDS	●		●	●
MauriceFlex cIEF	cIEF Fractionation				●
MauriceFlex Fractionation	cIEF Fractionation				●

You can run samples for a standard cIEF or CE-SDS batch in 96-well plates or in up to 48 sample vials with an integrated 0.2 mL insert. You can run your sample(s) for a MauriceFlex cIEF or MauriceFlex Fractionation batch in a 96-well plate on a MauriceFlex system. Fractions for a MauriceFlex Fractionation batch are collected in the 96-well plates.



Maurice C.



Maurice

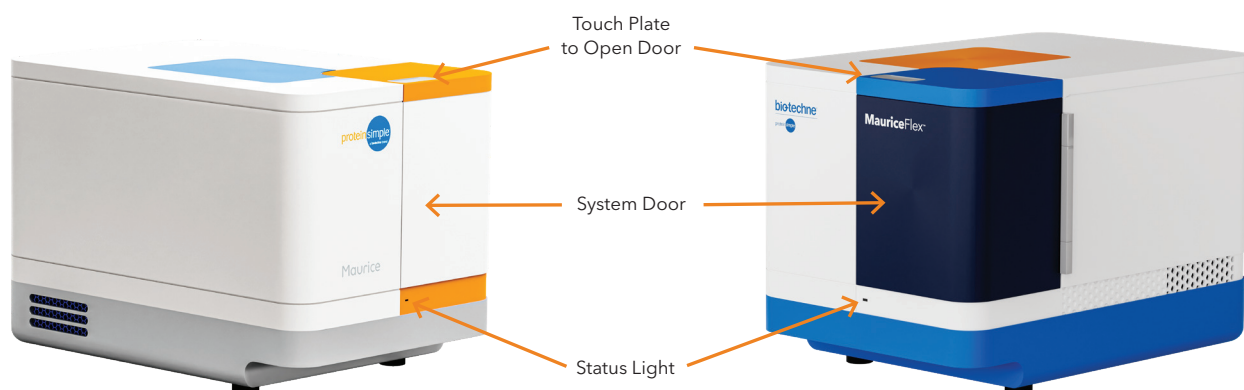


MauriceFlex



Maurice S.

External Components



!WARNING!

You can't replace or service any parts on Maurice systems except for the power entry fuse.

System Door

Maurice's door gives you access to the inside of the instrument to load cartridges, reagents and samples. To open the door, first make sure the status light is a steady blue. Then just touch the metal touch plate on the top. The blue status light will flash and the door will automatically open. Close it by pushing the door from the front until you hear the latch engage and see the status light turn solid.

NOTE: Maurice's door must be closed before starting a batch.

Status Light

The LED on Maurice's front panel tells you what he's doing. Here's what his different status lights mean:

- **Start-up (magenta):** You've just turned on the power and Maurice is warming up.
- **Ready (steady blue):** Maurice is powered on and ready to go.

NOTE: If after you power Maurice on the LED doesn't change from magenta to blue, the system didn't initialize. Please call ProteinSimple Technical Support.

- **Opening Door (long blue flash followed by blue pulses):** Maurice's door is opening.
- **Running (pulsing blue):** Maurice is running a batch.

- **Paused (blue pulse with magenta flashes):** The batch is paused. You can edit the batch and open Maurice's door to add new samples or reagents.
- **Trying to Open Door While Running (red flash):** Maurice's door can't be opened when he's running.
- **Error (steady red):** Maurice has detected an error. To get more information on the error, check the Status pane in the Run Summary Screen in Compass for iCE.



Internal Components

Cartridge Slot

The cartridge slot holds Maurice's ready-to-go application cartridges. The cartridge it holds depends on the system:

- **Maurice:** cIEF, CE-SDS PLUS and Turbo CE-SDS Cartridges only
- **Maurice C.:** cIEF Cartridges only
- **Maurice S.:** CE-SDS PLUS and Turbo CE-SDS Cartridges only
- **MauriceFlex:** cIEF, CE-SDS PLUS, Turbo CE-SDS and cIEF Fractionation Cartridges

The lights on either side of the cartridge slot will be **orange** after Maurice disengages the cartridge when the door is opened at the end of a batch, and whenever the slot is empty.



The lights change to **blue** once a cartridge is installed correctly.

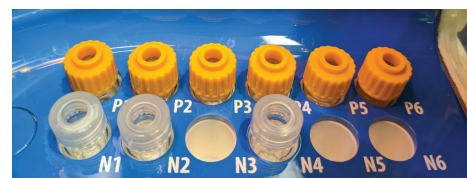
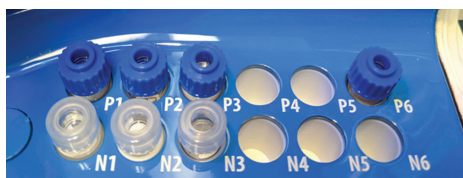


NOTE: You can find cartridge prep, installation and post-run procedures in Chapter 6: “Running cIEF Applications”, Chapter 8: “Running MauriceFlex cIEF Applications”, Chapter 10: “Running MauriceFlex Fractionation Applications”, Chapter 12: “Running CE-SDS PLUS Applications” and Chapter 14: “Running Turbo CE-SDS Applications”

Sample and Reagent Platform

Maurice’s sample and reagent platform for performing standard cIEF applications using a cIEF cartridge and CE-SDS applications using a CE-SDS PLUS or Turbo CE-SDS cartridge has two rows for batch reagents. These reagents are kept at room temperature.

- **Row P (top):** These reagents are loaded under pressure during the batch. Only use glass reagent vials with pressure caps in this row. Use **blue** pressure caps with cIEF reagents and **orange** pressure caps with CE-SDS reagents.
- **Row N (bottom):** Only use reagent vials with clear screw caps in this row.



Reagent platform for cIEF applications (left) and CE-SDS applications (right) on Maurice, Maurice C. and Maurice S.



Reagent platform for cIEF applications (left) and CE-SDS applications (right) on MauriceFlex.

IMPORTANT:

Reagent vials on MauriceFlex must be locked in place by sliding the locking mechanism from the left to the right before starting a batch.

Use glass reagent vials, 2 mL (PN 046-017) for standard cIEF and CE-SDS applications on Maurice, Maurice C., Maurice S. and MauriceFlex.

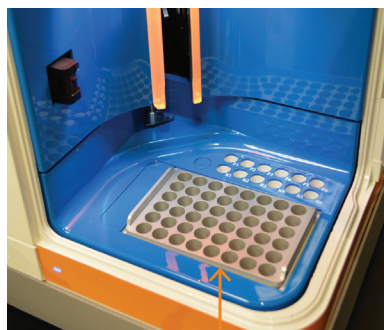
The sample block holds either a 96-well plate or 48-vial metal insert and is temperature-controlled. You can set it to 4 °C, 10 °C, 15 °C or turn the temperature control off in Compass for iCE software.

Sample cooling turns on when the run starts, and takes a few minutes to reach the temperature setting. After a run, the sample block stays at the set temperature until you open Maurice's door, then it shuts off until you start the next run. This prevents excess condensation.

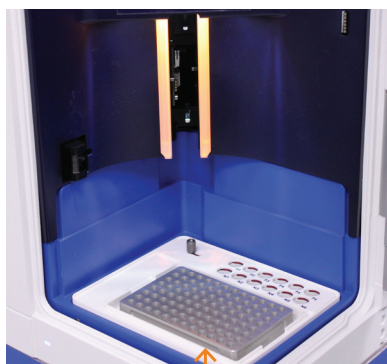
NOTE: Because Maurice holds the sample block temperature after a run until you open the door, samples are still viable for your next run and after overnight runs.



96-well Plate Insert for
Maurice, Maurice C. and Maurice S.



48-vial Insert for
Maurice, Maurice C. and Maurice S.



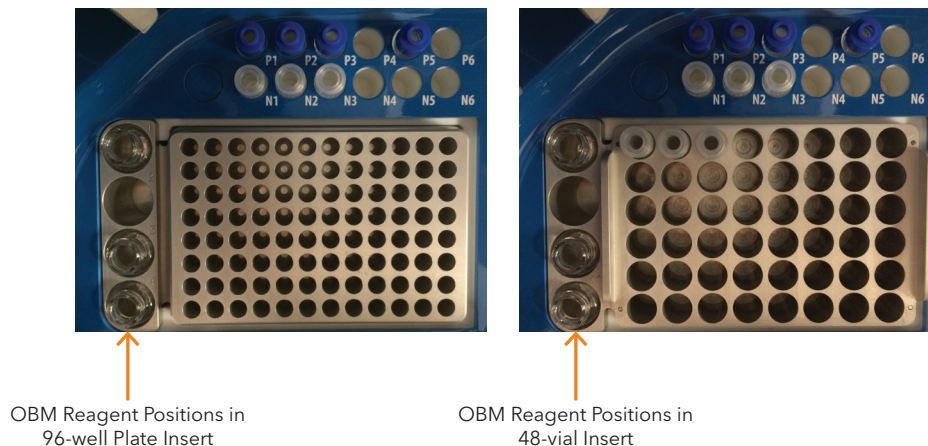
96-well Plate Insert for
MauriceFlex



48-vial Insert for
MauriceFlex

Maurice-OBM and Maurice C.-OBM instruments will also have four additional reagent locations in the 96-well plate and 48-vial metal inserts. These reagents are also temperature controlled at the same setting used for the samples.

- **Column M (left):** These reagents are used for on-board mixing. Only use 6 mL reagent vials in this row.



NOTES:

On-board mixing is only available on Maurice and Maurice C. systems.

When you're using a 96-well plate, well A1 should be in the top left corner of the insert.

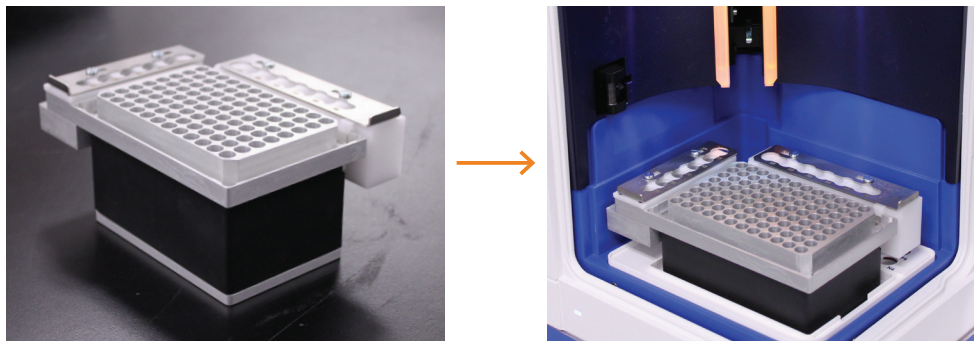
You can only use V-bottom plates with the 96-well plate insert. To ensure performance, we recommend only using 96-well plates from ProteinSimple (PN 046-021).

Remove plate lids before inserting a 96-well plate into Maurice.

You can find info on where to load reagents and samples for cIEF applications in "Step 4: Load Samples and Reagents" on page 99, for CE-SDS PLUS applications in "Step 4: Load Samples and Reagents" on page 299 and for Turbo CE-SDS application in "Step 4: Load Samples and Reagents" on page 368.

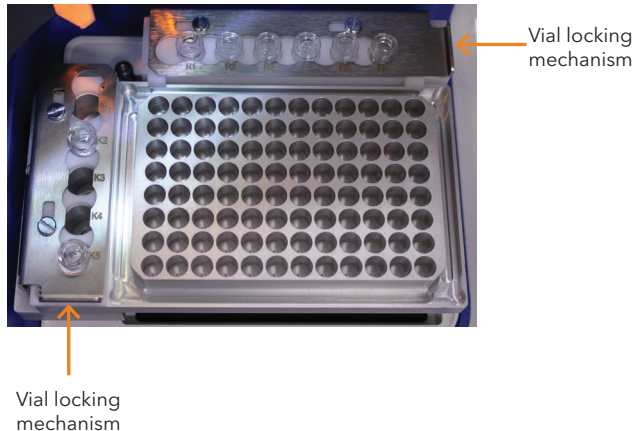
Fractionation Adapter

MauriceFlex uses a fractionation adapter for use with MauriceFlex cIEF or MauriceFlex Fractionation Applications.



The fractionation adapter has one row and one column for batch reagents.

- **Row R (top):** These reagents are kept at room temperature. Only use glass vials with insert, 0.3 mL (PN 110-0018) for the Fluorescence Calibration Standard and 2 mL crimp top glass vials (PN 110-0019) for reagents in this row. Caps are not required on the vials.
- **Column K (left):** These reagents are temperature controlled. Only use 2 mL crimp top glass vials (PN 110-0019) vials for reagents in this row. Caps are not required on the vials.



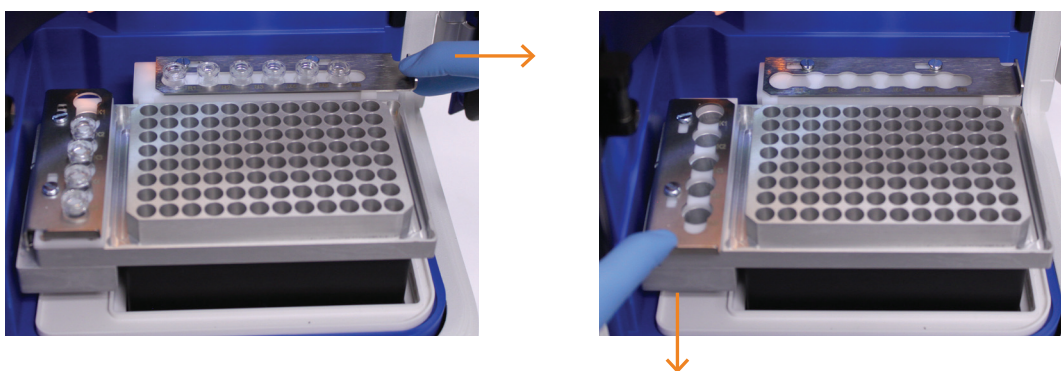
NOTES:

Remove all the vials in rows P and N of the reagent platform before installing the fractionation adapter if a CE-SDS or standard cIEF application was previously performed.

When installing the fractionation adapter, confirm the adapter is seated and sitting flat in MauriceFlex.

Reagent vials on MauriceFlex must be locked in place before starting a batch. The vial locking mechanism default state is locked. Pull the locking mechanism on Row R to the right and the locking mechanism on Column K down to unlock and place vials in the fractionation adapter. Release the locking mechanism to lock the vials in place.

If the fractionation adapter locking mechanism does not spring back upon release, check that all the vials are seated.



Unlocked Row

Unlocked Column

The fractionation adapter can hold a 96-well plate and is temperature-controlled. You can set it to 10 °C, 15 °C or turn the temperature control off in Compass for iCE software.

Sample cooling turns on when the run starts, and takes a few minutes to reach the temperature setting. After a run, the sample block stays at the set temperature until you open MauriceFlex's door, then it shuts off until you start the next run. This prevents excess condensation.

NOTES:

MauriceFlex holds the sample block temperature after a run until you open the door.

When placing the 96-well plate in MauriceFlex, well A1 should be in the top left corner of the insert.

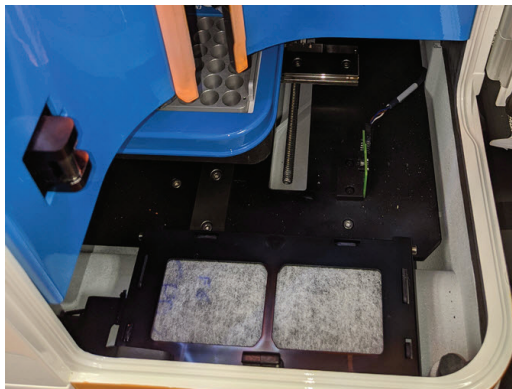
You can only use V-bottom plates with the 96-well plate insert. To ensure performance, we recommend only using 96-well plates from ProteinSimple (PN 046-021).

Remove plate lids before inserting a 96-well plate into MauriceFlex.

You can find info on where to load reagents and samples for MauriceFlex cIEF applications in "Step 2: Load Samples and Reagents" on page 161 and for MauriceFlex Fractionation applications in "Step 2: Load Samples and Reagents" on page 226.

Filter

The optional Maurice System Filter Upgrade adds a fan and filter assembly to the Maurice system to help eliminate odors or vapors that may be present in sample plates from substances such as beta-mercaptoethanol (β ME). The Maurice Filter Kit (PN 046-576) comes with replacement carbon filters and plate lids to cover sample plates when transporting them to and from Maurice.



For more information on filters see “Installing or Replacing the System Filter” on page 795.

Using the Plate Lid

The plate lid is provided for use with Maurice 96-well plates (PN 046-021). We recommend using the lid whenever you run reduced samples on a Maurice CE-SDS PLUS and Turbo CE-SDS batches. For best results, keep the lid on the plate whenever it's not in the Maurice sample block, including after samples are loaded in the wells, when transporting the plate around the lab, and immediately after a batch is completed and the door is opened.

IMPORTANT

Be sure to remove the plate lid before you start a batch! Leaving the lid on the plate will result in damage to the cartridge.

Rear Panel

Located on Maurice's rear panel is the power entry, power switch and network connector.



- **System Power** - The main system power components consist of the power input, fuse and power switch

!WARNING!

Only use the power supply cord provided with Maurice. If the cord is damaged, please contact ProteinSimple Technical Support.

!WARNING!

You can't replace or service any parts on Maurice except the power entry fuse.

!WARNING! SHOCK HAZARD

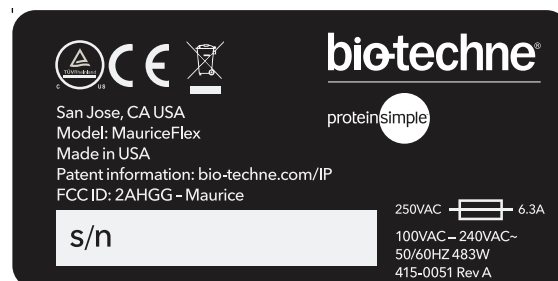
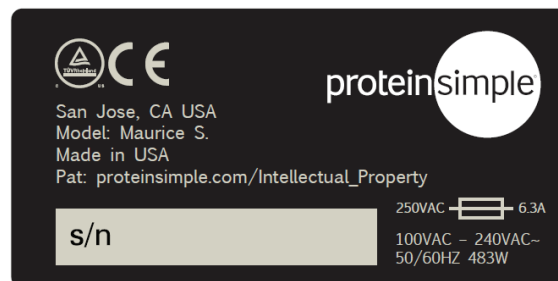
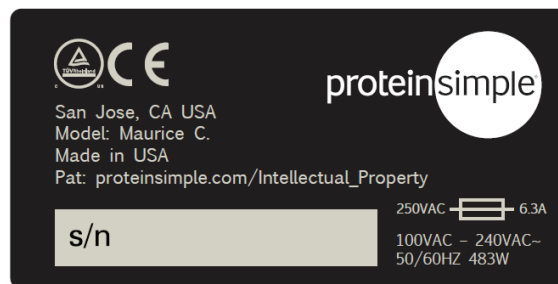
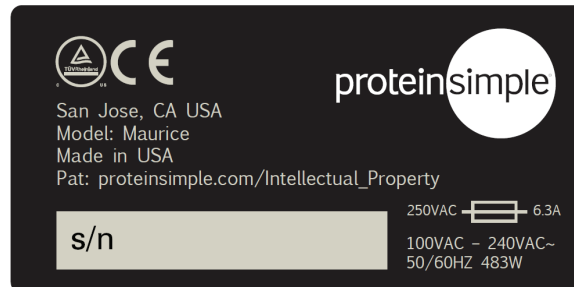
Disconnect the power cord from Maurice's power input to disconnect power to the instrument.

-
- **Network connection** - A 10/100/1000 Mbps Ethernet (RJ-45 connector) is used to connect Maurice to a computer or local network.

NOTE: Serial numbers are used to identify individual instruments.

System Labels

A full system label is located on the rear panel. It includes the ProteinSimple location, system model, power requirements, serial number and certification markings.

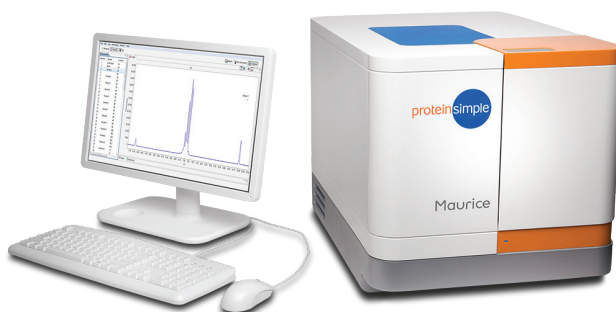


A serial number label is located on the Maurice system's front lower right side, on the system base.



Computer Workstation

The PC has two built-in Ethernet ports, one is used for Maurice and the other is available for your company's network. ProteinSimple configures one port to have a fixed IP for a local link connection to the instrument, the other is configurable by users and will typically use a DHCP for dynamic IP. Set the IP address to 172.30.1.2 subnet 255.255.255.0 to connect to Maurice.



Chapter 4:

Compass for iCE Overview

Chapter Overview

- Launching Compass for iCE
- Compass for iCE Overview
- Software Menus
- Changing the Compass for iCE Main Window Layout
- Viewing the User Guide
- Visiting the Bio-Techne Academy
- Checking for and Installing New Versions of Compass for iCE
- Viewing Release Notes
- Sending Run Files to Technical Support
- Compass for iCE Version Information
- Directory and File Information

Launching Compass for iCE



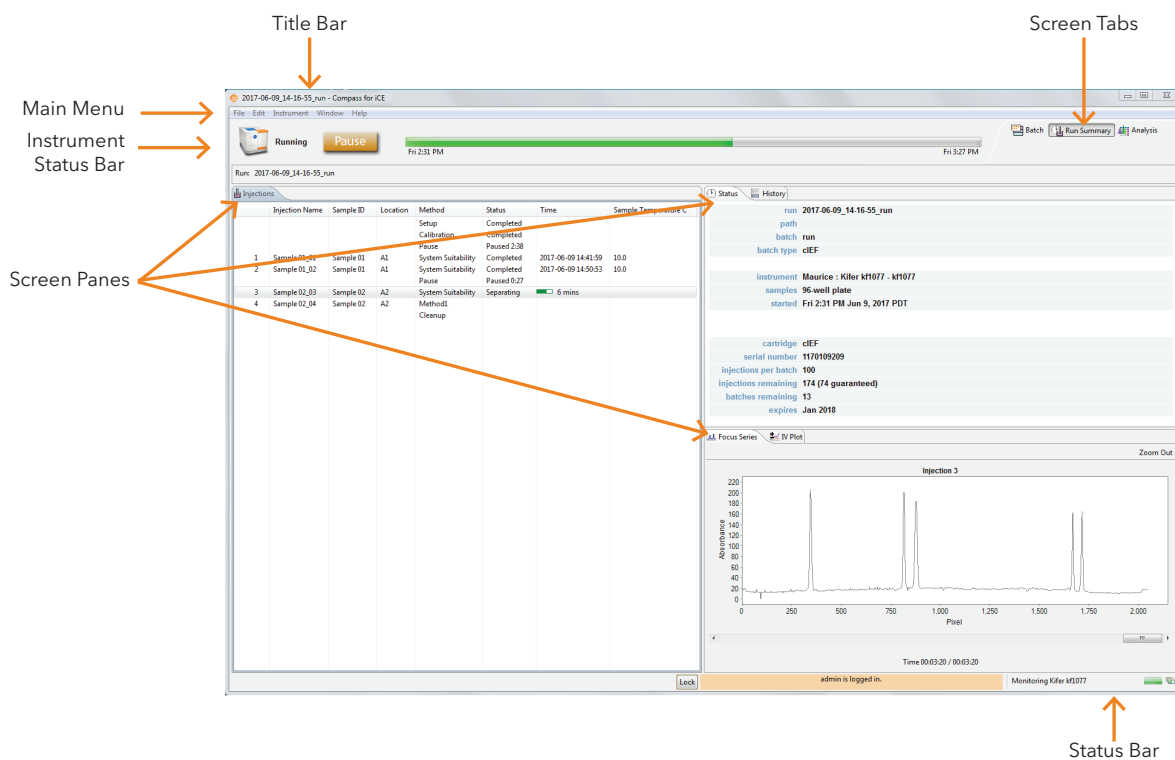
To open Compass for iCE, just double-click the icon on the computer desktop.

Compass for iCE Overview

Compass for iCE has three main screens:

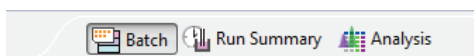
- **Batch** - You'll create and review your batch.
- **Run Summary** - Check out the status of your run.
- **Analysis** - Take a look at the data from your experiment.

Each screen has these components:



Changing the Screen View

To toggle between the Batch, Run Summary and Analysis screens, just click the button in the screen tab located in the upper right corner of the main window.



Batch Screen

The Batch screen is used to create, view, and edit batches. You can assign samples to 96-well plate wells or vials, create and modify methods, customize your injection list and assign methods to each of your injections. The Batch screen is also where you can define how many fractions to collect and which row to start collecting fractions when performing a MauriceFlex Fractionation batch.

Batch: Maurice_CE-SDS_TURBO_Reduced

Layout

10°C

Wash Run Run Run

Injections

Injection Name	Sample ID	Location	Method	Notes
1 Sample 01_01	Sample 01	A1	Reduced IgG	
2 Sample 01_02	Sample 01	A1	Reduced IgG	
3 Sample 01_03	Sample 01	A1	Reduced IgG	
4 Sample 01_04	Sample 01	A1	Reduced IgG	
5 Sample 01_05	Sample 01	A1	Reduced IgG	
6 Sample 01_06	Sample 01	A1	Reduced IgG	
7 Sample 01_07	Sample 01	A1	Reduced IgG	
8 Sample 01_08	Sample 01	A1	Reduced IgG	
9 Sample 01_09	Sample 01	A1	Reduced IgG	
10 Sample 01_10	Sample 01	A1	Reduced IgG	
11 Sample 01_11	Sample 01	A1	Reduced IgG	
12 Sample 01_12	Sample 01	A1	Reduced IgG	
13 Sample 02_13	Sample 02	A2	Reduced IgG	
14 Sample 02_14	Sample 02	A2	Reduced IgG	
15 Sample 02_15	Sample 02	A2	Reduced IgG	
16 Sample 02_16	Sample 02	A2	Reduced IgG	
17 Sample 02_17	Sample 02	A2	Reduced IgG	
18 Sample 02_18	Sample 02	A2	Reduced IgG	
19 Sample 02_19	Sample 02	A2	Reduced IgG	
20 Sample 02_20	Sample 02	A2	Reduced IgG	
21 Sample 02_21	Sample 02	A2	Reduced IgG	
22 Sample 02_22	Sample 02	A2	Reduced IgG	
23 Sample 02_23	Sample 02	A2	Reduced IgG	
24 Sample 02_24	Sample 02	A2	Reduced IgG	

Methods

Name	Sample Load	Separation
Reduced IgG	8 sec 3500 Volts	5.5 min 4200 Volts
Non-reduced IgG	8 sec 3500 Volts	8.0 min 4200 Volts
MW Markers	8 sec 3500 Volts	8.0 min 4200 Volts

DemoData_Maurice cIEF - Compass for iCE

File Edit Instrument Window Help

Batch: DemoData_Maurice cIEF

Layout

10°C

MC FI Cal Water Empty

Water

1 2 3 4 5 6 7 8

A B C D E F

Methods

Injection Name	Sample ID	Location	Method	Notes
1 System Suit_01	System Suit	A1	Method1	
2 Peptide Mix_02	Peptide Mix	A2	Method1	
3 mAb_03	mAb	A3	Method2	
4 System Suit_04	System Suit	A1	Method1	
5 Peptide Mix_05	Peptide Mix	A2	Method1	
6 mAb_06	mAb	A3	Method2	

Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	90	3.38, 10.17		
Method2	1.0 min 1500 Volts, 6.0 min 3000 Volts	6 Exposures	90	4.05, 9.99		

MauriceFlex cIEF - Compass for iCE

File Edit Instrument Window Help

Batch: MauriceFlex cIEF

Layout

10°C

MC FI Cal Water Water Water Empty

Sample Catholyte Water

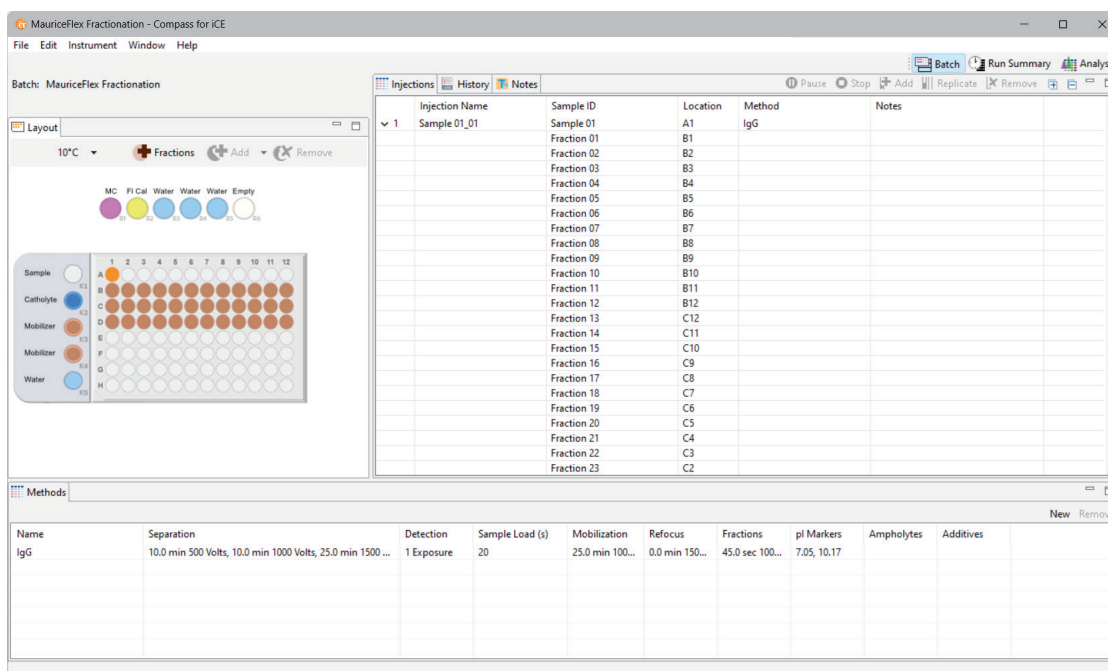
1 2 3 4 5 6 7 8 9 10 11 12

A B C D E F G H

Methods

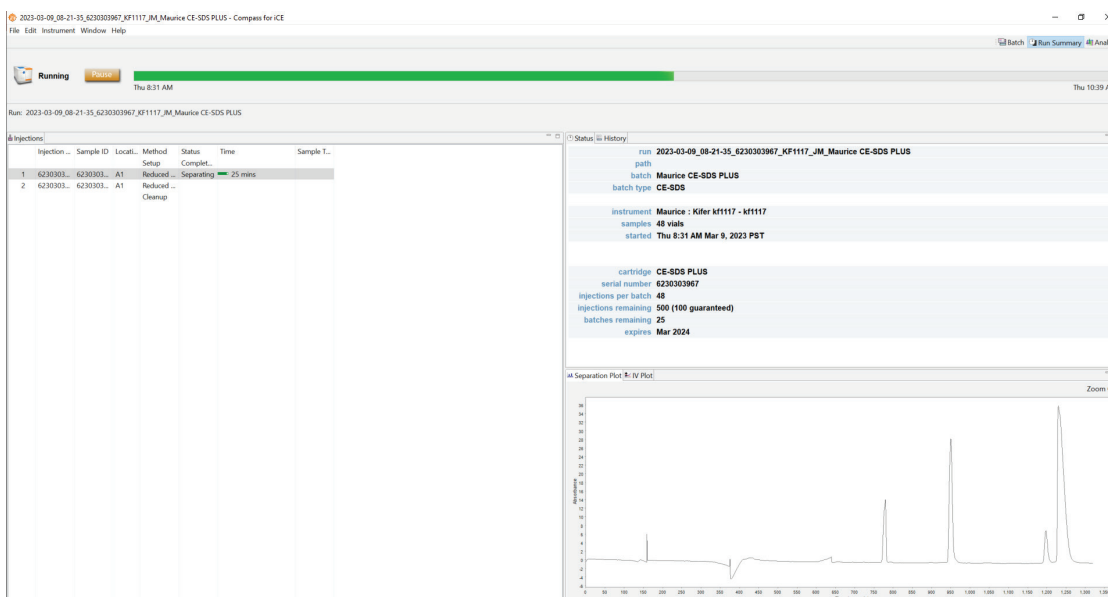
Injection Name	Sample ID	Location	Method	Notes
1 Sample 01_01	Sample 01	A1	IgG	
2 Sample 02_02	Sample 02	A2	IgG	
3 Sample 03_03	Sample 03	A3	IgG	
4 Sample 04_04	Sample 04	A4	IgG	

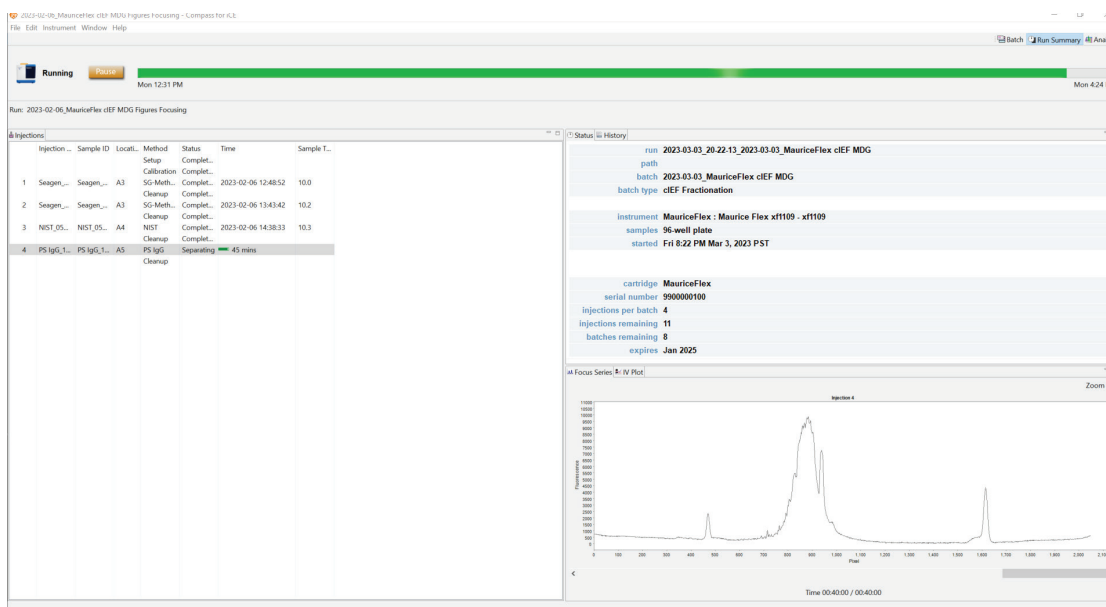
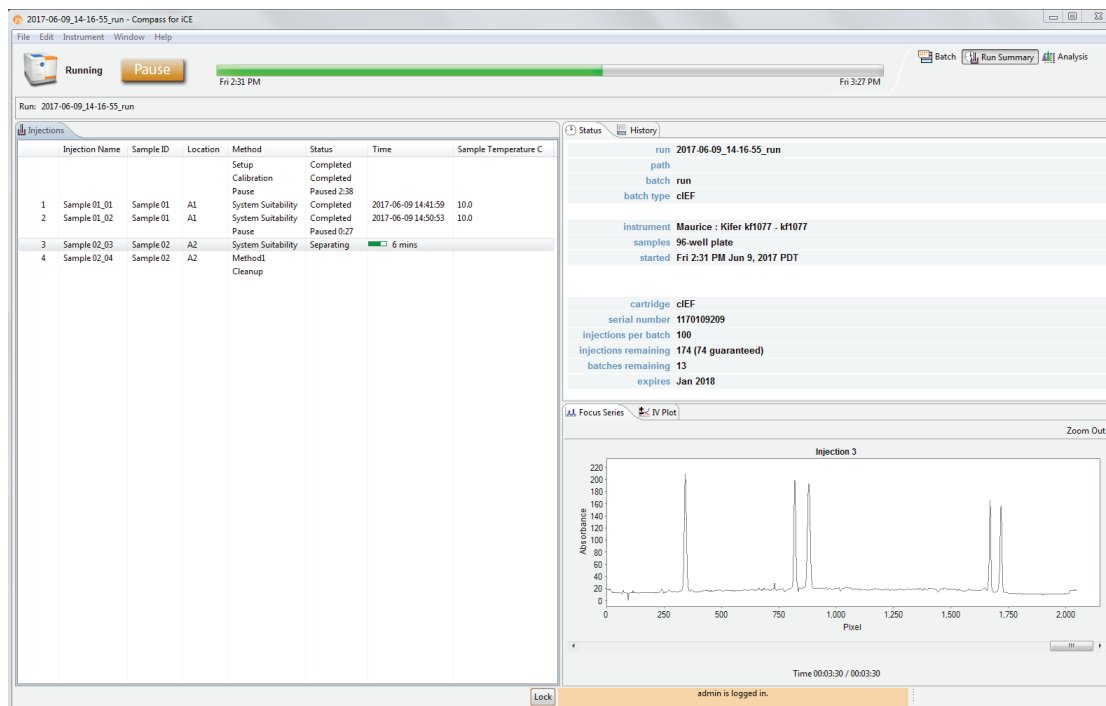
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25.0 min 1500 ...	1 Exposure	20	7.05, 10.17		

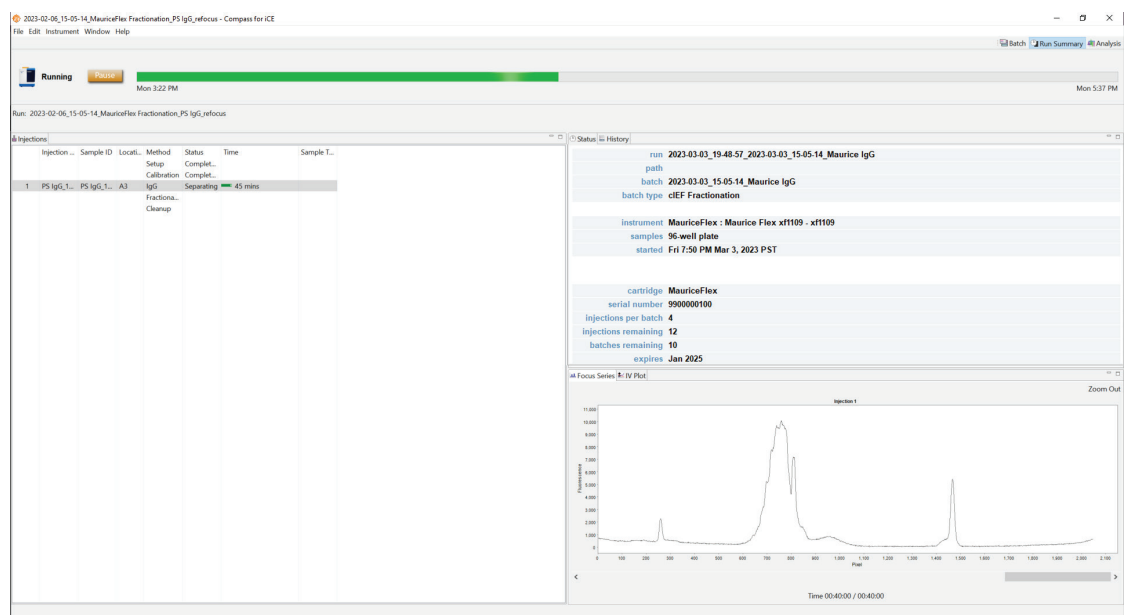


Run Summary Screen

The Run Summary screen is used to monitor status of a batch in progress, view the CE-SDS Separation plot or cIEF, MauriceFlex cIEF and MauriceFlex Fractionation Focus series for each injection and view the current and voltage plots for each injection.

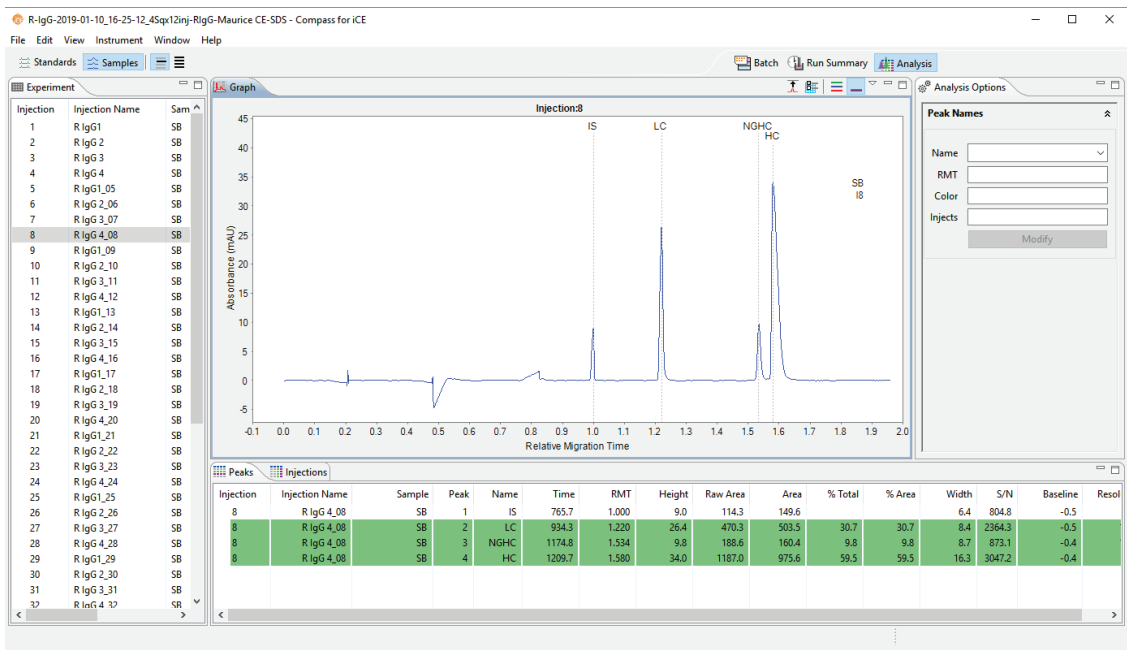


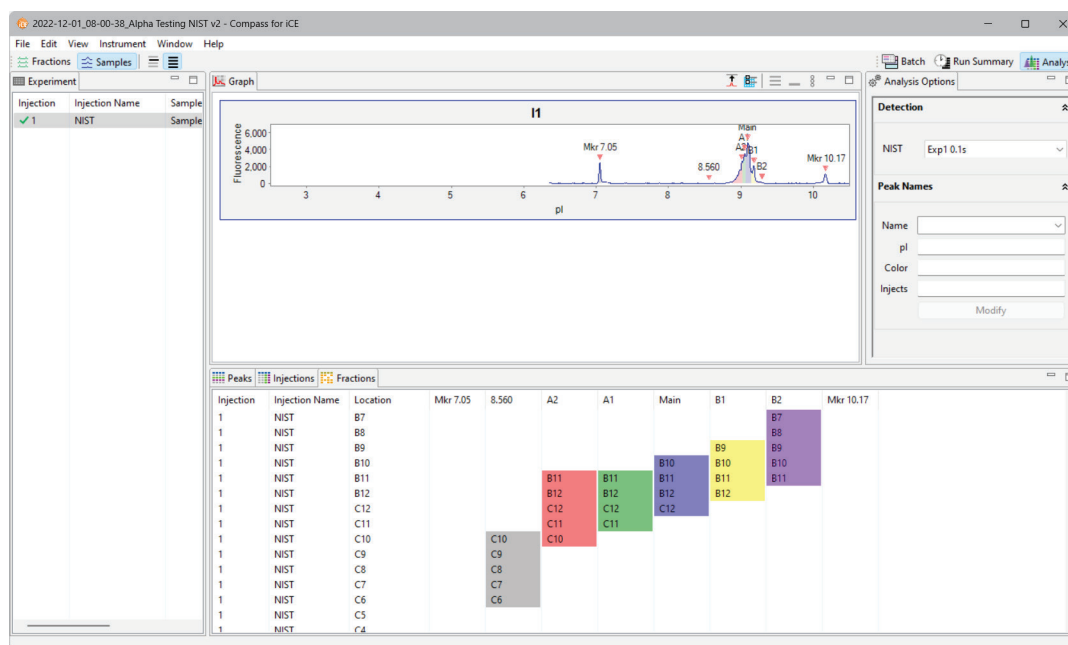
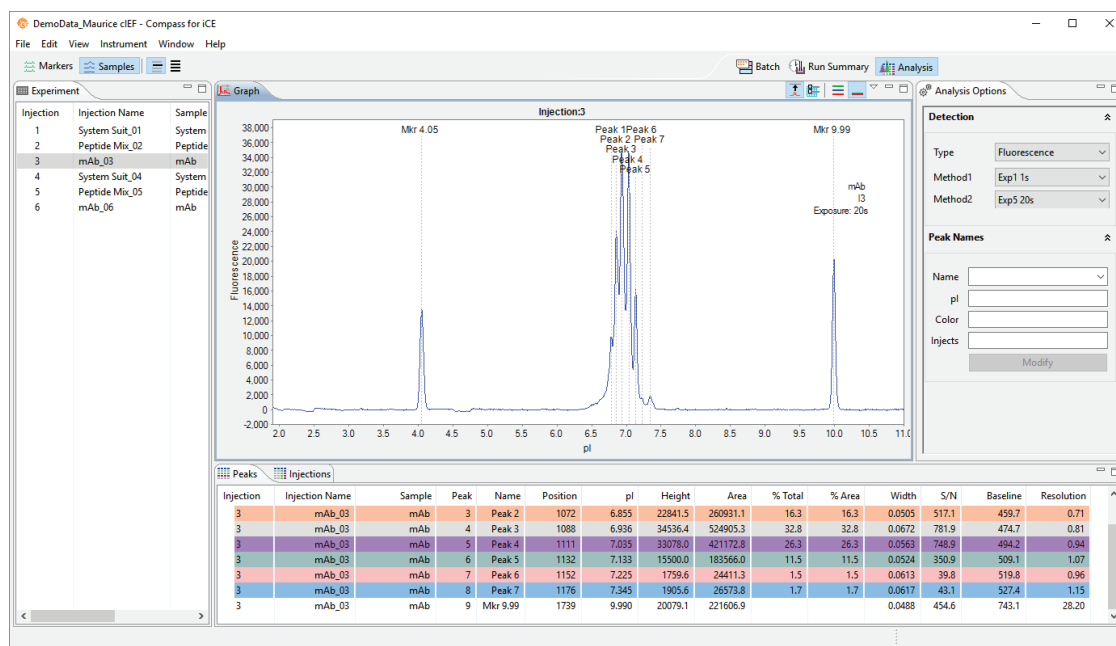




Analysis Screen

The Analysis screen is used to view data from your batch, including the graph view (electropherograms) and a table with your results, and analyze data for completed runs. For MauriceFlex Fractionation batches, you can also view a table predicting where each sample peak is collected into a 96-well plate.

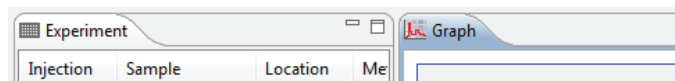




Screen Panes

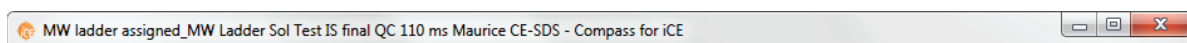
Each of the Batch, Run Summary and Analysis screens have multiple panes that let you view the individual components of a batch, method or data file. Each pane has a labeled tab and a unique icon. We'll describe panes specific to each screen later in the individual screen sections.

The active pane in a screen is blue. To view a pane, click in the pane or on its tab. The screen panes can be minimized and moved. To return to the default view, select **Default Layout** in the Window menu. The example below shows panes in the Batch screen, and the Graph pane is active:



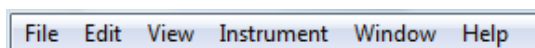
Title Bar

In the title bar you will see the batch file name and the icons that allow the main Compass for iCE window to be minimized, maximized or closed.



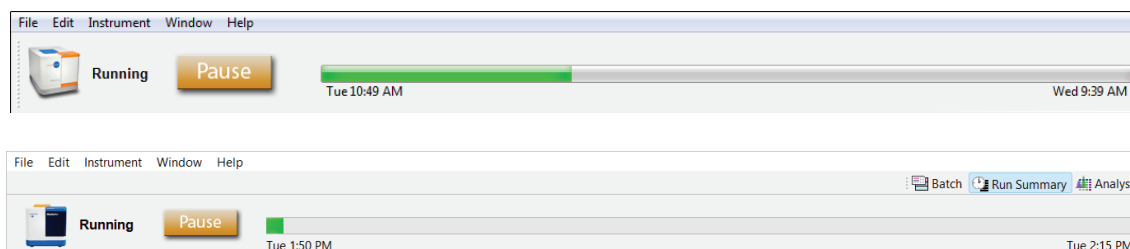
Main Menu

Access to various software, instrument and screen operations is available through the main menu. More details on menu commands can be found in “Software Menus” on page 44 .



Instrument Status Bar

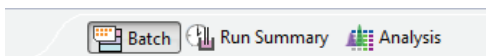
The instrument status bar is used to start batches and cleaning protocols, indicate system status and show run progress. More details on instrument control and status can be found in Chapter 17: “Controlling Maurice, Maurice C., Maurice S. and MauriceFlex”.



NOTE: You will only see the instrument status bar when Compass for iCE is connected to an instrument. There is no status bar on computer workstations that you’re only using for data analysis.

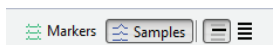
Screen Tab

The screen tab lets you move between Batch, Run Summary or Analysis screens and is located in the upper right corner of the main window. Just click a button to view a screen.



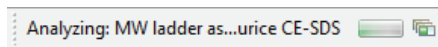
View Bar

The view bar is only displayed in the Analysis screen as part of the main menu, and allows you to switch between viewing standards or sample data for a CE-SDS batch, markers or sample data for a cIEF batch and fractions or sample for a MauriceFlex Fractionation batch. You can also switch between data for a single injection or all injections in the batch, or grouped injection data. View bar options are in “Analysis Options Pane” on page 483 for cIEF and MauriceFlex cIEF applications, page 671 for CE-SDS applications or page 580 for MauriceFlex fractionation application.



Compass for iCE Status Bar

The status bar is in the lower right corner of the main window. It displays active software processes and their progress.

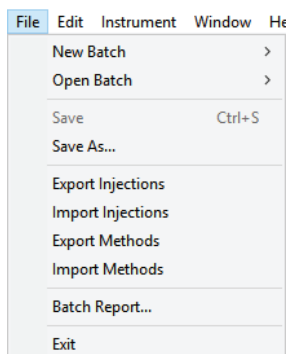


Software Menus

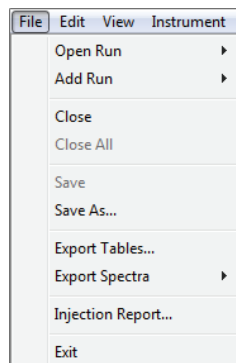
Some of the items in the Compass for iCE main menu are available in specific screens only, and menu commands change depending on which screen is active. You can find menus and commands available for each screen in the Chapter 5: “cIEF Batches”, Chapter 7: “MauriceFlex cIEF Batches”, Chapter 9: “MauriceFlex Fractionation Batches”, Chapter 11: “CE-SDS PLUS Batches”, Chapter 13: “Turbo CE-SDS Batches”, Chapter 16: “Run Status”, Chapter 18: “cIEF Data Analysis”, Chapter 20: “CE-SDS Data Analysis”, Chapter 19: “MauriceFlex Fractionation Data Analysis”.

File Menu

The File menu contains basic file commands. The file menu options change when you change between the Batch View or Run Summary/Analysis tab.



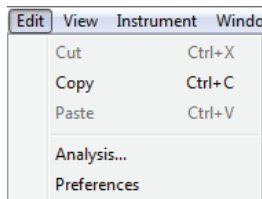
File Menu: Batch View



File Menu: Run Summary/Analysis

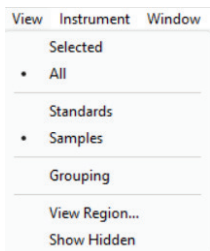
Edit Menu

The Edit menu contains basic editing commands, analysis and preferences options. Specific details on preferences are described in Chapter 21: “Setting Your Preferences”.

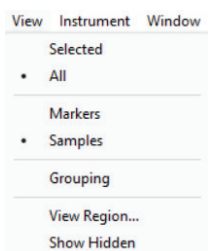
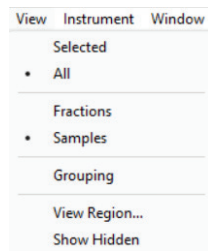


View Menu

The View menu is only available in the Analysis screen, and allows you to change how your data is displayed. For more info on view options check out “Analysis Options Pane” on page 483 for cIEF applications, page 671 for CE-SDS applications or page 580 for Maurice Fractionation and “Using Groups” on page 498 for cIEF applications or page 684 for CE-SDS applications.

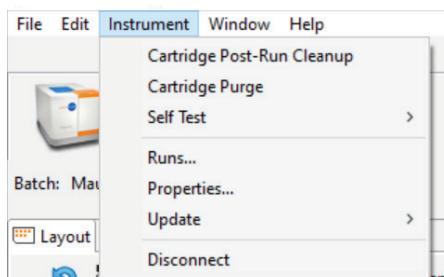


CE-SDS Application

Standard cIEF and
MauriceFlex cIEF ApplicationMauriceFlex Fractionation
Application

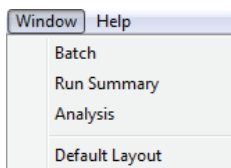
Instrument Menu

The Instrument menu is only available when the software is connected directly to your instrument. You can learn more about instrument control options in Chapter 17: “Controlling Maurice, Maurice C., Maurice S. and MauriceFlex”.



Window Menu

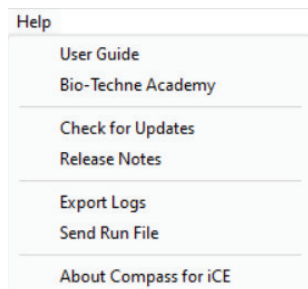
The Window menu lets you to switch between the Batch, Run Summary or Analysis screens, and restore screens to the default layout.



- **Batch** - Displays the Batch screen where you create, view, and edit batches.
- **Run Summary** - Displays the Run Summary screen which lets you view the status of a batch in progress.
- **Analysis** - Displays the Analysis screen that lets you view electropherograms and results and change analysis parameters
- **Default Layout** - Restores the individual panes in the current screen back to their default size and location.

Help Menu

The Help menu gives you access to Help, software updates, release notes and other software info.



- **User Guide** - Displays the User Guide for Maurice, Maurice C., Maurice S, and MauriceFlex.
- **Bio-Techne Academy** - Redirects you to the online Bio-Techne Academy where you can watch online training videos for Maurice, register for online workshops and access product collateral.
- **Check for Updates** - Automatically checks to see if a new version of Compass for iCE is available.
- **Release Notes** - Displays the software release notes for the current and prior versions.
- **Export Log** - Sends a zipped log file directly to ProteinSimple Technical Support.
- **Send Run File** - Sends a zipped run file directly to ProteinSimple Technical Support.
- **About Compass for iCE** - Displays the software version and build information.

Changing the Compass for iCE Main Window Layout

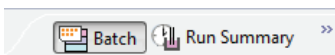
You can easily resize the main window and the individual panes in each screen. Screen panes can also be moved outside of the main window.

Resizing the Main Compass for iCE Window

To resize the main window, roll the mouse over a corner or border until the sizing arrow appears. Then just click and drag to resize.

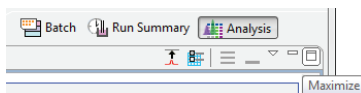
Resizing the Screen Tab

The screen tab can be sized to show all or just some of the screen buttons. To resize, roll the mouse over the left edge of the tab until the sizing arrow appears, then click and drag to resize. If a screen button is hidden, a double arrow will display in the tab. Just click to display and select the hidden screen.

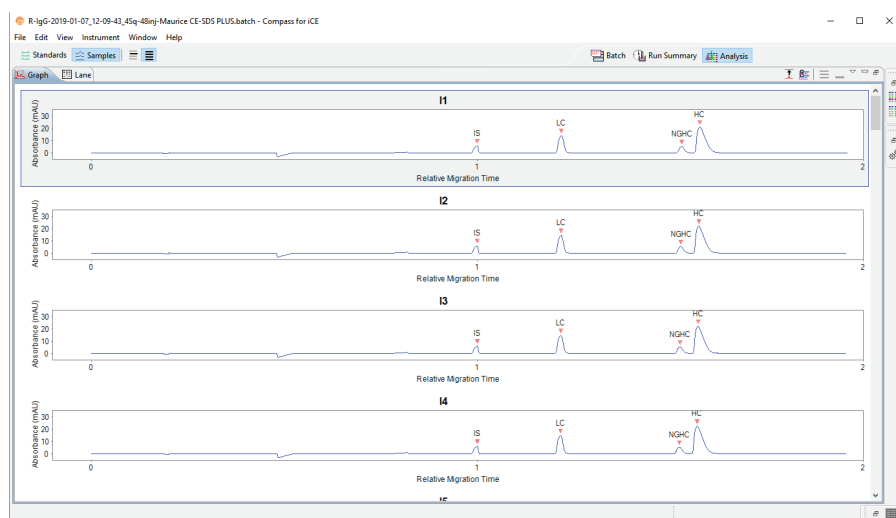


Resizing Screen Panes

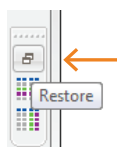
- **To resize a pane** - Roll the mouse over the pane border until the sizing arrow appears. Then just click and drag to resize.
- **To maximize a pane** - Click the maximize button in the upper right corner or double-click the tab.



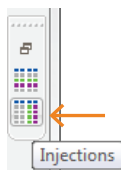
The other panes in the screen will automatically minimize to pane bars in the task area along the window border.



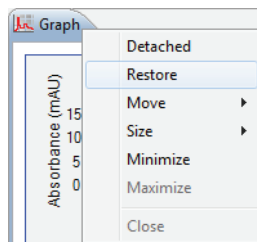
- **To restore all minimized panes** - Click Restore on the minimized pane bar.



- **To restore only one minimized pane** - Click the pane icon on the minimized pane bar.



- To restore a maximized pane to its original size - Double-click the tab or right click the tab and click **Restore**.



- To restore all panes to their original sizes - Select **Window** in the main menu and click **Default Layout**.

Changing the Location of Screen Panes

Panes can be moved to different locations within a screen.

- To move a pane - Click on its tab and drag it to the new location. As the pane is moved, area guides will display to assist you in choosing a drop location.



Area guides with a black arrow let you know that if the pane is dropped at that location, it will be resized and relocated as an individual pane in that area of the screen.

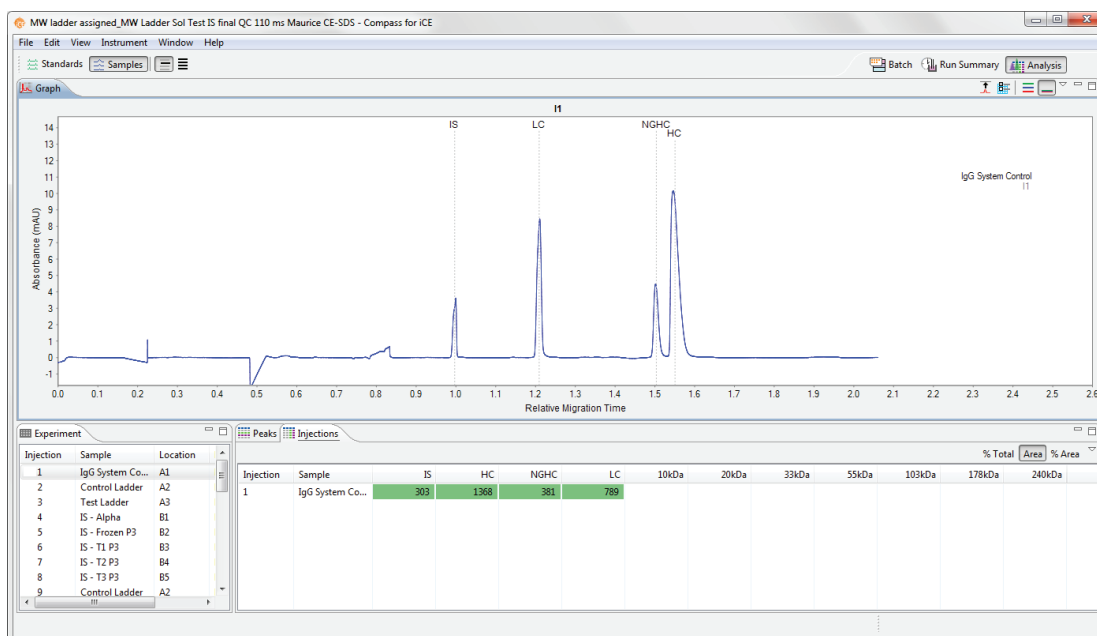


Area guides with a folder let you know that if the pane is dropped at that location, it will be added as a new tab in an area with one or more pane tabs.

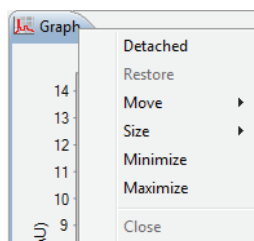


Area guides with a window let you know that if the pane is dropped at that location, it will be a separate window outside the Compass for iCE main window.

This example shows the Analysis screen after moving the Graph pane:



- **To detach a pane from the main window** - Click on its tab and drag it outside the main Compass for iCE window or right click the tab and click **Detached**.



- **To move a detached pane back inside the main window** - Right click the tab and deselect **Detached**.
- **To restore all panes to their original locations** - Select **Window** in the main menu and click **Default Layout**.

Restoring the Main Window to the Default Layout

To restore screen pane sizes and locations to the original Compass for iCE layout, select **Window** from the main menu and click **Default Layout**.

Viewing the User Guide

Select **Help** and click **User Guide** to view Maurice Systems User Guide.

- If the computer you're using has an internet connection, the latest online version of the User Guide PDF will display.
- If your computer doesn't have an internet connection, you'll first need to download the User Guide PDF from another computer or contact tech support for a copy. Upload the PDF file to the C:\Program Files\Compass for iCE folder to access the User Guide from the Help menu.

Visiting the Bio-Techne Academy

Select **Help** and click **Bio-Techne Academy** to get redirected to this online resource. Registered users can view online training videos, registers for online workshops or watch previous workshops and view instrument collateral. Training videos in Chinese, Korean, Japanese and French are also available. Enter Maurice in the search window to view Maurice-related content only.

Checking for and Installing New Versions of Compass for iCE

The software can automatically check to see if a newer version of software is available. To do this:

1. Make sure the computer being used has an active internet connection.
2. Select **Help** and click **Check for Updates**. If an update is found, a screen will display with the new version that's available.

3. Click **Finish** to start the download and install the update.
4. Follow the on-screen instructions to complete the software installation.
5. Reboot the computer before using the new version of software.

Viewing Release Notes

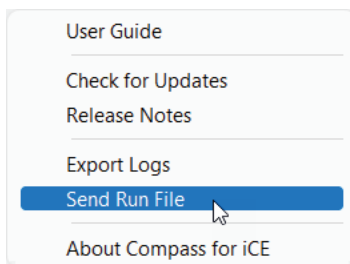
Select **Help** and click **Release Notes** to view feature updates and bug fixes for new and past versions of Compass for iCE. If the computer you're using has an internet connection, the latest online version of the release notes PDF will display. When an internet connection isn't available, the release notes PDF shipped with the original installer for the software will open instead. We recommend you review these notes whenever a software update is installed.

NOTE: You can contact ProteinSimple Technical Support to request the release notes for new versions of Compass for iCE before you install it.

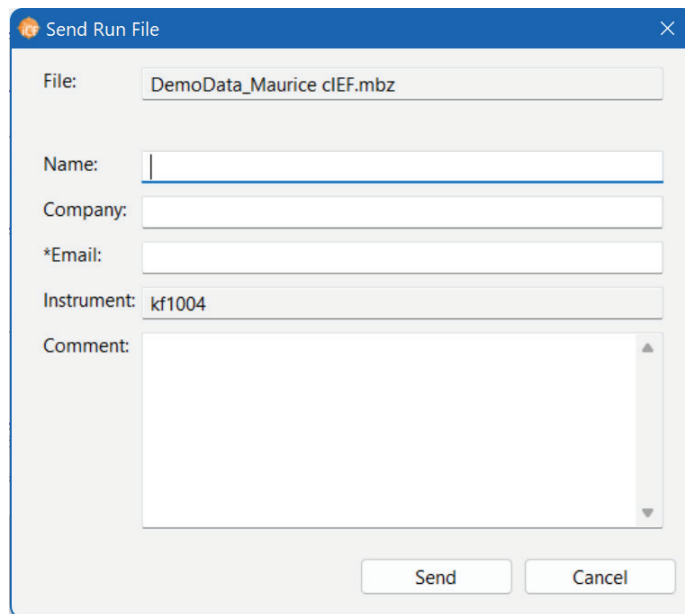
Sending Run Files to Technical Support

If the computer you're using has an internet connection, Compass for iCE can zip and send a run file directly to ProteinSimple Technical Support.

1. Open the run file in Compass for iCE.
2. Select **Help** in the main menu and click **Send Run File**:



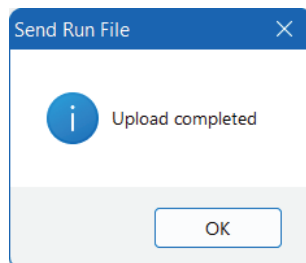
3. Enter your **Name**, **Company** and **E-mail** address, and any details in the **Comments** section:



The 'Send Run File' dialog box contains the following fields and controls:

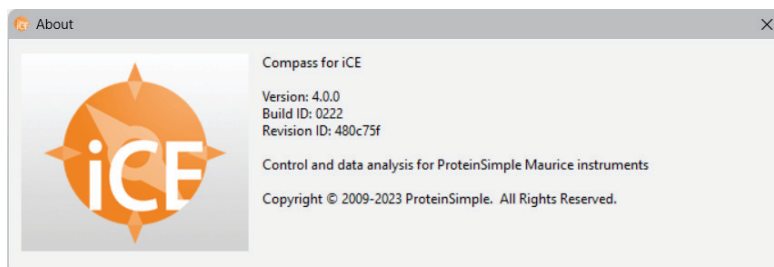
- File:** A text field containing 'DemoData_Maurice cIEF.mbz'.
- Name:** An empty text field.
- Company:** An empty text field.
- *Email:** An empty text field.
- Instrument:** A text field containing 'kf1004'.
- Comment:** A large, empty text area with a vertical scrollbar.
- Buttons:** 'Send' and 'Cancel' buttons at the bottom right.

4. Click **Send**. When the file upload is complete, the following message displays. Click **OK**.



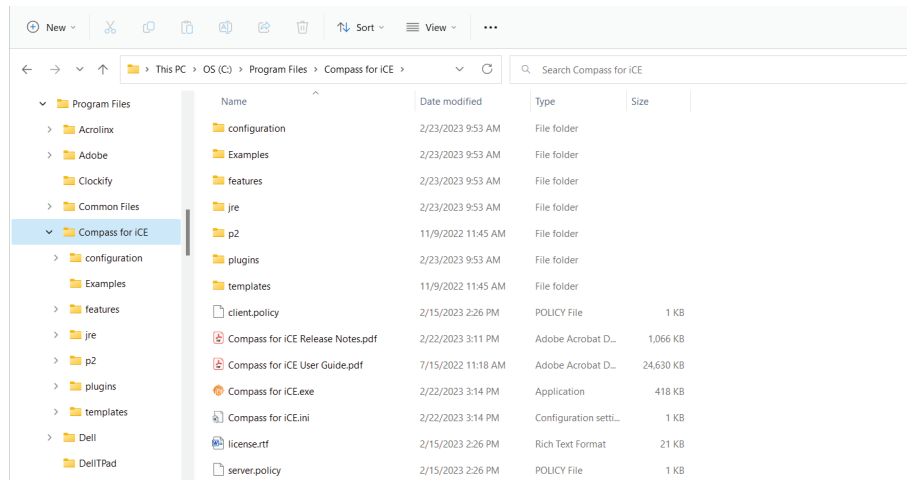
Compass for iCE Version Information

Select **Help** and click **About Compass for iCE** to view the software version and build number information.

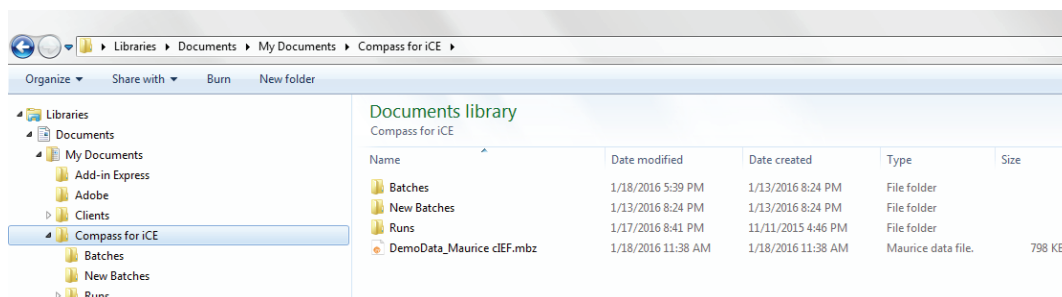


Directory and File Information

The main Compass for iCE directory is located in the **Program Files** folder, and also contains PDF files of the User Guide for Maurice, Maurice C., Maurice S. and MauriceFlex.



Batch and run files are located in the **Documents** folder in the User directory on your computer:



- **Batches Folder** - Contains all batch files that you've saved.
- **New Batches Folder** - Contains Maurice batch template files.
- **Runs Folder** - Contains all batch data files. Data is automatically written to this folder.

NOTE: When a Compass for iCE software update is performed, the template s in the New Batch folder are overwritten. If you have customized these batches, we recommend saving them in a unique subfolder prior to updating the software, then transferring them back to the New Batch folder after the update to avoid losing your customizations.

File Types

These file types are used by Compass for iCE:

- **Batch Files** - Use a *.batch file extension.
- **Run Files** - Use a *.mbz file extension. The default file format for run files is Date_Time_BatchName. An example run file name would be 2016-01-28_18-50-53_CE-SDS.mbz.
- **Analysis Settings Files** - Exported analysis settings files use a *.settings file extension.

Chapter 5:

cIEF Batches

Chapter Overview

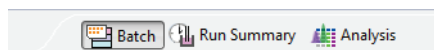
- Overview
- Batch Screen Overview
- Opening a Batch
- Creating a New Batch
- Viewing Replicate Injections
- On-board Mixing (Maurice and Maurice C.)
- Batch History
- Making Changes to a Batch
- Viewing and Editing Batches in Completed Runs
- Copying and Pasting Injection Names or Sample IDs from other Documents
- Importing and Exporting Injections
- Importing and Exporting Methods
- Batch Reports

Overview

Standard cIEF batches can be run on Maurice, Maurice C. or MauriceFlex using a cIEF cartridge.

Batch Screen Overview

You can use the Batch screen to create, view and edit batches. To get to this screen, click the **Batch** screen tab:



Batch Screen Panes

The Batch screen has five panes:

- **Layout** - Displays a map of either the 96-well plate or 48-vial tray of batch sample locations.
- **Injections** - Lists the injections, sample ID, sample locations and methods that Maurice, Maurice C. or MauriceFlex will execute for each sample in the batch.
- **History** - Lists all batch file events from initial creation to the most current update.
- **Notes** - Lets you enter specific information about your batch.
- **Methods** - Lets you create methods and enter method parameters used in the batch.

The screenshot shows the Batch screen interface with the following panes:

- Layout Pane:** Displays a 96-well plate map. The top row (A) is highlighted, showing wells 1 through 8. The temperature is set to 10°C. Buttons for 'Add' and 'Remove' are visible.
- Injections Pane:** A table listing injections with columns: Injection Name, Sample ID, Location, Method, and Notes.

Injection Name	Sample ID	Location	Method	Notes
System Suitability_01	System Suitability	A1	System Suitability	
mAb 11 Blank_02	mAb 11 Blank	A2	mAb Method	
mAb 11 Ref. Std_03	mAb 11 Ref. Std.	A3	mAb Method	
mAb 11 Prep 20160121_04	mAb 11 Prep 20160121	A4	mAb Method	
mAb 11 Prep 20160121_05	mAb 11 Prep 20160121	A4	mAb Method	
mAb 11 Prep 20160121_06	mAb 11 Prep 20160121	A4	mAb Method	
mAb 11 Ref. Std_07	mAb 11 Ref. Std.	A3	mAb Method	
mAb 11 Blank_08	mAb 11 Blank	A2	mAb Method	
- Methods Pane:** A table listing methods with columns: Name, Separation, Detection, Sample Load (s), pI Markers, Ampholytes, and Additives.

Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
System Suitability	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17		
mAb Method	1.0 min 1500 Volts, 6.0 min 3000 Volts	5 Exposures	55	4.05, 9.99		

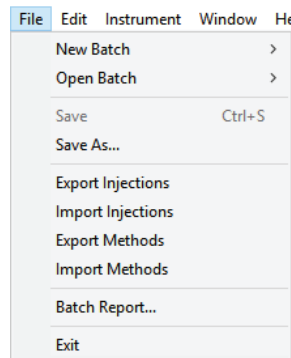
Software Menus Active in the Batch Screen

These main menu items are active in the Batch screen:

- File
- Edit
- Instrument (when the software is connected to Maurice, Maurice C. or MauriceFlex)
- Window
- Help

File Menu

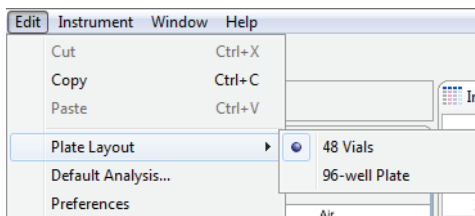
These File menu options are active:



- **New Batch** - Creates a new batch from a starter template.
- **Open Batch** - Opens an existing batch.
- **Save/Save As** - Saves the open batch.
- **Export Injections** - Exports injections from the current batch as a .csv file.
- **Import Injections** - Imports injections into the current batch from a .csv file.
- **Export Methods** - Exports method(s) from the current batch as separate files.
- **Import Methods** - Imports saved method(s) into the current batch.
- **Batch Report** - Exports a table of sample and method details for each injection in the batch as a PDF file.
- **Exit** - Closes Compass for iCE.

Edit Menu

The following Edit menu options are active in the Batch screen:



- **Cut** - Cuts the information currently selected.
- **Copy** - Copies the information currently selected.
- **Paste** - Pastes the copied information.

NOTE: Cut, Copy and Paste work differently depending on the pane you're in. How and when you can use them is described throughout the chapter.

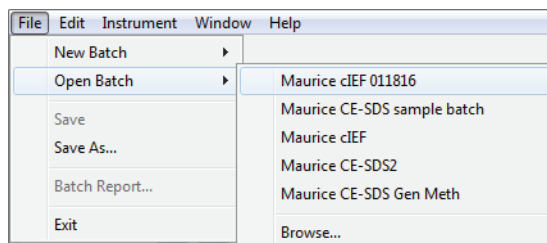
- **Plate Layout** - Lets you select between a 96-well plate or a 48-vial tray to run samples in.
- **Default Analysis** - Displays the default settings that will be used to analyze the data generated with your batch.
- **Default Analysis View** - Displays the default settings that will be used to view the data generated with your batch.
- **Preferences** - Lets you set and save your preferences for Access Control, data export, graph colors, grouped data and Twitter settings. See Chapter 13: "Setting Your Preferences" for more information.

Opening a Batch

NOTE: cIEF batches that use the FL458 nm fluorescence detection filter can only be opened in Compass for iCE 2.1 or later.

To open an existing batch:

1. Select **File** in the main menu and click **Open Batch**.



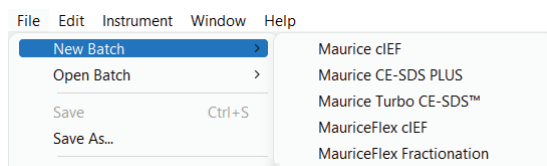
2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file. Select one of those or click **Browse** to open the Batches folder and select a different one.
3. To make changes to the batch, see the steps in “Creating a New Batch” on page 59. When you’re done, select **File** from the main menu and click **Save**.

Creating a New Batch

To create a new batch, we recommend that you use one of the template methods and make any necessary modifications from there.

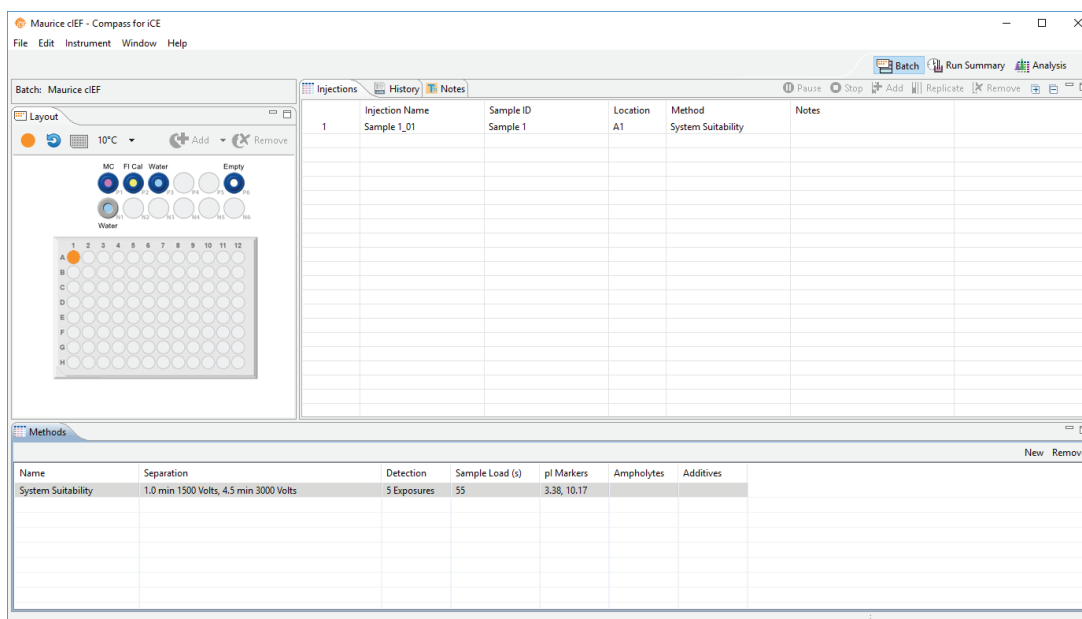
Step 1 - Open a Template Batch

1. Select **File** in the main menu and click **New Batch**:



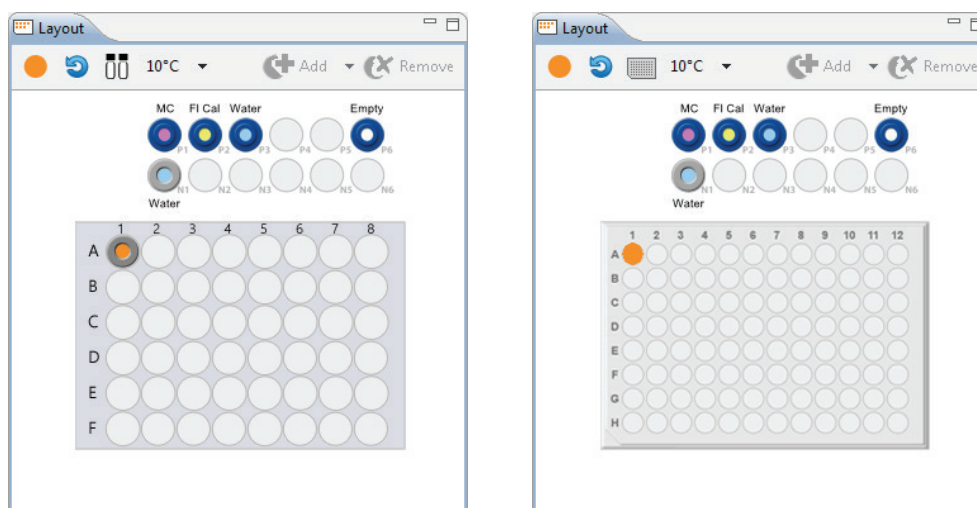
NOTE: Only template batches specific to your Maurice will be available when you are connected to the instrument. Template batches will not appear in the menu if that application type cannot be performed on your instrument.

2. Select **Maurice cIEF**. A batch using the default method will display.



Step 2 - Assign Your Samples

The Layout pane shows the default locations of where batch reagent should be placed in your Maurice system's samples and reagents platform, and a map of either the 96-well plate or the 48-vial tray used for samples.



The same reagent locations are used for every batch:

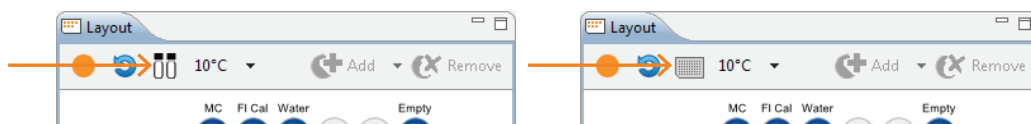
- **P1** - 0.5% Methyl Cellulose with **blue pressure cap**
- **P2** - Fluorescence Calibration Standard with **blue pressure cap**
- **P3** - Water vial with **blue pressure cap**
- **P6** - Empty vial (air) with **blue pressure cap**
- **N1** - Water vial with **clear screw cap**

If you are using Maurice-OBM or Maurice C.-OBM to do on-board mixing:

- **N2** - Water vial with **clear screw cap**
- **N3** - Empty vial with **clear screw cap**
- **M1** - 6 mL IEF Separation Mix vial
- **M2** - N/A for 48 samples or less, add a 6 mL IEF Separation Mix vial when running more than 48 samples
- **M3** - 6 mL water vial
- **M4** - 6 mL water vial

NOTE: On-board mixing is only available on Maurice and Maurice C. systems.

1. To assign samples, select 48 vials or a 96-well plate depending on what you're running. Clicking on the vial/plate icon toggles between formats.



2. Select your samples:

To import samples using a saved injections file:

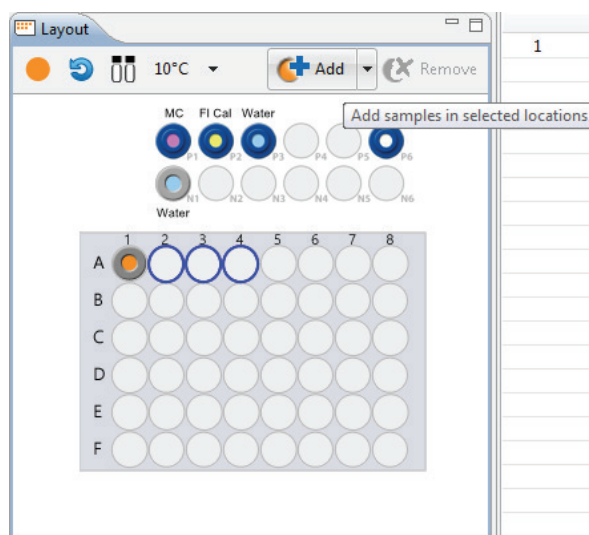
- a. Select **File** in the main menu and click **Import Injections**.
- b. Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.
- c. Skip to step 3 on page 63.

NOTE: For more information on setting up an injections file for import, see “Importing and Exporting Injections” on page 84

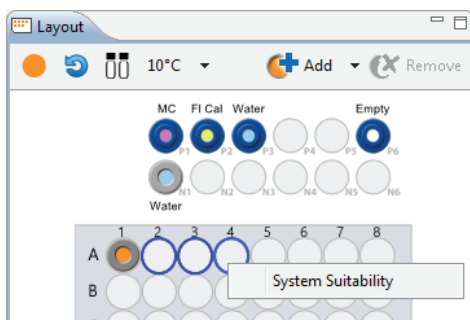
To select samples manually:

- **Add samples and select methods later:** Use your mouse to highlight the well or vial positions your samples are located in, then click **Add**. For this example we're using vials.

NOTE: The template batch automatically adds a sample in well or vial A1 by default, but if you don't have a sample in this position you can remove it after you've added new positions for your samples.



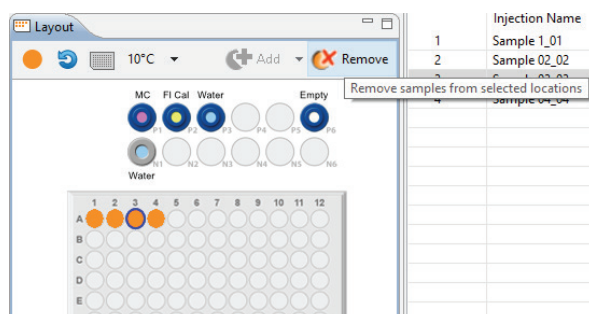
- **Add samples with preassigned methods:** Highlight the well or vial positions your samples are located in, then right-click and select a method.



Either option of adding samples populates the Injections table:

Injections History Notes						Pause	Stop	Add	Replicate	Remove		
	Injection Name	Sample ID	Location	Method	Notes							
1	Sample 1_01	Sample 1	A1	System Suitability								
2	Sample 02_02	Sample 02	A2	System Suitability								
3	Sample 03_03	Sample 03	A3	System Suitability								
4	Sample 04_04	Sample 04	A4	System Suitability								

To remove samples, just highlight the ones you want to remove and click **Remove**. They'll also be removed in the Injections table. You can also right-click well(s) or vial(s) in the Layout pane and click **Remove Sample(s)**.

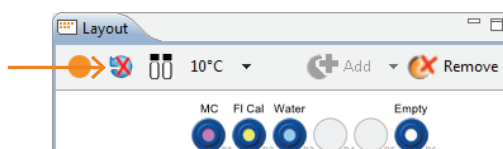


3. Compass for iCE can monitor the current during a separation for you, stop it if an abnormal current profile is detected and reinject the sample for you automatically. Reinjections are enabled by default, but if you don't want to use this feature just click the reinject icon to toggle it off.

NOTES:

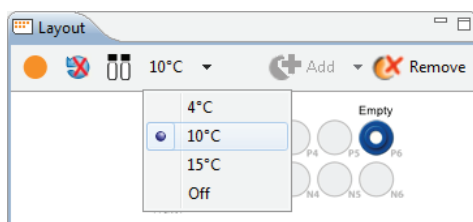
If a sample was reinjected, the injection row will be flagged in the Injections pane in the Run Summary screen. See “Injection Flags” on page 414 for more info.

A maximum of 10 reinjections are allowed per batch.



4. The sample platform is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The two reagent rows are kept at room temperature.



Step 3 - Assign Your Method Parameters

NOTE: We recommend using the default method parameters. Please see the [Maurice cIEF Method Development Guide](#) for more information on method optimization.

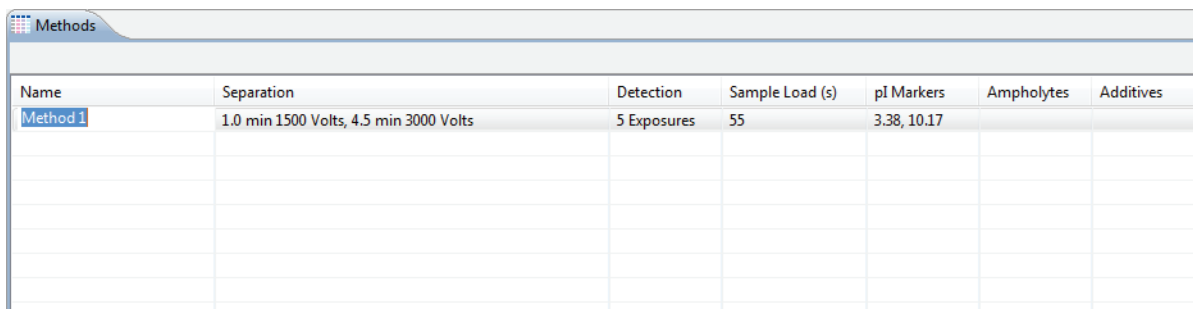
The Methods pane lists all of the methods used in a batch. You can add and remove methods here and change method parameters.

To import a saved method:

1. Select **File** in the main menu and click **Import Method**.
2. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

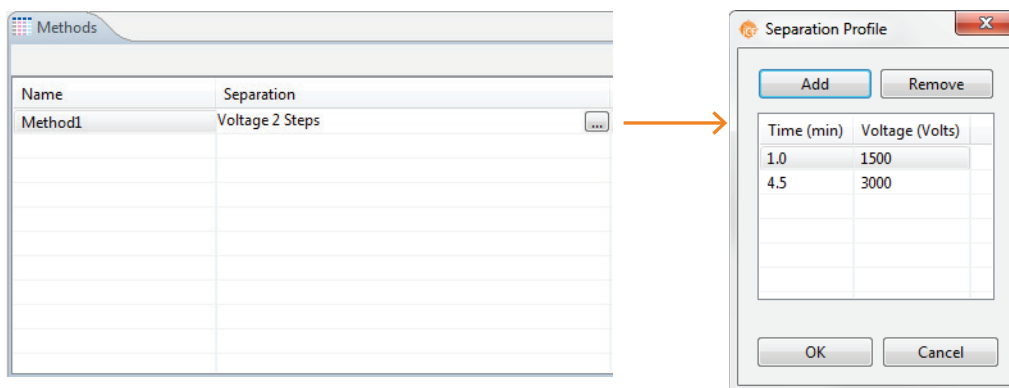
To create a new method or make changes to an existing one:

1. In the Methods table, click the first cell in the Name column and enter a new method name if needed.



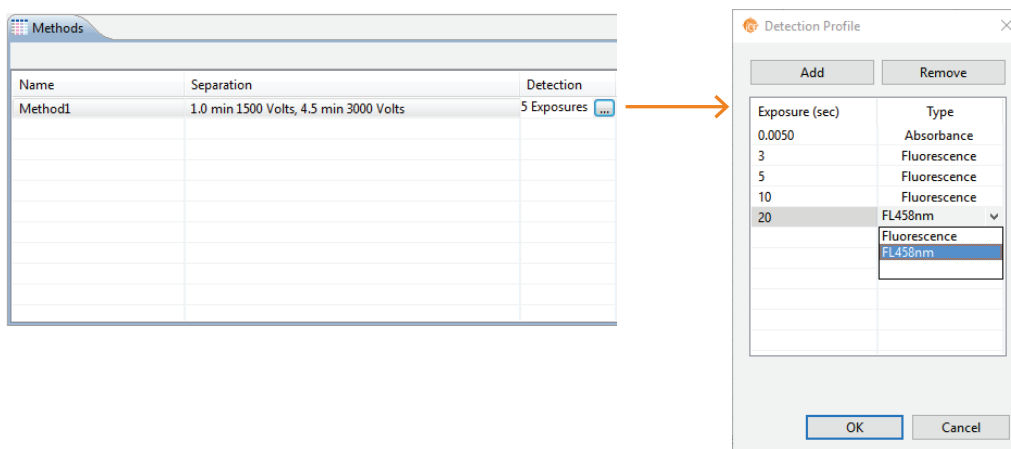
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17		

2. Click the first cell in the Separation column, then click the selection button [...] to set the pre-focusing and focusing time (in minutes) and voltage (V).



- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a load time (in minutes) and voltage value (in V).
- **To remove a profile step:** Select the row you want to remove and click **Remove**.

- Click the first cell in the Detection column the selection button [...] to set your exposure times for absorption and fluorescence detection modes.



- **To change the exposure time:** Just click in a cell under **Exposure** and type the new value(s) in seconds.

NOTES:

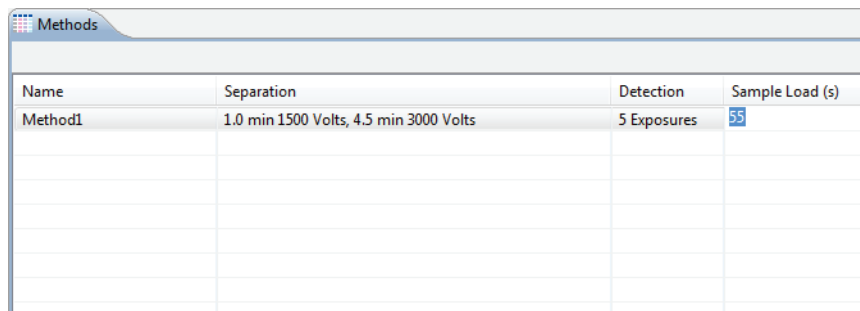
For standard Maurice, Maurice C. and MauriceFlex systems, the first exposure uses Absorbance detection and is a default setting that can't be changed or removed. Fluorescence detections (native fluorescence or FL458 nm) can be changed or removed.

If your Maurice system has the optional 458 nm filter installed, you can choose either the native fluorescence or the FL458 nm fluorescence filter for any of the remaining exposures. The FL458 nm filter enables detection of fluorescence emission at a longer wavelength to analyze molecules other than proteins, such as small molecule drugs in antibody-drug conjugates (ADCs). When used, ratiometric analysis of fluorescence and absorbance data can be applied to support applications including drug-antibody ratio (DAR) analysis.

- **To change the fluorescence detection for the exposure:** Click in a cell under **Type** and select an option.
- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in an exposure time (in seconds). **Optional:** If your Maurice has the FL458 nm filter, you can change the fluorescence detection for the exposure. Click in the **Type** cell to select an option.
- **To remove a profile step:** Select the row you want to remove and click **Remove**.

- Click the first cell in the Sample Load(s) column and set the load time in seconds.

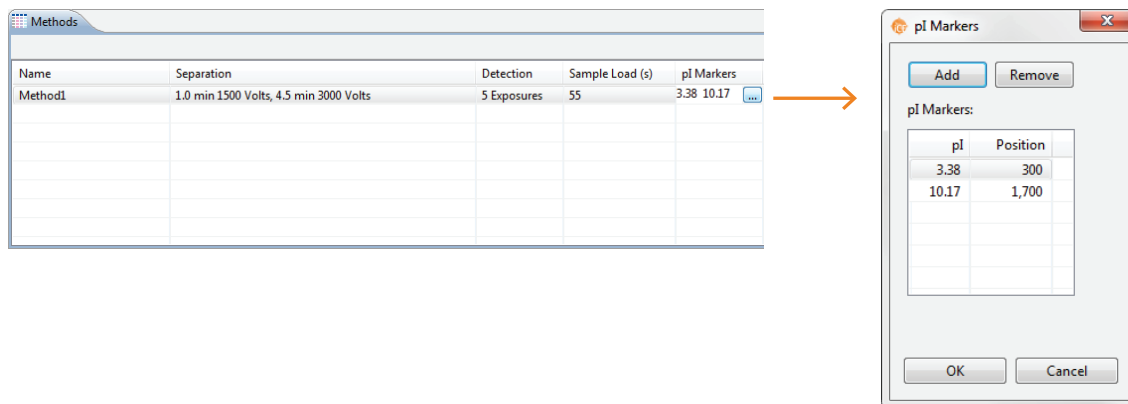
NOTE: We recommend using the default Sample Load time of 55 seconds. Please contact ProteinSimple Technical Support if you have questions on the Sample Load time to use for your application.



Name	Separation	Detection	Sample Load (s)
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55

- Click the first cell in the pI Markers column to select pI markers. Add new markers or remove existing ones then click OK.

NOTE: When you edit the pI markers in the method for a batch, Compass for iCE automatically creates a Markers group in the pI Markers Analysis settings for you.



- **To change a pI marker and position:** Just click in a cell under **pI** or **Position** and type the new value(s).
- **To add a pI marker:** Click **Add**. A new row will be added in the table. Then just type in a pI and a position (in pixels).
- **To remove a pI marker:** Select the row you want to remove and click **Remove**.

6. Optional: Click the first cell in the Ampholytes column and enter the ampholytes you're using.

Methods						
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17	Pharmalyte 3-10	

7. Optional: Click the first cell in the Additives column and enter any additives you're using.

Methods						
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17	Pharmalyte 3-10	Urea

8. You can now:

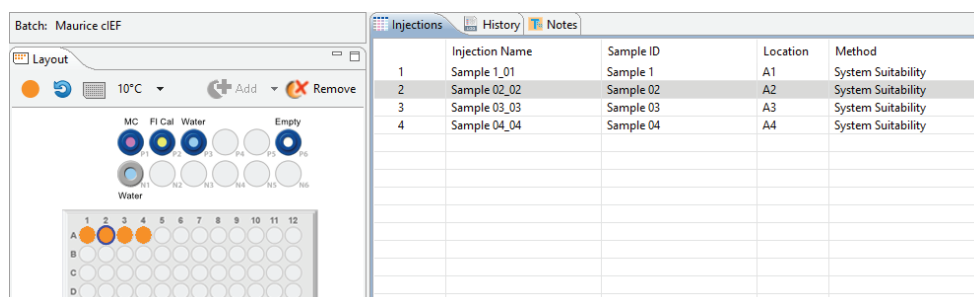
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.
- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.

Step 4 - Set Up Your Injections

The Injection pane lists all sample injections for the batch. Any samples that were added in “Step 2 - Assign Your Samples” are automatically added to this list.

NOTE: If you imported injections in Step 2, you can update the injections table and assign methods as needed using the information that follows. Otherwise skip to “Step 5 - Setup On-Board Mixing (Optional)” on page 70.

Selecting an injection in the table also selects the sample the injection will be made from in the plate or vial map in the Layout pane and vice-versa.



1. To add sample names, click the **Sample ID** cell for the injection and type a name.

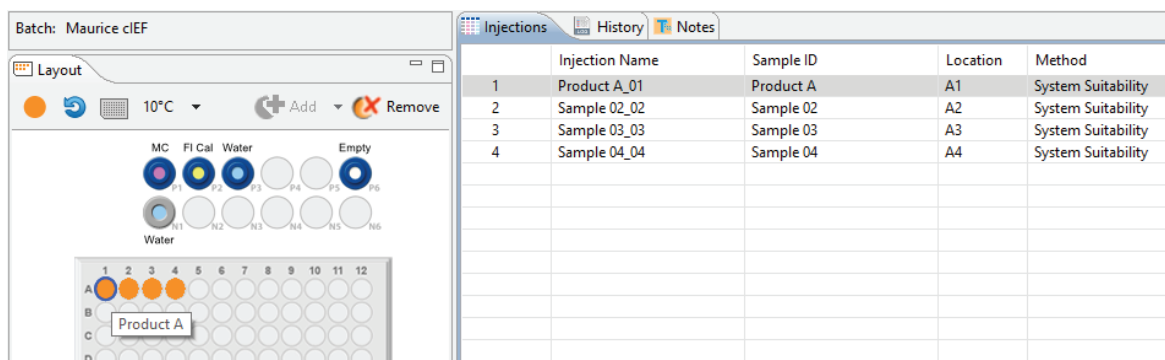
NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 82 for more info.

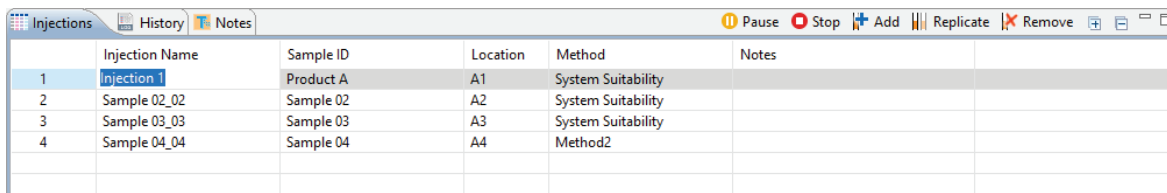
If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

Injections History Notes					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	System Suitability	

The sample name also displays when you hover the mouse over the sample in the plate or vial map:



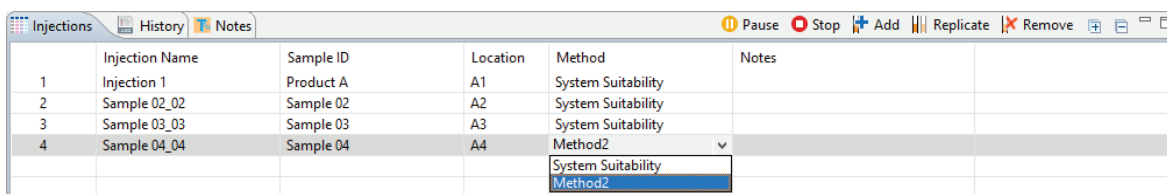
- Injection names are automatically set to 'Sample ID_injection number'. To change the name, click the **Injection Name** cell for the injection and type a name. Each injection name must be unique. If they aren't when you save the batch, Compass for ICE highlights any non-unique injection names so you can change them.



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	

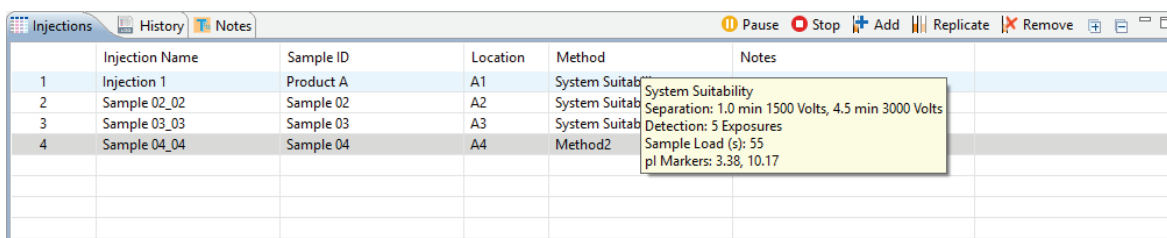
NOTE: Changing the injection name won't affect the sample ID.

- Assign your methods to each injection. Just click the **Method** cell for the injection and select a method from the drop down menu. If you added your samples with pre-assigned methods and don't need to make any changes, skip to the next step.



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	

Hovering over a method name displays the method parameters:

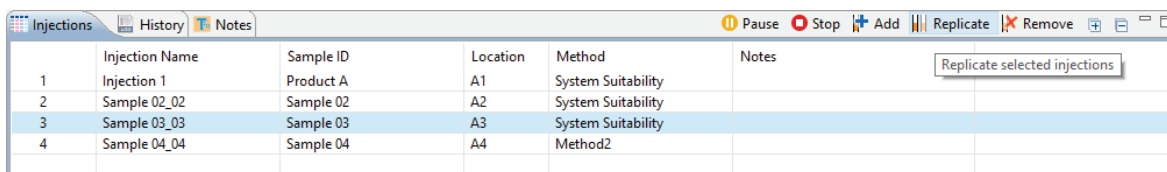


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	

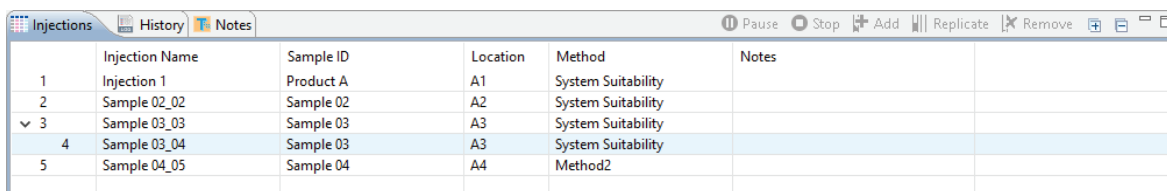
System Suitability
 Separation: 1.0 min 1500 Volts, 4.5 min 3000 Volts
 Detection: 5 Exposures
 Sample Load (s): 55
 pI Markers: 3.38, 10.17

4. You can add or remove injections as needed using a few different options. If you don't need to make any changes here, just skip to the next step.

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

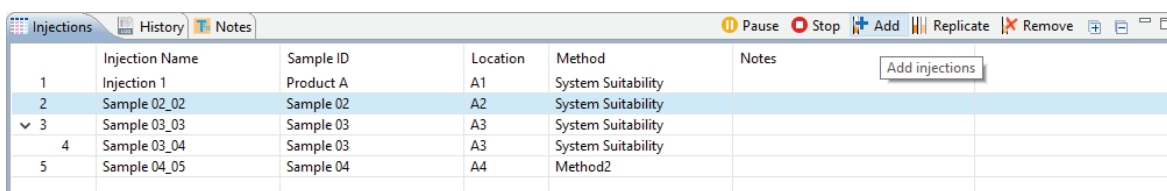


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 03_04	Sample 03	A3	System Suitability	
5	Sample 04_05	Sample 04	A4	Method2	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 03_04	Sample 03	A3	System Suitability	
5	Sample 04_05	Sample 04	A4	Method2	

- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

Step 5 - Setup On-Board Mixing (Optional)

If you're using Maurice-OBM or Maurice C.-OBM, you can tell Compass for iCE which samples to automatically mix for you. Up to 96 samples can be on-board mixed in a batch. You can also have a mix of premixed samples and samples to on-board mix in the same batch.

NOTES:

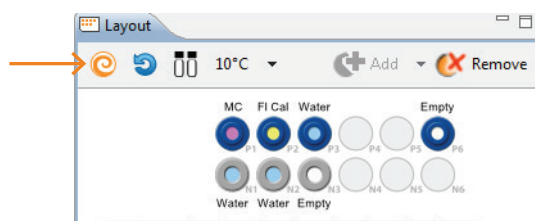
OBM is only available on Maurice and Maurice C. systems.

To use the OBM function, you must load 25 µL of sample in every sample vial or plate well that will be mixed on board. Maurice will automatically mix 25 µL of each sample with 100 µL of IEF Separation Mix prior to injection.

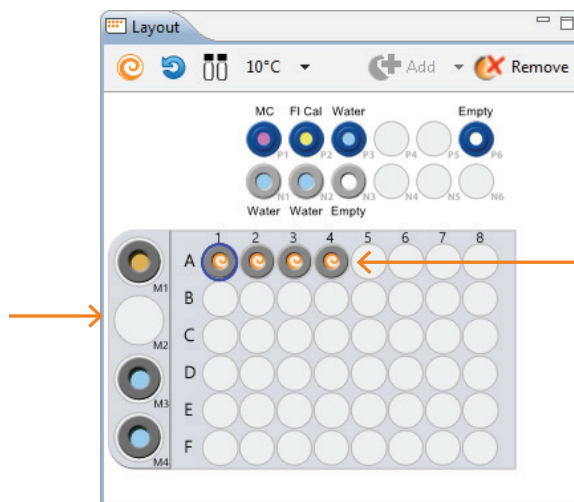
1. Make sure you've placed the following on-board mixing reagents in Maurice:

- **N2** - Water vial with **clear screw cap**
- **N3** - Empty vial with **clear screw cap**
- **M1** - 6 mL IEF Separation Mix vial
- **M2** - 6 mL IEF Separation Mix vial (use only when running more than 48 samples)
- **M3** - 6 mL water vial
- **M4** - 6 mL water vial

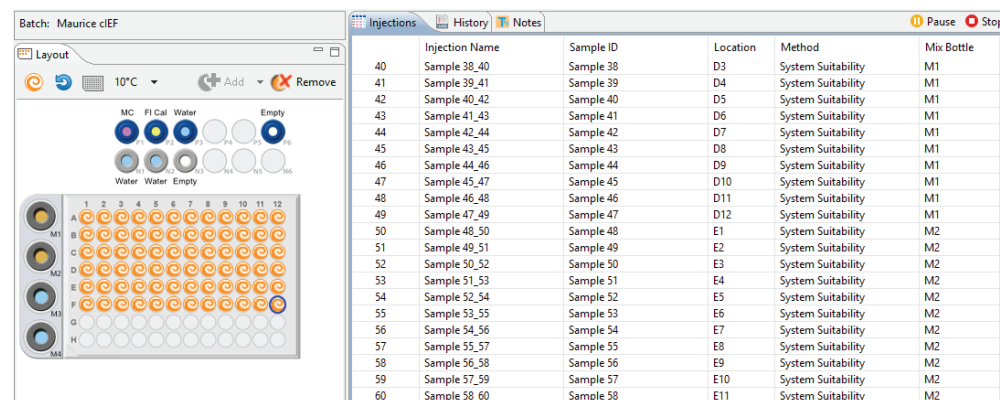
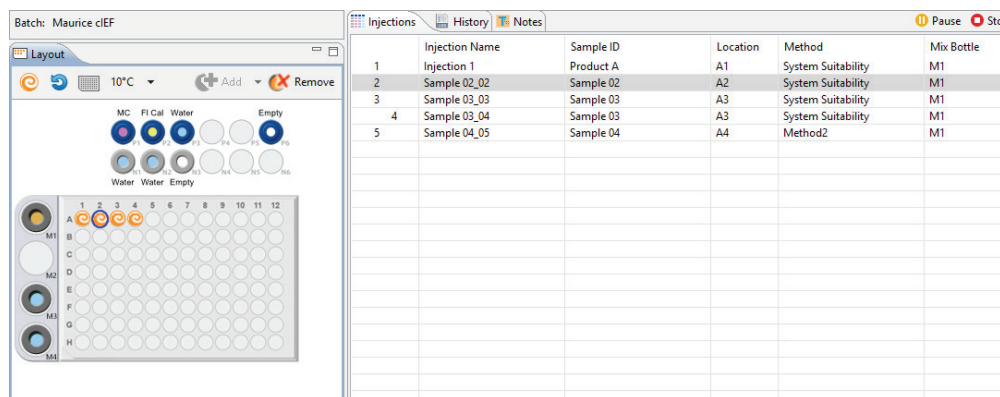
2. The on-board mixing function is disabled by default. Click the **on-board mixing icon** to toggle it on.



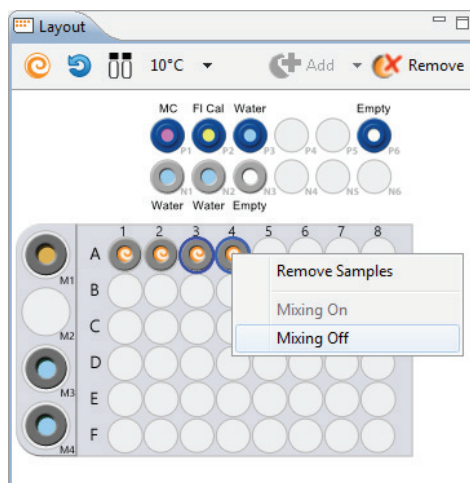
All sample positions will now display the mixing icon, and the on-board mixing reagent vial positions will also display on the left:



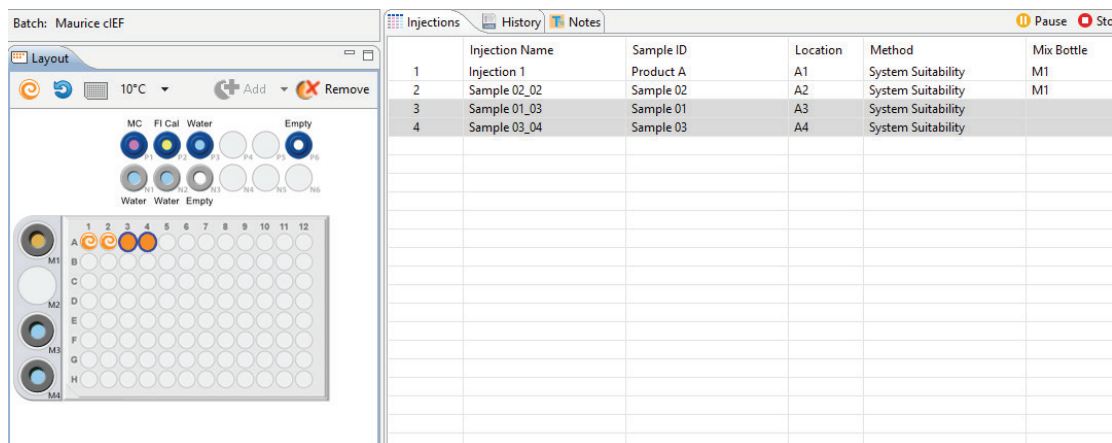
The Injection pane will also display a Mix Bottle column showing the position of the IEF Separation Mix vial that will be used for each sample. The vial in the M1 position is used for samples 1–48, and the vial in M2 will automatically display and be used for samples 49–96:



- To turn on-board mixing off for specific samples, select those well(s) or vial(s), right click and select **Mixing Off**.



The sample positions will display solid orange indicating mixing is off, and the Mix Bottle designation in the Injections pane is removed:



Step 6 - Add Programmed Pauses and Stops (Optional)

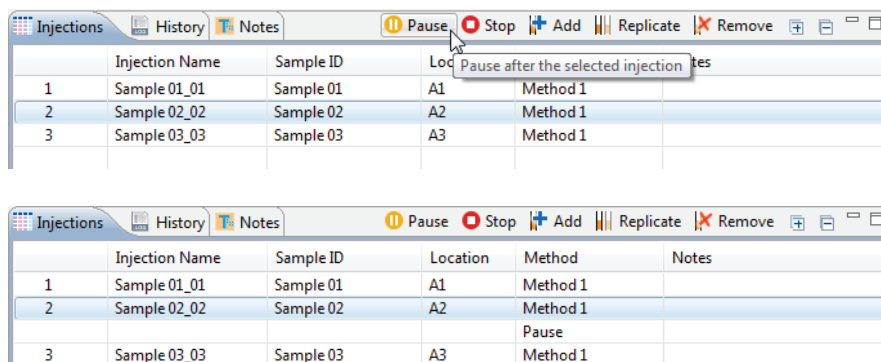
You can program pauses and stops in the batch to automatically pause or stop after injections even when you aren't in front of Maurice.

To program a pause:

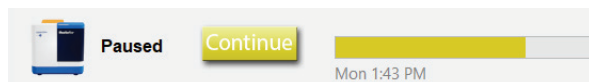
Pauses are helpful when you want to review injection results before continuing with the rest of the batch.

1. Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete.

NOTE: Maurice can tweet you when the batch pauses. See “Setting Up Maurice Systems to Send Tweets” on page 764.

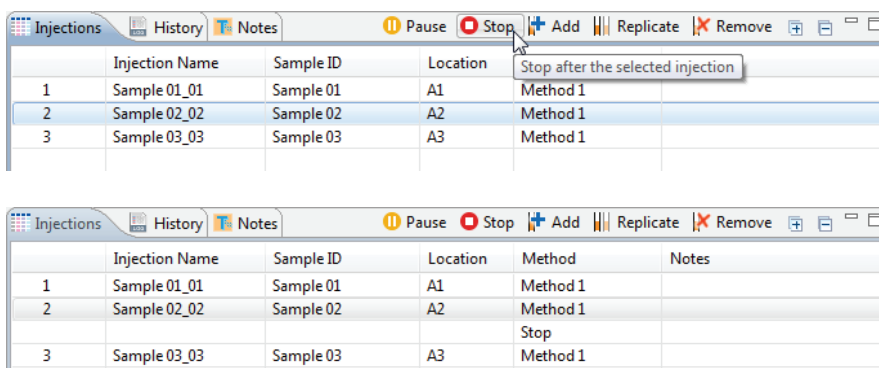


- To resume the batch, click **Continue** in the instrument status bar:



To stop the run after a specific injection:

- Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.

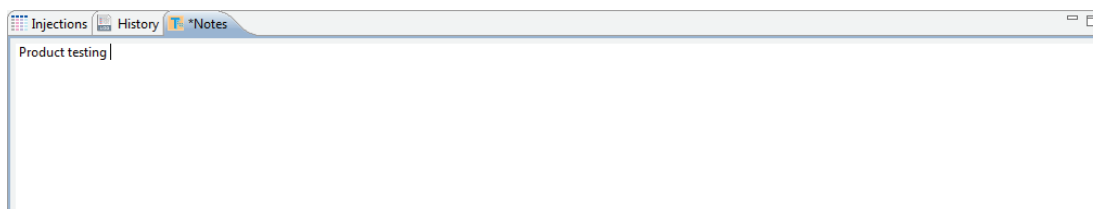


Data collected for injections completed up to the point where the batch was stopped are saved in the .mbz file.

- Be sure to complete your normal post-batch procedures.

Step 7 - Add Batch Notes (Optional)

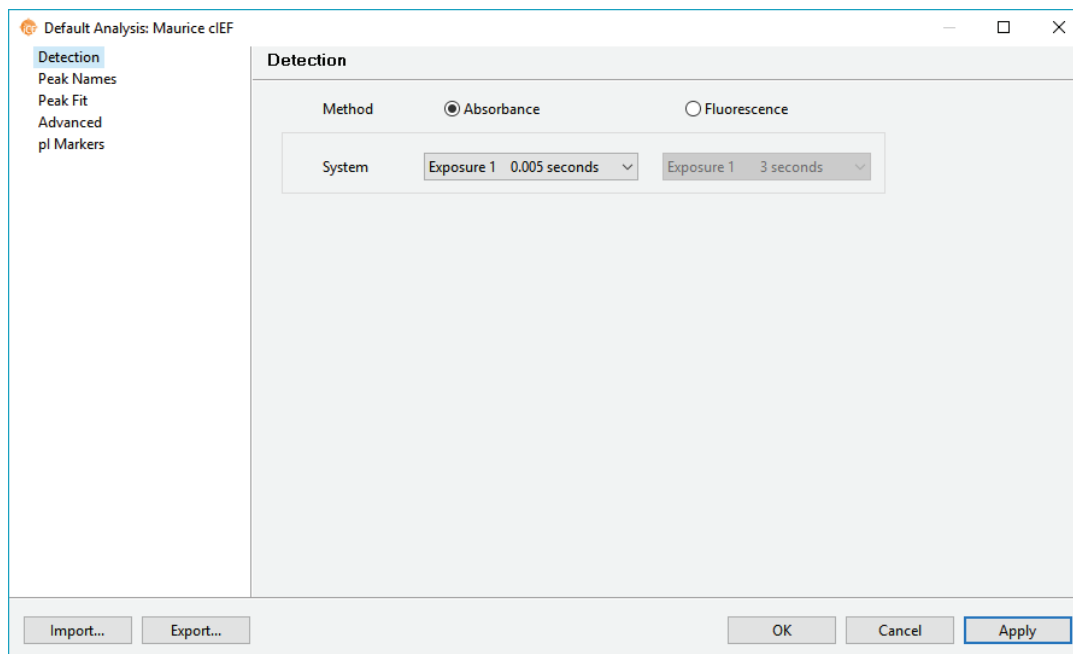
- Click on the **Notes** pane.
- Click in the notes area and type any information you want to add about your batch.



Step 8 - Modify Default Analysis Parameters (Optional)

You can preset the parameters used to analyze data generated with the batch. We recommend using the default parameters for cIEF applications, but if you need to modify parameters:

1. Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



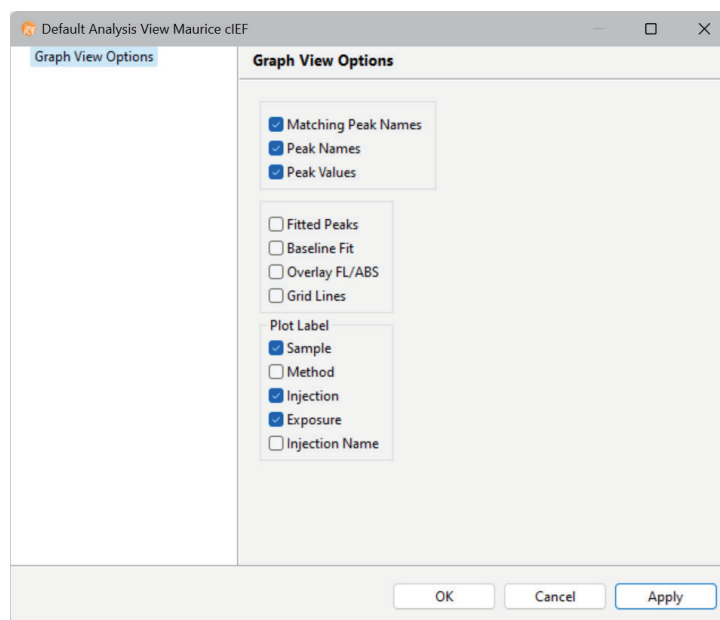
2. Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to "Analysis Settings Overview" on page 528.

Step 9 - Modify Default Analysis View Parameters (Optional)

You can preset the parameters used to display the electropherogram data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button.

To modify the parameters:

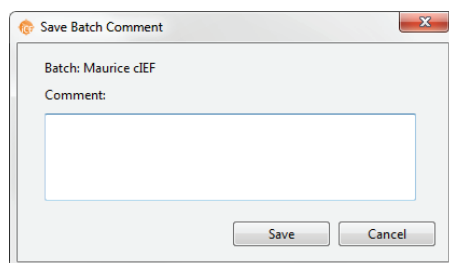
1. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



2. Change the parameters you want to, then click **Apply**. Click **OK** when you are done changing display parameters. For detailed information on graph view options, please refer to “Customizing the Data Display” on page 509.

Step 10 - Save Your Batch

1. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



2. Enter a name for your batch then click **Save**.

Viewing Replicate Injections

Injections with replicates have an arrow next to the injection number. You can expand or collapse the list of replicate injections by toggling the arrow:

The first screenshot shows the 'Injections' table with injection 1 expanded, displaying a list of 20 rows. The second screenshot shows the same table with injection 1 collapsed, displaying only the main injection row.

	Sample ID	Location	Method
▶ 1	Product A	A1	Method1
▶ 6	Product B	A2	Method1
11	Product C	A3	Method1
12	Product D	A4	Method1
13	Product E	A5	Method1
14	Product A	A6	Method2
15	Product B	B1	Method2
16	Product C	B2	Method2
17	Product D	B3	Method2
18	Product E	B4	Method2
19	Test	B5	Method2
20	Test	B6	Method2

	Sample ID	Location	Method
▶ 1	Product A	A1	Method1
2	Product A	A1	Method1
3	Product A	A1	Method1
4	Product A	A1	Method1
5	Product A	A1	Method1
▶ 6	Product B	A2	Method1
11	Product C	A3	Method1
12	Product D	A4	Method1
13	Product E	A5	Method1
14	Product A	A6	Method2
15	Product B	B1	Method2
16	Product C	B2	Method2

- To show all replicate injections in the batch, click the **Expand All Injections** button.

The screenshot shows the 'Injections' table with all replicate injections expanded. An orange arrow points to the 'Expand All Injections' button in the top right corner of the table.

	Sample ID	Location	Method	Notes
▶ 1	Product A	A1	Method1	
2	Product A	A1	Method1	
3	Product A	A1	Method1	
4	Product A	A1	Method1	
5	Product A	A1	Method1	
▶ 6	Product B	A2	Method1	
7	Product B	A2	Method1	
8	Product B	A2	Method1	
9	Product B	A2	Method1	
10	Product B	A2	Method1	
11	Product C	A3	Method1	
12	Product D	A4	Method1	
13	Product E	A5	Method1	
14	Product A	A6	Method2	
15	Product B	B1	Method2	
16	Product C	B2	Method2	
17	Product D	B3	Method2	
18	Product E	B4	Method2	
19	Test	B5	Method2	
20	Test	B6	Method2	

- To hide all replicate injections in the batch, click the **Collapse All Injections** button.

The screenshot shows the 'Injections' table with all replicate injections collapsed. An orange arrow points to the 'Collapse All Injections' button in the top right corner of the table.

	Sample ID	Location	Method	Notes
▶ 1	Product A	A1	Method1	
▶ 6	Product B	A2	Method1	
11	Product C	A3	Method1	
12	Product D	A4	Method1	
13	Product E	A5	Method1	
14	Product A	A6	Method2	
15	Product B	B1	Method2	
16	Product C	B2	Method2	
17	Product D	B3	Method2	
18	Product E	B4	Method2	
19	Test	B5	Method2	
20	Test	B6	Method2	

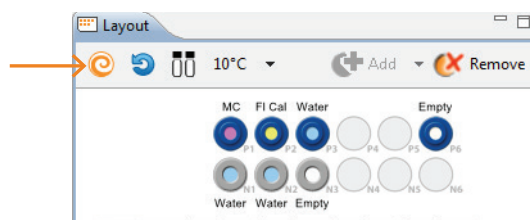
On-board Mixing (Maurice and Maurice C.)

Maurice's On-board Mixing function (OBM) can be used to automate preparation of cIEF samples by adding and aspirating the ampholyte mix into your protein samples. OBM is really useful when you need to analyze proteins using platform methods and when your proteins have a limited stability in their sample solution. It also minimizes assay complexity while maximizing mixing accuracy and consistency.

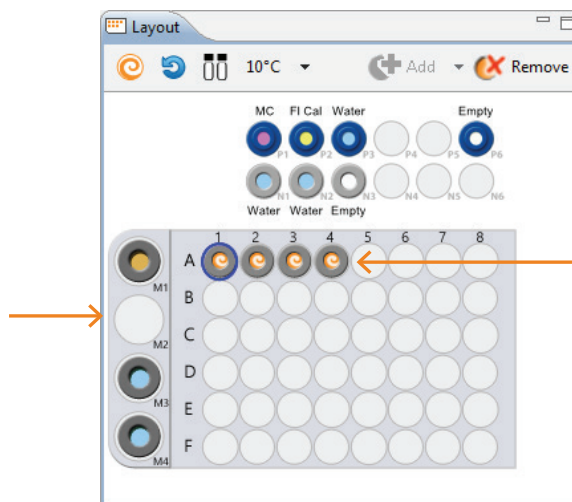
You can use OBM with samples in both 96-well plates and in vials, and on up to 96 samples in a single batch. The mixing protocol uses a fixed set of aspiration volume, speeds, and cycle parameters. The OBM protocol adds 100 μ L of IEF Separation Mix to 25 μ L of sample so the IEF Separation Mix components should be 1.25x that of the final desired concentration. Reagent positions M1–M4 on Maurice's sample and reagent platform all hold 6 mL vials that are used for on-board mixing.

- **M1 and M2:** Used for the IEF Separation Mix that contain the ampholytes, methyl cellulose, and excipient components of a cIEF sample solution. M1 is used for samples 1–48 and M2 is used for samples 49–96. The minimum IEF Separation Mix vial volume is 1.5 mL, and the maximum is 6.0 mL.
- **M3 and M4:** Used for DI water which is used in both aspirating the IEF Separation Mix and cleaning the OBM pipette between sample preparation cycles. Place two full, fresh 6 mL vials of DI water in positions M3 and M4 prior to running an OBM batch.

The on-board mixing function is disabled by default. Click the **on-board mixing icon** to toggle it on.



All sample positions will now display the mixing icon, and the on-board mixing reagent vial positions will also display on the left:



The Injection pane will also display a Mix Bottle column showing the position of the IEF Separation Mix vial that will be used for each sample. The vial in the M1 position is used for samples 1–48, and the vial in M2 will automatically display and be used for samples 49–96:

Batch: Maurice cIEF

Layout | 10°C | Add | Remove

MC FI Cal Water Empty
P1 P2 P3 P4 P5 P6
Water Water Empty
N1 N2 N3 N4 N5 N6

Injection Name	Sample ID	Location	Method	Mix Bottle
1 Injection 1	Product A	A1	System Suitability	M1
2 Sample 02_02	Sample 02	A2	System Suitability	M1
3 Sample 03_03	Sample 03	A3	System Suitability	M1
4 Sample 03_04	Sample 03	A3	System Suitability	M1
5 Sample 04_05	Sample 04	A4	Method2	M1

Batch: Maurice cIEF

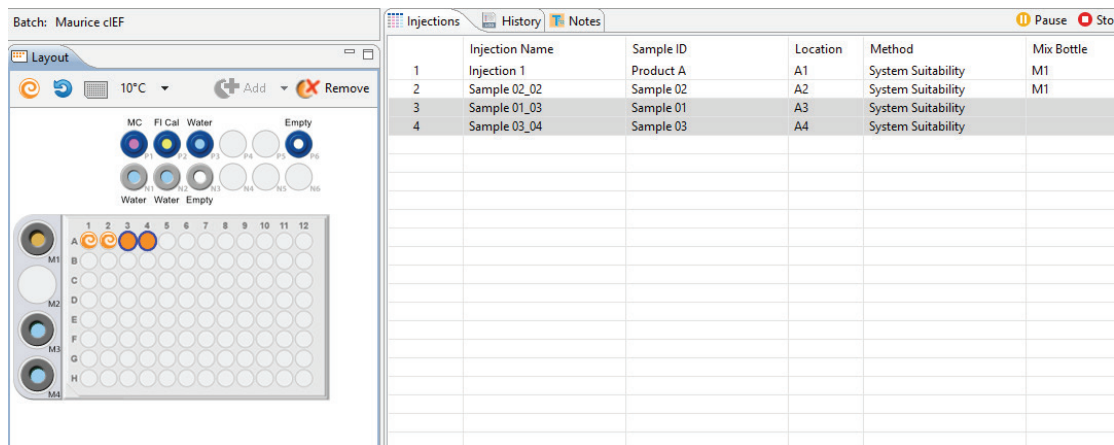
Layout | 10°C | Add | Remove

MC FI Cal Water Empty
P1 P2 P3 P4 P5 P6
Water Water Empty
N1 N2 N3 N4 N5 N6

Injection Name	Sample ID	Location	Method	Mix Bottle
40 Sample 38_40	Sample 38	D3	System Suitability	M1
41 Sample 39_41	Sample 39	D4	System Suitability	M1
42 Sample 40_42	Sample 40	D5	System Suitability	M1
43 Sample 41_43	Sample 41	D6	System Suitability	M1
44 Sample 42_44	Sample 42	D7	System Suitability	M1
45 Sample 43_45	Sample 43	D8	System Suitability	M1
46 Sample 44_46	Sample 44	D9	System Suitability	M1
47 Sample 45_47	Sample 45	D10	System Suitability	M1
48 Sample 46_48	Sample 46	D11	System Suitability	M1
49 Sample 47_49	Sample 47	D12	System Suitability	M1
50 Sample 48_50	Sample 48	E1	System Suitability	M2
51 Sample 49_51	Sample 49	E2	System Suitability	M2
52 Sample 50_52	Sample 50	E3	System Suitability	M2
53 Sample 51_53	Sample 51	E4	System Suitability	M2
54 Sample 52_54	Sample 52	E5	System Suitability	M2
55 Sample 53_55	Sample 53	E6	System Suitability	M2
56 Sample 54_56	Sample 54	E7	System Suitability	M2
57 Sample 55_57	Sample 55	E8	System Suitability	M2
58 Sample 56_58	Sample 56	E9	System Suitability	M2
59 Sample 57_59	Sample 57	E10	System Suitability	M2
60 Sample 58_60	Sample 58	E11	System Suitability	M2

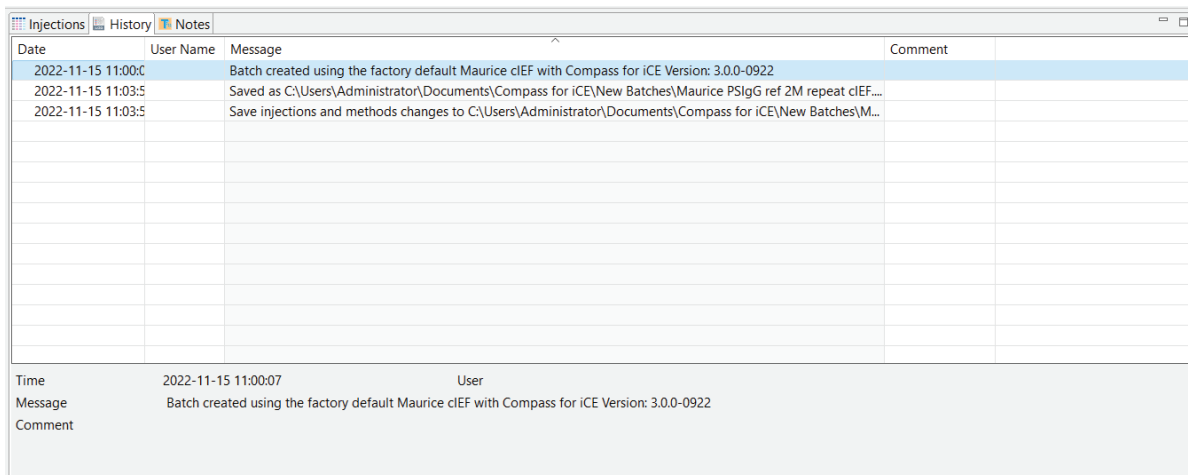
To turn on-board mixing off for specific samples, select those well(s) or vial(s), right click and select **Mixing Off**.

The sample positions will display solid orange indicating mixing is off, and the Mix Bottle designation in the Injections pane is removed:



Batch History

Clicking on the History pane shows the batch event history, starting with the date and time the batch was created through the most current event. Clicking on a row in the table displays the full event details in the box under the table.



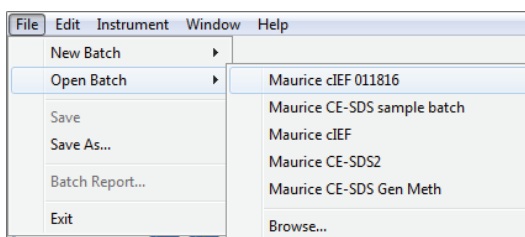
- **Date:** Date and time of the batch event.
- **User Name:** User that initiated the event. User names only display if you're using the Access Control feature to help satisfy the 21CFR Part 11 data security requirements. Go to "Enabling Access Control" on page 781 to learn how to set it up.
- **Message:** Description of the event that took place.
- **Comment:** Comments entered by the user when the batch was saved.

You can copy the information in the History pane to use in other documents:

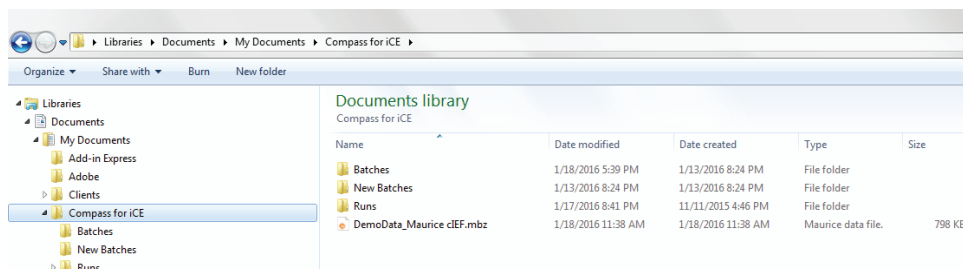
1. Click the **History** pane to make sure it's active.
2. Click **Edit** in the main menu and select **Copy**.
3. Open a document and click **Paste**.

Making Changes to a Batch

1. Select **File** in the main menu and click **Open Batch**.



2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file.
 - Select one of these files or click **Browse** to open the Batch folder and select a different one.
 - To open a batch used in a run file, click **Browse**. Go to the Run folder in the Compass for iCE directory and select the run file that uses the batch you want to edit.

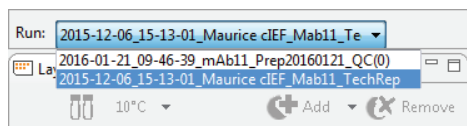


3. To make changes to the batch, see the steps in “Creating a New Batch” on page 59 “Creating a New Batch”. Then select **File** from the main menu and click **Save** or **Save As**.

Viewing and Editing Batches in Completed Runs

1. Click the **Analysis** screen and open your run file(s).
2. After the run opens, click the **Batch** screen to view the batch used with the run.

If you opened more than one run file, you can switch between viewing batches for each file. Just click the **down arrow** in the Run box and select the batch you want to view from the drop down list.



3. In completed run files, you can only make edits to the **sample** and **method** names in a batch, and these changes are saved to the run file only. Update sample and method names as needed.
4. Click the **Analysis** screen. Then select **File** from the main menu and click **Save** or **Save As** to save the new changes to the run file.

Copying and Pasting Injection Names or Sample IDs from other Documents

Injection names and Sample IDs can be copied from other documents and pasted into the Injection pane to make batch setup faster.

To paste a list of injection names into the batch:

1. Select the list of injection names in a document (Microsoft® Word®, Excel®, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 2	My Sample 2
4		My Injection 3	My Sample 3
5		My Injection 4	My Sample 4
6		My Injection 5	My Sample 5
7		My Injection 6	My Sample 6
8		My Injection 7	My Sample 7
9		My Injection 8	My Sample 8
10		My Injection 9	My Sample 9
11		My Injection 10	My Sample 10

2. Select an injection in the Injection pane, then hold the Ctrl+V keys to paste. You can also select the injection, right click and select **Paste**.

⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove ⚙ 📄					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	Method2	
3	Sample 03_03	Sample 03	A3	Method2	
4	Sample 04_04	Sample 04	A4	Method2	
5	Sample 05_05	Sample 05	A5	Method2	
6	Sample 06_06	Sample 06	A6	Method2	
7	Sample 07_07	Sample 07	A7	Method2	
8	Sample 08_08	Sample 08	A8	Method2	
9	Sample 09_09	Sample 09	A9	Method2	
10	Sample 10_10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

The injection names are pasted into the Injection pane:

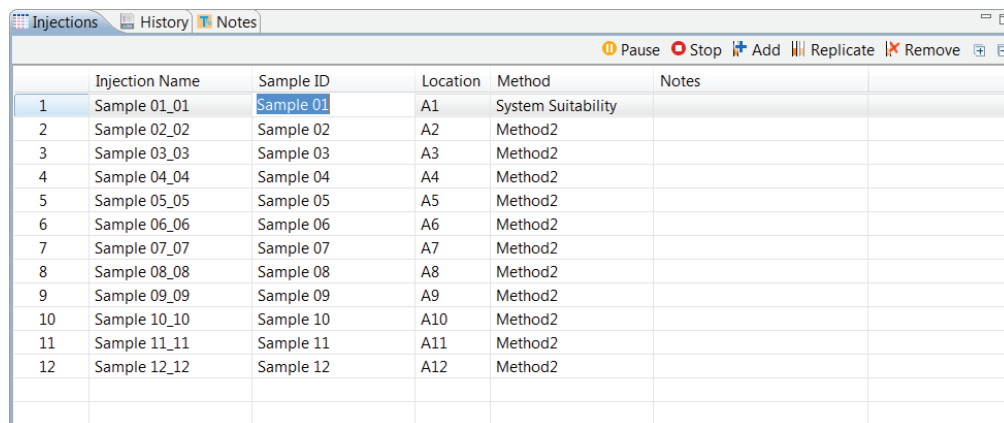
⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove ⚙ 📄					
	Injection Name	Sample ID	Location	Method	Notes
1	My Injection 1	Sample 01	A1	System Suitability	
2	My Injection 2	Sample 02	A2	Method2	
3	My Injection 3	Sample 03	A3	Method2	
4	My Injection 4	Sample 04	A4	Method2	
5	My Injection 5	Sample 05	A5	Method2	
6	My Injection 6	Sample 06	A6	Method2	
7	My Injection 7	Sample 07	A7	Method2	
8	My Injection 8	Sample 08	A8	Method2	
9	My Injection 9	Sample 09	A9	Method2	
10	My Injection 10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

To paste a list of Sample IDs into the batch:

1. Select the list of Sample IDs in a document (Microsoft Word, Excel, etc.) and select **Copy** or hold the Ctrl+C keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 2	My Sample 2
4		My Injection 3	My Sample 3
5		My Injection 4	My Sample 4
6		My Injection 5	My Sample 5
7		My Injection 6	My Sample 6
8		My Injection 7	My Sample 7
9		My Injection 8	My Sample 8
10		My Injection 9	My Sample 9
11		My Injection 10	My Sample 10
12			

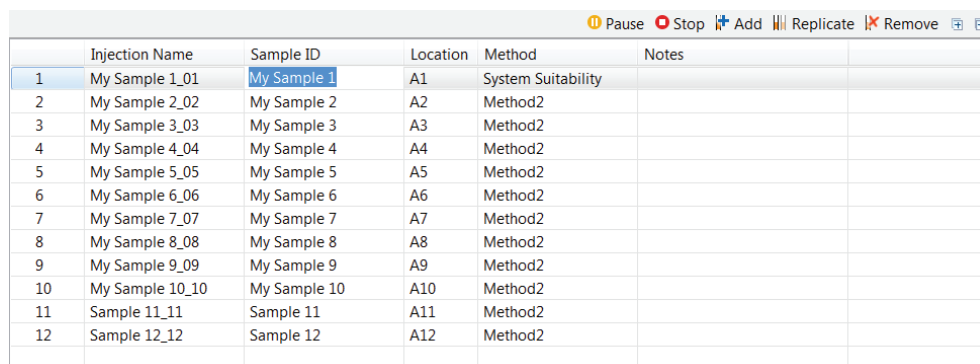
2. Select a Sample ID in the Injection pane, then hold the Ctrl+V keys to paste. You can also select the Sample ID, right click and select **Paste**.



	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	Method2	
3	Sample 03_03	Sample 03	A3	Method2	
4	Sample 04_04	Sample 04	A4	Method2	
5	Sample 05_05	Sample 05	A5	Method2	
6	Sample 06_06	Sample 06	A6	Method2	
7	Sample 07_07	Sample 07	A7	Method2	
8	Sample 08_08	Sample 08	A8	Method2	
9	Sample 09_09	Sample 09	A9	Method2	
10	Sample 10_10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

The Sample IDs are pasted into the Injection pane.

NOTE: If you paste in new sample IDs when the default injection names are still being used, the injection names will automatically update to match the sample ID. If you've already changed the injection names manually, updating the sample IDs will not update the injection names.



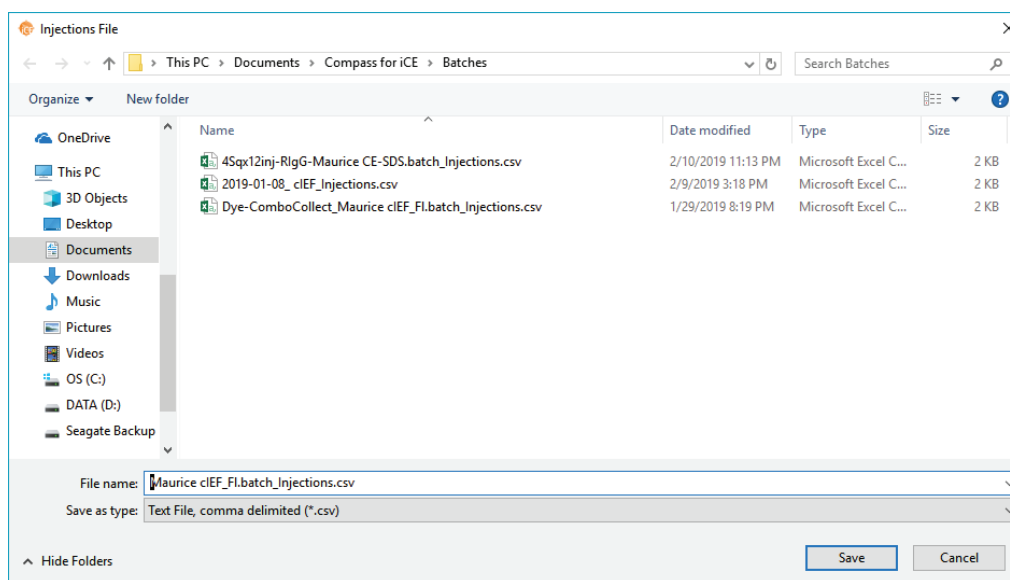
	Injection Name	Sample ID	Location	Method	Notes
1	My Sample 1_01	My Sample 1	A1	System Suitability	
2	My Sample 2_02	My Sample 2	A2	Method2	
3	My Sample 3_03	My Sample 3	A3	Method2	
4	My Sample 4_04	My Sample 4	A4	Method2	
5	My Sample 5_05	My Sample 5	A5	Method2	
6	My Sample 6_06	My Sample 6	A6	Method2	
7	My Sample 7_07	My Sample 7	A7	Method2	
8	My Sample 8_08	My Sample 8	A8	Method2	
9	My Sample 9_09	My Sample 9	A9	Method2	
10	My Sample 10_10	My Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

Importing and Exporting Injections

Injections in an open batch or run file can be exported as a separate file. This lets you import the same or a modified set of injections into another batch later, rather than having to enter the information manually.

Exporting Injections

1. Open the batch or run you want to export injections from.
2. In the Batch screen, select **File** in the main menu and click **Export Injections**. The following window displays:



3. The default directory is Compass for iCE/Batches. Change the directory if needed.
4. The default file name will be batchname.batch_Injections.csv. Change the name if needed and click **Save**.

Creating an Injections File

To create a new injections file .csv, it's easiest to export injections from an existing batch and use that as your template.

1. Follow the instructions in "Exporting Injections" above to export injections from an existing cIEF batch.
2. Open the .csv file in a program that provides a table/spreadsheet format.
3. To add or change injection information:
 - a. Type in a unique name in each cell in the Injection Name column. For replicate injections, type > in front of the injection name, for example: >Injection 1.
 - b. Type in a sample ID, location and method for each injection in their respective columns.
 - c. Optional: Enter a mix bottle (M1 or M2) if on-board mixing will be used for any of the injections. If you leave the mix bottle cell blank, mixing will not be used for that injection.
 - d. Optional: Type in notes if needed.

NOTE: Programmed pauses and stops can't be added in the injections file. They can be added in the Injections pane after importing your injections into Compass for iCE.

	A	B	C	D	E	F
1	Injection Name	Sample ID	Location	Method	Mix Bottle	Notes
2	Inj01	R01_SSPP	E4	System Suitability		
3	Inj02	S01_Her2-A1_None_0M Urea	D1	Method_M458		
4	>Inj02	S01_Her2-A1_None_0M Urea	D1	Method_M458		
5	>Inj02	S01_Her2-A1_None_0M Urea	D1	Method_M458		
6	Inj03	S02_Her2-B5_None_0M Urea	D2	Method_M458		
7	Inj04	S03_Her2-B7_None_0M Urea	D3	Method_M458		
8	Inj05	S04_Her2-A1_CPM_0M Urea	D4	Method_M458		
9	Inj06	S05_Her2-B5_CPM_0M Urea	D5	Method_M458		
10	Inj07	S06_Her2-B7_CPM_0M Urea	D6	Method_M458		
11	Inj08	S07_Her2-A1_CPM_4M Urea	D7	Method_M458		
12	Inj09	S08_Her2-B5_CPM_4M Urea	D8	Method_M458		
13	Inj10	S09_Her2-B7_CPM_4M Urea	D9	Method_M458		

4. Save the .csv file.

Importing Injections

1. Open the batch you want to import injections into, or open a new batch.
2. Select **File** in the main menu and click **Import Injections**.
3. Select an injections file (*.csv) and click **OK**. The imported information will display in the Layout and Injections panes.

NOTE: Imported injections will replace any injections already in the Injections table, they will not append an existing table.

Importing and Exporting Methods

Methods in an open batch or run file can be exported as a separate file. This lets you import the same method into another batch later, rather than having to re-enter method information manually.

Importing a Method

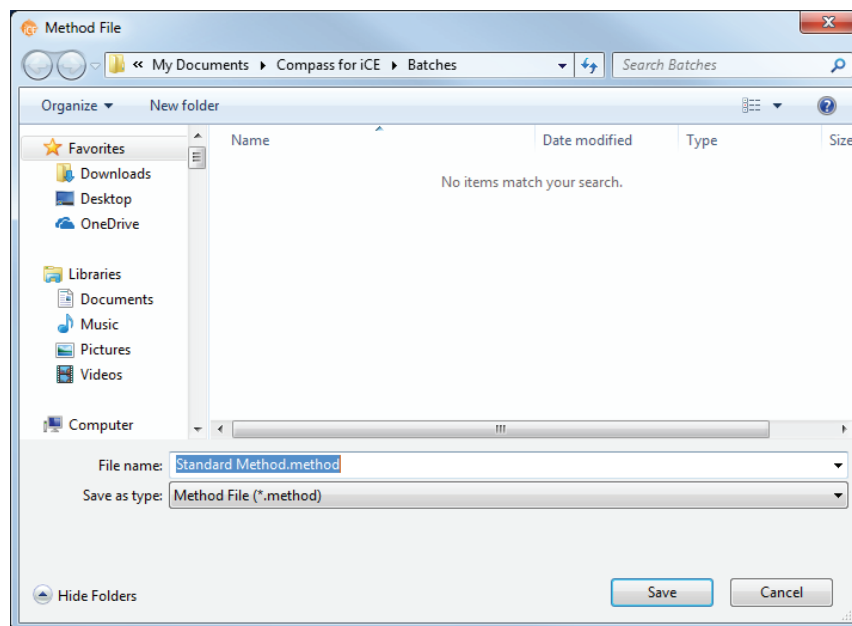
NOTE: Importing a method imports information into the Batch window's Method pane only.

1. Open the batch you want to import the method into.
2. Select **File** in the main menu and click **Import Method**.
3. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

Exporting a Method

NOTE: Exporting a method exports information in the Batch window's Method pane only.

1. Open the batch you want to export the method from.
2. In the Method pane, select the method you want to export. To export more than one method at the same time, hold the Ctrl key to select multiple methods.
3. Select **File** in the main menu and click **Export Method**. The following window displays:



4. The default directory is Compass for iCE/Batches. Change the directory if needed.
5. Enter a method file name and click **Save**. The settings will be saved as a *.method file.

Batch Reports

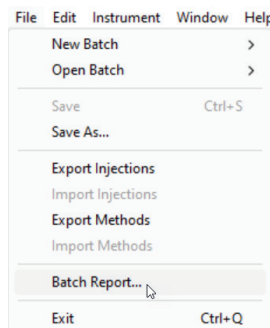
You can export a PDF file of sample and method details for each injection in the batch.

NOTES:

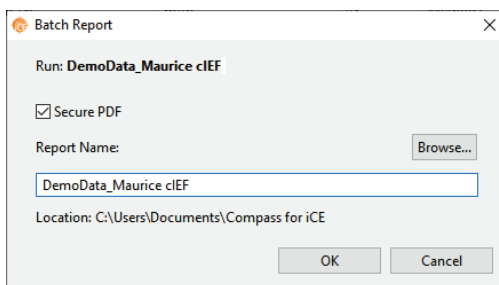
You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it's saved in a separate run folder. Existing reports aren't overwritten.

1. Go to the **Analysis** or **Run Summary** screen, then click **File > Open Run** and select a run file (if you don't have one open already).
2. After the run opens, go to the **Batch** screen.
3. Select **File** from the main menu and click **Batch Report**.

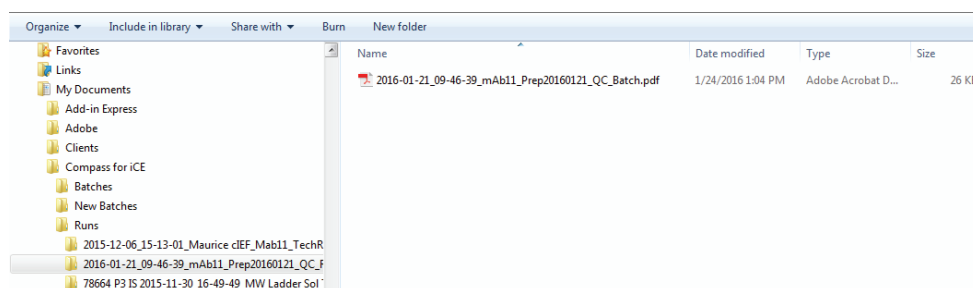


4. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.



Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don't want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.

5. The Batch Report PDF is exported to the **Runs** folder in the Compass for iCE directory. It'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Here's an example Batch Report:

cIEF Batch: Maurice cIEF

Injections

Injection	Injection Name	Sample ID	Location	Method	Notes
1	System Suit_01	System Suit	A1	Method1	
2	Peptide Mix_02	Peptide Mix	A2	Method1	
3	mAb_03	mAb	A3	Method2	
4	System Suit_04	System Suit	A1	Method1	
5	Peptide Mix_05	Peptide Mix	A2	Method1	
6	mAb_06	mAb	A3	Method2	

Methods

Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min, 1500 Volts 4.5 min, 3000 Volts	Absorbance, 0.005 sec Fluorescence, 1 sec Fluorescence, 3 sec Fluorescence, 5 sec Fluorescence, 10 sec	90	3.38 pI, 250 pixels 10.17 pI, 1700 pixels		
Method2	1.0 min, 1500 Volts 6.0 min, 3000 Volts	Absorbance, 0.005 sec Fluorescence, 1 sec Fluorescence, 5 sec Fluorescence, 10 sec Fluorescence, 15 sec Fluorescence, 20 sec	90	4.05 pI, 500 pixels 9.99 pI, 1700 pixels		

Batch Log

Date	User Name	Message	Comment
10/27/2015 12:01 PM		Batch created using the factory default Maurice cIEF	
10/27/2015 12:11 PM		Save protocol and template changes	Auto-saved

Created By: Andrea Sat 6:27 PM Mar 4, 2023 PST
 C:\Users\Andrea\Documents\Compass for iCE\Runs\DemoData_Maurice cIEF.mibz
 Computer: DESKTOP-1FMT605 Software Version: Compass for ICE 4.0.0, Build ID: 0222

Page 1 of 2



Chapter 6:

Running cIEF Applications

Chapter Overview

- Overview
- Before You Throw the Switch
- Power Up
- Running cIEF Applications
- Editing a Running Batch
- Post-batch Procedures
- Checking Your Data

Overview

Standard cIEF applications can be run on Maurice, Maurice C. or MauriceFlex systems using a cIEF cartridge.

Before You Throw the Switch

Ensure that everyone using Maurice have:

- Received instruction in general safety practices for laboratories.
- Received instruction in specific safety practices for Maurice.
- Received instruction on handling of biohazards (if biohazardous materials are to be used on Maurice).
- Read and understood all Product Inserts and related Safety Data Sheets (SDS).

Power Up

1. Turn on the computer connected to Maurice.
2. Turn on Maurice's main power switch.
3. Wait for Maurice to initialize. His indicator light will turn from magenta to blue when he's done.
4. Double-click the Compass for iCE software icon on the computer desktop to open the application.
5. Connect Maurice to Compass for iCE.

Running cIEF Applications

What You'll Need

- Maurice cIEF Cartridges
- Maurice cIEF Method Development Kit (optional)
- Maurice System Suitability Kit (optional)
- Maurice cIEF Fluorescence Calibration Standard
- Maurice cIEF pI Markers (3.38, 4.05, 5.85, 6.14, 7.05, 8.4, 9.50, 9.99, or 10.17)
- 0.5% Methyl Cellulose Solution
- 1% Methyl Cellulose Solution
- iCE Electrolyte Kit
- Deionized (DI) water

- Glass reagent vials, 2 mL



MauriceFlex crimp top glass vial (PN 110-0019, left) and glass reagent vial (PN 046-017, right)
Only glass reagent vials (PN 046-017) should be used to prepare cIEF batch reagents.

- 96-well plate or vials with integrated inserts for samples
- Clear screw caps for vials
- Blue pressure caps for vials
- P10, P20, P200, P1000 pipettes and tips
- Electrolyte pipette
- Vortexer
- Table-top centrifuge with plate adapter or vial adapter (12 mm, 2 mL vials)

IMPORTANT:

Use glass reagent vials, 2 mL (PN 046-017) to prepare cIEF batch reagents.

Step 1: Prep Your Markers, Samples and Reagents

NOTES:

You can prepare your samples to run either in 96-well plates or vials.

If you need to seal the 96-well plate during your run, we recommend the Slit Seal, 96 well plate seal from BioChromato (<https://biochromato.com/slit-seal/>). It can be used in both absorbance and native fluorescence modes. If you're currently using X-Pierce adhesive film (PN XP-100, Excel Scientific), we recommend using it in absorbance mode only.

System Suitability Peptide Panel (Optional)

NOTES:

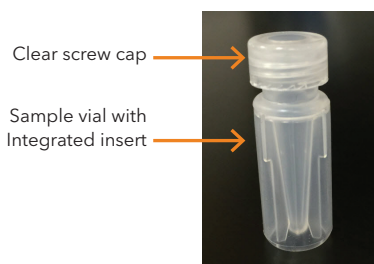
Run the System Suitability Peptide Panel when you need to confirm performance on Maurice.

The System Suitability Peptide Panel is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

1. Using scissors, carefully cut the top the package off leaving the sealing strip intact.
2. Take out the strip of tubes and cut one clear tube of lyophilized System Suitability Peptide Panel from the strip. Return the remaining tubes to the original package, reseal tightly and store at 2–8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Add 40 µL of DI water to the tube. Gently resuspend by pipetting up and down to mix.
5. Add 160 µL of the System Suitability Test Mix to the freshly reconstituted Peptide Panel. Gently mix by pipetting up and down. Transfer this solution to a 1.5 mL microcentrifuge tube.
6. Vortex the tube 3 times, 5 seconds each.
7. Centrifuge the tube at 10,000 xg for 3 minutes to sediment any particulates.
8. Carefully aspirate the top 160 µL of the solution and pipette it into a sample vial with integrated insert or well of a 96-well plate. You'll want to insert the pipette tip all the way to the bottom of the insert or well when you dispense the solution to avoid introducing bubbles.

NOTE: Make sure to check for and remove any bubbles at the bottom of the sample vial or well.

9. If you're using vials, close the sample vial with a clear screw cap.



Samples

1. In a microcentrifuge tube, prepare your sample at a concentration of 1 mg/mL in a final volume of 40 μ L in DI water.
2. In a separate tube, prepare IEF Separation Mix containing your chosen pI marker(s).

NOTE: Check out the Method Development Guide for suggested IEF Separation Mix recipes.

3. Add 160 μ L of IEF Separation Mix to the 40 μ L of your sample.
4. Vortex the tube 3 times, 5 seconds each.
5. Centrifuge the tube at 10,000 $\times g$ for 3 minutes to sediment any particulates
6. Carefully aspirate the top 160 μ L of the sample and pipette it into your sample vial with integrated insert or well of a 96-well plate by inserting the pipette tip all the way to the bottom to avoid introducing bubbles.

Note: Make sure to check for and remove any bubbles at the bottom of the sample vial or well.

7. If you're using vials, close the sample vial with a clear screw cap.
8. Spin the plate or sample vials at 1,000 $\times g$ for 5 minutes using the appropriate centrifuge adapter.

pI Markers

1. Open the vial of lyophilized pI marker by lifting the center tab and gently pulling it back to break the metal seal. Slowly remove the rubber stopper.
2. Add 210 μ L of DI water to the vial.
3. Put the rubber stopper back in the vial and vortex 3 times, 5 seconds each time to completely reconstitute the lyophilized cake. Repeat this step for all pI markers you're using.
4. Aliquot 20 μ L of each reconstituted pI marker into separate tubes for storage.

NOTES:

Keep your reconstituted pI markers on ice until you're ready to add them to your sample or IEF Separation Mix.

If you'll use the pI markers within a month, store the aliquots at 2-8 °C. Otherwise, store the aliquots at -20 °C. They'll be stable up to 6 months.

5. Use 2 μ L of each pI marker for every 200 μ L of sample.

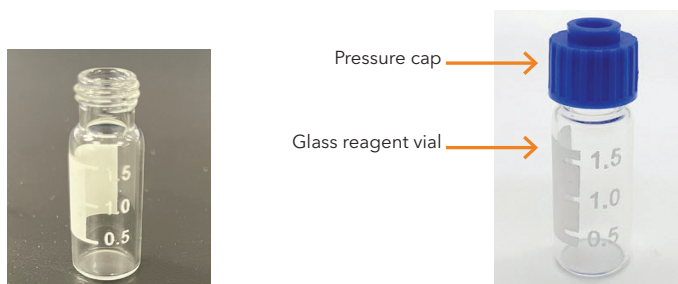
Reagents

IMPORTANT:

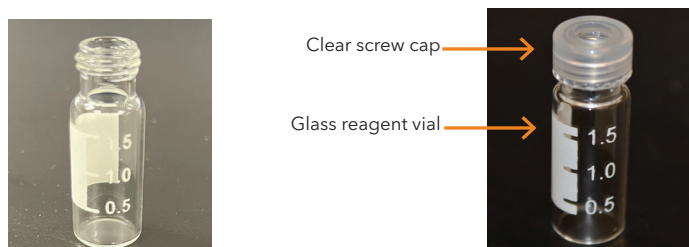
Use glass reagent vials, 2 mL (PN 046-017) to prepare cIEF batch reagents.

NOTE: Don't reuse reagents, vials or the clear screw caps. Always keep the blue pressure caps paired with their respective reagents. If you want to reuse pressure caps, first wash them thoroughly with DI water, soak overnight in DI water, rinse caps thoroughly with DI water again and air dry them before reusing.

1. Pipette 2 mL of 0.5% Methyl Cellulose into a glass reagent vial, label and close with a **blue pressure cap**.



2. Pipette 500 μ L of Fluorescence Calibration Standard in a glass reagent vial, label and close with a **blue pressure cap**.
3. Pipette 2 mL of DI water into a glass reagent vial, label and close with a **blue pressure cap**.
4. Close an empty glass reagent vial with a **blue pressure cap**.
5. Pipette 2 mL of DI water into a glass reagent vial, label and close with a **clear screw cap**.
6. If you will be doing on-board mixing, also prepare the following reagents. If not, skip to the next step.
 - a. Pipette 2 mL of DI water into a glass reagent vial, label and close with a **clear screw cap**.
 - b. Close an empty glass reagent vial with a **clear screw cap**.



- c. If you are running 48 samples or less, add 6 mL of IEF Separation Mix to one 6 mL glass reagent vial. For more than 48 samples, prepare two vials.
- d. Add 6 mL of DI water to two 6 mL glass reagent vials.

Step 2: Prep the cIEF Cartridge

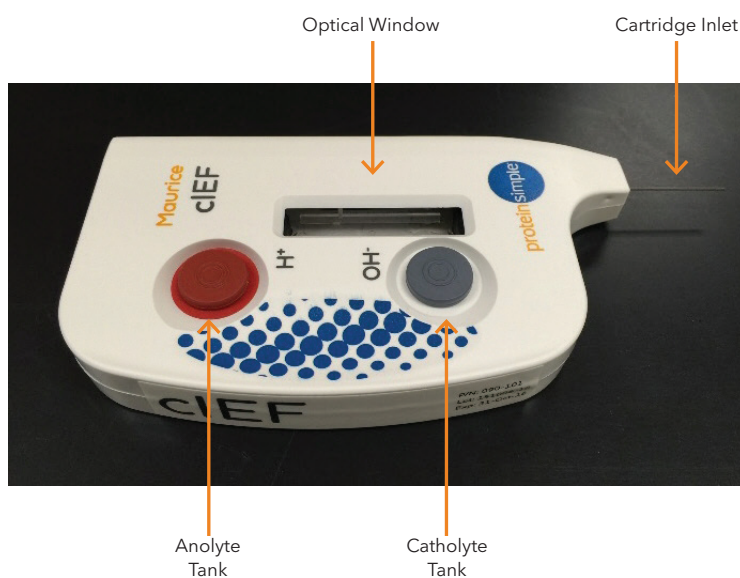
NOTE: A cIEF Cartridge is guaranteed for 100 injections and a maximum of 20 batches. Its RFID will keep track of how many are left for you. The absolute injection limit of the cartridge is 200.

1. Take the cIEF Cartridge out of its packaging. Save the packaging, you'll need it later to store the cartridge.

NOTE: When you remove the cartridge from its packaging, make sure the cartridge inlet doesn't come in contact with any surface.



2. Put the cartridge on a flat surface with its electrolyte tanks facing up.
3. Remove the stoppers from both electrolyte tanks.



4. Add 2 mL of Catholyte solution to the OH⁻ electrolyte tank (white port).
5. Add 2 mL Anolyte solution to the H⁺ electrolyte tank (red port).

NOTE: Make sure you don't overfill the electrolyte tanks.

6. Seal each tank with the rubber stoppers. Use the grey stopper for the OH⁻ tank and the red one for the H⁺ tank. If excess liquid comes out of the tank, make sure to wipe it with a laboratory wipe.



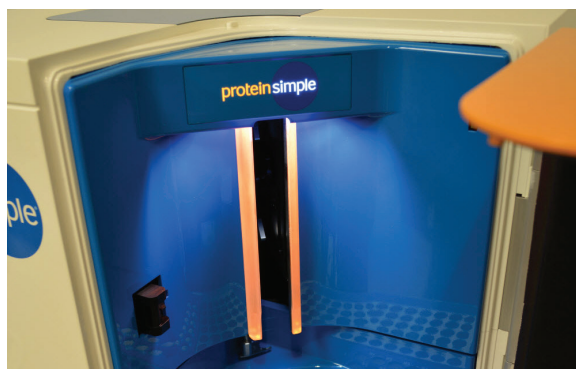
Step 3: Install the Cartridge

1. Open Maurice's door by touching the metal plate on top of the door.



NOTE: The indicator light on Maurice's front panel will blink rapidly as the door disengages.

- Swing the door open to access the cartridge slot and the reagents and samples platform. The lights on either side of the slot will be **orange**.



- Double check to make sure you've got electrolytes loaded and the tanks are properly sealed with the stoppers.
- Lift the cartridge and hold it vertically using the finger holds on either side, cartridge inlet down, with the cIEF label facing you.
- Gently insert it into the slot. You'll feel the alignment groove on the cartridge align inside the slot when you insert it.



- Continue to slide the cartridge into the slot until the locking mechanism engages. The lights on either side of the slot will change to **blue** once the cartridge is installed correctly.



Step 4: Load Samples and Reagents

1. Place the reagent vials into their respective positions on the sample and reagents platform:

NOTES:

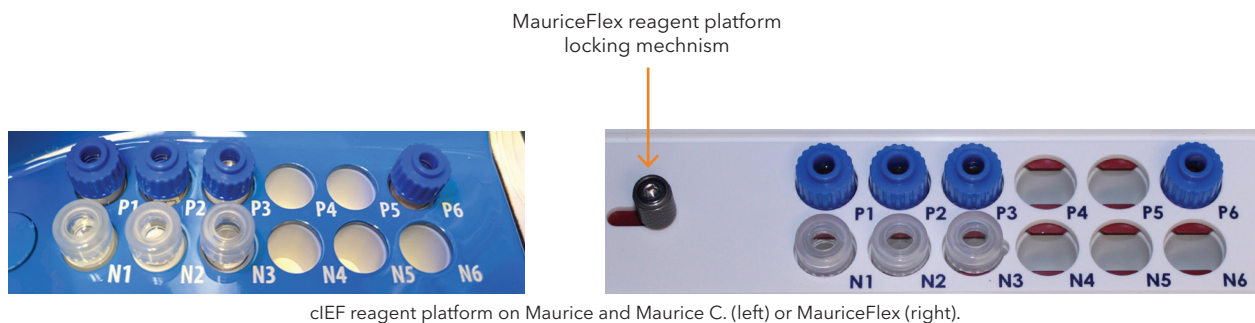
The two reagent rows in Maurice's reagent and sample platform are kept at room temperature.

Pressure caps are blue and have a raised, ringed surface. They should only be used in Reagent Row P. Use reagent vials with clear screw caps in Row N.

IMPORTANT:

Batch reagents should be prepared in glass reagent vials, 2 mL (PN 046-017).

- **P1** - 0.5% Methyl Cellulose with **blue pressure cap**
- **P2** - Fluorescence Calibration Standard with **blue pressure cap**
- **P3** - Water with **blue pressure cap**
- **P6** - Empty vial (air) with **blue pressure cap**
- **N1** - Water with **clear screw cap**



IMPORTANT:

Reagent vials on MauriceFlex must be locked in place by sliding the locking mechanism from the left to the right before starting a batch. A warning will appear in the Start Run window if they are not locked when you start the batch.

Optional: If you are doing on-board mixing, add these additional reagents:

- **N2** - Water with **clear screw cap**
- **N3** - Empty vial with **clear screw cap**
- **M1** - 6 mL IEF Separation Mix

- **M2** - N/A for 48 samples or less, add a 6 mL IEF Separation Mix vial when running more than 48 samples
- **M3** - 6 mL water
- **M4** - 6 mL water

NOTE: On-board mixing is only available on Maurice and Maurice C. systems.

Depending on what you prepared your samples in, put your 96-well sample plate in the metal plate insert or your sample vials in the metal vial insert.



OBM Reagents M1-M4

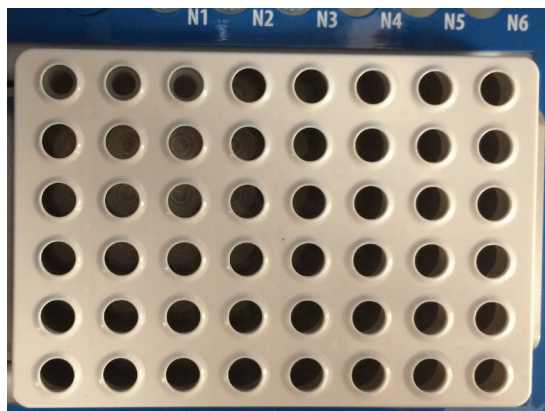
NOTES:

If you need to seal the 96-well plate during your run, we recommend the Slit Seal, 96 well plate seal from BioChromato (<https://biochromato.com/slit-seal/>). It can be used in both absorbance and native fluorescence modes. If you're currently using X-Pierce adhesive film (PN XP-100, Excel Scientific), we recommend using it in absorbance mode only.

Plate seals should NOT be used when running CE-SDS mode.

Well A1 on the 96-well plate should be in the top left corner of the insert.

2. If you are using a vial tray, place the condensation lid on top of the vials.



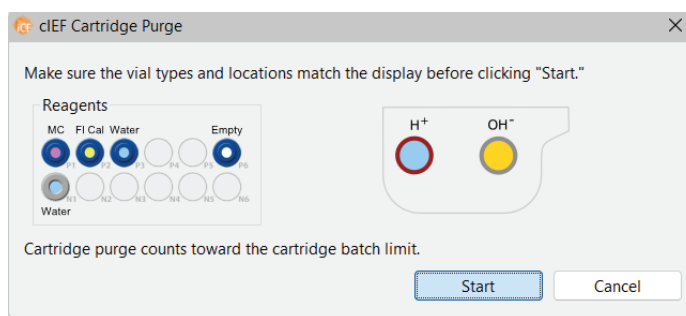
3. Close the instrument door. Maurice locks it automatically.

Step 5: Check for Cartridge Alerts

1. If your cartridge was last used in a run with an error you will need to perform a Cartridge Purge. The Maurice system will remind you if a Cartridge Purge is required.
 - a. To purge the cartridge, click the brown **Purge** button in the instrument status bar.



- b. Confirm that the required batch reagents are loaded and that the cartridge is prepped. Then click **Start**.



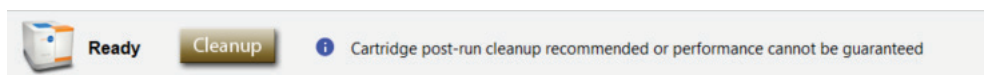
- c. Once the purge is done, perform a Cartridge Post-Run Cleanup by clicking on the brown **Cleanup** button that will appear in the instrument status bar. See step 2 for more info.

NOTE: If there's only one batch left on the cartridge, you'll get a message that it can't be used after purging.

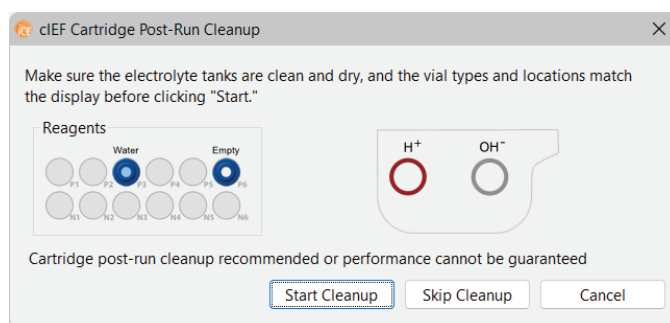
2. If you just completed a Cartridge Purge or if the cartridge you're using did not undergo a cartridge post-run cleanup after the last run, you will be asked if you want to perform a Cartridge Post-Run Cleanup.

To start the Post-Run Cleanup:

- Remove the cartridge from Maurice and remove the anolyte and catholyte from the cartridge. See "Post-batch Procedures" on page 120 for more info.
- Confirm there is a vial of Water (P3) and Air (P6) in place.
- Click the brown **Cleanup** button in the instrument status bar.



- Click **Start Cleanup** in the prompt that appears.

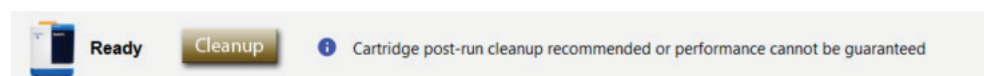


NOTE: You can also start a Post-Run Cleanup by selecting **Instrument > Cartridge Post-Run Cleanup**. When you start a Cartridge Post-Run Cleanup from the instrument menu, click **Start** in the prompt that appears to begin the cleanup.

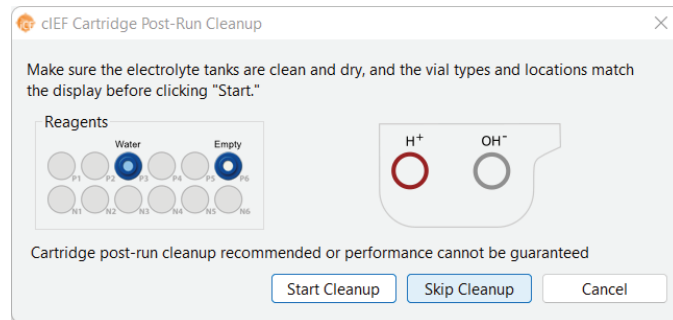
- Once the cartridge post-run cleanup is completed, re-prepare the cartridge and install the cartridge in Maurice. See "Step 2: Prep the cIEF Cartridge" on page 96 for more information.
- Click on the green **Start** button in the instrument status menu to start the batch.

To start the run without a Post-Run Cleanup:

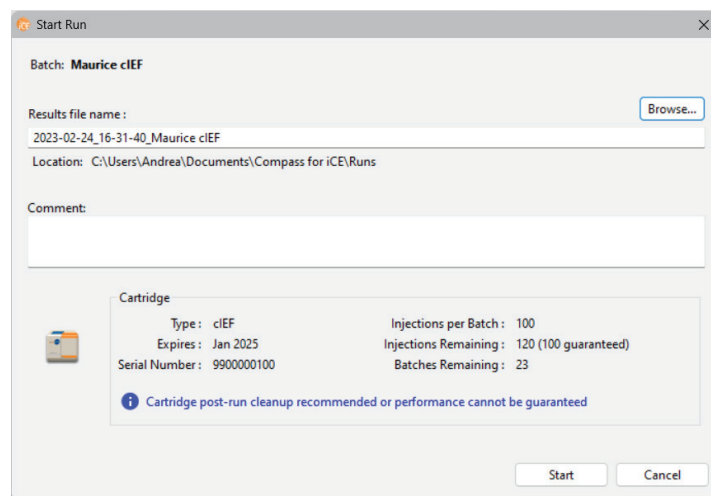
- Click the brown **Cleanup** button in the instrument status bar



- b. Click **Skip Cleanup** in the prompt that appears.



- c. Click **Start** in the Start Run window. The run summary and injection report will record that the cartridge did not undergo a post-run cleanup before starting the batch.

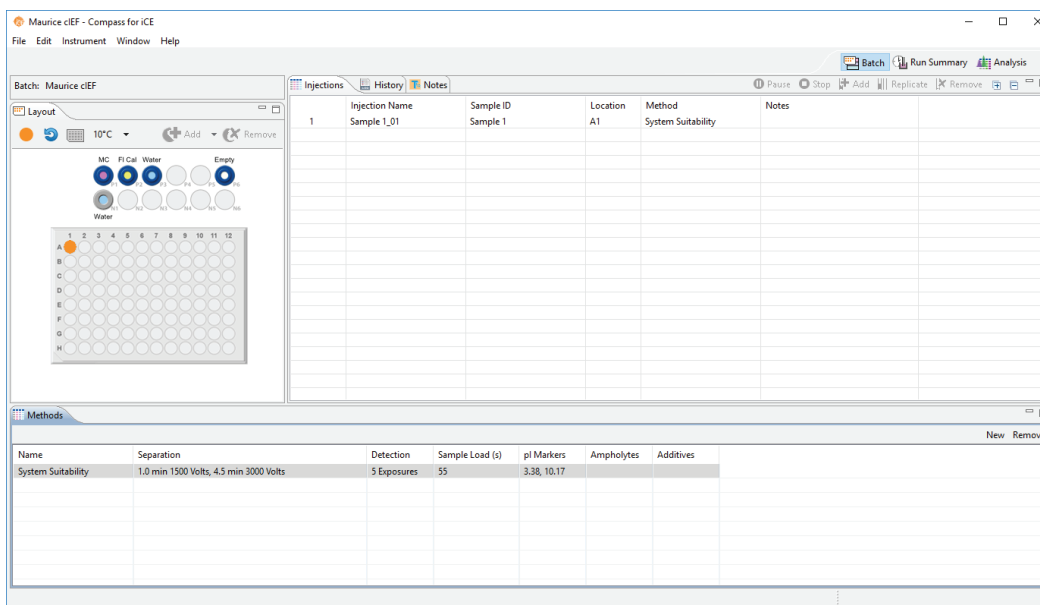


To start the run with a different cartridge:

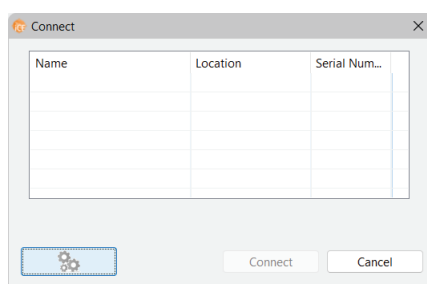
- If necessary, click **Cancel** in the cIEF Cartridge Post-Run Cleanup window.
- Open Maurice's door, remove the first cartridge from Maurice and prepare a second cartridge. See "Step 2: Prep the cIEF Cartridge" on page 96 for more information.

Step 6: Create a Batch

1. Launch Compass for iCE.
2. Select the **Batch** tab. This is where you'll enter sample/injection information, methods and batch parameters.



3. To create a batch, make sure your Maurice system is connected to Compass for iCE. Select **Instrument** and click **Connect**.
 - a. If your instrument is listed, select your Maurice system and click **Connect**.
 - b. If your instrument isn't listed, click on the **Settings** button and connect by typing in your instrument IP address.



To create a new batch:

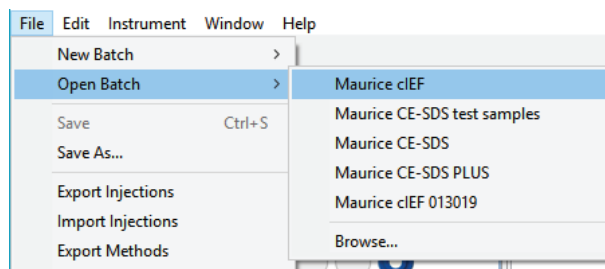
- On Maurice systems - in the main menu, select **File > New Batch > Maurice cIEF**
- On Maurice C. systems - in the main menu, select **File > New Batch**
- On MauriceFlex systems - in the main menu, select **File > New Batch > Maurice cIEF**

To use an existing batch: In the main menu, select **File > Open Batch**.

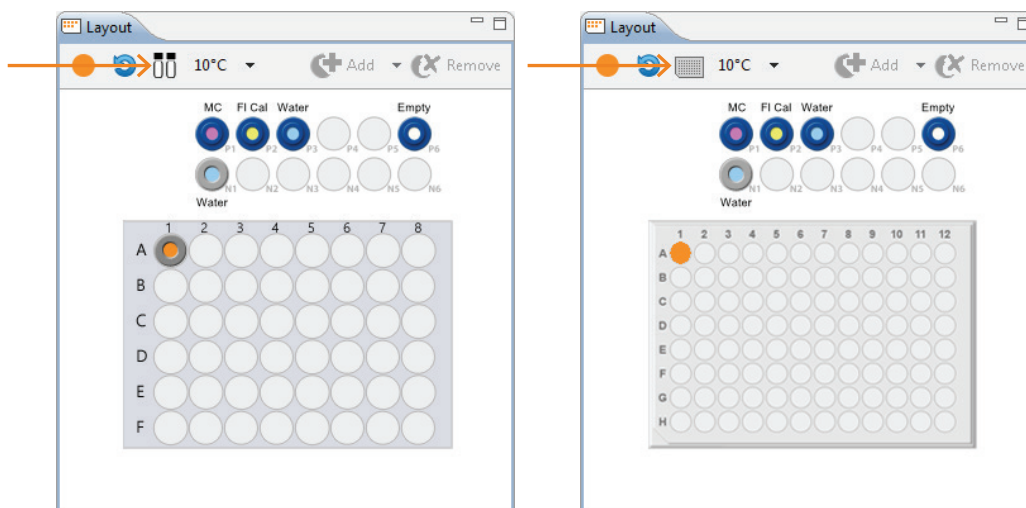
NOTES:

cIEF batches that use the FL458 nm fluorescence detection filter can only be opened in Compass for iCE 2.1 or later.

If you're making changes to an existing batch, follow the steps in this section. Otherwise skip to "Step 7: Start the Batch" on page 116.



4. In the Layout pane, clicking on the vial/plate icon toggles between formats. Select 48 vials or a 96-well plate depending on what you're running.



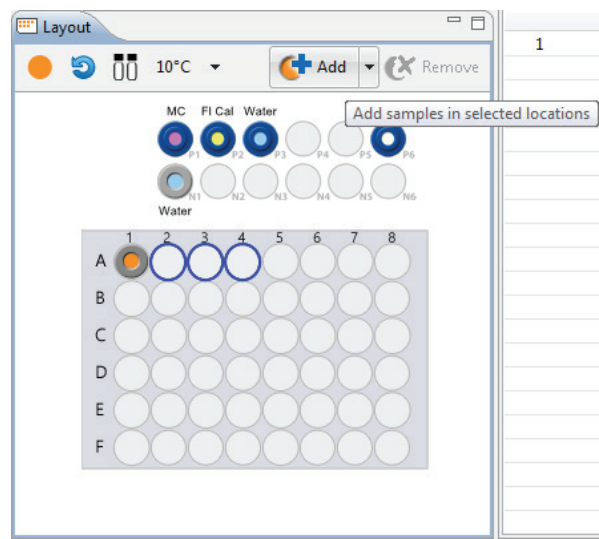
5. Add your samples:

To import samples using a saved injections file:

- a. Select **File** in the main menu and click **Import Injections**.
- b. Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.

NOTE: For more information on setting up an injections file for import, see "Importing and Exporting Injections" on page 84.

To select samples manually: Use your mouse to highlight the well positions for each of your samples, then click Add.

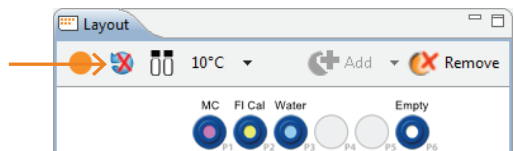


This populates the Injections table:

Injections						
History Notes						
	Injection Name	Sample ID	Location	Method	Notes	
1	Sample 1_01	Sample 1	A1	System Suitability		
2	Sample 02_02	Sample 02	A2	System Suitability		
3	Sample 03_03	Sample 03	A3	System Suitability		
4	Sample 04_04	Sample 04	A4	System Suitability		

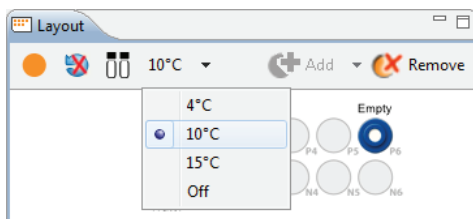
6. Compass for iCE can monitor the current during a separation for you, stop it if an abnormal current profile is detected and reinject the sample for you automatically. Reinjections are enabled by default, but if you don't want to use this feature just click the reinject icon to toggle it off.

NOTES:
If a sample was reinjected, the injection row will be flagged in the Injections pane in the Run Summary screen. See “Injection Flags” on page 414 for more info.
A maximum of 10 reinjections are allowed per batch.



7. The sample platform is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The two reagent rows are kept at room temperature.



8. Enter your method parameters in the Methods pane.

NOTE: We recommend using the default method parameters. Please see the Maurice cIEF Method Development Guide for more information on method optimization.

To import a saved method:

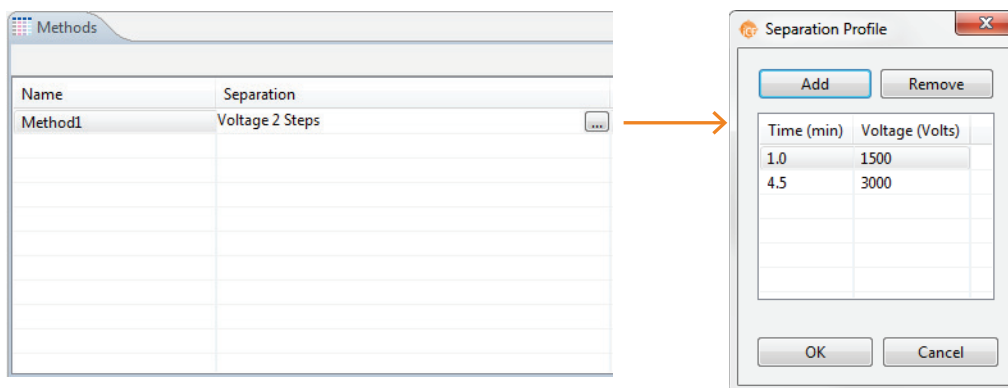
- a. Select **File** in the main menu and click **Import Method**.
- b. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

To create a new method or make changes to an existing one:

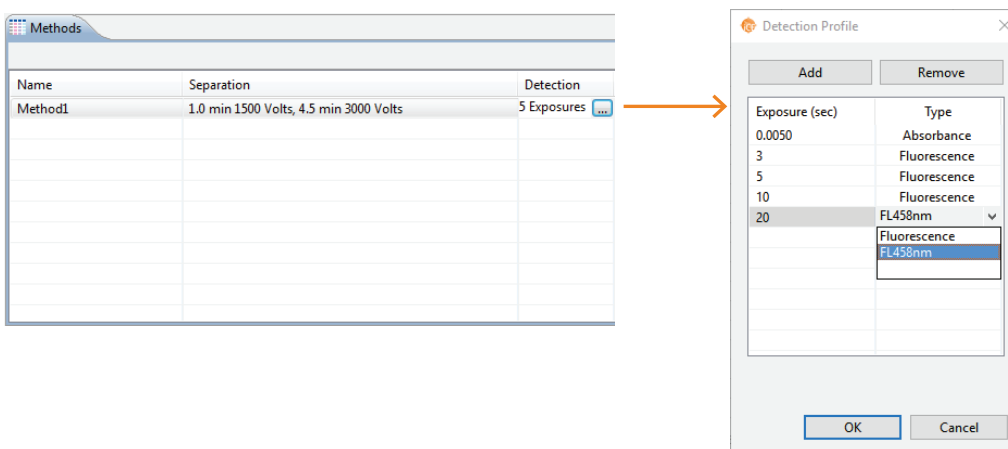
- a. Click the first cell in the Name column and enter a method name.

Methods						
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method 1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17		

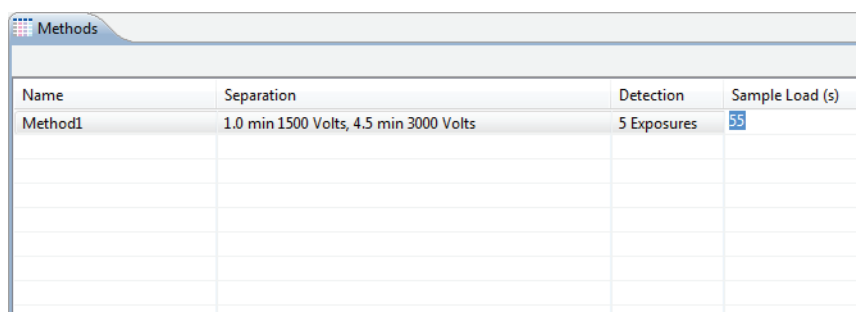
- b. Click the first cell in the Separation column, then click the selection button [...] to set the pre-focusing and focusing time (in minutes) and voltage.



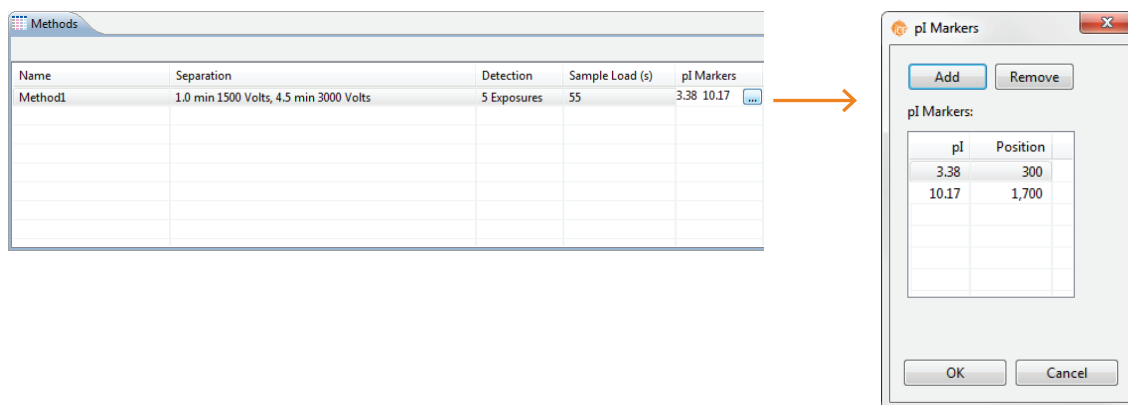
- c. Click the first cell in the Detection column then click the selection button [...] to set your exposure times for absorption and fluorescence detection modes. You can also select between native or FL458 nm fluorescence if the optional FL458 nm filter is installed on your system.



- d. Click the first cell in the Sample Load(s) column and set the load time in seconds.



- e. Click the first cell in the pI Markers column to select pI markers. Add new markers or remove existing ones then click **OK**.



- f. Optional: Click the first cell in the Ampholytes column and enter the ampholytes you're using.

Methods						
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17	Pharmalyte 3-10	

- g. Optional: Click the first cell in the Additives column and enter any additives you're using.

Methods						
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives
Method1	1.0 min 1500 Volts, 4.5 min 3000 Volts	5 Exposures	55	3.38, 10.17	Pharmalyte 3-10	Urea

9. You can now:

- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.

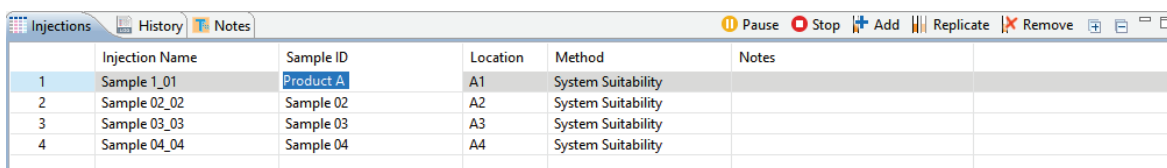
10. In the Injections pane:

- **To add or change sample names:** Click the **Sample ID** cell for the injection and type a name.

NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 82 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.



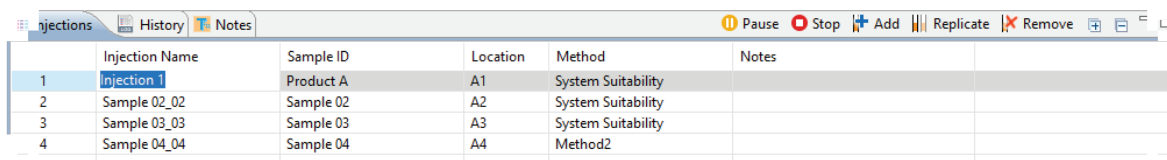
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	System Suitability	

- **To change injection names:** Click the **Injection Name** cell for the injection and type a name.

NOTES:

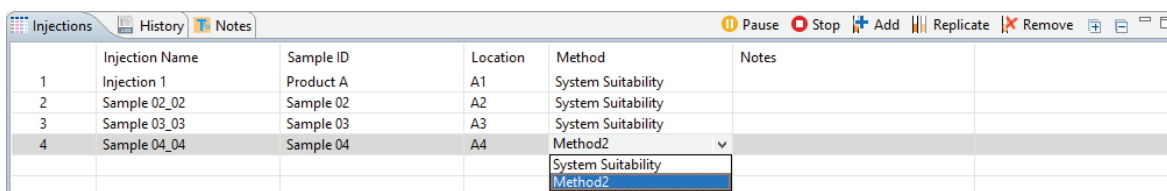
Each injection name must be unique.

Changing the injection name won’t affect the sample ID.



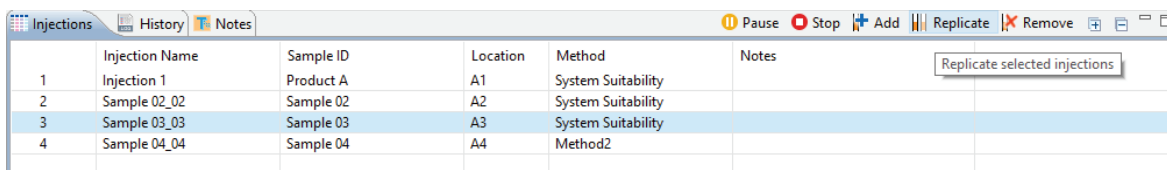
	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	

- **To assign methods for each injection:** Click the **Method** cell for the injection and select a method from the drop down menu.

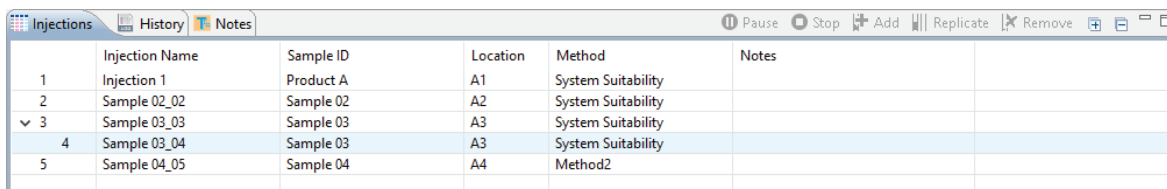


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

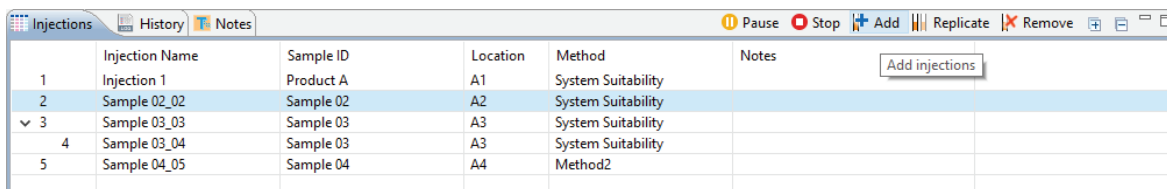


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 04_04	Sample 04	A4	Method2	



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 03_04	Sample 03	A3	System Suitability	
5	Sample 04_05	Sample 04	A4	Method2	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	System Suitability	
3	Sample 03_03	Sample 03	A3	System Suitability	
4	Sample 03_04	Sample 03	A3	System Suitability	
5	Sample 04_05	Sample 04	A4	Method2	

- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

11. If your Maurice has the on-board mixing (OBM) option, you can tell Compass for iCE which samples to automatically mix for you. Otherwise you can skip to the next step. Up to 96 samples can be on-board mixed in a batch. You can also have a mix of premixed samples and samples to on-board mix in the same batch.

NOTES:

OBM is only available on Maurice and Maurice C. systems.

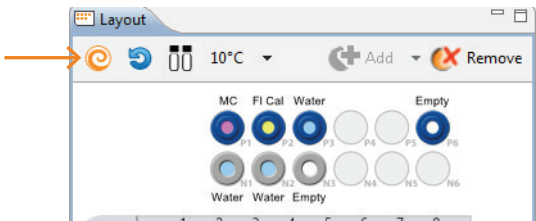
To use the OBM function, you must load 25 µL of sample in every sample vial or plate well that will be mixed on board. Maurice will automatically mix 25 µL of each sample with 100 µL of IEF Separation Mix prior to injection.

a. Make sure you've placed the following on-board mixing reagents in Maurice:

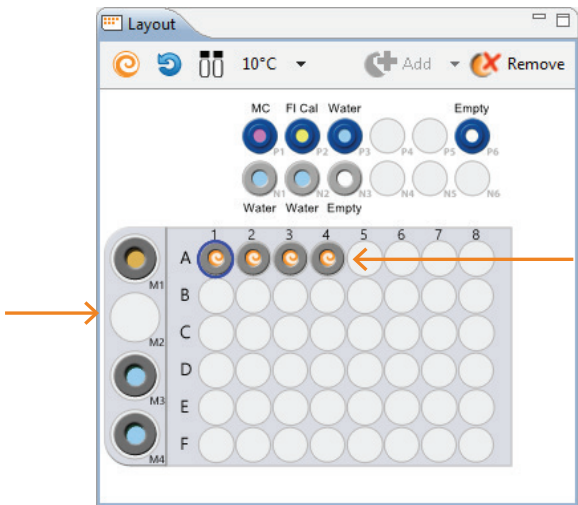
- **M1** - 6 mL IEF Separation Mix vial
- **M2** - 6 mL IEF Separation Mix vial (use only when running more than 48 samples)
- **M3** - 6 mL water vial

- **M4** - 6 mL water vial

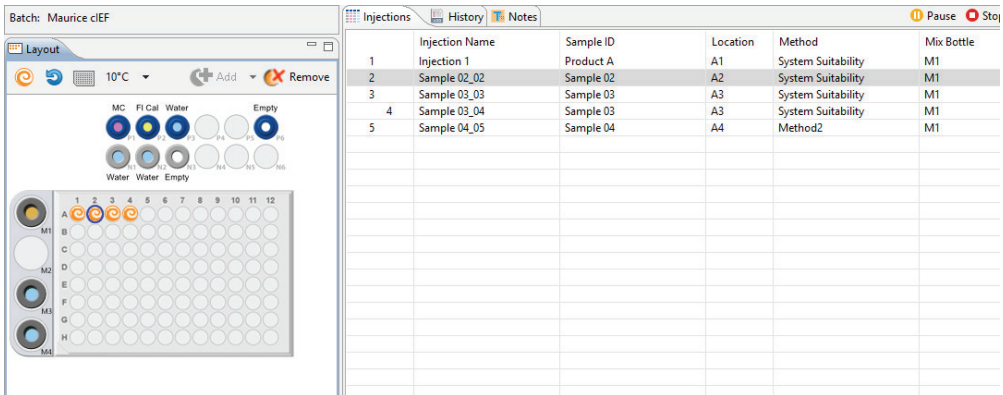
b. The on-board mixing function is disabled by default. Click the **on-board mixing icon** to toggle it on.



All sample positions will now display the mixing icon, and the on-board mixing reagent vial positions will also display on the left:



The Injection pane will also display a Mix Bottle column showing the position of the IEF Separation Mix vial that will be used for each sample. The vial in the M1 position is used for samples 1–48, and the vial in M2 will automatically display and be used for samples 49–96:



Batch: Maurice cIEF

Layout

10°C

Add Remove

MC FI Cal Water Empty

P1 P2 P3 P4 P5 P6

Water Water Empty

N1 N2 N3 N4 N5 N6

A 1 2 3 4 5 6 7 8 9 10 11 12

M1

B

M2

C

M3

D

M4

E

F

G

H

Injections History Notes

Pause Stop

	Injection Name	Sample ID	Location	Method	Mix Bottle
40	Sample 38_40	Sample 38	D3	System Suitability	Mix Bottle
41	Sample 39_41	Sample 39	D4	System Suitability	M1
42	Sample 40_42	Sample 40	D5	System Suitability	M1
43	Sample 41_43	Sample 41	D6	System Suitability	M1
44	Sample 42_44	Sample 42	D7	System Suitability	M1
45	Sample 43_45	Sample 43	D8	System Suitability	M1
46	Sample 44_46	Sample 44	D9	System Suitability	M1
47	Sample 45_47	Sample 45	D10	System Suitability	M1
48	Sample 46_48	Sample 46	D11	System Suitability	M1
49	Sample 47_49	Sample 47	D12	System Suitability	M1
50	Sample 48_50	Sample 48	E1	System Suitability	M2
51	Sample 49_51	Sample 49	E2	System Suitability	M2
52	Sample 50_52	Sample 50	E3	System Suitability	M2
53	Sample 51_53	Sample 51	E4	System Suitability	M2
54	Sample 52_54	Sample 52	E5	System Suitability	M2
55	Sample 53_55	Sample 53	E6	System Suitability	M2
56	Sample 54_56	Sample 54	E7	System Suitability	M2
57	Sample 55_57	Sample 55	E8	System Suitability	M2
58	Sample 56_58	Sample 56	E9	System Suitability	M2
59	Sample 57_59	Sample 57	E10	System Suitability	M2
60	Sample 58_60	Sample 58	E11	System Suitability	M2

c. To turn on-board mixing off for specific samples, select those well(s) or vial(s), right click and select **Mixing Off**.

Layout

10°C

Add Remove

MC FI Cal Water Empty

P1 P2 P3 P4 P5 P6

Water Water Empty

N1 N2 N3 N4 N5 N6

A 1 2 3 4 5 6 7 8

M1

B

M2

C

M3

D

M4

E

F

Remove Samples

Mixing On

Mixing Off

The sample positions will display solid orange indicating mixing is off, and the Mix Bottle designation in the Injections pane is removed:

Batch: Maurice cIEF

Layout

10°C

Add Remove

MC FI Cal Water Empty

P1 P2 P3 P4 P5 P6

Water Water Empty

N1 N2 N3 N4 N5 N6

A 1 2 3 4 5 6 7 8 9 10 11 12

M1

B

M2

C

M3

D

M4

E

F

G

H

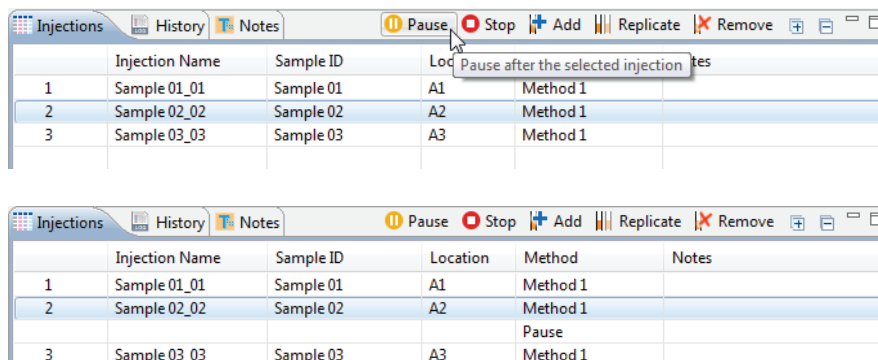
Injections History Notes

Pause Stop

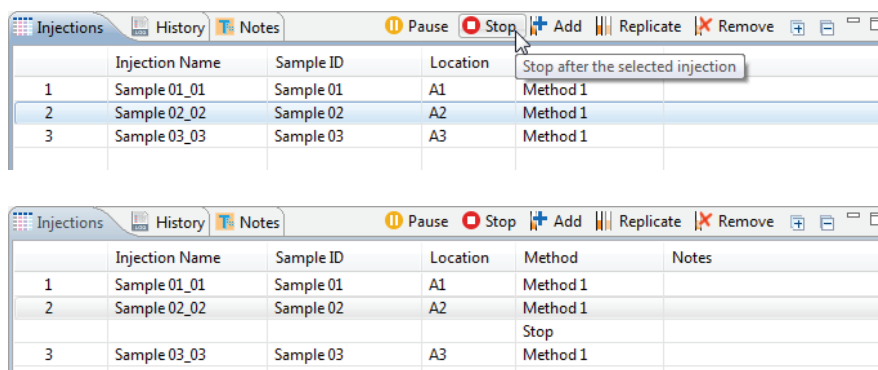
	Injection Name	Sample ID	Location	Method	Mix Bottle
1	Injection 1	Product A	A1	System Suitability	M1
2	Sample 02_02	Sample 02	A2	System Suitability	M1
3	Sample 01_03	Sample 01	A3	System Suitability	
4	Sample 03_04	Sample 03	A4	System Suitability	

12. Add programmed pauses or stops in the batch to automatically pause or stop after injections (optional).

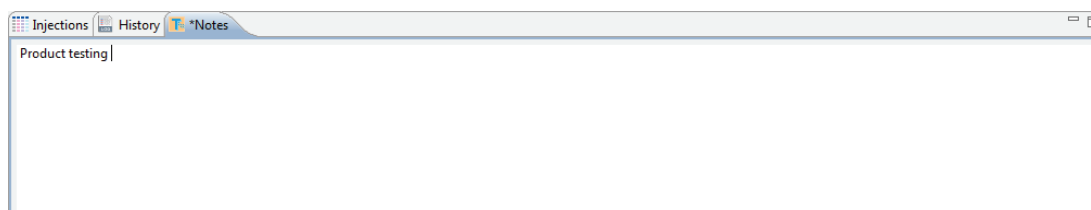
- **To program a pause:** Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete. To resume the batch, click **Continue** in the instrument status bar.



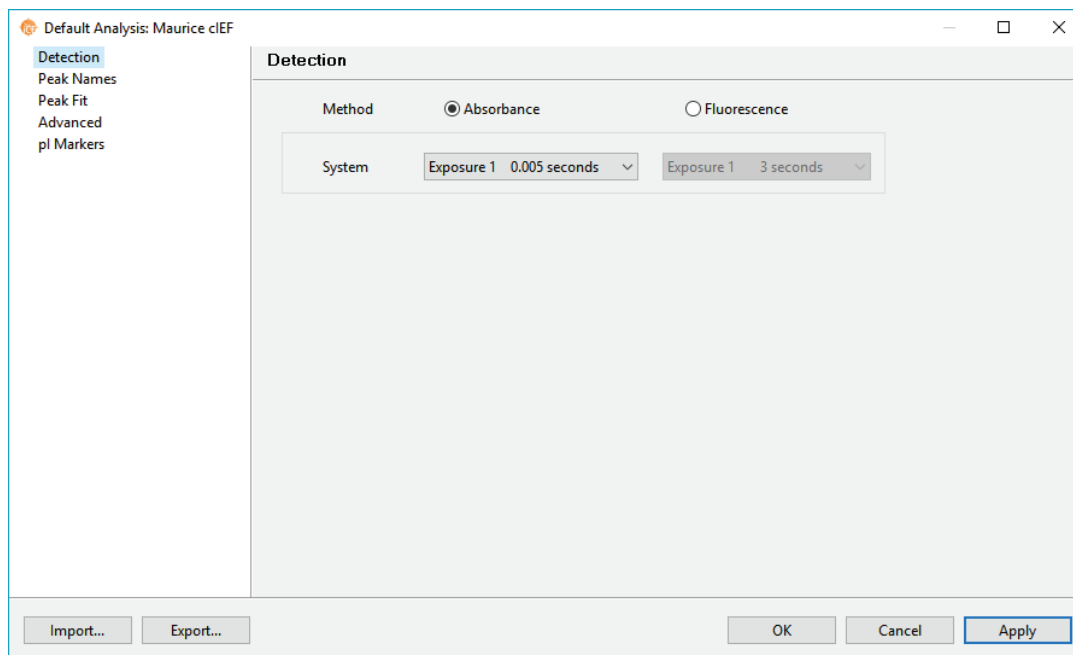
- **To stop the run after a specific injection:** Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.



13. Click on the **Notes** pane, then click in the notes area and type any information you want to add about your batch (optional).



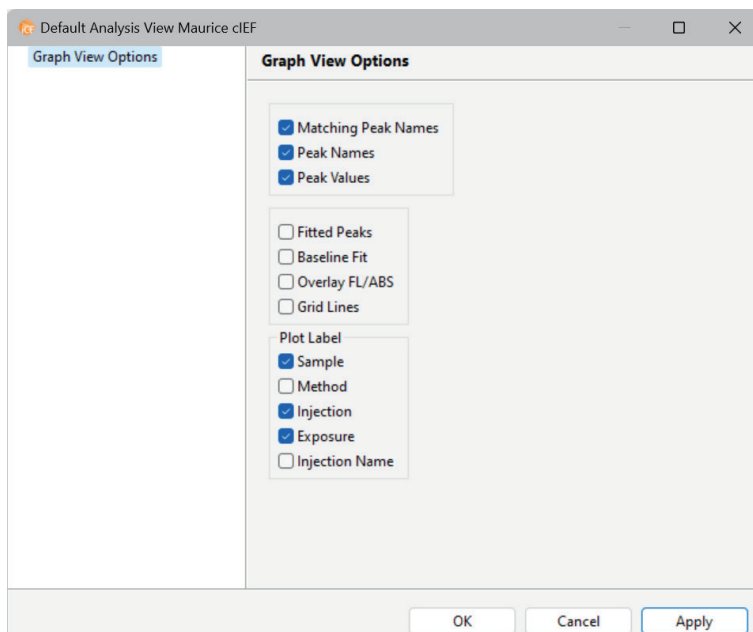
14. You can preset the parameters used to analyze data generated with the batch (optional). We recommend using the default parameters for cIEF applications, but if you want to modify parameters:
- Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



- Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to "Analysis Settings Overview" on page 528.

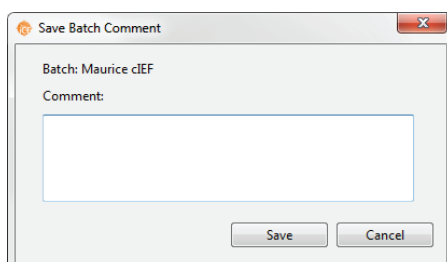
15. You can preset the parameters used to view data when the batch is complete (optional). Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button:

- a. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



- b. Change the parameters you want to, then click OK. For detailed graph view options, please refer to “Customizing the Data Display” on page 509.

16. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



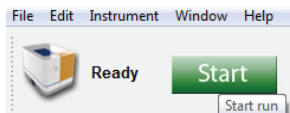
17. Enter a name for your batch then click **Save**.

Step 7: Start the Batch

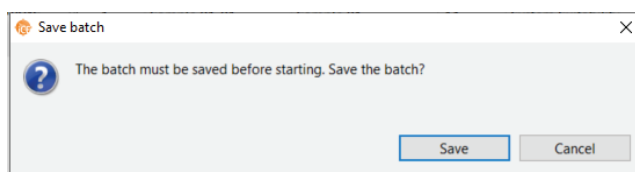
1. Make sure Maurice is connected to Compass for iCE. If he isn't, select **Instrument** and click **Connect**. Then select your Maurice system and click **Connect**.

- Click on the green **Start** button to start your batch.

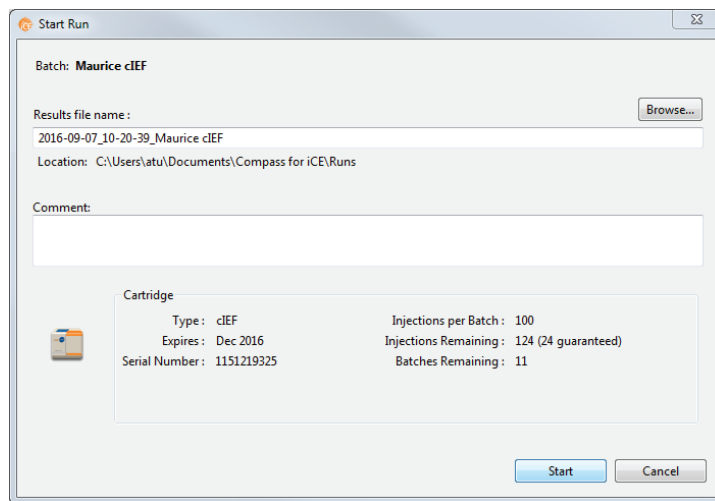
NOTE: An alert may appear if the wrong cartridge is installed or if a cartridge post-run cleanup or cartridge purge is recommended. See “Step 5: Check for Cartridge Alerts” on page 101 for more information.



- If you have made any changes, you will be asked to save your batch before starting the run. Click **Save**



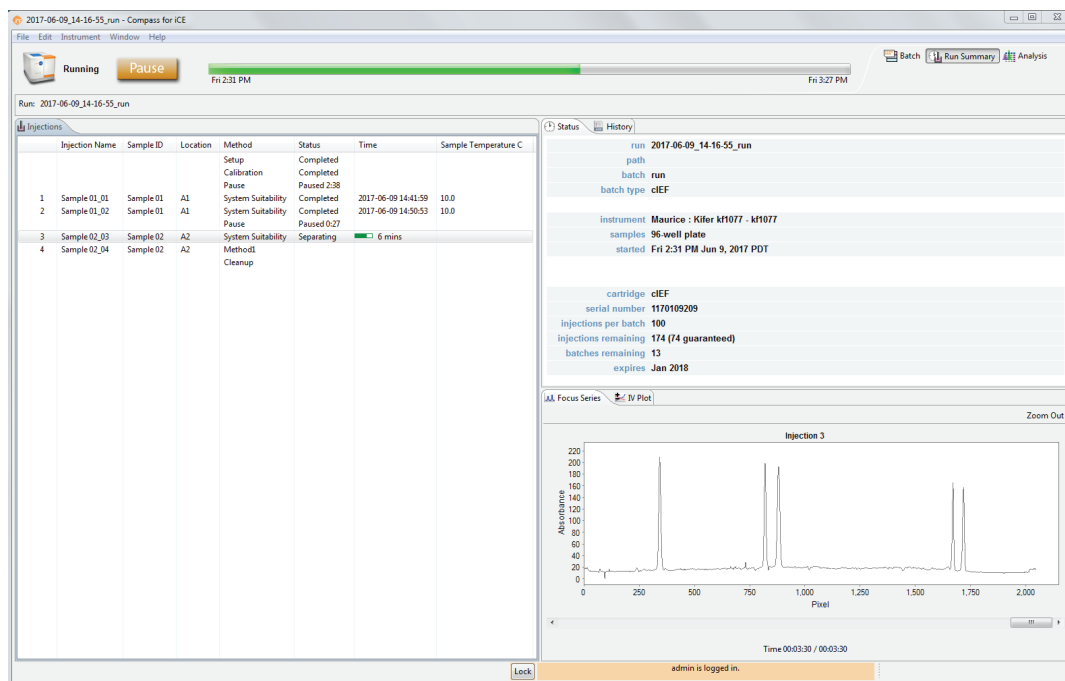
- The Start Run box shows you the number of injections and batches remaining on your cartridge. Make sure you still have enough injections left before you start the run. If not, prep a new cartridge.
- Click in the **Results File name** box if you want to change the default run file name. Otherwise just leave it as is.



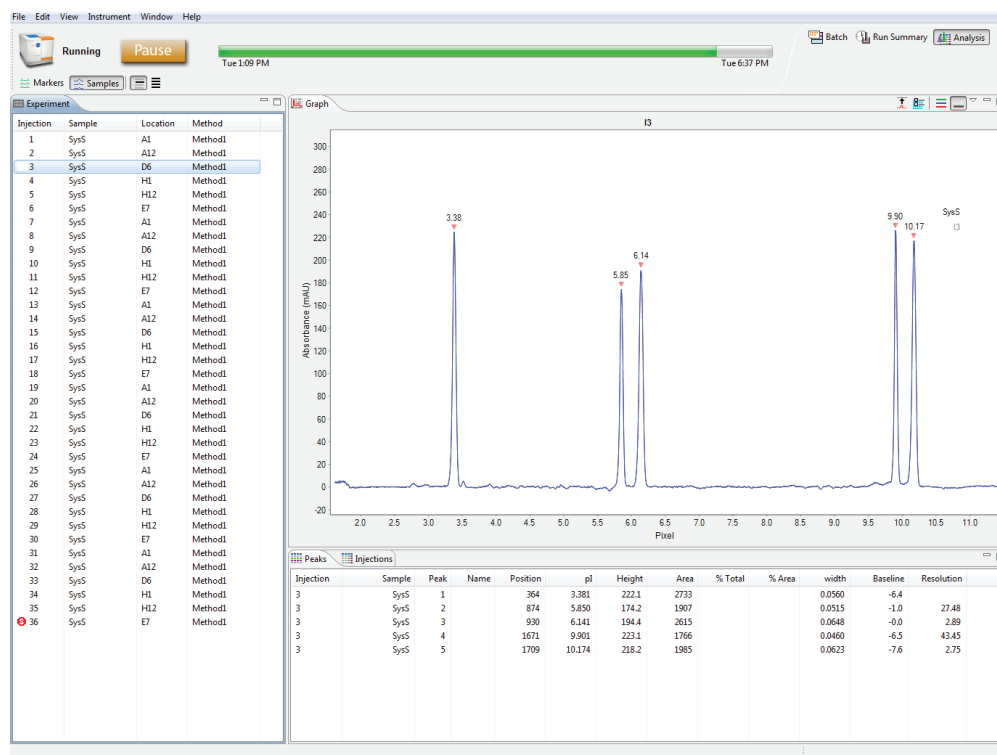
- If you don't want to save the file to the default Runs folder, click **Browse** to select a different location.
- Enter any run details you'd like in the Comments box (optional).
- Click **Start** to start the run.

NOTE: The indicator light on Maurice's front panel will pulse slowly while he's running the batch.

To monitor the progress of your batch, click the **Run Summary** tab. See Chapter 16: “Run Status” for more details.



To view results, analyze results or change analysis parameters on completed injections while the batch is still running, click the **Analysis** tab. See Chapter 18: “cIEF Data Analysis” for more details.



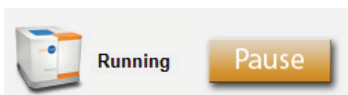
When your batch is complete, you can view electropherograms for all injections in the Analysis tab, and all batch and injection details in the Run Summary tab.

NOTE: If you'd like Maurice to let you know when your run is done, you can set him up to tweet you. Go to "Setting Up Maurice Systems to Send Tweets" on page 764 for more info.

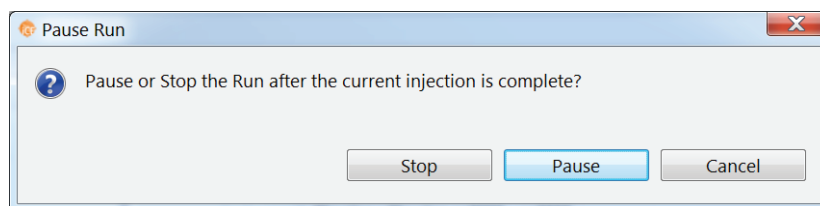
Editing a Running Batch

You can pause a batch after it's started to make updates or add samples and reagents.

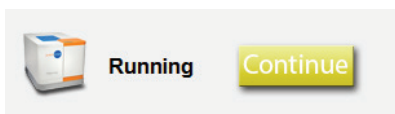
1. Select **Instrument > Pause** or select the **Pause** button.



2. Click **Pause** in the pop-up window.



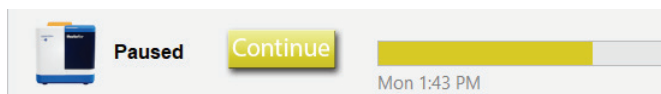
The Pause button will change to Continue but Maurice's status will stay at Running until Compass for iCE completes the current injection.



- You can now make these edits to the batch:
 - Add sample(s)
 - Add injection(s) to the end of the batch
 - Modify injections that haven't completed and aren't greyed out
 - Create a new method
 - Update the method. Once a method is run it can only be updated. Highlight the method you want to edit and click **Update** in the Methods pane. The software will copy the parameters into a new method that you can modify.

3. When Maurice completes the current injection, his status will change to Paused and the progress bar will turn yellow.

NOTE: Maurice's status light will change from a pulsing blue to a blue pulse with magenta flashes which means he's paused the run.



The status in the Injection pane also changes to Paused and the pause duration updates live. You can now open the instrument door to:

- Add new samples
- Add on-board mixing reagents if the paused batch was set up to do on-board mixing and your Maurice has this option.

IMPORTANT

The door must be closed within 15 minutes of opening it to avoid potential assay performance issues.

NOTE: If the door is open for more than 30 seconds, a white light near the cartridge will start to flash and a warning message displays in Compass for iCE until the door is closed.

4. To continue the batch, click **Continue**. Changes you made to the batch are saved and Maurice will continue with the next sample injection.

Post-batch Procedures

When the batch is done:

1. Open Maurice's door. The lights on either side of the cartridge slot will be **orange** as Maurice will have already disengaged the cartridge.
2. Remove your sample. Leave the Water (P3) and Air (P6) vials in place if your cartridge still has injections left since they will be needed for the post-run cleanup step. Remove the remaining reagent vials.

3. Gently pull out the cartridge making sure the cartridge inlet doesn't come into contact with any surface.



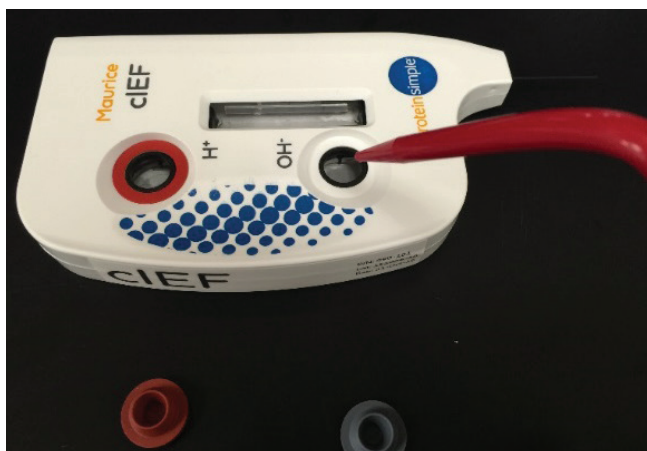
If you're at 100 injections, you've reached the limit of guaranteed performance for the cartridge. To dispose of a finished cartridge, put it in its original packing and discard it per your institution's safety and waste disposal guidelines. Discard any remaining reagent vials.

NOTE: Disposal of cartridges, sample plates and vials depends on the samples you ran. If you aren't sure what your sample origins are, we recommend you dispose of everything in biohazard waste.

If the cartridge will be used again, clean up and store the cartridge:

- a. Put the cartridge on a flat surface with its electrolyte tanks facing up.
- b. Remove the stoppers from both the electrolyte tanks.
- c. Using an electrolyte pipette or low vacuum, aspirate the solutions from each tank.
- d. Fill each tank with 2 mL DI water, then aspirate it out. Repeat this rinse 3 times.

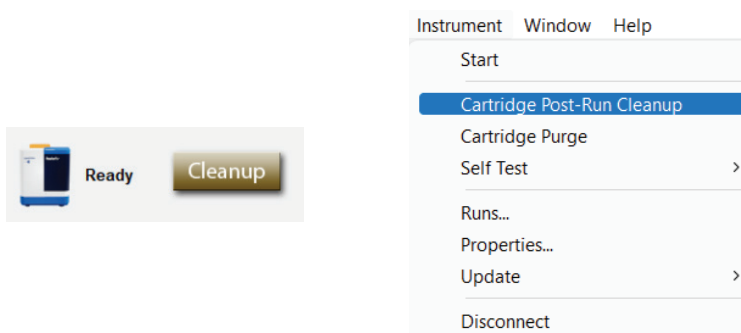
NOTE: Make sure not to get any liquid on the cartridge's optical window.



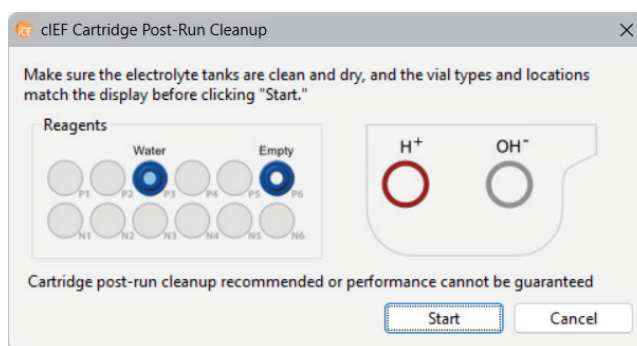
- e. Aspirate all the remaining liquid and make sure that the tanks are dry.

NOTE: If you are using Compass for iCE v3.0.0, go to step f. If you are using an earlier version of Compass for iCE, remove all the reagent vials and skip to step k.

- f. Put the stoppers back on the tanks and install the cartridge in Maurice.
- g. Verify there is at least 1.5 mL of DI water in the Water (P3) vial and air in the empty (P6) vial.
- h. Click on the brown **Cleanup** button in the Instrument Status Bar or select **Instrument** and **Cartridge Post-Run Cleanup** in the Compass main menu.



- i. You'll get the following message. Click **Start**. It'll only take 6 minutes.



- j. Open Maurice's door. Gently pull out the cartridge making sure the cartridge inlet doesn't come in contact with any surface.
- k. Remove the reagent vials.
- l. Leave the stoppers off to allow the tanks to air dry.
- m. Put the cartridge and stoppers back in the protective packaging and store at room temperature.

!WARNING! SHARPS HAZARD

The capillary inlet of the cartridge may present a potential sharps hazard. Dispose of used cartridges according to your organizations health and safety regulations.

**!WARNING! BIOHAZARD**

Cartridges, sample plates and vials should be handled by procedures recommended in the CDC/NIH manual: Biosafety in Microbiological and Biomedical Laboratories (BMBL). The manual is available from the U.S. Government Printing Office or online at <http://www.cdc.gov/biosafety/publications/bmbl5/>.

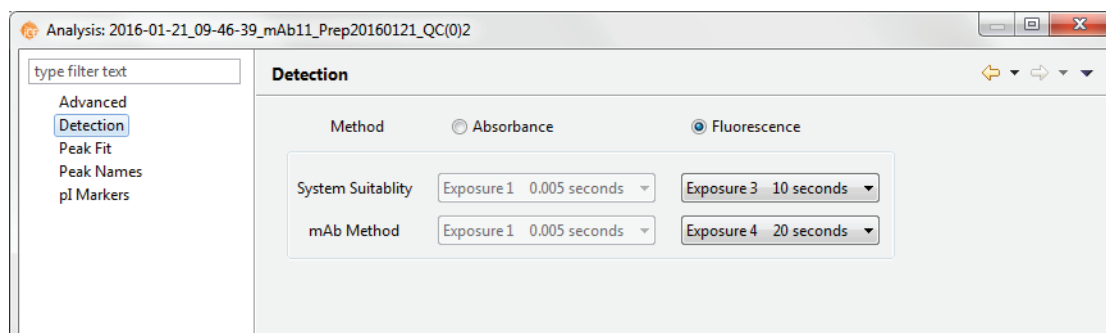
Depending on the samples used, the cartridges, plates and vials may constitute a chemical or a biohazard. Dispose of the cartridges, plates and vials in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations. Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste content before you store, handle, or dispose of chemical waste.

Checking Your Data

Compass for iCE detects your sample protein and pI marker peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

Step 1: Select Your Detection Mode

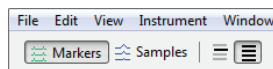
1. Go to the **Analysis** screen and open your run (if it isn't already open).
2. The data displays in absorbance mode by default. If you want to look at fluorescence data instead, select **Edit** from the main menu and click **Analysis**. In the Analysis window, select **Detection** in the left sidebar, then click **Fluorescence** in the Detection page.



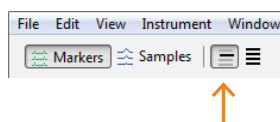
Step 2: Check Your pI Markers

To make sure your pI markers are identified correctly:

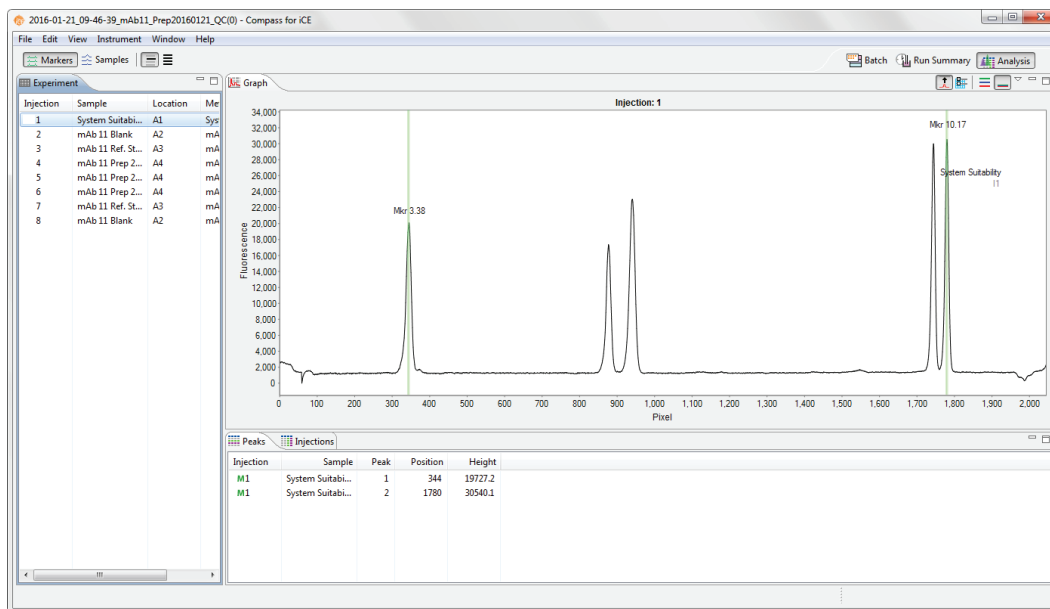
1. Go to the **Analysis** screen.
2. Click **Markers** in the View bar.



3. Click the **View Selected** icon in the View bar.

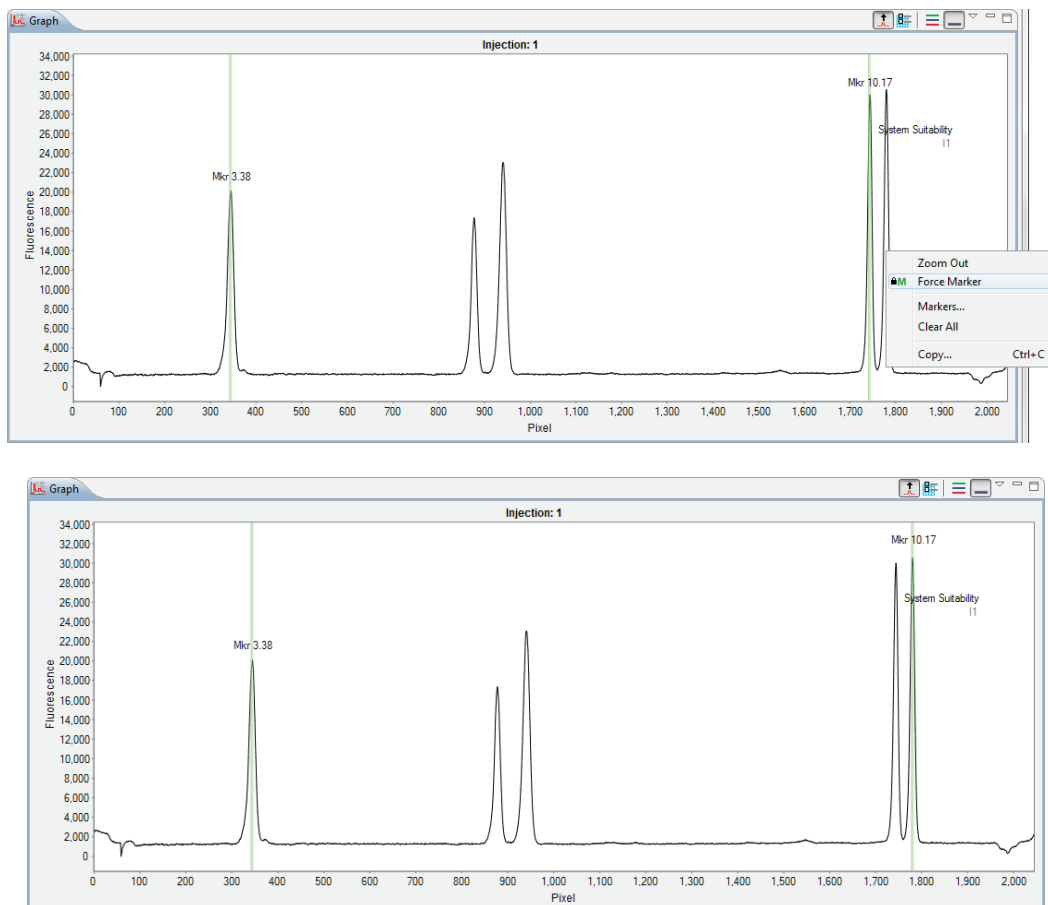


4. Click **Injection 1** in the Experiment pane.
5. Check that your pI markers in the electropherogram have been correctly identified. Each marker peak will have a green vertical line running through it and be labeled Mkr. They're also identified with an **M** in the Peaks table.



6. If your pI markers aren't identified correctly, here's how to manually correct them:

To set an unidentified peak as a pI marker: Right-click the peak in the electropherogram or Peaks table and select **Force Marker**. Compass will assign that peak as a pI marker, and correctly reassign the remaining pI marker peaks.



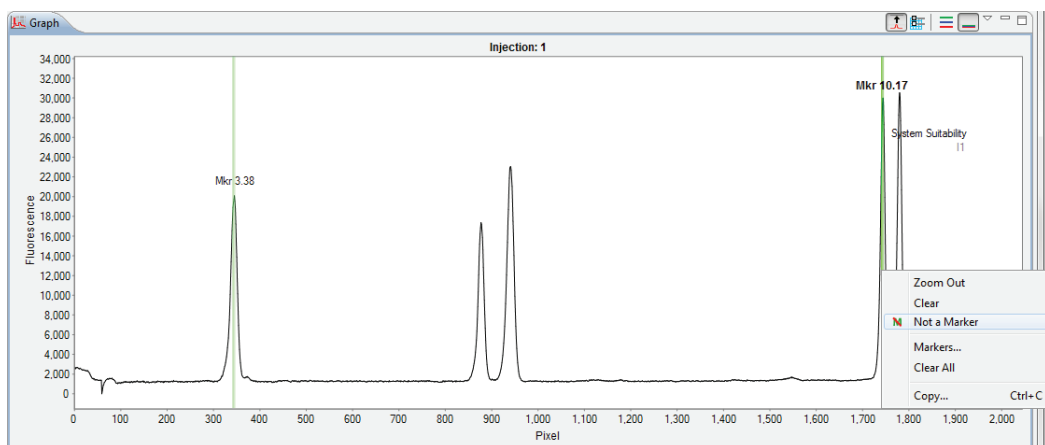
A lock icon indicating the pI marker was set manually will display next to the peak in the Peaks table and a check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Peaks					Injections			
Injection	Sample	Peak	Position	Height				
1	System Suitabi...	14	1315	1315.8				
1	System Suitabi...	15	1422	1433.6				
1	System Suitabi...	16	1549	1627.6				
1	System Suitabi...	17	1743	29396.8				
1	System Suitabi...	18	1780	30540.1				
1	System Suitabi...	19	1959	1399.5				
1	System Suitabi...	20	2018	1470.4				

Experiment			
Injection	Sample	Location	Me
1	System Suitabi...	A1	Sys
2	mAb 11 Blank	A2	mA
3	mAb 11 Ref. St...	A3	mA
4	mAb 11 Prep 2...	A4	mA
5	mAb 11 Prep 2...	A4	mA
6	mAb 11 Prep 2...	A4	mA
7	mAb 11 Ref. St...	A3	mA
8	mAb 11 Blank	A2	mA

NOTE: To remove pI marker peak assignments that were made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**. To clear all the manual settings for the injection, click **Clear All**.

If an incorrect peak is identified as a pI marker: Right-click the peak in the electropherogram or Peaks table and select **Not a Marker**. Compass should correctly reassign the remaining peaks as pI markers and update the Peaks table.



7. Repeat the previous steps for the remaining pI marker peaks as needed in the current injection and for all other injections to make sure all your pI markers are identified correctly.

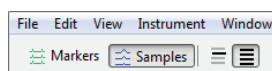
Step 3: Checking Sample Peaks

All detected peaks will be labeled automatically with the calculated protein pI.

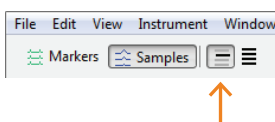
NOTE: The reported protein pI in Compass may vary slightly from predicted pIs based on sample, buffer, and method conditions.

To make sure your sample proteins are identified correctly:

1. Click **Samples** in the View bar.



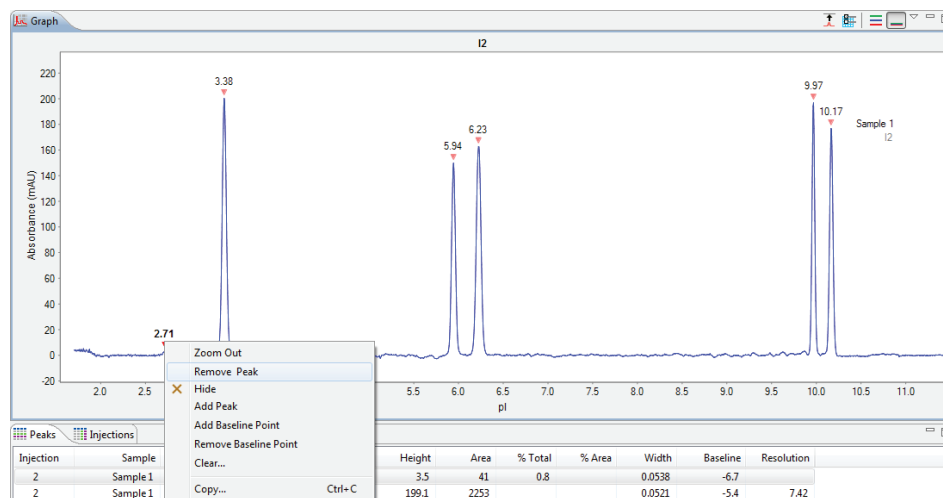
2. Click the **View Selected** icon in the View bar.



3. Click **Injection 1** in the Experiment pane.

4. If your sample peaks aren't identified correctly, here's how to manually correct them:

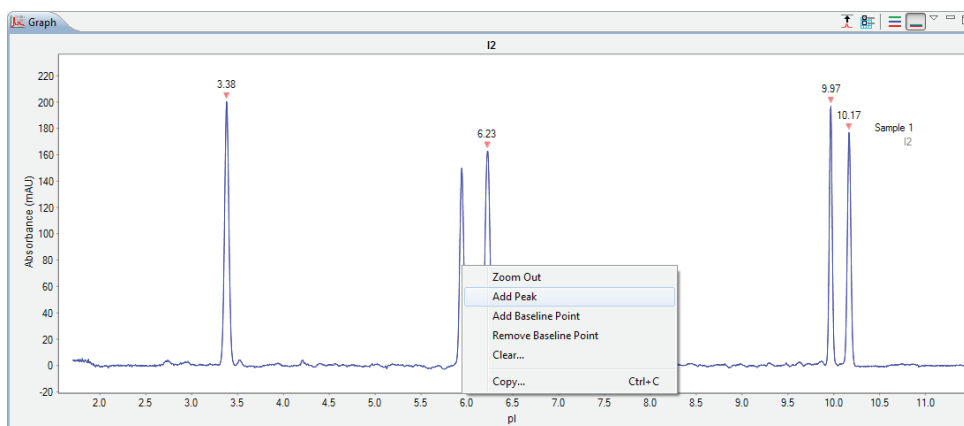
If a peak is incorrectly identified as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass will no longer identify it as a sample peak in the electropherogram and the peak data will be removed in the results table.



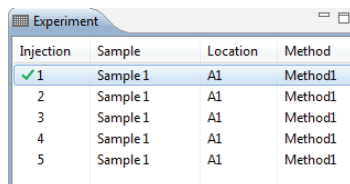
A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Injection	Sample	Location	Method
✓ 1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1

To identify a peak as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.



Injection	Sample	Location	Method
✓ 1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear**.

5. Repeat the previous steps for the remaining injections to make sure all sample peaks are correctly identified.

Step 4: Assigning Peak Names

Compass can also optionally identify and automatically name sample peaks using user-specified peak name settings. For more information on how to do this, see “Manual Peak Integration” on page 474.

Chapter 7:

MauriceFlex cIEF Batches

Chapter Overview

- Overview
- Batch Screen Overview
- Opening a Batch
- Creating a New Batch
- Viewing Replicate Injections
- Batch History
- Making Changes to a Batch
- Viewing and Editing Batches in Completed Runs
- Copying and Pasting Injection Names or Sample IDs from other Documents
- Importing and Exporting Injections
- Importing and Exporting Methods
- Batch Reports

Overview

MauriceFlex cIEF batches can be run on MauriceFlex systems using a cIEF Fractionation cartridge.

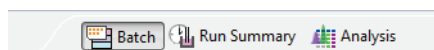
NOTES:

We recommend first optimizing methods with cIEF batches using a cIEF cartridge on Maurice, Maurice C. or MauriceFlex. Then, use recommendations in the MauriceFlex cIEF Fractionation Method Development Guide for tips to modify the method for a MauriceFlex cIEF batch using a cIEF Fractionation cartridge on MauriceFlex to confirm successful method transfer before performing a MauriceFlex Fractionation batch.

On-board mixing is not available on MauriceFlex.

Batch Screen Overview

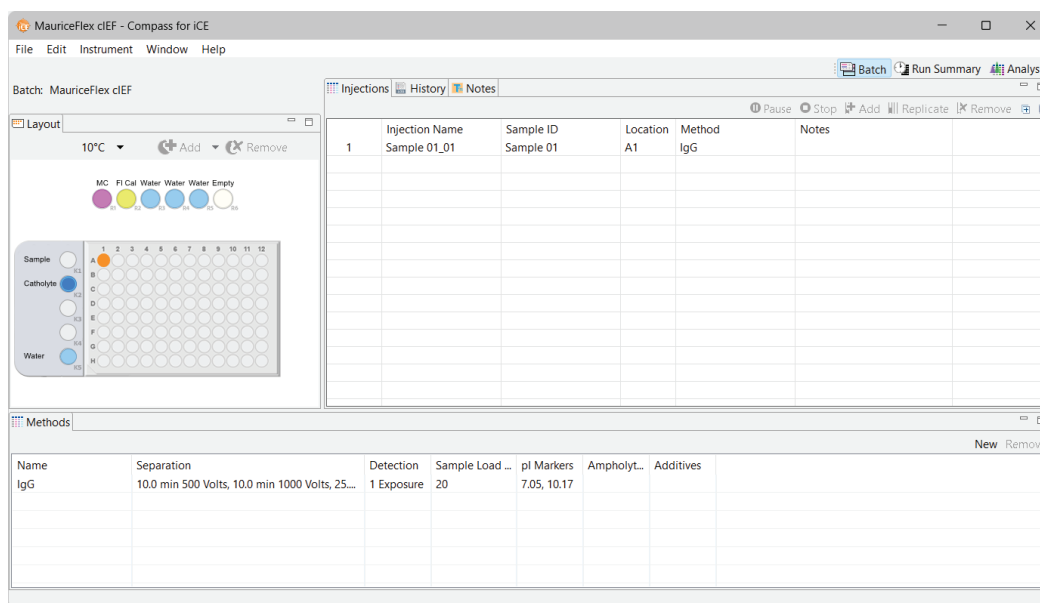
You can use the Batch screen to create, view and edit batches. To get to this screen, click the **Batch** screen tab:



Batch Screen Panes

The Batch screen has five panes:

- **Layout** - Displays a map of the 96-well plate for MauriceFlex cIEF batch sample locations. Batch reagent locations are also displayed.
- **Injections** - Lists the injections, sample ID, sample locations and methods that MauriceFlex will execute for each sample in the batch.
- **History** - Lists all batch file events from initial creation to the most current update.
- **Notes** - Lets you enter specific information about your batch.
- **Methods** - Lets you create methods and enter method parameters used in the batch.



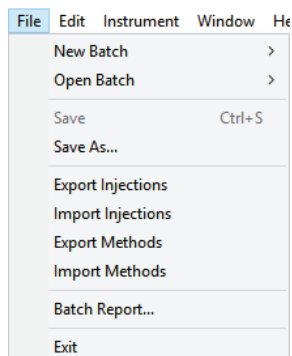
Software Menus Active in the Batch Screen

These main menu items are active in the Batch screen:

- File
- Edit
- Instrument (when the software is connected to MauriceFlex)
- Window
- Help

File Menu

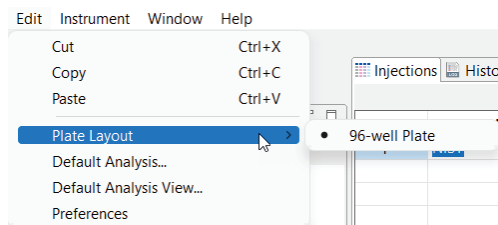
These File menu options are active:



- **New Batch** - Creates a new batch from a starter template.
- **Open Batch** - Opens an existing batch.
- **Save/Save As** - Saves the open batch.
- **Export Injections** - Exports injections from the current batch as a .csv file.
- **Import Injections** - Imports injections into the current batch from a .csv file.
- **Export Methods** - Exports method(s) from the current batch as separate files.
- **Import Methods** - Imports saved method(s) into the current batch.
- **Batch Report** - Exports a table of sample and method details for each injection in the batch as a PDF file.
- **Exit** - Closes Compass for iCE.

Edit Menu

The following Edit menu options are active in the Batch screen:



- **Cut** - Cuts the information currently selected.
- **Copy** - Copies the information currently selected.
- **Paste** - Pastes the copied information.

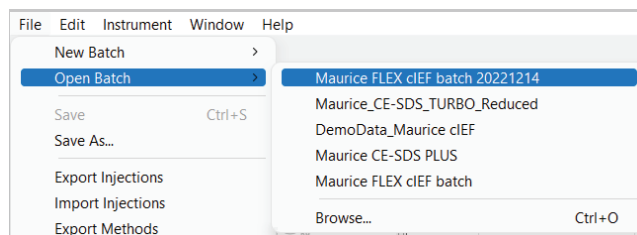
NOTE: Cut, Copy and Paste work differently depending on the pane you're in. How and when you can use them is described throughout the chapter.

- **Plate Layout** - Indicates a 96-well plate will be used to run samples.
- **Default Analysis** - Displays the default settings that will be used to analyze the data generated with your batch.
- **Default Analysis View** - Displays the default settings that will be used to view the data generated with your batch.
- **Preferences** - Lets you set and save your preferences for Access Control, data export, graph colors, grouped data and Twitter settings. See Chapter 21: "Setting Your Preferences" for more information.

Opening a Batch

To open an existing batch:

1. Select **File** in the main menu and click **Open Batch**.



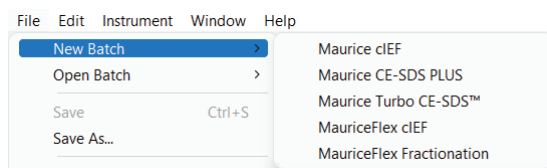
2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file. Select one of those or click **Browse** to open the Batches folder and select a different one.
3. To make changes to the batch, see the steps in “Creating a New Batch” on page 133. When you’re done, select **File** from the main menu and click **Save**.

Creating a New Batch

To create a new batch, we recommend that you use one of the template methods and make any necessary modifications from there.

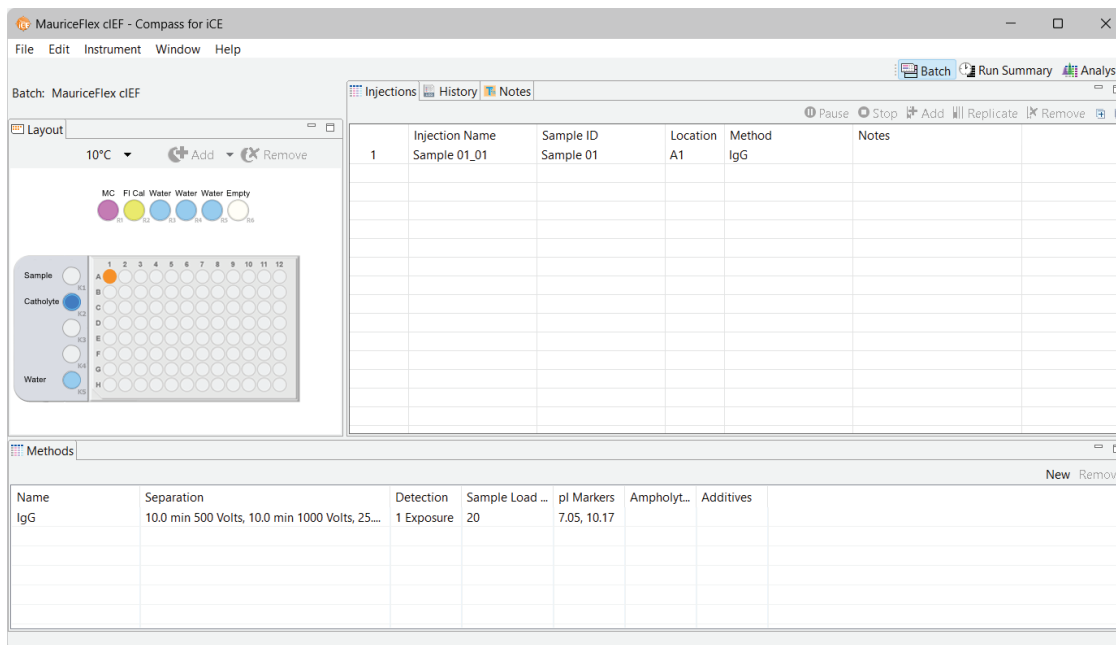
Step 1 - Open a Template Batch

1. Select **File** in the main menu and click **New Batch**:



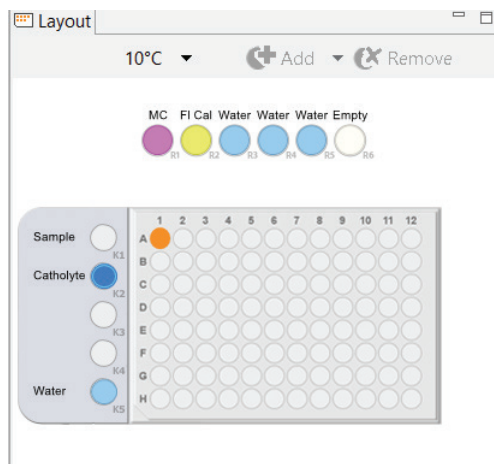
NOTE: Only template batches specific to your Maurice system will be available when you are connected to the instrument. Template batches will not appear in the menu if that application type cannot be performed on your instrument.

2. Select **MauriceFlex cIEF**. A batch using the default method will display.



Step 2 - Assign Your Samples

The Layout pane shows the default locations of where batch reagent should be placed in the fractionation adapter.



The same reagent locations are used for every batch:

- **R1** - 0.5% Methyl Cellulose
- **R2** - Fluorescence Calibration Standard
- **R3** - Water vial

- **R4** - Water vial
- **R5** - Water vial
- **R6** - Empty vial (air)
- **K2** - Catholyte Solution
- **K5** - Water vial

1. Select your samples:

To import samples using a saved injections file:

- Select **File** in the main menu and click **Import Injections**.
- Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.
- Skip to step 2 on page 136.

NOTE: For more information on setting up an injections file for import, see “Importing and Exporting Injections” on page 150.

To select samples manually:

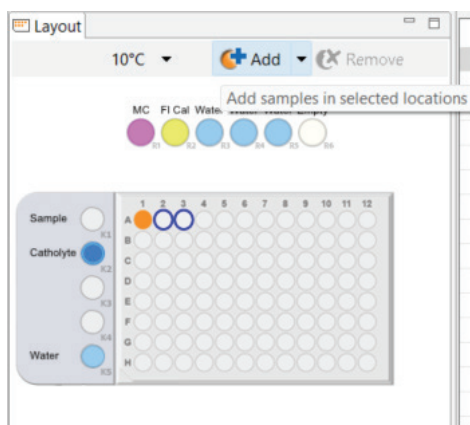
- **Add samples and select methods later:** Use your mouse to highlight the well or vial positions your samples are located in, then click **Add**.

NOTES:

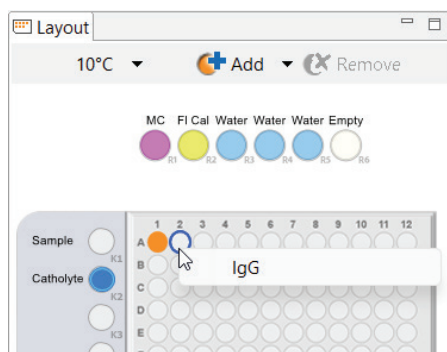
The template batch automatically adds a sample in well A1 by default.

Samples can be assigned to any well in the 96-well plate or in position K1 of the fractionation adapter if the sample is prepared in a vial.

The maximum number of injections per batch for a MauriceFlex cIEF batch is four.



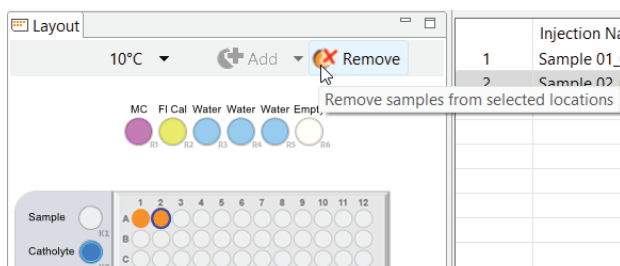
- **Add samples with preassigned methods:** Highlight the well or vial positions your samples are located in, then right-click and select a method.



Either option of adding samples populates the Injections table:

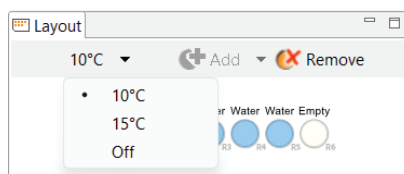
Injections History Notes					
Pause Stop Add Replicate Remove					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 02_02	Sample 02	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	

To remove samples, just highlight the ones you want to remove and click **Remove**. They'll also be removed in the Injections table. You can also right-click well(s) or vial(s) in the Layout pane and click **Remove Sample(s)**.



2. The 96-well plate on the fractionation adapter is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the down arrow next to the temperature and select an option.

NOTE: The reagent row (R) is kept at room temperature while the reagent column (K) and sample plate are temperature controlled.



Step 3 - Assign Your Method Parameters

NOTE: We recommend using the default method parameters. Please see the MauriceFlex cIEF Fractionation Method Development Guide for more information on method optimization.

The Methods pane lists all of the methods used in a batch. You can add and remove methods here and change method parameters.

To import a saved method:

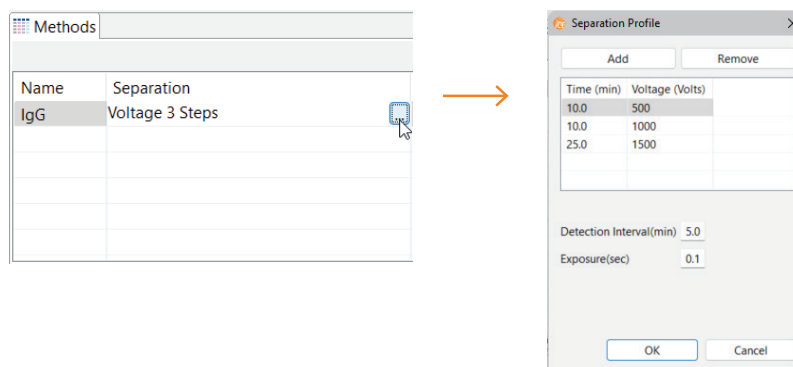
1. Select **File** in the main menu and click **Import Method**.
2. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

To create a new method or make changes to an existing one:

1. In the Methods table, click the first cell in the Name column and enter a new method name if needed.

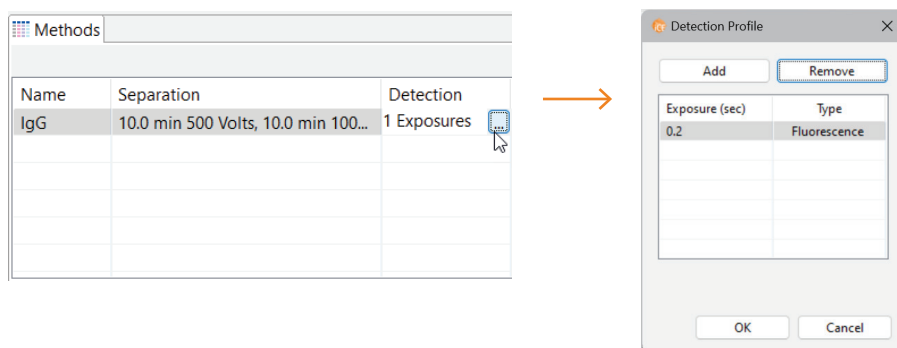
Name	Separation	Detection	Sample Load ...	pl Markers	Ampholyt...	Additives
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25...	1 Exposure	20	7.05, 10.17		

2. Click the first cell in the Separation column, then click the selection button [...] to set the pre-focusing and focusing time (in minutes) and voltage (V). You can also set the Detection Interval (in minutes) and Exposure (in seconds) to adjust how often an image is taken using the listed Exposure time.



- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
- **To change the detection interval or exposure parameters:** Click on the field and type the new value(s).

- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a separation time (in minutes) and voltage value (in V).
 - **To remove a profile step:** Select the row you want to remove and click **Remove**.
3. Click the first cell in the Detection column the selection button [...] to set your fluorescence detection time for the final focused image.



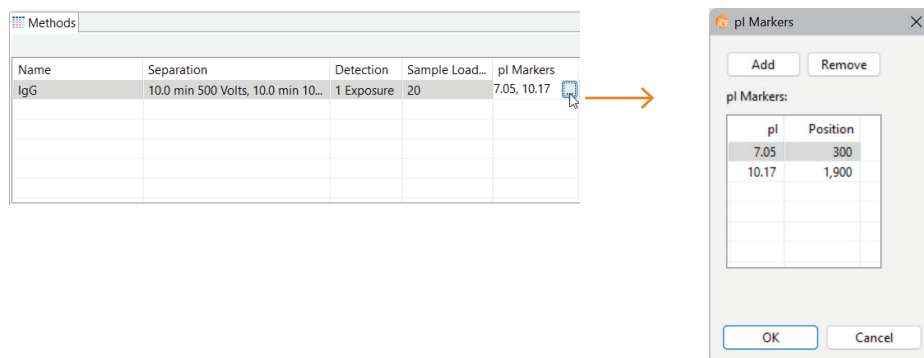
- **To change the exposure time:** Just click in a cell under **Exposure** and type the new value(s) in seconds.
 - **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in an exposure time (in seconds).
 - **To remove a profile step:** Select the row you want to remove and click **Remove**.
4. Click the first cell in the Sample Load(s) column and set the load time in seconds.

NOTE: We recommend using the default Sample Load time of 20 seconds. Please contact ProteinSimple Technical Support if you have questions on the Sample Load time to use for your application.

Methods			
Name	Separation	Detection	Sample Load ...
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25....	1 Exposure	20

5. Click the first cell in the pI Markers column to select pI markers. Add new markers or remove existing ones then click **OK**.

NOTE: When you edit the pI markers in the method for a batch, Compass for iCE automatically creates a Markers group in the pI Markers Analysis settings for you.



- **To change a pI marker and position:** Just click in a cell under **pI** or **Position** and type the new value(s).
 - **To add a pI marker:** Click **Add**. A new row will be added in the table. Then just type in a pI and a position (in pixels).
 - **To remove a pI marker:** Select the row you want to remove and click **Remove**.
6. Optional: Click the first cell in the Ampholytes column and enter the ampholytes you're using.

Methods							
							New Remove
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives	
IgG	10.0 min 500 Volts, 10.0 min 1000 Vol...	1 Exposure	20	7.05, 10.17	Pharmalyte 3-10	Urea	

7. Optional: Click the first cell in the Additives column and enter any additives you're using.

Methods							
							New Remove
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives	
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25.0 ...	1 Exposure	20	7.05, 10.17	Pharmalyte ...	Urea	

8. You can now:

- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.
- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.

Step 4 - Set Up Your Injections

The Injection pane lists all sample injections for the batch. Any samples that were added in “Step 2 - Assign Your Samples” are automatically added to this list.

NOTE: If you imported injections in Step 2, you can update the injections table and assign methods as needed using the information that follows.

Selecting an injection in the table also selects the sample the injection will be made from in the plate or vial map in the Layout pane and vice-versa.

The screenshot shows the MauriceFlex cIEF software interface. On the left is the 'Layout' pane, which includes a temperature control set to 10°C, 'Add' and 'Remove' buttons, and a plate/vial map with wells labeled A, B, C and 1 through 12. On the right is the 'Injections' table, which lists sample injections with columns for Injection Name, Sample ID, Location, Method, and Notes.

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 02_02	Sample 02	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	

1. To add sample names, click the **Sample ID** cell for the injection and type a name.

NOTES:

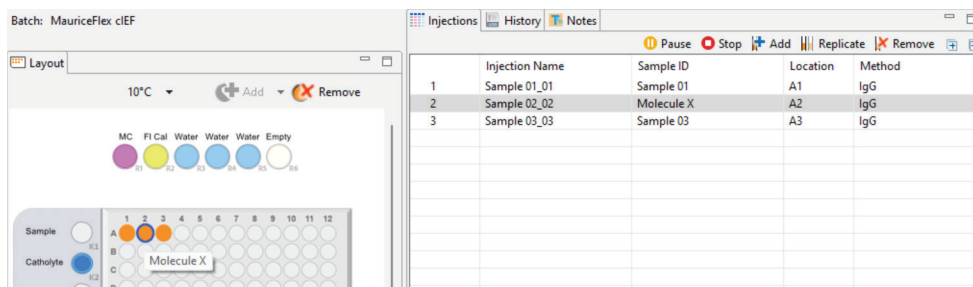
Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 149 “for more info.”

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

This close-up screenshot of the 'Injections' table shows the 'Sample ID' column for the second injection. The original 'Sample 02' has been manually edited to 'Molecule X', and the 'Injection Name' has automatically updated from 'Sample 02_02' to 'Molecule X_02'.

	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	Molecule X_02	Molecule X	A2	IgG
3	Sample 03_03	Sample 03	A3	IgG

The sample name also displays when you hover the mouse over the sample in the plate map:



- Injection names are automatically set to 'Sample ID_injection number'. To change the name, click the **Injection Name** cell for the injection and type a name. Each injection name must be unique. If they aren't when you save the batch, Compass for iCE highlights any non-unique injection names so you can change them.

The screenshot shows the 'Injections' table with the 'Injection Name' cell for row 2 selected. The table data is as follows:

	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	IgG	Sample 02	A2	IgG
3	Sample 03_03	Sample 03	A3	IgG

NOTE: Changing the injection name won't affect the sample ID.

- Assign your methods to each injection. Just click the **Method** cell for the injection and select a method from the drop down menu. If you added your samples with pre-assigned methods and don't need to make any changes, skip to the next step.

The screenshot shows the 'Injections' table with the 'Method' cell for row 3 open, displaying a dropdown menu. The table data is as follows:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	

The dropdown menu for row 3 shows 'Method2' selected.

Hovering over a method name displays the method parameters:

The screenshot shows the 'Injections' table with a tooltip displayed over the 'Method2' cell in row 2. The table data is as follows:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	Method2	
3	Sample 03_03	Sample 03	A3	IgG	

The tooltip for 'Method2' displays the following parameters:

- Separation: 10.0 min 500 Volts, 10.0 min 1000 Volts, 25.0 min 1500 Volts
- Detection: 1 Exposure
- Sample Load (s): 20
- pI Markers: 7.05, 10.17

4. You can add or remove injections as needed using a few different options. If you don't need to make any changes here, just skip to the next step.

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

The first screenshot shows the 'Injections' table with the following data:

	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	Molecule X_02	Molecule X	A2	Method2
3	Sample 03_03	Sample 03	A3	IgG

The second screenshot shows the 'Injections' table after replicating the second row. The data is now:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	Method2	
3	Molecule X_03	Molecule X	A2	Method2	
4	Sample 03_04	Sample 03	A3	IgG	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.

The screenshot shows the 'Injections' table with the following data:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	Method2	
3	Sample 03_03	Sample 03	A3	IgG	

- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

Step 5 - Add Programmed Pauses and Stops (Optional)

You can program pauses and stops in the batch to automatically pause or stop after injections even when you aren't in front of MauriceFlex.

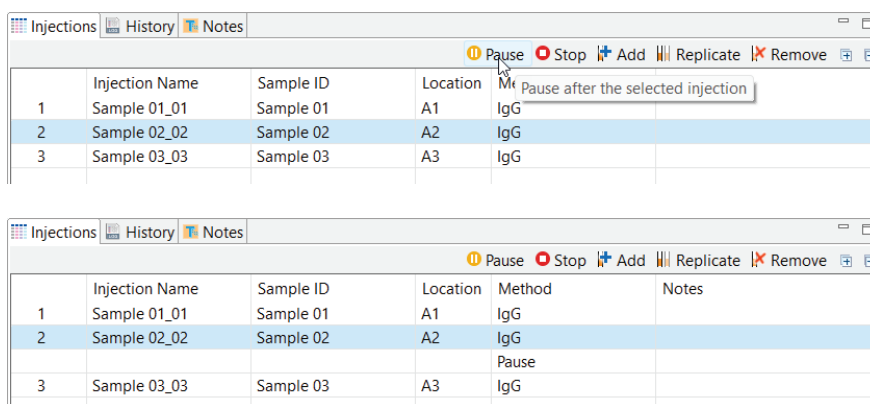
Note: The cartridge cleanup is performed at the end of the batch and after an injection. A programmed pause will occur after the injection before the cartridge cleanup starts.

To program a pause:

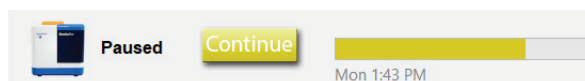
Pauses are helpful when you want to review injection results before continuing with the rest of the batch.

1. Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete.

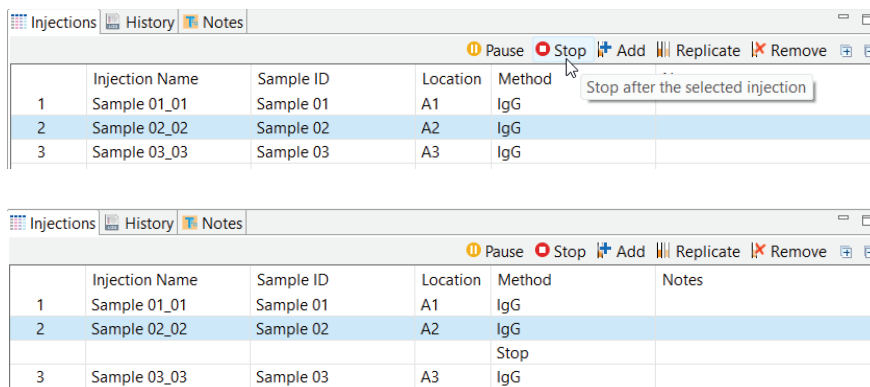
NOTE: MauriceFlex can tweet you when the batch pauses. See “Setting Up Maurice Systems to Send Tweets” on page 764.



2. To resume the batch, click **Continue** in the instrument status bar:

**To stop the run after a specific injection:**

1. Highlight the injection you want the batch to stop at and click **Stop**. MauriceFlex will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.

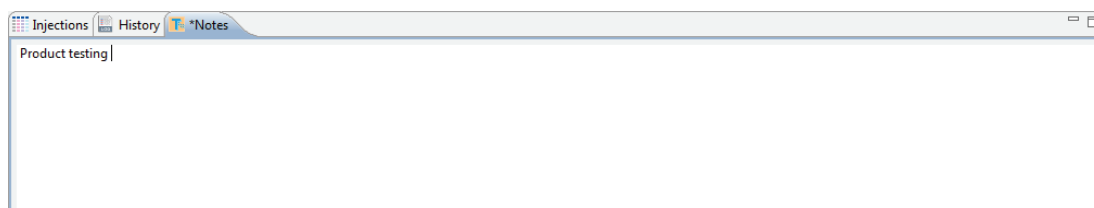


Data collected for injections completed up to the point where the batch was stopped are saved in the .mbz file.

2. Be sure to complete your normal post-batch procedures.

Step 6 - Add Batch Notes (Optional)

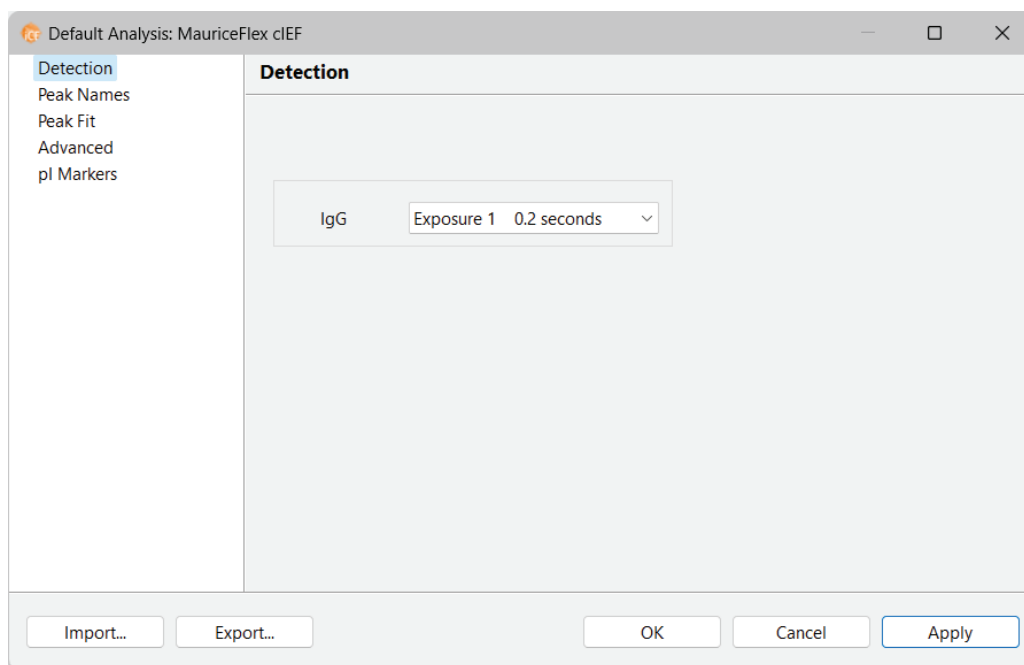
1. Click on the **Notes** pane.
2. Click in the notes area and type any information you want to add about your batch.



Step 7 - Modify Default Analysis Parameters (Optional)

You can preset the parameters used to analyze data generated with the batch. We recommend using the default parameters for MauriceFlex cIEF applications, but if you need to modify parameters:

1. Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



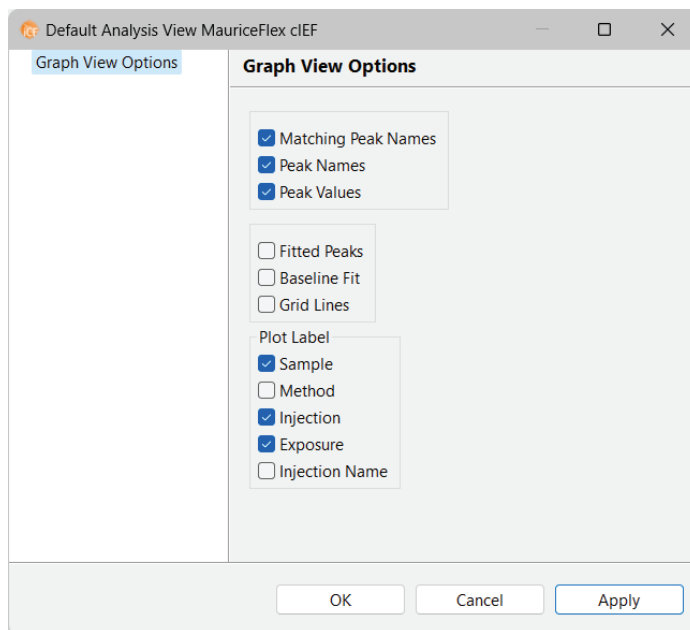
2. Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to "Analysis Settings Overview" on page 528.

Step 8 - Modify Default Analysis View Parameters (Optional)

You can present the parameters used to display the electropherogram data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button.

To modify the parameters:

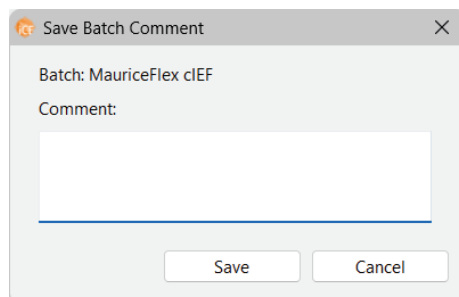
1. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



2. Change the parameters you want to, then click **Apply**. Click **OK** when you are done changing display parameters. For detailed information on graph view options, please refer to “Customizing the Data Display” on page 509.

Step 9 - Save Your Batch

1. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



2. Enter a name for your batch then click **Save**.

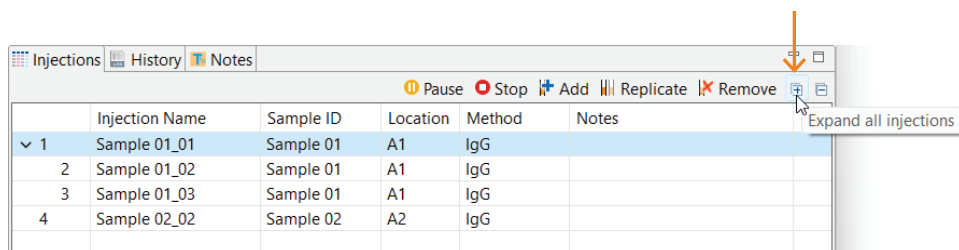
Viewing Replicate Injections

Injections with replicates have an arrow next to the injection number. You can expand or collapse the list of replicate injections by toggling the arrow:

Injections History Notes				
	Injection Name	Sample ID	Location	Method
> 1	Sample 01_01	Sample 01	A1	IgG
4	Sample 02_02	Sample 02	A2	IgG

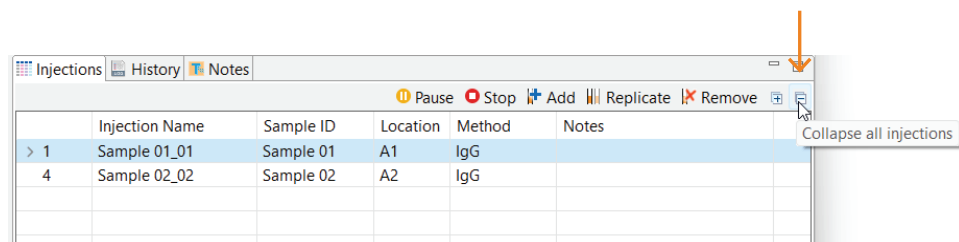
Injections History Notes				
	Injection Name	Sample ID	Location	Method
▼ 1	Sample 01_01	Sample 01	A1	IgG
2	Sample 01_02	Sample 01	A1	IgG
3	Sample 01_03	Sample 01	A1	IgG
4	Sample 02_02	Sample 02	A2	IgG

- To show all replicate injections in the batch, click the **Expand All Injections** button.



Injections History Notes					
	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 01_02	Sample 01	A1	IgG	
3	Sample 01_03	Sample 01	A1	IgG	
4	Sample 02_02	Sample 02	A2	IgG	

- To hide all replicate injections in the batch, click the **Collapse All Injections** button.



Injections History Notes					
	Injection Name	Sample ID	Location	Method	Notes
> 1	Sample 01_01	Sample 01	A1	IgG	
4	Sample 02_02	Sample 02	A2	IgG	

Batch History

Clicking on the History pane shows the batch event history, starting with the date and time the batch was created through the most current event. Clicking on a row in the table displays the full event details in the box under the table.

Date	User Name	Message	Comment
2023-01-13 14:49:08		Started run: 2023-01-13_14-49-17_MauriceFlex cIEF_user manual Batch: MauriceFlex cIEF.batch from Compass for iCE v4.0.0-0110	5221108010
> 2023-01-16 13:12:12		Saved analysis and methods changes from Compass for iCE v4.0.0-0113	

Time	2023-01-13 14:49:08	User
Message	Started run: 2023-01-13_14-49-17_MauriceFlex cIEF_user manual Batch: MauriceFlex cIEF.batch from Compass for iCE v4.0.0-0110	
Comment	5221108010	

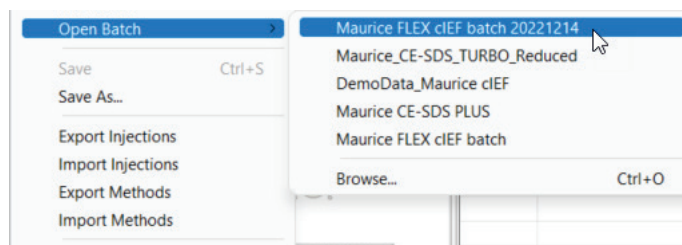
- **Date:** Date and time of the batch event.
- **User Name:** User that initiated the event. User names only display if you're using the Access Control feature to help satisfy the 21CFR Part 11 data security requirements. Go to "Enabling Access Control" on page 781 to learn how to set it up.
- **Message:** Description of the event that took place.
- **Comment:** Comments entered by the user when the batch was saved.

You can copy the information in the History pane to use in other documents:

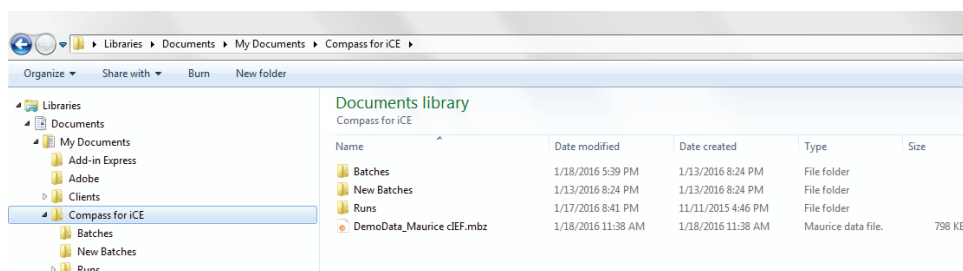
1. Click the **History** pane to make sure it's active.
2. Click **Edit** in the main menu and select **Copy**.
3. Open a document and click **Paste**.

Making Changes to a Batch

1. Select **File** in the main menu and click **Open Batch**.



2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file.
 - Select one of these files or click **Browse** to open the Batch folder and select a different one.
 - To open a batch used in a run file, click **Browse**. Go to the Run folder in the Compass for iCE directory and select the run file that uses the batch you want to edit.

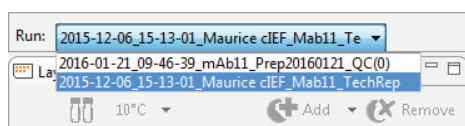


3. To make changes to the batch, see the steps in “Creating a New Batch” on page 133. Then select **File** from the main menu and click **Save** or **Save As**.

Viewing and Editing Batches in Completed Runs

1. Click the **Analysis** screen and open your run file(s).
2. After the run opens, click the **Batch** screen to view the batch used with the run.

If you opened more than one run file, you can switch between viewing batches for each file. Just click the **down arrow** in the Run box and select the batch you want to view from the drop down list.



3. In completed run files, you can only make edits to the **sample** and **method** names in a batch, and these changes are saved to the run file only. Update sample and method names as needed.
4. Click the **Analysis** screen. Then select **File** from the main menu and click **Save** or **Save As** to save the new changes to the run file.

Copying and Pasting Injection Names or Sample IDs from other Documents

Injection names and Sample IDs can be copied from other documents and pasted into the Injection pane to make batch setup faster.

To paste a list of injection names into the batch:

1. Select the list of injection names in a document (Microsoft® Word®, Excel®, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 1	My Sample 2
4		My Injection 1	My Sample 3
5		My Injection 1	My Sample 4

2. Select an injection in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the injection, right click and select **Paste**.

Injections History Notes					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 02_02	Sample 02	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	
4	Sample 04_04	Sample 04	A4	IgG	

The injection names are pasted into the Injection pane:

Injections History Notes					
	Injection Name	Sample ID	Location	Method	Notes
1	My Injection 1	Sample 01	A1	IgG	
2	My Injection 2	Sample 02	A2	IgG	
3	My Injection 3	Sample 03	A3	IgG	
4	My Injection 4	Sample 04	A4	IgG	

To paste a list of Sample IDs into the batch:

1. Select the list of Sample IDs in a document (Microsoft Word, Excel, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 1	My Sample 2
4		My Injection 1	My Sample 3
5		My Injection 1	My Sample 4

2. Select a Sample ID in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the Sample ID, right click and select **Paste**.

Injections History Notes Pause Stop Add Replicate Remove					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 02_02	Sample 02	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	
4	Sample 04_04	Sample 04	A4	IgG	

The Sample IDs are pasted into the Injection pane.

NOTE: If you paste in new sample IDs when the default injection names are still being used, the injection names will automatically update to match the sample ID. If you've already changed the injection names manually, updating the sample IDs will not update the injection names.

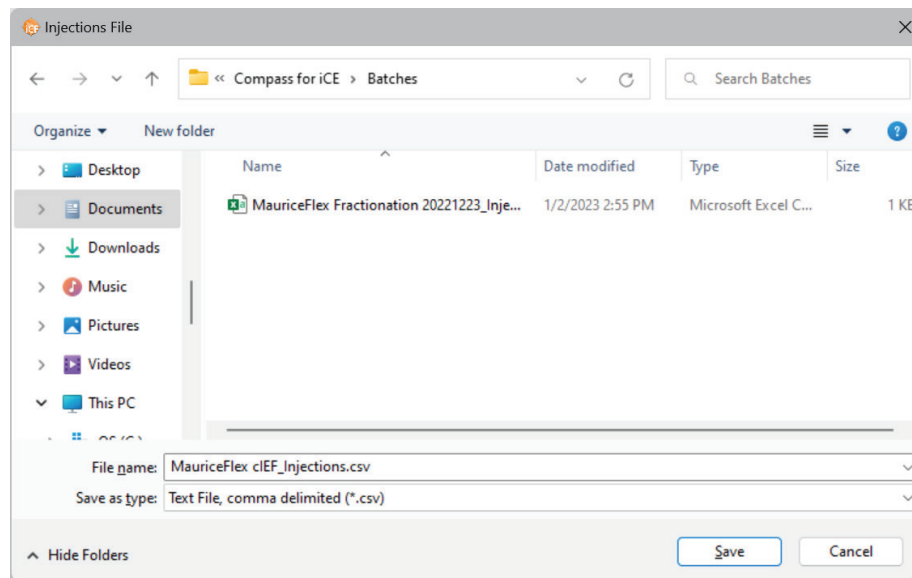
Injections History Notes Pause Stop Add Replicate Remove					
	Injection Name	Sample ID	Location	Method	Notes
1	My Sample 1_01	My Sample 1	A1	IgG	
2	My Sample 2_02	My Sample 2	A2	IgG	
3	My Sample 3_03	My Sample 3	A3	IgG	
4	My Sample 4_04	My Sample 4	A4	IgG	

Importing and Exporting Injections

Injections in an open batch or run file can be exported as a separate file. This lets you import the same or a modified set of injections into another batch later, rather than having to enter the information manually.

Exporting Injections

1. Open the batch or run you want to export injections from.
2. In the Batch screen, select **File** in the main menu and click **Export Injections**. The following window displays:



3. The default directory is Compass for iCE/Batches. Change the directory if needed.
4. The default file name will be batchname.batch_Injections.csv. Change the name if needed and click **Save**.

Creating an Injections File

To create a new injections file .csv, it's easiest to export injections from an existing batch and use that as your template.

1. Follow the instructions in "Exporting Injections" above to export injections from an existing cIEF batch.
2. Open the .csv file in a program that provides a table/spreadsheet format.
3. To add or change injection information:
 - a. Type in a unique name in each cell in the Injection Name column. For replicate injections, type > in front of the injection name, for example: >Injection 1.
 - b. Type in a sample ID, location and method for each injection in their respective columns.
 - c. Optional: Type in notes if needed.

NOTES:

Mix bottles are not used for MauriceFlex cIEF batches. Compass for iCE ignores this column when importing injections into a MauriceFlex cIEF batch.

Programmed pauses and stops can't be added in the injections file. They can be added in the Injections pane after importing your injections into Compass for iCE.

	A	B	C	D	E	F
1	Injection Name	Sample ID	Location	Method	Mix Bottle	Notes
2	Inj01	R01_SSPP	A1	Method_M458		
3	Inj02	S01_Her2_A1_None_0 M Urea	A2	Method_M458		
4	>Inj02	S01_Her2_A1_None_0 M Urea	A2	Method_M458		

4. Save the .csv file.

Importing Injections

1. Open the batch you want to import injections into, or open a new batch.
2. Select **File** in the main menu and click **Import Injections**.
3. Select an injections file (*.csv) and click **OK**. The imported information will display in the Layout and Injections panes.

NOTE: Imported injections will replace any injections already in the Injections table, they will not append an existing table.

Importing and Exporting Methods

Methods in an open batch or run file can be exported as a separate file. This lets you import the same method into another batch later, rather than having to re-enter method information manually.

Importing a Method

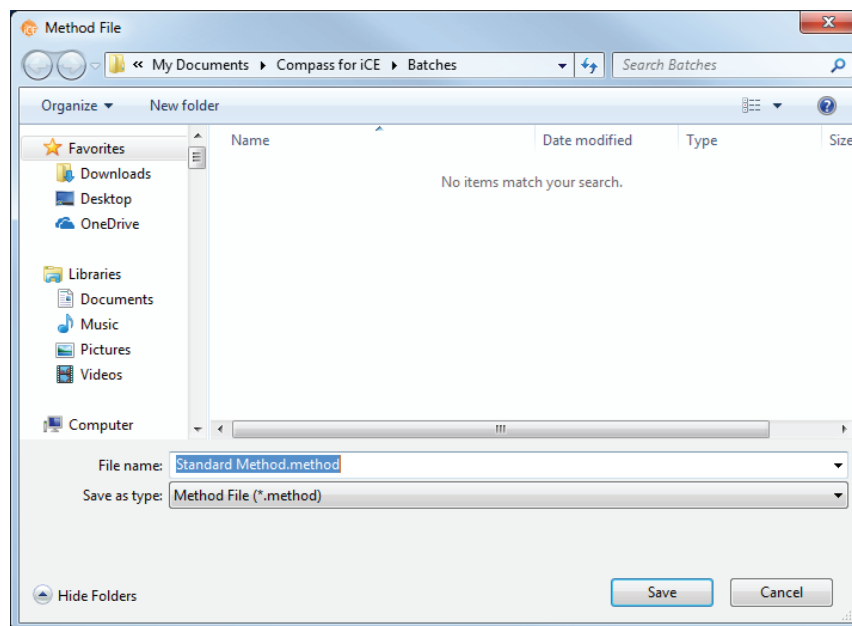
NOTE: Importing a method imports information into the Batch window's Method pane only.

1. Open the batch you want to import the method into.
2. Select **File** in the main menu and click **Import Method**.
3. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

Exporting a Method

NOTE: Exporting a method exports information in the Batch window's Method pane only.

1. Open the batch you want to export the method from.
2. In the Method pane, select the method you want to export. To export more than one method at the same time, hold the Ctrl key to select multiple methods.
3. Select **File** in the main menu and click **Export Method**. The following window displays:



4. The default directory is Compass for iCE/Batches. Change the directory if needed.
5. Enter a method file name and click **Save**. The settings will be saved as a *.method file.

Batch Reports

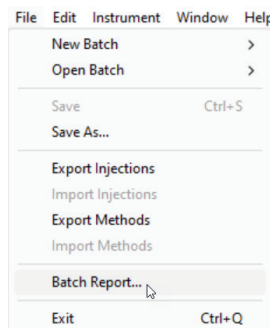
You can export a PDF file of sample and method details for each injection in the batch.

NOTES:

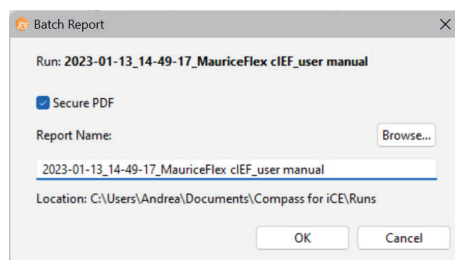
You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” for more information.

Any time a new report is run, it's saved in a separate run folder. Existing reports aren't overwritten.

1. Go to the **Analysis** or **Run Summary** screen, then click **File > Open Run** and select a run file (if you don't have one open already).
2. After the run opens, go to the **Batch** screen.
3. Select **File** from the main menu and click **Batch Report**.

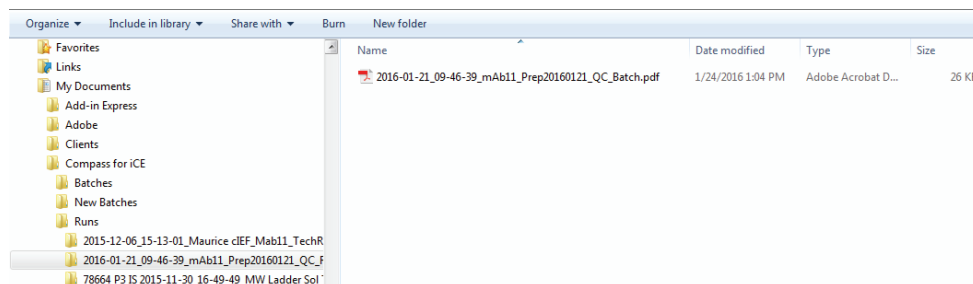


4. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.



Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don't want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.

5. The Batch Report PDF is exported to the **Runs** folder in the **Compass for iCE** directory. It'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Here's an example Batch Report:

cIEF Batch: MauriceFlex cIEF

Injections

Injection	Injection Name	Sample ID	Location	Method	Notes
1	sys suit_01	sys suit	A3	system suit_1	
2	sys suit_02	sys suit	A3	system suit_2	

Methods

Name	Separation	Detection	Sample Load (s)	pl Markers	Ampholytes	Additives
system suit_1	10.0 min, 500 Volts 10.0 min, 1000 Volts 10.0 min, 1500 Volts Detection Exposure: 0.2 sec Interval: 5.0 min	Fluorescence, 0.2 sec	20	3.38 pl, 350 pixels 10.17 pl, 1900 pixels		
system suit_2	10.0 min, 500 Volts 10.0 min, 1000 Volts 10.0 min, 2000 Volts Detection Exposure: 0.2 sec Interval: 5.0 min	Fluorescence, 0.2 sec	20	3.38 pl, 350 pixels 10.17 pl, 1900 pixels		

Batch Log

Date	User Name	Message	Comment
2023-01-13 14:07:35		Batch created using the factory default MauriceFlex cIEF with Compass for ICE Version: 4.0.0-0110	
2023-01-13 14:09:23		Saved as C:\Users\xiaojing.shen\Documents\Compass for ICE\Batches\MauriceFlex cIEF batch from Compass for ICE v4.0.0-0110	

Created By: Andrea Thu 3:56 PM Feb 23, 2023 PST (SECURED)
C:\Users\Andrea\Documents\Compass for ICE\Runs\2023-01-13_14-49-17_MauriceFlex cIEF_user manual.mbz
Computer: DESKTOP-1PM7G05 Software Version: Compass for ICE 4.0.0, Build ID: 0222

Page 1 of 2



cIEF Batch: MauriceFlex cIEF

Date	User Name	Message	Comment
2023-01-13 14:09:23		Save injections and methods changes to C:\Users\xiaojing.shen\Documents\Compass for ICE\Batches\MauriceFlex cIEF batch from Compass for ICE v4.0.0-0110	
2023-01-13 14:49:17		Save injections changes to C:\Users\xiaojing.shen\Documents\Compass for ICE\Batches\MauriceFlex cIEF batch from Compass for ICE v4.0.0-0110	
2023-01-13 14:56:45		Runtime injections changes	
2023-01-13 14:57:55		Runtime methods changes	

Created By: Andrea Thu 3:56 PM Feb 23, 2023 PST (SECURED)
C:\Users\Andrea\Documents\Compass for ICE\Runs\2023-01-13_14-49-17_MauriceFlex cIEF_user manual.mbz
Computer: DESKTOP-1PM7G05 Software Version: Compass for ICE 4.0.0, Build ID: 0222

Page 2 of 2



Chapter 8:

Running MauriceFlex cIEF Applications

Chapter Overview

- Overview
- Before You Throw the Switch
- Power Up
- Running MauriceFlex cIEF Applications
- Editing a Running Batch
- Post-batch Procedures
- Checking Your Data

Overview

MauriceFlex cIEF applications can be run on MauriceFlex system using a cIEF Fractionation cartridge.

NOTES:

We recommend first optimizing methods with cIEF batches using a cIEF cartridge on Maurice, Maurice C. or MauriceFlex. Then, use recommendations in the MauriceFlex cIEF Fractionation Method Development Guide for tips to modify the method for a MauriceFlex cIEF batch using a cIEF Fractionation cartridge on MauriceFlex to confirm successful method transfer before performing a MauriceFlex Fractionation batch.

cIEF Fractionation cartridges support 15 injections.

On-board mixing is not available on MauriceFlex.

Before You Throw the Switch

Ensure that everyone using MauriceFlex have:

- Received instruction in general safety practices for laboratories.
- Received instruction in specific safety practices for MauriceFlex.
- Received instruction on handling of biohazards (if biohazardous materials are to be used on MauriceFlex).
- Read and understood all Product Inserts and related Safety Data Sheets (SDS).

Power Up

1. Turn on the computer connected to MauriceFlex.
2. Turn on MauriceFlex's main power switch.
3. Wait for MauriceFlex to initialize. His indicator light will turn from magenta to blue when he's done.
4. Double-click the Compass for iCE software icon on the computer desktop to open the application.
5. Connect MauriceFlex to Compass for iCE.

Running MauriceFlex cIEF Applications

What You'll Need

- MauriceFlex cIEF Fractionation Cartridges
- MauriceFlex cIEF Fractionation Method Development Kit (optional)

- Maurice cIEF Fluorescence Calibration Standard
- Maurice cIEF pI Markers (3.38, 4.05, 5.85, 6.14, 7.05, 8.4, 9.50, 9.99, or 10.17)
- 0.5% Methyl Cellulose Solution
- 1% Methyl Cellulose Solution
- iCE Electrolyte Kit
- Deionized (DI) water
- MauriceFlex glass crimp top reagent vials, 2 mL



MauriceFlex glass crimp top reagent vial (PN 110-0019, left) and glass reagent vial (PN 046-017, right)
Only MauriceFlex glass crimp top reagent vials (PN 110-0019) should be used to prepare MauriceFlex cIEF batch reagents.

- MauriceFlex glass vials with insert, 0.3 mL for the Fluorescence Calibration Standard
- 96-well plate for samples
- P10, P20, P200, P1000 pipettes and tips
- Electrolyte pipette
- Vortexer
- Table-top centrifuge with plate adapter or vial adapter (12 mm, 2 mL vials)

IMPORTANT:

Use MauriceFlex glass crimp top reagent vials, 2 mL (PN 110-0019) and glass vials with insert, 0.3 mL (PN 110-0018) to prepare MauriceFlex cIEF batch reagents.

Step 1: Prep Your Markers, Samples and Reagents

NOTES:

You can prepare your samples in a 96-well plate or vial.

Samples

1. In a microcentrifuge tube, prepare your sample at a concentration of 2.5 – 10 mg/mL in a final volume of 25 μ L in DI water.
2. In a separate tube, prepare IEF Separation Mix containing your chosen pI marker(s).

NOTES:

Check out the MauriceFlex cIEF Fractionation Method Development Guide for suggested IEF Separation Mix recipes.

If you are running a method optimized for cIEF analysis, increase the concentration of arginine and both pI markers in your IEF Separation Mix. Check out the MauriceFlex cIEF Fractionation Method Development Guide for more information.

3. Add 100 μ L of IEF Separation Mix to the 25 μ L of your sample.
4. Thoroughly vortex the sample to mix completely.
5. Centrifuge the tube at 13,000 $\times g$ for 5 minutes to remove air bubbles and sediment any particulates.
6. Carefully aspirate the top 100 μ L of the sample and pipette it into a well in a 96-well plate by inserting the pipette tip all the way to the bottom to avoid introducing bubbles.
7. Spin the plate for 5 minutes at 1000 $\times g$ using a centrifuge plate adapter.

NOTES:

If you're preparing sample in a sample vial with insert, 0.2 mL (PN 046-083), increase the sample and IEF Separation Mix volumes in steps 1–3 by 3-fold and transfer a minimum of 300 μ L of sample to the vial.

Make sure to check for and remove any bubbles at the bottom of the well.

pI Markers

1. Open the vial of lyophilized pI marker by lifting the center tab and gently pulling it back to break the metal seal. Slowly remove the rubber stopper.
2. Add 210 μ L of DI water to the vial.
3. Put the rubber stopper back in the vial and vortex 3 times, 5 seconds each time to completely reconstitute the lyophilized cake. Repeat this step for all pI markers you're using.

4. Aliquot 20 μ L of each reconstituted pI marker into separate tubes for storage.

NOTES:

Keep your reconstituted pI markers on ice until you're ready to add them to your sample or IEF Separation Mix.

If you'll use the pI markers within a month, store the aliquots at 2-8 °C. Otherwise, store the aliquots at -10 °C to -30 °C. They'll be stable up to 6 months.

5. Use 4 μ L of each pI marker for every 200 μ L of sample.

Reagents

IMPORTANT:

Use MauriceFlex crimp top glass reagent vials, 2 mL (PN 110-0019) and glass vials with insert, 0.3 mL (PN 110-0018) to prepare MauriceFlex cIEF batch reagents. Glass reagent vials, 2 mL (PN 046-017) should only be used when running a standard cIEF or CE-SDS batch.

NOTE: Don't reuse reagents or vials

1. Pipette 350 μ L of Fluorescence Calibration Standard in a MauriceFlex glass vial with insert, 0.3 mL (PN 110-0018) and label.



2. Pipette 2 mL of 0.5% Methyl Cellulose into a MauriceFlex crimp top glass reagent vial and label.



3. Prepare four vials of water by pipetting 2 mL of DI water into a MauriceFlex crimp top glass reagent vial (PN 110-0019). Label the vials.
4. Pipette 2 mL of Catholyte solution to a MauriceFlex crimp top glass reagent vial and label.
5. Label an empty MauriceFlex crimp top glass reagent vial.

NOTES:

Make sure you don't overfill the vials, especially the methyl cellulose vial, to avoid introducing bubbles to your run. Wipe excess liquid at the mouth of the vial with a laboratory wipe.

Vials do not need to be capped.

Step 2: Load Samples and Reagents

1. Open MauriceFlex's door by touching the metal plate on top of the door.



NOTE: The indicator light on MauriceFlex's front panel will blink rapidly as the door disengages.

2. Swing the door open to access the cartridge slot and the reagents and samples platform. The lights on either side of the slot will be **orange**.

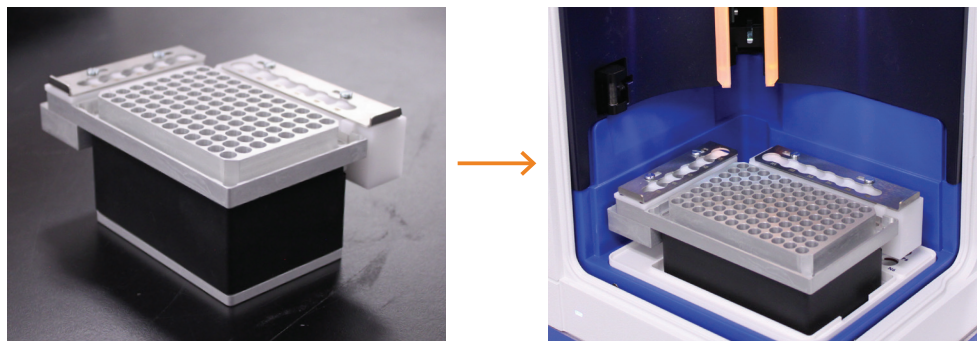


3. Install the fractionation adapter in MauriceFlex.

NOTES:

Remove all the vials in rows P and N of the reagent platform before installing the fractionation adapter if a CE-SDS or standard cIEF application was previously performed.

When installing the fractionation adapter, confirm the adapter is seated and sitting flat in MauriceFlex.



4. Place the reagent vials into their respective positions on the fractionation adapter.

IMPORTANT:

Reagent vials on MauriceFlex must be locked in place before starting a batch. The vial locking mechanism default state is locked. Pull the locking mechanism on Row R to the right and the locking mechanism on Column K down to unlock and place vials in the fractionation adapter. Release the locking mechanism to lock the vials in place.

If the fractionation adapter locking mechanism does not spring back upon release, check that all the vials are seated.

NOTES:

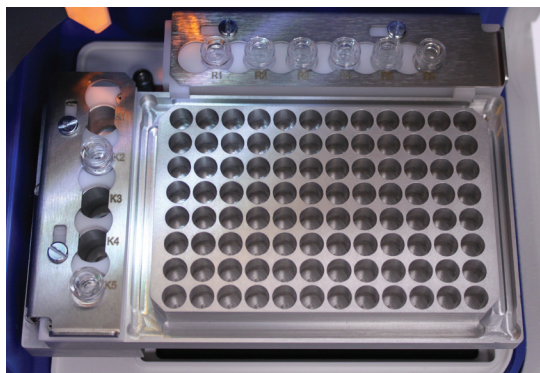
The reagent row (R) on the fractionation adapter is kept at room temperature. The reagent column (K) and sample plate are temperature controlled.

MauriceFlex cIEF batch reagents should be prepared in 2 mL crimp top glass vials (PN 110-0019) and glass vials with insert, 0.3 mL (110-0018).

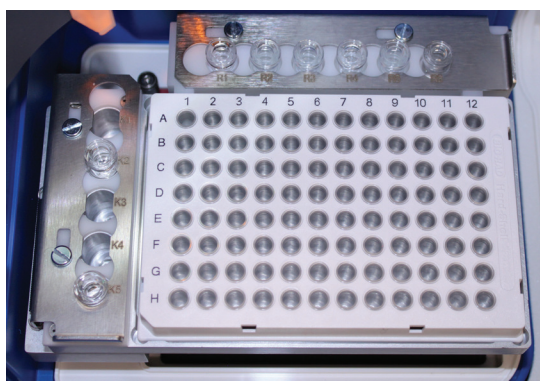
Vials do not need to be capped before placing them on the fractionation adapter.

- **R1** - 0.5% Methyl Cellulose
- **R2** - Fluorescence Calibration Standard
- **R3** - Water
- **R4** - Water

- R5 - Water
- R6 - Empty vial (air)
- K2 - Catholyte Solution
- K5 - Water



5. Place your 96-well sample plate on the Fractionation Adapter. Well A1 should be in the top left corner of the adapter.



Step 3: Prep the cIEF Fractionation Cartridge

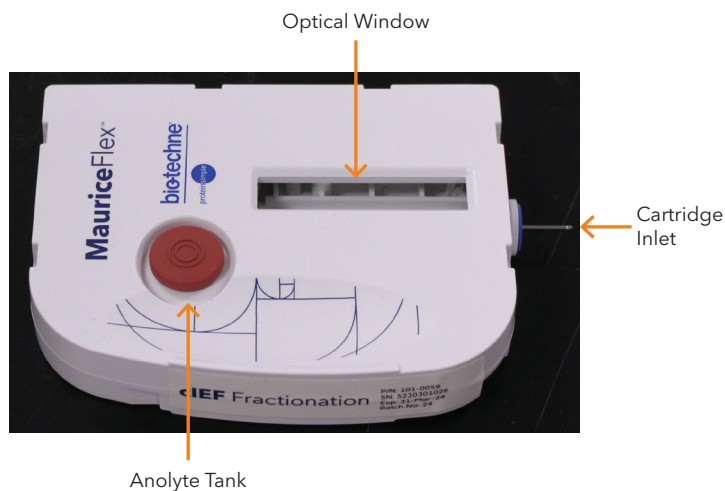
NOTE: A cIEF Fractionation Cartridge supports a maximum of 15 injections and a maximum of 15 batches. Its RFID will keep track of how many are left for you.

1. Take the cIEF Fractionation Cartridge out of its packaging. Save the packaging, you'll need it later to store the cartridge.

NOTE: When you remove the cartridge from its packaging, make sure the cartridge inlet doesn't come in contact with any surface.



2. Put the cartridge on a flat surface with the electrolyte tank facing up.
3. Remove the stopper from the anolyte tank.



4. Add 2 mL Anolyte Solution to the H⁺ electrolyte tank (red port).

NOTE: Make sure you don't overfill the electrolyte tanks.

5. Seal the Anolyte tank with the rubber stopper. If excess liquid comes out of the tank, make sure to wipe it with a laboratory wipe.



Step 4: Install the Cartridge

1. Double check to make sure you've got anolyte loaded and the tank is properly sealed with the stopper.
2. Lift the cartridge and hold it vertically using the finger holds on either side, cartridge inlet down, with the cIEF Fractionation label facing you.
3. Gently insert it into the slot. You'll feel the alignment groove on the cartridge align inside the slot when you insert it.



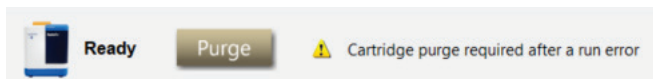
4. Continue to slide the cartridge into the slot until the locking mechanism engages. The lights on either side of the slot will change to **blue** once the cartridge is installed correctly.



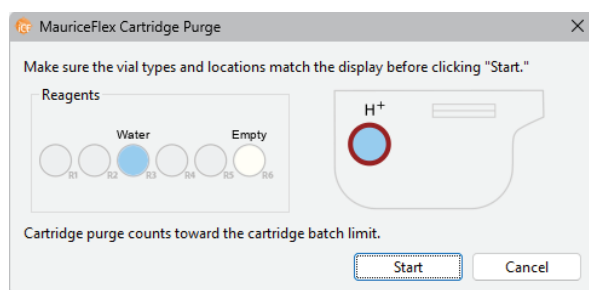
5. Close the instrument door. MauriceFlex locks it automatically.

Step 5: Check for Cartridge Alerts

1. If your cartridge was last used in a run with an error, you will need to perform a Cartridge Purge. MauriceFlex will remind you if a Cartridge Purge is required.
- a. To purge the cartridge, click the brown **Purge** button in the instrument status bar.



- b. Confirm that the required batch reagents are loaded and that the cartridge is prepped. Then click **Start**.



- c. Once the purge is done, perform a Cartridge Post-Run Cleanup by clicking on the brown **Cleanup** button that will appear in the instrument status bar. See step 2 for more information.

NOTE: If there's only one batch left on the cartridge, you'll get a message that it can't be used after purging.

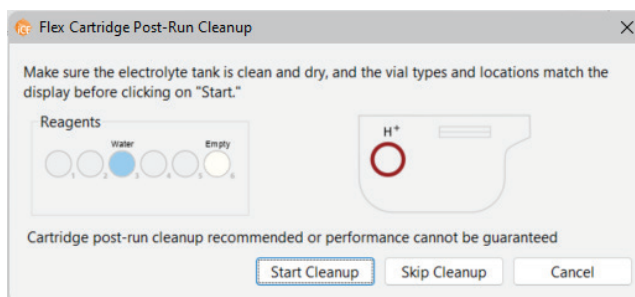
2. If you just completed a Cartridge Purge or if the cartridge you're using did not undergo a cartridge post-run cleanup after the last run, you will be asked if you want to perform a Cartridge Post-Run Cleanup.

To start the Post-Run Cleanup:

- Remove the cartridge from MauriceFlex and remove the analyte. See "Post-batch Procedures" on page 182 for more information.
- Confirm there is a vial of Water (P3) and Air (P6) in place
- Click the brown **Cleanup** button in the instrument status bar.



- Click **Start Cleanup** in the prompt that appears.

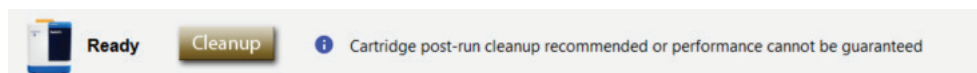


NOTE: You can also start a Post-Run Cleanup by selecting **Instrument > Cartridge Post-Run Cleanup**. When you start a Cartridge Post-Run Cleanup from the instrument menu, click **Start** in the prompt that appears to begin the cleanup.

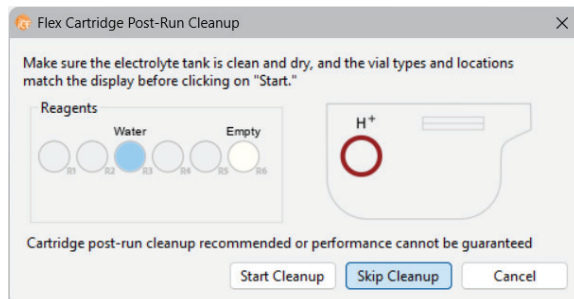
- Once the cartridge post-run cleanup is completed, re-prepare the cartridge and install the cartridge in MauriceFlex. See "Step 3: Prep the cIEF Fractionation Cartridge" on page 163 for more information.
- Click on the green **Start** button in the instrument status menu to start the batch.

To start the run without a Post-Run Cleanup:

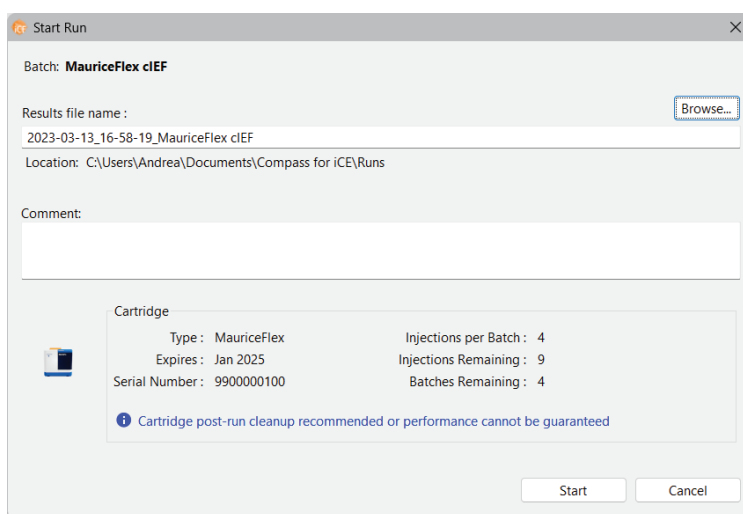
- Click the brown **Cleanup** button in the instrument status bar



- b. Click **Skip Cleanup** in the prompt that appears.



- c. Click **Start** in the Start Run window. The run summary and injection report will record that the cartridge did not undergo a post-run cleanup before starting the batch.



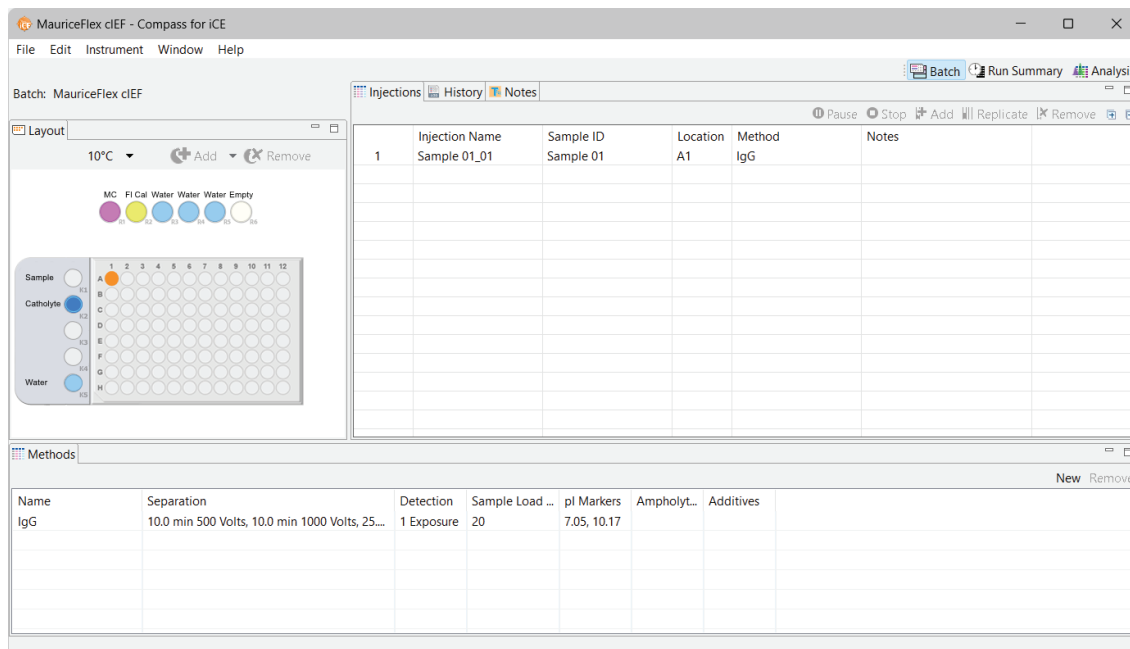
To start the run with a different cartridge:

- If necessary, click **Cancel** in the cIEF Fractionation Cartridge Post-Run Cleanup window.
- Open MauriceFlex's door, remove the first cartridge from MauriceFlex and prepare a second cartridge. See "Step 3: Prep the cIEF Fractionation Cartridge" on page 163 for more information.

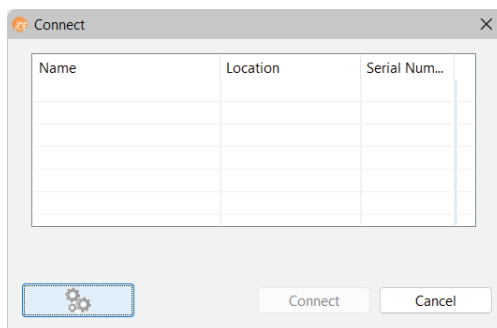
Step 6: Create a Batch

- Launch Compass for iCE.

2. Select the **Batch** tab. This is where you'll enter sample/injection information, methods and batch parameters.

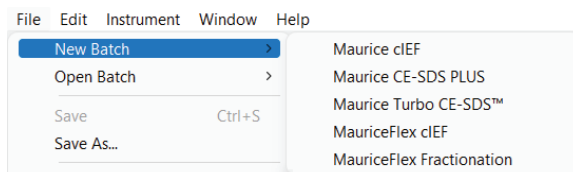


3. To create a batch, make sure MauriceFlex is connected to Compass for iCE. Select **Instrument** and click **Connect**.
 - a. If your instrument is listed, select your MauriceFlex system and click **Connect**.
 - b. If your instrument isn't listed, click on the **Settings** button and connect by typing in your instrument IP address.



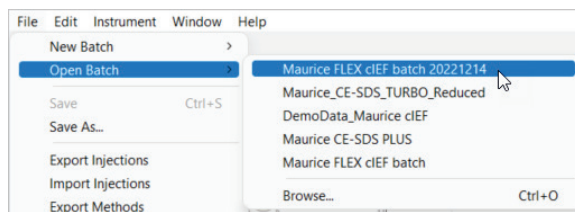
To create a new batch:

- In the main menu, select **File > New Batch > MauriceFlex cIEF**



To use an existing batch: In the main menu, select **File > Open Batch**.

NOTE: If you're making changes to an existing batch, follow the steps in this section. Otherwise skip to "Step 7: Start the Batch" on page 178.



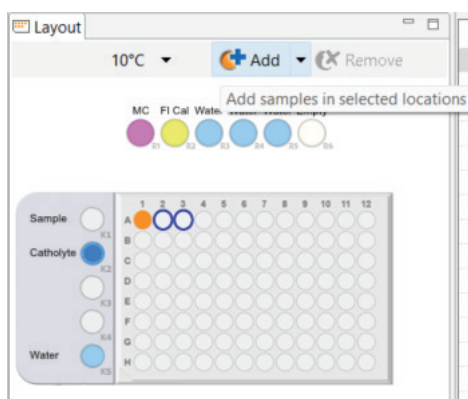
4. Add your samples:

To import samples using a saved injections file:

- a. Select **File** in the main menu and click **Import Injections**.
- b. Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.

NOTE: For more information on setting up an injections file for import, see "Importing and Exporting Injections" on page 150.

To select samples manually: Use your mouse to highlight the well positions for each of your samples, then click **Add**.

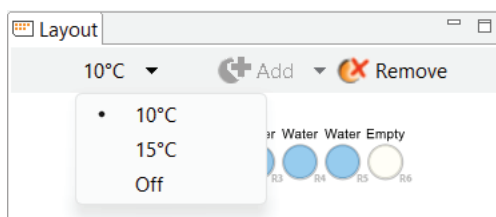


This populates the Injections table:

Injections					
History Notes Pause Stop Add Replicate Remove					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 02_02	Sample 02	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	

- The fractionation adapter is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The reagent row (R) is kept at room temperature. The reagent column (K) and sample plate are temperature controlled.



- Enter your method parameters in the Methods pane.

NOTE: We recommend using the default method parameters. Please see the MauriceFlex cIEF Fractionation Method Development Guide for more information on method optimization.

To import a saved method:

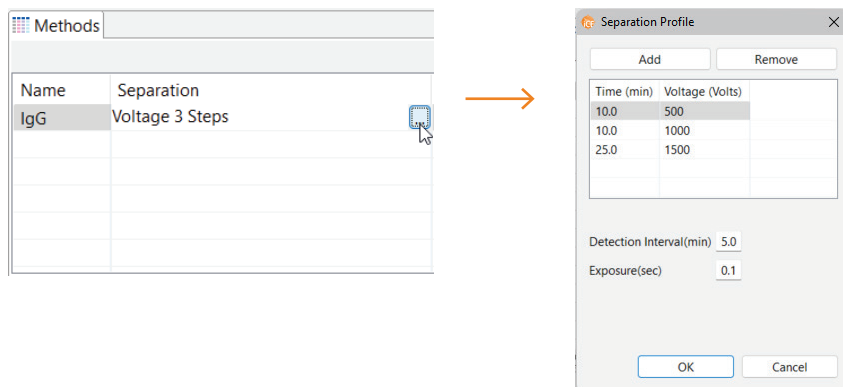
- Select **File** in the main menu and click **Import Method**.
- Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

To create a new method or make changes to an existing one:

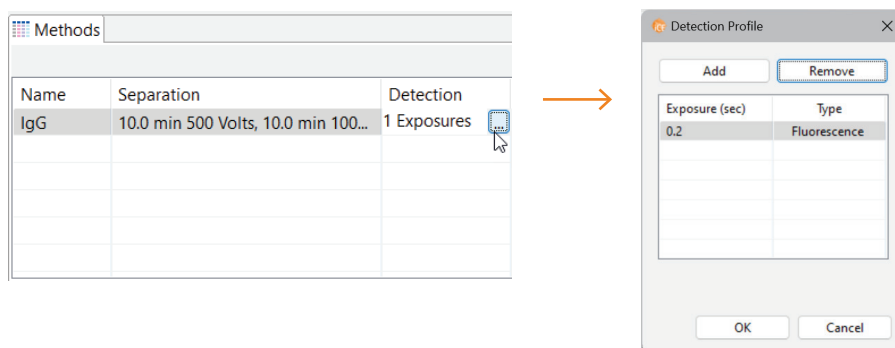
- Click the first cell in the Name column and enter a method name.

Methods						
New Remove						
Name	Separation	Detection	Sample Load ...	pI Markers	Ampholyt...	Additives
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25....	1 Exposure	20	7.05, 10.17		

- b. Click the first cell in the Separation column, then click the selection button [...] to set the pre-focusing and focusing time (in minutes) and voltage (V). You can also set the Detection Interval (in minutes) and Exposure (in seconds) to adjust how often an image is taken using the listed Exposure time.



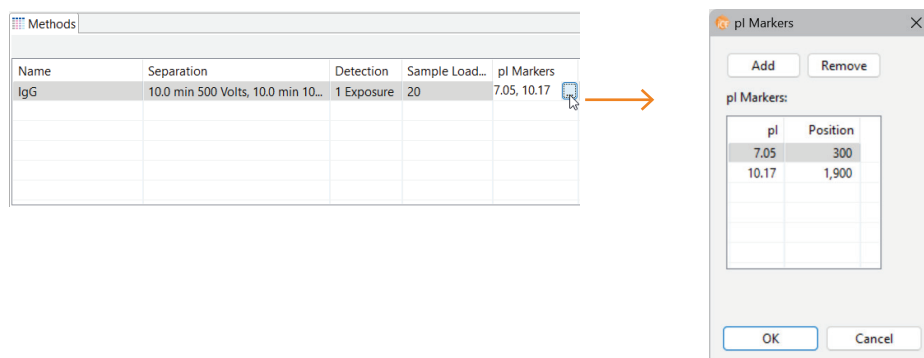
- c. Click the first cell in the Detection column the selection button [...] to set your fluoresences detection time for the final focused image.



- d. Click the first cell in the Sample Load(s) column and set the load time in seconds.

Methods			
Name	Separation	Detection	Sample Load ...
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25...	1 Exposure	20

- e. Click the first cell in the pI Markers column to select pI markers. Add new markers or remove existing ones then click **OK**.



- f. Optional: Click the first cell in the Ampholytes column and enter the ampholytes you're using.

Methods							
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives	
IgG	10.0 min 500 Volts, 10.0 min 1000 Vol...	1 Exposure	20	7.05, 10.17	Pharmalyte 3-10	Urea	

- g. Optional: Click the first cell in the Additives column and enter any additives you're using.

Methods							
Name	Separation	Detection	Sample Load (s)	pI Markers	Ampholytes	Additives	
IgG	10.0 min 500 Volts, 10.0 min 1000 Volts, 25.0 ...	1 Exposure	20	7.05, 10.17	Pharmalyte ...	Urea	

7. You can now:

- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.

8. In the Injections pane:

- **To add or change sample names:** Click the **Sample ID** cell for the injection and type a name.

NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 149 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	Sample 02_02	Molecule X	A2	IgG
3	Sample 03_03	Sample 03	A3	IgG

- **To change injection names:** Click the Injection Name cell for the injection and type a name.

NOTES:

Each injection name must be unique.

Changing the injection name won’t affect the sample ID.

	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	IgG	Sample 02	A2	IgG
3	Sample 03_03	Sample 03	A3	IgG

- **To assign methods for each injection:** Click the Method cell for the injection and select a method from the drop down menu.

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

The first screenshot shows the 'Injections' table with the following data:

	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	Molecule X_02	Molecule X	A2	Method2
3	Sample 03_03	Sample 03	A3	IgG

The second screenshot shows the 'Injections' table after replicating the second row:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	Method2	
3	Molecule X_03	Molecule X	A2	Method2	
4	Sample 03_04	Sample 03	A3	IgG	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.

The screenshot shows the 'Injections' table with the following data:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Molecule X_02	Molecule X	A2	Method2	
3	Sample 03_03	Sample 03	A3	IgG	

- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

9. Add programmed pauses or stops in the batch to automatically pause or stop after injections (optional).

- **To program a pause:** Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete. To resume the batch, click **Continue** in the instrument status bar.

The first screenshot shows the 'Injections' table with the following data:

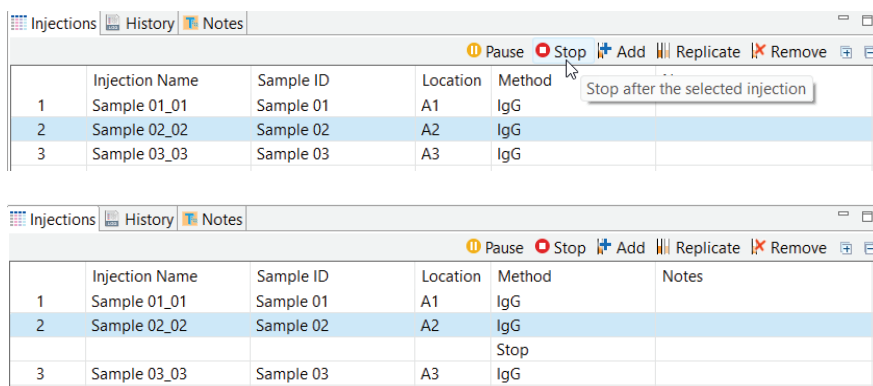
	Injection Name	Sample ID	Location	Method
1	Sample 01_01	Sample 01	A1	IgG
2	Sample 02_02	Sample 02	A2	IgG
3	Sample 03_03	Sample 03	A3	IgG

The second screenshot shows the 'Injections' table after pausing at the second row:

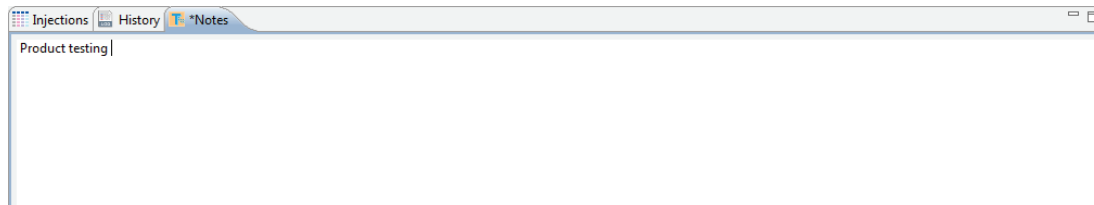
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
2	Sample 02_02	Sample 02	A2	IgG	
3	Sample 03_03	Sample 03	A3	IgG	

NOTE: A cleanup step will occur after each injection on the MauriceFlex cartridge. A programmed pause will occur after the injection completes but before the cleanup step begins.

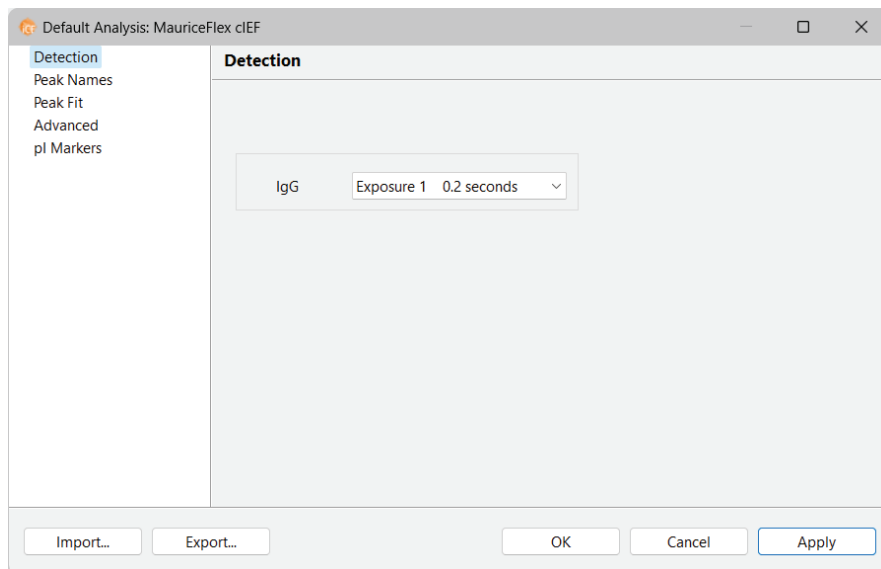
- **To stop the run after a specific injection:** Highlight the injection you want the batch to stop at and click **Stop**. MauriceFlex will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.



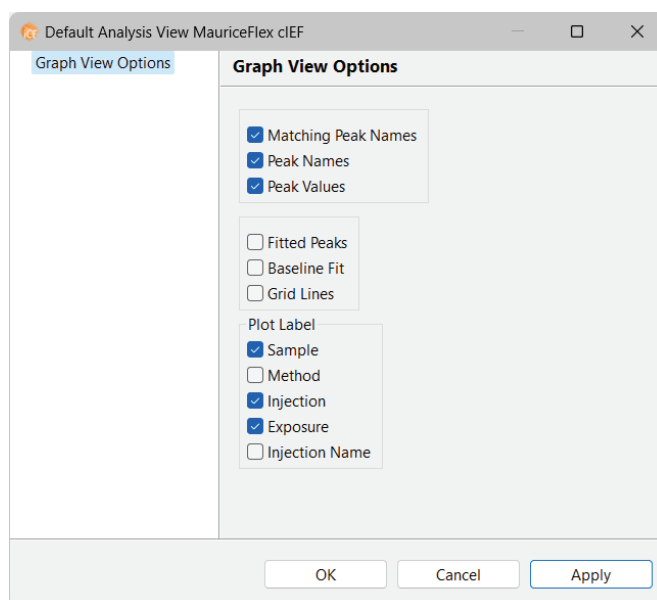
10. Click on the **Notes** pane, then click in the notes area and type any information you want to add about your batch (optional).



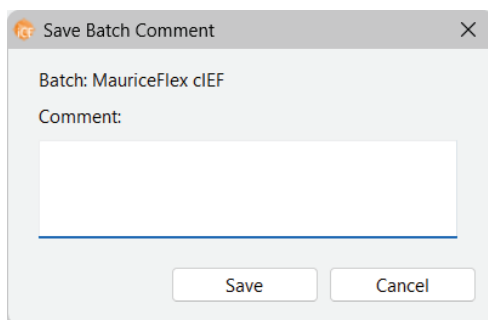
11. You can preset the parameters used to analyze data generated with the batch (optional). We recommend using the default parameters for MauriceFlex cIEF applications, but if you want to modify parameters:
- Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



- Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to "Analysis Settings Overview" on page 528.
12. You can preset the parameters used to view data when the batch is complete (optional). Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button:
- Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



- b. Change the parameters you want to, then click OK. For detailed information on graph view options, please refer to “Customizing the Data Display” on page 509.
13. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



14. Enter a name for your batch then click **Save**.

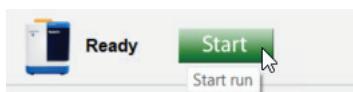
Step 7: Start the Batch

1. Make sure MauriceFlex is connected to Compass for iCE. If he isn't, select **Instrument** and click **Connect**. Then select your MauriceFlex system and click **Connect**.
2. Click on the green **Start** button to start your batch.

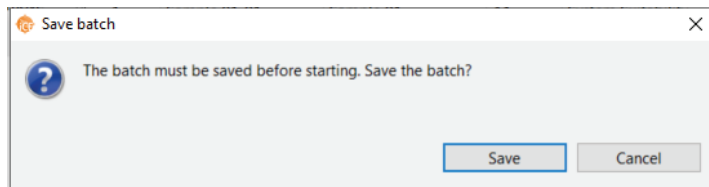
NOTES:

An alert may appear if the wrong cartridge is installed or if a cartridge post-run cleanup or cartridge purge is recommended. See “Step 5: Check for Cartridge Alerts” for more information.

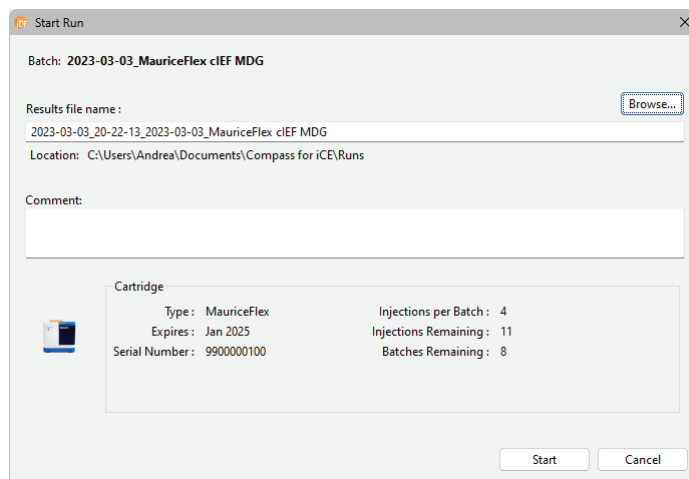
An alert may appear if you are starting a MauriceFlex cIEF batch and have not installed the fractionation adapter. See “Adapter and Insert Alerts” on page 403 for more information.



3. If you have made any changes, you will be asked to save your batch before starting the run. Click **Save**.



- The Start Run box shows you the number of injections and batches remaining on your cartridge. Make sure you still have enough injections left before you start the run. If not, prep a new cartridge.
- Click in the **Results File name** box if you want to change the default run file name. Otherwise just leave it as is.



Start Run

Batch: 2023-03-03_MauriceFlex cIEF MDG

Results file name: Browse...

Location: C:\Users\Andrea\Documents\Compass for ICE\Runs

Comment:

Cartridge

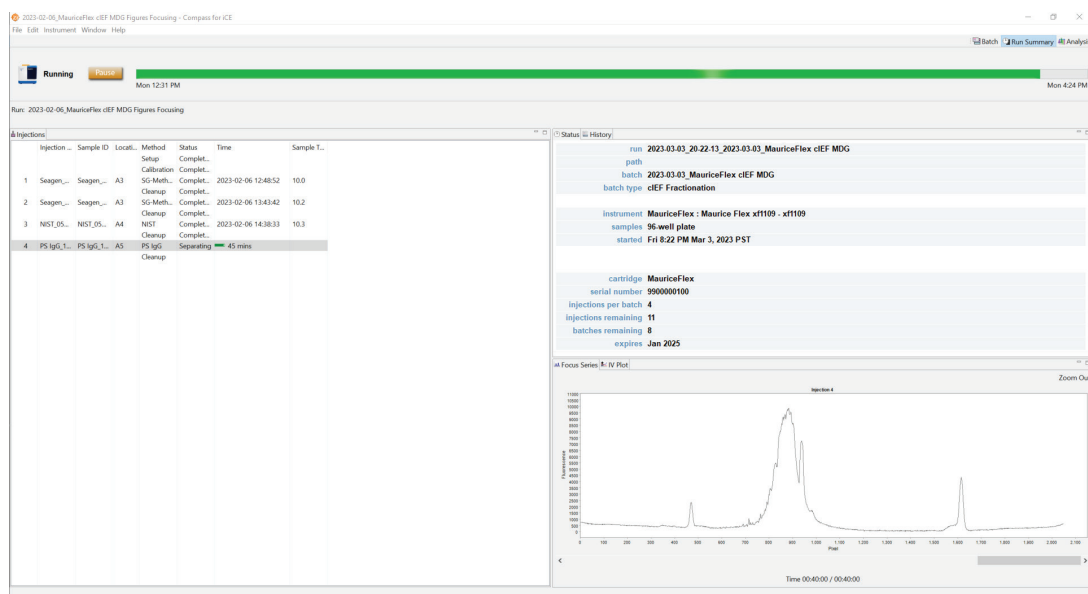
Type: MauriceFlex	Injections per Batch: 4
Expires: Jan 2025	Injections Remaining: 11
Serial Number: 9900000100	Batches Remaining: 8

Start Cancel

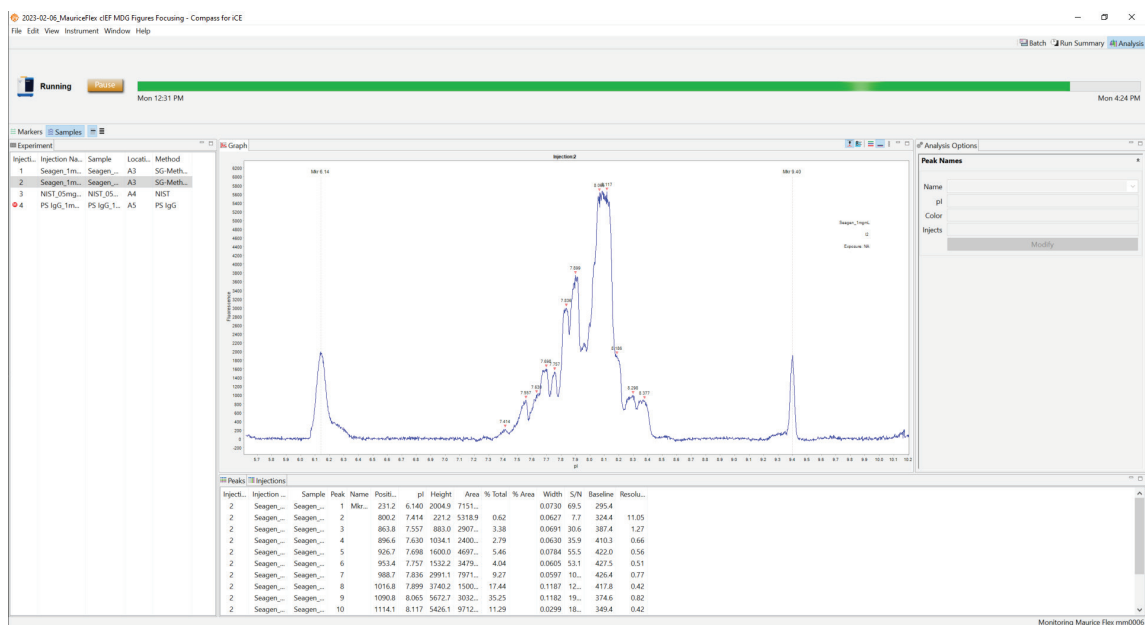
- If you don't want to save the file to the default Runs folder, click **Browse** to select a different location.
- Enter any run details you'd like in the Comments box (optional).
- Click **Start** to start the run.

NOTE: The indicator light on MauriceFlex's front panel will pulse slowly while he's running the batch.

To monitor the progress of your batch, click the **Run Summary** tab. See Chapter 16: "Run Status" for more details.



To view results, analyze results or change analysis parameters on completed injections while the batch is still running, click the **Analysis** tab. See Chapter 18: “cIEF Data Analysis” for more details.



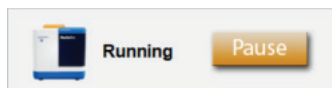
When your batch is complete, you can view electropherograms for all injections in the Analysis tab, and all batch and injection details in the Run Summary tab.

NOTE: If you'd like MauriceFlex to let you know when your run is done, you can set him up to tweet you. Go to “Setting Up Maurice Systems to Send Tweets” on page 764 for more info.

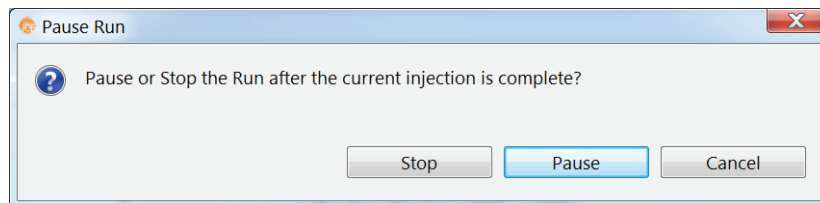
Editing a Running Batch

You can pause a batch after it's started to make updates or add samples and reagents.

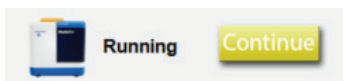
1. Select **Instrument > Pause** or select the **Pause** button.



2. Click **Pause** in the pop-up window.



The Pause button will change to Continue but MauriceFlex's status will stay at Running until Compass for iCE completes the current injection.



- You can now make these edits to the batch:
 - Add sample(s)
 - Add injection(s) to the end of the batch
 - Modify injections that haven't completed and aren't greyed out
 - Create a new method
 - Update the method. Once a method is run it can only be updated. Highlight the method you want to edit and click **Update** in the Methods pane. The software will copy the parameters into a new method that you can modify.
3. When MauriceFlex completes the current injection, his status will change to Paused and the progress bar will turn yellow.

NOTE: MauriceFlex's status light will change from a pulsing blue to a blue pulse with magenta flashes which means he's paused the run.



The status in the Injection pane also changes to Paused and the pause duration updates live. You can now open the instrument door to:

- Add new samples

IMPORTANT

The door must be closed within 15 minutes of opening it to avoid potential assay performance issues.

NOTE: If the door is open for more than 30 seconds, a white light near the cartridge will start to flash and a warning message displays in Compass for iCE until the door is closed.

4. To continue the batch, click **Continue**. Changes you made to the batch are saved and MauriceFlex will continue with the next sample injection.

Post-batch Procedures

When the batch is done:

1. Open MauriceFlex's door. The lights on either side of the cartridge slot will be **orange** as MauriceFlex will have already disengaged the cartridge.
2. Remove your sample. Leave the Water (P3) and Air (P6) vials in place if your cartridge still has injections left since they will be needed for the post-run cleanup step. Remove the remaining reagent vials.

NOTE: You can remove reagent vials from the MauriceFlex fractionation adapter when the locking mechanism is pulled to the right (row R) or pulled down (column K). Release the mechanism to lock remaining vials in place before starting the Post-Run Cleanup.

3. Gently pull out the cartridge making sure the cartridge inlet doesn't come into contact with any surface.



If you're at 15 injections, you've reached the limit of supported injections for the cartridge. To dispose of a finished cartridge, put it in its original packing and discard it per your institution's safety and waste disposal guidelines. Discard any remaining reagent vials.

NOTE: Disposal of cartridges, sample plates and vials depends on the samples you ran. If you aren't sure what your sample origins are, we recommend you dispose of everything in biohazard waste.

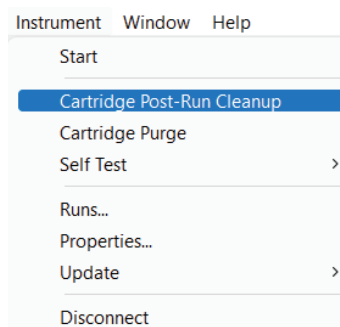
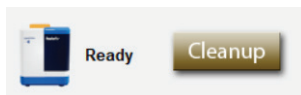
If the cartridge will be used again, clean up and store the cartridge:

- a. Put the cartridge on a flat surface with its electrolyte tank facing up.
- b. Remove the stoppers from the Anolyte tank.
- c. Using an electrolyte pipette or low vacuum, aspirate the solution from the tank.
- d. Fill the tank with 3 mL DI water, then aspirate it out. Repeat this rinse 2 more times.

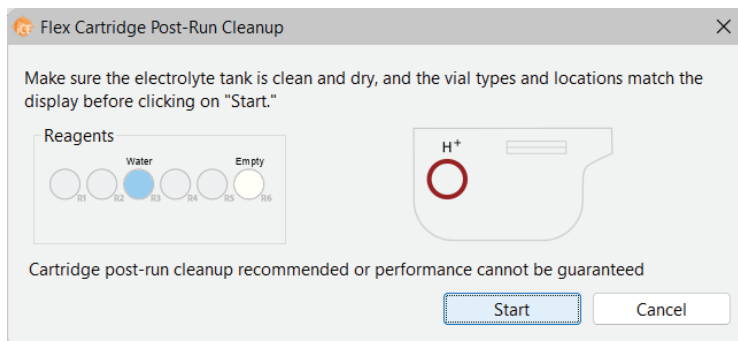
NOTE: Make sure not to get any liquid on the cartridge's optical window.



- e. Aspirate all the remaining liquid and make sure that the tank is dry.
- f. Put the stopper back on the tank and install the cartridge in MauriceFlex.
- g. Verify there is at least 2 mL of DI water in the Water (P3) vial and air in the empty (P6) vial.
- h. Click on the brown **Cleanup** button in the Instrument Status Bar or select **Instrument** and **Cartridge Post-Run Cleanup** in the Compass main menu.



- i. You'll get the following message. Click **Start**. It'll only take 6 minutes.



- j. Open MauriceFlex's door. Gently pull out the cartridge making sure the cartridge inlet doesn't come in contact with any surface.
- k. Remove the reagent vials.
- l. Leave the stopper off to allow the tank to air dry.
- m. Put the cartridge and stopper back in the protective packaging and store at room temperature.

!WARNING! SHARPS HAZARD

The capillary inlet of the cartridge may present a potential sharps hazard. Dispose of used cartridges according to your organizations health and safety regulations.



!WARNING! BIOHAZARD

Cartridges, sample plates and vials should be handled by procedures recommended in the CDC/NIH manual: Biosafety in Microbiological and Biomedical Laboratories (BMBL). The manual is available from the U.S. Government Printing Office or online at <http://www.cdc.gov/biosafety/publications/bmbl5/>.

Depending on the samples used, the cartridges, plates and vials may constitute a chemical or a biohazard. Dispose of the cartridges, plates and vials in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations. Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste content before you store, handle, or dispose of chemical waste.

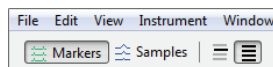
Checking Your Data

Compass for iCE detects your sample protein and pI marker peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

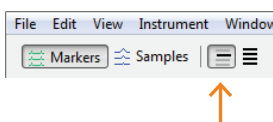
Step 1: Check Your pI Markers

To make sure your pI markers are identified correctly:

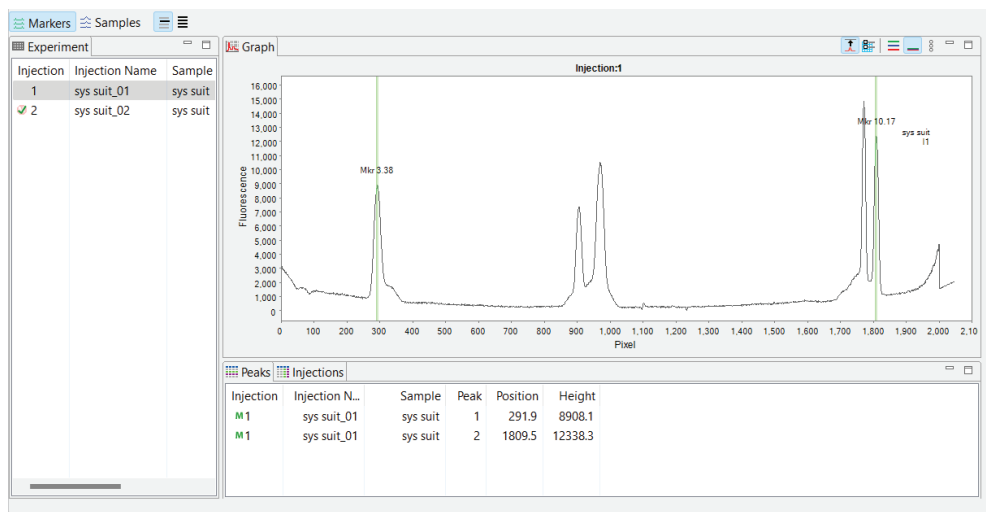
1. Go to the **Analysis** screen.
2. Click **Markers** in the View bar.



3. Click the **View Selected** icon in the View bar.

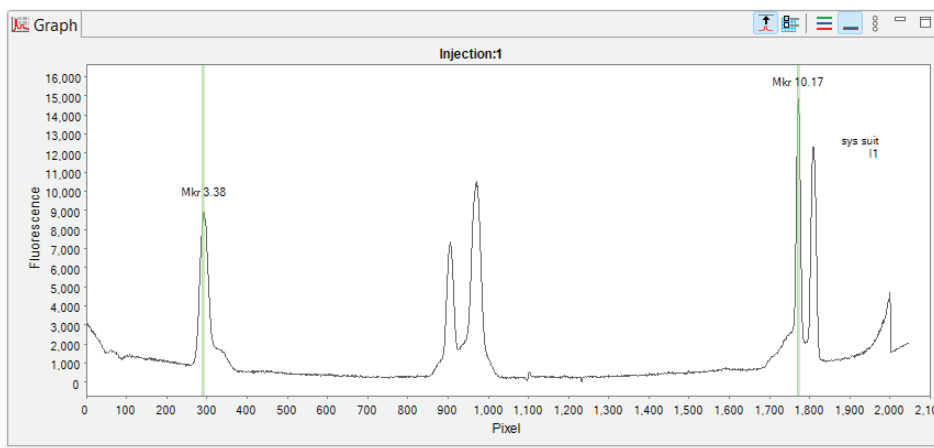
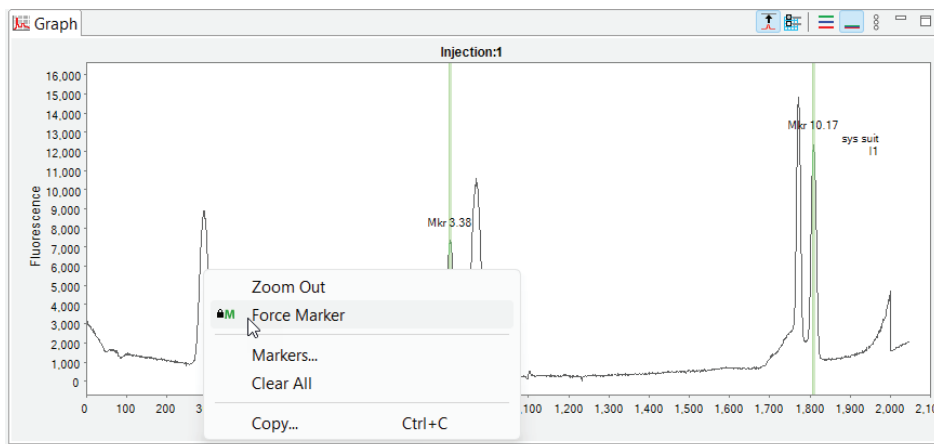


4. Click **Injection 1** in the Experiment pane.
5. Check that your pI markers in the electropherogram have been correctly identified. Each marker peak will have a green vertical line running through it and be labeled Mkr. They're also identified with an **M** in the Peaks table.



6. If your pI markers aren't identified correctly, here's how to manually correct them:

To set an unidentified peak as a pI marker: Right-click the peak in the electropherogram or Peaks table and select **Force Marker**. Compass will assign that peak as a pI marker, and correctly reassign the remaining pI marker peaks.



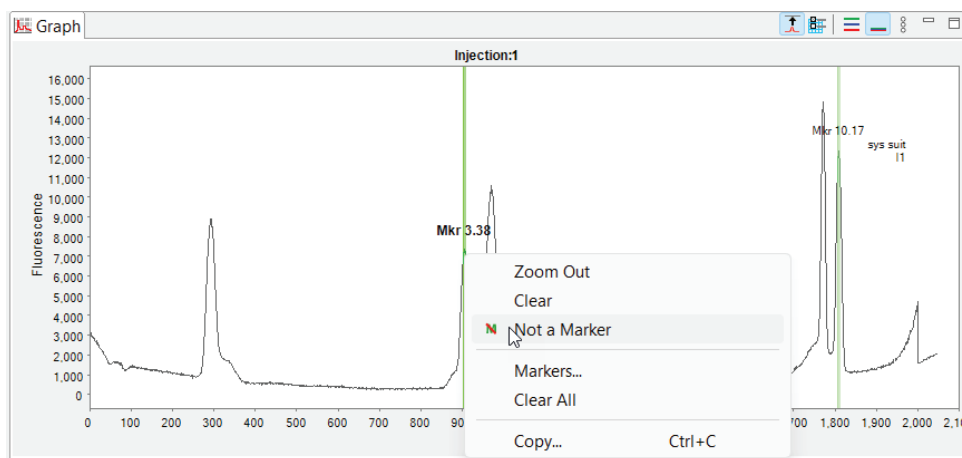
A lock icon indicating the pI marker was set manually will display next to the peak in the Peaks table and a check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Peaks		
Injection	Injection Name	Sample
M 2	sys suit_02	sys suit
🔒 M 2	sys suit_02	sys suit

Experiment		
Injection	Injection Name	Sample
1	sys suit_01	sys suit
✓ 2	sys suit_02	sys suit

NOTE: To remove pI marker peak assignments that were made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**. To clear all the manual settings for the injection, click **Clear All**.

If an incorrect peak is identified as a pI marker: Right-click the peak in the electropherogram or Peaks table and select **Not a Marker**. Compass should correctly reassign the remaining peaks as pI markers and update the Peaks table.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment		
Injection	Injection Name	Sample
✓ 1	sys suit_01	sys suit
2	sys suit_02	sys suit

- Repeat the previous steps for the remaining pI marker peaks as needed in the current injection and for all other injections to make sure all your pI markers are identified correctly.

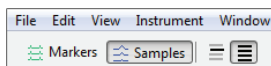
Step 2: Checking Sample Peaks

All detected peaks will be labeled automatically with the calculated protein pI.

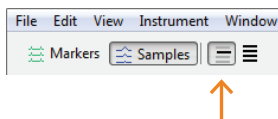
NOTE: The reported protein pI in Compass may vary slightly from predicted pIs based on sample, buffer, and method conditions.

To make sure your sample proteins are identified correctly:

- Click **Samples** in the View bar.



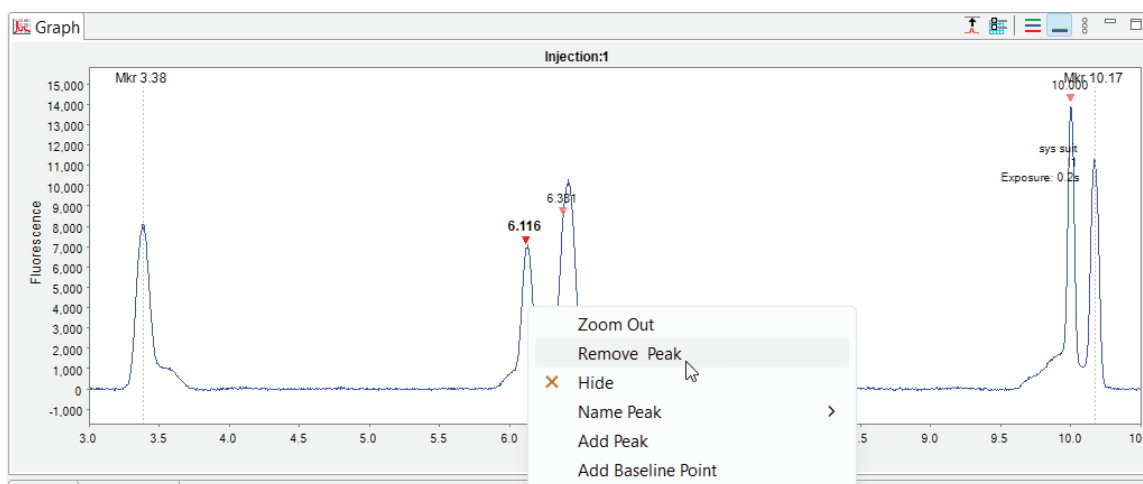
2. Click the **View Selected** icon in the View bar.



3. Click **Injection 1** in the Experiment pane.

4. If your sample peaks aren't identified correctly, here's how to manually correct them:

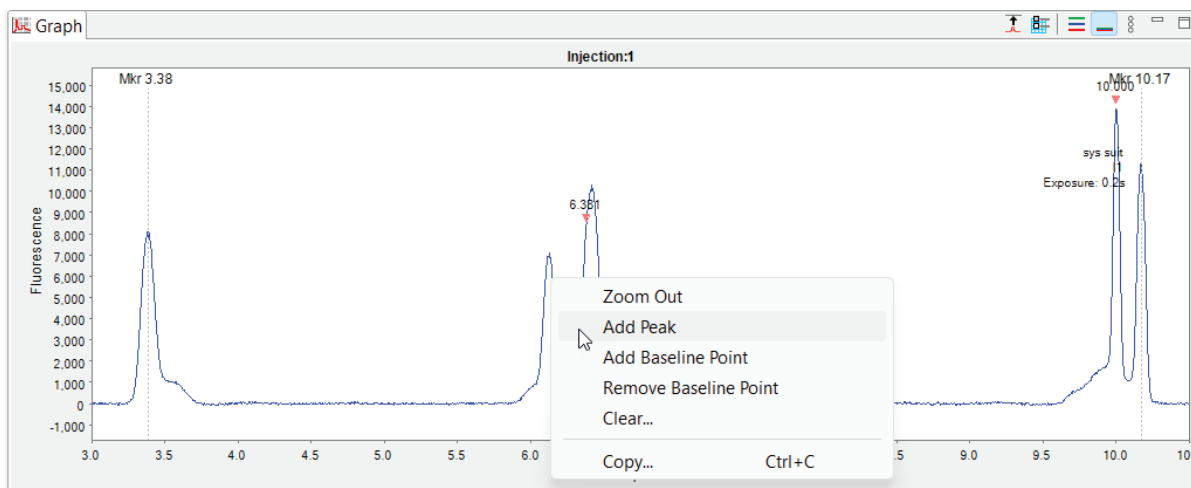
If a peak is incorrectly identified as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass will no longer identify it as a sample peak in the electropherogram and the peak data will be removed in the results table.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment		
Injection	Injection Name	San
✓ 1	sys suit_01	sys
2	sys suit_02	sys

To identify a peak as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment		
Injection	Injection Name	San
✓ 1	sys suit_01	sys
2	sys suit_02	sys

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear**.

- Repeat the previous steps for the remaining injections to make sure all sample peaks are correctly identified.

Step 3: Assigning Peak Names

Compass can also optionally identify and automatically name sample peaks using user-specified peak name settings. For more information on how to do this, see “Manual Peak Integration” on page 547.

Chapter 9:

MauriceFlex Fractionation Batches

Chapter Overview

- Overview
- Batch Screen Overview
- Opening a Batch
- Creating a New Batch
- Viewing Fractions To Be Collected
- Batch History
- Making Changes to a Batch
- Viewing and Editing Batches in Completed Runs
- Copying and Pasting an Injection Name or Sample ID from other Documents
- Importing and Exporting an Injection
- Importing and Exporting Methods
- Batch Reports

Overview

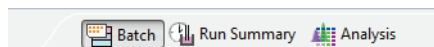
MauriceFlex Fractionation batches can be run on MauriceFlex systems using a cIEF Fractionation cartridge.

NOTES:

We recommend first optimizing methods with cIEF batches using a cIEF cartridge on Maurice, Maurice C. or MauriceFlex. Then, use recommendations in the MauriceFlex cIEF Fractionation Method Development Guide to modify the method for a MauriceFlex cIEF batch using a cIEF Fractionation cartridge on MauriceFlex to confirm successful method transfer before performing a MauriceFlex Fractionation batch.

Batch Screen Overview

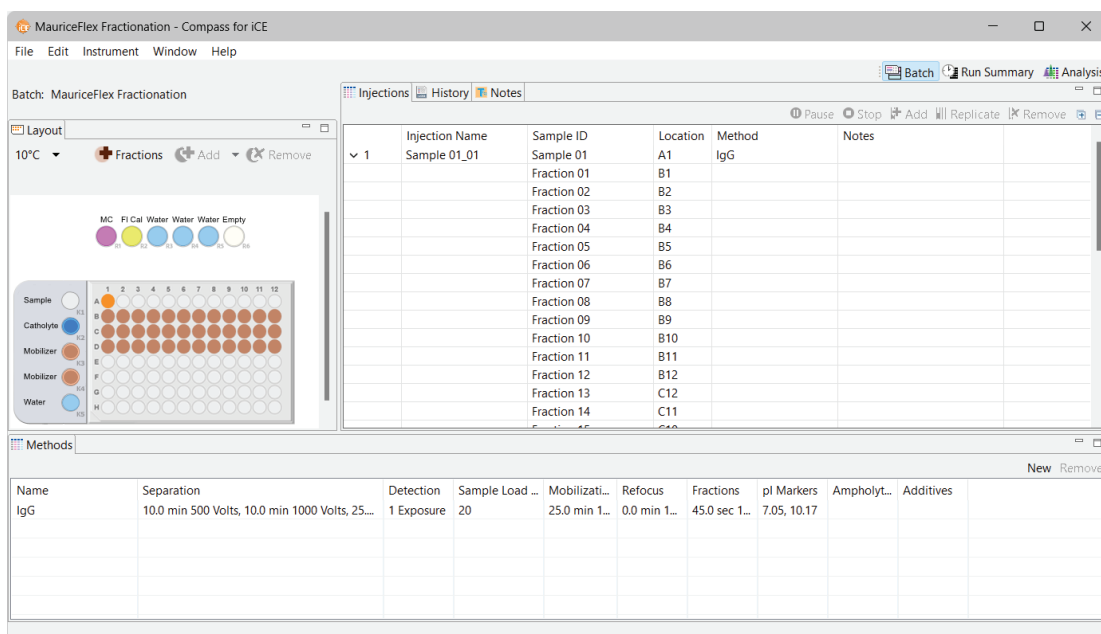
You can use the Batch screen to create, view and edit batches. To get to this screen, click the **Batch** screen tab:



Batch Screen Panes

The Batch screen has five panes:

- **Layout** - Displays a map of the 96-well plate for MauriceFlex Fractionation batch sample locations. Batch reagent locations are also displayed.
- **Injections** - Lists the injections, sample ID, sample locations and methods that MauriceFlex will execute for each sample in the batch.
- **History** - Lists all batch file events from initial creation to the most current update.
- **Notes** - Lets you enter specific information about your batch.
- **Methods** - Lets you create methods and enter method parameters used in the batch.



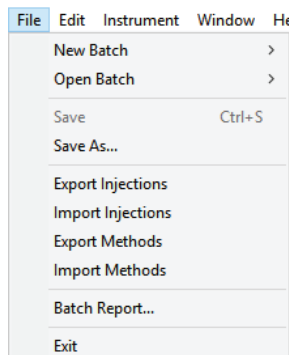
Software Menus Active in the Batch Screen

These main menu items are active in the Batch screen:

- File
- Edit
- Instrument (when the software is connected to MauriceFlex)
- Window
- Help

File Menu

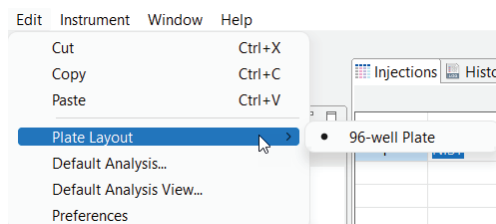
These File menu options are active:



- **New Batch** - Creates a new batch from a starter template.
- **Open Batch** - Opens an existing batch.
- **Save/Save As** - Saves the open batch.
- **Export Injections** - Exports injections from the current batch as a .csv file.
- **Import Injections** - Imports injections into the current batch from a .csv file.
- **Export Methods** - Exports method(s) from the current batch as separate files.
- **Import Methods** - Imports saved method(s) into the current batch.
- **Batch Report** - Exports a table of sample and method details for each injection in the batch as a PDF file.
- **Exit** - Closes Compass for iCE.

Edit Menu

The following Edit menu options are active in the Batch screen:



- **Cut** - Cuts the information currently selected.
- **Copy** - Copies the information currently selected.
- **Paste** - Pastes the copied information.

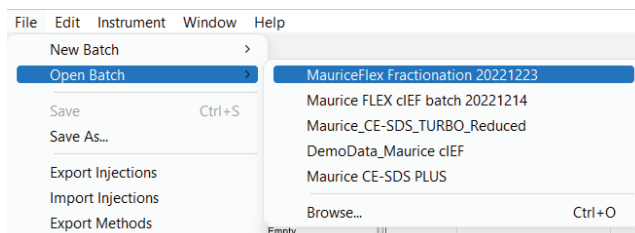
NOTE: Cut, Copy and Paste work differently depending on the pane you're in. How and when you can use them is described throughout the chapter.

- **Plate Layout** - Indicates a 96-well plate will be used to run samples.
- **Default Analysis** - Displays the default settings that will be used to analyze the data generated with your batch.
- **Default Analysis View** - Displays the default settings that will be used to view the data generated with your batch.
- **Preferences** - Lets you set and save your preferences for Access Control, data export, graph colors, grouped data and Twitter settings. See Chapter 21: "Setting Your Preferences" for more information.

Opening a Batch

To open an existing batch:

1. Select **File** in the main menu and click **Open Batch**.



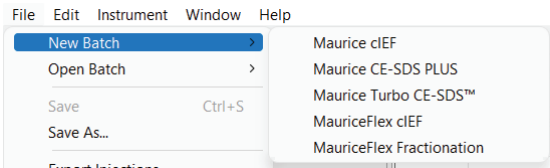
2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file. Select one of those or click **Browse** to open the Batches folder and select a different one.
3. To make changes to the batch, see the steps in "Creating a New Batch" on page 194. When you're done, select **File** from the main menu and click **Save**.

Creating a New Batch

To create a new batch, we recommend that you use one of the template methods and make any necessary modifications from there.

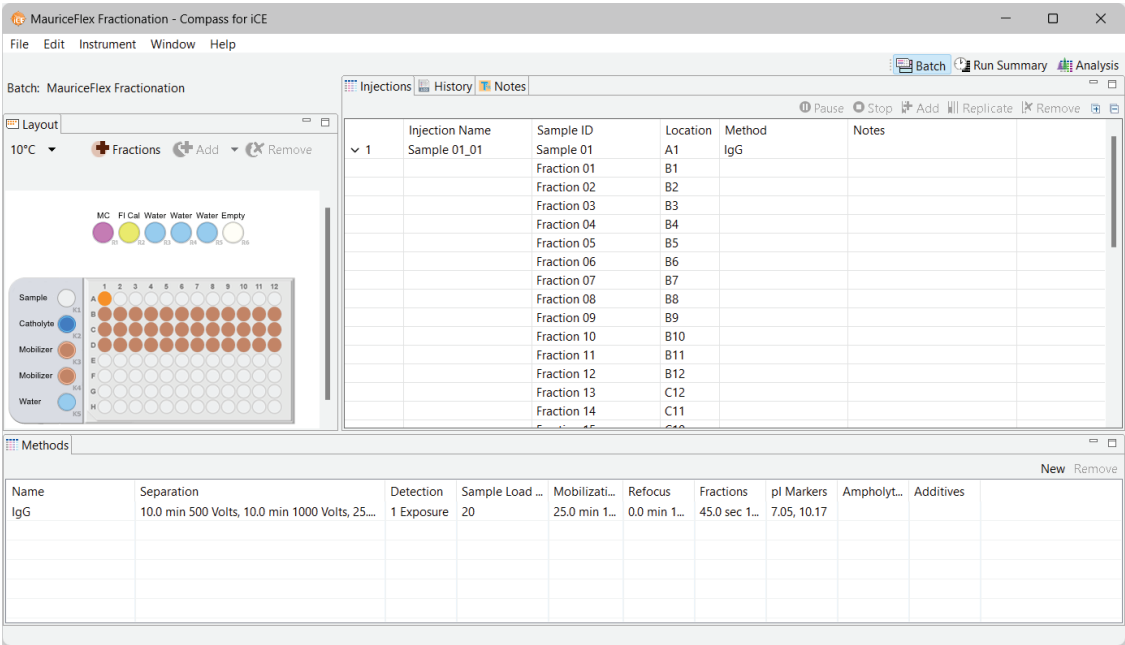
Step 1 - Open a Template Batch

1. Select **File** in the main menu and click **New Batch**:



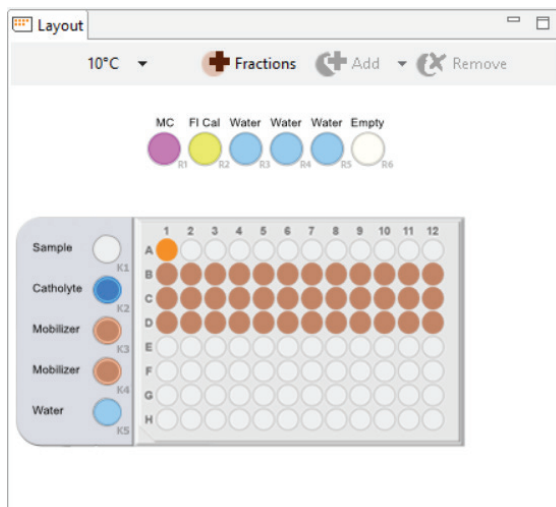
NOTE: Only template batches specific to your Maurice will be available when you are connected to the instrument. Template batches will not appear in the menu if that application type cannot be performed on your instrument.

2. Select **MauriceFlex Fractionation** A batch using the default method will display.



Step 2 - Assign Your Samples

The Layout pane shows the default locations of where batch reagent should be placed in the fractionation adapter.



The same reagent locations are used for every batch:

- **R1** - 0.5% Methyl Cellulose
- **R2** - Fluorescence Calibration Standard
- **R3** - Water vial
- **R4** - Water vial
- **R5** - Water vial
- **R6** - Empty vial (air)
- **K2** - Catholyte Solution
- **K3** - Mobilizer Solution (5 mM ammonium acetate)
- **K4** - Mobilizer Solution (5 mM ammonium acetate)
- **K5** - Water vial

1. Select your sample:

To import a sample using a saved injections file:

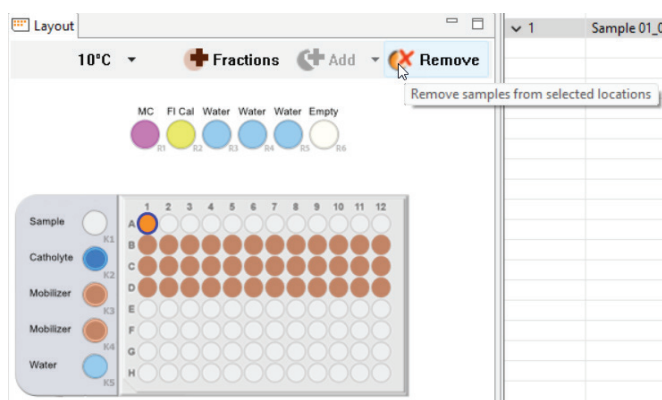
- Select **File** in the main menu and click **Import Injections**.
- Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.
- Skip to step 3 on page 199.

NOTE: For more information on setting up an injections file for import, see “Importing and Exporting an Injection” on page 214.

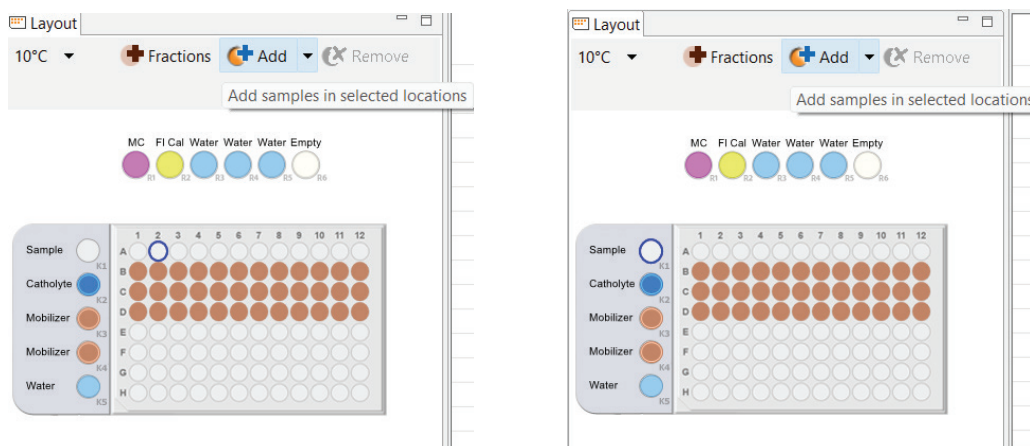
To change the sample location manually and select the method later:

NOTE: The template batch automatically adds the sample in well A1 by default. You can change the sample location to any well in row A or to position K1 if your sample is in a Maurice sample vial with integrated insert.

- a. In the Layout pane, highlight well A1 and click **Remove**. The sample will also be removed from the Injections table. You can also right-click on the well and click **Remove Sample**.

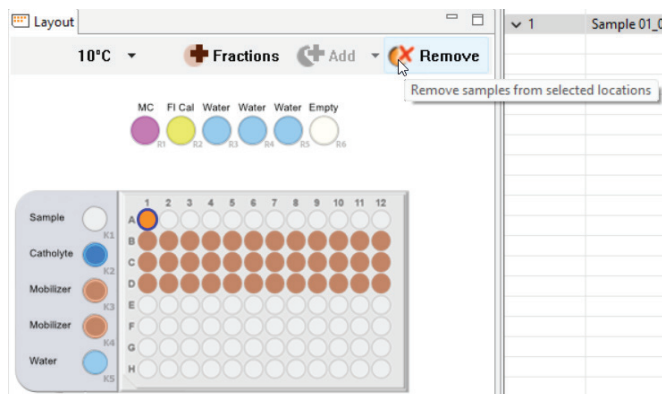


- b. Use your mouse to highlight the well in Row A that your sample is located in or position K1, then click **Add**. The Injections table will automatically update

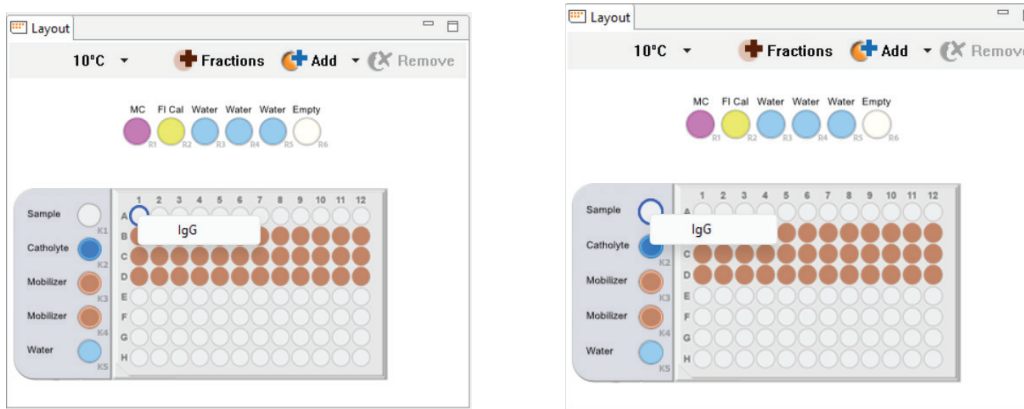


To change the sample location with preassigned methods:

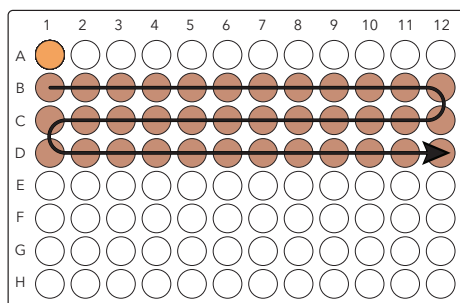
- Highlight well A1 and click **Remove**. The sample will also be removed from the Injections table. You can also right-click on the well and click **Remove Sample**.



- Highlight the well your sample is located in or position K1, then right-click and select a method. The Injections table will automatically update.



- Select where sample peak fractions will be collected in the 96-well plate. All fraction collection starts in column 1 and will be collected in a serpentine pattern. You can define which row to start collecting fractions and how many fractions to collect.



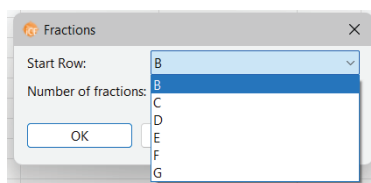
NOTES:

The sample location must first be defined before fraction locations can be assigned.

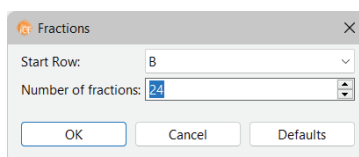
A minimum of 24 fractions must be collected.

To assign fraction collection locations:

- a. In the Layout pane, click **Fractions**.
- b. A prompt will appear. In the **Start Row** pulldown, select the row to start collecting fractions.



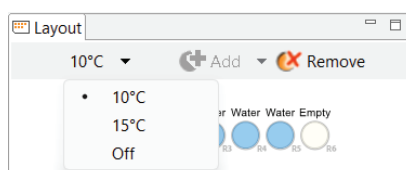
- c. In the **Number of fractions** field, type the number of fractions to collect or use the arrows to increase or decrease the number of fractions.



NOTE: The number entered will be highlighted in red if the number is below the minimum (24) or higher than the available wells on the plate based on where you start collecting fractions.

- d. Click **Defaults** to reset fields to default settings.
 - e. Click **OK**. The plate map in the Layout pane and the Injections list will automatically update.
3. The 96-well plate on the fractionation adapter is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The reagent row (R) is kept at room temperature while the reagent column (K) and sample plate are temperature controlled.



Step 3 - Assign Your Method Parameters

NOTE: We recommend using the default method parameters. Please see the Maurice cIEF Fractionation Method Development Guide for more information on method optimization.

The Methods pane lists all of the methods used in a batch. You can add and remove methods here and change method parameters.

To import a saved method:

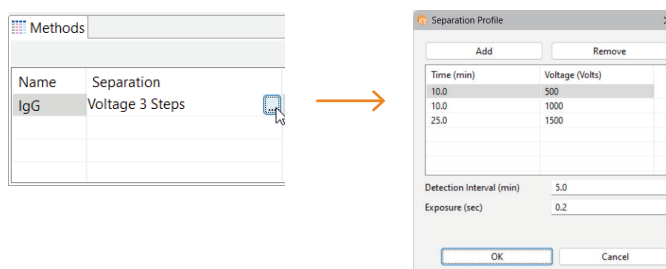
1. Select **File** in the main menu and click **Import Method**.
2. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

To create a new method or make changes to an existing one:

1. In the Methods table, click the first cell in the Name column and enter a new method name if needed.

Name	Separation	Detection	Sample Load ...	Mobilizati...	Refocus	Fractions	pl Markers	Ampholyt...	Additives
IgG	10.0 min 500 Volts, 1...	1 Exposure	20	25.0 min 1...	0.0 min 150...	45.0 sec 1...	7.05, 10.17		

2. Click the first cell in the Separation column, then click the selection button [...] to set the pre-focusing and focusing time (in minutes) and voltage (V). You can also set the Detection Interval (in minutes) and Exposure (in seconds) to adjust how often an image is taken using the listed Exposure time.



- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
- **To change the detection interval or exposure parameters:** Click on the field and type the new value(s).
- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a load time (in minutes) and voltage value (in V).
- **To remove a profile step:** Select the row you want to remove and click **Remove**.

3. The Detection Profile exposure setting is linked to and defined by the Separation step and cannot be changed.
4. Click the first cell in the Sample Load(s) column and set the load time in seconds.

NOTE: We recommend using the default Sample Load time of 20 seconds. Please contact ProteinSimple Technical Support if you have questions on the Sample Load time to use for your application.

Methods			
Name	Separation	Detection	Sample Load (s)
IgG	10.0 min 500 Volts, 1...	1 Exposure	20

5. Click the first cell in the Mobilization column, then click selection button [...] to set the mobilization time (in minutes) and voltage (V). You can also set the Detection Interval (in minutes) and Exposure (in seconds) to adjust how often an image is taken using the listed Exposure time.

Methods				
Name	Separation	Detection	Sample Load ...	Mobilization
IgG	10.0 min 500 Volts, ...	1 Exposure	20	25.0 min 1000 Vo... ...

→

Time (min)
Voltage (Volts)

Detection Interval (min)
Exposure (sec)

- To change time and voltage parameters: Just click in a cell under **Time** or **Voltage** and type the new value(s).
- To change the detection interval or exposure parameters: Click on the field and type the new value(s).

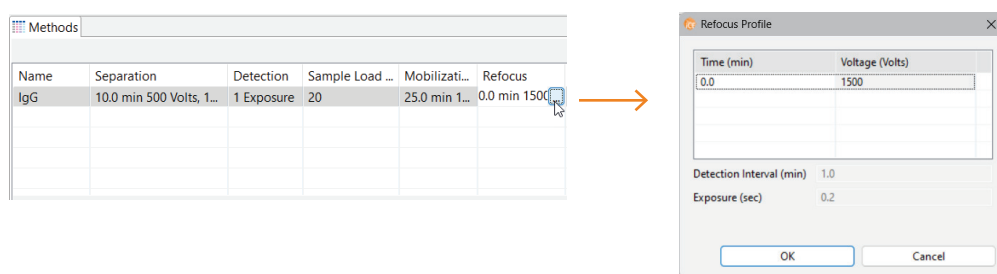
6. An optional refocusing step can be added right before fraction collection to improve fraction resolution. Click the first cell in the Refocus column, then click the selection button [...] to set the refocusing time (in minutes).

NOTES:

The Compass for iCE fraction peak predictor will not be available if a refocusing step is added to the method. A warning will display in the refocus profile and Start Run window and will be recorded in the run summary when the time parameter is changed.

The refocusing step voltage is linked and defined by the Separation step and cannot be changed.

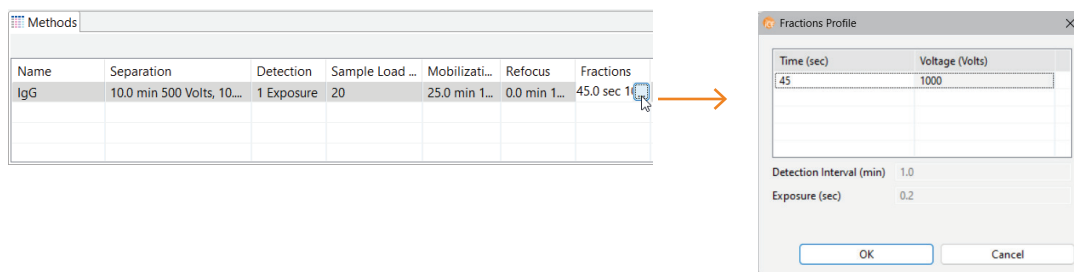
Refer to the MauriceFlex cIEF Fractionation Method Development Guide for refocusing times and voltage setting tips.



- **To change time:** Just click in a cell under **Time** and type the new value(s).

7. Click the first cell in the Fractions column, then click the selection button [...] to set the fraction collection time (in seconds) and voltage (V).

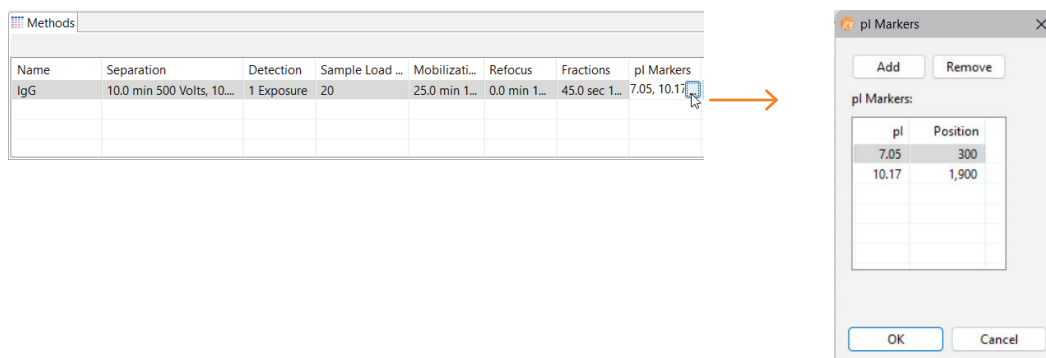
NOTE: The voltage, detection interval and exposure settings are linked to and defined by the Mobilization step settings and cannot be changed.



- **To change time parameters:** Just click in a cell under **Time** and type the new value(s).

8. Click the first cell in the pI Markers column to select pI markers. Add new markers or remove existing ones then click **OK**.

NOTE: When you edit the pI markers in the method for a batch, Compass for iCE automatically creates a Markers group in the pI Markers Analysis settings for you.



- **To change a pI marker and position:** Just click in a cell under **pI** or **Position** and type the new value(s).
 - **To add a pI marker:** Click **Add**. A new row will be added in the table. Then just type in a pI and a position (in pixels).
 - **To remove a pI marker:** Select the row you want to remove and click **Remove**.
9. Optional: Click the first cell in the Ampholytes column and enter the ampholytes you're using.

Methods									
Name	Separation	Detection	Sample Load ...	Mobilizati...	Refocus	Fractions	pI Markers	Ampholytes	
IgG	10.0 min 500 Volts...	1 Exposure	20	25.0 min 1...	0.0 min 1...	45.0 sec 1...	7.05, 10.17	Pharmalyte 3-10	

10. Optional: Click the first cell in the Additives column and enter any additives you're using.

Methods									
									New Remove
Name	Separation	Detection	Sample Load ...	Mobilizati...	Refocus	Fractions	pI Markers	Ampholytes	Additives
IgG	10.0 min 500 Volts, 10.0...	1 Exposure	20	25.0 min 1...	0.0 min 1...	45.0 sec 1...	7.05, 10.17	Pharmalyte 3-10	Urea

11. You can now:

- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.
- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.

Step 4 - Set Up Your Injection

The Injection pane lists information for the batch sample injection and lists fractionation collection locations in the 96-well plate.

Injection Name	Sample ID	Location	Method
▼ 1 Sample 01_01	Sample 01	A1	IgG
	Fraction 01	B1	
	Fraction 02	B2	
	Fraction 03	B3	
	Fraction 04	B4	
	Fraction 05	B5	
	Fraction 06	B6	
	Fraction 07	B7	
	Fraction 08	B8	
	Fraction 09	B9	
	Fraction 10	B10	
	Fraction 11	B11	
	Fraction 12	B12	
	Fraction 13	C12	
	Fraction 14	C11	
	Fraction 15	C10	
	Fraction 16	C9	
	Fraction 17	C8	
	Fraction 18	C7	
	Fraction 19	C6	
	Fraction 20	C5	
	Fraction 21	C4	
	Fraction 22	C3	
	Fraction 23	C2	
	Fraction 24	C1	
	Fraction 25	D1	

NOTE: If you imported the injection in Step 2, you can update the injections table and assign a method as needed using the information that follows.

1. To change the sample name, click the **Sample ID** cell and type a name.

NOTES:

A Sample ID can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting an Injection Name or Sample ID from other Documents” on page 213 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		
		Fraction 02	B2		
		Fraction 03	B3		

The sample name also displays when you hover the mouse over the sample in the plate map:

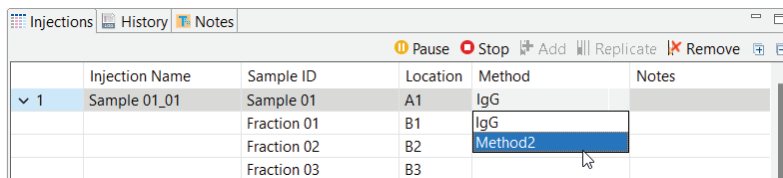
	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		
		Fraction 02	B2		
		Fraction 03	B3		
		Fraction 04	B4		
		Fraction 05	B5		
		Fraction 06	B6		
		Fraction 07	B7		
		Fraction 08	B8		
		Fraction 09	B9		
		Fraction 10	B10		
		Fraction 11	B11		

2. The Injection name is automatically set to ‘Sample ID_injection number’. To change the name, click the **Injection Name** cell and type a name. The injection name must be unique. If it isn’t when you save the batch, Compass for iCE highlights any non-unique injection names so you can change them.

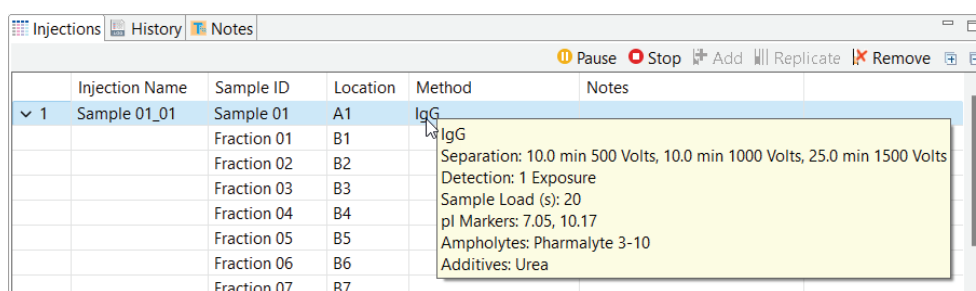
	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		
		Fraction 02	B2		

NOTE: Changing the injection name won’t affect the sample ID.

- Assign your methods to each injection. Just click the **Method** cell for the injection and select a method from the drop down menu. If you added your samples with pre-assigned methods and don't need to make any changes, skip to the next step.



Hovering over a method name displays the method parameters:



Step 5 - Add Programmed Pauses and Stops (Optional)

You can program pauses and stops in the batch to automatically pause or stop after injections even when you aren't in front of MauriceFlex.

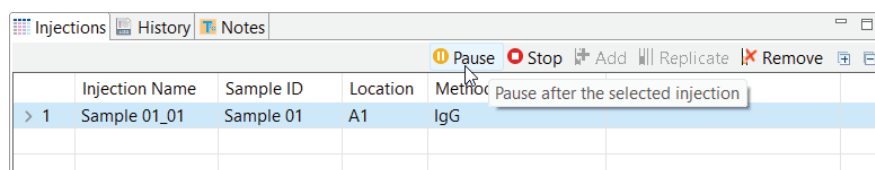
Note: The cartridge cleanup is performed at the end of the batch and after an injection. A programmed pause will occur after the injection before the cartridge cleanup starts.

To program a pause:

Pauses are helpful when you want to review injection results before continuing with the rest of the batch.

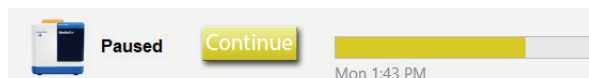
- Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete.

NOTE: MauriceFlex can tweet you when the batch pauses. See “Setting Up Maurice Systems to Send Tweets” on page 764.



Injections History Notes						
<div> Pause Stop Add Replicate Remove </div>						
	Injection Name	Sample ID	Location	Method	Notes	
> 1	Sample 01_01	Sample 01	A1	IgG		
				Pause		

- To resume the batch, click **Continue** in the instrument status bar:



To stop the run after a specific injection:

- Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.

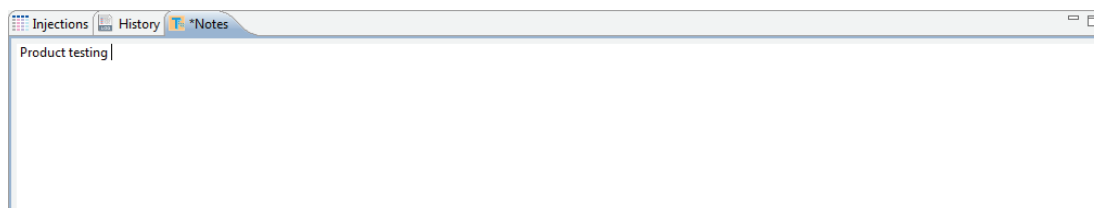
Injections History Notes						
<div> Pause Stop Add Replicate Remove </div>						
	Injection Name	Sample ID	Location	Method	Notes	
> 1	Sample 01_01	Sample 01	A1	IgG		
				Stop		

Data collected for injections completed up to the point where the batch was stopped are saved in the .mbz file.

- Be sure to complete your normal post-batch procedures.

Step 6 - Add Batch Notes (Optional)

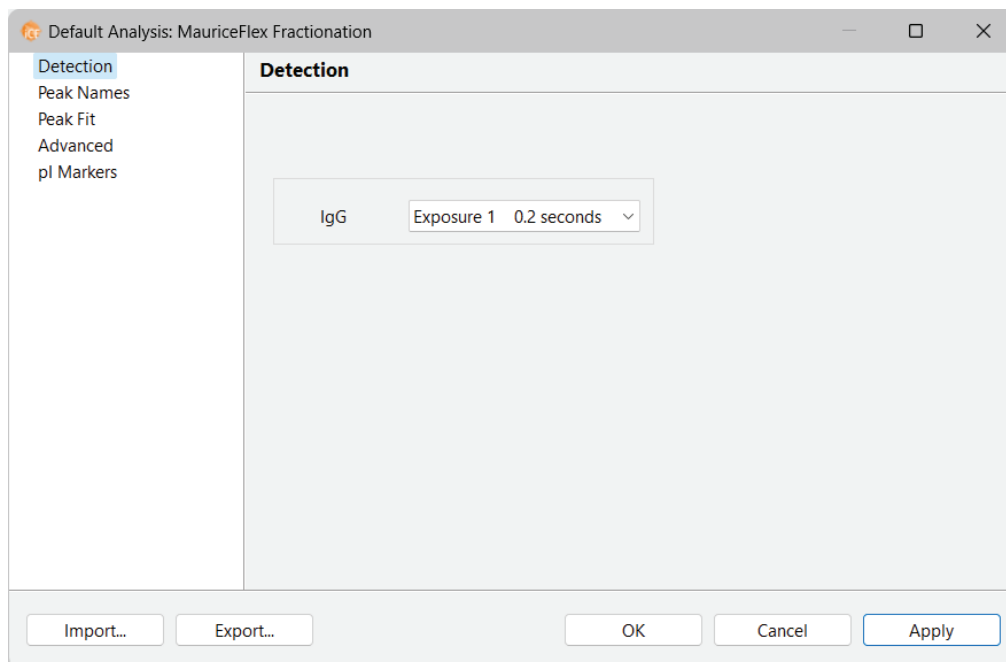
- Click on the **Notes** pane.
- Click in the notes area and type any information you want to add about your batch.



Step 7 - Modify Default Analysis Parameters (Optional)

You can preset the parameters used to analyze data generated with the batch. We recommend using the default parameters for MauriceFlex Fractionation applications, but if you need to modify parameters:

1. Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



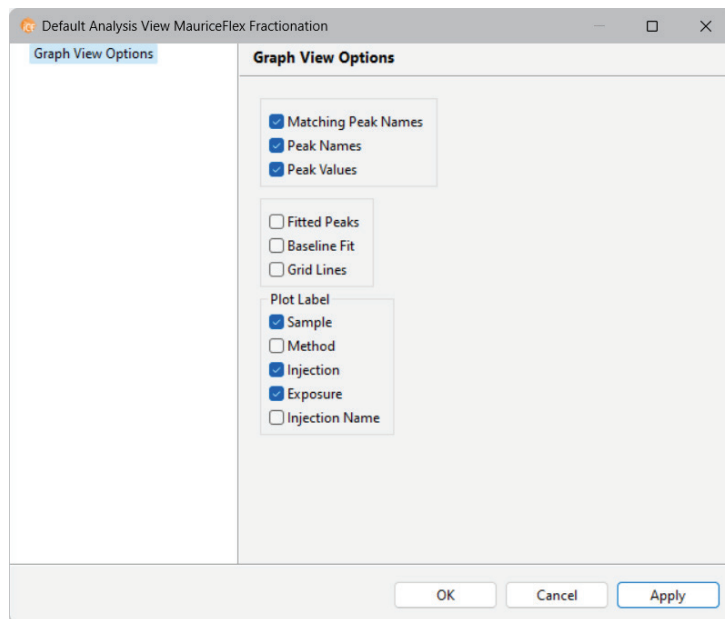
2. Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to "Analysis Settings Overview" on page 620.

Step 8 - Modify Default Analysis View Parameters (Optional)

You can preset the parameters used to display the electropherogram data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button.

To modify the parameters:

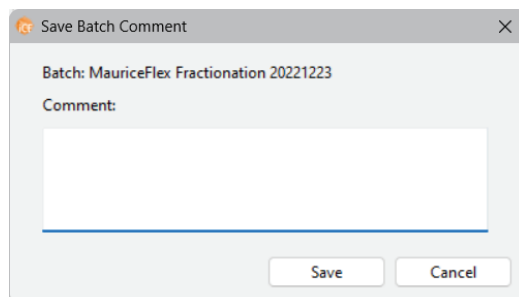
1. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



2. Change the parameters you want to, then click **Apply**. Click **OK** when you are done changing display parameters. For detailed information on on graph view options, please refer to “Customizing the Data Display” on page 603.

Step 9 - Save Your Batch

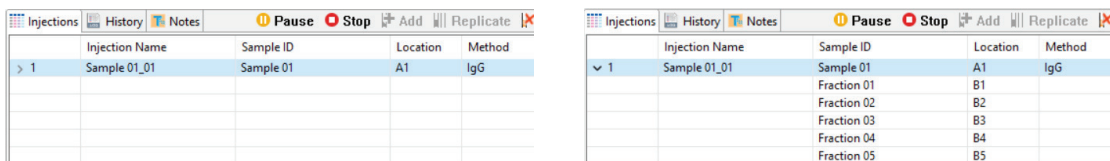
1. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



2. Enter a name for your batch then click **Save**.

Viewing Fractions To Be Collected

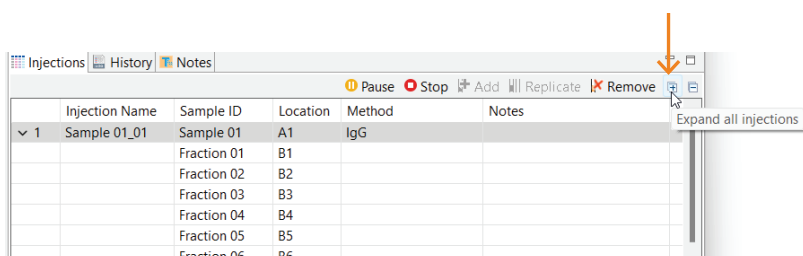
An arrow will appear next to the injection number. Toggle the arrow to expand or collapse the list of fractions that will be collected.



	Injection Name	Sample ID	Location	Method
> 1	Sample 01_01	Sample 01	A1	IgG

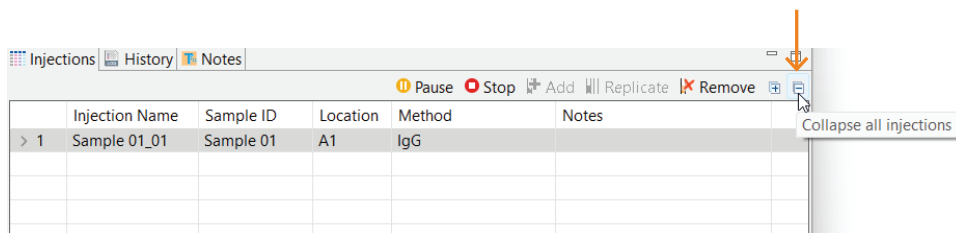
	Injection Name	Sample ID	Location	Method
▼ 1	Sample 01_01	Sample 01	A1	IgG
		Fraction 01	B1	
		Fraction 02	B2	
		Fraction 03	B3	
		Fraction 04	B4	
		Fraction 05	B5	

- To show all fractions collected in the batch, click the **Expand All Injections** button.



	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		
		Fraction 02	B2		
		Fraction 03	B3		
		Fraction 04	B4		
		Fraction 05	B5		
		Fraction 06	B6		

- To hide all fractions collected in the batch, click the **Collapse All Injections** button.



	Injection Name	Sample ID	Location	Method	Notes
> 1	Sample 01_01	Sample 01	A1	IgG	

Batch History

Clicking on the History pane shows the batch event history, starting with the date and time the batch was created through the most current event. Clicking on a row in the table displays the full event details in the box under the table.

[illegible]

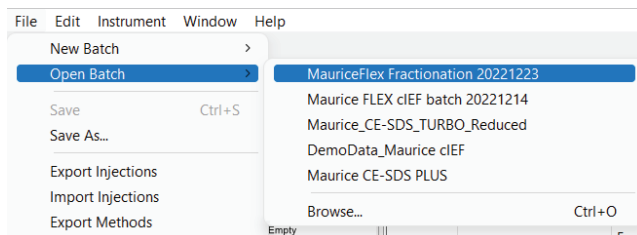
- **Date:** Date and time of the batch event.
- **User Name:** User that initiated the event. User names only display if you're using the Access Control feature to help satisfy the 21CFR Part 11 data security requirements. Go to "Enabling Access Control" on page 781 to learn how to set it up.
- **Message:** Description of the event that took place.
- **Comment:** Comments entered by the user when the batch was saved.

You can copy the information in the History pane to use in other documents:

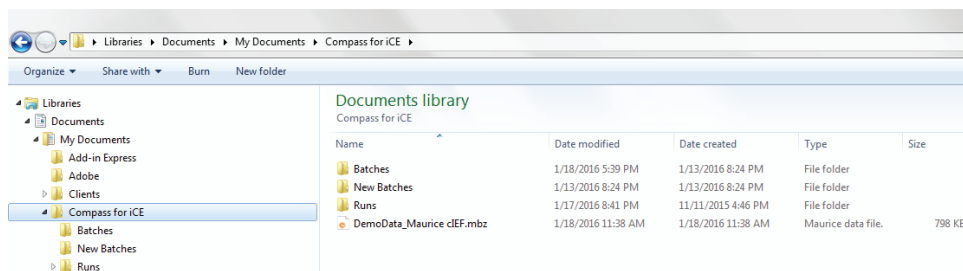
1. Click the **History** pane to make sure it's active.
2. Click **Edit** in the main menu and select **Copy**.
3. Open a document and click **Paste**.

Making Changes to a Batch

1. Select **File** in the main menu and click **Open Batch**.



2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file.
 - Select one of these files or click **Browse** to open the Batch folder and select a different one.
 - To open a batch used in a run file, click **Browse**. Go to the Run folder in the Compass for iCE directory and select the run file that uses the batch you want to edit.

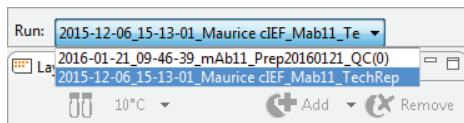


3. To make changes to the batch, see the steps in “Creating a New Batch” on page 194. Then select **File** from the main menu and click **Save** or **Save As**.

Viewing and Editing Batches in Completed Runs

1. Click the **Analysis** screen and open your run file(s).
2. After the run opens, click the **Batch** screen to view the batch used with the run.

If you opened more than one run file, you can switch between viewing batches for each file. Just click the **down arrow** in the Run box and select the batch you want to view from the drop down list.



3. In completed run files, you can only make edits to the **sample** and **method** names in a batch, and these changes are saved to the run file only. Update sample and method names as needed.

- Click the **Analysis** screen. Then select **File** from the main menu and click **Save** or **Save As** to save the new changes to the run file.

Copying and Pasting an Injection Name or Sample ID from other Documents

An injection name and Sample ID can be copied from other documents and pasted into the Injection pane to make batch setup faster.

To paste an injection name into the batch:

- Select the injection name in a document (Microsoft®, Word®, Excel®, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3			

- Select the injection cell in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the injection, right click and select **Paste**.

Injections	History	Notes	Pause	Stop	Add	Replicate	Remove			
1	Injection Name	Sample ID	Location	Method	Notes					
1	Sample 01_01	Sample 01	A1	IgG						
		Fraction 01	B1							
		Fraction 02	B2							

The injection name is pasted into the Injection pane:

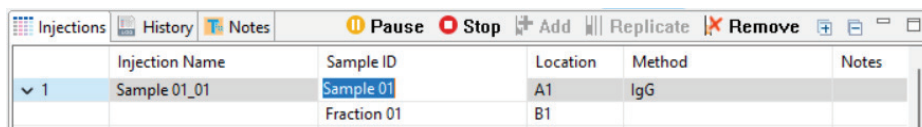
Injections	History	Notes	Pause	Stop	Add	Replicate	Remove			
1	Injection Name	Sample ID	Location	Method	Notes					
1	My Injection 1	Sample 01	A1	IgG						
		Fraction 01	B1							
		Fraction 02	B2							

To paste a Sample ID into the batch:

- Select the Sample ID in a document (Microsoft®, Word®, Excel®, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3			

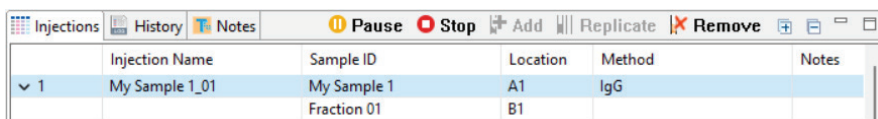
2. Select a Sample ID in the Injection pane, then hold the Ctrl+V keys to paste. You can also select the Sample ID, right click and select **Paste**.



	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		

The Sample ID is pasted into the Injection pane.

NOTE: If you paste in a new sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you've already changed the injection name manually, updating the sample ID will not update the injection names.



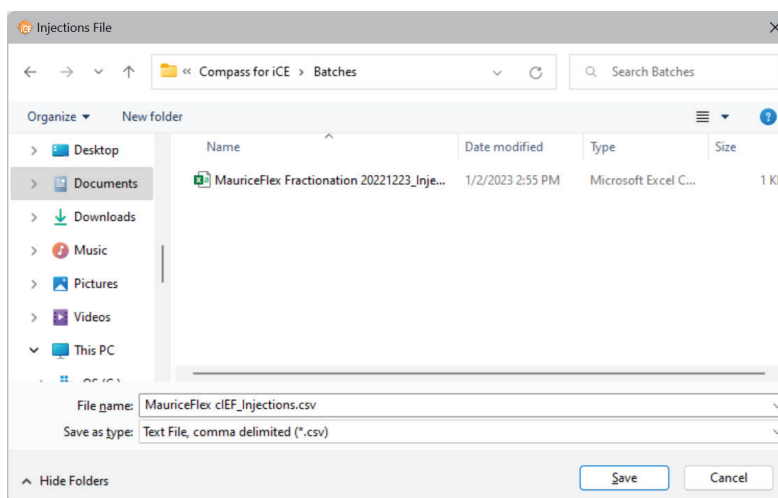
	Injection Name	Sample ID	Location	Method	Notes
1	My Sample 1_01	My Sample 1	A1	IgG	
		Fraction 01	B1		

Importing and Exporting an Injection

An injections in an open batch or run file can be exported as a separate file. This lets you import the same or modified injection into another batch later, rather than having to enter the information manually.

Exporting an Injection

1. Open the batch or run you want to export the injection from.
2. In the Batch screen, select **File** in the main menu and click **Export Injections**. The following window displays:



3. The default directory is Compass for iCE/Batches. Change the directory if needed.
4. The default file name will be batchname.batch_Injections.csv. Change the name if needed and click **Save**.

Creating an Injection File

To create a new injection file .csv, it's easiest to export the injection from an existing batch and use that as your template.

1. Follow the instructions in "Exporting an Injection" above to export the injection from an existing cIEF batch.
2. Open the .csv file in a program that provides a table/spreadsheet format.
3. To add or change injection information:
 - a. Type in a unique name in each cell in the Injection Name column.
 - b. Type in a sample ID, location and method for each injection in their respective columns.
 - c. Optional: Type in notes if needed.

NOTES:

Mix bottles are not used for MauriceFlex cIEF batches. Compass for iCE ignores this column when importing injections into a MauriceFlex cIEF batch.

Programmed pauses and stops can't be added in the injections file. They can be added in the Injections pane after importing your injections into Compass for iCE.

	A	B	C	D	E	F
1	Injection Name	Sample ID	Location	Method	Mix Bottle	Notes
2	Inj01	R01_SSPP	E4	System Suitability		

4. Save the .csv file.

Importing an Injection

1. Open the batch you want to import the injection into, or open a new batch.
2. Select **File** in the main menu and click **Import Injections**.
3. Select an injections file (*.csv) and click **OK**. The imported information will display in the Layout and Injections panes.

NOTE: The imported injection will replace any injections already in the Injections table, they will not append an existing table.

Importing and Exporting Methods

Methods in an open batch or run file can be exported as a separate file. This lets you import the same method into another batch later, rather than having to re-enter method information manually.

Importing a Method

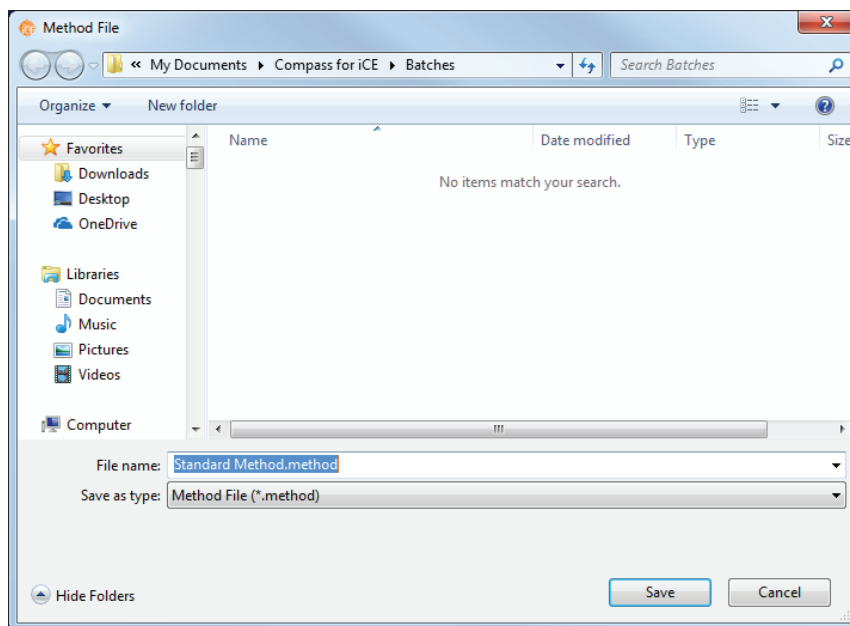
NOTE: Importing a method imports information into the Batch window's Method pane only.

1. Open the batch you want to import the method into.
2. Select **File** in the main menu and click **Import Method**.
3. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

Exporting a Method

NOTE: Exporting a method exports information in the Batch window's Method pane only.

1. Open the batch you want to export the method from.
2. In the Method pane, select the method you want to export. To export more than one method at the same time, hold the **Ctrl** key to select multiple methods.
3. Select **File** in the main menu and click **Export Method**. The following window displays:



4. The default directory is Compass for iCE/Batches. Change the directory if needed.
5. Enter a method file name and click **Save**. The settings will be saved as a *.method file.

Batch Reports

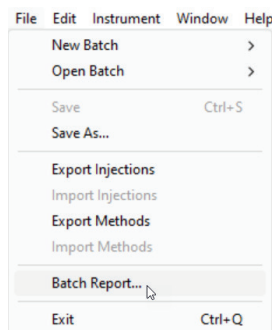
You can export a PDF file of sample and method details for each injection in the batch.

NOTES:

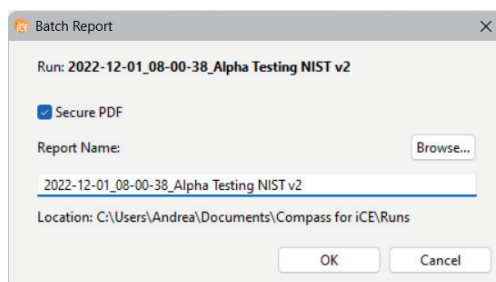
You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it’s saved in a separate run folder. Existing reports aren’t overwritten.

1. Go to the **Analysis** or **Run Summary** screen, then click **File > Open Run** and select a run file (if you don’t have one open already).
2. After the run opens, go to the **Batch** screen.
3. Select **File** from the main menu and click **Batch Report**.

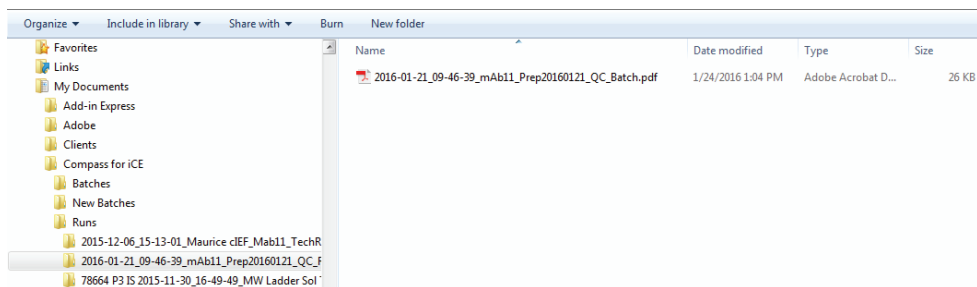


4. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.



Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don’t want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.

5. The Batch Report PDF is exported to the Runs folder in the Compass for iCE directory. It'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Here's an example Batch Report:

MauriceFlex Fractionation Batch: Alpha Testing NIST

Injections

Injection	Injection Name	Sample ID	Location	Method	Notes
1	NIST	Sample 01	A1	NIST	
		Fraction 01	B1		
		Fraction 02	B2		
		Fraction 03	B3		
		Fraction 04	B4		
		Fraction 05	B5		
		Fraction 06	B6		
		Fraction 07	B7		
		Fraction 08	B8		
		Fraction 09	B9		
		Fraction 10	B10		
		Fraction 11	B11		
		Fraction 12	B12		
		Fraction 13	C12		
		Fraction 14	C11		
		Fraction 15	C10		
		Fraction 16	C9		
		Fraction 17	C8		
		Fraction 18	C7		
		Fraction 19	C6		
		Fraction 20	C5		
		Fraction 21	C4		
		Fraction 22	C3		
		Fraction 23	C2		
		Fraction 24	C1		

Created By: Andrea Fri 9:24 AM Feb 24, 2023 PST (SECURED)
 C:\Users\Andrea\Documents\Compass for iCE\Runs\2022-12-01_08-00-38_Alpha Testing NIST v2.mtz
 Computer: DESKTOP-IPMT002 Software Version: Compass for iCE 4.0.0, Build ID: 0222

Page 1 of 2



MauriceFlex Fractionation Batch: Alpha Testing NIST

Methods

Name	Separation	Detection	Sample Load (s)	Mobilization	Refocus	Fractions	pl Markers
NIST	10.0 min, 500 Volts 10.0 min, 1000 Volts 25.0 min, 1500 Volts Detection Exposure: 0.1 sec Interval: 5.0 min	Fluorescence, 0.1 sec	30	25.0 min 1000 Volts Detection Exposure: 0.1 sec Interval: 1.0 min	0.0 min 1500 Volts Detection Exposure: 0.1 sec Interval: 1.0 min	45.0 sec 1000 Volts Detection Exposure: 0.1 sec Interval: 1.0 min	7.05 pl, 300 pixels 10.17 pl, 1800 pixels

Methods (continued)

Name	Ampholytes	Additives
NIST		

Batch Log

Date	User Name	Message	Comment
2022-11-04 08:07:40		Batch created using the factory default Maurice Flex Fractions with Compass for iCE Version: 4.0.0-1102	
2022-11-04 08:33:39		Saved as Z:\shared\ppd\science\rd science\team members\mason\Flex Cartridge Validation - Backup\Alpha Testing\Alpha Testing NIST.batch from Compass for iCE v4.0.0-1102	
2022-11-04 08:33:39		Save injections changes to Z:\shared\ppd\science\rd science\team members\mason\Flex Cartridge Validation - Backup\Alpha Testing\Alpha Testing NIST.batch from Compass for iCE v4.0.0-1102	

Created By: Andrea Fri 6:24 AM Feb 24, 2023 PST (SECURED)
 C:\Users\Andrea\Documents\Compass for iCE\Runs\2022-12-01_08-00-38_Alpha Testing NIST v2.mbz
 Computer: DESKTOP-1FM7005 Software Version: Compass for iCE 4.0.0, Build ID: 0222

Page 2 of 2



Chapter 10:

Running MauriceFlex Fractionation Applications

Chapter Overview

- Overview
- Before You Throw the Switch
- Power Up
- MauriceFlex Fractionation Workflow
- Running MauriceFlex Fractionation Applications
- Editing a Running Batch
- Post-batch Procedures
- Checking Your Data

Overview

MauriceFlex Fractionation applications can be run on MauriceFlex system using a cIEF Fractionation cartridge. You can then confirm which fractions contain the desired peak by running a standard cIEF batch on a Maurice, Maurice C. or MauriceFlex system with a cIEF cartridge before mass spectrometry analysis.

NOTES:

We recommend first optimizing methods with cIEF batches using a cIEF cartridge on Maurice, Maurice C. or MauriceFlex. Then, use recommendations in the MauriceFlex cIEF Fractionation Method Development Guide for tips to modify the method for a MauriceFlex cIEF batch using a cIEF Fractionation cartridge on MauriceFlex to confirm successful method transfer before performing a MauriceFlex Fractionation batch.

cIEF Fractionation cartridges support 15 injections.

On-board mixing is not available on MauriceFlex.

Before You Throw the Switch

Ensure that everyone using MauriceFlex have:

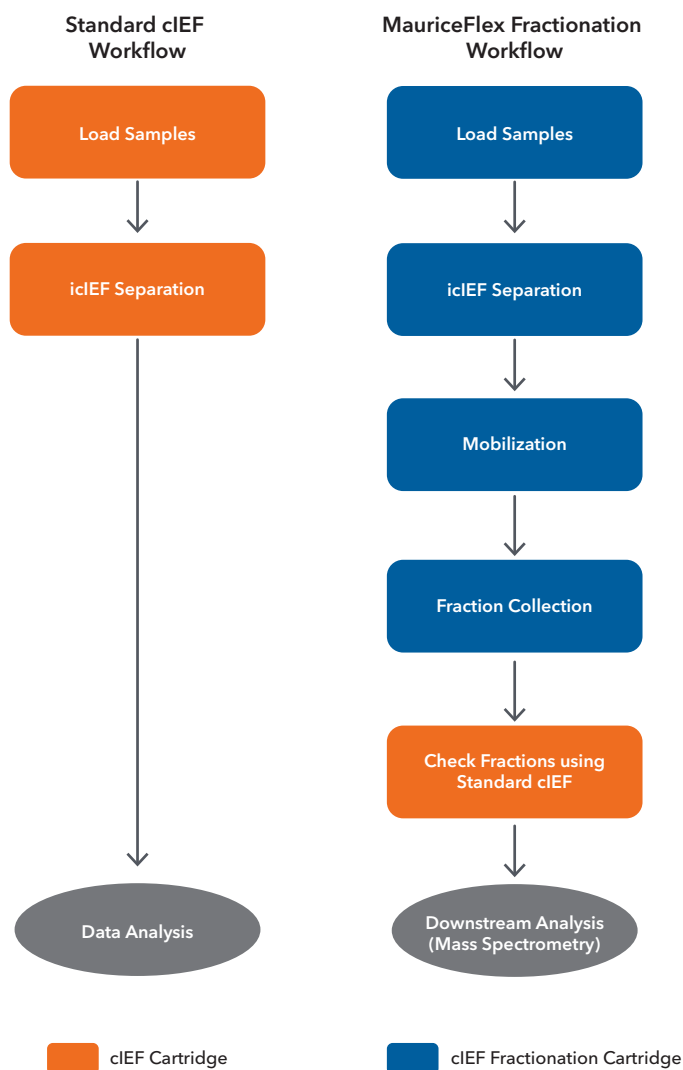
- Received instruction in general safety practices for laboratories.
- Received instruction in specific safety practices for MauriceFlex.
- Received instruction on handling of biohazards (if biohazardous materials are to be used on MauriceFlex).
- Read and understood all Product Inserts and related Safety Data Sheets (SDS).

Power Up

1. Turn on the computer connected to MauriceFlex.
2. Turn on MauriceFlex's main power switch.
3. Wait for MauriceFlex to initialize. His indicator light will turn from magenta to blue when he's done.
4. Double-click the Compass for iCE software icon on the computer desktop to open the application.
5. Connect MauriceFlex to Compass for iCE.

MauriceFlex Fractionation Workflow

Perform a MauriceFlex Fractionation batch to isolate specific peaks of interest from your Maurice cIEF electropherogram. Then run your fractions using a cIEF batch to confirm the peak identity and purity in your fraction before mass spectrometry analysis.



Comparison of standard cIEF and MauriceFlex Fractionation applications.

Running MauriceFlex Fractionation Applications

What You'll Need

- MauriceFlex cIEF Fractionation Cartridges
- MauriceFlex cIEF Fractionation Method Development Kit (optional)
- Maurice cIEF Fluorescence Calibration Standard
- Maurice cIEF pI Markers (3.38, 4.05, 5.85, 6.14, 7.05, 8.4, 9.50, 9.99, or 10.17)
- 0.5% Methyl Cellulose Solution
- 1% Methyl Cellulose Solution
- iCE Electrolyte Kit
- Deionized (DI) water
- MauriceFlex glass crimp top reagent vials, 2 mL



MauriceFlex glass crimp top reagent vial (PN 110-0019, left) and glass reagent vial (PN 046-017, right)
Only MauriceFlex glass crimp top reagent vials (PN 110-0019) should be used to prepare MauriceFlex Fractionation batch reagents.

- Glass vials with insert, 0.3 mL for the Fluorescence Calibration Standard 96-well plate for samples
- P10, P20, P200, P1000 pipettes and tips
- Electrolyte pipette
- Vortexer
- Table-top centrifuge with plate adapter or vial adapter (12 mm, 2 mL vials)

IMPORTANT:

Use MauriceFlex glass crimp top reagent vials, 2 mL (PN 110-0019) and glass vials with insert, 0.3 mL (PN 110-0018) to prepare MauriceFlex Fractionation batch reagents.

Step 1: Prep Your Markers, Samples and Reagents

NOTE: You can prepare your samples in a 96-well plate or vial.

Mobilizer Solution

1. Add 20 μ L of 2 M Ammonium Acetate to 8 mL of DI or LC-MS grade water and mix thoroughly.

NOTE: Use LC-MS grade water if fractions will be directly analyzed by mass spectrometry.

2. Store the prepared Mobilizer Solution on ice until ready to dispense in the 96-well plate.

Samples

1. In a microcentrifuge tube, prepare your sample at a concentration of 2.5–10 mg/mL in a final volume of 25 μ L in DI water.
2. In a separate tube, prepare IEF Separation Mix containing your chosen pI marker(s).

NOTES:

Check out the MauriceFlex cIEF Fractionation Method Development Guide for suggested IEF Separation Mix recipes.

If you are running a method optimized for cIEF analysis, increase the concentration of arginine and both pI markers in your IEF Separation Mix. Check out the MauriceFlex cIEF Fractionation Method Development Guide for more information.

3. Add 100 μ L of IEF Separation Mix to the 25 μ L of your sample.
4. Thoroughly vortex the sample to mix completely.
5. Centrifuge the tube at 13,000 xg for 5 minutes to remove air bubbles and sediment any particulates.
6. Carefully aspirate the top 100 μ L of the sample and pipette it into a well in a 96-well plate by inserting the pipette tip all the way to the bottom to avoid introducing bubbles.
7. Pipette 30 μ L of Mobilizer Solution (5 mM ammonium acetate) to each fraction well in a 96-well plate as defined by your MauriceFlex Fractionation batch. The number of wells will depend on the number of fractions being collected.
8. Spin the plate for 5 minutes at 1000 xg using a centrifuge plate adapter.

NOTES:

If you're preparing sample in a sample vial with insert, 0.2 mL (PN 046-083), increase the sample and IEF Separation Mix volumes in steps 1–3 by 3-fold and transfer a minimum of 300 μ L of sample to the vial.

Make sure to check for and remove any bubbles at the bottom of the well.

pI Markers

1. Open the vial of lyophilized pI marker by lifting the center tab and gently pulling it back to break the metal seal. Slowly remove the rubber stopper.
2. Add 210 μL of DI water to the vial.
3. Put the rubber stopper back in the vial and vortex 3 times, 5 seconds each time to completely reconstitute the lyophilized cake. Repeat this step for all pI markers you're using.
4. Aliquot 20 μL of each reconstituted pI marker into separate tubes for storage.

NOTES:

Keep your reconstituted pI markers on ice until you're ready to add them to your sample or IEF Separation Mix.

If you'll use the pI markers within a month, store the aliquots at 2-8 °C. Otherwise, store the aliquots at -10 °C to -30 °C. They'll be stable up to 6 months.

5. Use 4 μL of each pI marker for every 200 μL of sample.

Reagents

IMPORTANT:

Use MauriceFlex crimp top glass reagent vials, 2 mL (PN 110-0019) and glass vials with insert, 0.3 mL (PN 110-0018) to prepare MauriceFlex Fractionation batch reagents. Glass reagent vials, 2 mL (PN 046-017) should only be used when running a standard cIEF or CE-SDS batch.

NOTE: Don't reuse reagents or vials

1. Pipette 350 μL of Fluorescence Calibration Standard in a MauriceFlex glass vial with insert, 3 mL (PN 110-0018) and label.



2. Pipette 2 mL of 0.5% Methyl Cellulose into a MauriceFlex crimp top glass reagent vial and label.



3. Pipette 2 mL of DI water into four MauriceFlex crimp top glass reagent vials. Label the vials.
4. Pipette 2 mL of Catholyte solution to a MauriceFlex crimp top glass reagent vial and label.
5. Pipette 2 mL of Mobilizer Solution (5 mM ammonium acetate) into two MauriceFlex crimp top glass reagent vials. Label the vials.
6. Label an empty MauriceFlex crimp top glass reagent vial.

NOTES:

Make sure you don't overfill the vials, especially the methyl cellulose vial, to avoid introducing bubbles to your run. Wipe excess liquid at the mouth of the vial with a laboratory wipe.

Vials do not need to be capped.

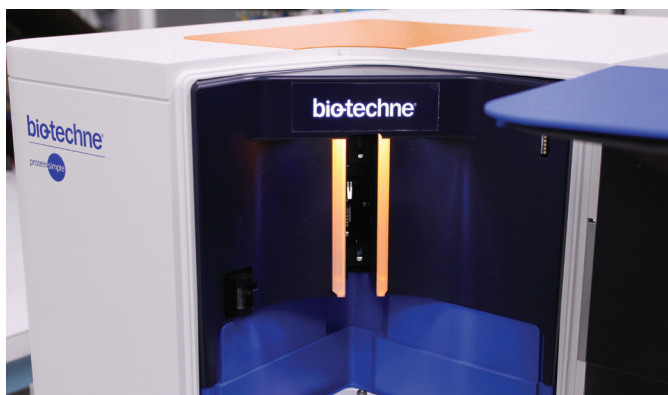
Step 2: Load Samples and Reagents

1. Open MauriceFlex's door by touching the metal plate on top of the door.



NOTE: The indicator light on MauriceFlex's front panel will blink rapidly as the door disengages.

- Swing the door open to access the cartridge slot and the reagents and samples platform. The lights on either side of the slot will be **orange**.

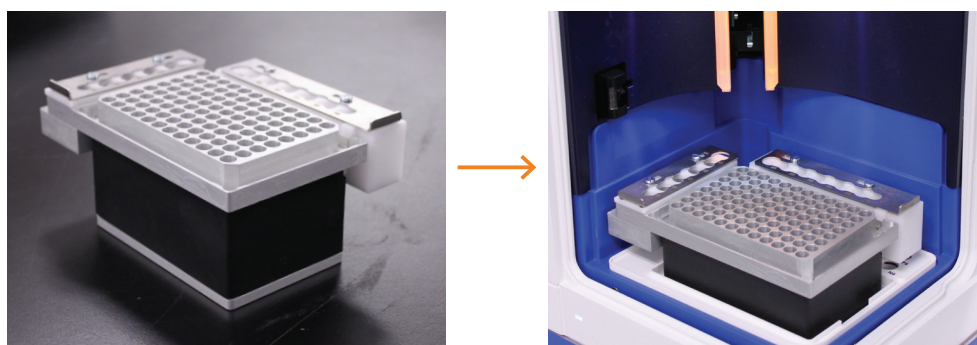


- Install the fractionation adapter in MauriceFlex.

NOTES:

Remove all the vials in rows P and N of the reagent platform before installing the fractionation adapter if a CE-SDS or standard cIEF application was previously performed.

When installing the fractionation adapter, confirm the adapter is seated and sitting flat in MauriceFlex.



- Place the reagent vials into their respective positions on the fractionation adapter.

IMPORTANT:

Reagent vials on MauriceFlex must be locked in place before starting a batch. The vial locking mechanism default state is locked. Pull the locking mechanism on Row R to the right and the locking mechanism on Column K down to unlock and place vials in the fractionation adapter. Release the locking mechanism to lock the vials in place.

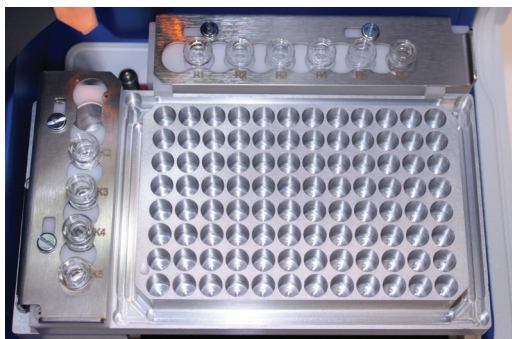
NOTES:

The reagent row (R) on the fractionation adapter is kept at room temperature. The reagent column (K) and sample plate are temperature controlled.

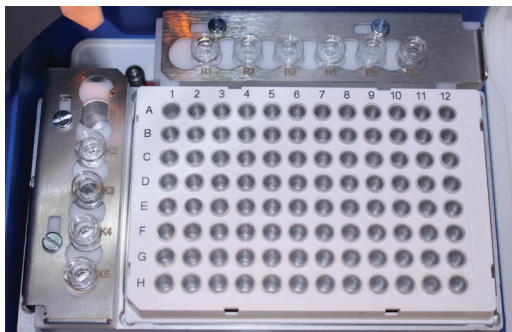
MauriceFlex Fractionation batch reagents should be prepared in 2 mL crimp top glass vials (PN 110-0019) and glass vials with insert, 0.3 mL (110-0018).

Vials do not need to be capped before placing them on the fractionation adapter.

- **R1** - 0.5% Methyl Cellulose
- **R2** - Fluorescence Calibration Standard
- **R3** - Water
- **R4** - Water
- **R5** - Water
- **R6** - Empty vial (air)
- **K2** - Catholyte Solution
- **K3** - Mobilizer Solution (5 mM Ammonium Acetate)
- **K4** - Mobilizer Solution (5 mM Ammonium Acetate)
- **K5** - Water



- Place your 96-well sample plate on the fractionation adapter. Well A1 should be in the top left corner of the adapter.



Step 3: Prep the cIEF Fractionation Cartridge

NOTE: A cIEF Fractionation Cartridge supports a maximum of 15 injections and a maximum of 15 batches. Its RFID will keep track of how many are left for you.

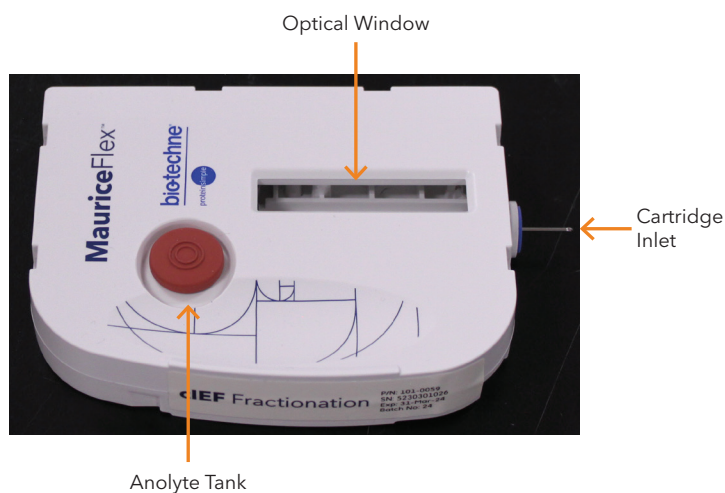
- Take the cIEF Fractionation Cartridge out of its packaging. Save the packaging, you'll need it later to store the cartridge.

NOTE: When you remove the cartridge from its packaging, make sure the cartridge inlet doesn't come in contact with any surface.



- Put the cartridge on a flat surface with the electrolyte tank facing up.

3. Remove the stopper from the Anolyte tank.



4. Add 2 mL Anolyte Solution to the H⁺ electrolyte tank (red port).

NOTE: Make sure you don't overfill the electrolyte tanks.

5. Seal the Anolyte tank with the rubber stopper. If excess liquid comes out of the tank, make sure to wipe it with a laboratory wipe.



Step 4: Install the Cartridge

1. Double check to make sure you've got analyte loaded and the tank is properly sealed with the stoppers.
2. Lift the cartridge and hold it vertically using the finger holds on either side, cartridge inlet down, with the cIEF Fractionation label facing you.
3. Gently insert it into the slot. You'll feel the alignment groove on the cartridge align inside the slot when you insert it.



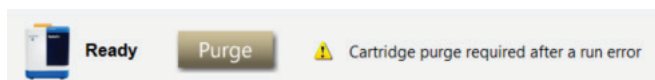
4. Continue to slide the cartridge into the slot until the locking mechanism engages. The lights on either side of the slot will change to **blue** once the cartridge is installed correctly.



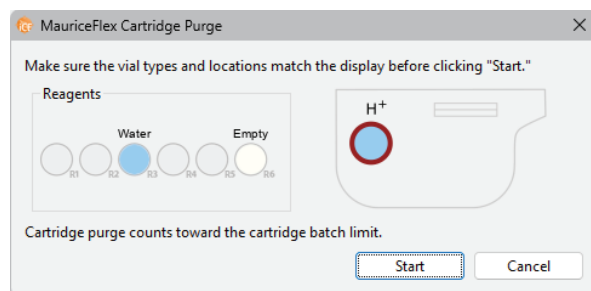
5. Close the instrument door. MauriceFlex locks it automatically.

Step 5: Check for Cartridge Alerts

1. If your cartridge was last used in a run with an error, you will need to perform a Cartridge Purge. The Maurice system will remind you if a Cartridge Purge is required.
- a. To purge the cartridge, click the brown **Purge** button in the instrument status bar.



- b. Confirm that the required reagents are loaded and that the cartridge is prepped. Then click **Start**.



- c. Once the purge is done, perform a Cartridge Post-Run Cleanup by clicking on the brown **Cleanup** button that will appear in the instrument status bar. See step 2 for more info.

NOTE: If there's only one batch left on the cartridge, you'll get a message that it can't be used after purging.

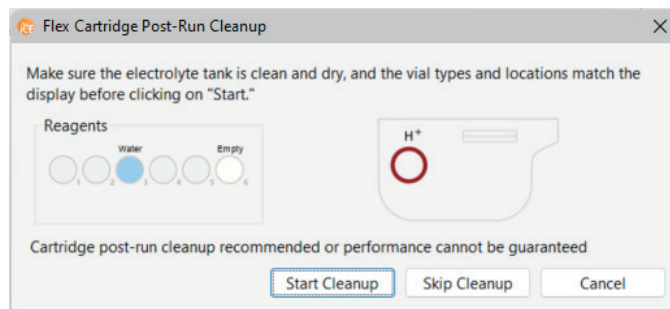
2. If you just completed a Cartridge Purge or if the cartridge you're using did not undergo a cartridge post-run cleanup after the last run, you will be asked if you want to perform a Cartridge Post-Run Cleanup.

To start the Post-Run Cleanup:

- a. Remove the cartridge from MauriceFlex and remove the Analyte from the cartridge. See "Post-batch Procedures" on page 248 for more information.
- b. Confirm there is a vial of Water (P3) and Air (P6) in place.
- c. Click the brown **Cleanup** button in the instrument status bar.



- d. Click **Start Cleanup** in the prompt that appears.

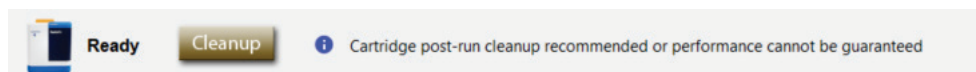


NOTE: You can also start a Post-Run Cleanup by selecting **Instrument > Cartridge Post-Run Cleanup**. When you start a Cartridge Post-Run Cleanup from the instrument menu, click **Start** in the prompt that appears to begin the cleanup.

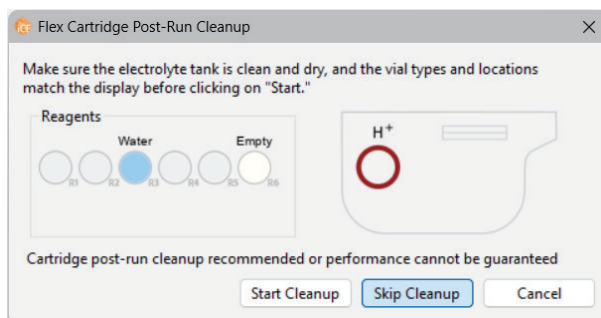
- e. Once the cartridge post-run cleanup is completed, re-prepare the cartridge and install the cartridge in MauriceFlex. See “Step 3: Prep the cIEF Fractionation Cartridge” on page 229 for more information.
- f. Click on the green **Start** button in the instrument status menu to start the batch.

To start the run without a Post-Run Cleanup:

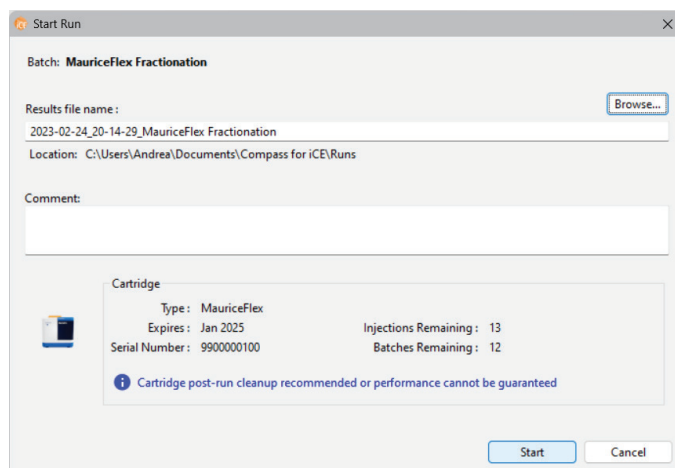
- a. Click the brown **Cleanup** button in the instrument status bar



- b. Click **Skip Cleanup** in the prompt that appears.



- c. Click **Start** in the Start Run window. The run summary and injection report will record that the cartridge did not undergo a post-run cleanup before starting the batch.

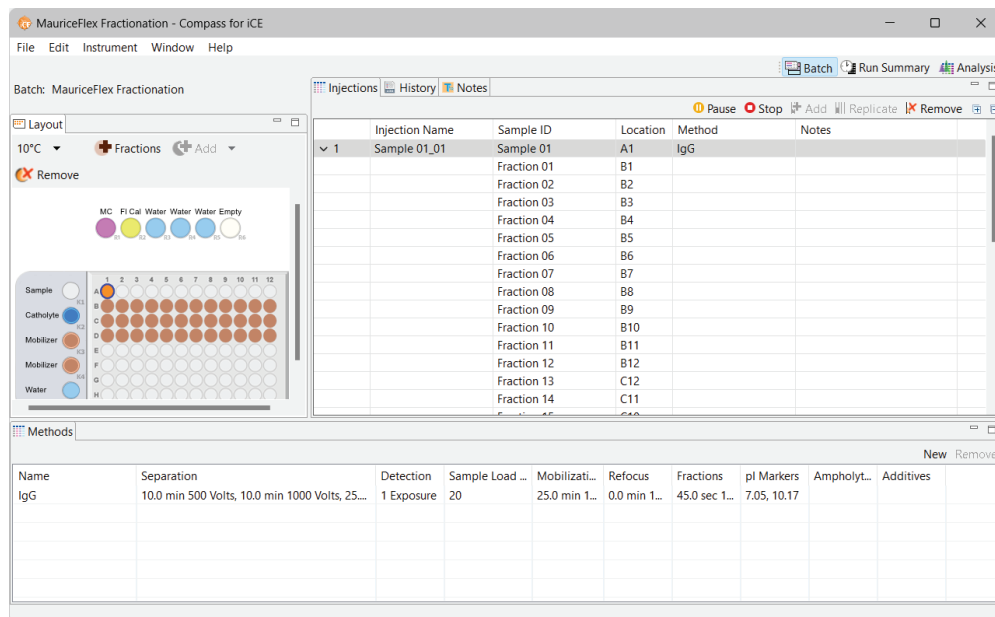


To start the run with a different cartridge:

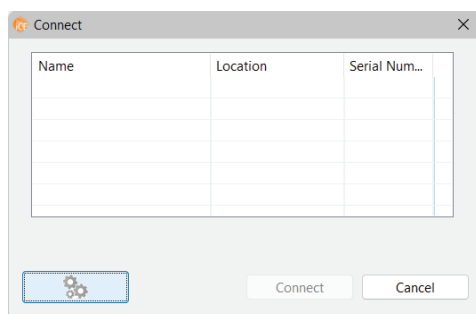
- If necessary, click **Cancel** in the cIEF Fractionation Cartridge Post-Run Cleanup window.
- Open MauriceFlex's door, remove the first cartridge from MauriceFlex and prepare a second cartridge. See "Step 3: Prep the cIEF Fractionation Cartridge" on page 229 for more information.

Step 6: Create a Batch

- Launch Compass for iCE.
- Select the **Batch** tab. This is where you'll enter sample/injection information, methods and batch parameters.

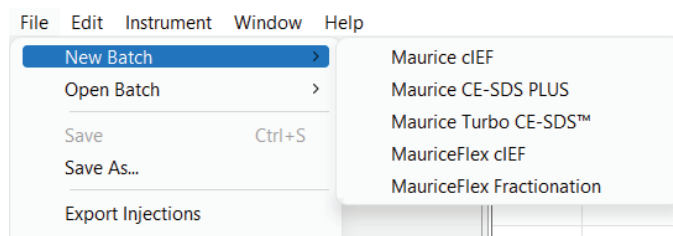


3. To create a batch, make sure MauriceFlex is connected to Compass for iCE. Select **Instrument** and click **Connect**.
 - a. If your instrument is listed, select your MauriceFlex system and click **Connect**.
 - b. If your instrument isn't listed, click on the **Settings** button and connect by typing in your instrument IP address.



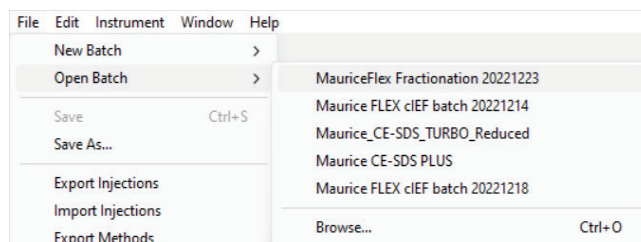
To create a new batch:

- In the main menu, select **File > New Batch > MauriceFlex Fractionation**



To use an existing batch: In the main menu, select **File > Open Batch**.

NOTE: If you're making changes to an existing batch, follow the steps in this section. Otherwise skip to "Step 7: Start the Batch" on page 244.



4. Add your sample:

To import a sample using a saved injection file:

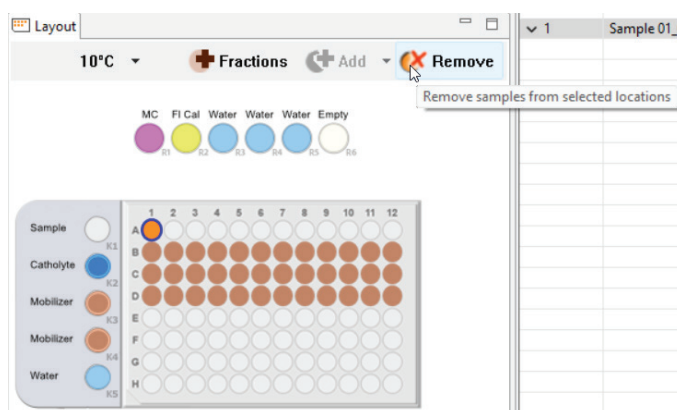
- a. Select **File** in the main menu and click **Import Injections**.

- b. Select an injection .csv file and click **OK**. The imported information will display in the Injections pane.

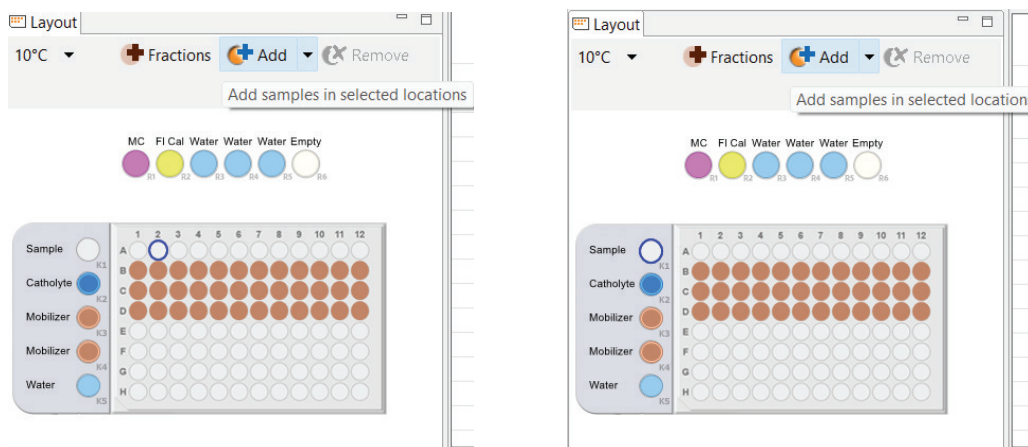
NOTE: For more information on setting up an injections file for import, see “Importing and Exporting an Injection” on page 214.

To change the sample location manually:

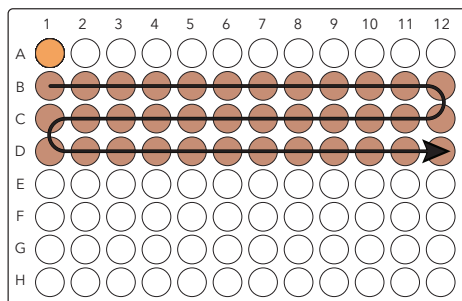
- a. In the Layout pane, highlight well A1 and click **Remove**. The sample will also be removed from the Injections table. You can also right-click on the well and click **Remove Sample**.



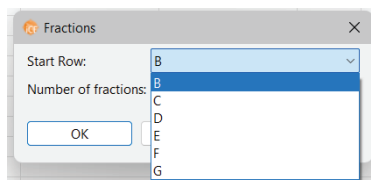
- b. Use your mouse to highlight the well your sample is located in or position K1, then click **Add**. The Injections table will automatically update.



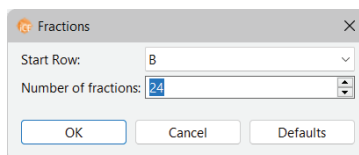
5. Assign fraction collection locations. Fraction collection starts in column 1 of the Start Row and will be collected in a serpentine pattern.



- a. In the Layout pane, click **Fractions**.
- b. A prompt will appear. In the **Start Row** pulldown, select the row to start collecting fractions.



- c. In the **Number of fractions** field, type the number of fractions to collect or use the arrows to increase or decrease the number of fractions.

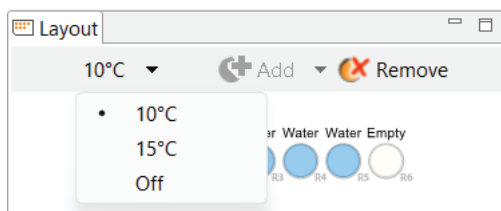


NOTE: The number entered will be highlighted in red if it is below the minimum (24) or higher than the available wells on the plate based on where you start collecting fractions.

- d. Click **Defaults** to reset fields to default settings.
- e. Click **OK**. The plate map in the Layout pane and the Injections list will automatically update.

6. The fractionation adapter is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The reagent row (R) is kept at room temperature. The reagent column (K) and sample plate are temperature controlled.



7. Enter your method parameters in the Methods pane.

NOTE: We recommend using the default method parameters. Please see the MauriceFlex cIEF Fractionation Method Development Guide for more information on method optimization.

To import a saved method:

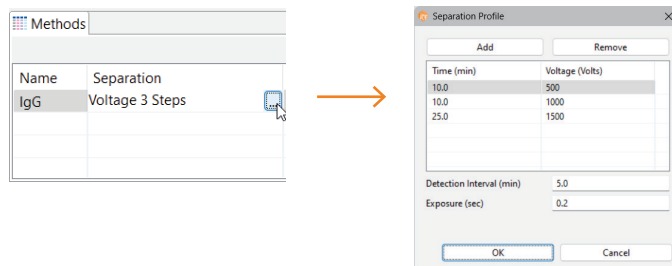
- Select **File** in the main menu and click **Import Method**.
- Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

To create a new method or make changes to an existing one:

- Click the first cell in the Name column and enter a method name.

Methods										New	Remove
Name	Separation	Detection	Sample Load ...	Mobilizati...	Refocus	Fractions	pI Markers	Ampholyt...	Additives		
10°C	10.0 min 500 Volts, 1...	1 Exposure	20	25.0 min 1...	0.0 min 150...	45.0 sec 1...	7.05, 10.17				

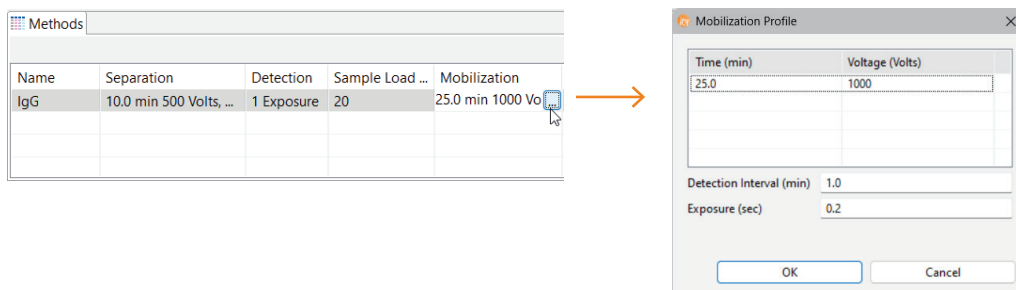
- b. Click the first cell in the Separation column, then click the selection button [...] to set the pre-focusing and focusing time (in minutes) and voltage (V). You can also set the Detection Interval (in minutes) and Exposure (in seconds) to adjust how often an image is taken using the listed Exposure time.



- c. The Detection Profile exposure setting is linked to and defined by the Separation step and cannot be changed.
- d. Click the first cell in the Sample Load(s) column and set the load time in seconds.

Methods			
Name	Separation	Detection	Sample Load (s)
IgG	10.0 min 500 Volts, 1...	1 Exposure	20

- e. Click the first cell in the Mobilization column, then click the selection button [...] to set the mobilization time (in minutes) and voltage. You can also set the Detection Interval (in minutes) and Exposure (in seconds) to adjust how often an image is taken using the listed Exposure time.



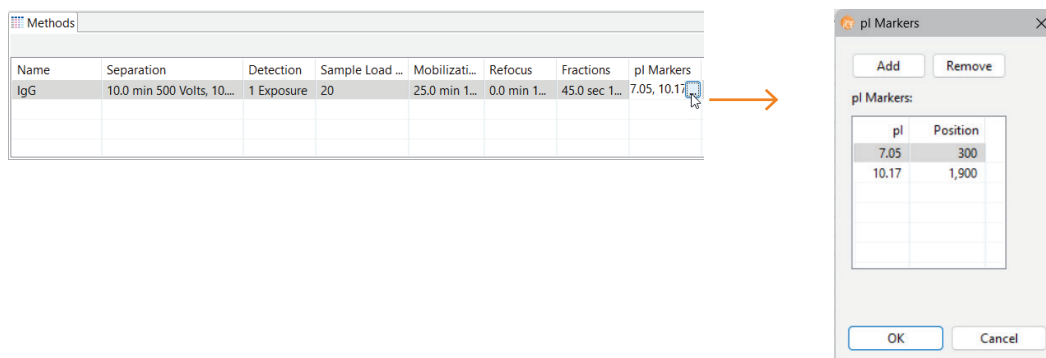
- f. Optional: Click on the first cell on the Refocus column, then click the selection button [...] to set the refocusing time (in minutes) and voltage.

NOTES:

The Compass for iCE fraction peak predictor is available but not applicable if a refocusing step is added to the method. A warning will display in the refocus profile and Start Run window and will be recorded in the run summary when then time parameter is changed.

Refer to the MauriceFlex cIEF Fractionation Method Development Guide for refocusing times and voltage setting tips.

- g. Click the first cell in the pI Markers column to select pI markers. Add new markers or remove existing ones then click **OK**.



- h. Optional: Click the first cell in the Ampholytes column and enter the ampholytes you're using.

Methods									
Name	Separation	Detection	Sample Load ...	Mobilizati...	Refocus	Fractions	pI Markers	Ampholytes	
IgG	10.0 min 500 Volts, 10.0...	1 Exposure	20	25.0 min 1...	0.0 min 1...	45.0 sec 1...	7.05, 10.17	Pharmalyte 3-10	

- i. Optional: Click the first cell in the Additives column and enter any additives you're using.

Methods									
									New Remove
Name	Separation	Detection	Sample Load ...	Mobilizati...	Refocus	Fractions	pI Markers	Ampholytes	Additives
IgG	10.0 min 500 Volts, 10.0...	1 Exposure	20	25.0 min 1...	0.0 min 1...	45.0 sec 1...	7.05, 10.17	Pharmalyte 3-10	Urea

8. You can now:

- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.

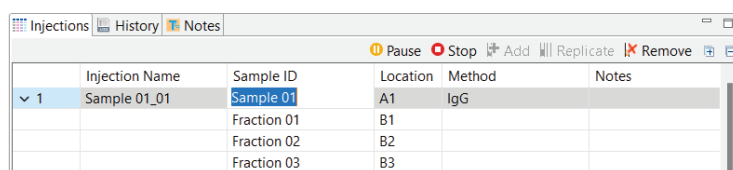
9. In the Injections pane:

- **To change the sample name:** Click the **Sample ID** cell and type a name.

NOTES:

The Sample ID can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting an Injection Name or Sample ID from other Documents” on page 213 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.



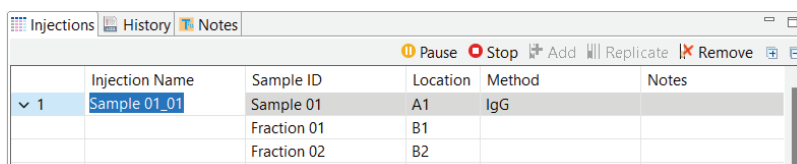
	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		
		Fraction 02	B2		
		Fraction 03	B3		

- **To change the injection name:** Click the **Injection Name** cell and type a name.

NOTES:

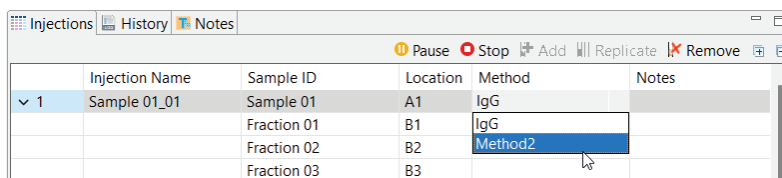
The injection name must be unique.

Changing the injection name won’t affect the sample ID.



	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1		
		Fraction 02	B2		

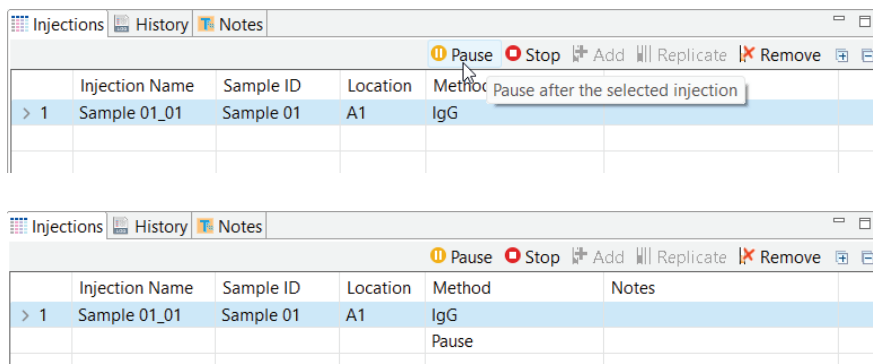
- **To assign a methods for the injection:** Click the **Method** cell and select a method from the drop down menu.



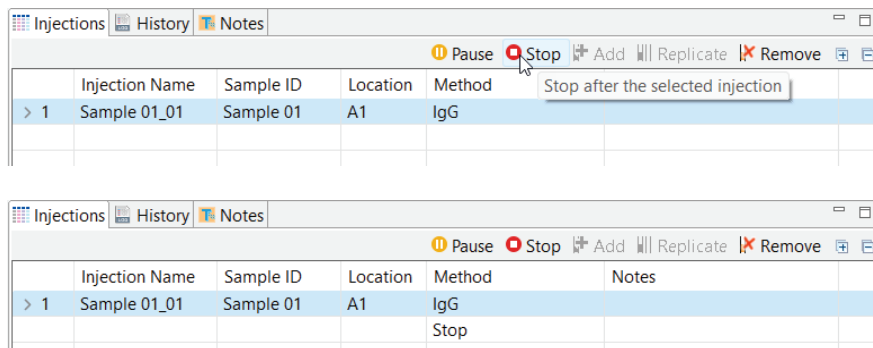
	Injection Name	Sample ID	Location	Method	Notes
▼ 1	Sample 01_01	Sample 01	A1	IgG	
		Fraction 01	B1	IgG	
		Fraction 02	B2	Method2	
		Fraction 03	B3		

Add a programmed pause or stop in the batch to automatically pause or stop after fractions have been collected (optional).

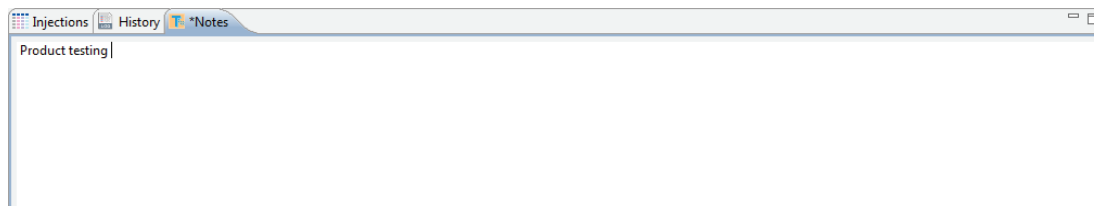
- **To program a pause:** Highlight the injection and click **Pause**. The batch will pause after fractions for that injection have been collected. To resume the batch, click **Continue** in the instrument status bar.



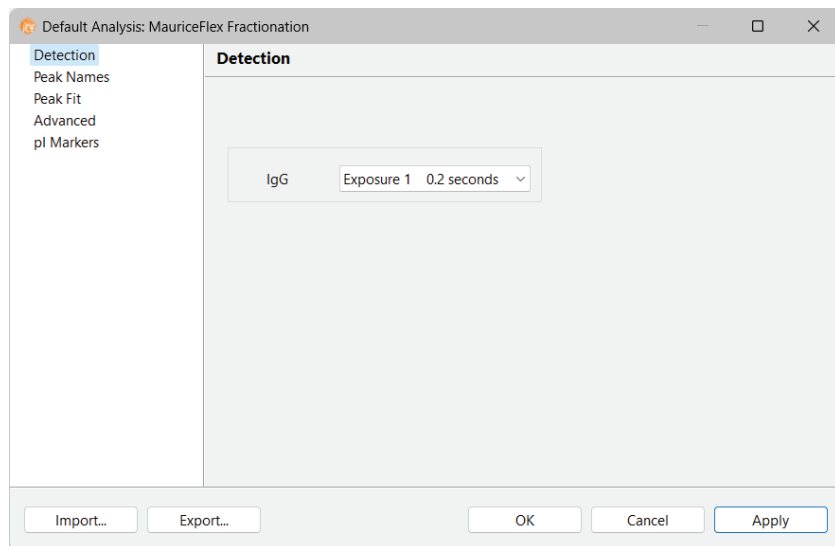
- **To program a stop:** Highlight the injection and click **Stop**. MauriceFlex will continue the batch through fraction collection, then stop the batch and perform end-run cleanup steps.



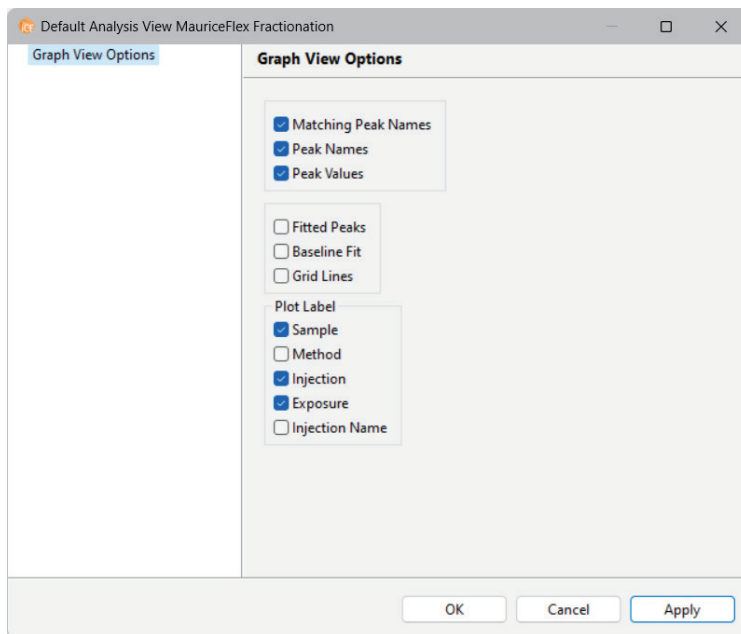
10. Click on the **Notes** pane, then click in the notes area and type any information you want to add about your batch (optional).



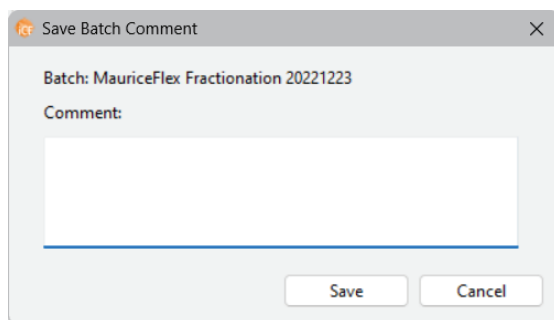
11. You can preset the parameters used to analyze data generated with the batch (optional). We recommend using the default parameters for MauriceFlex Fractionation applications, but if you want to modify parameters:
- Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



- Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to "Analysis Settings Overview" on page 620.
12. You can preset the parameters used to display the electropherogram data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button: Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



- a. Change the parameters you want to, then click **OK**. For detailed information on graph view options, please refer to “Customizing the Data Display” on page 603.
13. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



14. Enter a name for your batch then click **Save**.

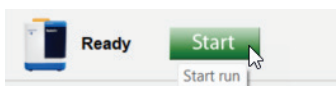
Step 7: Start the Batch

1. Make sure MauriceFlex is connected to Compass for iCE. If he isn't, select **Instrument** and click **Connect**. Then select your MauriceFlex system and click **Connect**.
2. Click on the green **Start** button to start your batch.

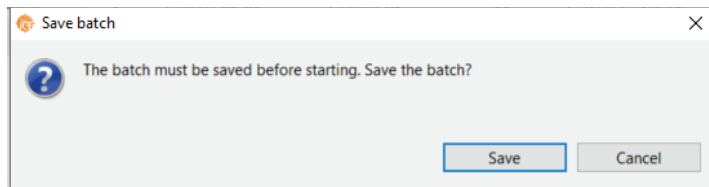
NOTES:

An alert may appear if the wrong cartridge is installed or if a cartridge post-run cleanup or cartridge purge is recommended. See “Step 5: Check for Cartridge Alerts” on page 232 for more information.

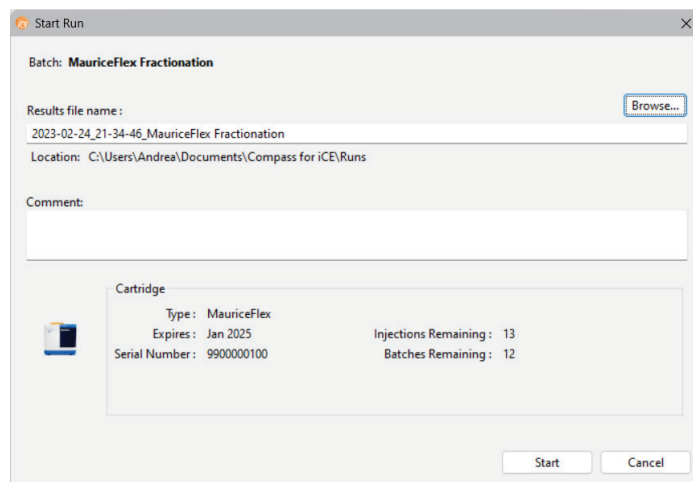
An alert may appear if you are starting a MauriceFlex Fractionation batch and have not installed the fractionation adapter. See “Adapter and Insert Alerts” on page 403 for more information.



3. If you have made any changes, you will be asked to save your batch before starting the run. Click **Save**.



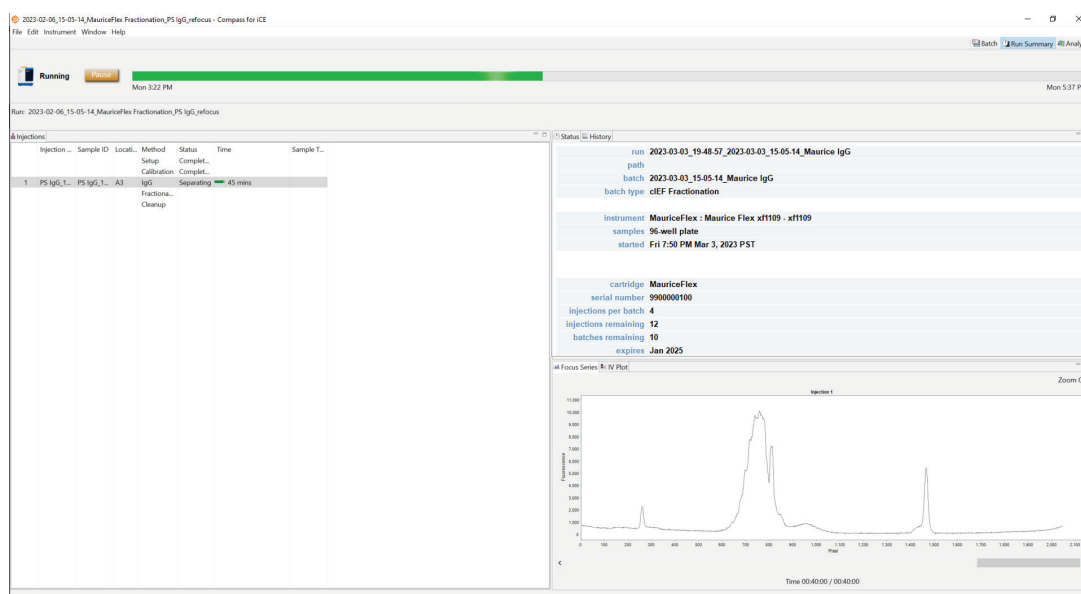
- The Start Run box shows you the number of injections and batches remaining on your cartridge. Make sure you still have enough injections left before you start the run. If not, prep a new cartridge.
- Click in the **Results File name** box if you want to change the default run file name. Otherwise just leave it as is.



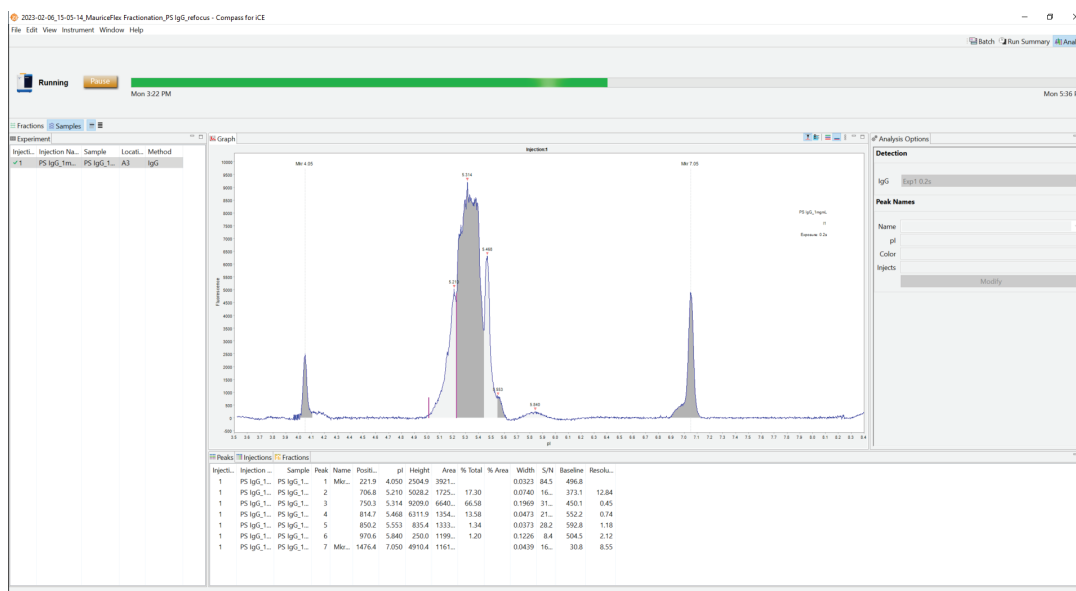
- If you don't want to save the file to the default Runs folder, click **Browse** to select a different location.
- Enter any run details you'd like in the Comments box (optional).
- Click **Start** to start the run.

NOTE: The indicator light on MauriceFlex's front panel will pulse slowly while he's running the batch.

To monitor the progress of your batch, click the **Run Summary** tab. See Chapter 16: "Run Status" for more details.



To view results, analyze results or change analysis parameters on completed injections while the batch is still running, click the **Analysis** tab. See Chapter 19: “MauriceFlex Fractionation Data Analysis” for more details.



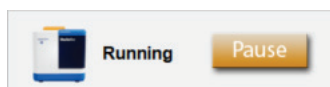
When your batch is complete, you can view electropherograms for the sample injection and mobilization in the Analysis tab, and all batch and injection details in the Run Summary tab.

NOTE: If you'd like MauriceFlex to let you know when your run is done, you can set him up to tweet you. Go to “Setting Up Maurice Systems to Send Tweets” on page 764 for more info.

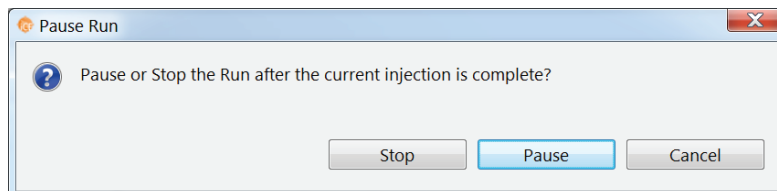
Editing a Running Batch

You can pause a batch after it's started to make updates or add samples and reagents.

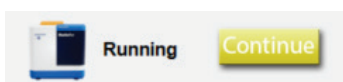
1. Select **Instrument** > **Pause** or select the **Pause** button.



NOTE: MauriceFlex can only be paused during a run after the calibration step or after the injection has completed.

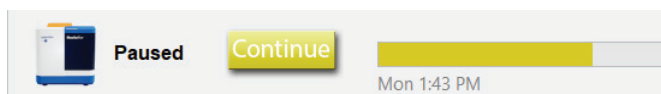
2. Click **Pause** in the pop-up window.

The Pause button will change to Continue but MauriceFlex's status will stay at Running until Compass for iCE completes the current injection.



- You can now make these edits to the batch:
 - Change the Sample Name
 - Change the Injection Name
 - Modify the injection if it isn't greyed out
 - Create a new method
 - Update the method.
3. When MauriceFlex completes the current injection, his status will change to Paused and the progress bar will turn yellow.

NOTE: MauriceFlex's status light will change from a pulsing blue to a blue pulse with magenta flashes which means he's paused the run.



The status in the Injection pane also changes to Paused and the pause duration updates live. You can now open the instrument door.

IMPORTANT

The door must be closed within 15 minutes of opening it to avoid potential assay performance issues.

NOTE: If the door is open for more than 30 seconds, a white light near the cartridge will start to flash and a warning message displays in Compass for iCE until the door is closed.

4. To continue the batch, click **Continue**. Changes you made to the batch are saved and MauriceFlex will continue with the batch.

Post-batch Procedures

When the batch is done:

1. Open MauriceFlex's door. The lights on either side of the cartridge slot will be **orange** as MauriceFlex will have already disengaged the cartridge.
2. Remove and save the 96-well plate. Collected fractions will be used to for further analysis.

IMPORTANT:

We recommend that you seal the plate to prevent evaporation if the fractions will not be used immediately.

NOTE: Check out the MauriceFlex cIEF Fractionation Method Development guide for more info on how confirm the identity and purity of peaks collected in your fractions using a Maurice cIEF batch.

3. If the sample was prepared in a vial, remove your sample. Leave the Water (P3) and Air (P6) vials in place if your cartridge still has injections left since they will be needed for the post-run cleanup step. Remove the remaining reagent vials.

NOTE: You can remove reagent vials from the MauriceFlex fractionation adapter when the locking mechanism is pulled to the right (row R) or pulled down (column K). Release the mechanism to lock remaining vials in place before starting the Post-Run Cleanup.

4. Gently pull out the cartridge making sure the cartridge inlet doesn't come into contact with any surface.



If you're at 15 injections, you've reached the limit of supported injections for the cartridge. To dispose of a finished cartridge, put it in its original packing and discard it per your institution's safety and waste disposal guidelines. Discard any remaining reagent vials.

NOTE: Disposal of cartridges, sample plates and vials depends on the samples you ran. If you aren't sure what your sample origins are, we recommend you dispose of everything in biohazard waste.

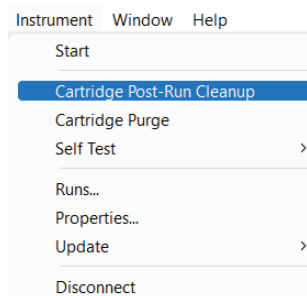
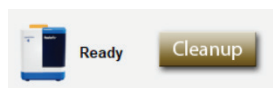
If the cartridge will be used again, clean up and store the cartridge:

- a. Put the cartridge on a flat surface with its electrolyte tank facing up.
- b. Remove the stoppers from the Anolyte tank.
- c. Using an electrolyte pipette or low vacuum, aspirate the solutions from the tank.
- d. Fill the tank with 3 mL DI water, then aspirate it out. Repeat this rinse 2 more times.

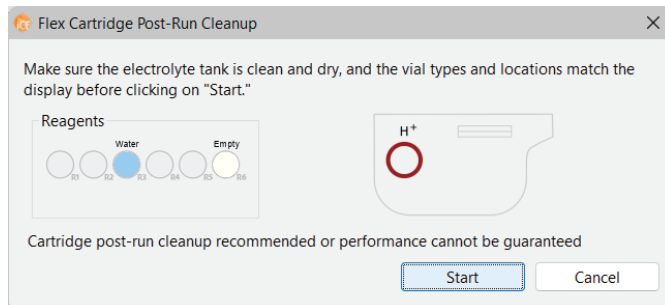
NOTE: Make sure not to get any liquid on the cartridge's optical window.



- e. Aspirate all the remaining liquid and make sure that the tank is dry.
- f. Put the stopper back on the tank and install the cartridge in MauriceFlex.
- g. Verify there is at least 2 mL of DI water in the Water (P3) vial and air in the empty (P6) vial.
- h. Click on the brown **Cleanup** button in the Instrument Status Bar or select **Instrument** and **Cartridge Post-Run Cleanup** in the Compass main menu.



- i. You'll get the following message. Click **Start**. It'll only take 6 minutes.



- j. Open MauriceFlex's door. Gently pull out the cartridge making sure the cartridge inlet doesn't come in contact with any surface.
- k. Remove the reagent vials.
- l. Leave the stopper off to allow the tank to air dry.
- m. Put the cartridge and stopper back in the protective packaging and store at room temperature.

!WARNING! SHARPS HAZARD

The capillary inlet of the cartridge may present a potential sharps hazard. Dispose of used cartridges according to your organizations health and safety regulations.



!WARNING! BIOHAZARD

Cartridges, sample plates and vials should be handled by procedures recommended in the CDC/NIH manual: Biosafety in Microbiological and Biomedical Laboratories (BMBL). The manual is available from the U.S. Government Printing Office or online at <http://www.cdc.gov/biosafety/publications/bmbl5/>.

Depending on the samples used, the cartridges, plates and vials may constitute a chemical or a biohazard. Dispose of the cartridges, plates and vials in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations. Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste content before you store, handle, or dispose of chemical waste.

Checking Your Data

Compass for iCE detects your sample protein, fraction and pI marker peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

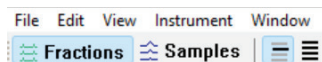
Check your data in the Analysis screen.

- When you select **Fractions** in the View bar, you'll see:
 - The focused injected sample (I1) - where you can make sure the pI markers for your sample are identified correctly.
 - Mobilization electropherograms (T0, T1, T2, etc) that show the sample moving off of the capillary- where you can confirm all the peaks in the mobilization electropherograms are correctly identified.
- When you select **Samples** in the view bar, you'll see:
 - The focused, injection sample (I1) - where you can make sure peaks in your sample injection are identified correctly and where you can name the peaks. Named peaks will automatically be updated in the mobilization electropherogram in the Fractions view bar.

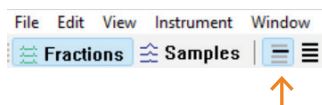
Step 1: Check the pI Markers for the Injected Sample

To make sure your pI markers are identified correctly:

1. Go to the **Analysis** screen.
2. Click **Fractions** in the View bar.

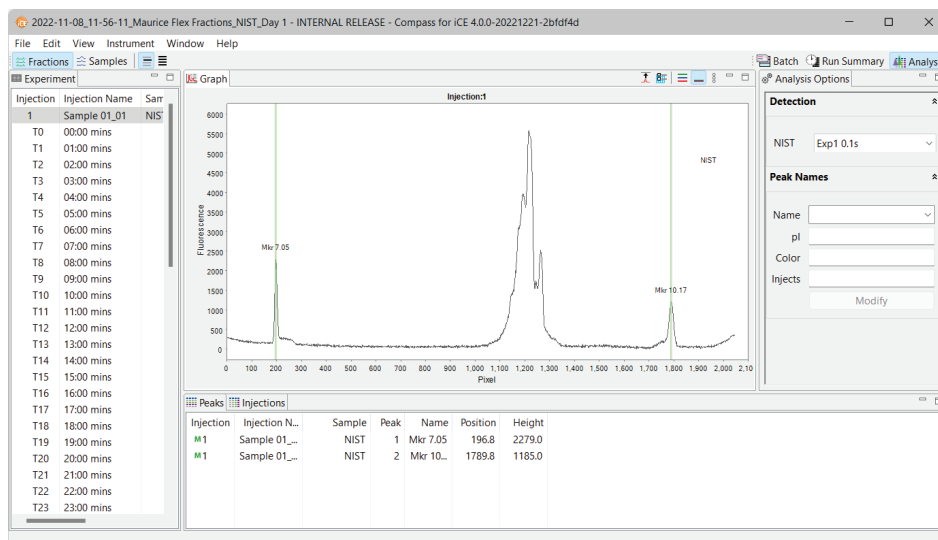


3. Click the **View Selected** icon in the View bar.



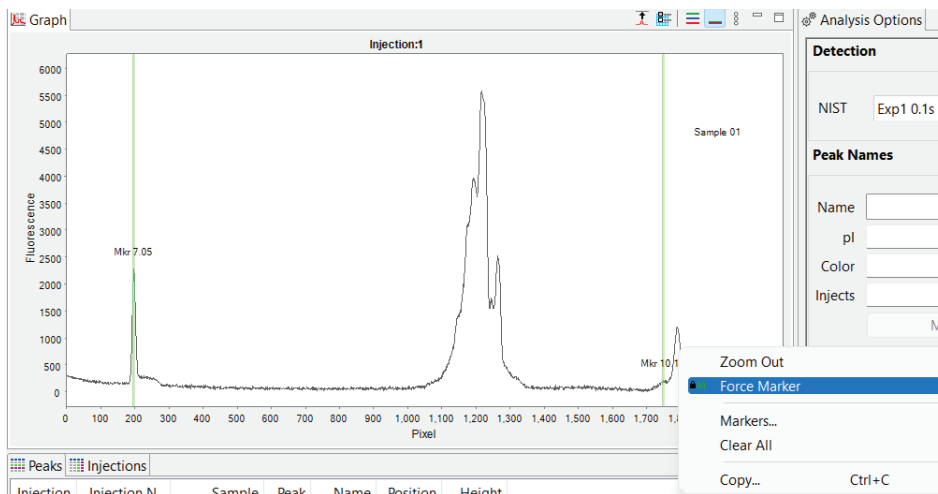
4. Click **Injection 1** in the Experiment pane.

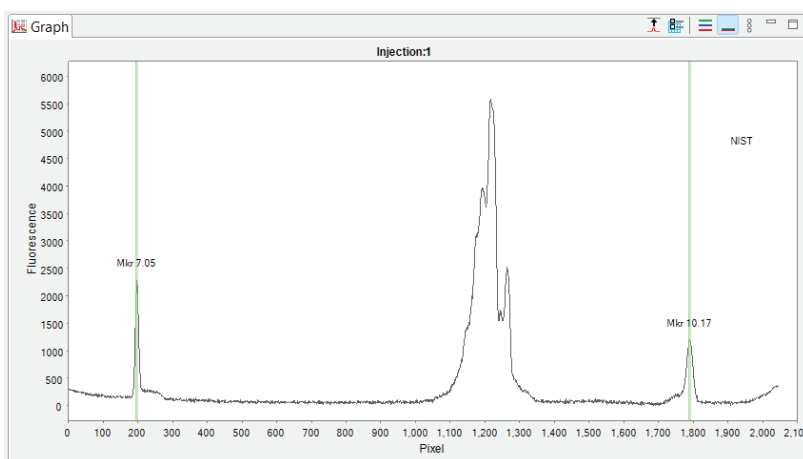
5. Check that your pI markers in the electropherogram have been correctly identified. Each marker peak will have a green vertical line running through it and be labeled Mkr. They're also identified with an **M** in the Peaks table.



6. If your pI markers aren't identified correctly, here's how to manually correct them:

To set an unidentified peak as a pI marker: Right-click the peak in the electropherogram or Peaks table and select **Force Marker**. Compass will assign that peak as a pI marker, and correctly reassign the remaining pI marker peaks.



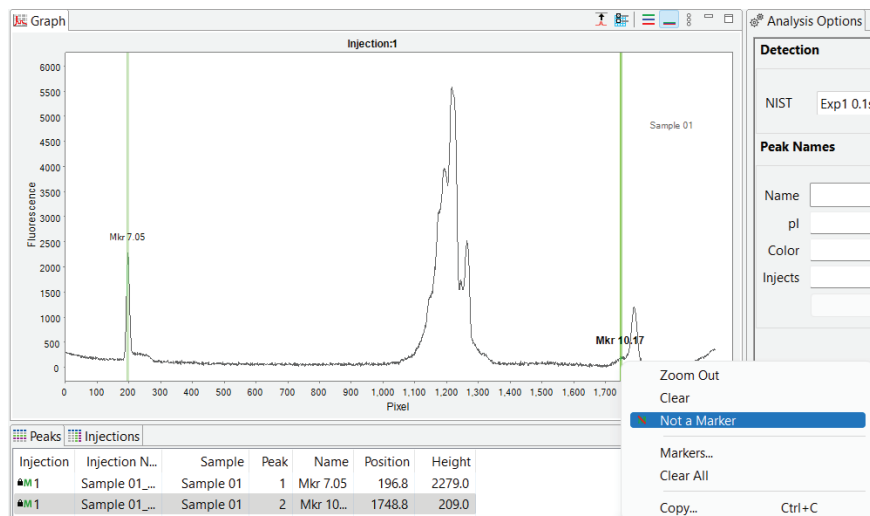


A lock icon indicating the pI marker was set manually will display next to the peak in the Peaks table and a check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

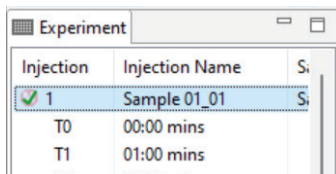
Peaks			Injections		
Injection	Injection Name	Sample	Injection	Injection Name	Sample
M 1	Sample 01_01	Sample 01	✓ 1	Sample 01_01	S
M 1	Sample 01_01	Sample 01	T0	00:00 mins	
			T1	01:00 mins	

NOTE: To remove pI marker peak assignments that were made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**. To clear all the manual settings for the injection, click **Clear All**.

If an incorrect peak is identified as a pI marker: Right-click the peak in the electropherogram or Peaks table and select **Not a Marker**. Compass should correctly reassign the remaining peaks as pI markers and update the Peaks table.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.



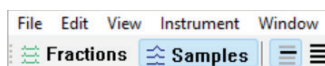
Step 2: Checking Sample Peaks

All detected peaks in your injected sample will be labeled automatically with the calculated protein pI.

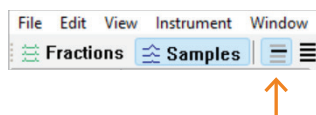
NOTE: The reported protein pI in Compass may vary slightly from predicted pIs based on sample, buffer, and method conditions.

To make sure peaks in your sample injection are identified correctly:

1. Click **Samples** in the View bar.



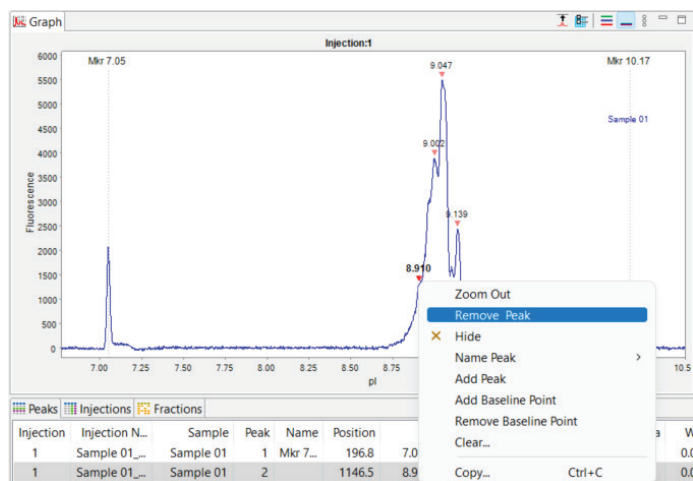
2. Click the **View Selected** icon in the View bar.



3. Click **Injection 1** in the Experiment pane.

4. If your sample peaks aren't identified correctly, here's how to manually correct them:

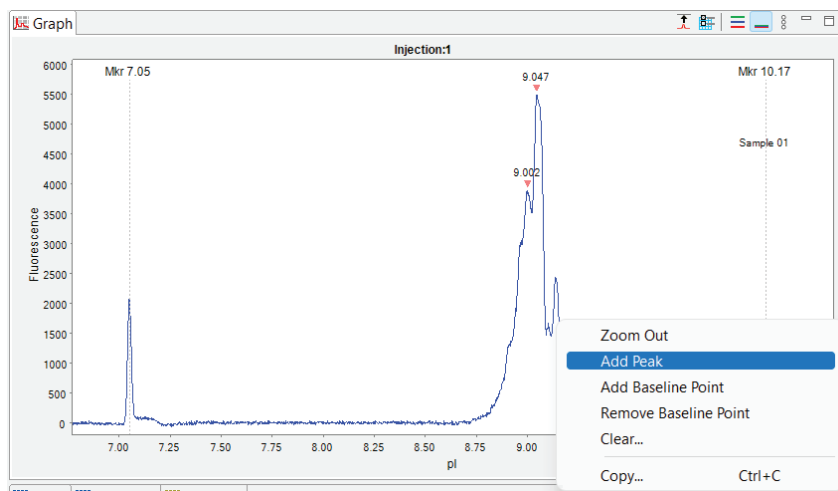
If a peak is incorrectly identified as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass will no longer identify it as a sample peak in the electropherogram and the peak data will be removed in the results table.



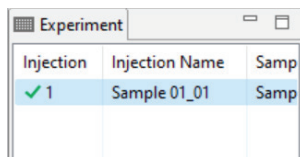
A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment		
Injection	Injection Name	Samp
✓ 1	Sample 01_01	Samp

To identify a peak as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.



Injection	Injection Name	Sample
✓ 1	Sample 01_01	Sample

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear**.

- Repeat the previous steps to make sure all sample peaks are correctly identified.

Step 3: Assigning Sample Peak Names

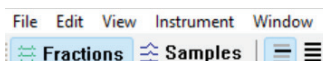
Compass can also optionally identify and name sample peaks in the focused injection sample using user-specified peak name settings. For more information on how to do this, see “Manual Peak Integration” on page 637.

Step 4: Checking Peaks in the Mobilization Electropherograms

All detected peaks in the mobilization electropherograms will be labeled automatically with either the pixel position or peak name as defined in Peak Names.

To make sure peaks in the mobilization electropherograms are identified correctly:

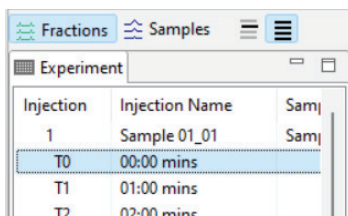
- Click **Fractions** in the View bar.



- Click the **View Selected** icon in the View bar.



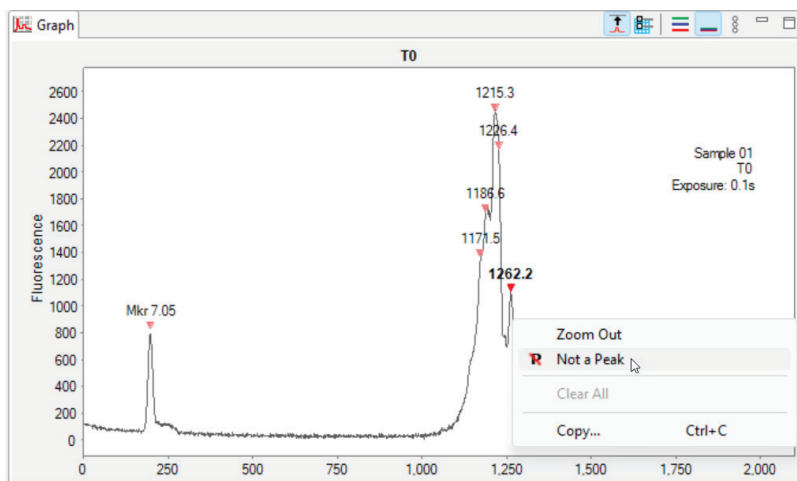
- Click **T0** in the Experiment pane.



Injection	Injection Name	Sample
1	Sample 01_01	Sample
T0	00:00 mins	
T1	01:00 mins	
T2	02:00 mins	

4. If the peaks aren't identified correctly, here's how to manually correct them:

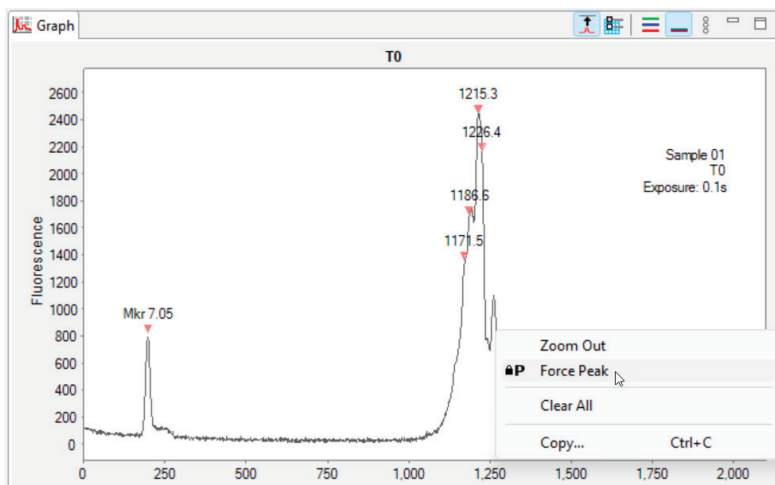
If a peak is incorrectly identified: Right-click the peak in the electropherogram or Peaks table and select **Not a Peak**. Compass will no longer identify it as a peak in the mobilization electropherogram and the peak data will be removed in the results table.



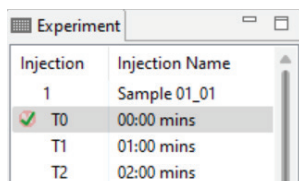
A check mark will appear next to the mobilization electropherogram in the Experiment pane to indicate a manual correction was made.

Experiment	
Injection	Injection Name
1	Sample 01_01
✓ T0	00:00 mins
T1	01:00 mins
T2	02:00 mins

To identify a peak: Right-click the peak in the electropherogram and select **Force Peak**. Compass will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A check mark will appear next to the mobilization electropherogram in the Experiment pane to indicate a manual correction was made.



Injection	Injection Name
1	Sample 01_01
✓ T0	00:00 mins
T1	01:00 mins
T2	02:00 mins

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear**.

5. Repeat the previous steps to make sure all peaks are correctly identified.

Step 5: Assigning Peak Names in the Mobilization Electropherograms

Existing peak names assigned to the focused injection sample are automatically updated in the mobilization electropherogram and can be manually adjusted. For more information on how to do this, see page 601.

Chapter 11:

CE-SDS PLUS Batches

Chapter Overview

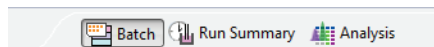
- Overview
- Batch Screen Overview
- Opening a Batch
- Creating a New Batch
- Viewing Replicate Injections
- Batch History
- Making Changes to a Batch
- Viewing and Editing Batches in Completed Runs
- Copying and Pasting Injection Names or Sample IDs from other Documents
- Importing and Exporting Injections
- Importing and Exporting Methods
- Batch Reports

Overview

Maurice CE-SDS PLUS batches can be run on Maurice, Maurice S., or MauriceFlex systems using a CE-SDS PLUS cartridge.

Batch Screen Overview

You can use the Batch screen to create, view, and edit batches. To get to this screen, click the **Batch** screen tab:



Batch Screen Panes

The Batch screen has five panes:

- **Layout** - Displays a map of either the 96-well plate or 48-vial tray of batch sample locations.
- **Injections** - Lists the injections, sample ID, sample locations and methods that Maurice or Maurice S. will execute for each sample in the batch.
- **History** - Lists all batch file events from initial creation to the most current update.
- **Notes** - Lets you enter specific information about your batch.
- **Methods** - Lets you create methods and enter method parameters used in the batch.

The screenshot shows the Maurice CE-SDS PLUS Batch screen. The top menu bar includes File, Edit, Instrument, Window, and Help. The main window has tabs for Batch, Run Summary, and Analysis. The Batch tab is active, showing a sub-menu with Injections, History, and Notes. The Layout pane on the left shows a 96-well plate map with columns 1-8 and rows A-F. The Injections pane shows a list of 15 injections with columns for Injection Name, Sample ID, Location, Method, and Notes. The Methods pane at the bottom shows a table with columns for Name, Sample Load, and Separation.

Injection Name	Sample ID	Location	Method	Notes
1 IgG System Control_01	Sample 01	A1	Method 1	
2 Control Ladder_02	Sample 02	A2	Method 2	
3 Test Ladder_03	Sample 03	A3	Method 2	
4 IS - Alpha_04	Sample 04	B1	Method 1	
5 IS - Alpha_05	Sample 08	B2	Method 2	
6 IS - Beta_06	Sample 05	B3	Method 1	
7 IS - Beta_07	Sample 06	B4	Method 2	
8 IS - Gamma_08	Sample 07	B5	Method 1	
9 IS - Gamma_09	Sample 09	B6	Method 2	
10 IgG System Control_10	Sample 01	A1	Method 1	
11 Control Ladder_11	Sample 02	A2	Method 2	
12 Test Ladder_12	Sample 03	A3	Method 2	
13 IS - Alpha_13	Sample 04	B1	Method 1	
14 IS - Alpha_14	Sample 08	B2	Method 2	
15 IS - Beta_15	Sample 05	B3	Method 1	

Name	Sample Load	Separation
Method 1	20 sec 4600 Volts	0.1 min 1150 Volts, 0.1 min 3450 Volts, 25.0 ...
Method 2	20 sec 4600 Volts	0.1 min 1150 Volts, 0.1 min 3450 Volts, 30.0 ...

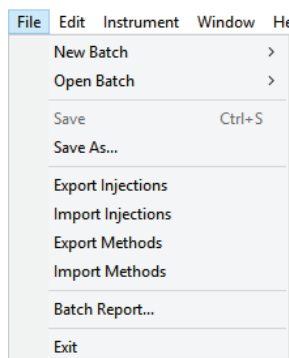
Software Menus Active in the Batch Screen

These main menu items are active in the Batch screen:

- File
- Edit
- Instrument (when the software is connected to Maurice, Maurice S. or MauriceFlex)
- Window
- Help

File Menu

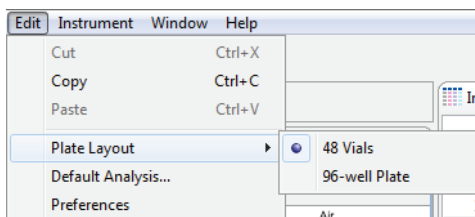
These File menu options are active:



- **New Batch** - Creates a new batch from a starter template.
- **Open Batch** - Opens an existing batch.
- **Save/Save As** - Saves the open batch.
- **Export Injections** - Exports injections from the current batch as a .csv file.
- **Import Injections** - Imports injections into the current batch from a .csv file.
- **Export Methods** - Exports method(s) from the current batch as separate files.
- **Import Methods** - Imports saved method(s) into the current batch.
- **Batch Report** - Exports a table of sample and method details for each injection in the batch as a PDF file.
- **Exit** - Closes Compass for iCE.

Edit Menu

The following Edit menu options are active in the Batch screen:



- **Cut** - Cuts the information currently selected.
- **Copy** - Copies the information currently selected.
- **Paste** - Pastes the copied information.

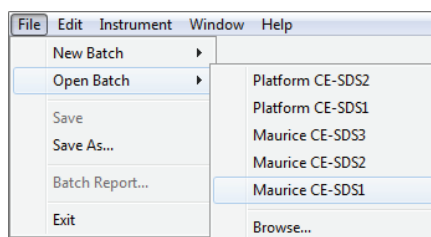
NOTE: Cut, Copy and Paste work differently depending on the pane you're in. How and when you can use them is described throughout the chapter.

- **Plate Layout** - Lets you select between a 96-well plate or a 48-vial tray to run samples in.
- **Default Analysis** - Displays the default settings that will be used to analyze the data generated with your batch.
- **Default Analysis View** - Displays the default settings that will be used to view the data generated with your batch.
- **Preferences** - Lets you set and save your preferences for Access Control, data export, graph colors, grouped data and Twitter settings. See Chapter 21, "Setting Your Preferences" for more information.

Opening a Batch

To open an existing batch:

1. Select **File** in the main menu and click **Open Batch**.



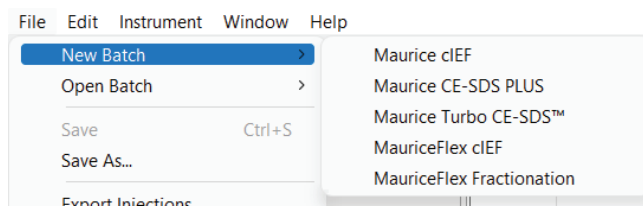
2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file. Select one of those or click **Browse** to open the Batches folder and select a different one.
3. To make changes to the batch, see the steps in “Creating a New Batch”. When you’re done, select **File** from the main menu and click **Save**.

Creating a New Batch

To create a new batch, we recommend that you use one of the template methods and make any necessary modifications from there.

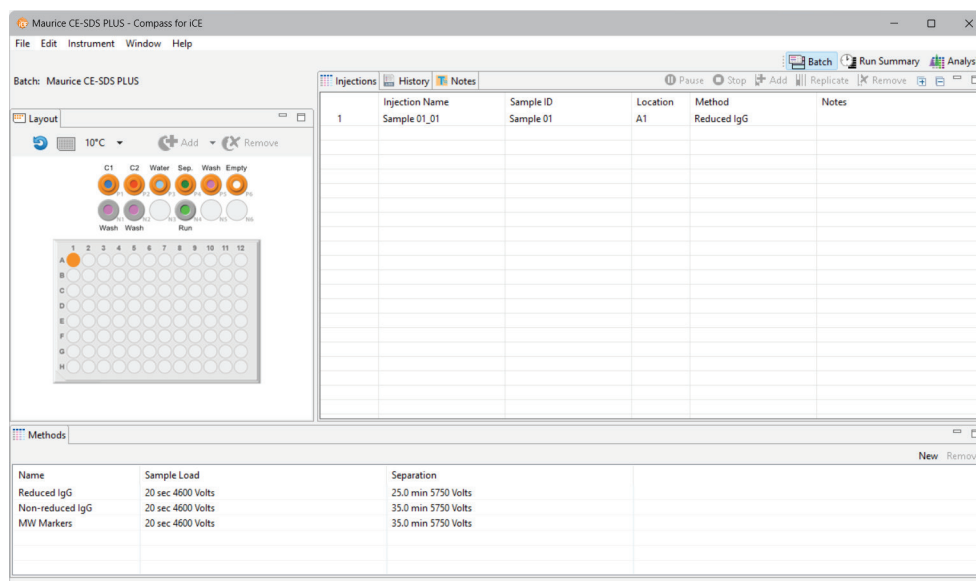
Step 1 - Open a Template Batch

1. Select **File** in the main menu and click **New Batch**:



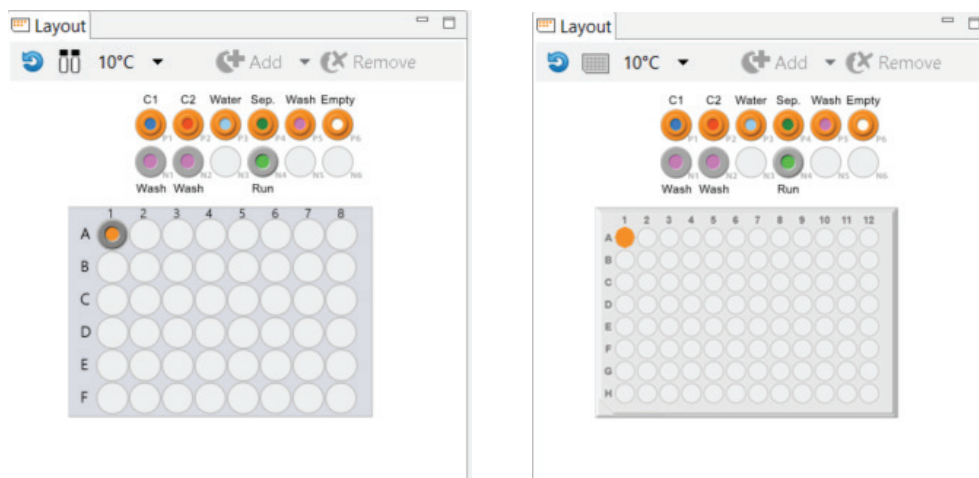
NOTE: Only template batches specific to your Maurice will be available when you are connected to the instrument. Template batches will not appear in the menu if that application type cannot be performed on your instrument.

2. Select **Maurice CE-SDS PLUS**. A batch using the default method will display.



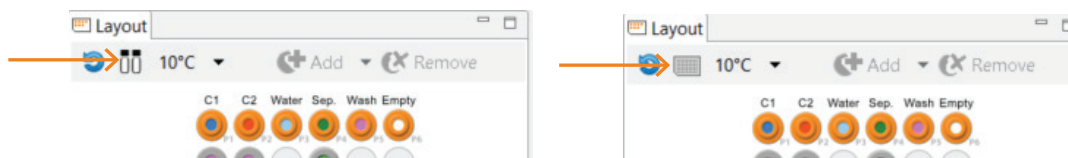
Step 2 - Assign Your Samples

The Layout pane shows the default locations of where batch reagent should be placed in your Maurice system's samples and reagents platform, and a map of either the 96-well plate or the 48-vial tray used for samples.



The same reagent locations are used for every batch:

- P1 - Conditioning Solution 1 with **orange pressure cap**
 - P2 - Conditioning Solution 2 with **orange pressure cap**
 - P3 - DI water with **orange pressure cap**
 - P4 - Separation Matrix with **orange pressure cap**
 - P5 - Wash Solution vial with **orange pressure cap**
 - P6 - Empty vial (air) with **orange pressure cap**
 - N1 - Wash Solution vial with **clear screw cap**
 - N2 - Wash Solution vial with **clear screw cap**
 - N4 - Running Buffer - Bottom with **clear screw cap**
1. To assign samples, select a 96-well plate or 48 vials depending on what you're running. Clicking on the vial/plate icon toggles between formats.



2. Select your samples:

To import samples using a saved injections file:

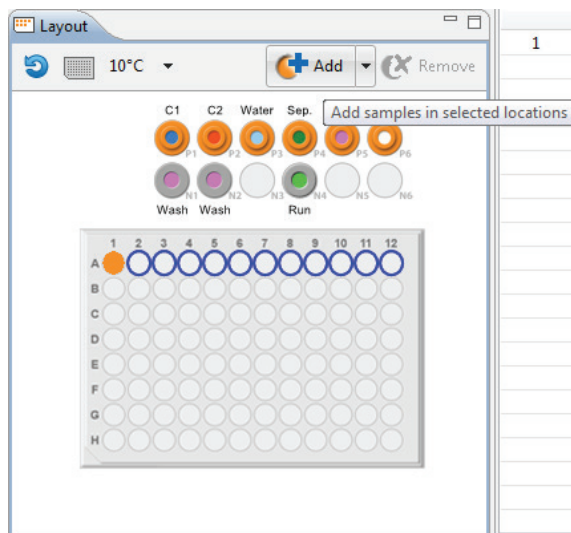
- Select **File** in the main menu and click **Import Injections**.
- Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.
- Skip to step 3 page 266.

NOTE: For more information on setting up an injections file for import, see “Importing and Exporting Injections” on page 281.

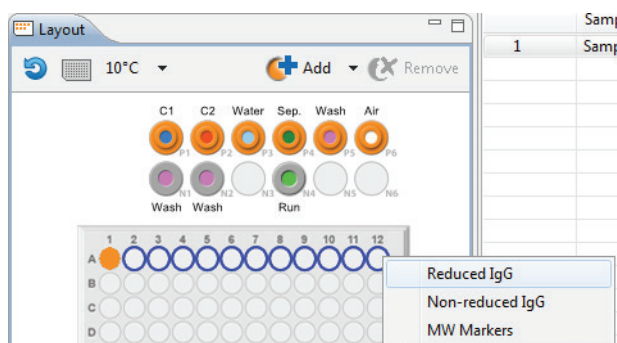
To select samples manually:

- **Add samples and select methods later:** Use your mouse to highlight the well or vial positions your samples are located in, then click **Add**. For this example we’re using a 96-well plate.

NOTE: The template batch automatically adds a sample in well or vial A1 by default, but if you don’t have a sample in this position you can remove it after you’ve added new positions for your samples.



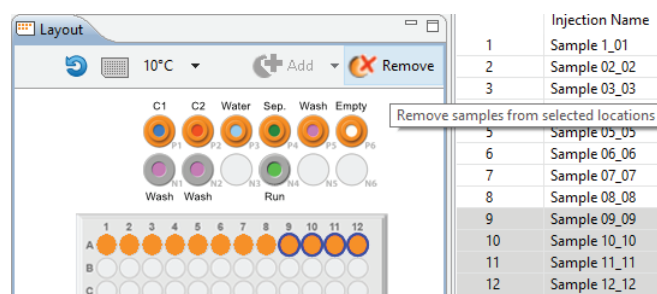
- **Add samples with preassigned methods:** Highlight the well or vial positions your samples are located in, then right-click and select a method.



Either option of adding samples populates the Injections table:

Injections					
History Notes					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Sample 1	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

To remove samples, just highlight the ones you want to remove and click **Remove**. They'll also be removed in the Injections table. You can also right-click well(s) or vial(s) in the Layout pane and click **Remove Sample(s)**.

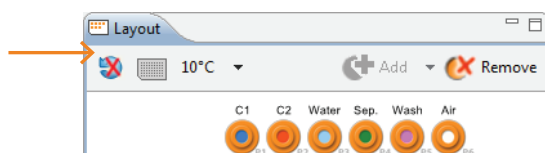


3. Compass for iCE can monitor the current during a separation for you, stop it if an abnormal current profile is detected and reinject the sample for you automatically. Reinjections are enabled by default, but if you don't want to use this feature just click the reinject icon to toggle it off.

NOTES:

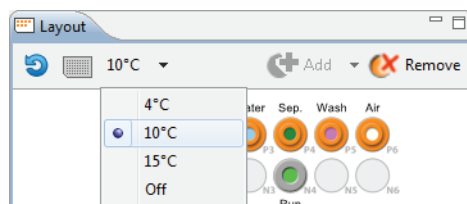
If a sample was reinjected, the injection row will be flagged in the Injections pane in the Run Summary screen. See “Injection Flags” on page 414 for more info.

A maximum of 10 reinjections are allowed per batch.



4. The sample platform is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The two reagent rows are kept at room temperature.



Step 3 - Assign Your Method Parameters

NOTE: There are three default methods. We recommend using the default method parameters for the listed sample types. Please see the [Maurice CE-SDS Application Guide](#) for more information on method optimization.

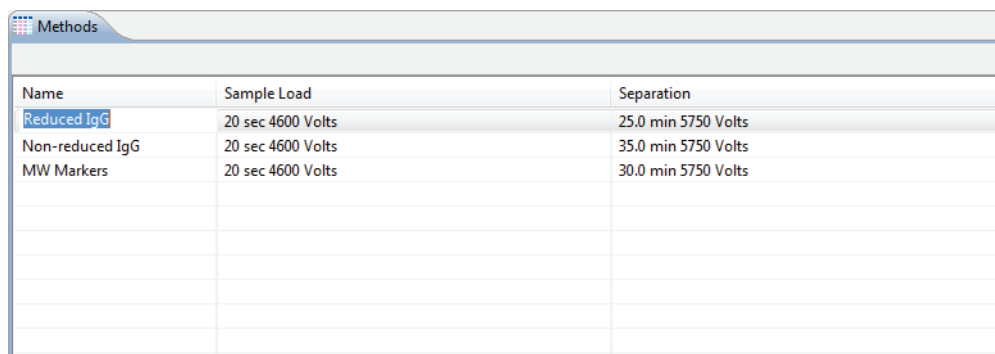
The Methods pane lists all of the methods used in a batch. You can add and remove methods here and change method parameters.

To import a saved method:

1. Select **File** in the main menu and click **Import Method**.
2. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

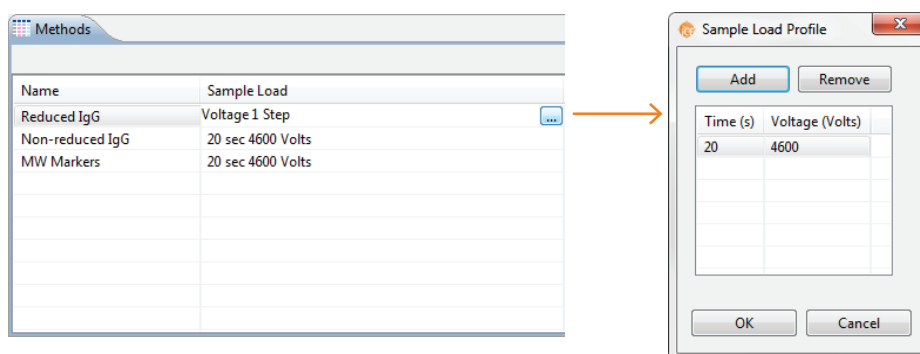
To create a new method or make changes to an existing one:

1. In the Methods table, click the first cell in the Name column and enter a new method name if needed.



Name	Sample Load	Separation
Reduced IgG	20 sec 4600 Volts	25.0 min 5750 Volts
Non-reduced IgG	20 sec 4600 Volts	35.0 min 5750 Volts
MW Markers	20 sec 4600 Volts	30.0 min 5750 Volts

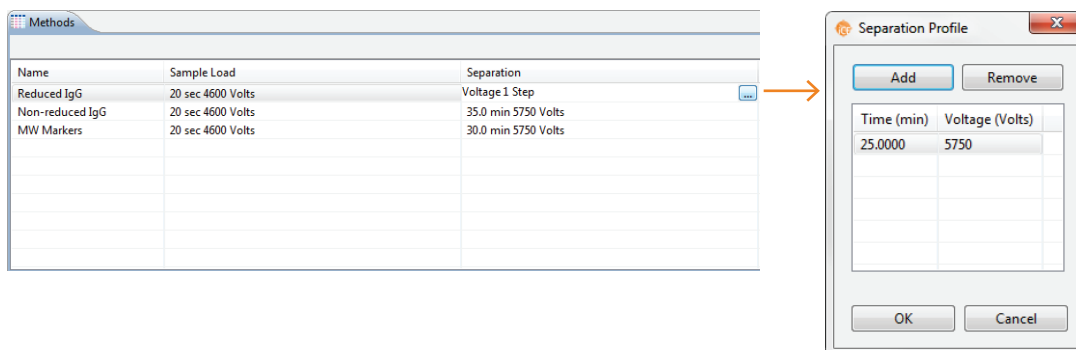
2. Click the first cell in the Sample Load column, then click the selection button [...] to set your sample load profile time (in seconds) and voltage.



- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a load time (in seconds) and voltage value.
- **To remove a profile step:** Select the row you want to remove and click **Remove**.

- Click the first cell in the Separation column the selection button [...] to set your separation profile parameters (in minutes) and voltage.

NOTE: Run your reduced IgG samples and IgG Standard for 25 minutes and the CE-SDS MW Markers for 35 minutes. Run your non-reduced IgG samples and IgG Standard for 35 minutes. The default separation voltage for all sample types is 5750 volts.



- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a load time (in seconds) and voltage value.
- **To remove a profile step:** Select the row you want to remove and click **Remove**.

- You can now:

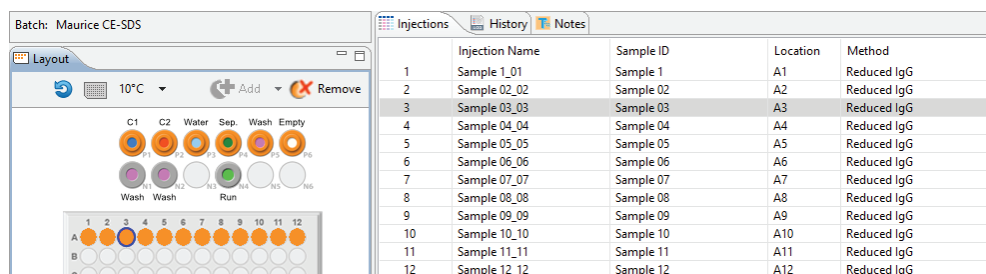
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.
- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.

Step 4 - Set Up Your Injections

The Injection pane lists all sample injections for the batch. Any samples that were added in Step 2 - Assign Your Samples are automatically added to this list.

NOTE: If you imported injections in Step 2, you can update the injections table and assign methods as needed using the information that follows. Otherwise skip to “Step 5 - Add Programmed Pauses and Stops (Optional)” on page 272.

Selecting an injection in the table also selects the sample the injection will be made from in the plate or vial map in the Layout pane and vice-versa.



1. To add sample names, click the **Sample ID** cell for the injection and type a name.

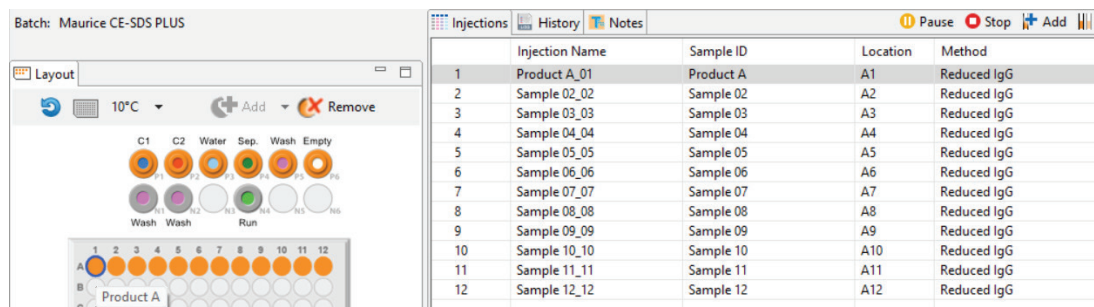
NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 279 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

Injections History Notes					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

The sample name also displays when you hover the mouse over the sample in the plate or vial map:



- Injection names are automatically set to 'Sample ID_injection number'. To change the name, click the **Injection Name** cell for the injection and type a name. Each injection name must be unique. If they aren't when you save the batch, Compass for iCE highlights any non-unique injection names so you can change them.

	Injection Name	Sample ID	Location	Method	Notes
1	Injection	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

NOTE: Changing the injection name won't affect the sample ID.

- Assign your methods to each injection. Just click the **Method** cell for the injection and select a method from the drop down menu. If you added your samples with pre-assigned methods and don't need to make any changes, skip to the next step.

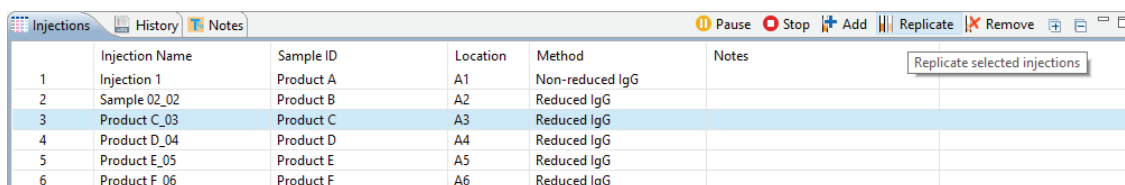
	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	

Hovering over a method name displays the method parameters:

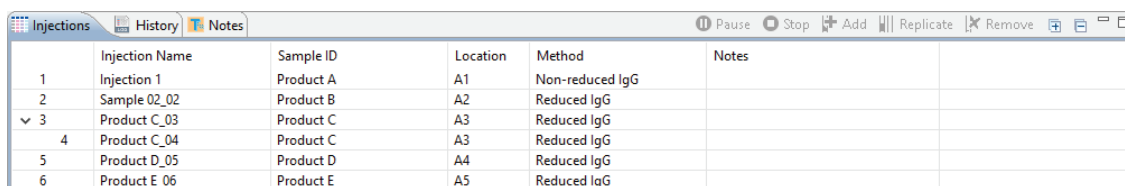
	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	

4. You can add or remove injections as needed using a few different options. If you don't need to make any changes here, just skip to the next step.

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

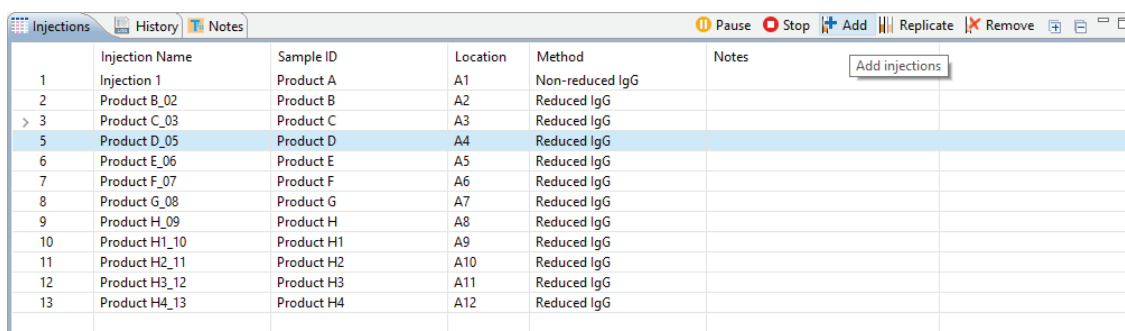


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product E_05	Product E	A5	Reduced IgG	
6	Product F_06	Product F	A6	Reduced IgG	



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product C_04	Product C	A3	Reduced IgG	
5	Product D_05	Product D	A4	Reduced IgG	
6	Product E_06	Product E	A5	Reduced IgG	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Product B_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product E_05	Product E	A5	Reduced IgG	
6	Product F_06	Product F	A6	Reduced IgG	
7	Product G_07	Product G	A7	Reduced IgG	
8	Product H_08	Product H	A8	Reduced IgG	
9	Product I_09	Product I	A9	Reduced IgG	
10	Product J_10	Product J	A10	Reduced IgG	
11	Product K_11	Product K	A11	Reduced IgG	
12	Product L_12	Product L	A12	Reduced IgG	
13	Product M_13	Product M	A13	Reduced IgG	

- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

Step 5 - Add Programmed Pauses and Stops (Optional)

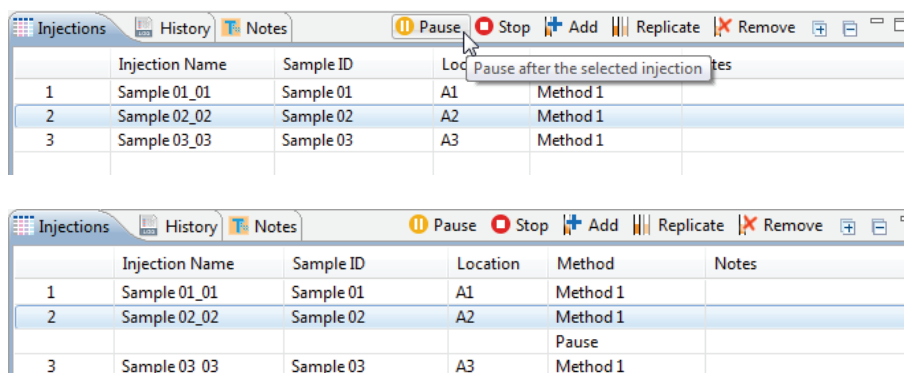
You can program pauses and stops in the batch to automatically pause or stop after injections even when you aren't in front of Maurice.

To program a pause:

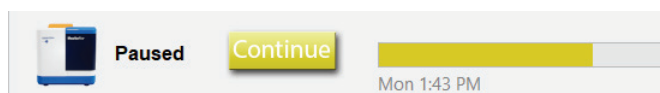
Pauses are helpful when you want to review injection results before continuing with the rest of the batch.

1. Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete.

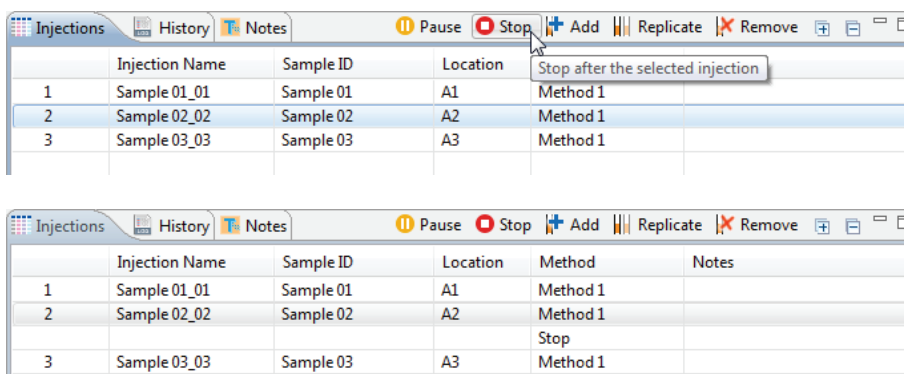
NOTE: Maurice can tweet you when the batch pauses. See “Setting Up Maurice Systems to Send Tweets” on page 764.



2. To resume the batch, click **Continue** in the instrument status bar:

**To stop the run after a specific injection:**

1. Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.

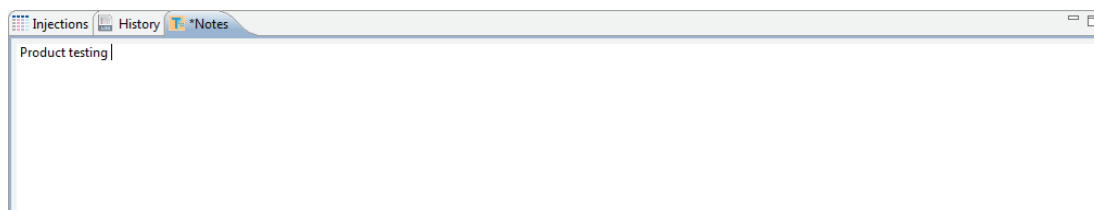


Data collected for injections completed up to the point where the batch was stopped are saved in the .mbz file.

2. Be sure to complete your normal post-batch procedures.

Step 6 - Add Batch Notes (Optional)

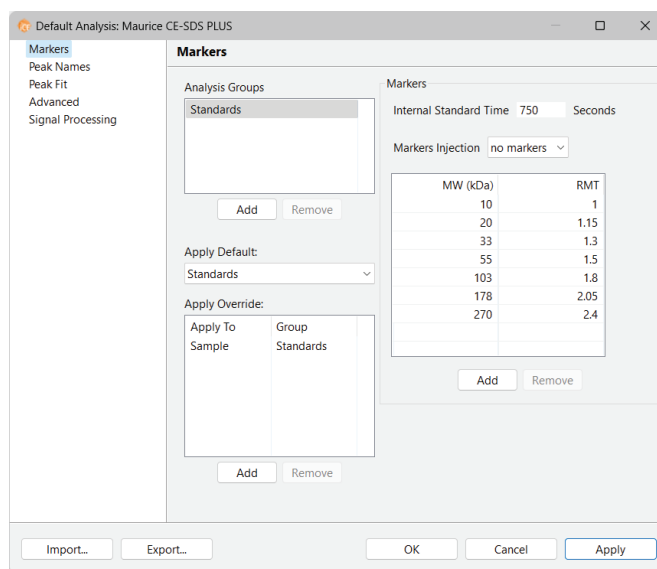
1. Click on the **Notes** pane.
2. Click in the notes area and type any information you want to add about your batch.



Step 7 - Modify Default Analysis Parameters (Optional)

You can preset the parameters used to analyze data generated with the batch. We recommend using the default parameters for CE-SDS PLUS applications, but if you need to modify parameters:

1. Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



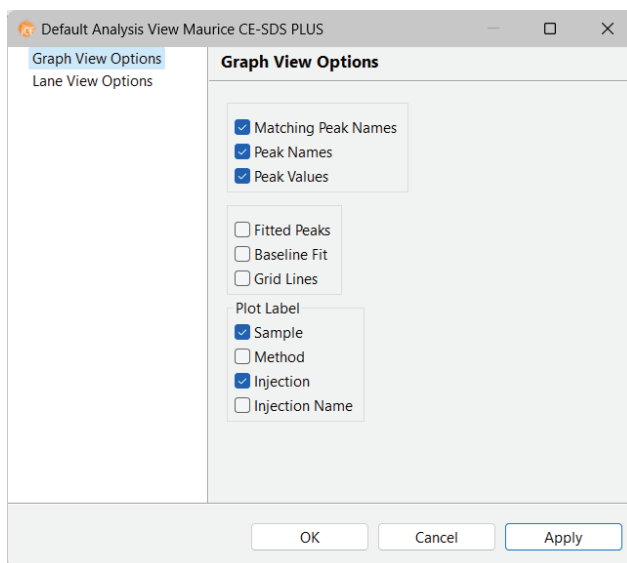
2. Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to “Analysis Settings Overview” on page 713

Step 8 - Modify Default Analysis View Parameters (Optional)

You can preset the parameters used to display the graph and lane view data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph View Options and Lane View Options buttons.

To modify the parameters:

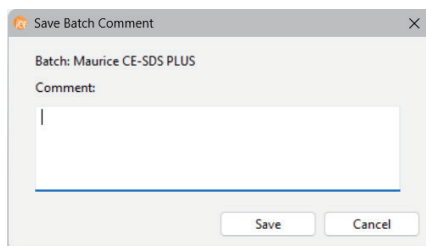
1. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



2. Change the parameters you want to, then click **Apply**. Click **OK** when you are done changing display parameters. For detailed information on graph and lane view options, please refer to “Customizing the Data Display” on page 695.

Step 9 - Save Your Batch

1. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



2. Enter a name for your batch then click **Save**.

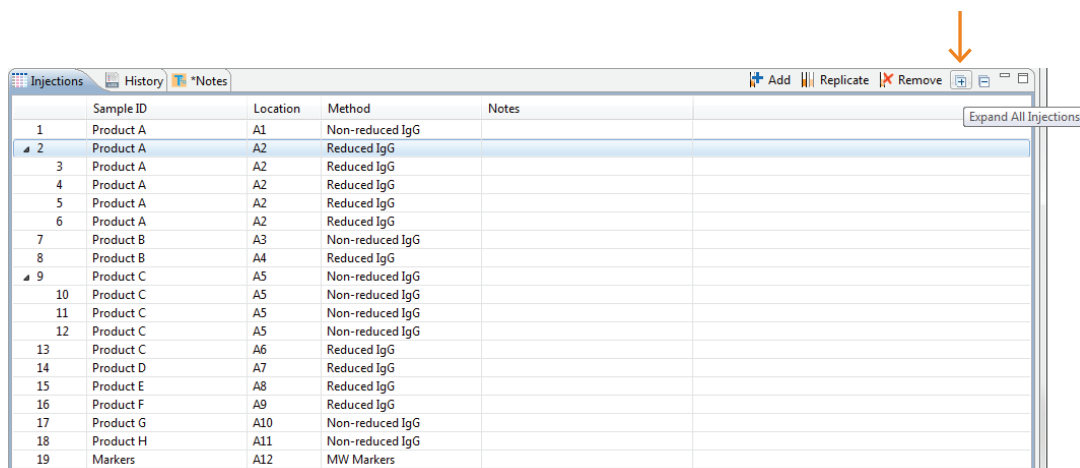
Viewing Replicate Injections

Injections with replicates have an arrow next to the injection number. You can expand or collapse the list of replicate injections by toggling the arrow:

	Sample ID	Location	Method
1	Product A	A1	Non-reduced IgG
▶ 2	Product A	A2	Reduced IgG
7	Product B	A3	Non-reduced IgG
8	Product B	A4	Reduced IgG
▶ 9	Product C	A5	Non-reduced IgG
13	Product C	A6	Reduced IgG
14	Product D	A7	Reduced IgG
15	Product E	A8	Reduced IgG
16	Product F	A9	Reduced IgG
17	Product G	A10	Non-reduced IgG
18	Product H	A11	Non-reduced IgG
19	Markers	A12	MW Markers

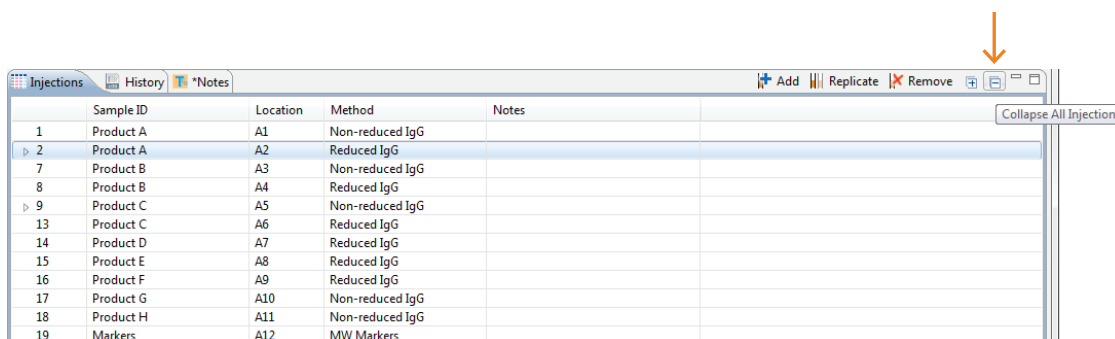
	Sample ID	Location	Method
1	Product A	A1	Non-reduced IgG
▶ 2	Product A	A2	Reduced IgG
3	Product A	A2	Reduced IgG
4	Product A	A2	Reduced IgG
5	Product A	A2	Reduced IgG
6	Product A	A2	Reduced IgG
7	Product B	A3	Non-reduced IgG
8	Product B	A4	Reduced IgG
▶ 9	Product C	A5	Non-reduced IgG
13	Product C	A6	Reduced IgG
14	Product D	A7	Reduced IgG
15	Product E	A8	Reduced IgG

- To show all replicate injections in the batch, click the **Expand All Injections** button.



	Sample ID	Location	Method	Notes
1	Product A	A1	Non-reduced IgG	
▶ 2	Product A	A2	Reduced IgG	
3	Product A	A2	Reduced IgG	
4	Product A	A2	Reduced IgG	
5	Product A	A2	Reduced IgG	
6	Product A	A2	Reduced IgG	
7	Product B	A3	Non-reduced IgG	
8	Product B	A4	Reduced IgG	
▶ 9	Product C	A5	Non-reduced IgG	
10	Product C	A5	Non-reduced IgG	
11	Product C	A5	Non-reduced IgG	
12	Product C	A5	Non-reduced IgG	
13	Product C	A6	Reduced IgG	
14	Product D	A7	Reduced IgG	
15	Product E	A8	Reduced IgG	
16	Product F	A9	Reduced IgG	
17	Product G	A10	Non-reduced IgG	
18	Product H	A11	Non-reduced IgG	
19	Markers	A12	MW Markers	

- To hide all replicate injections in the batch, click the **Collapse All Injections** button.



	Sample ID	Location	Method	Notes
1	Product A	A1	Non-reduced IgG	
▶ 2	Product A	A2	Reduced IgG	
7	Product B	A3	Non-reduced IgG	
8	Product B	A4	Reduced IgG	
▶ 9	Product C	A5	Non-reduced IgG	
13	Product C	A6	Reduced IgG	
14	Product D	A7	Reduced IgG	
15	Product E	A8	Reduced IgG	
16	Product F	A9	Reduced IgG	
17	Product G	A10	Non-reduced IgG	
18	Product H	A11	Non-reduced IgG	
19	Markers	A12	MW Markers	

Batch History

Clicking on the History pane shows the batch event history, starting with the date and time the batch was created through the most current event. Clicking on a row in the table displays the full event details in the box under the table.

[illegible]

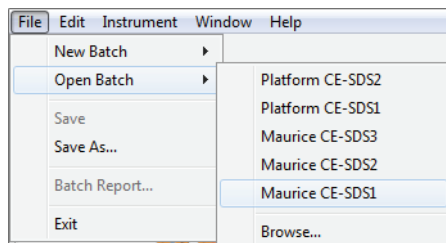
- **Date:** Date and time of the batch event.
- **User Name:** User that initiated the event. User names only display if you're using the Access Control feature to help satisfy the 21CFR Part 11 data security requirements. Go to "Enabling Access Control" on page 781 to learn how to set it up.
- **Message:** Description of the event that took place.
- **Comment:** Comments entered by the user when the batch was saved.

You can copy the information in the History pane to use in other documents:

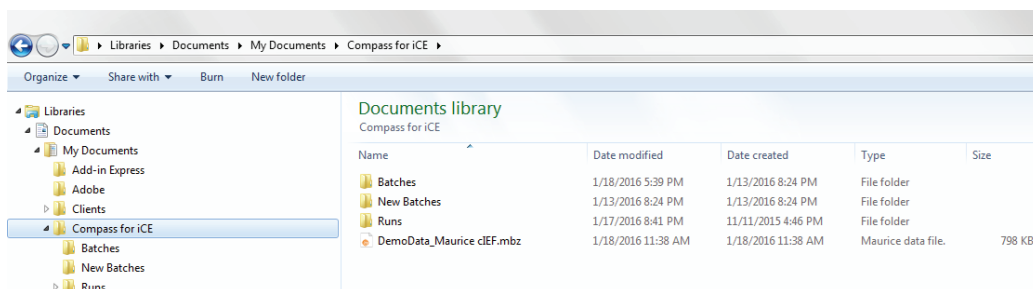
1. Click the **History** pane to make sure it's active.
2. Click **Edit** in the main menu and select **Copy**.
3. Open a document and click **Paste**.

Making Changes to a Batch

1. Select **File** in the main menu and click **Open Batch**.



2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file.
 - Select one of these files or click **Browse** to open the Batch folder and select a different one.
 - To open a batch used in a run file, click **Browse**. Go to the Run folder in the Compass for iCE directory and select the run file that uses the batch you want to edit.

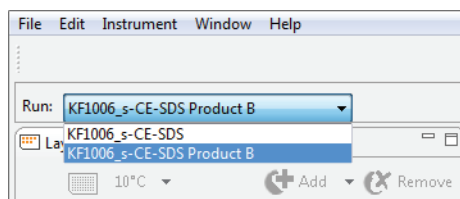


3. To make changes to the batch, see the steps in “Creating a New Batch” on page 84. Then select **File** from the main menu and click **Save** or **Save As**.

Viewing and Editing Batches in Completed Runs

1. Click the **Analysis** screen and open your run file(s).
2. After the run opens, click the **Batch** screen to view the batch used with the run.

If you opened more than one run file, you can switch between viewing batches for each file. Just click the **down arrow** in the Run box and select the batch you want to view from the drop down list.



- In completed run files, you can only make edits to the **sample** and **method** names in a batch, and these changes are saved to the run file only. Update sample and method names as needed.
- Click the **Analysis** screen. Then select **File** from the main menu and click **Save** or **Save As** to save the new changes to the run file.

Copying and Pasting Injection Names or Sample IDs from other Documents

Injection names and Sample IDs can be copied from other documents and pasted into the Injection pane to make batch setup faster.

To paste a list of injection names into the batch:

- Select the list of injection names in a document (Microsoft® Word®, Excel®, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 2	My Sample 2
4		My Injection 3	My Sample 3
5		My Injection 4	My Sample 4
6		My Injection 5	My Sample 5
7		My Injection 6	My Sample 6
8		My Injection 7	My Sample 7
9		My Injection 8	My Sample 8
10		My Injection 9	My Sample 9
11		My Injection 10	My Sample 10

- Select an injection in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the injection, right click and select **Paste**.

⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove 📄					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	Method2	
3	Sample 03_03	Sample 03	A3	Method2	
4	Sample 04_04	Sample 04	A4	Method2	
5	Sample 05_05	Sample 05	A5	Method2	
6	Sample 06_06	Sample 06	A6	Method2	
7	Sample 07_07	Sample 07	A7	Method2	
8	Sample 08_08	Sample 08	A8	Method2	
9	Sample 09_09	Sample 09	A9	Method2	
10	Sample 10_10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

The injection names are pasted into the Injection pane:

⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove ⌵ 📄					
	Injection Name	Sample ID	Location	Method	Notes
1	My Injection 1	Sample 01	A1	System Suitability	
2	My Injection 2	Sample 02	A2	Method2	
3	My Injection 3	Sample 03	A3	Method2	
4	My Injection 4	Sample 04	A4	Method2	
5	My Injection 5	Sample 05	A5	Method2	
6	My Injection 6	Sample 06	A6	Method2	
7	My Injection 7	Sample 07	A7	Method2	
8	My Injection 8	Sample 08	A8	Method2	
9	My Injection 9	Sample 09	A9	Method2	
10	My Injection 10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

To paste a list of Sample IDs into the batch:

1. Select the list of Sample IDs in a document (Microsoft Word, Excel, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 2	My Sample 2
4		My Injection 3	My Sample 3
5		My Injection 4	My Sample 4
6		My Injection 5	My Sample 5
7		My Injection 6	My Sample 6
8		My Injection 7	My Sample 7
9		My Injection 8	My Sample 8
10		My Injection 9	My Sample 9
11		My Injection 10	My Sample 10
12			

2. Select a Sample ID in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the Sample ID, right click and select **Paste**.

📄 Injections 📄 History 📄 Notes ⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove ⌵ 📄					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	Method2	
3	Sample 03_03	Sample 03	A3	Method2	
4	Sample 04_04	Sample 04	A4	Method2	
5	Sample 05_05	Sample 05	A5	Method2	
6	Sample 06_06	Sample 06	A6	Method2	
7	Sample 07_07	Sample 07	A7	Method2	
8	Sample 08_08	Sample 08	A8	Method2	
9	Sample 09_09	Sample 09	A9	Method2	
10	Sample 10_10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

The Sample IDs are pasted into the Injection pane.

NOTE: If you paste in new sample IDs when the default injection names are still being used, the injection names will automatically update to match the sample ID. If you've already changed the injection names manually, updating the sample IDs will not update the injection names.

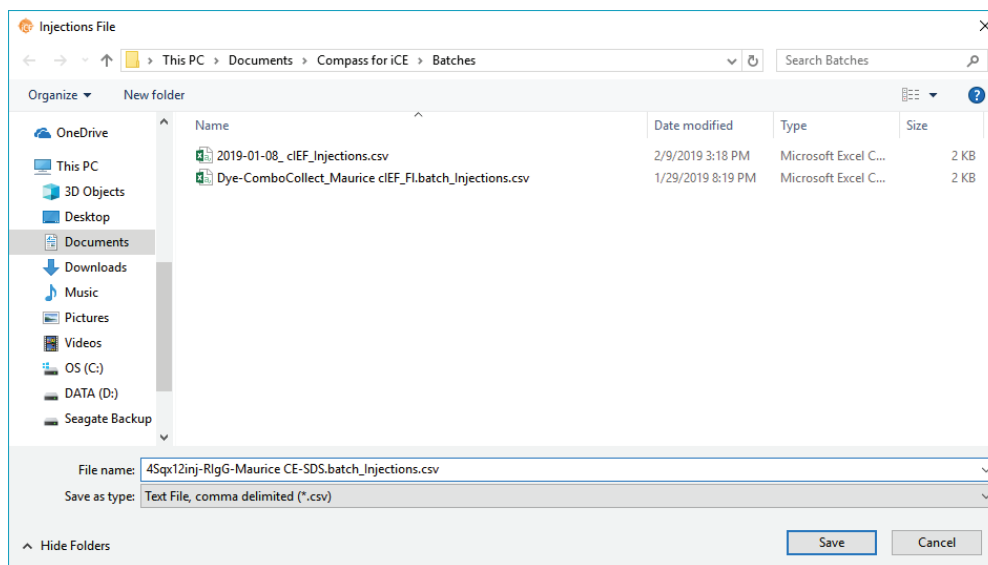
⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove ⚙ 🗑					
	Injection Name	Sample ID	Location	Method	Notes
1	My Sample 1_01	My Sample 1	A1	System Suitability	
2	My Sample 2_02	My Sample 2	A2	Method2	
3	My Sample 3_03	My Sample 3	A3	Method2	
4	My Sample 4_04	My Sample 4	A4	Method2	
5	My Sample 5_05	My Sample 5	A5	Method2	
6	My Sample 6_06	My Sample 6	A6	Method2	
7	My Sample 7_07	My Sample 7	A7	Method2	
8	My Sample 8_08	My Sample 8	A8	Method2	
9	My Sample 9_09	My Sample 9	A9	Method2	
10	My Sample 10_10	My Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

Importing and Exporting Injections

Injections in an open batch or run file can be exported as a separate file. This lets you import the same or a modified set of injections into another batch later, rather than having to enter the information manually.

Exporting Injections

1. Open the batch or run you want to export injections from.
2. In the Batch screen, select **File** in the main menu and click **Export Injections**. The following window displays:



3. The default directory is Compass for iCE/Batches. Change the directory if needed.
4. The default file name will be batchname.batch_Injections.csv. Change the name if needed and click **Save**.

Creating an Injections File

To create a new injections file .csv, it's easiest to export injections from an existing batch and use that as your template.

1. Follow the instructions in Exporting Injections above to export injections from an existing CE-SDS PLUS batch.
2. Open the .csv file in a program that provides a table/spreadsheet format.
3. To add or change injection information:
 - a. Type in a unique name in each cell in the Injection Name column. For replicate injections, type > in front of the injection name, for example: >Injection 1.
 - b. Type in a sample ID, location and method for each injection in their respective columns.
 - c. Optional: Type in notes if needed.

NOTES:

Mix bottle is not used for CE-SDS PLUS batches. Compass for iCE ignores this column when importing injections into CE-SDS PLUS batches.

Programmed pauses and stops can't be added in the injections file. They can be added in the Injections pane after importing your injections.

	A	B	C	D	E	F
1	Injection Name	Sample ID	Location	Method	Mix Bottle	Notes
2	R IgG1	SB	B2	Reduced IgG		
3	R IgG 2	SB	B3	Reduced IgG		
4	>R IgG 2	SB	B3	Reduced IgG		
5	>R IgG 2	SB	B3	Reduced IgG		
6	R IgG 3	SB	C2	Reduced IgG		
7	R IgG 4	SB	C3	Reduced IgG		
8	R IgG1_05	SB	B2	Reduced IgG		
9	R IgG 2_06	SB	B3	Reduced IgG		
10	R IgG 3_07	SB	C2	Reduced IgG		

4. Save the .csv file.

Importing Injections

1. Open the batch you want to import injections into, or open a new batch.
2. Select **File** in the main menu and click **Import Injections**.
3. Select an injections file (*.csv) and click **OK**. The imported information will display in the Layout and Injections panes.

NOTE: Imported injections will replace any injections already in the Injections table, they will not append an existing table.

Importing and Exporting Methods

Methods in an open batch or run file can be exported as a separate file. This lets you import the same method into another batch later, rather than having to re-enter method information manually.

Importing a Method

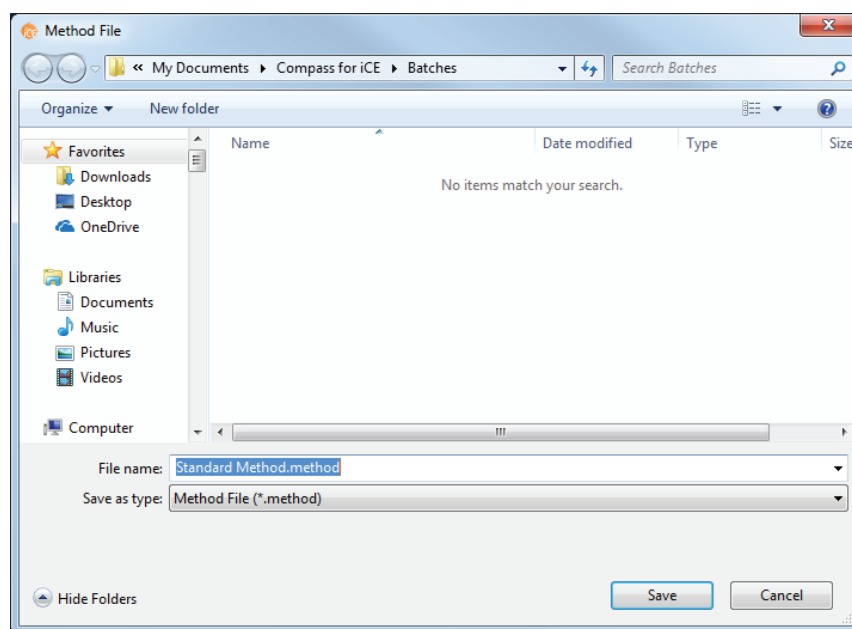
NOTE: Importing a method imports information into the Batch window's Method pane only.

1. Open the batch you want to import the method into.
2. Select **File** in the main menu and click **Import Method**.
3. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

Exporting a Method

NOTE: Exporting a method exports information in the Batch window's Method pane only.

1. Open the batch you want to export the method from.
2. In the Method pane, select the method you want to export. To export more than one method at the same time, hold the **Ctrl** key to select multiple methods.
3. Select **File** in the main menu and click **Export Method**. The following window displays:



4. The default directory is Compass for iCE/Batches. Change the directory if needed.
5. Enter a method file name and click **Save**. The settings will be saved as a *.method file.

Batch Reports

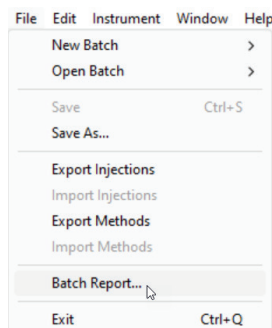
You can export a PDF file of sample and method details for each injection in the batch.

NOTES:

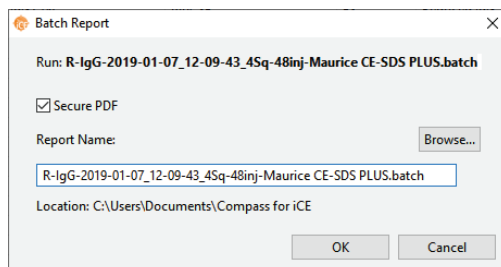
You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it’s saved in a separate run folder. Existing reports aren’t overwritten.

1. Go to the **Analysis** or **Run Summary** screen, then click **File > Open Run** and select a run file (if you don’t have one open already).
2. After the run opens, go to the **Batch** screen.
3. Select **File** from the main menu and click **Batch Report**.

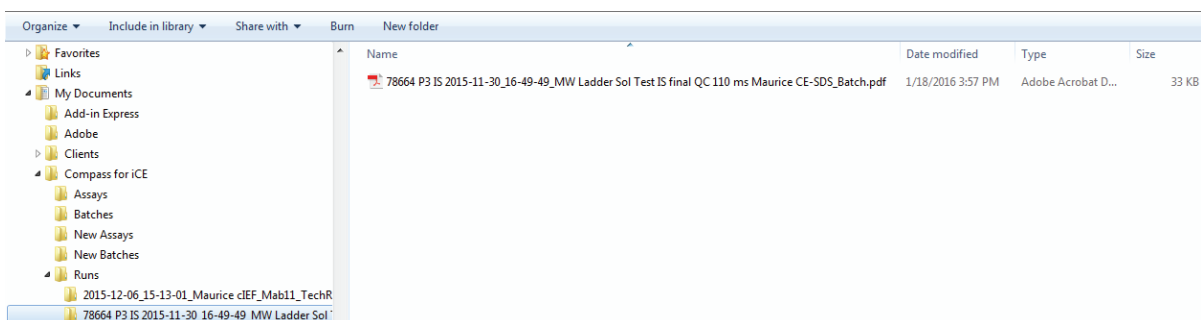


4. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.



Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don’t want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.

5. The Batch Report PDF is exported to the Runs folder in the Compass for iCE directory. It'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Here's an example Batch Report:

CE-SDS Batch: 4Sqx12inj-RlgG-Maurice CE-SDS

Injection	Injection Name	Sample ID	Location	Method	Sample Load	Separation
1	R IgG1	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
2	R IgG 2	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
3	R IgG 3	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
4	R IgG 4	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
5	R IgG1_05	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
6	R IgG 2_06	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
7	R IgG 3_07	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
8	R IgG 4_08	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
9	R IgG1_09	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
10	R IgG 2_10	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
11	R IgG 3_11	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
12	R IgG 4_12	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
13	R IgG1_13	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
14	R IgG 2_14	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
15	R IgG 3_15	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
16	R IgG 4_16	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
17	R IgG1_17	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
18	R IgG 2_18	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
19	R IgG 3_19	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
20	R IgG 4_20	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
21	R IgG1_21	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
22	R IgG 2_22	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
23	R IgG 3_23	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
24	R IgG 4_24	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
25	R IgG1_25	SB	B2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
26	R IgG 2_26	SB	B3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
27	R IgG 3_27	SB	C2	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts
28	R IgG 4_28	SB	C3	Reduced IgG	20 sec 4600 Volts	25.0 min, 5750 Volts

Created By: Jacquelyn Sat 4:47 PM Feb 2, 2019 CST
 C:\Users\Jacquelyn\Documents\ProteinSimple\Maurice\User Guide\Rev 12 edits\Data from Andrea - no edits\R-IgG-2019-01-10_16-25-12_45qx12inj-RlgG-Maurice CE-SDS.mbz
 Computer: DESKTOP-C7FPQ08

Page 1 of 3



CE-SDS Batch: 4Sqx12inj-RlgG-Maurice CE-SDS

Batch Log

Date	User Name	Message	Comment
2019-01-10 16:22:44		Batch created using the factory default Maurice CE-SDS with Compass for iCE Version: 2.1.0-1219	
2019-01-10 16:25:08		Saved as C:\Users\jkazakova\Documents\Compass for iCE\Batches\4Sqx12inj-RlgG-Maurice CE-SDS.batch from Compass for iCE v2.1.0-1219	
2019-01-10 16:25:08		Save injections changes from Compass for iCE v2.1.0-1219	

Created By: Jacquelyn Sat 4:47 PM Feb 2, 2019 CST Software Version: 2.1.0, Build ID: 0130
C:\Users\Jacquelyn\Documents\Clients\ProteinSimple\Maurice\User Guide\Rev 12 edits\Data from Andrea - no edits\RlgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS.mbz
Computer: DESKTOP-CTFPQ08

Page 3 of 3



Chapter 12:

Running CE-SDS PLUS Applications

Chapter Overview

- Overview
- Before You Throw the Switch
- Power Up
- Running CE-SDS PLUS Applications
- Editing a Running Batch
- Post-batch Procedures
- Checking Your Data

Overview

CE-SDS PLUS applications can be run on Maurice, Maurice S. and MauriceFlex systems with a CE-SDS PLUS cartridge.

Before You Throw the Switch

Ensure that everyone using Maurice have:

- Received instruction in general safety practices for laboratories.
- Received instruction in specific safety practices for Maurice.
- Received instruction on handling of biohazards (if biohazardous materials are to be used on Maurice).
- Read and understood all Product Inserts and related Safety Data Sheets (SDS).

Power Up

1. Turn on the computer connected to Maurice.
2. Turn on Maurice's main power switch.
3. Wait for Maurice to initialize. His indicator light will turn from magenta to blue when he's done.
4. Double-click the Compass for iCE software icon on the computer desktop to open the application.
5. Connect Maurice to Compass for iCE.

Running CE-SDS PLUS Applications

What You'll Need

- Maurice CE-SDS PLUS Application Kit which include:
 - Maurice CE-SDS PLUS Cartridges
 - Cartridge Cleaning Vials
 - Separation Matrix
 - Running Buffer (Top and Bottom)
 - CE-SDS PLUS 1X Sample Buffer
 - Wash Solution
 - Conditioning Solutions (1 and 2)
 - 25X Internal Standard

- Glass reagent vials, 2 mL
- 96-well plates
- Clear screw caps for vials
- Orange pressure caps for vials
- Maurice CE-SDS IgG Standard (optional)
- Maurice CE-SDS MW Markers (optional)
- β -mercaptoethanol (β ME, >98% = 14.2 M) for reducing conditions
- Iodoacetamide (IAM, 250 mM) for alkylation at non-reducing conditions
- Deionized (DI) water
- Sample vials with integrated inserts for samples (optional)
- P10, P20, P200, P1000 and pipette tips
- Water bath or thermocycler
- Vortexer
- Table-top centrifuge with plate adapter or vial adapter (12 mm, 2 mL vials)

Step 1: Prep Your Internal Standard, Samples and Reagents

NOTE: You can prepare your samples to run either in 96-well plates or vials. Using 96-well plates is the default method.

Internal Standard

NOTES:

The Internal Standard is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Aliquot the reconstituted solution into appropriately sized vials and store at -70 to -90 °C for long term storage. For short-term storage (< 1 week), the solution can be stored at 2-8 °C.

Prepare the Internal Standard in the same buffer as your sample.

1. Open the vial of lyophilized 25X Internal Standard by lifting the center tab and gently pulling it back to break the metal seal. Then slowly remove the rubber stopper.
2. Reconstitute by adding 240 μ L of 1X Sample Buffer. Pipette up and down a few times to mix thoroughly. This results in a 25X Internal Standard solution.

NOTE: Don't vortex the reconstituted Internal Standard during prep.

Sample Prep Under Reducing Conditions

NOTES:

Prepare a minimum of 50 μL of sample for a CE-SDS PLUS batch.

Reduced IgG Sample

1. In a microcentrifuge tube, dilute your IgG sample with 1X Sample Buffer to a concentration of 0.25-1.00 mg/mL in a final volume of 50 μL .
-

NOTE: Dilute at least 1:1 with CE-SDS PLUS 1X Sample Buffer.

2. Add 2 μL of reconstituted 25X Internal Standard for every 50 μL of sample volume.
 3. Add 2.5 μL of 14.2 M β -mercaptoethanol for every 50 μL of sample volume.
 4. Mix thoroughly.
-

NOTE: Heating in the presence of SDS linearizes the protein and allows for SDS binding. The reducing agents break up inter- and intra-molecular disulfide bonds.

5. Centrifuge the tube and heat the mixture in a water bath or thermocycler at 70 °C for 10 minutes.
 6. Put the tube on ice for 5 minutes.
 7. Vortex briefly and spin down.
-

Reduced IgG Standard (Optional)

NOTES:

The IgG Standard is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Prepare the reduced IgG Standard in CE-SDS PLUS 1X Sample Buffer.

1. Using scissors, carefully cut the top of the foil package, leaving the sealing strip intact.

2. Take out the strip of tubes and carefully cut one pink tube of lyophilized IgG Standard from the strip. Put the unopened tubes back in the package, seal tightly and store at 2-8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Reconstitute the IgG Standard with 50 µL of 1X Sample Buffer. Gently resuspend by pipetting the solution up and down. Transfer the solution to a microcentrifuge tube.
5. Add 2 µL of reconstituted 25X Internal Standard.
6. Add 2.5 µL of 14.2 M β-mercaptoethanol.
7. Mix thoroughly by vortex.
8. Centrifuge the tube and heat mixture in a water bath or thermocycler at 70 °C for 10 minutes.
9. Put the tube on ice for 5 minutes.
10. Vortex briefly and spin down.

CE-SDS Molecular Weight (MW) Markers (Optional)

NOTES:

The CE-SDS MW Markers are lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Prepare the MW Markers in CE-SDS PLUS 1X Sample Buffer.

1. Using scissors, carefully cut the top of the foil package leaving the sealing strip intact.
2. Take out the strip of tubes and carefully cut one green tube of lyophilized CE-SDS MW Markers from the strip. Put the unopened tubes back in the package, seal tightly and store at 2-8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Reconstitute the CE-SDS MW Markers with 50 µL of 1X Sample Buffer. Gently resuspend by pipetting the solution up and down. Transfer the solution to a microcentrifuge tube.
5. Add 2 µL of reconstituted 25X Internal Standard.
6. Add 2.5 µL of 14.2 M β-mercaptoethanol.
7. Mix thoroughly.
8. Centrifuge the tube and heat the mixture in a water bath or thermocycler at 70 °C for 10 minutes.
9. Put the tube on ice for 5 minutes.
10. Vortex briefly and spin down.

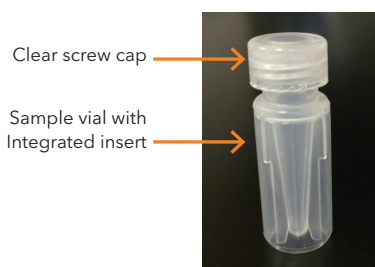
Spin Samples, Standards and CE-SDS MW Markers

If you're using a 96-well plate:

1. Transfer 50 μ L of each of your samples, IgG Standard and CE-SDS MW Markers to their designated wells in a 96-well plate.
2. Cover the plate with a lid and spin for 10 minutes at 1000 xg in a centrifuge with a plate adapter.
3. Pop any bubbles in the samples with a clean pipette tip.

If you're using vials:

1. Transfer 50 μ L of each of your samples, IgG Standard and CE-SDS MW Markers to their designated sample vials with insert, 0.2 mL.
2. Close the vials with a clear screw cap.
3. Place the vials in a centrifuge using a vial adapter (12 mm, 2 mL vials) and spin for 10 minutes at 1000 xg.



Sample Prep Under Non-reducing Conditions

Alkylation Reagent

NOTES:

We use a 250 mM solution of iodoacetamide (IAM) as an alkylating reagent.

Prepare a fresh 250 mM solution of iodoacetamide in DI water before use.

1. Weigh out 46 mg of IAM directly into a 1.5 mL microcentrifuge tube.
2. Add 1 mL of DI water to the tube and mix thoroughly.

Non-reduced IgG Sample

NOTE: Prepare a minimum of 50 μ L of sample for a CE-SDS PLUS batch.

1. In a microcentrifuge tube, dilute your IgG sample with 1X Sample Buffer to a concentration of 0.25-1.0 mg/mL in a final volume of 50 μ L.

NOTE: Dilute at least 1:1 with CE-SDS PLUS 1X Sample Buffer.

2. Add 2 μ L of reconstituted 25X Internal Standard for every 50 μ L of sample volume.
3. Add 2.5 μ L of 250 mM IAM for every 50 μ L of sample volume.
4. Mix thoroughly.
5. Centrifuge the tube and heat the mixture in a water bath or thermocycler at 70 °C for 10 minutes.
6. Put the tube on ice for 5 minutes.
7. Vortex briefly and spin down.

Non-reduced IgG Standard (Optional)

NOTES:

The IgG Standard is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Prepare the non-reduced IgG Standard in CE-SDS PLUS 1X Sample Buffer.

1. Using scissors, carefully cut the top of the foil package, leaving the sealing strip intact.
2. Take out the strip of tubes and carefully cut one pink tube of lyophilized IgG Standard from the strip. Put the unopened tubes back in the package, seal tightly and store at 2-8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Reconstitute the IgG Standard with 50 μ L of 1X Sample Buffer. Gently resuspend by pipetting the solution up and down. Transfer the solution to a microcentrifuge tube.
5. Add 2 μ L of reconstituted 25X Internal Standard.
6. Add 2.5 μ L of 250 mM IAM.
7. Mix thoroughly by vortex.

NOTE: Heating in the presence of SDS linearizes the protein and allows for SDS binding. The alkylating agents prevent disulfide-bond scrambling catalyzed by free sulfhydryl groups. This minimizes the appearance of fragments under non-reducing conditions.

8. Centrifuge the tube and heat mixture in a water bath or thermocycler at 70 °C for 10 minutes.

9. Put the tube on ice for 5 minutes.
10. Vortex briefly and spin down.

Spin Samples and Standards

If you're using a 96-well plate:

1. Transfer 50 μ L of each of your samples and IgG Standard to their designated wells in a 96-well plate.
2. Cover the plate with a lid and spin for 10 minutes at 1000 xg in a centrifuge with a plate adapter.
3. Pop any bubbles in the samples with a clean pipette tip.

If you're using vials:

1. Transfer 50 μ L of your samples and IgG Standard to their designated sample vials with insert, 0.2 mL.
2. Close the vials with a clear screw cap.
3. Place the vials in a centrifuge using vial adapter (12 mm, 2 mL vials) and spin for 10 minutes at 1000 xg.

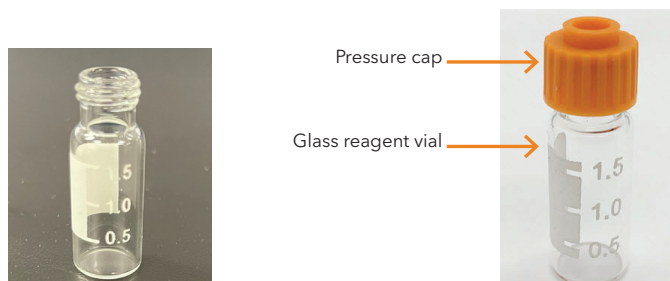
Reagents

IMPORTANT:

Use glass reagent vials, 2 mL (PN 046-017) to prepare CE-SDS PLUS batch reagents.

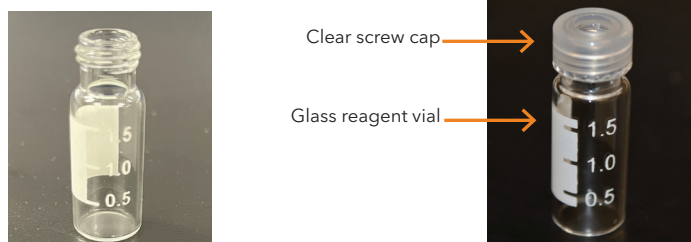
NOTE: Don't reuse reagents, vials or the clear screw caps. Always keep the orange pressure caps paired with their respective reagents. If you want to reuse pressure caps, first wash them thoroughly with DI water, soak overnight in DI water, rinse caps thoroughly with DI water again and air dry them before reusing.

1. Pipette 1.5 mL of Conditioning Solution 1 into a glass reagent vial, label and close with an **orange pressure cap**.



2. Pipette 1.5 mL of Conditioning Solution 2 into a glass reagent vial, label and close with an **orange pressure cap**.
3. Pipette 1.0 mL of Wash Solution into a glass reagent vial, label each and close with an **orange pressure cap**.

- Pipette 1.5 mL of Wash Solution into two glass reagent vials, label each and close the vials with **clear screw caps**.



- Pipette 1 mL of Separation Matrix into a glass reagent vial, label and close with an **orange pressure cap**.
- Pipette 1 mL of Running Buffer - Bottom into one glass reagent vial, label and close with a **clear screw cap**.
- Pipette 1.5 mL of DI water into a glass reagent vials, label and close with an **orange pressure cap**.
- Close an empty glass reagent vial with an **orange pressure cap**.

Step 2: Prep the Cartridge

NOTES:

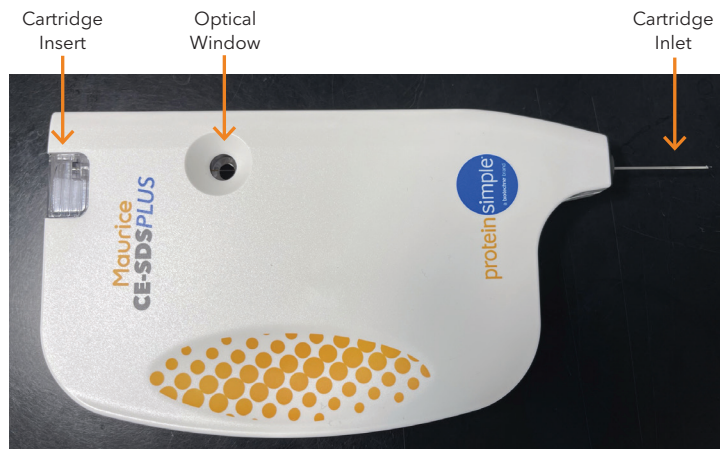
A CE-SDS PLUS Cartridge is guaranteed for 100 injections, with a maximum of 48 injections per batch and a maximum of 25 batches. The absolute injection limit of the cartridge is 500.

- Take the CE-SDS PLUS Cartridge out of its packaging. Save the packaging, you'll need it later to store the cartridge.

NOTE: When you remove the cartridge from its packaging, make sure the cartridge inlet doesn't come in contact with any surface.



2. Pull the cartridge insert out of the cartridge.



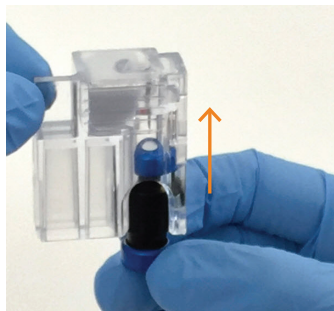
3. Grab a fresh vial of Top Running Buffer from 2-8 °C storage and slide it into the cartridge insert so that the metal pin on the side of the vial is facing out. Press the vial up until it is completely inside the cartridge insert.

NOTES:

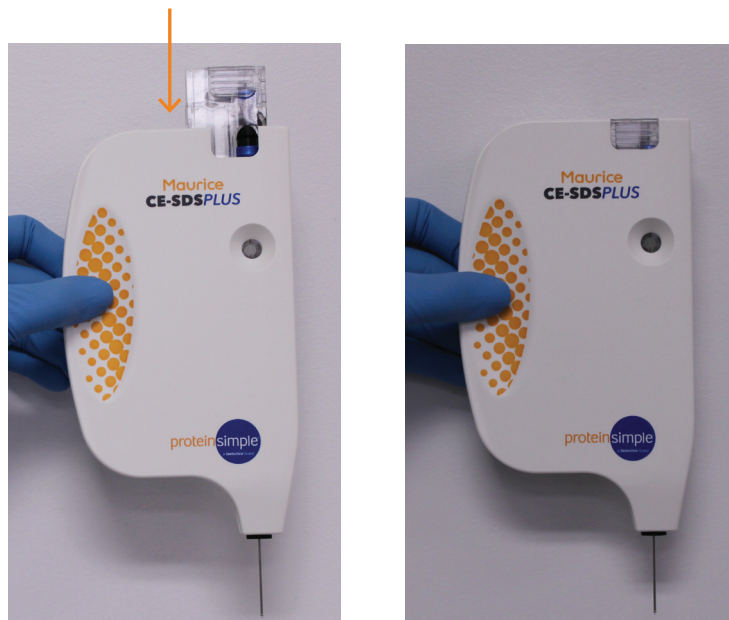
The Top Running Buffer vial should be stored at 2-8 °C when not in use.

The Top Running Buffer vial has metal pins on either side, so no specific orientation is necessary.

Once you've inserted the Top Running Buffer vial, the cartridge insert and the cartridge **must** be kept in an upright position at all times.



4. Slide the cartridge insert back into the cartridge.



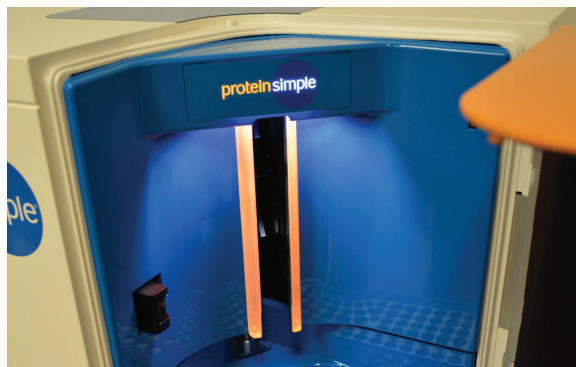
Step 3: Install the Cartridge

1. Open Maurice's door by touching the metal plate on top of the door.



NOTE: The indicator light on Maurice's front panel will blink rapidly as the door disengages.

- Swing the door open to access the cartridge slot and the reagents and samples platform. The lights on either side of the slot will be **orange**.



- Lift the cartridge and hold it vertically using the finger holds on either side, cartridge inlet down, with the CE-SDS PLUS label facing you.
- Gently insert it into the slot. You'll feel the alignment groove on the cartridge align inside the slot when you insert it.



- Continue to slide the cartridge into the slot until the locking mechanism engages. The lights on either side of the slot will change to **blue** once the cartridge is installed correctly.



Step 4: Load Samples and Reagents

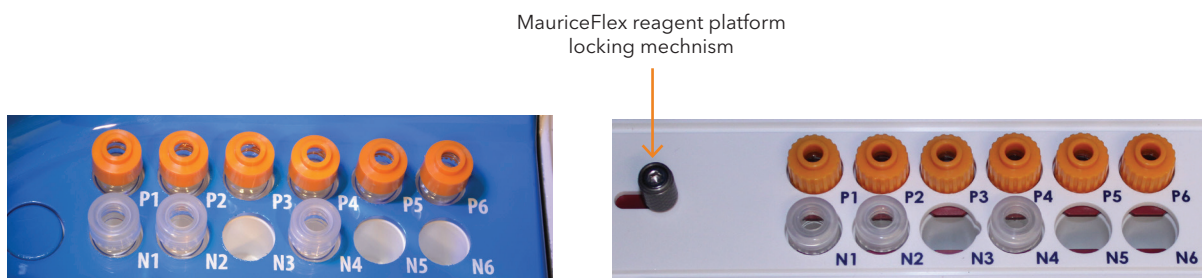
1. Place the reagent vials into their respective positions in the sample and reagents platform:

NOTES:

The two reagent rows in Maurice's reagent and sample platform are kept at room temperature.

Pressure caps are orange and have a raised, ringed surface. They should only be used in Reagent Row P. Use reagent vials with clear screw caps in Row N.

- P1 - Conditioning Solution 1 with **orange pressure cap**
- P2 - Conditioning Solution 2 with **orange pressure cap**
- P3 - DI water with **orange pressure cap**
- P4 - Separation Matrix with **orange pressure cap**
- P5 - Wash Solution with **orange pressure cap**
- P6 - Empty vial (air) with **orange pressure cap**
- N1 - Wash Solution with **clear screw cap**
- N2 - Wash Solution with **clear screw cap**
- N4 - Running Buffer - Bottom with **clear screw cap**



CE-SDS PLUS reagent platform on Maurice and Maurice S. (left) or MauriceFlex (right).

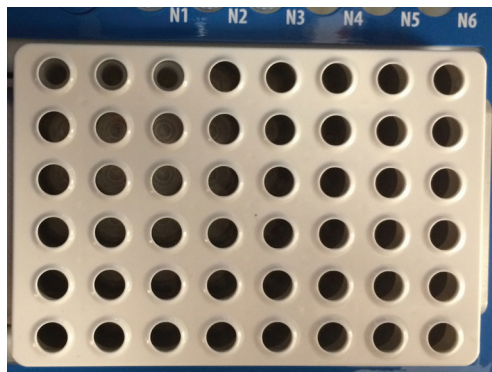
IMPORTANT:

Reagent vials on MauriceFlex must be locked in place by sliding the locking mechanism from the left to the right before starting a batch. A warning will appear in the Start Run window if they are not locked when you start the batch.

- Depending on what you prepared your samples in, put your 96-well sample plate in the metal plate insert or your sample vials in the metal vial insert. If you have a lid on the 96-well sample plate, be sure to remove it before closing the instrument door!

NOTE: Well A1 on the 96-well plate should be in the top left corner of the insert.

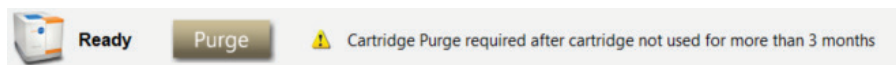
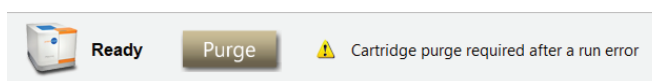
- If you are using a vial tray, place the condensation lid on top of the vials.



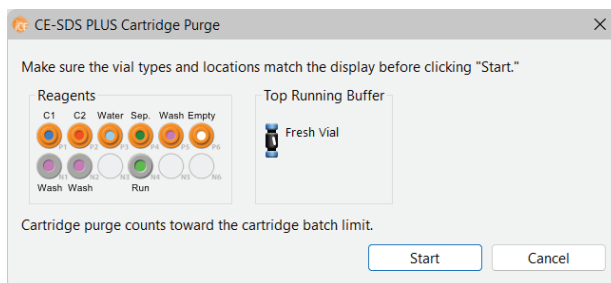
- Close the instrument door. Maurice locks it automatically.

Step 5: Check for Cartridge Alerts

- If your cartridge was last used in a run with an error or if it has not been used in the last three months, you will need to perform a Cartridge Purge. The Maurice system will remind you if a Cartridge Purge is required.
 - To purge the cartridge, click the brown **Purge** button in the instrument status bar.



- Load the reagents required for the cartridge purge and ensure the cartridge is prepped. Then click **Start**.



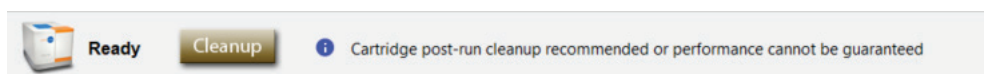
- c. Once the purge is done, perform a Cartridge Post-Run Cleanup by clicking on the brown **Cleanup** button that will appear in the instrument status bar. See step 2 for more info.

NOTE: If there's only one batch left on the cartridge, you'll get a message that it can't be used after purging.

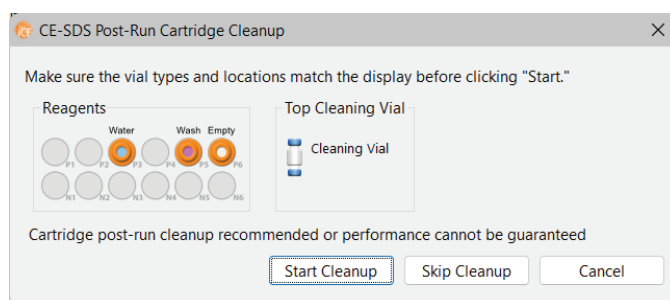
2. If you just completed a Cartridge Purge or if the cartridge you're using did not undergo a cartridge post-run cleanup after the last run, you will be asked if you want to perform a Cartridge Post-Run Cleanup.

To start the Post-Run Cleanup:

- a. Remove the cartridge from Maurice, remove the cartridge insert and check the saturation sensor in the back of the cartridge insert. See "Post-batch Procedures" on page 316 for more information.
- b. Replace the Top Running Buffer vial with a Cleaning Vial in the cartridge insert.
- c. Slide the cartridge insert back into the cartridge.
- d. Re-install the cartridge in Maurice.
- e. Confirm there is a vial of Water (P3), Wash Solution (P5) and Air (P6) in place.
- f. Click the brown **Cleanup** button in the instrument status bar.



- g. Click **Start Cleanup** in the prompt that appears.

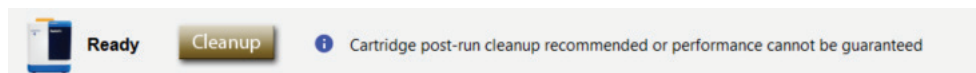


NOTE: You can also start a Post-Run Cleanup by selecting **Instrument > Cartridge Post-Run Cleanup**. When you start a Cartridge Post-Run Cleanup from the instrument menu, click **Start** in the prompt that appears to begin the cleanup.

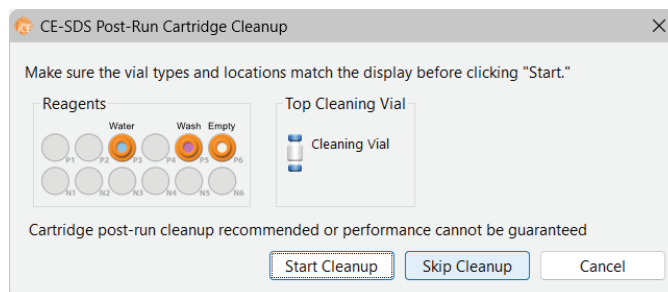
- h. Once the cartridge post-run cleanup is completed, re-prepare the cartridge and install the cartridge in Maurice. See "Step 2: Prep the Cartridge" on page 295 for more information.
- i. Click on the green **Start** button in the instrument status menu to start the batch.

To start the run without a Post-Run Cleanup:

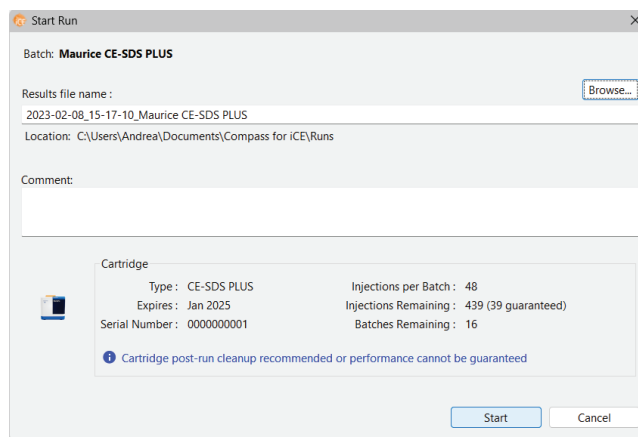
- a. Click the brown **Cleanup** button in the instrument status bar



- b. Click **Skip Cleanup** in the prompt that appears.



- c. Click **Start** in the Start Run window. The run summary and injection report will record that the cartridge did not undergo a post-run cleanup before starting the batch.

**To start the run with a different cartridge:**

- a. If necessary, click **Cancel** in the CE-SDS PLUS Cartridge Post-Run Cleanup window.
- b. Open Maurice's door, remove the first cartridge from Maurice and prepare a second cartridge. See "Step 2: Prep the Cartridge" on page 295 for more information.

Step 6: Create a Batch

1. Launch Compass for iCE.
2. Select the **Batch** tab. This is where you'll enter sample/injection information, methods and batch parameters.

The screenshot shows the 'Batch: Maurice CE-SDS PLUS' window. The 'Layout' tab on the left shows a 12-well plate layout with columns labeled C1, C2, Water, Sep, Wash, Empty and rows labeled A through H. The 'Injections' tab is active, displaying a table with columns: Injection Name, Sample ID, Location, Method, and Notes. The first row contains 'Sample 01_01', 'Sample 01', 'A1', 'Reduced IgG', and an empty 'Notes' field. The 'Methods' tab at the bottom shows a table with columns: Name, Sample Load, and Separation. It lists three methods: 'Reduced IgG' (20 sec 4600 Volts, 25.0 min 5750 Volts), 'Non-reduced IgG' (20 sec 4600 Volts, 35.0 min 5750 Volts), and 'MW Markers' (20 sec 4600 Volts, 35.0 min 5750 Volts).

Injection Name	Sample ID	Location	Method	Notes
Sample 01_01	Sample 01	A1	Reduced IgG	

Name	Sample Load	Separation
Reduced IgG	20 sec 4600 Volts	25.0 min 5750 Volts
Non-reduced IgG	20 sec 4600 Volts	35.0 min 5750 Volts
MW Markers	20 sec 4600 Volts	35.0 min 5750 Volts

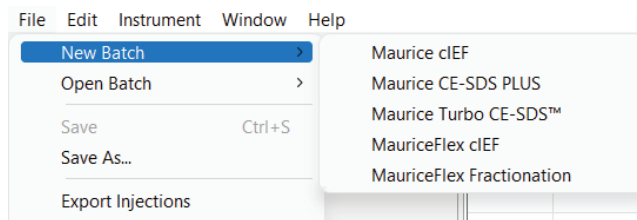
3. To create a batch, make sure your Maurice system is connected to Compass for iCE. Select **Instrument** and click **Connect**.
 - a. If your instrument is listed, select your Maurice system and click **Connect**.
 - b. If your instrument isn't listed, click on the **Settings** button and connect by typing in your instrument IP address.

The 'Connect' dialog box has a table with columns: Name, Location, and Serial Num... Below the table is a 'Settings' button (represented by a gear icon) and 'Connect' and 'Cancel' buttons.

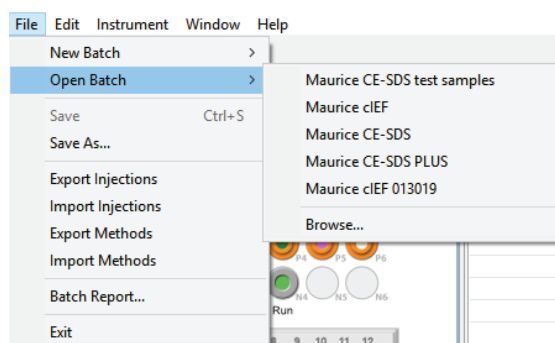
Name	Location	Serial Num...

To create a new batch:

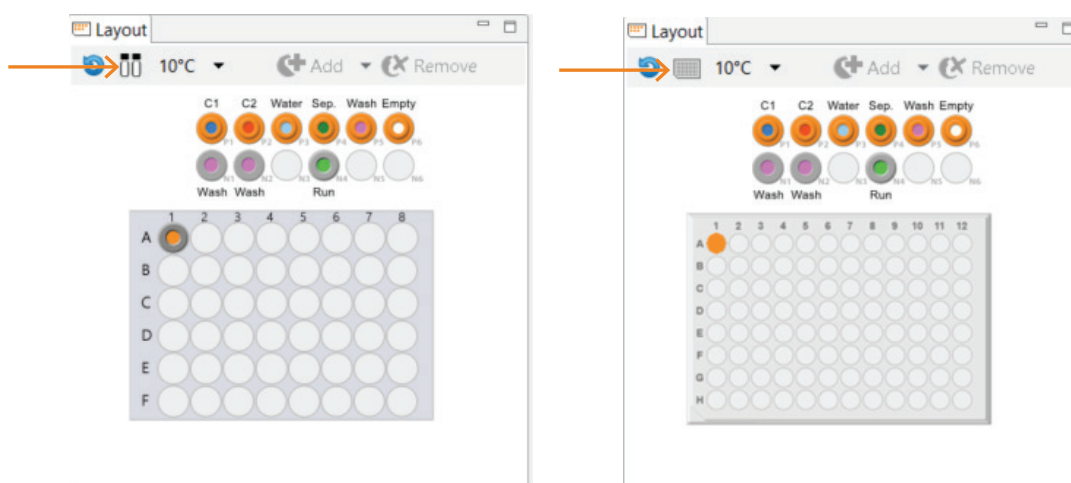
- In the main menu, select **File > New Batch > Maurice CE-SDS Plus**.

**To use an existing batch:** In the main menu, select **File > Open Batch**.

NOTE: If you're making changes to an existing batch, follow the steps in this section. Otherwise skip to "Step 7: Start the Batch" on page 312.



4. In the Layout pane, clicking on the vial/plate icon toggles between formats. Select a 96-well plate or 48 vials depending on what you're running.



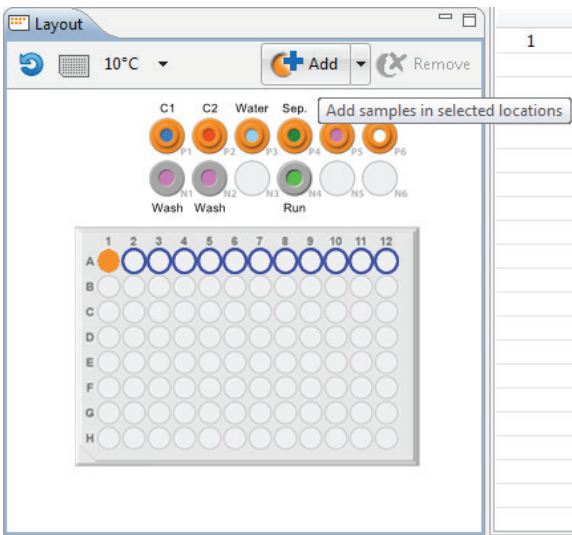
5. Add your samples:

To import samples using a saved injections file:

- a. Select **File** in the main menu and click **Import Injections**.
- b. Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.

NOTE: For more information on setting up an injections file for import, see “Importing and Exporting Injections” on page 281.

To select samples manually: Use your mouse to highlight the well positions for each of your samples, then click **Add**.



This populates the Injections table:

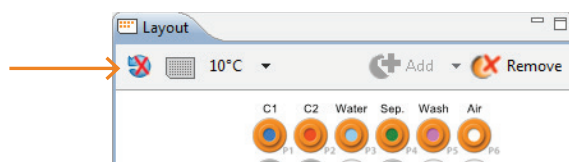
	Injection Name	Sample ID	Location	Method	Notes	
1	Sample 1_01	Sample 1	A1	Reduced IgG		
2	Sample 02_02	Sample 02	A2	Reduced IgG		
3	Sample 03_03	Sample 03	A3	Reduced IgG		
4	Sample 04_04	Sample 04	A4	Reduced IgG		
5	Sample 05_05	Sample 05	A5	Reduced IgG		
6	Sample 06_06	Sample 06	A6	Reduced IgG		
7	Sample 07_07	Sample 07	A7	Reduced IgG		
8	Sample 08_08	Sample 08	A8	Reduced IgG		
9	Sample 09_09	Sample 09	A9	Reduced IgG		
10	Sample 10_10	Sample 10	A10	Reduced IgG		
11	Sample 11_11	Sample 11	A11	Reduced IgG		
12	Sample 12_12	Sample 12	A12	Reduced IgG		

6. Compass for iCE can monitor the current during a separation for you, stop it if an abnormal current profile is detected and reinject the sample for you automatically. Reinjections are enabled by default, but if you don't want to use this feature just click the reinject icon to toggle it off.

NOTES:

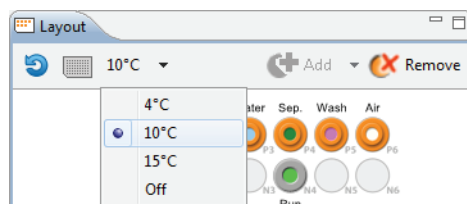
If a sample was reinjected, the injection row will be flagged in the Injections pane in the Run Summary screen. See "Injection Flags" on page 414 for more info.

A maximum of 10 reinjections are allowed per batch.



7. The sample platform is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The two reagent rows are kept at room temperature.



8. Enter your method parameters in the pane.

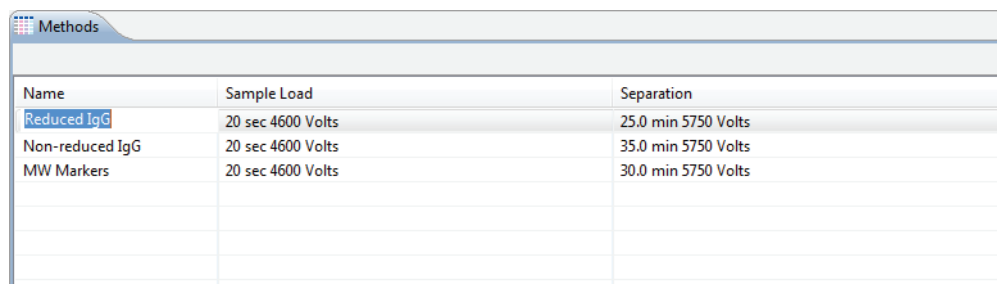
NOTE: There are three default methods. We recommend using the default method parameters for the listed samples. Please see the [Maurice CE-SDS Application Guide](#) for more information on method optimization.

To import a saved method:

- a. Select **File** in the main menu and click **Import Method**.
- b. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

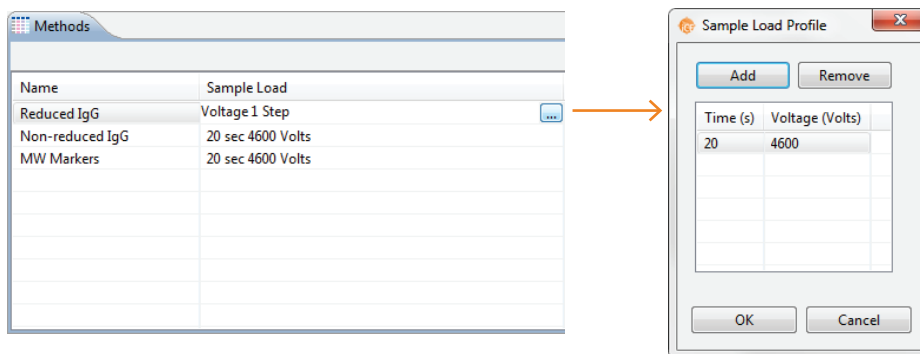
To create a new method or make changes to an existing one:

- a. Click the first cell in the Name column and enter a new method name if needed.



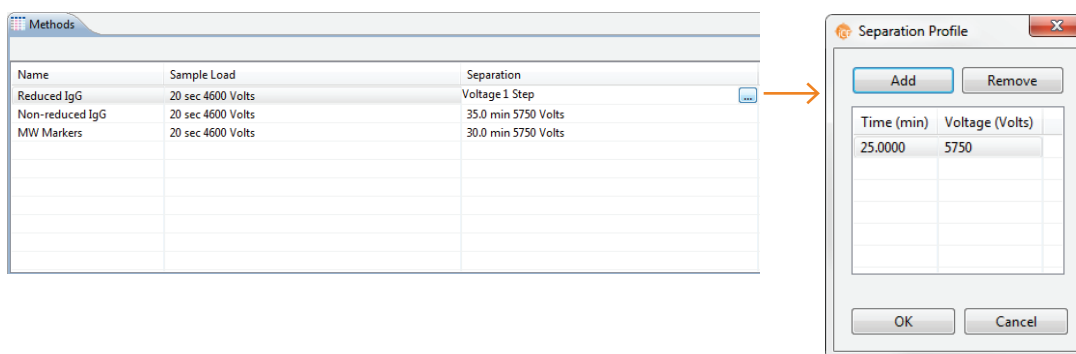
Name	Sample Load	Separation
Reduced IgG	20 sec 4600 Volts	25.0 min 5750 Volts
Non-reduced IgG	20 sec 4600 Volts	35.0 min 5750 Volts
MW Markers	20 sec 4600 Volts	30.0 min 5750 Volts

- b. Click the first cell in the Sample Load column, then click then click the selection button [...] to set your load profile time (in seconds) and voltage.



- c. Click the first cell in the Separation column then click the selection button [...] to set your separation time (in minutes) and voltage.

NOTE: Run your reduced IgG samples and IgG Standard for 25 minutes and the CE-SDS MW Markers for 35 minutes. Run your non-reduced IgG samples and IgG Standard for 35 minutes. The default separation voltage for all sample types is 5750 volts.



9. You can now:

- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.

10. In the Injections pane:

- **To add or change sample names:** Click the **Sample ID** cell for the injection and type a name.

NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 103 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

- **To change injection names:** Click the **Injection Name** cell for the injection and type a name.

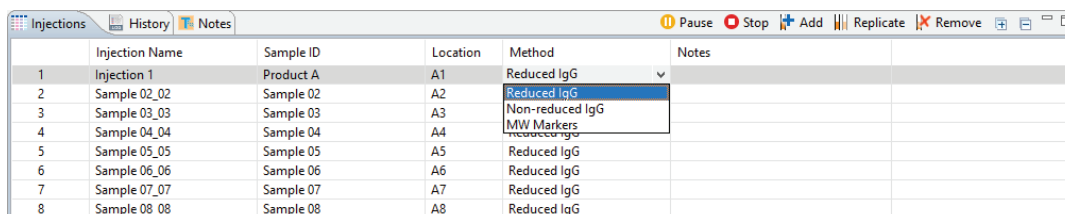
NOTES:

Each injection name must be unique.

Changing the injection name won’t affect the sample ID.

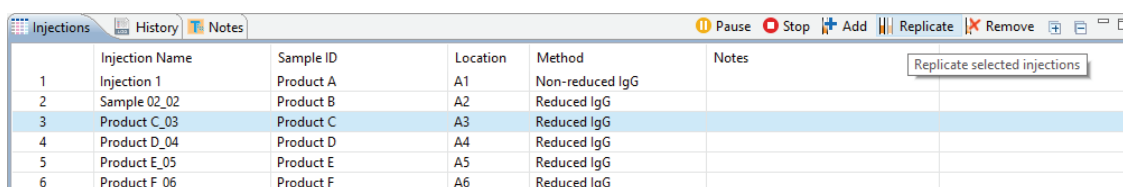
	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

- **To assign methods for each injection:** Click the **Method** cell for the injection and select a method from the drop down menu.

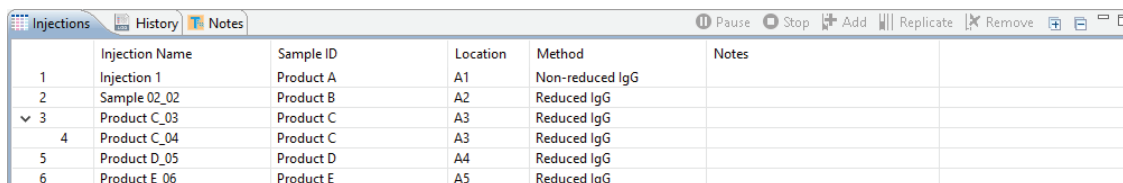


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Non-reduced IgG	
4	Sample 04_04	Sample 04	A4	MW Markers	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

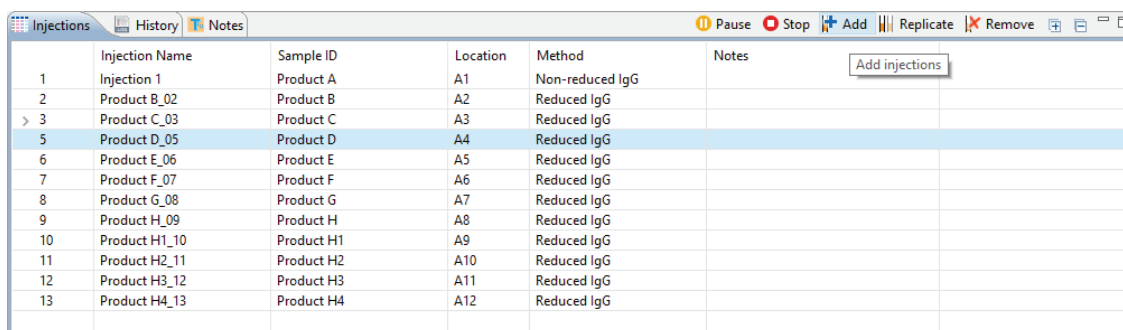


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product E_05	Product E	A5	Reduced IgG	
6	Product F_06	Product F	A6	Reduced IgG	



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product C_04	Product C	A3	Reduced IgG	
5	Product D_05	Product D	A4	Reduced IgG	
6	Product E_06	Product E	A5	Reduced IgG	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.

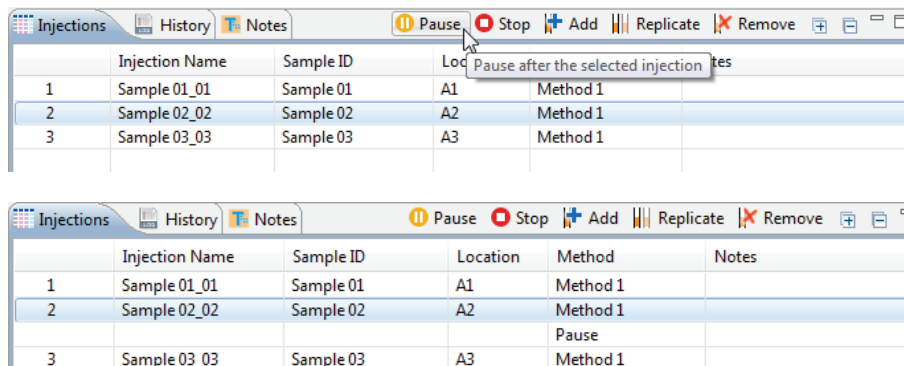


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Product B_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product D_05	Product D	A4	Reduced IgG	
6	Product E_06	Product E	A5	Reduced IgG	
7	Product F_07	Product F	A6	Reduced IgG	
8	Product G_08	Product G	A7	Reduced IgG	
9	Product H_09	Product H	A8	Reduced IgG	
10	Product H1_10	Product H1	A9	Reduced IgG	
11	Product H2_11	Product H2	A10	Reduced IgG	
12	Product H3_12	Product H3	A11	Reduced IgG	
13	Product H4_13	Product H4	A12	Reduced IgG	

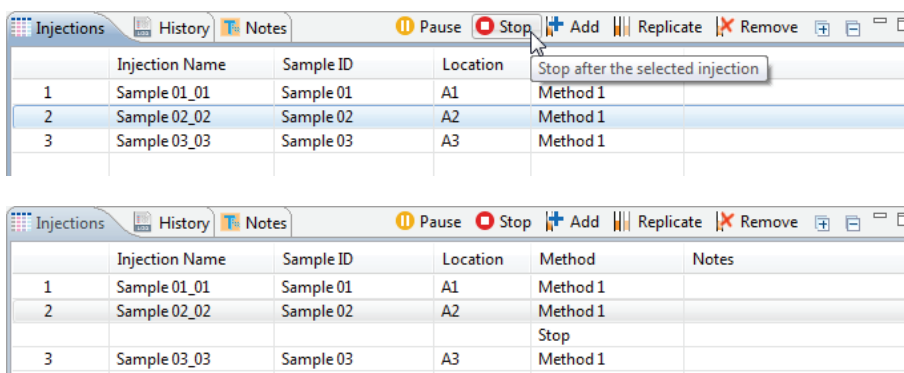
- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

11. Add programmed pauses or stops in the batch to automatically pause or stop after injections (optional).

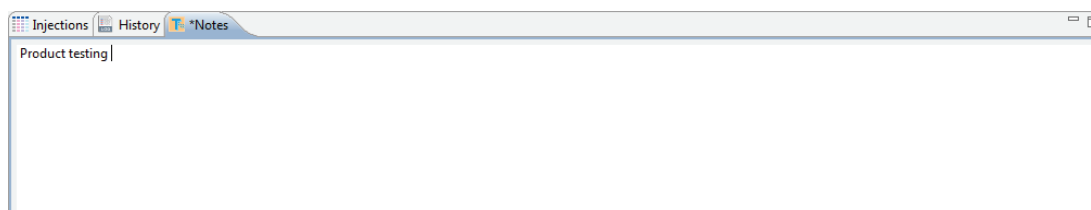
- **To program a pause:** Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete. To resume the batch, click **Continue** in the instrument status bar.



- **To stop the run after a specific injection:** Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.



12. Click on the **Notes** pane, then click in the notes area and type any information you want to add about your batch (optional).



13. You can preset the parameters used to analyze data generated with the batch (optional). We recommend using the default parameters for CE-SDS PLUS applications, but if you want to modify parameters:
- Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:

Default Analysis: Maurice CE-SDS PLUS

Markers

Peak Names
Peak Fit
Advanced
Signal Processing

Analysis Groups

Standards

Add Remove

Apply Default:
Standards

Apply Override:

Apply To	Group
Sample	Standards

Add Remove

Markers

Internal Standard Time 750 Seconds

Markers Injection no markers

MW (kDa)	RMT
10	1
20	1.15
33	1.3
55	1.5
103	1.8
178	2.05
270	2.4

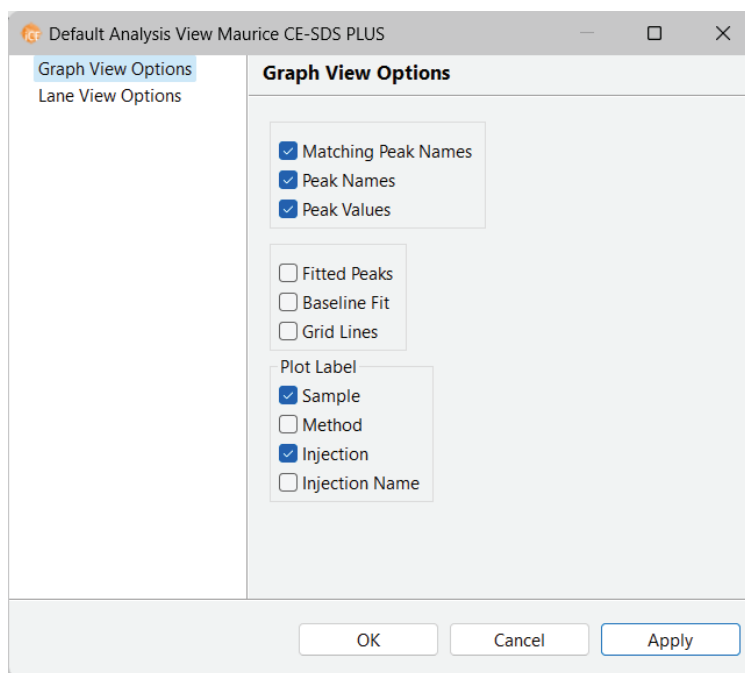
Add Remove

Import... Export... OK Cancel Apply

- Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to “Analysis Settings Overview” on page 713.

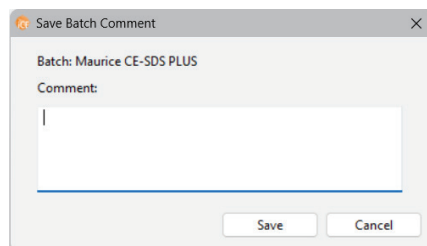
14. You can preset the parameters used to view data when the batch is complete (optional). Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options button:

- a. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



- b. Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to “Customizing the Data Display” on page 695.

15. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



16. Enter a name for your batch then click **Save**.

Step 7: Start the Batch

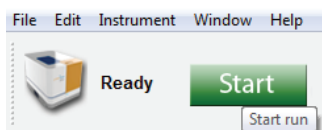
1. Make sure Maurice is connected to Compass for iCE. If he isn't, select **Instrument** and click **Connect**. Then select your Maurice system and click **Connect**.

2. Click **Start** to start your batch.

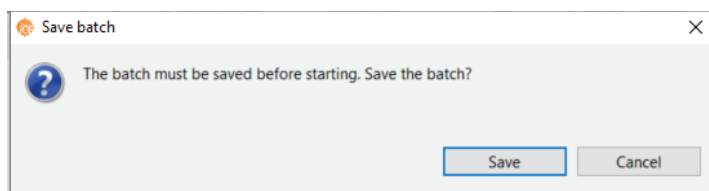
NOTES:

An alert may appear if the wrong cartridge is installed or if a cartridge post-run cleanup or cartridge purge is recommended. See page “Step 5: Check for Cartridge Alerts” on page 300 for more information.

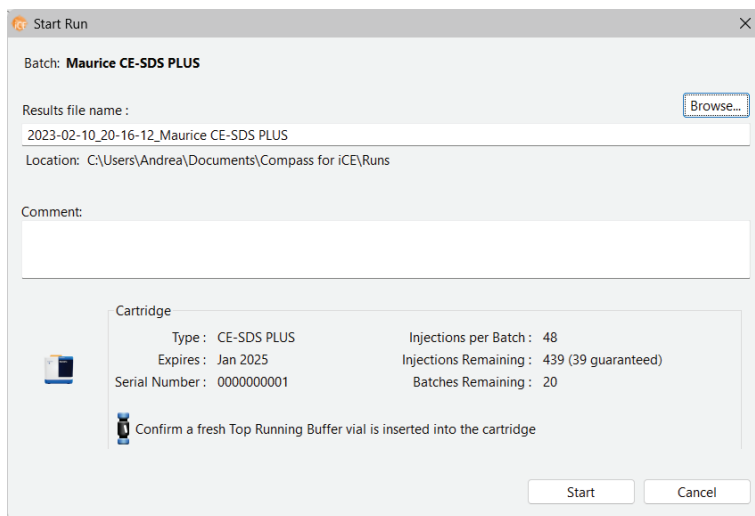
An alert may appear if the wrong adapter is installed or if you are starting a CE-SDS PLUS batch on MauriceFlex and have not locked the reagent vials in place. See “Adapter and Insert Alerts” on page 403 for more information.



3. You will be asked to save your batch before starting the run. Click **Save**.



4. The Start Run box shows you the number of injections and batches remaining on your cartridge. Make sure you still have enough injections left before you start the run. If not, prep a new cartridge.
5. Click in the **Results File name** box if you want to change the default run file name. Otherwise just leave it as is.

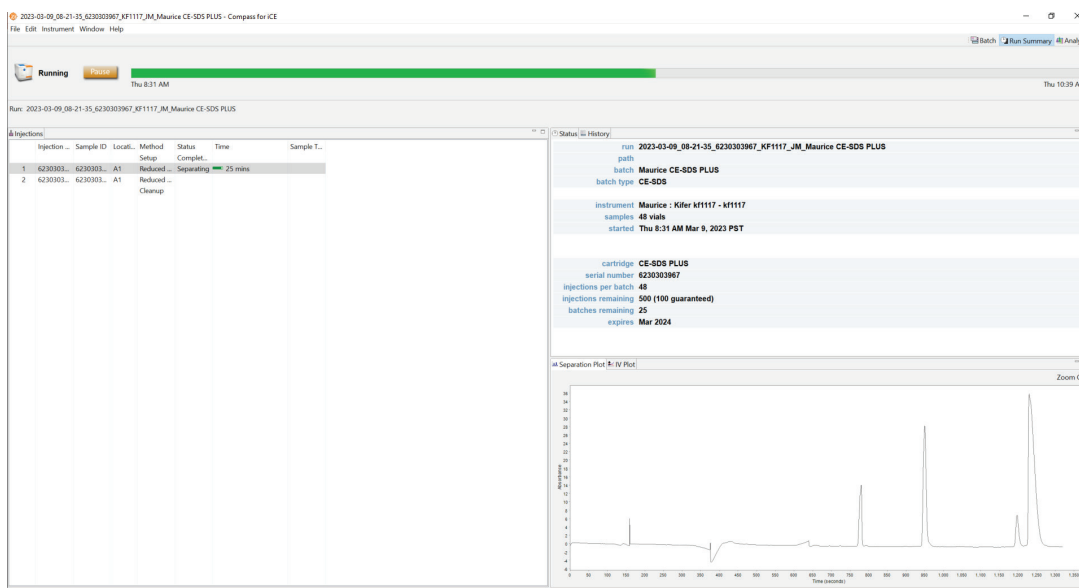


6. If you don't want to save the file to the default Runs folder, click **Browse** to select a different location.

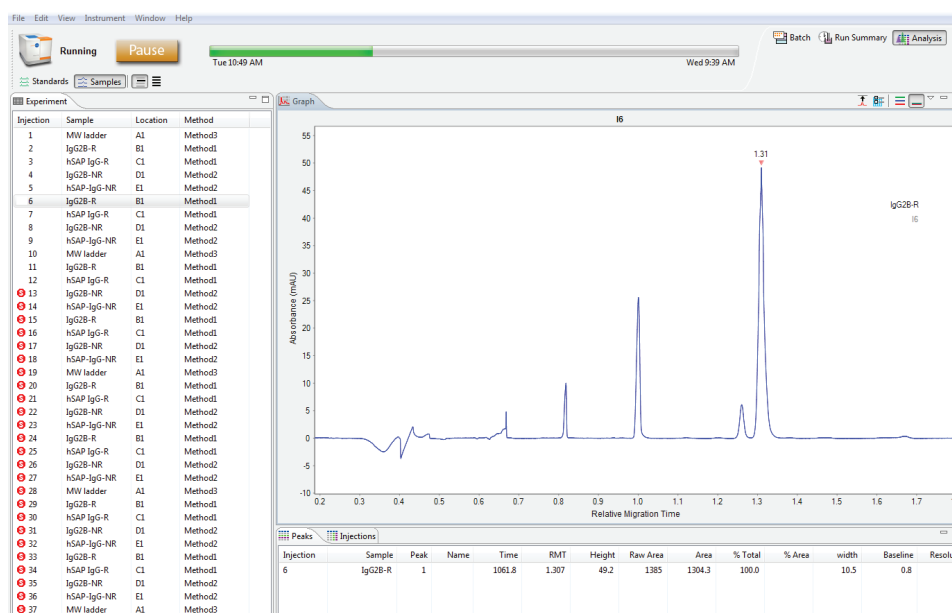
- Enter any run details you'd like in the Comments box (optional).
- Click **Start** to start the run.

NOTE: The indicator light on Maurice's front panel will pulse slowly while he's running the batch.

To monitor the progress of your batch, click the **Run Summary** tab. See Chapter 16: "Run Status" for more details.



To view results, analyze results or change analysis parameters on completed injections while the batch is still running, click the **Analysis** tab. See Chapter 20: "CE-SDS Data Analysis" for more details.



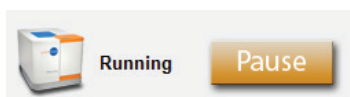
When your batch is complete, you can view electropherograms for all injections in the Analysis tab, and all batch and injection details in the Run Summary tab.

NOTE: If you'd like Maurice to let you know when your run is done, you can set him up to tweet you. Go to "Setting Up Maurice Systems to Send Tweets" on page 764 for more info.

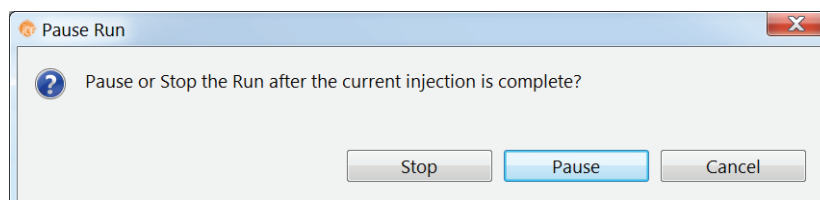
Editing a Running Batch

You can pause a batch after it's started to make updates or add samples and reagents.

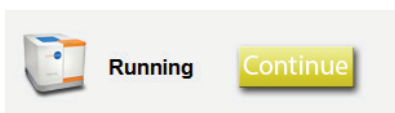
1. Select **Instrument > Pause** or select the **Pause** button.



2. Click **Pause** in the pop-up window.



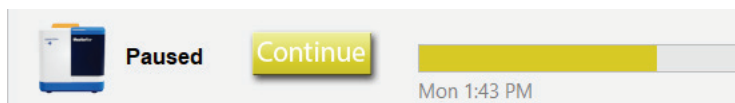
The Pause button will change to Continue but Maurice's status will stay at Running until Compass for iCE completes the current injection or calibration.



- You can now make these edits to the batch:
 - Add sample(s)
 - Add injection(s) to the end of the batch
 - Modify injections that haven't completed and aren't greyed out
 - Create a new method
 - Update the method. Once a method is run it can only be updated. Highlight the method you want to edit and click **Update** in the Methods pane. The software will copy the parameters into a new method that you can modify.

3. When Maurice completes the current injection, his status will change to Paused and the progress bar will turn yellow.

NOTE: Maurice's status light will change from a pulsing blue to a blue pulse with magenta flashes which means he's paused the run.



The status in the Injection pane also changes to Paused and the pause duration updates live. You can now open the instrument door to:

- Add new samples

IMPORTANT

The door must be closed within 15 minutes of opening it to avoid potential assay performance issues.

NOTE: If the door is open for more than 30 seconds, a white light near the cartridge will start to flash and a warning message displays in Compass for iCE until the door is closed.

4. To continue the batch, click **Continue**. Changes you made to the batch are saved and Maurice will continue with the next sample injection.

NOTE: The D2 lamp is turned off for safety reasons when the door is opened. Once the batch is continued, Maurice will pause for seven additional minutes before the next injection to let the lamp re-stabilize. During that time, the batch status in Compass for iCE will display injection loading.

Post-batch Procedures

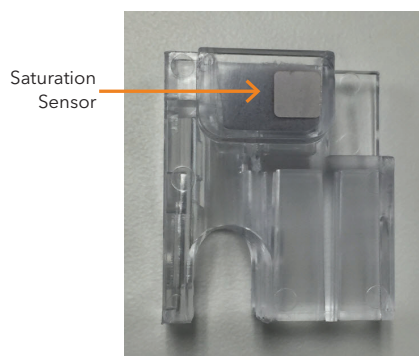
When the batch is done:

1. Open Maurice's door. The lights on either side of the cartridge slot will be **orange** as Maurice will have already disengaged the cartridge.
2. If you're using the optional Maurice Filter Kit to contain β ME odor, cover the 96-well sample plate immediately with a plate lid.
3. Remove your samples. Leave the Water (P3), Wash Solution (P5), and Air (P6) vials in place if your cartridge still has injections left since they will be needed for the cartridge post-run cleanup step. Discard the remaining reagent vials.
4. Gently pull out the cartridge making sure the cartridge inlet doesn't come into contact with any surface.

NOTE: If you see any separation matrix sticking to the capillary inlet, soak it in DI water for 5 minutes. Then wipe it using a lint-free laboratory wipe that's been moistened with DI water.



5. Pull the cartridge insert out.
6. Remove the Top Running Buffer vial and dispose of it according to your institution's safety and waste disposal guidelines.
7. Check the saturation sensor on the back of the cartridge insert. If it's red, you'll need to use a new cartridge insert for your next batch. If the saturation sensor isn't red, you can keep using the current cartridge insert with that cartridge.



NOTE: Don't dispose of the cartridge insert unless the saturation sensor is red.

If you're at 100 injections, you've reached the limit of guaranteed performance for the CE-SDS PLUS cartridge. To dispose of a finished cartridge, put it in its original packing and discard it along with the cartridge insert and the Top Running Buffer vial per your institution's safety and waste disposal guidelines. Discard the cleaning vial you've used with that cartridge too.

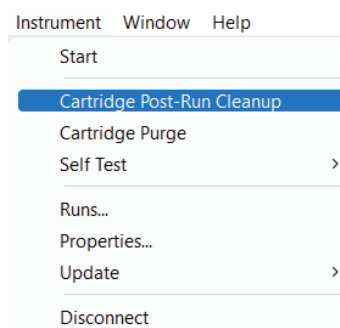
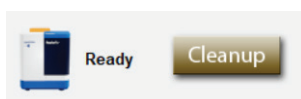
NOTE: Disposal of cartridges, sample plates and vials depends on the samples you ran. If you aren't sure what your sample origins are, we recommend you dispose of everything in biohazard waste.

If the cartridge will be used again, clean up and store the cartridge:

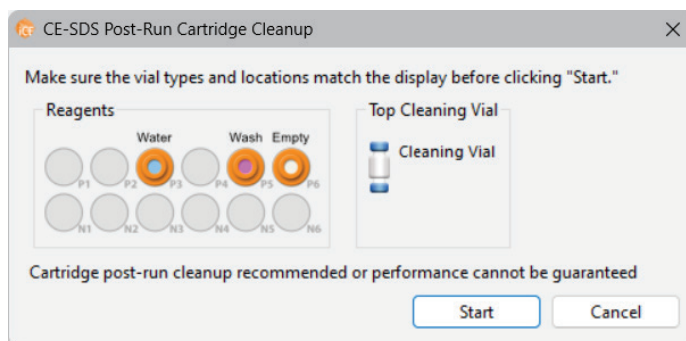
- Verify that there is 1.5 mL of DI water in the Water (P3) vial, 1.5 mL of Wash Solution in the Wash Solution (P5) vial, and air in the empty (P6) vial.
- Insert a Cleaning Vial into the cartridge insert.



- Slide the cartridge insert back into the cartridge.
- Insert the cartridge in Maurice.
- Click on the brown **Cleanup** button in the Instrument Status Bar or select **Instrument** and **Cartridge Post-Run Cleanup** in the Compass main menu.



- You'll get the following message. Click **Start**. It'll only take six minutes.



- g. Once the cleanup procedure is done, discard the reagent vials and remove the cartridge.
- h. Pull the insert from the cartridge.
- i. Remove the Cleaning Vial and push the empty insert back into the cartridge.

NOTE: The cleaning vial is paired with the cartridge and can be used for a maximum of five Cartridge Cleanup cycles of that cartridge. Dispose of the cleaning vial when you dispose of the cartridge. Don't use it with other cartridges.

- j. Put the cartridge back in its protective packaging and store it at room temperature.

!WARNING! SHARPS HAZARD

The capillary inlet of the cartridge may present a potential sharps hazard. Dispose of used cartridges according to your organizations health and safety regulations.



!WARNING! BIOHAZARD

Cartridges, sample plates and vials should be handled by procedures recommended in the CDC/NIH manual: Biosafety in Microbiological and Biomedical Laboratories (BMBL). The manual is available from the U.S. Government Printing Office or online at <http://www.cdc.gov/biosafety/publications/bmbl5/>.

Depending on the samples used, the cartridges, plates and vials may constitute a chemical or a biohazard. Dispose of the cartridges, plates and vials in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations. Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste content before you store, handle, or dispose of chemical waste.

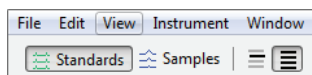
Checking Your Data

Compass for iCE detects your sample proteins, CE-SDS MW Markers and Internal Standard peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

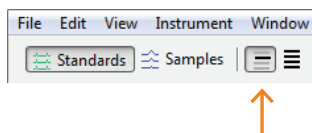
Step 1: Check Your Internal Standard

To make sure your Internal Standard is identified correctly:

1. Go to the **Analysis** screen and open your run (if it isn't already open).
2. Click **Standards** in the View bar.

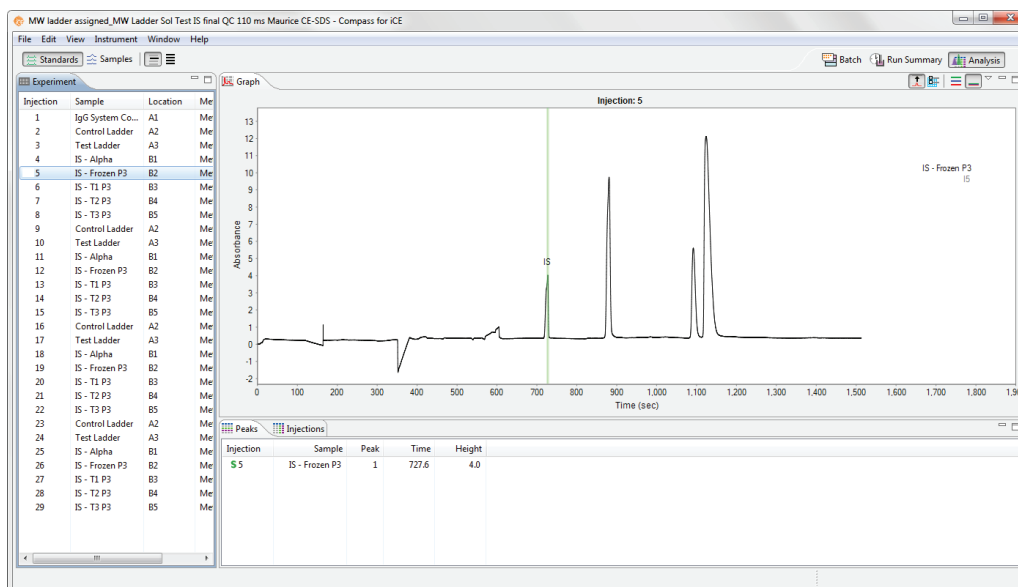


3. Click the **View Selected** icon in the View bar.



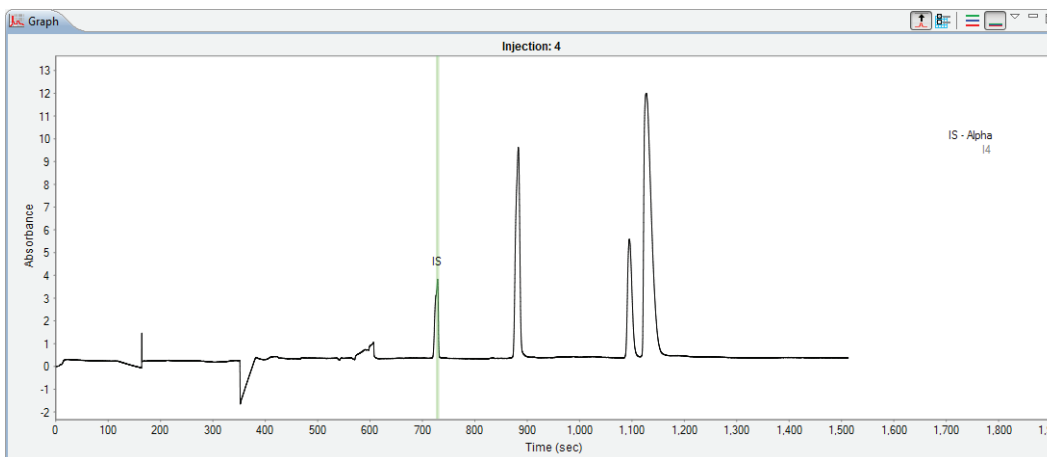
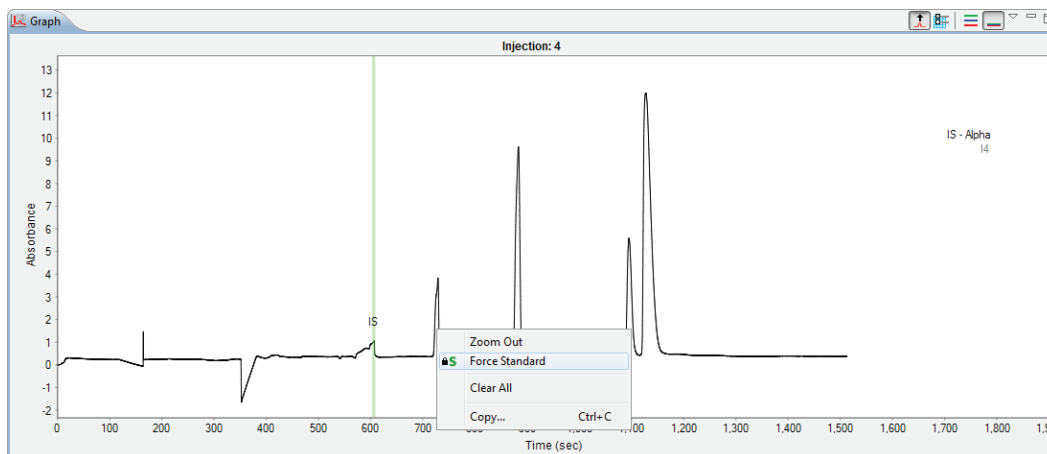
4. Click **Injection 1** in the Experiment pane.

5. Check that your Internal Standard in the electropherogram has been correctly identified. It'll be labeled **IS** and will have a green vertical line running through it. The Internal Standard is also identified with an **S** in the Peaks table.



6. If your Internal Standard isn't identified correctly, here's how to manually correct it:

To set an unidentified peak as the Internal Standard: Right-click the peak in the electropherogram or Peaks table and select **Force Standard**. Compass will assign that peak as the Internal Standard.



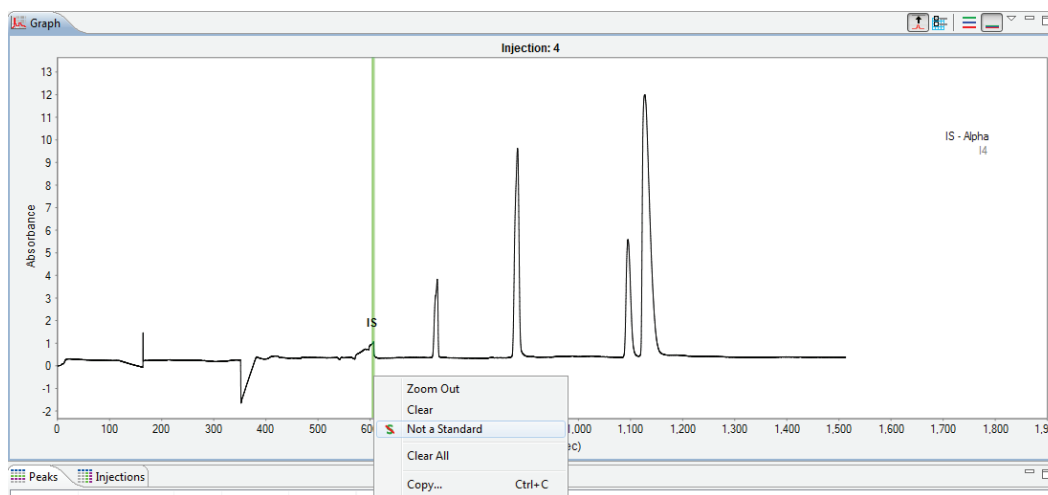
A lock icon indicating the Internal Standard was set manually will display next to the peak in the Peaks table and a check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Peaks				
Injection	Sample	Peak	Time	Height
1	Sample 1	11	548.2	115.1
1	Sample 1	12	557.7	115.2
1	Sample 1	13	572.0	134.0
1	Sample 1	14	583.5	149.7
1	Sample 1	15	590.6	190.0
1	Sample 1	16	710.8	230.3
1	Sample 1	17	714.6	278.5

Experiment			
Injection	Sample	Location	Method
1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1

NOTE: To remove an Internal Standard peak assignment that was made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**. To clear all the manual settings for the injection, click **Clear All**.

If an incorrect peak is identified as the Internal Standard: Right-click the peak in the electropherogram or Peaks table and select **Not a Standard**.



A check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Experiment			
Injection	Sample	Location	Method
✓ 1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1

- Repeat the previous steps for all other injections to make sure your Internal Standard is identified correctly.

Step 2: Set Your Molecular Weight (MW) Markers

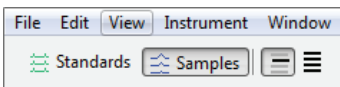
NOTE: You'll only need to do this if you ran the CE-SDS MW Markers. If you didn't, you can skip to the next section.

Compass reports the relative migration time (RMT) of your sample in the Peaks table. If you also want to know the relative molecular weight of your sample, you can run the CE-SDS MW Markers as one of your injections.

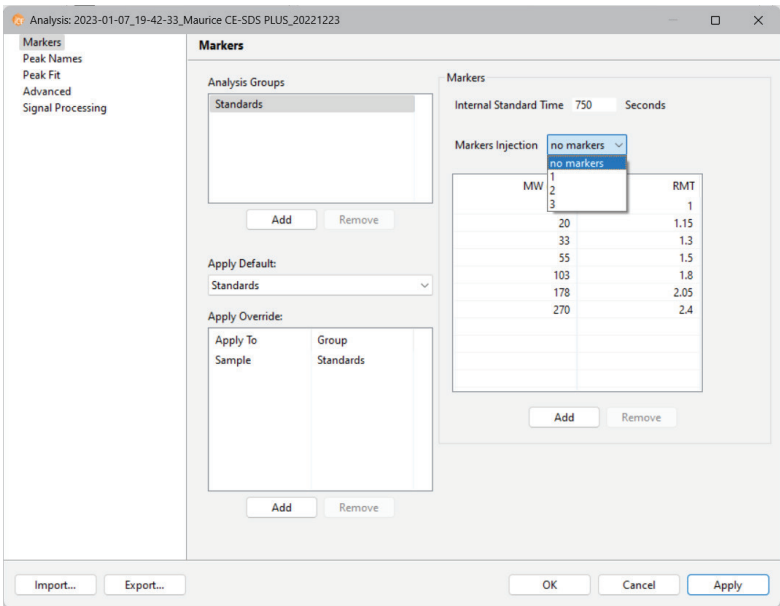
You'll see these sizing markers when you run the CE-SDS MW Markers: 10, 20, 33, 55, 103, 178, and 270 kDa.

To get MW data:

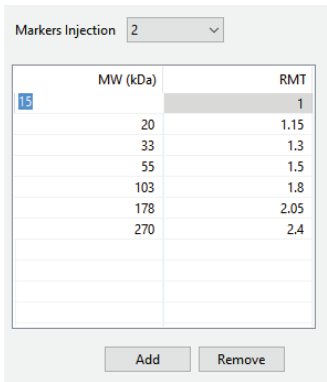
1. Click **Samples** in the View bar.



2. Select **Edit** from the main menu and click **Analysis**. In the Analysis window, select **Markers** in the left sidebar. Then click the **Markers Injection** drop down menu to select the injection you ran your CE-SDS MW Markers in.

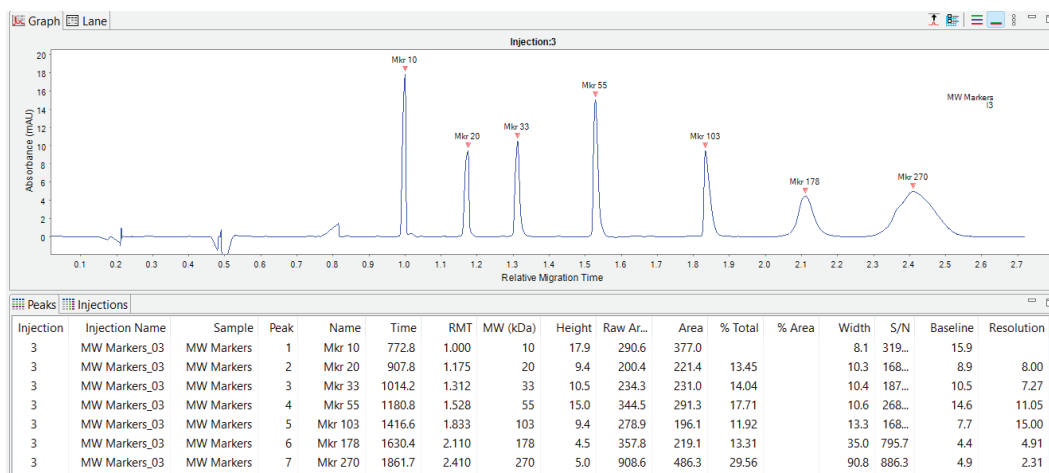


3. The default Maurice CE-SDS MW Markers molecular weights and relative migration time (RMT) values are already populated in the table. If you'd like to use these values, skip to the next step. If you're using different markers, click in the MW and RMT cells to type new values, click a row and select **Remove** to delete, or click **Add** to add a new one.



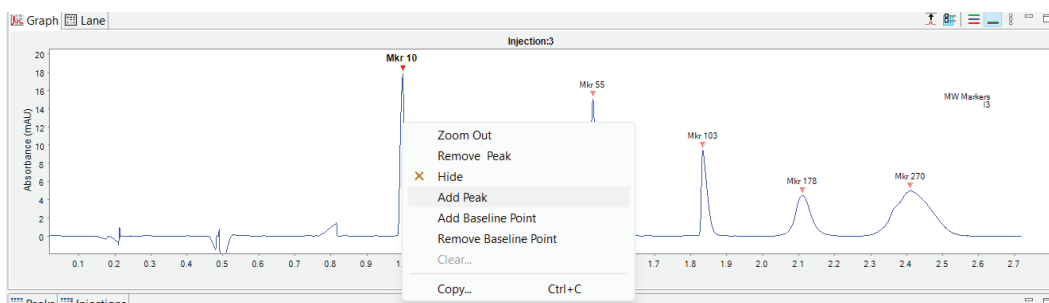
- Click **OK** to close the Analysis window. Compass will automatically assign the molecular weights to your markers and label them Mkr. A MW (kDa) column will also now display in the Peaks table.

NOTE: The Mkr 10 peak is also the Internal Standard in every sample.



- It's always a good idea to verify that all your CE-SDS MW Markers are identified correctly. Here's how to manually correct them:

To set an unidentified peak as a MW Marker: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will assign that peak as a MW Marker, and correctly reassign the remaining marker peaks.

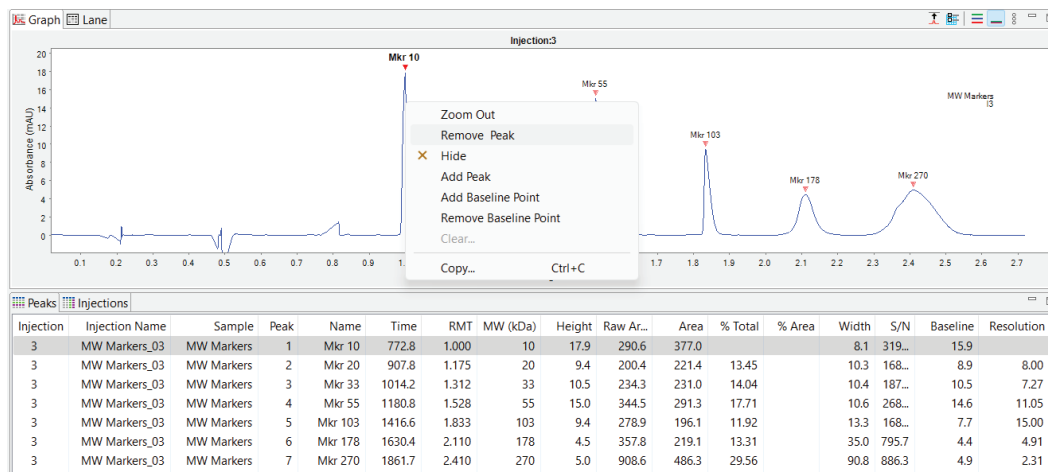


A check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Experiment			
Injection	Sample	Location	Me
1	IgG System Co...	A1	Me
✓ 2	Control Ladder	A2	Me
3	Test Ladder	A3	Me
4	IS - Alpha	B1	Me

NOTE: To remove MW Marker peak assignments that were made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**.

If an incorrect peak is identified as a MW Marker: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass should correctly reassign the remaining peaks as markers and update the Peaks table.



A check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

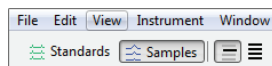
Experiment			
Injection	Sample	Location	Me
1	IgG System Co...	A1	Me
✓ 2	Control Ladder	A2	Me
3	Test Ladder	A3	Me
4	IS - Alpha	B1	Me

Step 3: Checking Sample Peaks

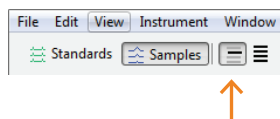
All detected peaks will be labeled automatically with the RMT (default) or apparent MW (if the CE-SDS MW Markers were run).

To make sure your sample proteins are identified correctly:

1. Click **Samples** in the View bar.



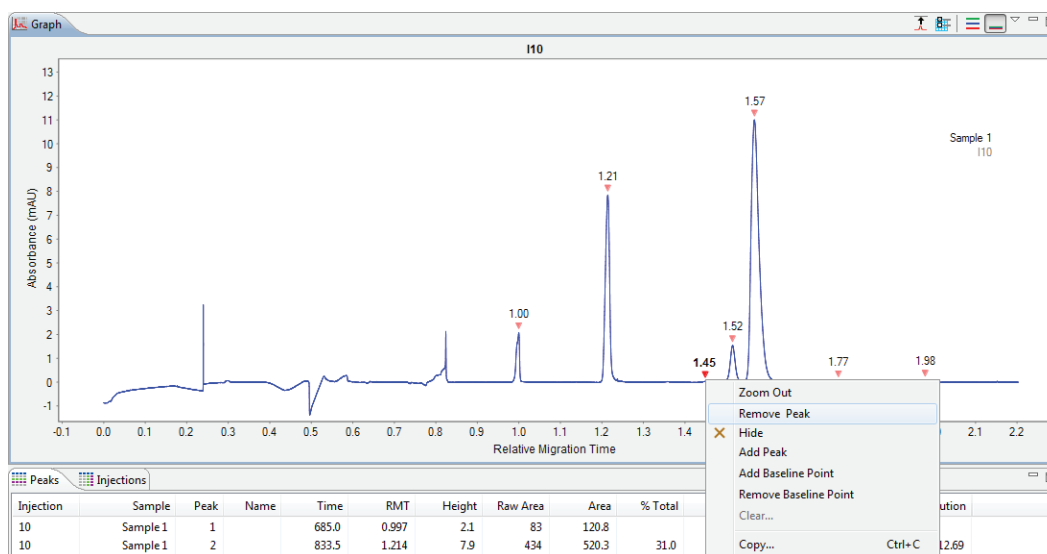
2. Click the **View Selected** icon in the View bar.



3. Click **Injection 1** in the Experiment pane.

4. If your sample peaks aren't identified correctly, here's how to manually correct them:

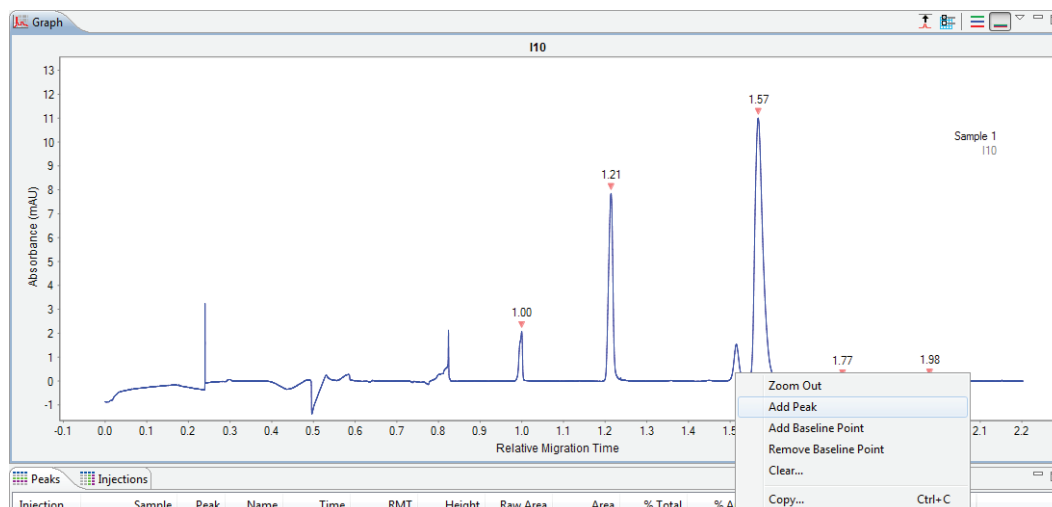
If a peak is incorrectly identified as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass will no longer identify it as a sample peak in the electropherogram and the peak data will be removed in the results table.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment			
Injection	Sample	Location	Method
1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1
8	Sample 1	A1	Method1
9	Sample 1	A1	Method1
✓ 10	Sample 1	A1	Method1
11	Sample 1	A1	Method1
12	Sample 1	A1	Method1

To identify a peak as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Injection	Sample	Location	Method
1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1
8	Sample 1	A1	Method1
9	Sample 1	A1	Method1
✓ 10	Sample 1	A1	Method1
11	Sample 1	A1	Method1
12	Sample 1	A1	Method1

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear**.

- Repeat the previous steps for the remaining injections to make sure all sample peaks are correctly identified.

Step 4: Assigning Peak Names

Compass can also optionally identify and automatically name sample peaks using user-specified peak name settings. For more information on how to do this, see “Manual Peak Integration” on page 739.

Chapter 13:

Turbo CE-SDS Batches

Chapter Overview

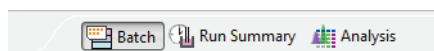
- Overview
- Batch Screen Overview
- Opening a Batch
- Creating a New Batch
- Viewing Replicate Injections
- Batch History
- Making Changes to a Batch
- Viewing and Editing Batches in Completed Runs
- Copying and Pasting Injection Names or Sample IDs from other Documents
- Importing and Exporting Injections
- Importing and Exporting Methods
- Batch Reports

Overview

Maurice Turbo CE-SDS batches can be run on Maurice, Maurice S., or MauriceFlex systems using a Turbo CE-SDS cartridge.

Batch Screen Overview

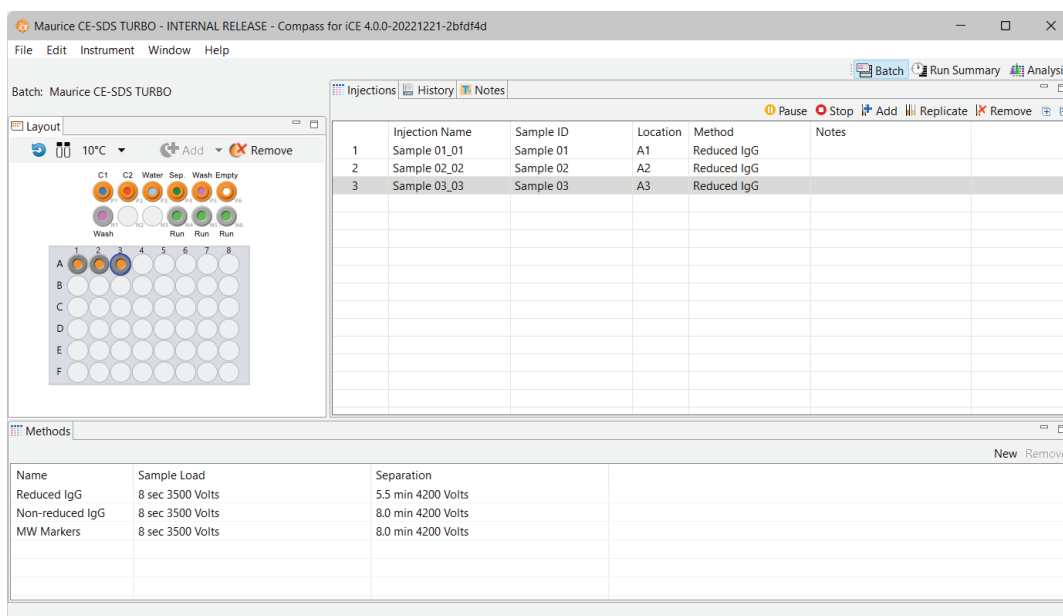
You can use the Batch screen to create, view, and edit batches. To get to this screen, click the **Batch** screen tab:



Batch Screen Panes

The Batch screen has five panes:

- **Layout** - Displays a map of either the 96-well plate or 48-vial tray of batch sample locations.
- **Injections** - Lists the injections, sample ID, sample locations and methods that Maurice or Maurice S. will execute for each sample in the batch.
- **History** - Lists all batch file events from initial creation to the most current update.
- **Notes** - Lets you enter specific information about your batch.
- **Methods** - Lets you create methods and enter method parameters used in the batch.



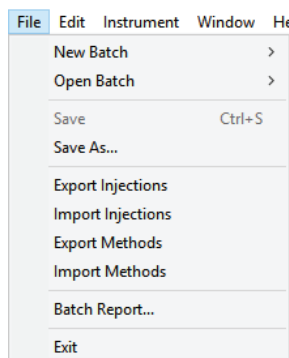
Software Menus Active in the Batch Screen

These main menu items are active in the Batch screen:

- File
- Edit
- Instrument (when the software is connected to Maurice, Maurice S. or MauriceFlex)
- Window
- Help

File Menu

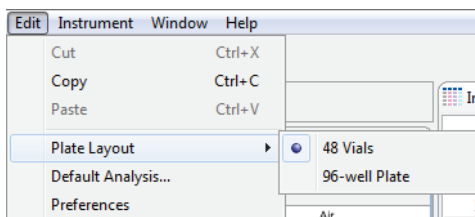
These File menu options are active:



- **New Batch** - Creates a new batch from a starter template.
- **Open Batch** - Opens an existing batch.
- **Save/Save As** - Saves the open batch.
- **Export Injections** - Exports injections from the current batch as a .csv file.
- **Import Injections** - Imports injections into the current batch from a .csv file.
- **Export Methods** - Exports method(s) from the current batch as separate files.
- **Import Methods** - Imports saved method(s) into the current batch.
- **Batch Report** - Exports a table of sample and method details for each injection in the batch as a PDF file.
- **Exit** - Closes Compass for iCE.

Edit Menu

The following Edit menu options are active in the Batch screen:



- **Cut** - Cuts the information currently selected.
- **Copy** - Copies the information currently selected.
- **Paste** - Pastes the copied information.

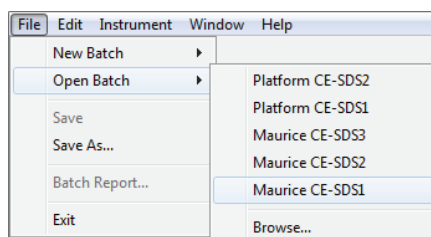
NOTE: Cut, Copy and Paste work differently depending on the pane you're in. How and when you can use them is described throughout the chapter.

- **Plate Layout** - Lets you select between a 96-well plate or a 48-vial tray to run samples in.
- **Default Analysis** - Displays the default settings that will be used to analyze the data generated with your batch.
- **Default Analysis View** - Displays the default settings that will be used to view the data generated with your batch.
- **Preferences** - Lets you set and save your preferences for Access Control, data export, graph colors, grouped data and Twitter settings. See Chapter 21, "Setting Your Preferences" for more information.

Opening a Batch

To open an existing batch:

1. Select **File** in the main menu and click **Open Batch**.



2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file. Select one of those or click **Browse** to open the Batches folder and select a different one.

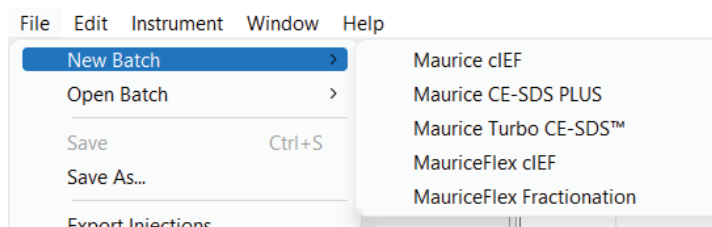
3. To make changes to the batch, see the steps in “Creating a New Batch”. When you’re done, select File from the main menu and click Save.

Creating a New Batch

To create a new batch, we recommend that you use one of the template methods and make any necessary modifications from there.

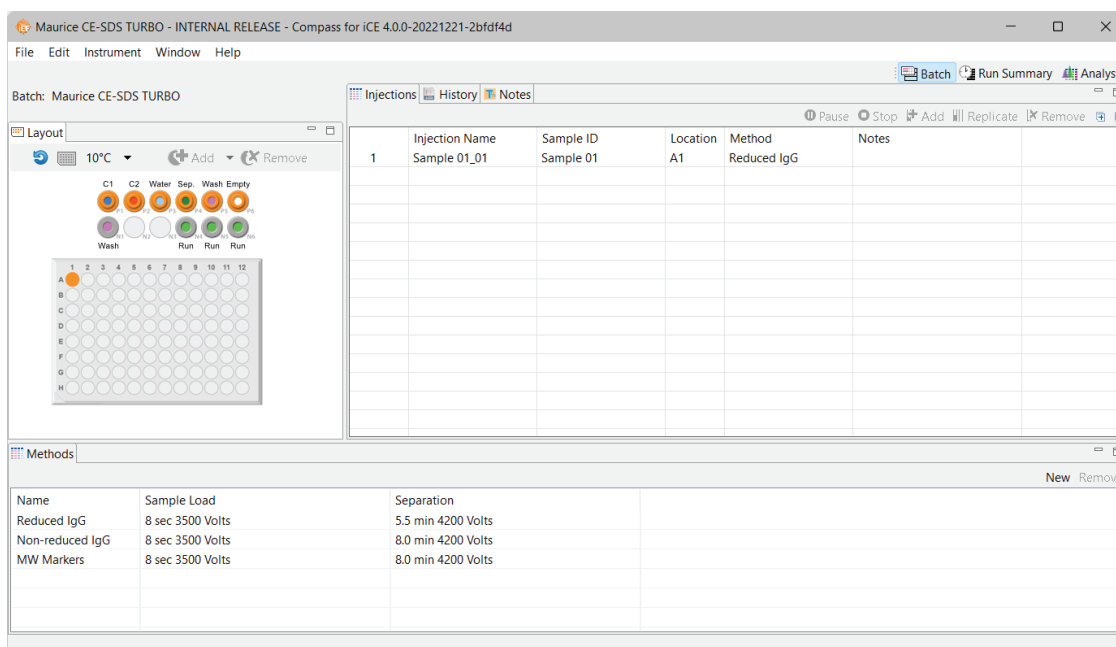
Step 1 - Open a Template Batch

1. Select **File** in the main menu and click **New Batch**:



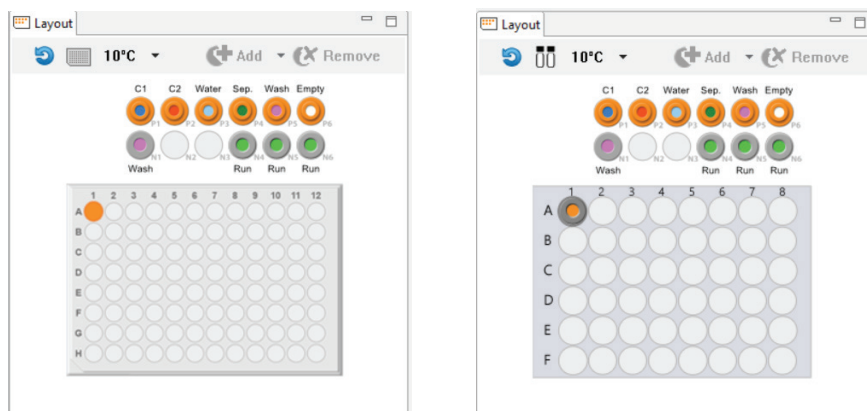
NOTE: Only template batches specific to your Maurice will be available when you are connected to the instrument. Template batches will not appear in the menu if that application type cannot be performed on your instrument.

2. Select **Maurice Turbo CE-SDS**. A batch using the default method will display.



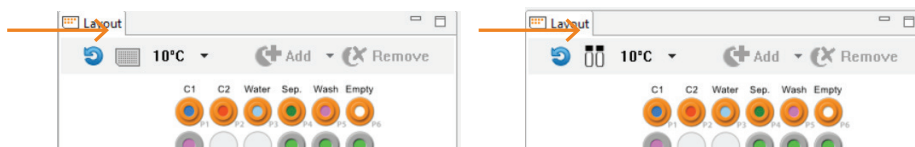
Step 2 - Assign Your Samples

The Layout pane shows the default locations of where batch reagent should be placed in your Maurice system's samples and reagents platform, and a map of either the 96-well plate or the 48-vial tray used for samples.



The same reagent locations are used for every batch:

- **P1** - Conditioning Solution 1 with **orange pressure cap**
 - **P2** - Conditioning Solution 2 with **orange pressure cap**
 - **P3** - DI water with **orange pressure cap**
 - **P4** - Separation Matrix with **orange pressure cap**
 - **P5** - Wash Solution vial with **orange pressure cap**
 - **P6** - Empty vial (air) with **orange pressure cap**
 - **N1** - Wash Solution vial with **clear screw cap**
 - **N4** - Running Buffer - Bottom with **clear screw cap**
 - **N5** - Running Buffer - Bottom with **clear screw cap**
 - **N6** - Running Buffer - Bottom with **clear screw cap**
1. To assign samples, select a 96-well plate or 48 vials depending on what you're running. Clicking on the vial/plate icon toggles between formats.



2. Select your samples:

To import samples using a saved injections file:

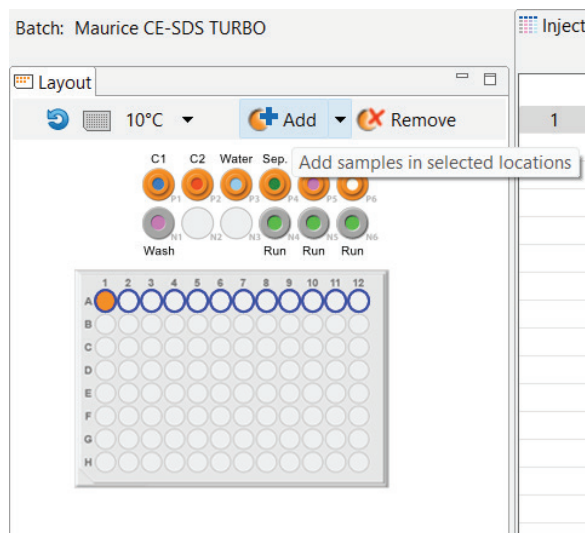
- Select **File** in the main menu and click **Import Injections**.
- Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.
- Skip to step 3 on page 336.

NOTE: For more information on setting up an injections file for import, see “Importing and Exporting Injections” on page 350.

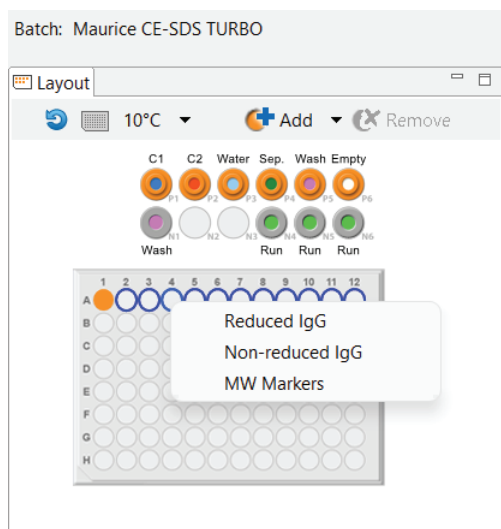
To select samples manually:

- **Add samples and select methods later:** Use your mouse to highlight the well or vial positions your samples are located in, then click **Add**. For this example we’re using a 96-well plate.

NOTE: The template batch automatically adds a sample in well or vial A1 by default, but if you don’t have a sample in this position you can remove it after you’ve added new positions for your samples.



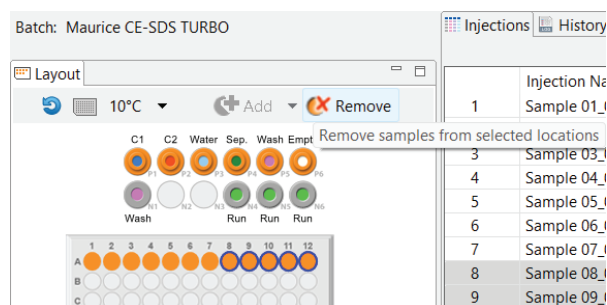
- **Add samples with preassigned methods:** Highlight the well or vial positions your samples are located in, then right-click and select a method.



Either option of adding samples populates the Injections table:

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Sample 1	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

To remove samples, just highlight the ones you want to remove and click **Remove**. They'll also be removed in the Injections table. You can also right-click well(s) or vial(s) in the Layout pane and click **Remove Sample(s)**.

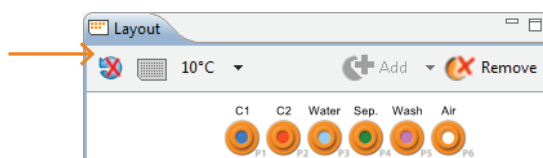


3. Compass for iCE can monitor the current during a separation for you, stop it if an abnormal current profile is detected and reinject the sample for you automatically. Reinjections are enabled by default, but if you don't want to use this feature just click the reinject icon to toggle it off.

NOTES:

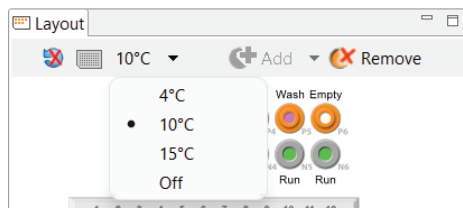
If a sample was reinjected, the injection row will be flagged in the Injections pane in the Run Summary screen. See "Injection Flags" on page 414 for more info.

A maximum of 10 reinjections are allowed per batch.



4. The sample platform is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The two reagent rows are kept at room temperature.



Step 3 - Assign Your Method Parameters

NOTE: There are three default methods. We recommend using the default method parameters for the listed sample types. Please see the [Maurice CE-SDS Application Guide](#) for more information on method optimization.

The Methods pane lists all of the methods used in a batch. You can add and remove methods here and change method parameters.

To import a saved method:

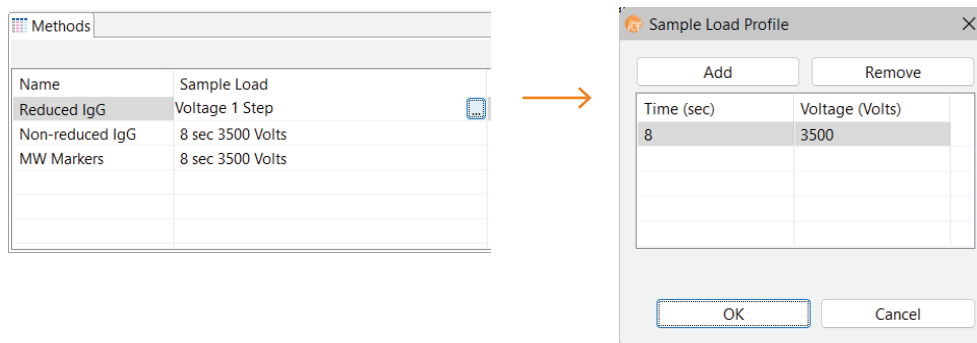
1. Select **File** in the main menu and click **Import Method**.
2. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

To create a new method or make changes to an existing one:

1. In the Methods table, click the first cell in the Name column and enter a new method name if needed.

Methods			
Name	Sample Load	Separation	
Reduced IgG	8 sec 3500 Volts	5.5 min 4200 Volts	
Non-reduced IgG	8 sec 3500 Volts	8.0 min 4200 Volts	
MW Markers	8 sec 3500 Volts	8.0 min 4200 Volts	

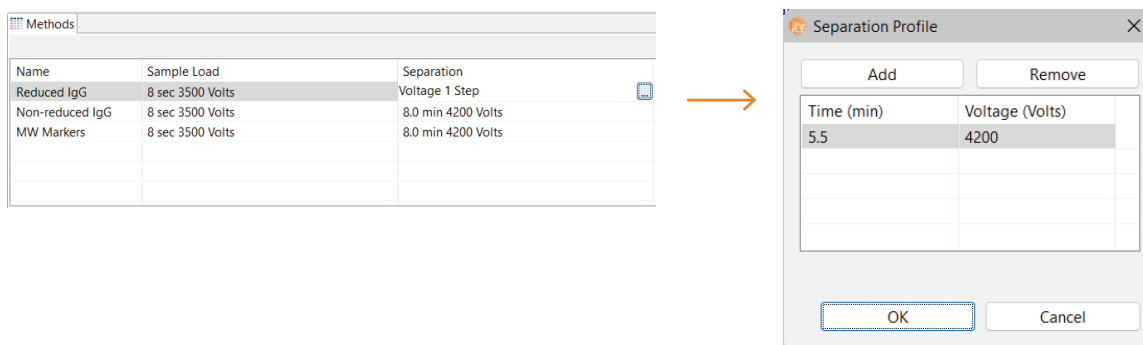
2. Click the first cell in the Sample Load column, then click the selection button [...] to set your sample load profile time (in seconds) and voltage.



- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
- **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a load time (in seconds) and voltage value.
- **To remove a profile step:** Select the row you want to remove and click **Remove**.

- Click the first cell in the Separation column the selection button [...] to set your separation profile parameters (in minutes) and voltage.

NOTE: Run your reduced IgG samples for 5.5 minutes and your non-reduced IgG samples, IgG Standard and CE-SDS MW Markers for 8.0 minutes. The default separation voltage for all sample types is 4200 volts.



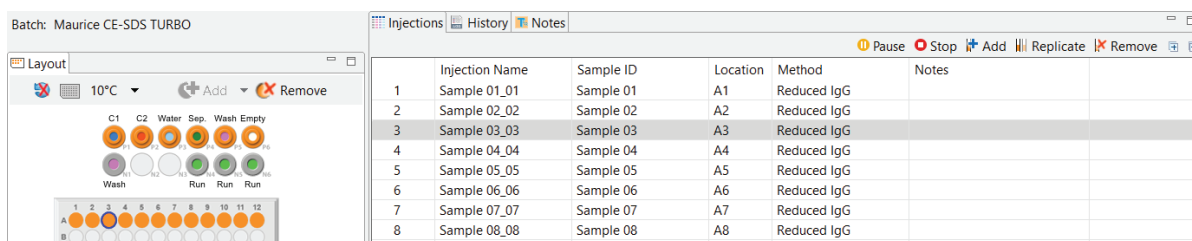
- **To change time and voltage parameters:** Just click in a cell under **Time** or **Voltage** and type the new value(s).
 - **To add a profile step:** Click **Add**. A new row will be added in the table. Then just type in a load time (in seconds) and voltage value.
 - **To remove a profile step:** Select the row you want to remove and click **Remove**.
- You can now:
 - Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.
 - Make updates to the remaining methods by repeating the prior steps.
 - Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.

Step 4 - Set Up Your Injections

The Injection pane lists all sample injections for the batch. Any samples that were added in *Step 2 - Assign Your Samples* are automatically added to this list.

NOTE: If you imported injections in Step 2, you can update the injections table and assign methods as needed using the information that follows. Otherwise skip to “Step 5 - Add Programmed Pauses and Stops (Optional)” on page 341.

Selecting an injection in the table also selects the sample the injection will be made from in the plate or vial map in the Layout pane and vice-versa.



1. To add sample names, click the **Sample ID** cell for the injection and type a name.

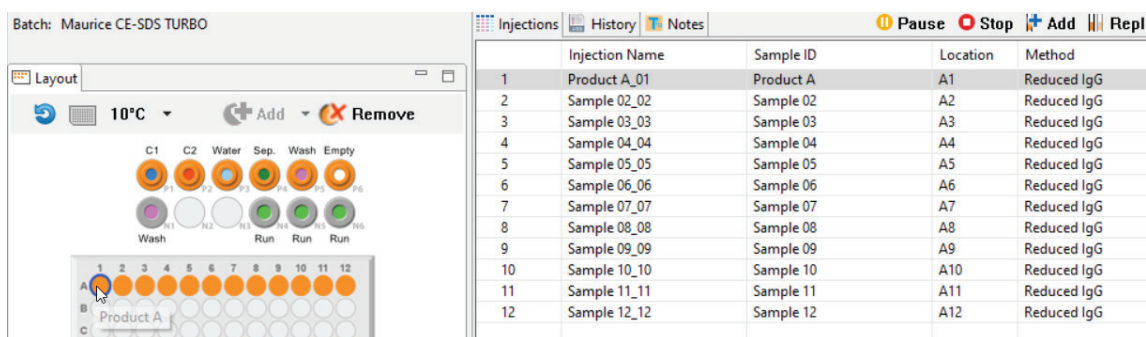
NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 348 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.

	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

The sample name also displays when you hover the mouse over the sample in the plate or vial map:



- Injection names are automatically set to 'Sample ID_injection number'. To change the name, click the **Injection Name** cell for the injection and type a name. Each injection name must be unique. If they aren't when you save the batch, Compass for iCE highlights any non-unique injection names so you can change them.

	Injection Name	Sample ID	Location	Method	Notes
1	Injection	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

NOTE: Changing the injection name won't affect the sample ID.

- Assign your methods to each injection. Just click the **Method** cell for the injection and select a method from the drop down menu. If you added your samples with pre-assigned methods and don't need to make any changes, skip to the next step.

	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	

Hovering over a method name displays the method parameters:

	Injection Name	Sample ID	Location	Method	Notes
1	Product A_01	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	

4. You can add or remove injections as needed using a few different options. If you don't need to make any changes here, just skip to the next step.

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product E_05	Product E	A5	Reduced IgG	
6	Product F_06	Product F	A6	Reduced IgG	

	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product C_04	Product C	A3	Reduced IgG	
5	Product D_05	Product D	A4	Reduced IgG	
6	Product E_06	Product E	A5	Reduced IgG	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.

	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Product B_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product E_05	Product E	A5	Reduced IgG	
6	Product F_06	Product F	A6	Reduced IgG	
7	Product G_07	Product G	A7	Reduced IgG	
8	Product H_08	Product H	A8	Reduced IgG	
9	Product H1_09	Product H1	A9	Reduced IgG	
10	Product H2_10	Product H2	A10	Reduced IgG	
11	Product H3_11	Product H3	A11	Reduced IgG	
12	Product H4_12	Product H4	A12	Reduced IgG	
13	Product H4_13	Product H4	A12	Reduced IgG	

- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

Step 5 - Add Programmed Pauses and Stops (Optional)

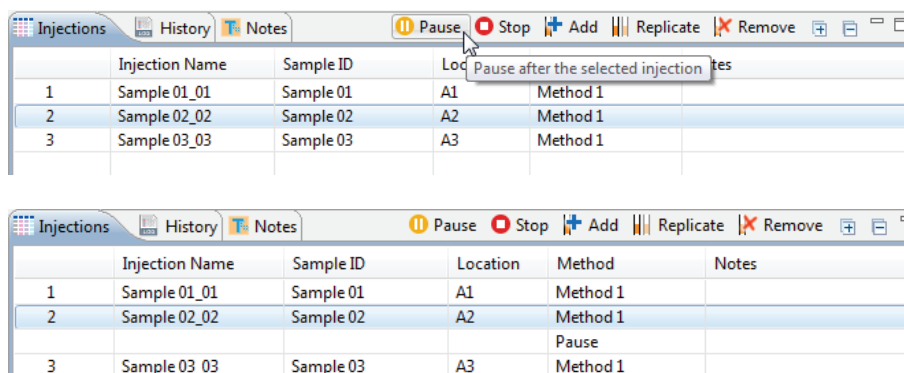
You can program pauses and stops in the batch to automatically pause or stop after injections even when you aren't in front of Maurice.

To program a pause:

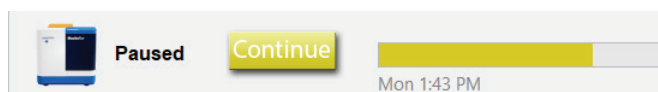
Pauses are helpful when you want to review injection results before continuing with the rest of the batch.

1. Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete.

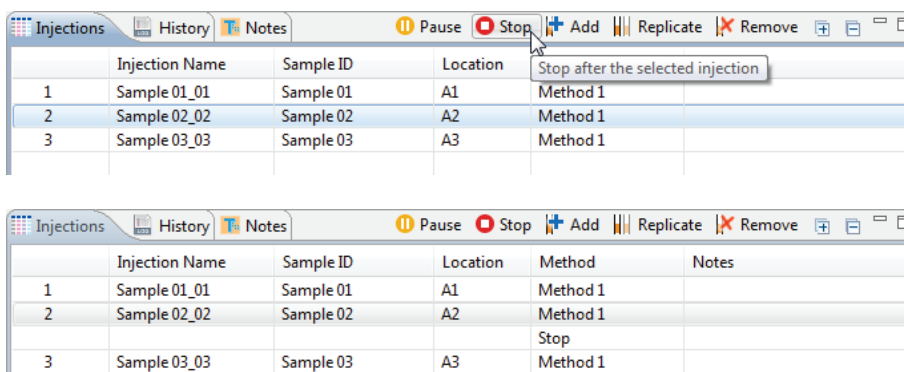
NOTE: Maurice can tweet you when the batch pauses. See “Setting Up Maurice Systems to Send Tweets” on page 764.



2. To resume the batch, click **Continue** in the instrument status bar:

**To stop the run after a specific injection:**

1. Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.

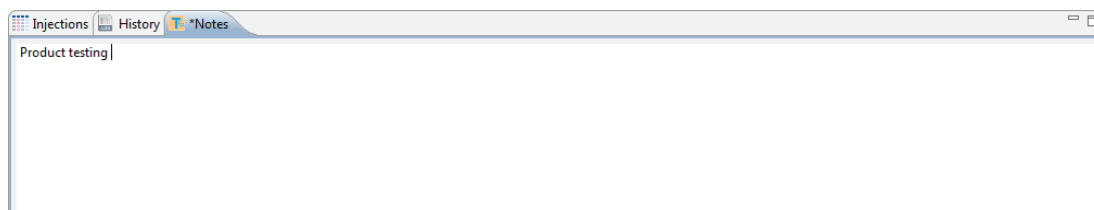


Data collected for injections completed up to the point where the batch was stopped are saved in the .mbz file.

2. Be sure to complete your normal post-batch procedures.

Step 6 - Add Batch Notes (Optional)

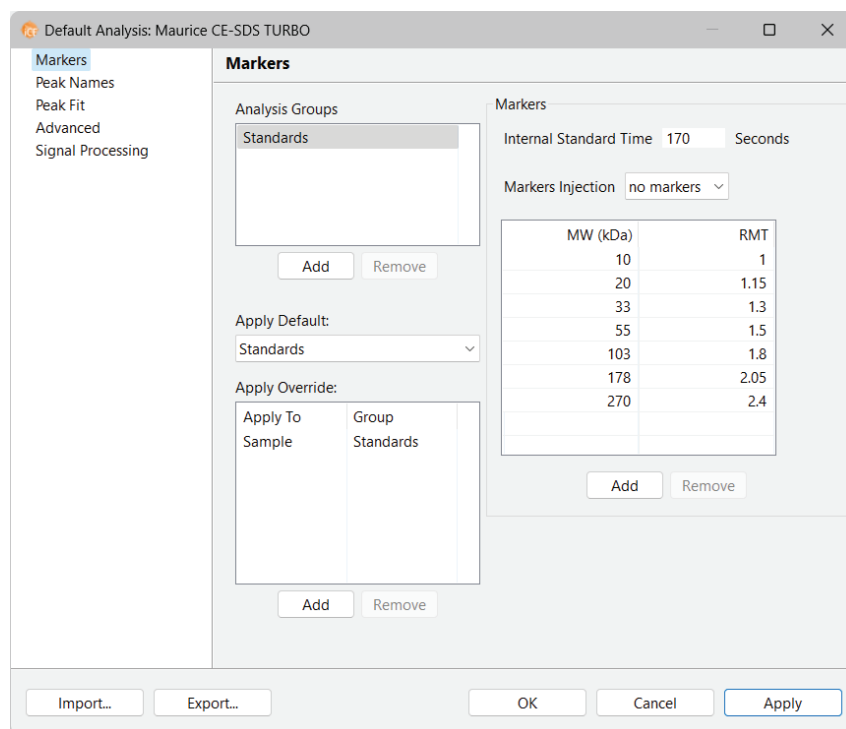
1. Click on the **Notes** pane.
2. Click in the notes area and type any information you want to add about your batch.



Step 7 - Modify Default Analysis Parameters (Optional)

You can preset the parameters used to analyze data generated with the batch. We recommend using the default parameters for Turbo CE-SDS applications, but if you need to modify parameters:

1. Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:



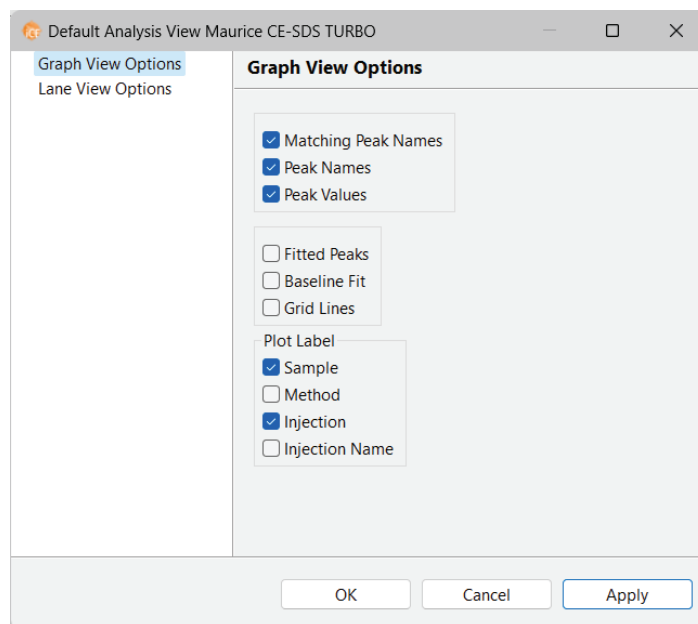
2. Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to “Analysis Settings Overview” on page 713.

Step 8 - Modify Default Analysis View Parameters (Optional)

You can preset the parameters used to display the graph and lane view data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph View Options and Lane View Options buttons.

To modify the parameters:

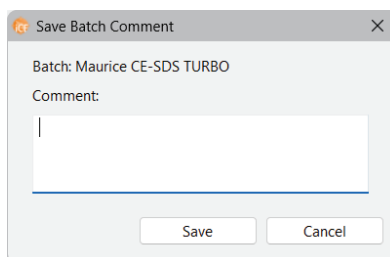
1. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



2. Change the parameters you want to, then click **Apply**. Click **OK** when you are done changing display parameters. For detailed information on data display parameters, please refer to “Customizing the Data Display” on page 695.

Step 9 - Save Your Batch

1. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



2. Enter a name for your batch then click **Save**.

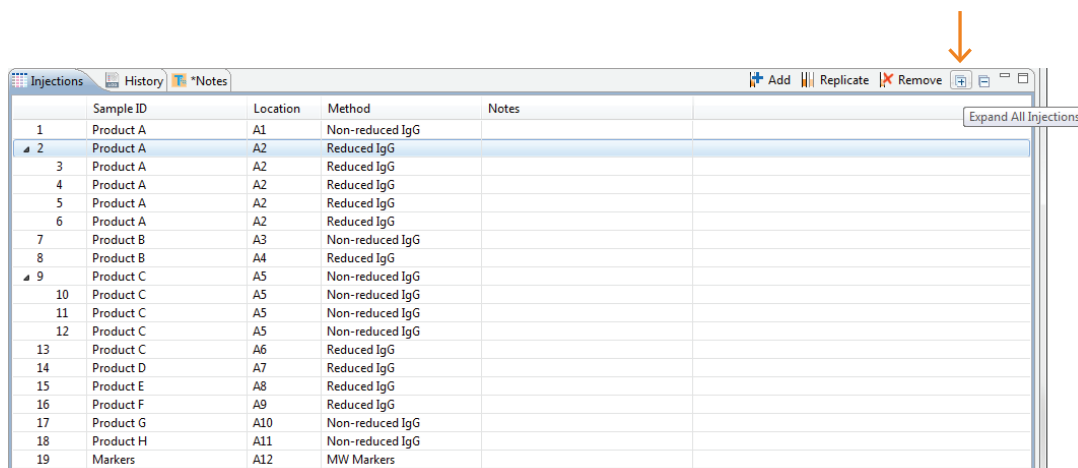
Viewing Replicate Injections

Injections with replicates have an arrow next to the injection number. You can expand or collapse the list of replicate injections by toggling the arrow:

	Sample ID	Location	Method
1	Product A	A1	Non-reduced IgG
▶ 2	Product A	A2	Reduced IgG
7	Product B	A3	Non-reduced IgG
8	Product B	A4	Reduced IgG
▶ 9	Product C	A5	Non-reduced IgG
13	Product C	A6	Reduced IgG
14	Product D	A7	Reduced IgG
15	Product E	A8	Reduced IgG
16	Product F	A9	Reduced IgG
17	Product G	A10	Non-reduced IgG
18	Product H	A11	Non-reduced IgG
19	Markers	A12	MW Markers

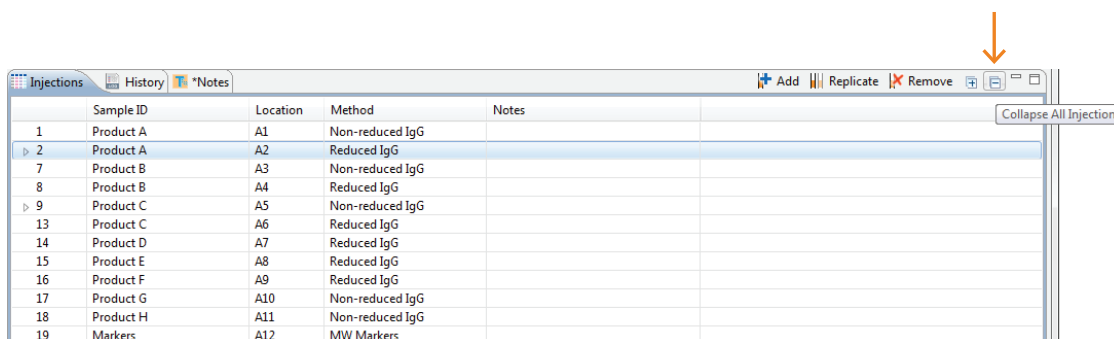
	Sample ID	Location	Method
1	Product A	A1	Non-reduced IgG
▶ 2	Product A	A2	Reduced IgG
3	Product A	A2	Reduced IgG
4	Product A	A2	Reduced IgG
5	Product A	A2	Reduced IgG
6	Product A	A2	Reduced IgG
7	Product B	A3	Non-reduced IgG
8	Product B	A4	Reduced IgG
▶ 9	Product C	A5	Non-reduced IgG
13	Product C	A6	Reduced IgG
14	Product D	A7	Reduced IgG
15	Product E	A8	Reduced IgG

- To show all replicate injections in the batch, click the **Expand All Injections** button.



	Sample ID	Location	Method	Notes
1	Product A	A1	Non-reduced IgG	
▶ 2	Product A	A2	Reduced IgG	
3	Product A	A2	Reduced IgG	
4	Product A	A2	Reduced IgG	
5	Product A	A2	Reduced IgG	
6	Product A	A2	Reduced IgG	
7	Product B	A3	Non-reduced IgG	
8	Product B	A4	Reduced IgG	
▶ 9	Product C	A5	Non-reduced IgG	
10	Product C	A5	Non-reduced IgG	
11	Product C	A5	Non-reduced IgG	
12	Product C	A5	Non-reduced IgG	
13	Product C	A6	Reduced IgG	
14	Product D	A7	Reduced IgG	
15	Product E	A8	Reduced IgG	
16	Product F	A9	Reduced IgG	
17	Product G	A10	Non-reduced IgG	
18	Product H	A11	Non-reduced IgG	
19	Markers	A12	MW Markers	

- To hide all replicate injections in the batch, click the **Collapse All Injections** button.



	Sample ID	Location	Method	Notes
1	Product A	A1	Non-reduced IgG	
▶ 2	Product A	A2	Reduced IgG	
7	Product B	A3	Non-reduced IgG	
8	Product B	A4	Reduced IgG	
▶ 9	Product C	A5	Non-reduced IgG	
13	Product C	A6	Reduced IgG	
14	Product D	A7	Reduced IgG	
15	Product E	A8	Reduced IgG	
16	Product F	A9	Reduced IgG	
17	Product G	A10	Non-reduced IgG	
18	Product H	A11	Non-reduced IgG	
19	Markers	A12	MW Markers	

Batch History

Clicking on the History pane shows the batch event history, starting with the date and time the batch was created through the most current event. Clicking on a row in the table displays the full event details in the box under the table.

Date	User Name	Message	Comment
2022-01-17 14:28:1		Batch created using the factory default Maurice CE-SDS TURBO with Compass for iCE Version: 3.0.0-0114	
2022-01-17 14:29:4		Saved as C:\Users\hxu\Documents\Compass for iCE\Batches\Maurice CE-SDS TURBO_D1-8_96x.batch fro...	
2022-01-17 14:29:4		Save injections changes to C:\Users\hxu\Documents\Compass for iCE\Batches\Maurice CE-SDS TURBO_...	

Time	User
2022-01-17 14:28:14	

Message
Batch created using the factory default Maurice CE-SDS TURBO with Compass for iCE Version: 3.0.0-0114

Comment

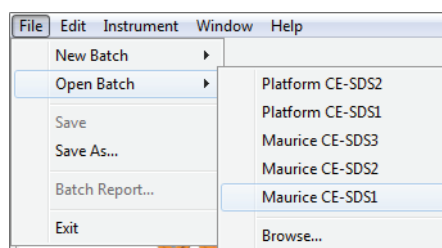
- **Date:** Date and time of the batch event.
- **User Name:** User that initiated the event. User names only display if you're using the Access Control feature to help satisfy the 21CFR Part 11 data security requirements. Go to "Enabling Access Control" on page 781 to learn how to set it up.
- **Message:** Description of the event that took place.
- **Comment:** Comments entered by the user when the batch was saved.

You can copy the information in the History pane to use in other documents:

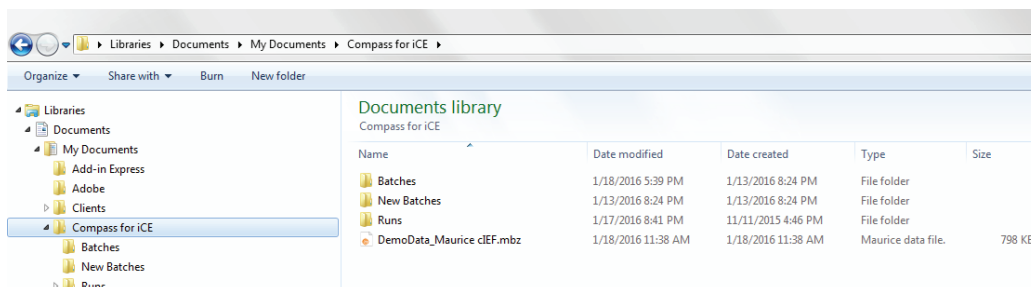
1. Click the **History** pane to make sure it's active.
2. Click **Edit** in the main menu and select **Copy**.
3. Open a document and click **Paste**.

Making Changes to a Batch

1. Select **File** in the main menu and click **Open Batch**.



2. A list of the last five batches opened will display based on the last location found. If the location of the file has changed, you may need to navigate to find the file.
 - Select one of these files or click **Browse** to open the Batch folder and select a different one.
 - To open a batch used in a run file, click **Browse**. Go to the Run folder in the Compass for iCE directory and select the run file that uses the batch you want to edit.

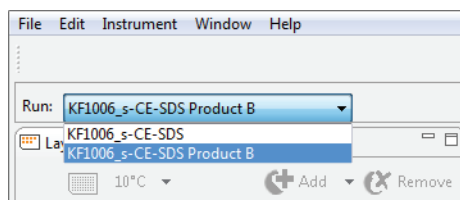


3. To make changes to the batch, see the steps in “Creating a New Batch” on page 332. Then select **File** from the main menu and click **Save** or **Save As**.

Viewing and Editing Batches in Completed Runs

1. Click the **Analysis** screen and open your run file(s).
2. After the run opens, click the **Batch** screen to view the batch used with the run.

If you opened more than one run file, you can switch between viewing batches for each file. Just click the **down arrow** in the Run box and select the batch you want to view from the drop down list.



3. In completed run files, you can only make edits to the **sample** and **method** names in a batch, and these changes are saved to the run file only. Update sample and method names as needed.
4. Click the **Analysis** screen. Then select **File** from the main menu and click **Save** or **Save As** to save the new changes to the run file.

Copying and Pasting Injection Names or Sample IDs from other Documents

Injection names and Sample IDs can be copied from other documents and pasted into the Injection pane to make batch setup faster.

To paste a list of injection names into the batch:

1. Select the list of injection names in a document (Microsoft® Word®, Excel®, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 2	My Sample 2
4		My Injection 3	My Sample 3
5		My Injection 4	My Sample 4
6		My Injection 5	My Sample 5
7		My Injection 6	My Sample 6
8		My Injection 7	My Sample 7
9		My Injection 8	My Sample 8
10		My Injection 9	My Sample 9
11		My Injection 10	My Sample 10

2. Select an injection in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the injection, right click and select **Paste**.

⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove 🔍 📄						
	Injection Name	Sample ID	Location	Method	Notes	
1	Sample 01_01	Sample 01	A1	System Suitability		
2	Sample 02_02	Sample 02	A2	Method2		
3	Sample 03_03	Sample 03	A3	Method2		
4	Sample 04_04	Sample 04	A4	Method2		
5	Sample 05_05	Sample 05	A5	Method2		
6	Sample 06_06	Sample 06	A6	Method2		
7	Sample 07_07	Sample 07	A7	Method2		
8	Sample 08_08	Sample 08	A8	Method2		
9	Sample 09_09	Sample 09	A9	Method2		
10	Sample 10_10	Sample 10	A10	Method2		
11	Sample 11_11	Sample 11	A11	Method2		
12	Sample 12_12	Sample 12	A12	Method2		

The injection names are pasted into the Injection pane:

⏸ Pause ⏹ Stop ➕ Add 🔄 Replicate ✖ Remove 🔍 📄						
	Injection Name	Sample ID	Location	Method	Notes	
1	My Injection 1	Sample 01	A1	System Suitability		
2	My Injection 2	Sample 02	A2	Method2		
3	My Injection 3	Sample 03	A3	Method2		
4	My Injection 4	Sample 04	A4	Method2		
5	My Injection 5	Sample 05	A5	Method2		
6	My Injection 6	Sample 06	A6	Method2		
7	My Injection 7	Sample 07	A7	Method2		
8	My Injection 8	Sample 08	A8	Method2		
9	My Injection 9	Sample 09	A9	Method2		
10	My Injection 10	Sample 10	A10	Method2		
11	Sample 11_11	Sample 11	A11	Method2		
12	Sample 12_12	Sample 12	A12	Method2		

To paste a list of Sample IDs into the batch:

1. Select the list of Sample IDs in a document (Microsoft Word, Excel, etc.) and select **Copy** or hold the **Ctrl+C** keys to copy.

	A	B	C
1			
2		My Injection 1	My Sample 1
3		My Injection 2	My Sample 2
4		My Injection 3	My Sample 3
5		My Injection 4	My Sample 4
6		My Injection 5	My Sample 5
7		My Injection 6	My Sample 6
8		My Injection 7	My Sample 7
9		My Injection 8	My Sample 8
10		My Injection 9	My Sample 9
11		My Injection 10	My Sample 10
12			

2. Select a Sample ID in the Injection pane, then hold the **Ctrl+V** keys to paste. You can also select the Sample ID, right click and select **Paste**.

Injections History Notes					
Pause Stop Add Replicate Remove					
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 01_01	Sample 01	A1	System Suitability	
2	Sample 02_02	Sample 02	A2	Method2	
3	Sample 03_03	Sample 03	A3	Method2	
4	Sample 04_04	Sample 04	A4	Method2	
5	Sample 05_05	Sample 05	A5	Method2	
6	Sample 06_06	Sample 06	A6	Method2	
7	Sample 07_07	Sample 07	A7	Method2	
8	Sample 08_08	Sample 08	A8	Method2	
9	Sample 09_09	Sample 09	A9	Method2	
10	Sample 10_10	Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

The Sample IDs are pasted into the Injection pane.

NOTE: If you paste in new sample IDs when the default injection names are still being used, the injection names will automatically update to match the sample ID. If you've already changed the injection names manually, updating the sample IDs will not update the injection names.

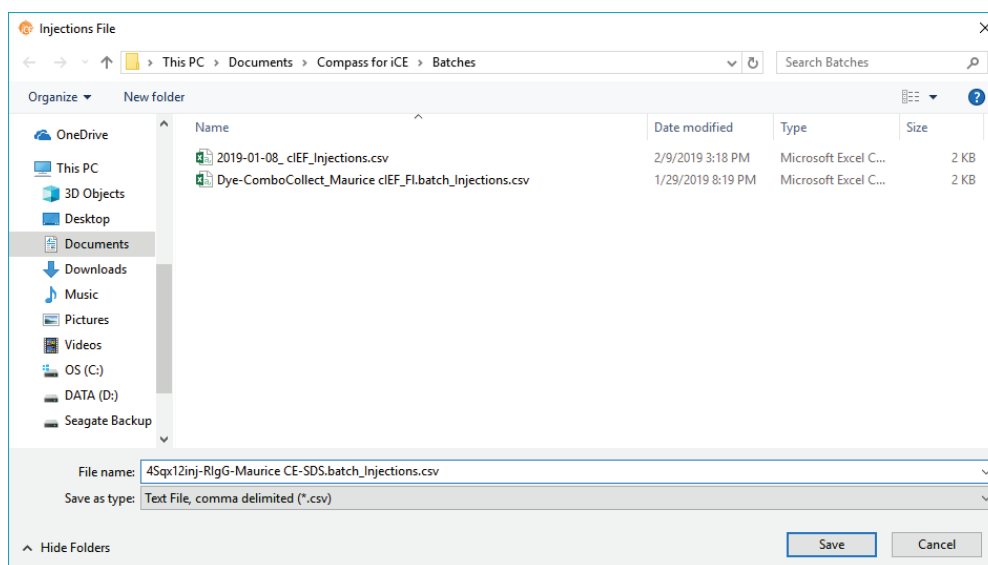
Pause Stop Add Replicate Remove					
	Injection Name	Sample ID	Location	Method	Notes
1	My Sample 1_01	My Sample 1	A1	System Suitability	
2	My Sample 2_02	My Sample 2	A2	Method2	
3	My Sample 3_03	My Sample 3	A3	Method2	
4	My Sample 4_04	My Sample 4	A4	Method2	
5	My Sample 5_05	My Sample 5	A5	Method2	
6	My Sample 6_06	My Sample 6	A6	Method2	
7	My Sample 7_07	My Sample 7	A7	Method2	
8	My Sample 8_08	My Sample 8	A8	Method2	
9	My Sample 9_09	My Sample 9	A9	Method2	
10	My Sample 10_10	My Sample 10	A10	Method2	
11	Sample 11_11	Sample 11	A11	Method2	
12	Sample 12_12	Sample 12	A12	Method2	

Importing and Exporting Injections

Injections in an open batch or run file can be exported as a separate file. This lets you import the same or a modified set of injections into another batch later, rather than having to enter the information manually.

Exporting Injections

1. Open the batch or run you want to export injections from.
2. In the Batch screen, select **File** in the main menu and click **Export Injections**. The following window displays:



3. The default directory is Compass for iCE/Batches. Change the directory if needed.
4. The default file name will be batchname.batch_Injections.csv. Change the name if needed and click **Save**.

Creating an Injections File

To create a new injections file .csv, it's easiest to export injections from an existing batch and use that as your template.

1. Follow the instructions in Exporting Injections above to export injections from an existing Turbo CE-SDS batch.
2. Open the .csv file in a program that provides a table/spreadsheet format.
3. To add or change injection information:
 - a. Type in a unique name in each cell in the Injection Name column. For replicate injections, type > in front of the injection name, for example: >Injection 1.
 - b. Type in a sample ID, location and method for each injection in their respective columns.

- c. Optional: Type in notes if needed.

NOTES:

Mix bottle is not used for Turbo CE-SDS batches. Compass for iCE ignores this column when importing injections into Turbo CE-SDS batches.

Programmed pauses and stops can't be added in the injections file. They can be added in the Injections pane after importing your injections.

	A	B	C	D	E	F
1	Injection Name	Sample ID	Location	Method	Mix Bottle	Notes
2	R IgG1	SB	B2	Reduced IgG		
3	R IgG 2	SB	B3	Reduced IgG		
4	>R IgG 2	SB	B3	Reduced IgG		
5	>R IgG 2	SB	B3	Reduced IgG		
6	R IgG 3	SB	C2	Reduced IgG		
7	R IgG 4	SB	C3	Reduced IgG		
8	R IgG1_05	SB	B2	Reduced IgG		
9	R IgG 2_06	SB	B3	Reduced IgG		
10	R IgG 3_07	SB	C2	Reduced IgG		

4. Save the .csv file.

Importing Injections

1. Open the batch you want to import injections into, or open a new batch.
2. Select **File** in the main menu and click **Import Injections**.
3. Select an injections file (*.csv) and click **OK**. The imported information will display in the Layout and Injections panes.

NOTE: Imported injections will replace any injections already in the Injections table, they will not append an existing table.

Importing and Exporting Methods

Methods in an open batch or run file can be exported as a separate file. This lets you import the same method into another batch later, rather than having to re-enter method information manually.

Importing a Method

NOTE: Importing a method imports information into the Batch window's Method pane only.

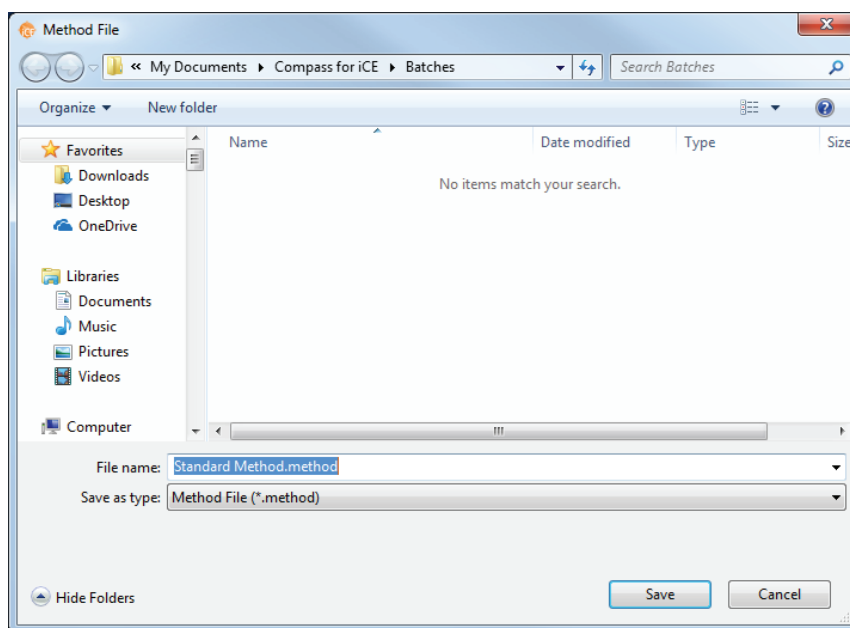
1. Open the batch you want to import the method into.

2. Select **File** in the main menu and click **Import Method**.
3. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

Exporting a Method

NOTE: Exporting a method exports information in the Batch window's Method pane only.

1. Open the batch you want to export the method from.
2. In the Method pane, select the method you want to export. To export more than one method at the same time, hold the **Ctrl** key to select multiple methods.
3. Select **File** in the main menu and click **Export Method**. The following window displays:



4. The default directory is Compass for iCE/Batches. Change the directory if needed.
5. Enter a method file name and click **Save**. The settings will be saved as a *.method file.

Batch Reports

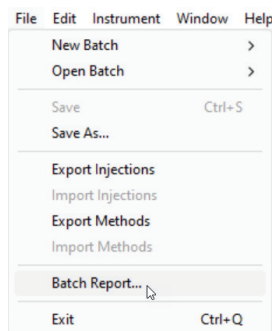
You can export a PDF file of sample and method details for each injection in the batch.

NOTES:

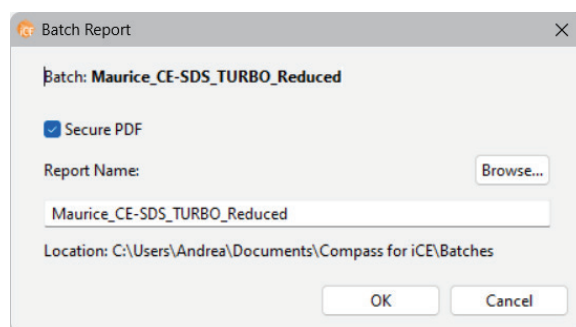
You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it’s saved in a separate run folder. Existing reports aren’t overwritten.

1. Go to the **Analysis** or **Run Summary** screen, then click **File > Open Run** and select a run file (if you don’t have one open already).
2. After the run opens, go to the **Batch** screen.
3. Select **File** from the main menu and click **Batch Report**.

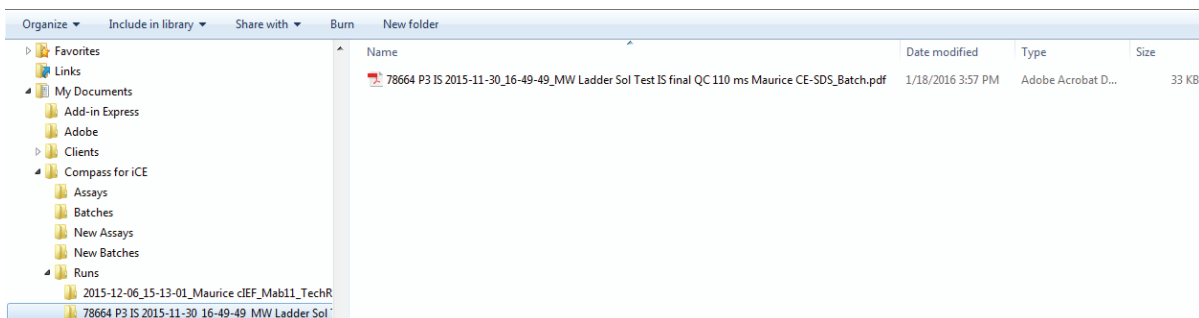


4. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.



Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don’t want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.

5. The Batch Report PDF is exported to the Runs folder in the Compass for iCE directory. It'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Here's an example Batch Report:

CE-SDS Batch: Maurice_CE-SDS_TURBO_Reduced

Injection	Injection Name	Sample ID	Location	Method	Sample Load	Separation
1	Sample 01_01	Sample 01	A1	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
13	Sample 02_13	Sample 02	A2	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
25	Sample 03_25	Sample 03	A3	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
37	Sample 04_37	Sample 04	A4	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
49	Sample 05_49	Sample 05	A5	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
61	Sample 06_61	Sample 06	A6	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
73	Sample 07_73	Sample 07	A7	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts
85	Sample 08_85	Sample 08	A8	Reduced IgG	8 sec 3500 Volts	5.5 min, 4200 Volts

Created By: Andrea Sat 11:56 PM Feb 25, 2023 PST (SECURED)
 C:\Users\Andrea\Documents\Compass for iCE\Batches\Maurice_CE-SDS_TURBO_Reduced\batch
 Computer: DESKTOP-1FM7G05 Software Version: Compass for iCE 4.0.0, Build ID: 0222

Page 1 of 2



CE-SDS Batch: Maurice_CE-SDS_TURBO_Reduced

Batch Log

Date	User Name	Message	Comment
2022-12-14 10:22:35		Saved as C:\Users\Andrea\Documents\Compass for iCE\Batches\Maurice_CE-SDS_TURBO_Reduced.batch from Compass for iCE v4.0.0-1102	
2022-12-14 10:22:35		Save Batch changes to C:\Users\Andrea\Documents\Compass for iCE\Batches\Maurice_CE-SDS_TURBO_Reduced.batch from Compass for iCE v4.0.0-1102	

Created By: Andrea Sat 11:59 PM Feb 25, 2023 PST (SECURED)
C:\Users\Andrea\Documents\Compass for iCE\Batches\Maurice_CE-SDS_TURBO_Reduced.batch
Computer: DESKTOP-1FM7005 Software Version: Compass for iCE 4.0.0, Build ID: 0222

Page 2 of 2



Chapter 14:

Running Turbo CE-SDS Applications

Chapter Overview

- Overview
- Before You Throw the Switch
- Power Up
- Running Turbo CE-SDS Applications
- Editing a Running Batch
- Post-batch Procedures
- Checking Your Data

Overview

Turbo CE-SDS applications can be run on Maurice, Maurice S. or MauriceFlex systems with a Turbo CE-SDS cartridge.

Before You Throw the Switch

Ensure that everyone using Maurice have:

- Received instruction in general safety practices for laboratories.
- Received instruction in specific safety practices for Maurice.
- Received instruction on handling of biohazards (if biohazardous materials are to be used on Maurice).
- Read and understood all Product Inserts and related Safety Data Sheets (SDS).

Power Up

1. Turn on the computer connected to Maurice.
2. Turn on Maurice's main power switch.
3. Wait for Maurice to initialize. His indicator light will turn from magenta to blue when he's done.
4. Double-click the Compass for iCE software icon on the computer desktop to open the application.
5. Connect Maurice to Compass for iCE.

Running Turbo CE-SDS Applications

What You'll Need

- Maurice Turbo CE-SDS Application Kit which include:
 - Maurice Turbo CE-SDS Cartridges
 - Separation Matrix
 - Turbo CE-SDS Running Buffer - Bottom
 - CE-SDS or CE-SDS PLUS 1X Sample Buffer
 - Wash Solution
 - Conditioning Solutions (1 and 2)
 - 25X Internal Standard
 - Glass reagent vials, 2 mL

- 96-well plates
- Clear screw caps for vials
- Orange pressure caps for vials
- Maurice CE-SDS IgG Standard (optional)
- Maurice CE-SDS MW Markers (optional)
- β -mercaptoethanol (β ME, >98% = 14.2 M) for reducing conditions
- Optional: Sodium hypochlorite solution (10-15%), for neutralizing β -mercaptoethanol (Sigma PN 425044)
- Iodoacetamide (IAM, 250 mM) for alkylation at non-reducing conditions
- Deionized (DI) water
- Sample vials with integrated inserts for samples (optional)
- P10, P20, P200, P1000 and pipette tips
- Water bath or thermocycler
- Vortexer
- Table-top centrifuge with plate adapter or vial adapter (12 mm, 2 mL vials)

Step 1: Prep Your Internal Standard, Samples and Reagents

NOTE: You can prepare your samples to run either in 96-well plates or vials. Using 96-well plates is the default method.

Internal Standard

NOTES:

The Internal Standard is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Aliquot the reconstituted solution into appropriately sized vials and store at -70 to -90 °C for long term storage. For short-term storage (< 1 week), the solution can be stored at 2-8 °C.

Prepare the Internal Standard in the same buffer as your sample.

1. Open the vial of lyophilized 25X Internal Standard by lifting the center tab and gently pulling it back to break the metal seal. Then slowly remove the rubber stopper.
2. Reconstitute by adding 240 μ L of 1X Sample Buffer. Pipette up and down a few times to mix thoroughly. This results in a 25X Internal Standard solution.

NOTE: Don't vortex the reconstituted Internal Standard during prep.

Sample Prep Under Reducing Conditions

NOTES:

Prepare a minimum of 100 μ L of sample for a Turbo CE-SDS batch.

Reduced IgG Sample

NOTE: If you are sample limited and running a Turbo CE-SDS batch, prepare your sample in a final volume of 50 μ L and dilute the denatured sample 1:1 in DI water to bring the volume to 100 μ L.

1. In a microcentrifuge tube, dilute your IgG sample with 1X Sample Buffer to a concentration of 0.25-1.00 mg/mL in a final volume of 100 μ L.
-

NOTE: Dilute at least 1:1 with either CE-SDS 1X Sample Buffer or CE-SDS PLUS 1X Sample Buffer.

2. Add 2 μ L of reconstituted 25X Internal Standard for every 50 μ L of sample volume.
 3. Add 2.5 μ L of 14.2 M β -mercaptoethanol for every 50 μ L of sample volume.
 4. Mix thoroughly.
-

NOTE: Heating in the presence of SDS linearizes the protein and allows for SDS binding. The reducing agents break up inter- and intra-molecular disulfide bonds.

5. Centrifuge the tube and heat the mixture in a water bath or thermocycler at 70 °C for 10 minutes.
 6. Put the tube on ice for 5 minutes.
 7. Vortex briefly and spin down.
-

Reduced IgG Standard (Optional)

NOTES:

The IgG Standard is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Prepare the reduced IgG Standard in either CE-SDS 1X Sample Buffer or CE-SDS PLUS 1X Sample Buffer.

1. Using scissors, carefully cut the top of the foil package, leaving the sealing strip intact.
2. Take out the strip of tubes and carefully cut one pink tube of lyophilized IgG Standard from the strip. Put the unopened tubes back in the package, seal tightly and store at 2-8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Reconstitute the IgG Standard with 50 µL of 1X Sample Buffer. Gently resuspend by pipetting the solution up and down. Transfer the solution to a microcentrifuge tube.
5. Add 2 µL of reconstituted 25X Internal Standard.
6. Add 2.5 µL of 14.2 M β-mercaptoethanol.
7. Mix thoroughly by vortex.
8. Centrifuge the tube and heat mixture in a water bath or thermocycler at 70 °C for 10 minutes.
9. Put the tube on ice for 5 minutes.
10. Vortex briefly and spin down.

CE-SDS Molecular Weight (MW) Markers (Optional)

NOTES:

The CE-SDS MW Markers are lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Prepare the MW Markers in either CE-SDS 1X Sample Buffer or CE-SDS PLUS 1X Sample Buffer.

1. Using scissors, carefully cut the top of the foil package leaving the sealing strip intact.
2. Take out the strip of tubes and carefully cut one green tube of lyophilized CE-SDS MW Markers from the strip. Put the unopened tubes back in the package, seal tightly and store at 2-8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Reconstitute the CE-SDS MW Markers with 50 µL of 1X Sample Buffer. Gently resuspend by pipetting the solution up and down. Transfer the solution to a microcentrifuge tube.
5. Add 2 µL of reconstituted 25X Internal Standard.
6. Add 2.5 µL of 14.2 M β-mercaptoethanol.
7. Mix thoroughly.
8. Centrifuge the tube and heat the mixture in a water bath or thermocycler at 70 °C for 10 minutes.
9. Put the tube on ice for 5 minutes.
10. Vortex briefly and spin down.

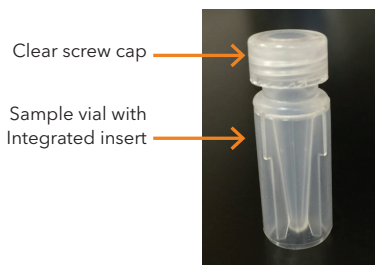
Spin Samples, Standards and CE-SDS MW Markers

If you're using a 96-well plate:

1. Transfer 100 μ L of each of your samples, IgG Standard and CE-SDS MW Markers to their designated wells in a 96-well plate.
2. Cover the plate with a lid and spin for 10 minutes at 1000 xg in a centrifuge with a plate adapter.
3. Pop any bubbles in the samples with a clean pipette tip.

If you're using vials:

1. Transfer 100 μ L of each of your samples, IgG Standard and CE-SDS MW Markers to their designated sample vials with insert, 0.2 mL.
2. Close the vials with a clear screw cap.
3. Place the vials in a centrifuge using a vial adapter (12 mm, 2 mL vials) and spin for 10 minutes at 1000 xg.



Sample Prep Under Non-reducing Conditions

Alkylation Reagent

NOTES:

We use a 250 mM solution of iodoacetamide (IAM) as an alkylating reagent.

Prepare a fresh 250 mM solution of iodoacetamide in DI water before use.

1. Weigh out 46 mg of IAM directly into a 1.5 mL microcentrifuge tube.
2. Add 1 mL of DI water to the tube and mix thoroughly.

Non-reduced IgG Sample**NOTES:**

Prepare a minimum of 100 μ L of sample for a Turbo CE-SDS batch.

If you are sample limited and running a Turbo CE-SDS batch, prepare your sample in a final volume of 50 μ L and dilute the denatured sample 1:1 in DI water to bring the volume to 100 μ L.

1. In a microcentrifuge tube, dilute your IgG sample with 1X Sample Buffer to a concentration of 0.25-1.0 mg/mL in a final volume of 50 μ L.

NOTE: Dilute at least 1:1 with CE-SDS 1X Sample Buffer or CE-SDS PLUS 1X Sample Buffer.

2. Add 2 μ L of reconstituted 25X Internal Standard for every 50 μ L of sample volume.
3. Add 2.5 μ L of 250 mM IAM for every 50 μ L of sample volume.
4. Mix thoroughly.
5. Centrifuge the tube and heat the mixture in a water bath or thermocycler at 70 °C for 10 minutes.
6. Put the tube on ice for 5 minutes.
7. Vortex briefly and spin down.

Non-reduced IgG Standard (Optional)**NOTES:**

The IgG Standard is lyophilized. Always reseal the foil bag containing the unopened tubes with desiccant to prevent moisture absorption.

Prepare the non-reduced IgG Standard in either CE-SDS 1X Sample Buffer or CE-SDS PLUS 1X Sample Buffer.

1. Using scissors, carefully cut the top of the foil package, leaving the sealing strip intact.
2. Take out the strip of tubes and carefully cut one pink tube of lyophilized IgG Standard from the strip. Put the unopened tubes back in the package, seal tightly and store at 2-8 °C.
3. Pierce the foil on the tube with a clean pipette tip.
4. Reconstitute the IgG Standard with 50 μ L of 1X Sample Buffer. Gently resuspend by pipetting the solution up and down. Transfer the solution to a microcentrifuge tube.
5. Add 2 μ L of reconstituted 25X Internal Standard.
6. Add 2.5 μ L of 250 mM IAM.

7. Mix thoroughly by vortex.

NOTE: Heating in the presence of SDS linearizes the protein and allows for SDS binding. The alkylating agents prevent disulfide-bond scrambling catalyzed by free sulfhydryl groups. This minimizes the appearance of fragments under non-reducing conditions.

8. Centrifuge the tube and heat mixture in a water bath or thermocycler at 70 °C for 10 minutes.
9. Put the tube on ice for 5 minutes.
10. Vortex briefly and spin down.

Spin Samples and Standards

If you're using a 96-well plate:

1. Transfer 100 µL of each of your samples and IgG Standard to their designated wells in a 96-well plate.
2. Cover the plate with a lid and spin for 10 minutes at 1000 xg in a centrifuge with a plate adapter.
3. Pop any bubbles in the samples with a clean pipette tip.

If you're using vials:

1. Transfer 100 µL of your samples and IgG Standard to their designated sample vials with insert, 0.2 mL.
2. Close the vials with a clear screw cap.
3. Place the vials in a centrifuge using vial adapter (12 mm, 2 mL vials) and spin for 10 minutes at 1000 xg.

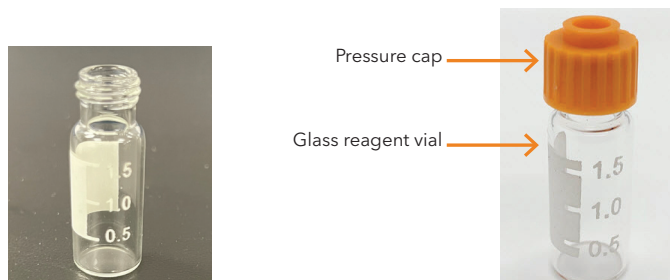
Reagents

IMPORTANT:

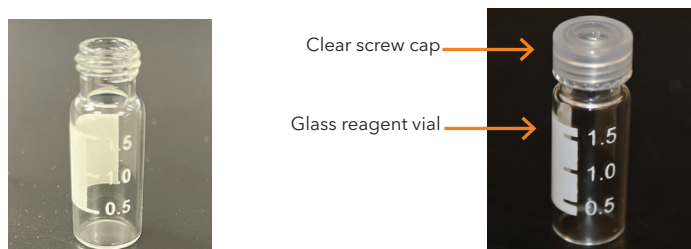
Use glass reagent vials, 2 mL (PN 046-017) to prepare Turbo CE-SDS batch reagents.

NOTE: Don't reuse reagents, vials or the clear screw caps. Always keep the orange pressure caps paired with their respective reagents. If you want to reuse pressure caps, first wash them thoroughly with DI water, soak overnight in DI water, rinse caps thoroughly with DI water again and air dry them before reusing.

1. Pipette 1.5 mL of Conditioning Solution 1 into a glass reagent vial, label and close with an **orange pressure cap**.



2. Pipette 1.5 mL of Conditioning Solution 2 into a glass reagent vial, label and close with an **orange pressure cap**.
3. Pipette 1.0 mL of Wash Solution into a glass reagent vial, label each and close with an **orange pressure cap**.
4. Pipette 1.5 mL of Wash Solution into a glass reagent vial, label each and close with a **clear screw caps**.



5. Pipette 1 mL of Separation Matrix into a glass reagent vial, label and close with an **orange pressure cap**.
6. Pipette 1 mL of Running Buffer - Bottom into three glass reagent vials, label each and close with a **clear screw cap**.
7. Pipette 1.5 mL of DI water into a glass reagent vials, label and close with an **orange pressure cap**.
8. Close an empty glass reagent vial with an **orange pressure cap**.

Step 2: Prep the Cartridge

NOTES:

A Turbo CE-SDS cartridge is guaranteed for 100 injections, with a maximum of 96 injections per batch and a maximum of 25 batches.

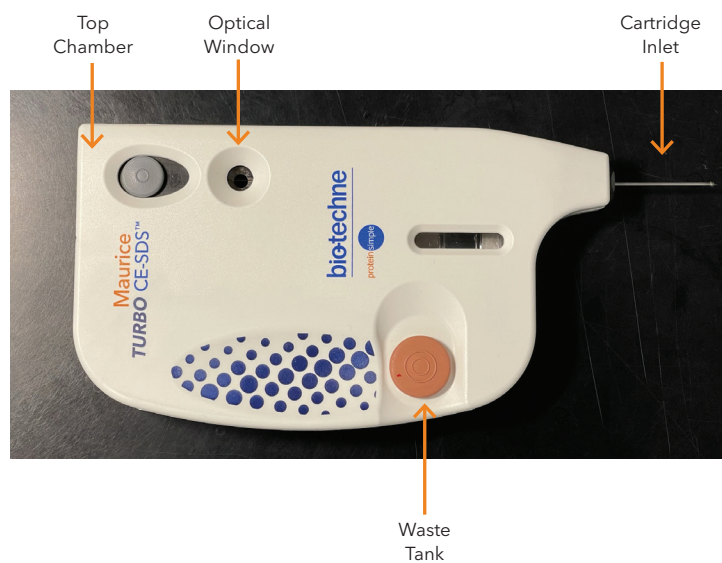
Prepare your samples and reagents before preparing the cartridge. Allowing the Separation Matrix to sit in the cartridge too long may result in cartridge clogs.

1. Take the Turbo CE-SDS Cartridge out of its packaging. Save the packaging, you'll need it later to store the cartridge.

NOTE: When you remove the cartridge from its packaging, make sure the cartridge inlet doesn't come in contact with any surface.



2. Place the cartridge on a flat surface and remove the stopper from the Top Chamber.



3. Add 750 μ L of Separation Matrix to the Top Chamber.



4. Pop any visible bubbles with a clean pipet tip.
5. Firmly close the chamber with the stopper.
6. Ensure the Waste Tank is firmly closed with the stopper.

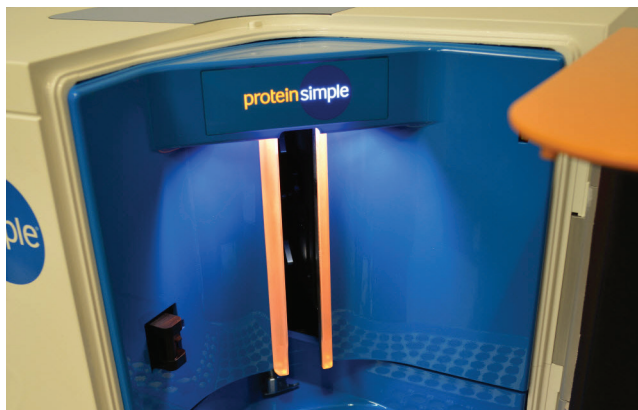
Step 3: Install the Cartridge

1. Open Maurice's door by touching the metal plate on top of the door.



NOTE: The indicator light on Maurice's front panel will blink rapidly as the door disengages.

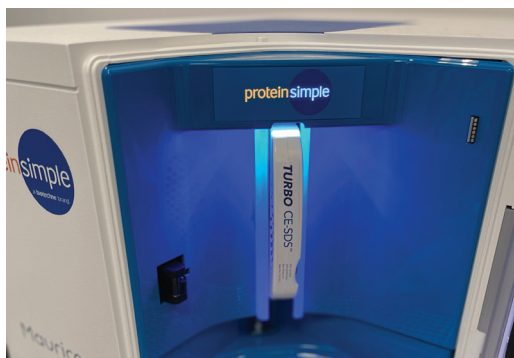
2. Swing the door open to access the cartridge slot and the reagents and samples platform. The lights on either side of the slot will be **orange**.



3. Lift the cartridge and hold it vertically using the finger holds on either side, cartridge inlet down, with the Turbo CE-SDS label facing you.
4. Gently insert it into the slot. You'll feel the alignment groove on the cartridge align inside the slot when you insert it.



5. Continue to slide the cartridge into the slot until the locking mechanism engages. The lights on either side of the slot will change to **blue** once the cartridge is installed correctly.



Step 4: Load Samples and Reagents

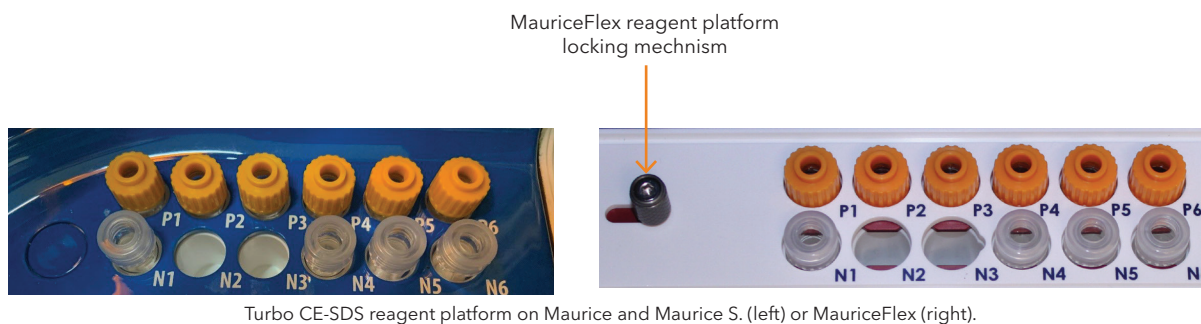
1. Place the reagent vials into their respective positions in the sample and reagents platform:

NOTES:

The two reagent rows in Maurice's reagent and sample platform are kept at room temperature.

Pressure caps are orange and have a raised, ringed surface. They should only be used in Reagent Row P. Use reagent vials with clear screw caps in Row N.

- P1 - Conditioning Solution 1 with **orange pressure cap**
- P2 - Conditioning Solution 2 with **orange pressure cap**
- P3 - DI water with **orange pressure cap**
- P4 - Separation Matrix with **orange pressure cap**
- P5 - Wash Solution with **orange pressure cap**
- P6 - Empty vial (air) with **orange pressure cap**
- N1 - Wash Solution with **clear screw cap**
- N4 - Turbo Running Buffer - Bottom with **clear screw cap**
- N5 - Turbo Running Buffer - Bottom with **clear screw cap**
- N6 - Turbo Running Buffer - Bottom with **clear screw cap**



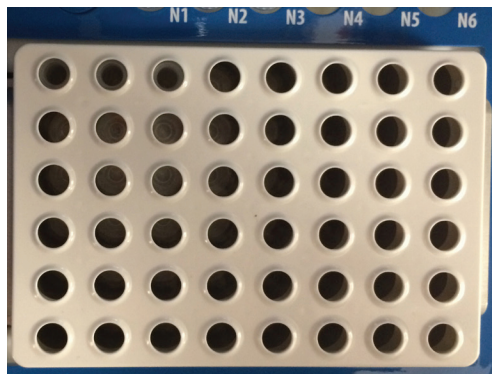
IMPORTANT:

Reagent vials on MauriceFlex must be locked in place by sliding the locking mechanism from the left to the right before starting a batch. A warning will appear in the Start Run window if they are not locked when you start the batch.

2. Depending on what you prepared your samples in, put your 96-well sample plate in the metal plate insert or your sample vials in the metal vial insert. If you have a lid on the 96-well sample plate, be sure to remove it before closing the instrument door!

NOTE: Well A1 on the 96-well plate should be in the top left corner of the insert.

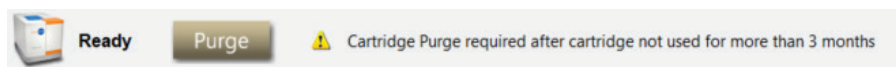
3. If you are using a vial tray, place the condensation lid on top of the vials.



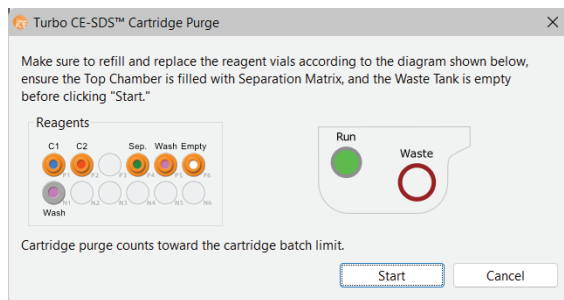
4. Close the instrument door. Maurice locks it automatically.

Step 5: Check for Cartridge Alerts

1. If your cartridge was last used in a run with an error or if it has not been used in the last three months, you will need to perform a Cartridge Purge. The Maurice system will remind you if a Cartridge Purge is required.
 - a. To purge the cartridge, click the brown **Purge** button in the instrument status bar.



- b. Confirm that the required batch reagents are loaded and that the the cartridge is prepped. Then click **Start**.



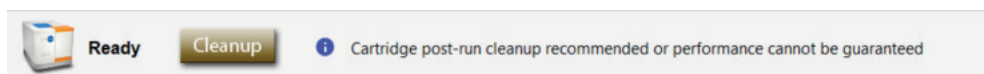
- c. Once the purge is done, perform a Cartridge Post-Run Cleanup by clicking on the brown **Cleanup** button that will appear in the instrument status bar. See step 2 for more info.

NOTE: If there's only one batch left on the cartridge, you'll get a message that it can't be used after purging.

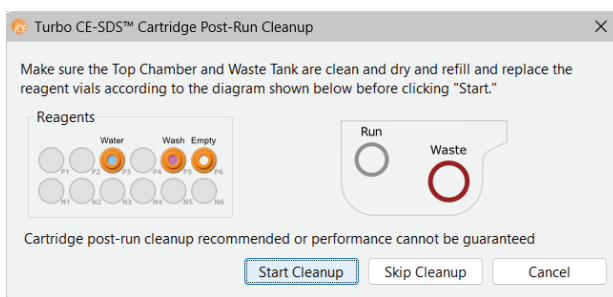
2. If you just completed a Cartridge Purge or if the cartridge you're using did not undergo a cartridge post-run cleanup after the last run, you will be asked if you want to perform a Cartridge Post-Run Cleanup.

To start the Post-Run Cleanup:

- a. Remove the cartridge from Maurice and remove the Separation Matrix from the Top Chamber and any liquid from the Waste tank from the cartridge. See "Post-batch Procedures" on page 385 for more info.
- b. Confirm there is a vial of Water (P3), Wash Solution (P5) and Air (P6) in place,
- c. Click the brown **Cleanup** button in the instrument status bar.



- d. Re-install the cartridge in Maurice.
- e. Click **Start Cleanup** in the prompt that appears.

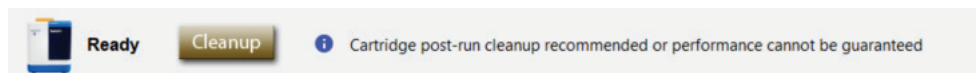


NOTE: You can also start a Post-Run Cleanup by selecting **Instrument > Cartridge Post-Run Cleanup**. When you start a Cartridge Post-Run Cleanup from the instrument menu, click **Start** in the prompt that appears to begin the cleanup.

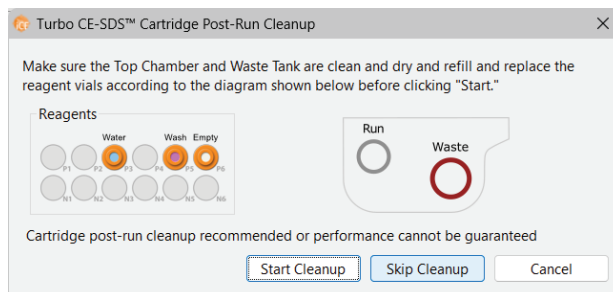
- f. Once the cartridge post-run cleanup is completed, re-prepare the cartridge and install the cartridge in Maurice. See "Step 2: Prep the Cartridge" on page 364 for more information.
- g. Click on the green **Start** button in the instrument status menu to start the batch.

To start the run without a Post-Run Cleanup:

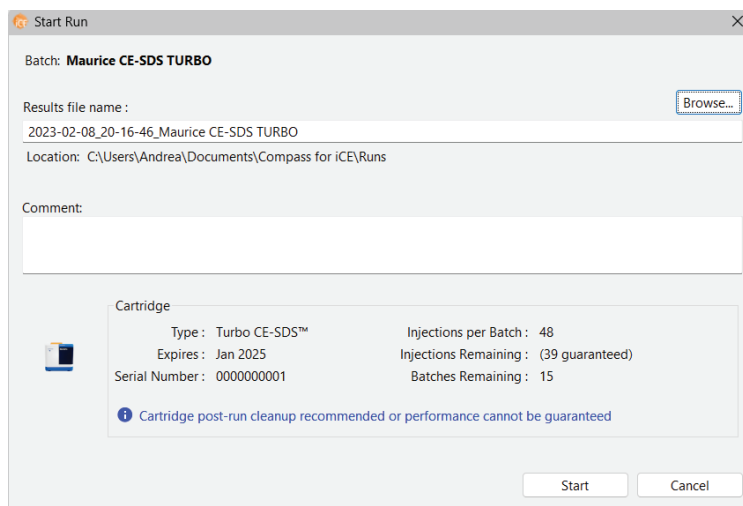
- a. Click the brown **Cleanup** button in the instrument status bar



- b. Click **Skip Cleanup** in the prompt that appears.



- c. Click **Start** in the Start Run window. The run summary and injection report will record that the cartridge did not undergo a post-run cleanup before starting the batch.

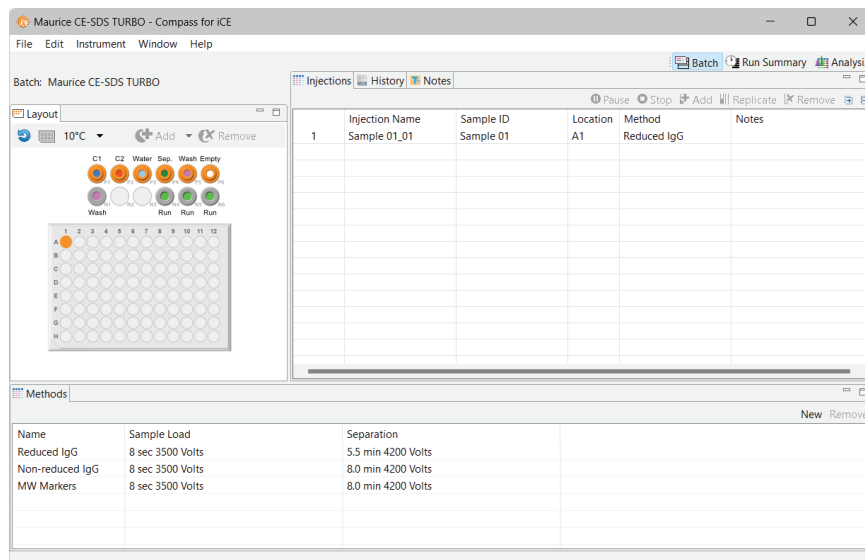
**To start the run with a different cartridge:**

- a. If necessary, click **Cancel** in the Turbo CE-SDS Cartridge Post-Run Cleanup window.
- b. Open Maurice's door, remove the first cartridge from Maurice and prepare a second cartridge. See "Step 2: Prep the Cartridge" on page 364 for more information.

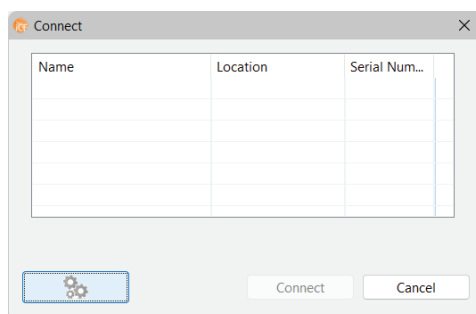
Step 6: Create a Batch

1. Launch Compass for iCE.

2. Select the **Batch** tab. This is where you'll enter sample/injection information, methods and batch parameters.

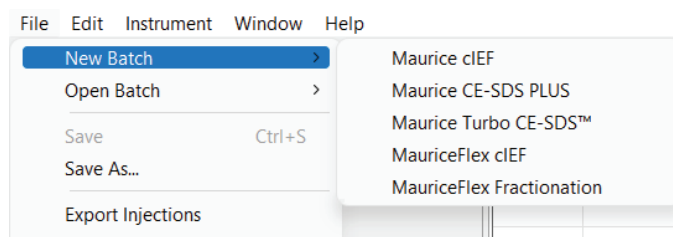


3. To create a batch, make sure your Maurice system is connected to Compass for iCE. Select **Instrument** and click **Connect**.
 - a. If your instrument is listed, select your Maurice system and click **Connect**.
 - b. If your instrument isn't listed, click on the **Settings** button and connect by typing in your instrument IP address.



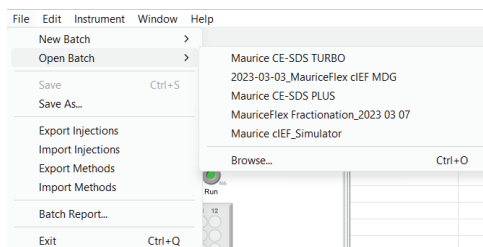
To create a new batch:

- In the main menu, select **File > New Batch > Maurice Turbo CE-SDS**.

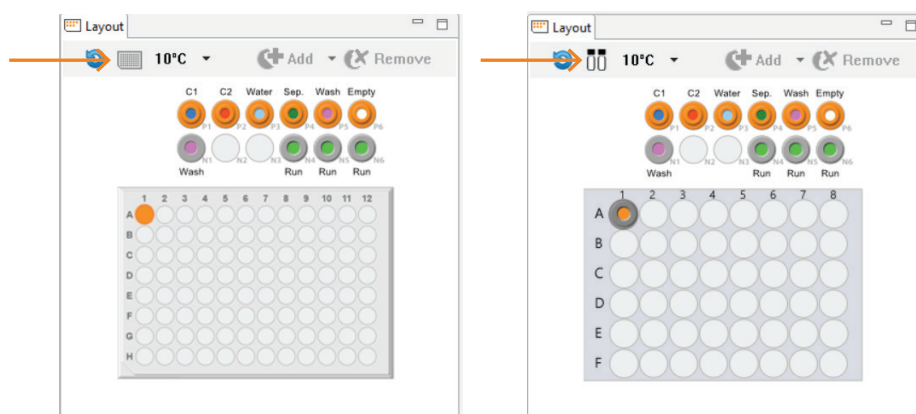


To use an existing batch: In the main menu, select **File > Open Batch**.

NOTE: If you're making changes to an existing batch, follow the steps in this section. Otherwise skip to "Step 7: Start the Batch" on page 381.



4. In the Layout pane, clicking on the vial/plate icon toggles between formats. Select a 96-well plate or 48 vials depending on what you're running.



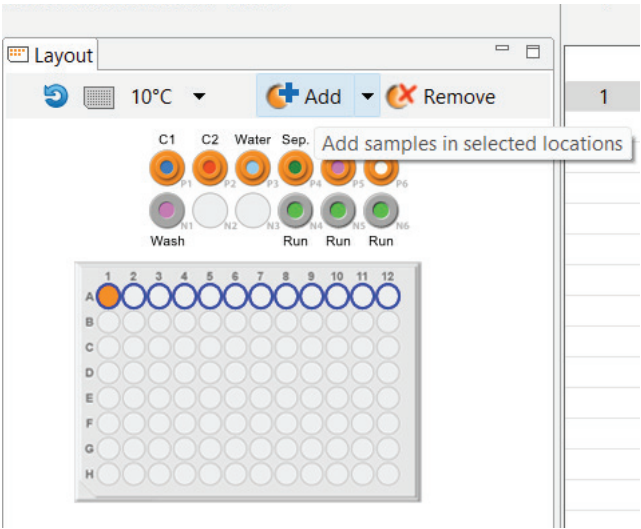
5. Add your samples:

To import samples using a saved injections file:

- Select **File** in the main menu and click **Import Injections**.
- Select an injections .csv file and click **OK**. The imported information will display in the Injections pane.

NOTE: For more information on setting up an injections file for import, see "Importing and Exporting Injections" on page 350.

To select samples manually: Use your mouse to highlight the well positions for each of your samples, then click **Add**.



This populates the Injections table:

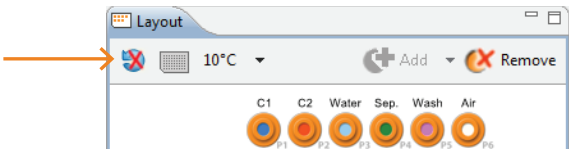
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Sample 1	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

6. Compass for iCE can monitor the current during a separation for you, stop it if an abnormal current profile is detected and reinject the sample for you automatically. Reinjects are enabled by default, but if you don't want to use this feature just click the reinject icon to toggle it off.

NOTES:

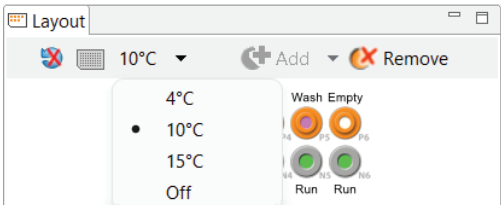
If a sample was reinjected, the injection row will be flagged in the Injections pane in the Run Summary screen. See “Injection Flags” on page 414 for more info.

A maximum of 10 reinjections are allowed per batch.



7. The sample platform is temperature controlled with a default setting of 10 °C. To change the temperature or turn it off, click the **down arrow** next to the temperature and select an option.

NOTE: The two reagent rows are kept at room temperature.



8. Enter your method parameters in the pane.

NOTE: There are three default methods. We recommend using the default method parameters for the listed samples. Please see the Maurice CE-SDS Application Guide for more information on method optimization.

To import a saved method:

- a. Select **File** in the main menu and click **Import Method**.
- b. Select a method file (*.method) and click **OK**. The imported information will display in the Method pane.

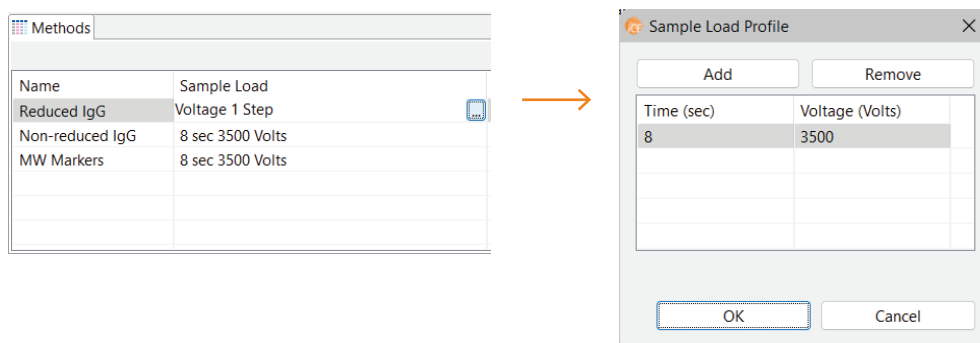
To create a new method or make changes to an existing one:

- a. Click the first cell in the Name column and enter a new method name if needed.

A screenshot of a software window titled "Methods". It contains a table with four columns. The first column is "Name", the second is "Sample Load", the third is "Separation", and the fourth is empty. The table has three rows of data. The first row is "Reduced IgG", "8 sec 3500 Volts", "5.5 min 4200 Volts". The second row is "Non-reduced IgG", "8 sec 3500 Volts", "8.0 min 4200 Volts". The third row is "MW Markers", "8 sec 3500 Volts", "8.0 min 4200 Volts". In the top right corner of the window, there are buttons labeled "New" and "Remove".

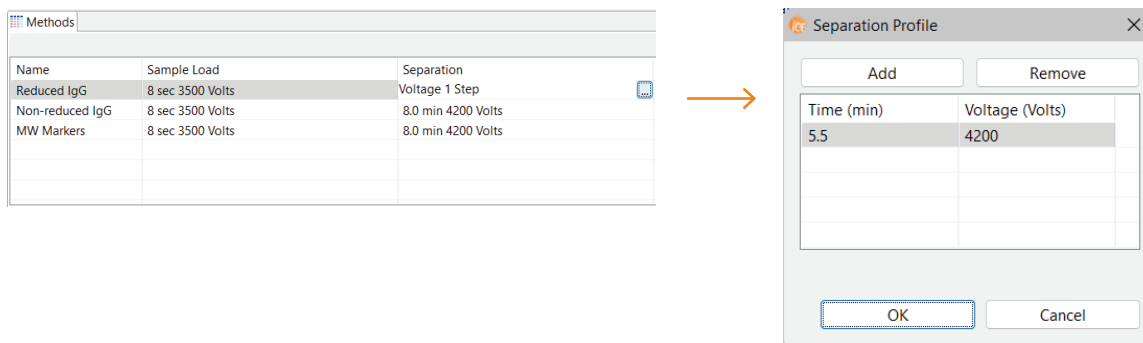
Name	Sample Load	Separation	
Reduced IgG	8 sec 3500 Volts	5.5 min 4200 Volts	
Non-reduced IgG	8 sec 3500 Volts	8.0 min 4200 Volts	
MW Markers	8 sec 3500 Volts	8.0 min 4200 Volts	

- b. Click the first cell in the Sample Load column, then click then click the selection button [...] to set your load profile time (in seconds) and voltage.



- c. Click the first cell in the Separation column then click the selection button [...] to set your separation time (in minutes) and voltage.

NOTE: Run your reduced IgG samples for 5.5 minutes and your non-reduced IgG samples, IgG Standard and CE-SDS MW Markers for 8.0 minutes. The default separation voltage for all sample types is 4200 volts.



9. You can now:

- Make updates to the remaining methods by repeating the prior steps.
- Select a method and click **Remove** in the upper right corner of the Methods pane to delete it.
- Click **New** in the upper right corner of the Methods pane to add a new one if you want to use different methods for different samples. Then just repeat the steps above for the new method.

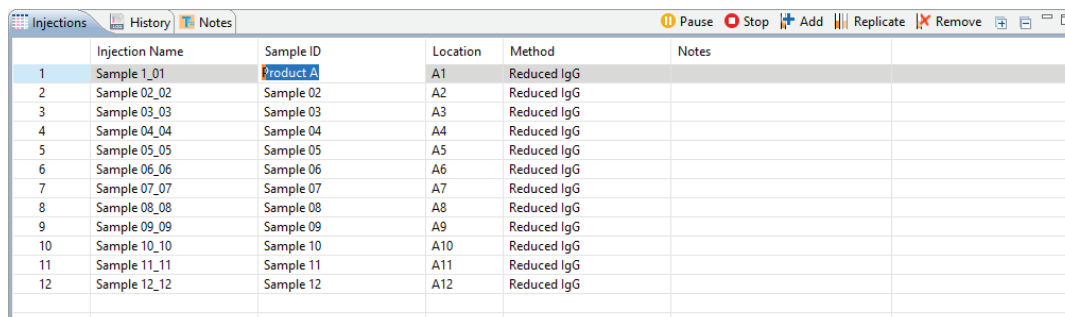
10. In the Injections pane:

- **To add or change sample names:** Click the **Sample ID** cell for the injection and type a name.

NOTES:

Sample IDs can be copied from other documents and pasted into Compass for iCE to make batch setup faster. Check out “Copying and Pasting Injection Names or Sample IDs from other Documents” on page 103 for more info.

If you change the sample ID when the default injection name is still being used, the injection name will automatically update to match the sample ID. If you’ve already changed the injection name manually, updating the sample ID will not update the injection name.



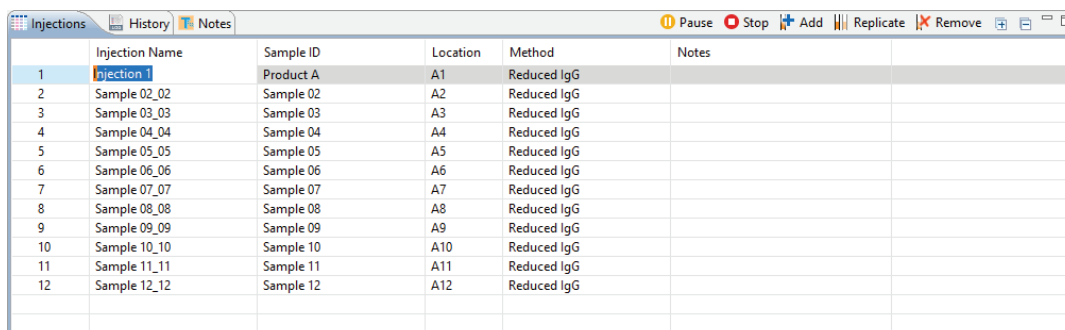
	Injection Name	Sample ID	Location	Method	Notes
1	Sample 1_01	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

- **To change injection names:** Click the **Injection Name** cell for the injection and type a name.

NOTES:

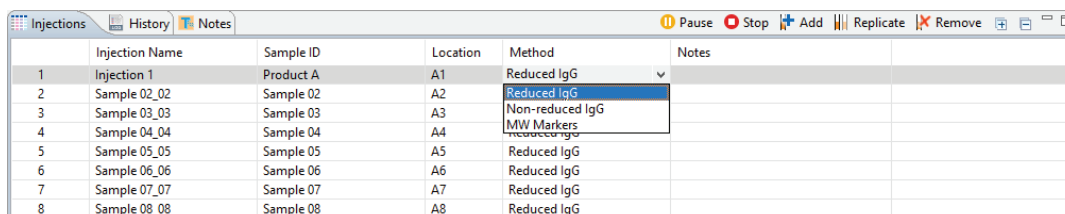
Each injection name must be unique.

Changing the injection name won’t affect the sample ID.



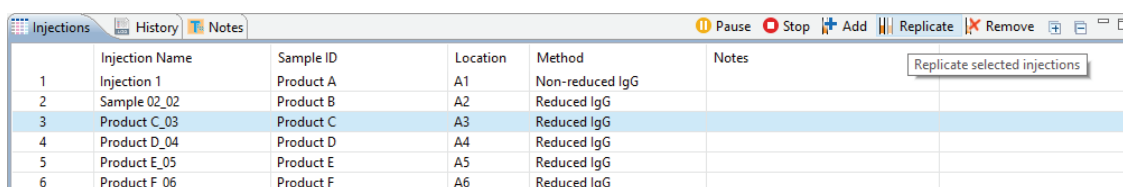
	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Reduced IgG	
4	Sample 04_04	Sample 04	A4	Reduced IgG	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	
9	Sample 09_09	Sample 09	A9	Reduced IgG	
10	Sample 10_10	Sample 10	A10	Reduced IgG	
11	Sample 11_11	Sample 11	A11	Reduced IgG	
12	Sample 12_12	Sample 12	A12	Reduced IgG	

- **To assign methods for each injection:** Click the **Method** cell for the injection and select a method from the drop down menu.

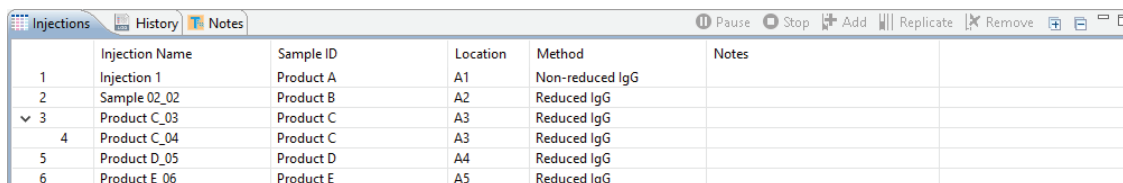


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Reduced IgG	
2	Sample 02_02	Sample 02	A2	Reduced IgG	
3	Sample 03_03	Sample 03	A3	Non-reduced IgG	
4	Sample 04_04	Sample 04	A4	MW Markers	
5	Sample 05_05	Sample 05	A5	Reduced IgG	
6	Sample 06_06	Sample 06	A6	Reduced IgG	
7	Sample 07_07	Sample 07	A7	Reduced IgG	
8	Sample 08_08	Sample 08	A8	Reduced IgG	

- **To add sequential replicate injections:** Highlight an injection and click **Replicate**. A new injection will be added under the row you selected

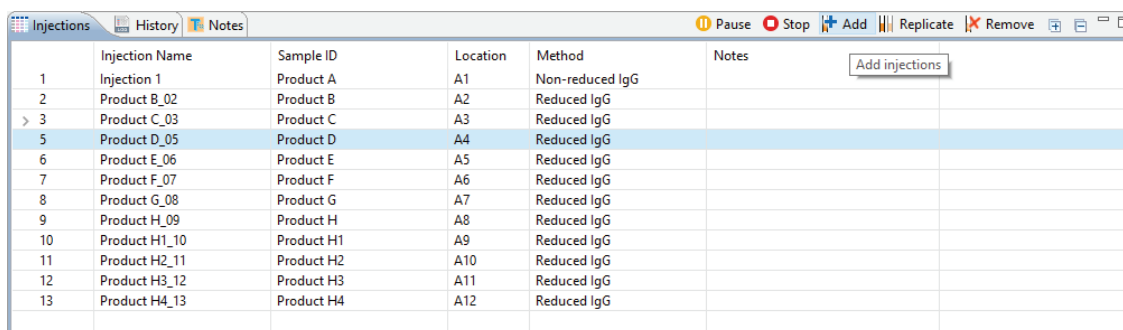


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product E_05	Product E	A5	Reduced IgG	
6	Product F_06	Product F	A6	Reduced IgG	



	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Sample 02_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product C_04	Product C	A3	Reduced IgG	
5	Product D_05	Product D	A4	Reduced IgG	
6	Product E_06	Product E	A5	Reduced IgG	

- **To add injections at the end of the batch:** Click on any sample position in the Layout pane or sample injection in the Injection table and click **Add**. A new injection using the same sample will be added to the end of the table. Assign the method if needed.

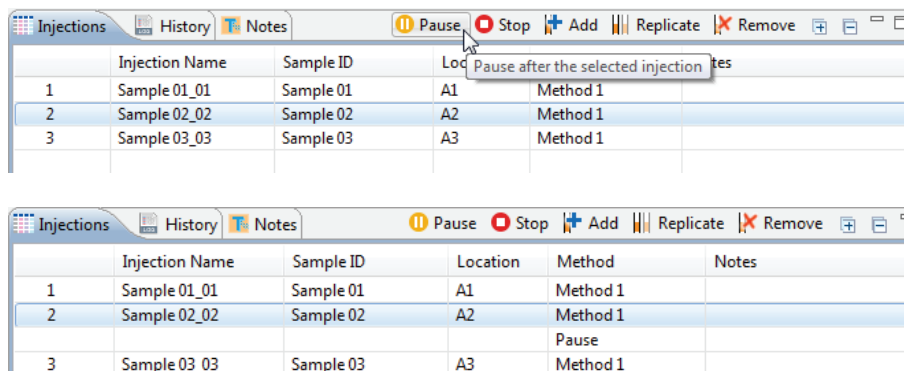


	Injection Name	Sample ID	Location	Method	Notes
1	Injection 1	Product A	A1	Non-reduced IgG	
2	Product B_02	Product B	A2	Reduced IgG	
3	Product C_03	Product C	A3	Reduced IgG	
4	Product D_04	Product D	A4	Reduced IgG	
5	Product D_05	Product D	A4	Reduced IgG	
6	Product E_06	Product E	A5	Reduced IgG	
7	Product F_07	Product F	A6	Reduced IgG	
8	Product G_08	Product G	A7	Reduced IgG	
9	Product H_09	Product H	A8	Reduced IgG	
10	Product H1_10	Product H1	A9	Reduced IgG	
11	Product H2_11	Product H2	A10	Reduced IgG	
12	Product H3_12	Product H3	A11	Reduced IgG	
13	Product H4_13	Product H4	A12	Reduced IgG	

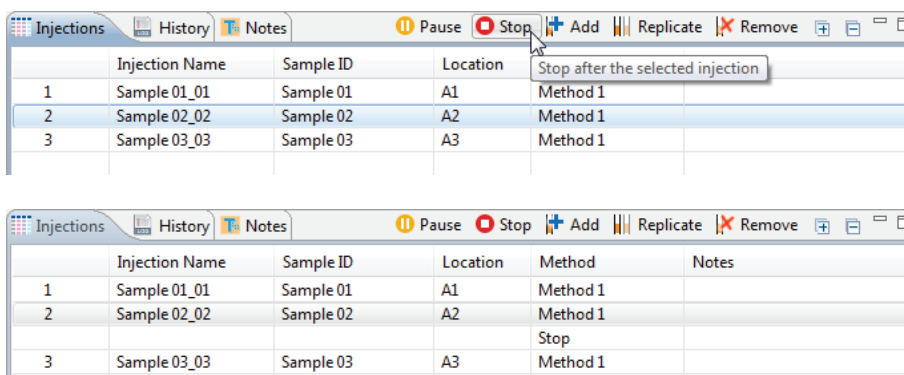
- **To copy and paste injections:** Right-click on an injection in the Injection table and select **Copy**. Click on the injection number above where you want to insert the copied injection, right click and select **Paste**. The copied injection will be inserted under the row you selected.

11. Add programmed pauses or stops in the batch to automatically pause or stop after injections (optional).

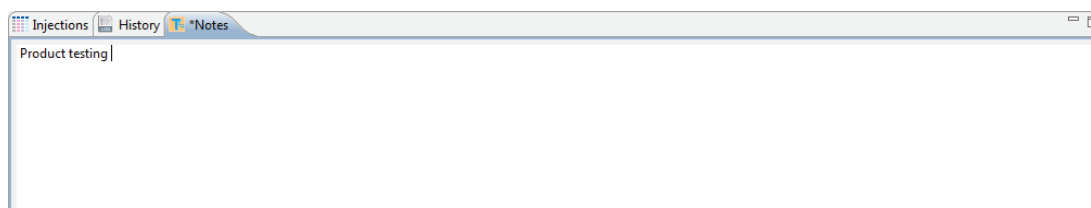
- **To program a pause:** Highlight the injection you want to pause at and click **Pause**. The batch will pause after that injection is complete. To resume the batch, click **Continue** in the instrument status bar.



- **To stop the run after a specific injection:** Highlight the injection you want the batch to stop at and click **Stop**. Maurice will continue the batch through the selected injection, then stop the batch and perform end-run cleanup steps.



12. Click on the **Notes** pane, then click in the notes area and type any information you want to add about your batch (optional).



13. You can preset the parameters used to analyze data generated with the batch (optional). We recommend using the default parameters for Turbo CE-SDS applications, but if you want to modify parameters:
- Select **Edit** from the main menu and click **Default Analysis**. The following screen will display:

Default Analysis: Maurice CE-SDS TURBO

Markers

Peak Names
Peak Fit
Advanced
Signal Processing

Analysis Groups

Standards

Add Remove

Apply Default:
Standards

Apply Override:

Apply To	Group
Sample	Standards

Add Remove

Markers

Internal Standard Time 170 Seconds

Markers Injection no markers

MW (kDa)	RMT
10	1
20	1.15
33	1.3
55	1.5
103	1.8
178	2.05
270	2.4

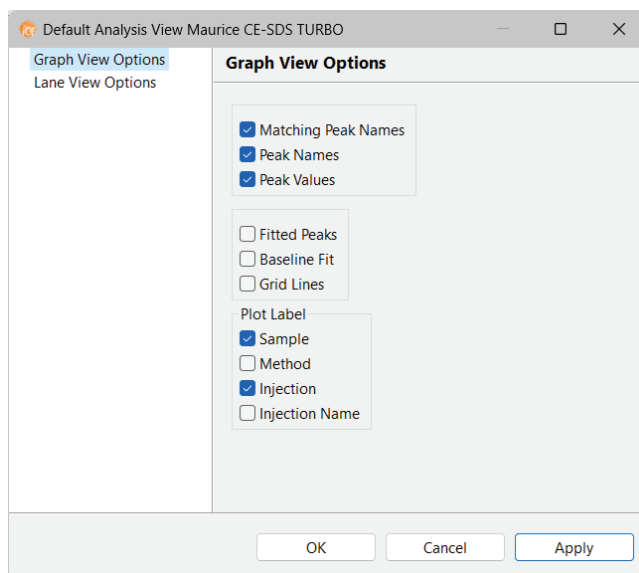
Add Remove

Import... Export... OK Cancel Apply

- Change the parameters you want to, then click **OK**. For detailed information on analysis parameters, please refer to “Analysis Settings Overview” on page 713.

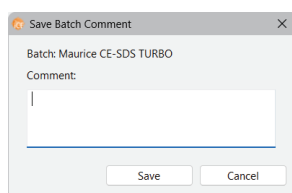
14. You can preset the parameters used to display the graph and lane view data when the batch is complete. Default Analysis View Parameters must be set before you start the batch and cannot be modified once the batch starts. However, you can still manually change parameters in the Analysis pane using the Graph Options and Lane Options buttons:

- a. Select **Edit** from the main menu and click **Default Analysis View**. The following screen will display:



- b. Change the parameters you want to, then click **OK**. For detailed information on graph and lane view options, please refer to “Customizing the Data Display” on page 695.

15. Once all of your sample, method and injection info is entered, select **File > Save**. Enter any comments on the batch if you want, then click **Save**.



16. Enter a name for your batch then click **Save**.

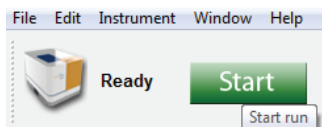
Step 7: Start the Batch

1. Make sure Maurice is connected to Compass for iCE. If he isn't, select **Instrument** and click **Connect**. Then select your Maurice system and click **Connect**.
2. Click **Start** to start your batch.

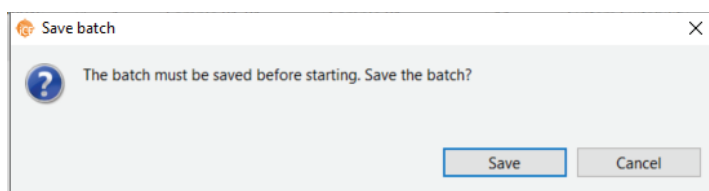
NOTES:

An alert may appear if the wrong cartridge is installed or if a cartridge post-run cleanup or cartridge purge is recommended. See “Step 5: Check for Cartridge Alerts” on page 369 for more information.

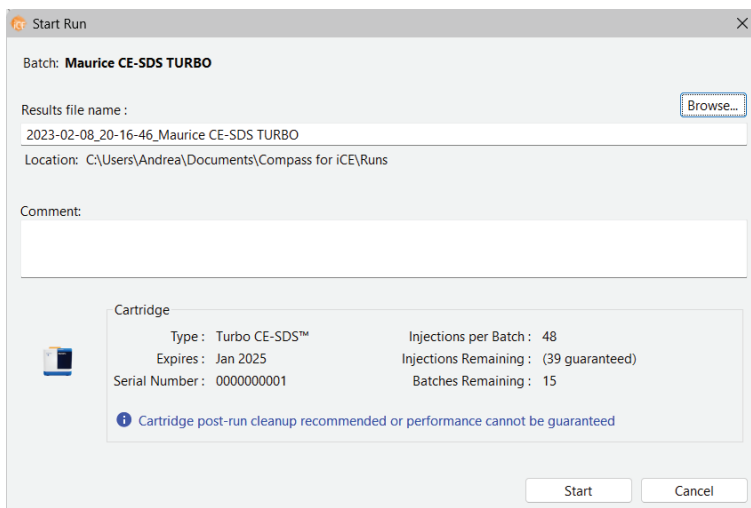
An alert may appear if the wrong adapter is installed or if you are starting a Turbo CE-SDS batch on MauriceFlex and have not locked the reagent vials in place. See “Adapter and Insert Alerts” on page 403 for more information.



3. You will be asked to save your batch before starting the run. Click **Save**.



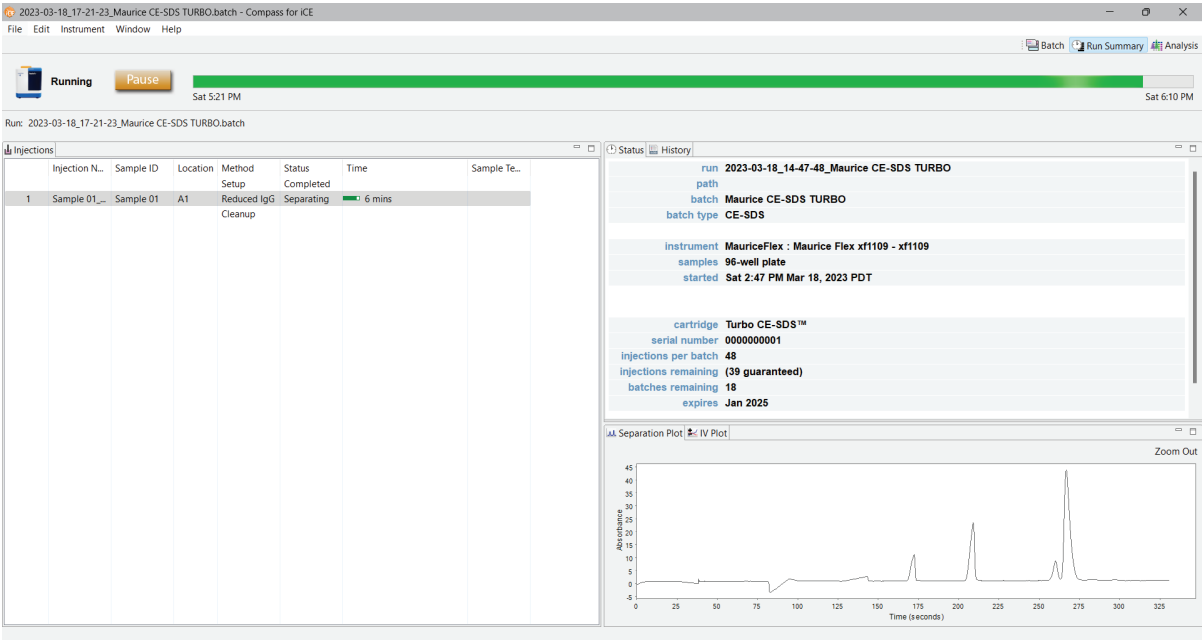
4. The Start Run box shows you the number of injections and batches remaining on your cartridge. Make sure you still have enough injections left before you start the run. If not, prep a new cartridge.
5. Click in the **Results File name** box if you want to change the default run file name. Otherwise just leave it as is.



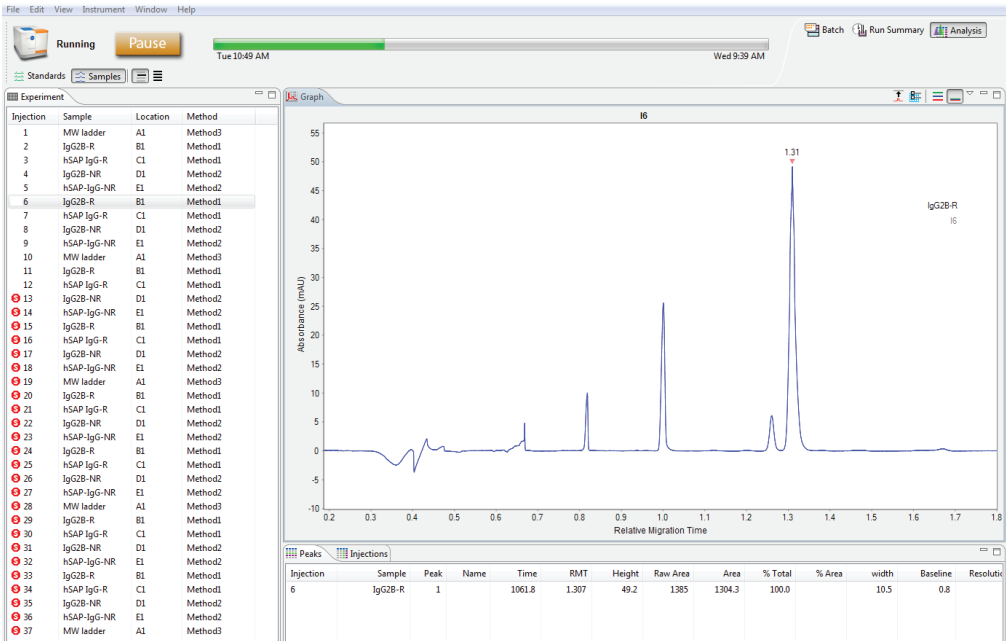
6. If you don't want to save the file to the default Runs folder, click **Browse** to select a different location.
7. Enter any run details you'd like in the Comments box (optional).
8. Click **Start** to start the run.

NOTE: The indicator light on Maurice’s front panel will pulse slowly while he’s running the batch.

To monitor the progress of your batch, click the **Run Summary** tab. See Chapter 16: “Run Status”for more details.



To view results, analyze results or change analysis parameters on completed injections while the batch is still running, click the **Analysis** tab. See Chapter 20: “CE-SDS Data Analysis” for more details.



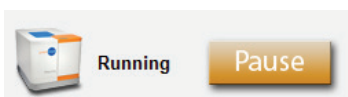
When your batch is complete, you can view electropherograms for all injections in the Analysis tab, and all batch and injection details in the Run Summary tab.

NOTE: If you'd like Maurice to let you know when your run is done, you can set him up to tweet you. Go to "Setting Up Maurice Systems to Send Tweets" on page 764 for more info.

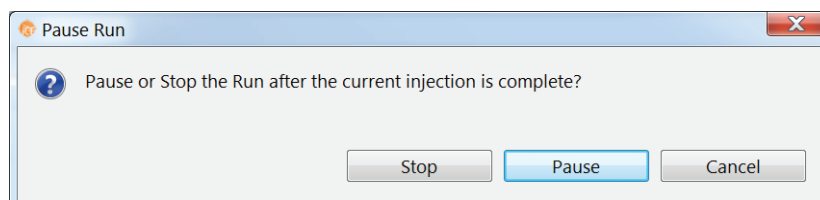
Editing a Running Batch

You can pause a batch after it's started to make updates or add samples and reagents.

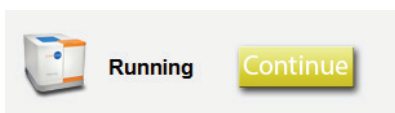
1. Select **Instrument** > **Pause** or select the **Pause** button.



2. Click **Pause** in the pop-up window.



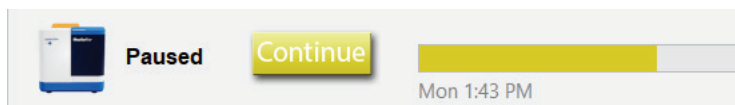
The Pause button will change to Continue but Maurice's status will stay at Running until Compass for iCE completes the current injection or calibration.



- You can now make these edits to the batch:
 - Add sample(s)
 - Add injection(s) to the end of the batch
 - Modify injections that haven't completed and aren't greyed out
 - Create a new method
 - Update the method. Once a method is run it can only be updated. Highlight the method you want to edit and click **Update** in the Methods pane. The software will copy the parameters into a new method that you can modify.

- When Maurice completes the current injection, his status will change to Paused and the progress bar will turn grey.

NOTE: Maurice's status light will change from a pulsing blue to a blue pulse with magenta flashes which means he's paused the run.



The status in the Injection pane also changes to Paused and the pause duration updates live. You can now open the instrument door to:

- Add new samples

IMPORTANT

The door must be closed within 15 minutes of opening it to avoid potential assay performance issues.

NOTE: If the door is open for more than 30 seconds, a white light near the cartridge will start to flash and a warning message displays in Compass for iCE until the door is closed.

- To continue the batch, click **Continue**. Changes you made to the batch are saved and Maurice will continue with the next sample injection.

NOTE: The D2 lamp is turned off for safety reasons when the door is opened. Once the batch is continued, Maurice will pause for 7 additional minutes before the next injection to let the lamp re-stabilize. During that time, the batch status in Compass for iCE will display injection loading.

Post-batch Procedures

When the batch is done:

- Open Maurice's door. The lights on either side of the cartridge slot will be **orange** as Maurice will have already disengaged the cartridge.
- If you're using the optional Maurice Filter Kit to contain β ME odor, cover the 96-well sample plate immediately with a plate lid.
- Remove your samples. Leave the Water (P3), Wash Solution (P5), and Air (P6) vials in place if your cartridge still has injection left as they will be needed for the cartridge post-run cleanup step. Discard the remaining reagent vials.
- Gently pull out the cartridge making sure the cartridge inlet doesn't come into contact with any surface.

NOTE: If you see any separation matrix sticking to the capillary inlet, soak it in DI water for 5 minutes. Then wipe it using a lint-free laboratory wipe that's been moistened with DI water.



If you're at 100 injections, you've reached the limit of guaranteed performance for the Turbo CE-SDS cartridge (note this is not necessarily the maximum injection limit). The cartridge will not be usable after 25 batches. To dispose of a finished cartridge, put it in its original packing and discard it per your institution's safety and waste disposal guidelines.

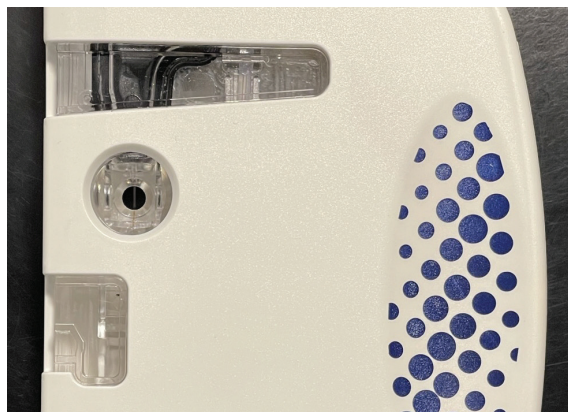
NOTE: Disposal of cartridges, sample plates and vials depends on the samples you ran. If you aren't sure what your sample origins are, we recommend you dispose of everything in biohazard waste.

If the cartridge will be used again, clean up and store the cartridge:

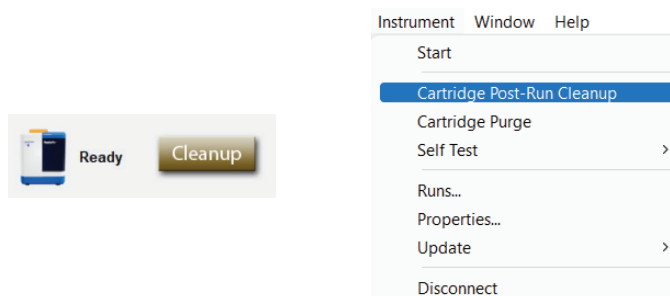
- Put the cartridge on a flat surface and remove the stopper from the Waste Tank.
- Tilt the cartridge so that all the liquid flows toward the Waste Tank opening, and aspirate out all of the liquid.
Optional: Before aspirating out the liquid, add 200 μ L of sodium hypochlorite solution (10-15%) to neutralize β -mercaptoethanol, then aspirate out all the liquid.
- Remove the stopper from the Top Chamber and aspirate out all the liquid.
- Dispense 2.7 mL of DI water into the Top Chamber to fill it. Aspirate out all the liquid. Repeat 2 more times.

NOTE: Tilt the cartridge so that all the liquid flows toward the Top Chamber opening while aspirating.

- e. Aspirate out all the remaining liquid. Use the transparent back of the Top Chamber to check if any liquid remains visible. Residual liquid can be removed at this point by tilting the cartridge.



- f. Verify that there is 1.5 mL of DI water in the Water (P3) vial, 1.5 mL of Wash Solution in the Wash Solution (P5) vial and air in the empty (P6) vial.
- g. Place the stoppers on the empty Waste Tank and Top Chamber. They should be firmly closed.
- h. Insert the cartridge in Maurice
- i. Click on the brown **Cleanup** button in the Instrument Status Bar or select **Instrument** and **Cartridge Post-Run Cleanup** in the Compass main menu.



- j. You'll get the following message. Click **Start**. It'll only take 7 minutes.



- k. Once the cleanup procedure is done, discard the reagent vials and remove the cartridge.

NOTES:

Keep cartridge upright while transporting.

See page 401 for information on what to do if a cartridge clog is detected during the Post-Run Cleanup.

- l. Place the cartridge on a flat surface and remove the stopper from the Waste Tank.
- m. Tilt the cartridge so that all the liquid flows toward the Waste Tank opening, and aspirate out. Leave the stopper off to allow to air dry.
- n. Remove the stopper from the Top Chamber and aspirate out all visible liquid. Leave the stopper off to allow to air dry.
- o. Put the cartridge and stoppers in its protective packaging and store at room temperature.

!WARNING! SHARPS HAZARD

The capillary inlet of the cartridge may present a potential sharps hazard. Dispose of used cartridges according to your organizations health and safety regulations.



!WARNING! BIOHAZARD

Cartridges, sample plates and vials should be handled by procedures recommended in the CDC/NIH manual: Biosafety in Microbiological and Biomedical Laboratories (BMBL). The manual is available from the U.S. Government Printing Office or online at <http://www.cdc.gov/biosafety/publications/bmbl5/>.

Depending on the samples used, the cartridges, plates and vials may constitute a chemical or a biohazard. Dispose of the cartridges, plates and vials in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations. Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste content before you store, handle, or dispose of chemical waste.

Checking Your Data

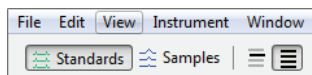
Compass for iCE detects your sample proteins, CE-SDS MW Markers and Internal Standard peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

Step 1: Check Your Internal Standard

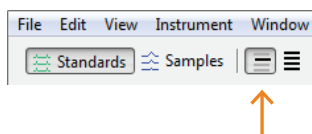
To make sure your Internal Standard is identified correctly:

1. Go to the **Analysis** screen and open your run (if it isn't already open).

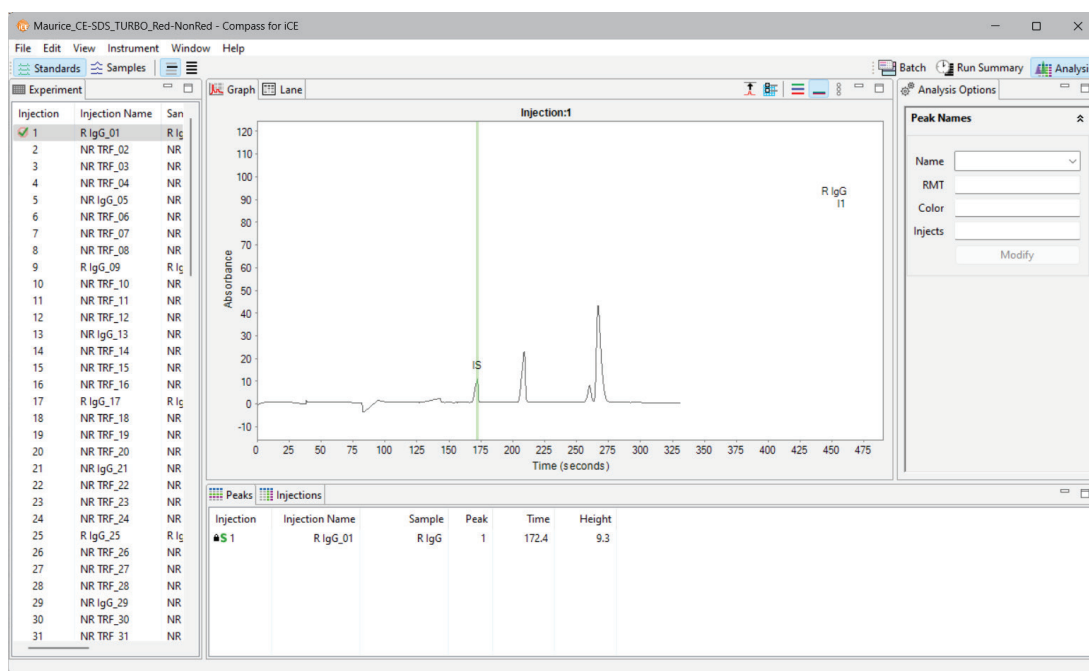
- Click **Standards** in the View bar.



- Click the **View Selected** icon in the View bar.

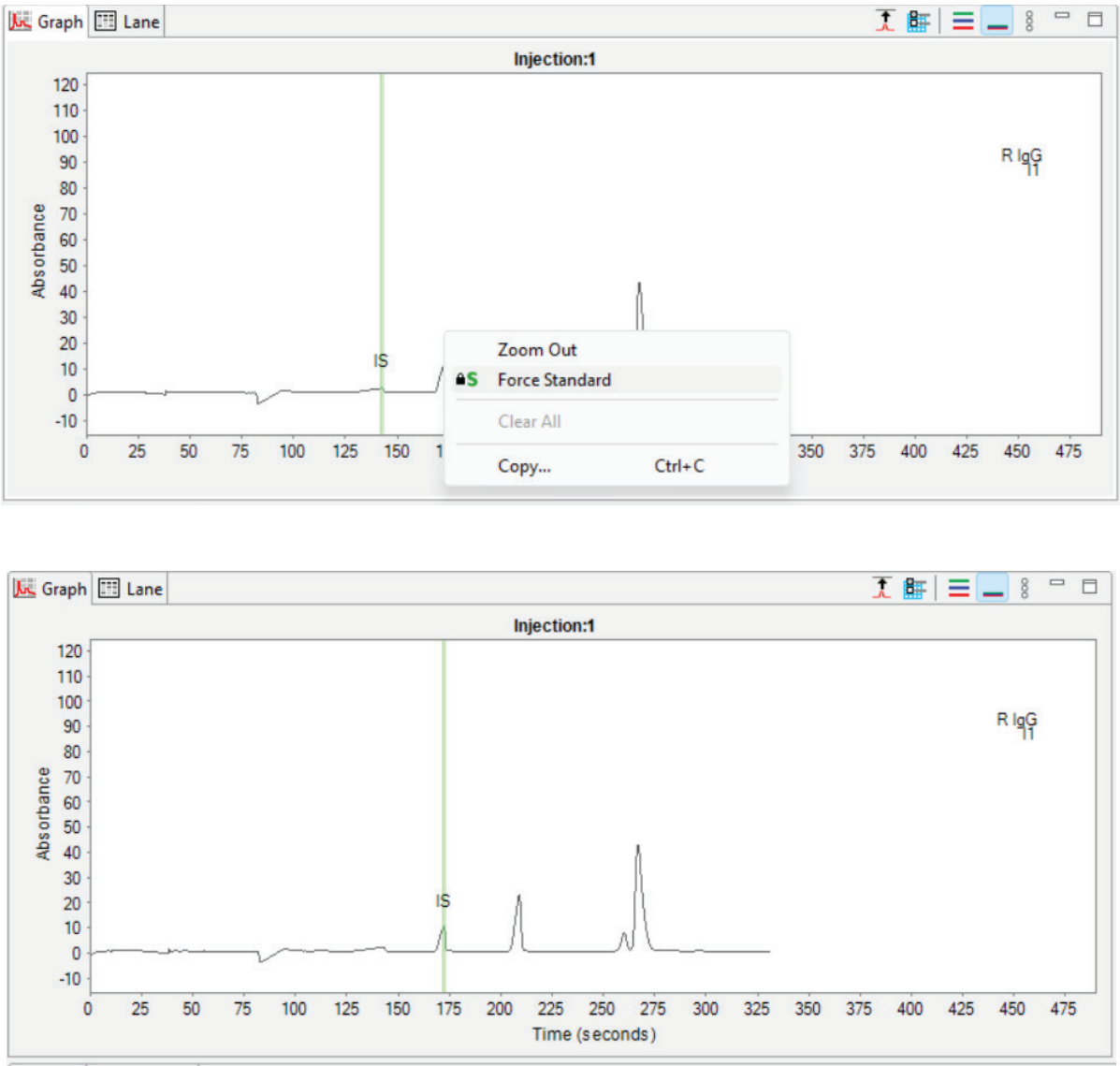


- Click **Injection 1** in the Experiment pane.
- Check that your Internal Standard in the electropherogram has been correctly identified. It'll be labeled IS and will have a green vertical line running through it. The Internal Standard is also identified with an **S** in the Peaks table.



6. If your Internal Standard isn't identified correctly, here's how to manually correct it:

To set an unidentified peak as the Internal Standard: Right-click the peak in the electropherogram or Peaks table and select **Force Standard**. Compass will assign that peak as the Internal Standard.



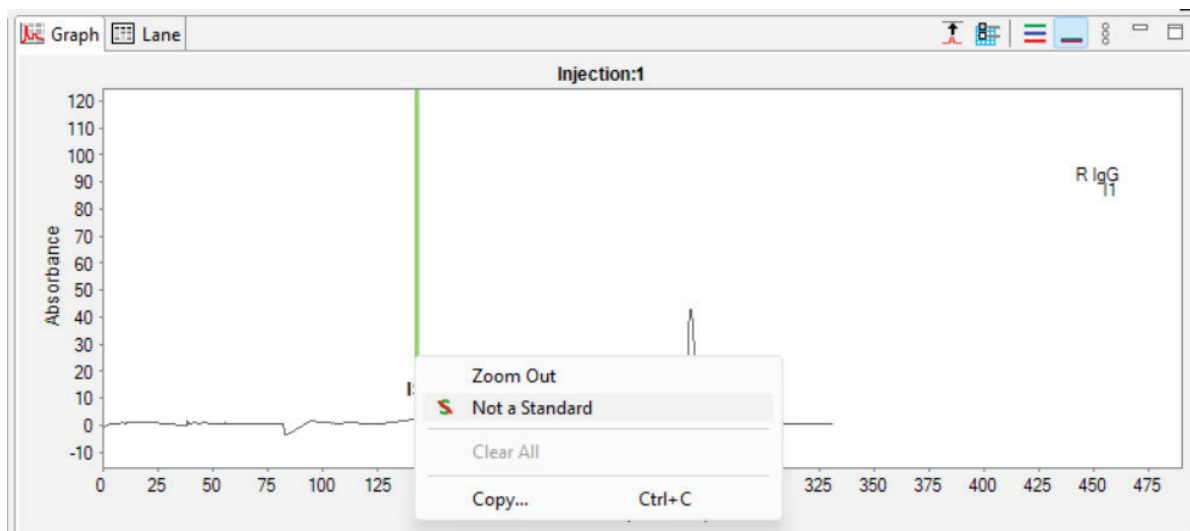
A lock icon indicating the Internal Standard was set manually will display next to the peak in the Peaks table and a check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Peaks			Injections		
Injection	Sample	Peak	Injection	Sample	Peak
1	Sample 1	11	1	Sample 1	12
1	Sample 1	13	1	Sample 1	14
1	Sample 1	15	1	Sample 1	16
1	Sample 1	17			

Experiment			
Injection	Sample	Location	Method
✓ 1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1

NOTE: To remove an Internal Standard peak assignment that was made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**. To clear all the manual settings for the injection, click **Clear All**.

If an incorrect peak is identified as the Internal Standard: Right-click the peak in the electropherogram or Peaks table and select **Not a Standard**.



A check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Experiment			
Injection	Sample	Location	Method
✓ 1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1

- Repeat the previous steps for all other injections to make sure your Internal Standard is identified correctly.

Step 2: Set Your Molecular Weight (MW) Markers

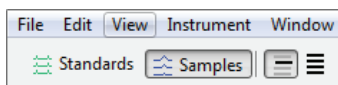
NOTE: You'll only need to do this if you ran the CE-SDS MW Markers. If you didn't, you can skip to the next section.

Compass reports the relative migration time (RMT) of your sample in the Peaks table. If you also want to know the relative molecular weight of your sample, you can run the CE-SDS MW Markers as one of your injections.

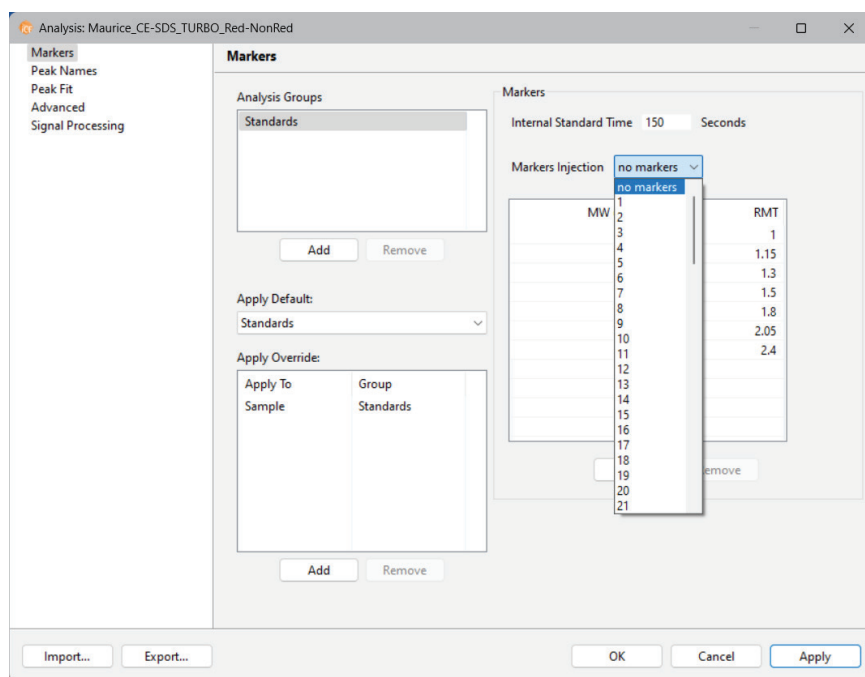
You'll see these sizing markers when you run the CE-SDS MW Markers: 10, 20, 33, 55, 103, 178, and 270 kDa.

To get MW data:

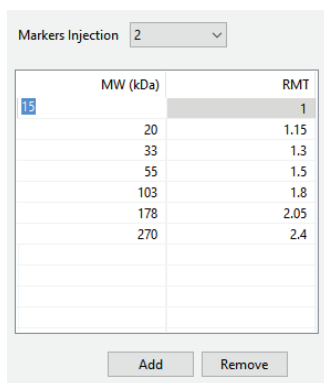
1. Click **Samples** in the View bar.



2. Select **Edit** from the main menu and click **Analysis**. In the Analysis window, select **Markers** in the left sidebar. Then click the **Markers Injection** drop down menu to select the injection you ran your CE-SDS MW Markers in.

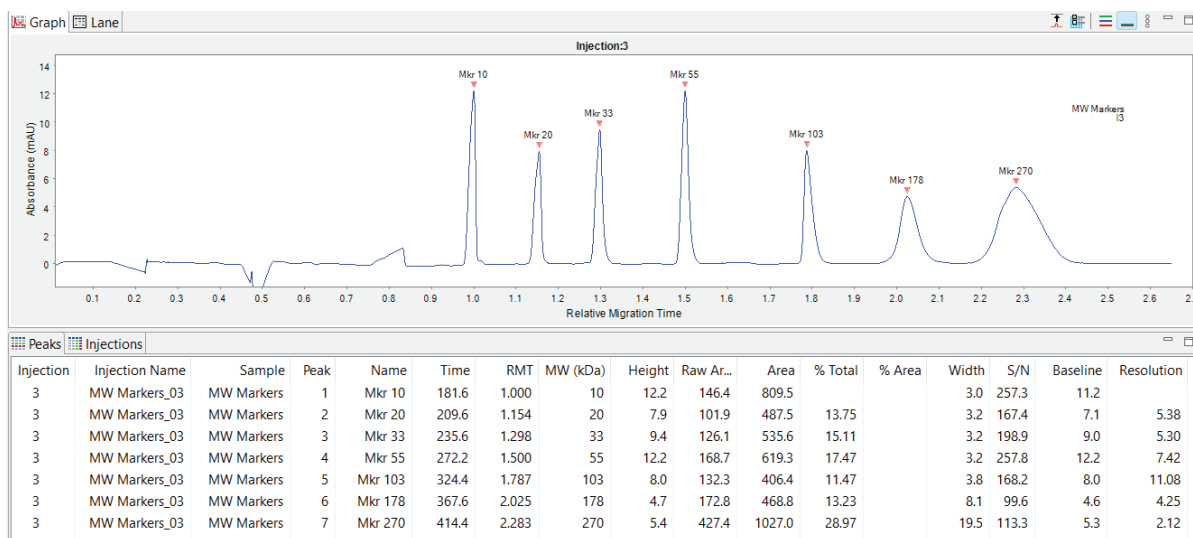


3. The default Maurice CE-SDS MW Markers molecular weights and relative migration time (RMT) values are already populated in the table. If you'd like to use these values, skip to the next step. If you're using different markers, click in the MW and RMT cells to type new values, click a row and select **Remove** to delete, or click **Add** to add a new one.



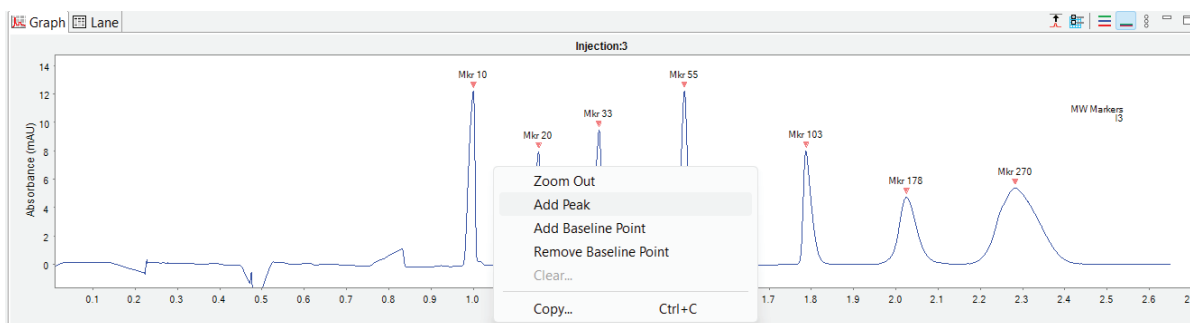
- Click **OK** to close the Analysis window. Compass will automatically assign the molecular weights to your markers and label them Mkr. A MW (kDa) column will also now display in the Peaks table.

NOTE: The Mkr 10 peak is also the Internal Standard in every sample.



- It's always a good idea to verify that all your CE-SDS MW Markers are identified correctly. Here's how to manually correct them:

To set an unidentified peak as a MW Marker: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will assign that peak as a MW Marker, and correctly reassign the remaining marker peaks.

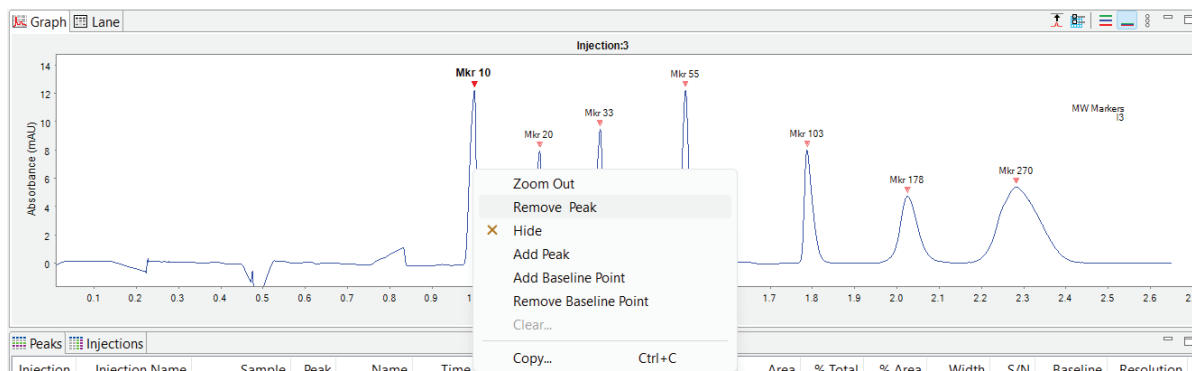


A check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Experiment			
Injection	Sample	Location	Me
1	IgG System Co...	A1	Me
✓ 2	Control Ladder	A2	Me
3	Test Ladder	A3	Me
4	IS - Alpha	B1	Me

NOTE: To remove MW Marker peak assignments that were made manually, right click on the peak in the electropherogram or Peaks table and click **Clear**.

If an incorrect peak is identified as a MW Marker: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass should correctly reassign the remaining peaks as markers and update the Peaks table.



A check mark will appear next to the injection in the Experiment pane to show a manual correction was made.

Experiment			
Injection	Sample	Location	Me
1	IgG System Co...	A1	Me
✓ 2	Control Ladder	A2	Me
3	Test Ladder	A3	Me
4	IS - Alpha	B1	Me

Step 3: Checking Sample Peaks

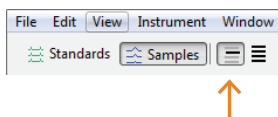
All detected peaks will be labeled automatically with the RMT (default) or apparent MW (if the CE-SDS MW Markers were run).

To make sure your sample proteins are identified correctly:

1. Click **Samples** in the View bar.

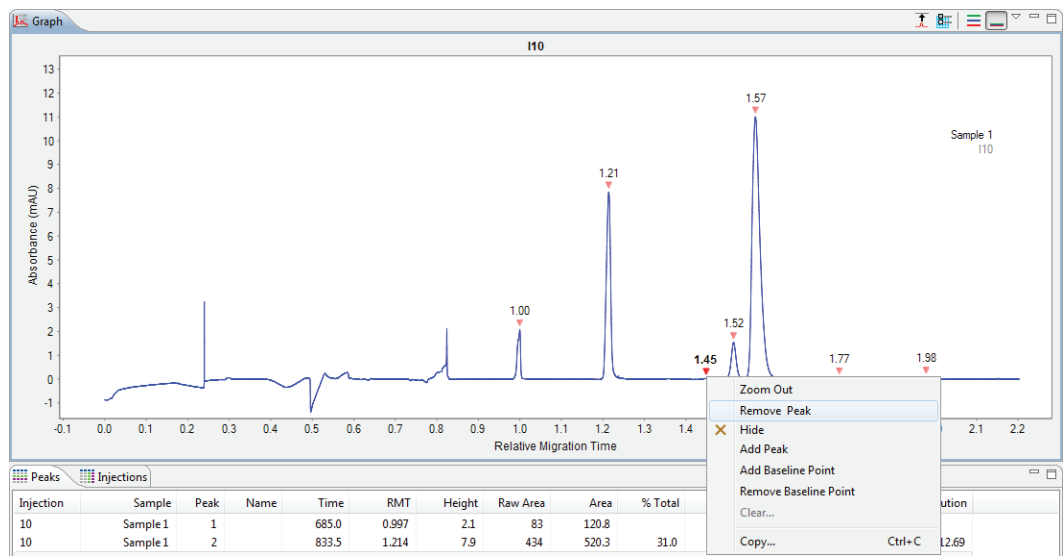


2. Click the **View Selected** icon in the View bar.



- 3. Click **Injection 1** in the Experiment pane.
- 4. If your sample peaks aren't identified correctly, here's how to manually correct them:

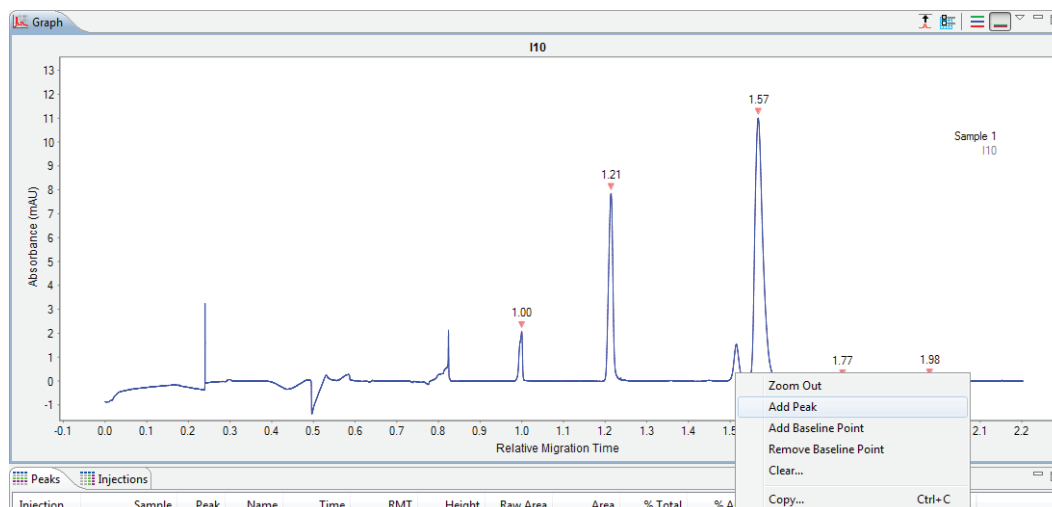
If a peak is incorrectly identified as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Remove Peak**. Compass will no longer identify it as a sample peak in the electropherogram and the peak data will be removed in the results table.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment			
Injection	Sample	Location	Method
1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1
8	Sample 1	A1	Method1
9	Sample 1	A1	Method1
✓ 10	Sample 1	A1	Method1
11	Sample 1	A1	Method1
12	Sample 1	A1	Method1

To identify a peak as a sample peak: Right-click the peak in the electropherogram or Peaks table and select **Add Peak**. Compass will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Injection	Sample	Location	Method
1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1
8	Sample 1	A1	Method1
9	Sample 1	A1	Method1
✓ 10	Sample 1	A1	Method1
11	Sample 1	A1	Method1
12	Sample 1	A1	Method1

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear**.

- Repeat the previous steps for the remaining injections to make sure all sample peaks are correctly identified.

Step 4: Assigning Peak Names

Compass can also optionally identify and automatically name sample peaks using user-specified peak name settings. For more information on how to do this, see “Manual Peak Integration” on page 739.

Chapter 15:

Cartridge and Adapter Alerts

Chapter Overview

- Overview
- Cartridge Alerts
- Adapter and Insert Alerts
- Compass Alerts

Overview

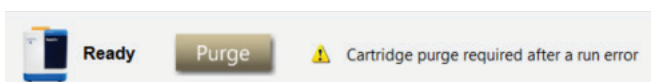
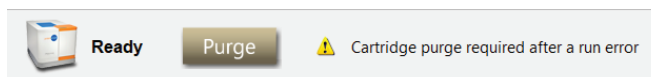
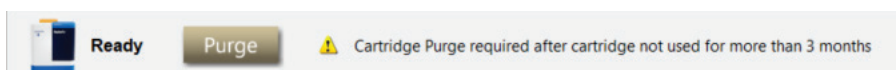
Maurice will double check that everything is set up properly when you install your cartridge, start your batch or pause your run.

Cartridge Alerts

Cartridge Maintenance

An alert may appear when you install your cartridge in Maurice if a cartridge purge or cartridge post-run cleanup is recommended for your cartridge.

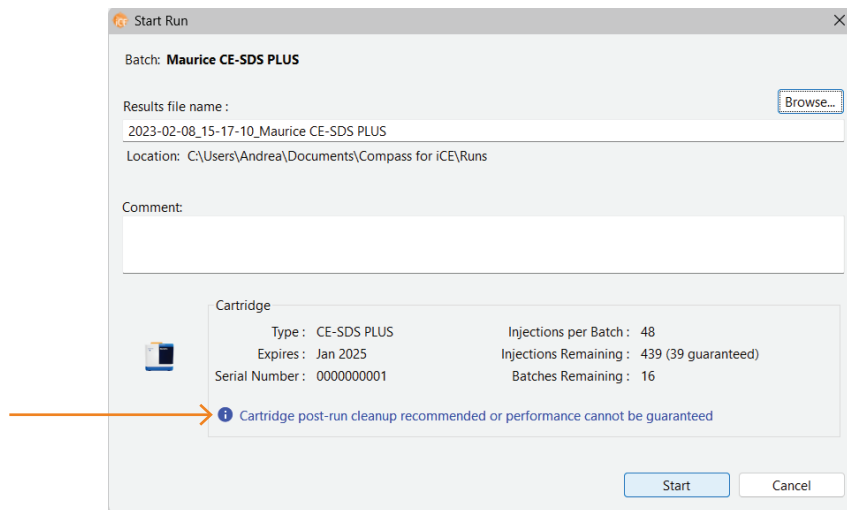
- If you install a cIEF or cIEF Fractionation cartridge that was used previously in a run that had an error, or a CE-SDS PLUS or Turbo CE-SDS cartridge that has not been used in 3 months or was used previously in a run that had an error, an alert that a cartridge purge is required will appear next to a brown Purge button. Run a cartridge post-run cleanup after performing a cartridge purge.
- See “Step 5: Check for Cartridge Alerts” for cIEF applications, page 166 for MauriceFlex cIEF applications, page 232 for MauriceFlex Fractionation applications, page 300 for CE-SDS PLUS applications, and page 369 for Turbo CE-SDS applications for more information.



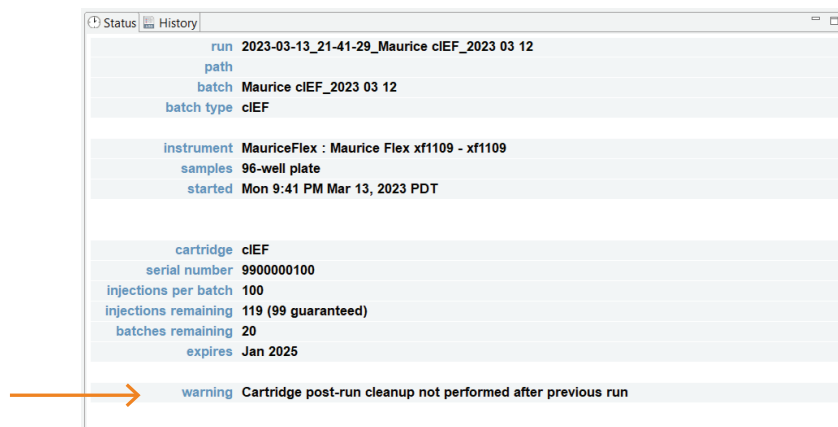
- If you install a cartridge that did not undergo a post-run cleanup after the last batch, an alert will appear next to a brown Cleanup button recommending you perform a cartridge post-run cleanup. The green Start button will appear after successful cleanup.
- See page 120 for cIEF applications, page 182 for MauriceFlex cIEF applications, page 248 for MauriceFlex Fractionation applications, page 316 for CE-SDS PLUS applications, and page 385 for Turbo CE-SDS applications for more information.



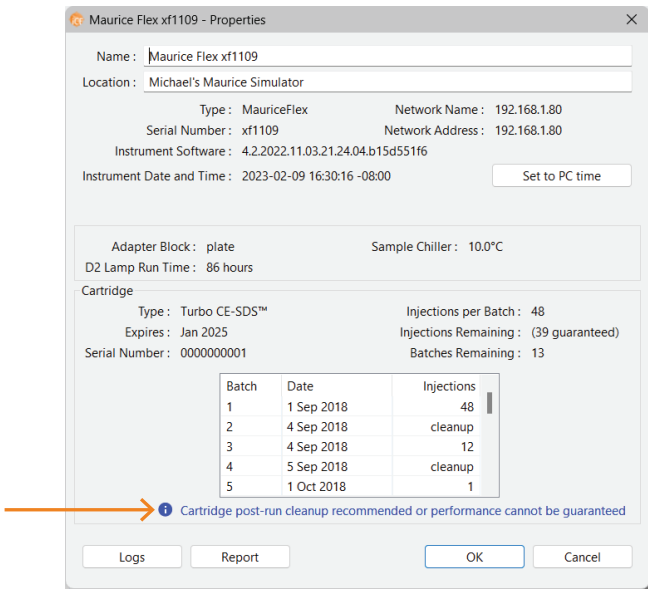
If you start the batch without performing the post-run clean up, a warning will appear in the Start Run window.



A warning that the post-run cleanup was not performed will appear at the bottom of the Run Summary pane if you proceed with the batch without running the post-run cleanup.

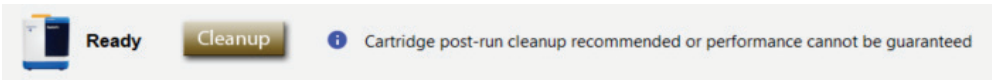


The warning will also be recorded in the Cartridge Properties and Injection Report.

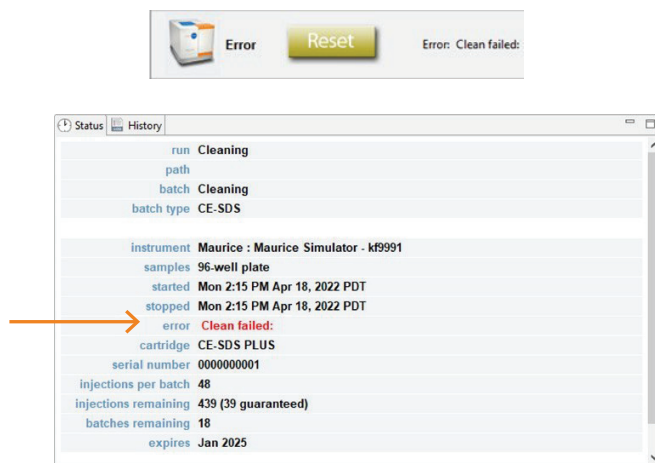


Model	Maurice OBM
Instrument S/N	kf1147
Software Version	Compass for iCE 3.0.0, Build ID: 0124
Firmware Version	4.1.2021.12.09.23.23.49.55b8e524f
Tray Type	96-well plate
Cartridge Type	Turbo CE-SDS™
Cartridge S/N	8211213001
Cartridge Expiration	Dec 2022
Injections Remaining	(4 guaranteed)
Batches Remaining	20
Warning	Cartridge post-run cleanup not performed after previous run

At the end of your batch, an alert will appear next to a brown Cleanup button reminding you to perform a cartridge post-run cleanup. The green start button will appear after successful cleanup.



- If the cartridge post-run cleanup was stopped during the cleanup or did not complete successfully, and error will appear next to a yellow Reset button.
- Click the yellow **Reset** button to clear the error.
- You can try running the post-run cleanup again, start a run without the cleanup or prepare another cartridge to use for your run.
- The error will be recorded in the Run Summary Status pane and Cartridge properties.



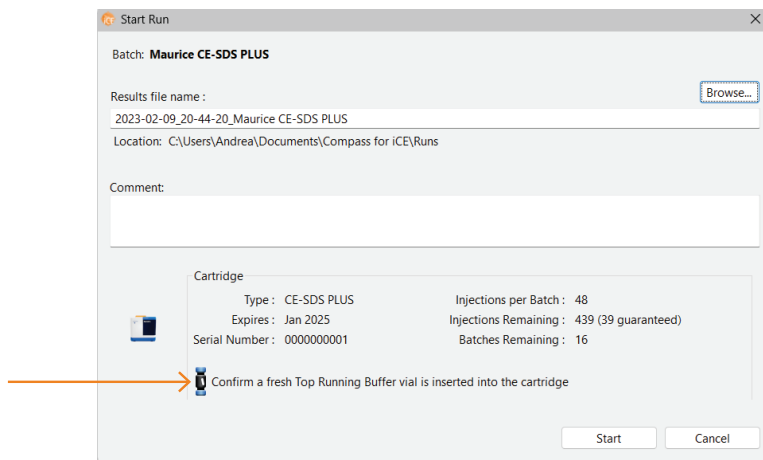
- If the cartridge post-run cleanup for your Turbo CE-SDS cartridge completed successfully but a clog is detected, an error will appear next to the Reset button and in the Run Summary Status pane.

When this happens

- Purge the cartridge. See page 369 for more information
- Perform a post-run cleanup and store the cartridge. See page 385 for more information on performing a post-run cleanup and storing the cartridge.
- If the purge or the post-run cleanup is not successful, contact Technical Support.

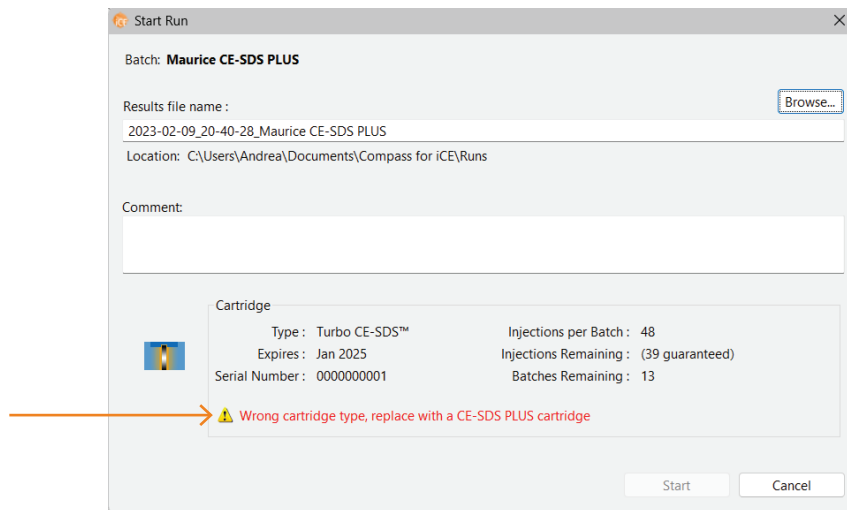
Cartridge Prep

An alert will appear in the Start Run window to ensure the cartridge has been correctly prepped.



Cartridge Type

An alert will appear in the Start Run window if the cartridge type that is installed in Maurice does not match your batch type. You will not be able to start the batch until the correct cartridge is installed.



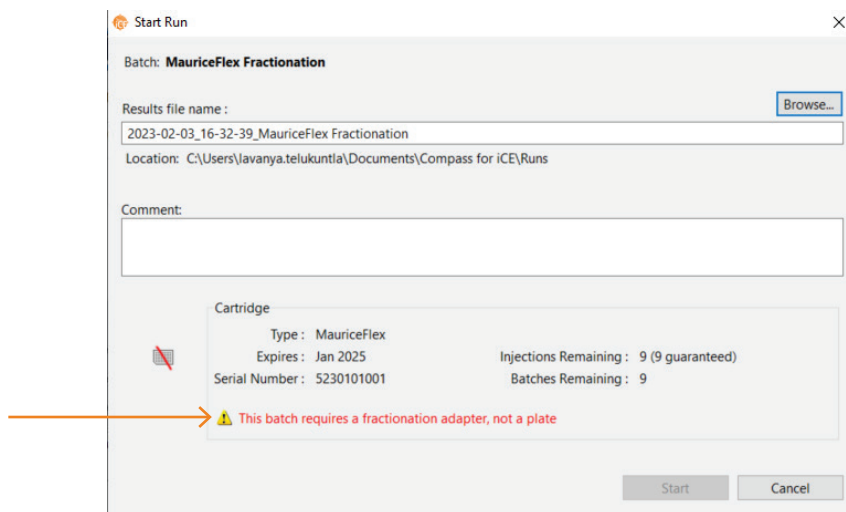
You will also see an alert in the Start Run window if no cartridge is detected. You will not be able to start the batch until a cartridge is installed.

Adapter and Insert Alerts

Maurice will check to ensure the correct sample or vial insert or adapter is installed before you start a batch or if you re-start a paused batch. .

Alerts when Starting a Batch

Alerts will appear in the Start Run window.



If you are running a standard cIEF or CE-SDS application on Maurice, Maurice C. and Maurice S., Compass for iCE will check to make sure that:

- The metal plate insert is installed if your sample is prepared in a 96-well plate.
- The metal vial insert is installed if your samples are prepared in vials.

If you are running a standard cIEF or CE-SDS application on MauriceFlex, Compass for iCE will check to make sure that:

- The fractionation adapter is not installed.
- The metal plate insert is installed if your sample is prepared in a 96-well plate.
- The metal vial insert is installed if your samples are prepared in vials.
- The reagent vials are locked in the reagent platform.

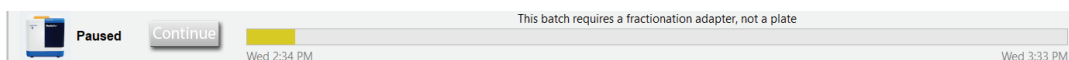
If you are running a MauriceFlex cIEF or MauriceFlex Fractionation application on MauriceFlex, Compass for iCE will check to make sure that:

- The fractionation adapter is installed.

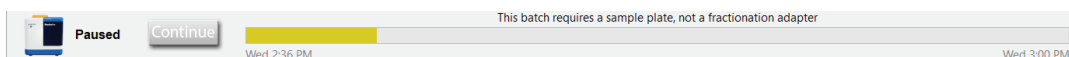
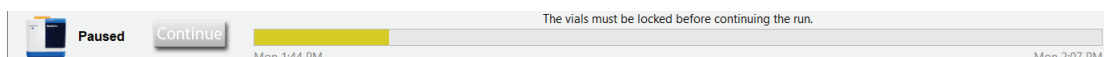
Alerts when Pausing a Batch

Alerts may appear when your batch is paused and click the yellow **Continue** button. .

- When you pause a standard cIEF or CE-SDS batch, an alert may appear if correct metal insert is not detected.



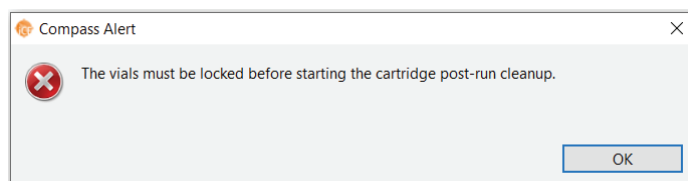
- When you pause a running MauriceFlex batch, an alert may also appear if the batch reagent vials are not locked or if the fractionation adapter is not detected.



Compass Alerts

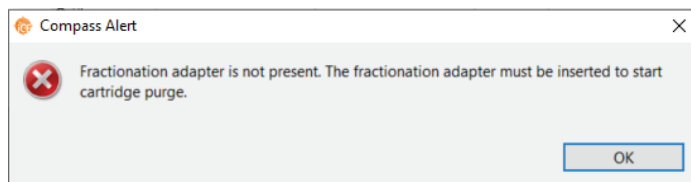
A Compass Alert window may appear during the Cartridge Post-Run Clean, Cartridge Purge, Cartridge Self-Test and Instrument Self-Test if you have the wrong cartridge, insert or adapter installed.

When performing a Cartridge Post-Run Cleanup, Compass Alerts will remind you to:



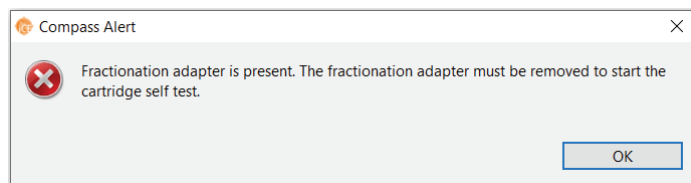
- Install a cartridge
- Lock reagent vials when cleaning a cIEF, CE-SDS PLUS or Turbo CE-SDS cartridge on MauriceFlex
- Install the fractionation adapter when cleaning a cIEF Fractionation cartridge on MauriceFlex
- Remove the fractionation adapter when cleaning a cIEF, CE-SDS PLUS or Turbo CE-SDS cartridge on MauriceFlex

When performing a Cartridge Purge, Compass Alerts will remind you to:



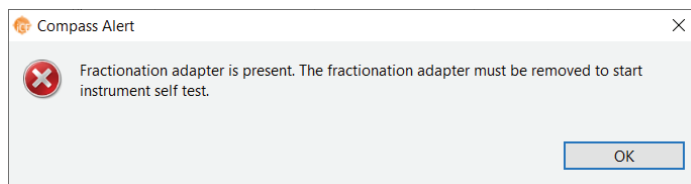
- Install a cartridge
- Lock reagent vials when purging a cIEF, CE-SDS PLUS or Turbo CE-SDS cartridge on MauriceFlex
- Install the fractionation adapter when purging a cIEF Fractionation cartridge on MauriceFlex
- Remove the fractionation adapter when purging a cIEF, CE-SDS PLUS or Turbo CE-SDS cartridge on MauriceFlex

When performing a Cartridge Self-Test, Compass Alerts will remind you to:



- Install a cartridge
- Lock reagent vials when testing a cIEF, CE-SDS PLUS or Turbo CE-SDS cartridge on MauriceFlex
- Remove the fractionation adapter when testing cartridges on MauriceFlex

When performing an Instrument Self-Test, Compass Alerts will remind you to:



- Remove the cartridge
- Remove the fractionation adapter when testing MauriceFlex

Chapter 16:

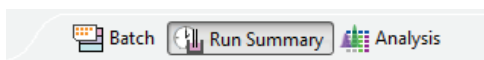
Run Status

Chapter Overview

- Run Summary Screen Overview
- Opening Run Files
- Batch Injection Information
- Run Status Information
- Viewing the Injection Focus Series (cIEF, MauriceFlex cIEF, MauriceFlex Fractionation Only)
- Viewing the Fractionation Focus Series (MauriceFlex Fractionation Only)
- Viewing the Separation Plot (CE-SDS PLUS and Turbo CE-SDS Only)
- Current and Voltage Plots
- Run History
- Viewing Multiple Events
- Viewing Run Errors or Warnings
- Injection Reports
- Switching Between Open Run Files
- Closing Run Files

Run Summary Screen Overview

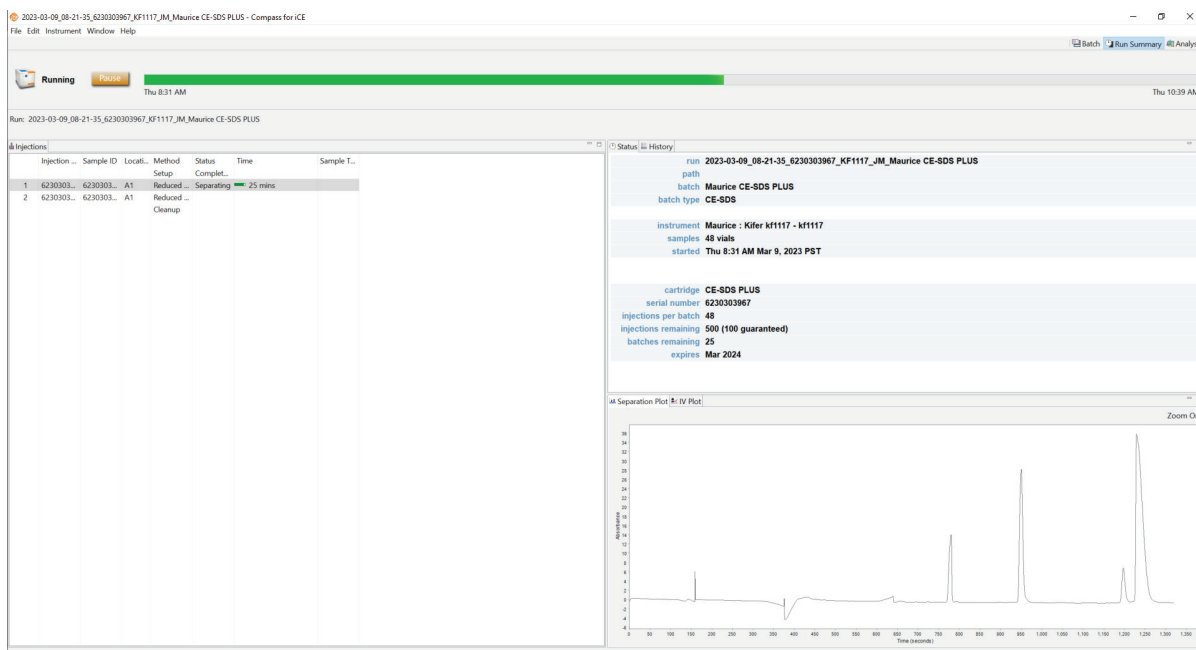
You can use the Run Summary screen to monitor the status of a batch in progress, see the CE-SDS PLUS and Turbo CE-SDS separation or cIEF, MauriceFlex cIEF and MauriceFlex Fractionation Focus series for your injections or the current and voltage plots for each injection. To get to this screen, click the **Run Summary** screen tab:

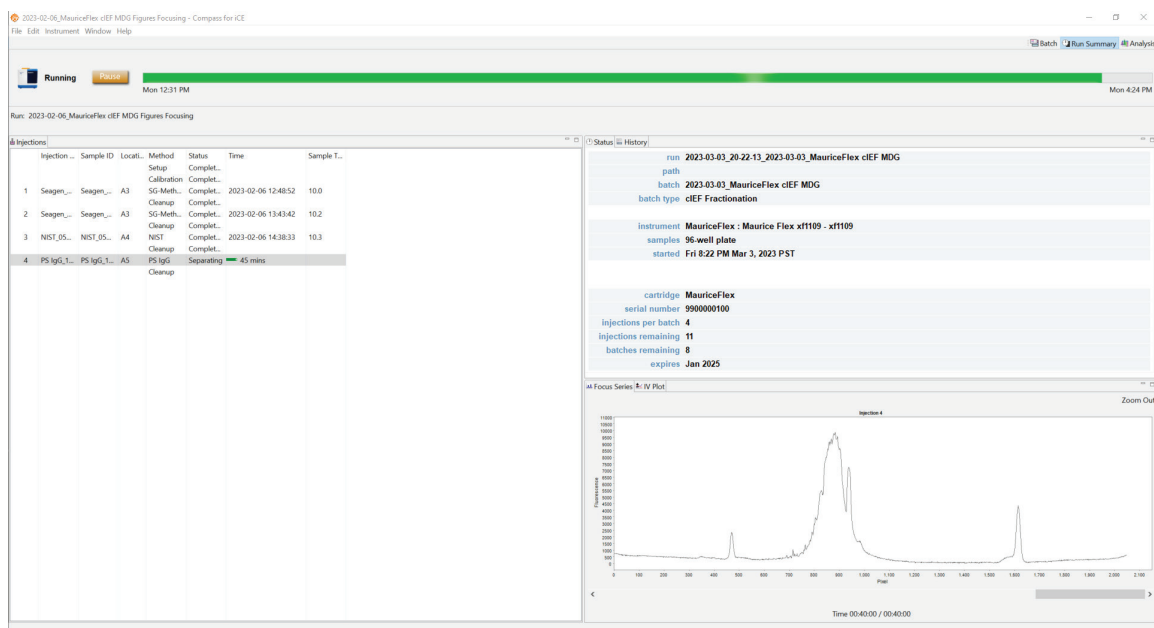
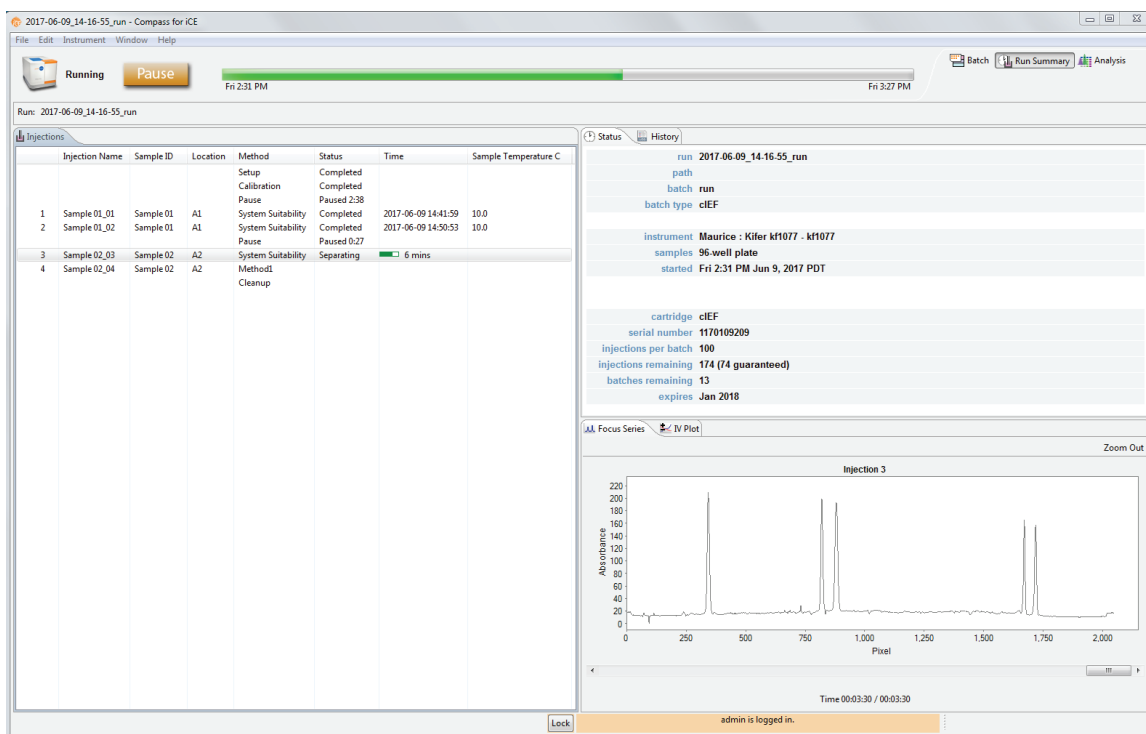


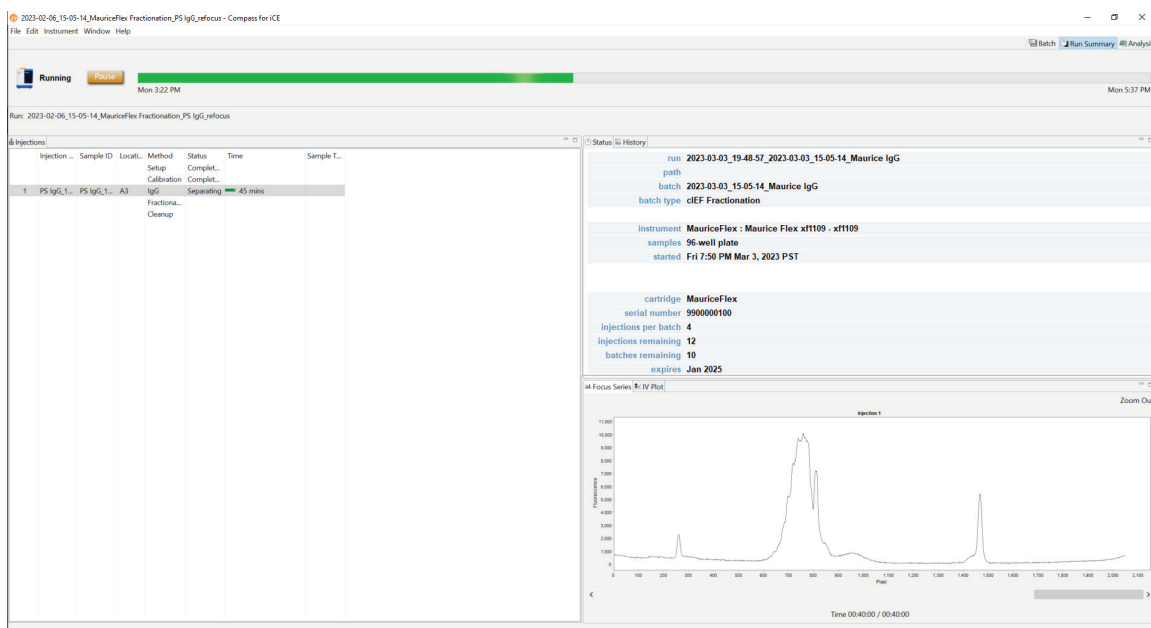
Run Summary Screen Panes

The Run Summary screen has five panes:

- **Injections** - Lists the sample IDs, sample locations and methods used for each injection in the run. It also shows the status of the current injection if a run is in progress.
- **Status** - Displays run file information and the current status of a run if one's in progress.
- **History** - Running history of all run file events from when the run was first started to the most current analysis update.
- **Separation Plot (CE-SDS PLUS and Turbo CE-SDS only)**- Lets you view the raw protein separation in the capillary for each injection.
- **Focus Series (cIEF, MauriceFlex cIEF and MauriceFlex Fractionation only)** - Lets you view the recorded focusing of proteins along the pH gradient in the capillary for each injection.
- **IV Plot** - Lets you view plots of the total current and voltage measured during the separation for each injection.







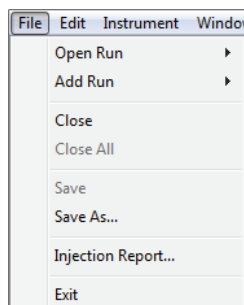
Software Menus Active in the Run Summary Screen

These main menu items are active in the Run Summary screen:

- File
- Edit
- Instrument (when the software is connected to an instrument)
- Window
- Help

File Menu

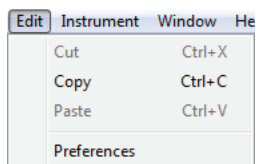
These File menu options are active:



- **Open Run** - Opens a run file.
- **Add Run** - Lets you open and view other run files besides the one that's already open.
- **Close** - Closes the run file currently being viewed.
- **Close All** - Closes all open run files.
- **Save/Save As** - If you made changes in the Analysis screen before you went to the Run Summary screen, this saves your changes to the run file.
- **Injection Report** - Exports the raw and analyzed data, IV plot, peaks table, sample and system info for individual injections as PDF files. You can also export the run history with all analysis events.
- **Exit** - Closes Compass for iCE.

Edit Menu

These Edit menu options are active:



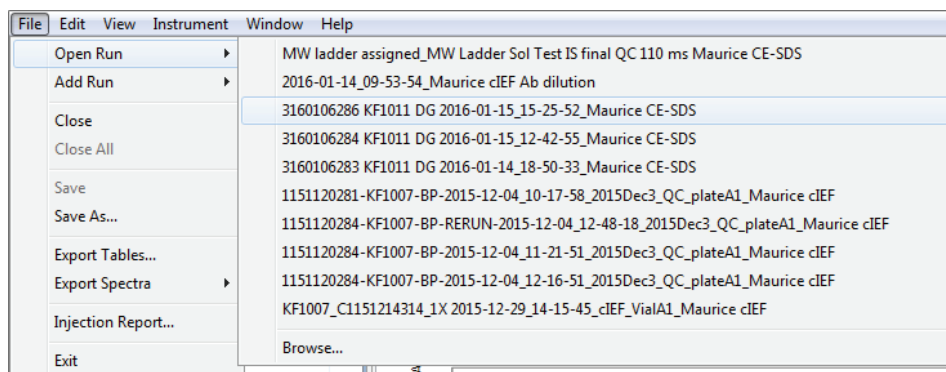
- **Copy** - Copies the information in the History pane so you can paste it into other documents.
- **Preferences** - Lets you set and save your preferences for Access Control, data export, graph colors, grouped data and Twitter settings. See Chapter 21: "Setting Your Preferences" for more information.

Opening Run Files

You can open one run file or multiple files at the same time to compare information between runs.

Opening One Run File

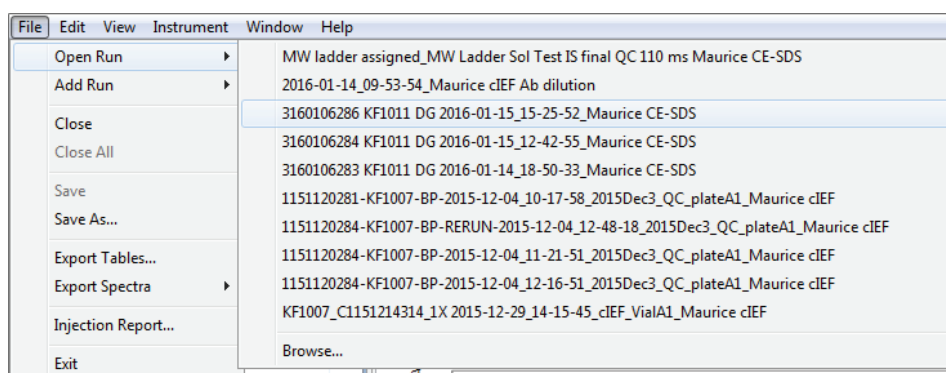
1. Select **File** in the main menu and click **Open Run**.



2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.

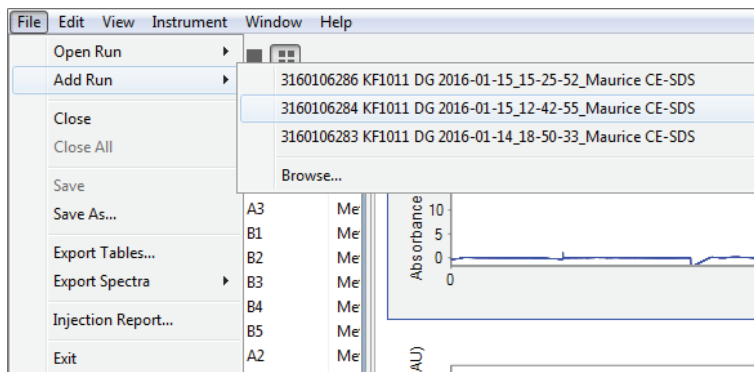
Opening Multiple Run Files

1. To open the first run file, select **File** in the main menu and click **Open Run**.



2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.

3. To open another run file, select **File** in the main menu and click **Add Run**.



4. A list of runs will display. You can only open a run that uses a similar application as the run that's already open (cIEF with MauriceFlex cIEF, MauriceFlex Fractionation only, or CE-SDS PLUS with Turbo CE-SDS), so the run files displayed are only for that application. Select one of these runs or click **Browse** to open the Runs folder and select a different file.
5. Repeat the last two steps to open additional runs.

Batch Injection Information

The Injections pane lists the system protocols (Setup and Cleanup) and injections performed during the run.

	Injection Name	Sample ID	Location	Method	Status
1	IgG System Co...	IgG System Co...	A1	Setup	Completed
2	Control Ladder...	Control Ladder	A2	Method1	Completed
3	Test Ladder_03	Test Ladder	A3	Method2	Completed
4	IS - Alpha_04	IS - Alpha	B1	Method1	Completed
5	IS - Frozen P3_...	IS - Frozen P3	B2	Method1	Completed
6	IS - T1 P3_06	IS - T1 P3	B3	Method1	Completed
7	IS - T2 P3_07	IS - T2 P3	B4	Method1	Completed
8	IS - T3 P3_08	IS - T3 P3	B5	Method1	Completed
9	Control Ladder...	Control Ladder	A2	Method2	Completed
10	Test Ladder_10	Test Ladder	A3	Method2	Completed
11	IS - Alpha_11	IS - Alpha	B1	Method1	Completed
12	IS - Frozen P3_...	IS - Frozen P3	B2	Method1	Completed
13	IS - T1 P3_13	IS - T1 P3	B3	Method1	Completed
14	IS - T2 P3_14	IS - T2 P3	B4	Method1	Completed
15	IS - T3 P3_15	IS - T3 P3	B5	Method1	Completed
16	Control Ladder...	Control Ladder	A2	Method2	Completed
17	Test Ladder_17	Test Ladder	A3	Method2	Completed
18	IS - Alpha_18	IS - Alpha	B1	Method1	Completed
19	IS - Frozen P3_...	IS - Frozen P3	B2	Method1	Completed
20	IS - T1 P3_20	IS - T1 P3	B3	Method1	Completed
21	IS - T2 P3_21	IS - T2 P3	B4	Method1	Completed
22	IS - T3 P3_22	IS - T3 P3	B5	Method1	Completed
23	Control Ladder...	Control Ladder	A2	Method2	Completed
24	Test Ladder_24	Test Ladder	A3	Method2	Completed
25	IS - Alpha_25	IS - Alpha	B1	Method1	Completed
26	IS - Frozen P3_...	IS - Frozen P3	B2	Method1	Completed
27	IS - T1 P3_27	IS - T1 P3	B3	Method1	Completed

CE-SDS

	Injection Name	Sample ID	Location	Method	Status
				Setup	Completed
				Calibration	Completed
1	System Suitabi...	System Suitabi...	A1	System S...	Completed
2	mAb 11 Blank...	mAb 11 Blank	A2	mAb Me...	Completed
3	mAb 11 Ref. St...	mAb 11 Ref. St...	A3	mAb Me...	Completed
4	mAb 11 Prep 2...	mAb 11 Prep 2...	A4	mAb Me...	Completed
5	mAb 11 Prep 2...	mAb 11 Prep 2...	A4	mAb Me...	Completed
6	mAb 11 Prep 2...	mAb 11 Prep 2...	A4	mAb Me...	Completed
7	mAb 11 Ref. St...	mAb 11 Ref. St...	A3	mAb Me...	Completed
8	mAb 11 Blank...	mAb 11 Blank	A2	Cleanup	Completed

Standard cIEF

	Injection Name	Sample ID	Location	Method	Status
1	sys suit_01	sys suit	A3	Setup	Completed
2	sys suit_02	sys suit	A3	Calibration	Completed
				system suit_1	Completed
				Cleanup	Completed
				system suit_2	Completed
				Cleanup	Completed

MauriceFlex cIEF

	Injection Name	Sample ID	Location	Method	Status
1	Sample 01_01	Sample 01	A1	Setup	Completed
				Calibration	Completed
				NIST	Completed
				Fractionati...	Completed
				Cleanup	Completed

MauriceFlex Fractionation

- Each injection includes information on:
 - Injection name
 - Sample ID
 - Sample location
 - Method
 - Run status
 - Time of injection
 - Sample temperature
- Clicking on an injection displays its data in the Focus Series (cIEF, MauriceFlex cIEF and MauriceFlex Fractionation) or Separation (CE-SDS PLUS and Turbo CE-SDS) and IV Plot panes. Hovering over a method name displays the method parameters:

	Sample ID	Location	Method	Status
			Setup	Completed
1	IgG System Co...	A1	Method1	Completed
2	Control Ladder	A2	Method2	
3	Test Ladder	A3	Method2	

Method1
 Sample Load: 20 sec 4600 Volts
 Separation: 0.1 min 1150 Volts, 0.1 min 3450 Volts, 25.0 min 5750 Volts

Injections				
	Sample ID	Location	Method	Status
			Setup	Completed
			Calibration	Completed
1	mAb11 Sample 1	A1	Method1	Completed
2	mAb11 Sample 2	A2	Method1 Separation: 1.0 min 1500 Volts, 6.0 min 3000 Volts Detection: 6 Exposures Sample Load: 90 seconds pI Markers: 4.05, 9.99	
3	mAb11 Sample 3	A3		
4	mAb11 Sample 4	A4		
5	mAb11 Sample 5	A5		

For runs in progress, the Status column displays:

- **Running** for Setup, Conditioning (CE-SDS only) and Cleanup protocols that are in progress
- **Loading** or **Separating** for injections in progress. Once the separation starts, a status bar displays next to the injection so you know when the separation will be done. Hovering your mouse over the progress bar tells you the time left for the injection.
- **Completed** for Setup, Conditioning and Cleanup protocols and injections that are done.

	Sample ID	Locat...	Method	Status	
			Setup	Completed	
1	MW Ladder	A1	MW Markers	Separating	31 mins
2	IgG-R	A8	Reduced IgG		5 mins remaining

For MauriceFlex Fractionation runs in progress, the fractionation step Status column additionally displays:

- **Mobilizing** for when Mobilization solution is being applied to the cartridge.
- **Refocusing** for when the optional refocusing step is being applied
- **Eluting (Fractionation number)** for when protein is mobilizing off of the capillary and being collected into fraction wells on the 96-well plate. The fraction being collected in the series will be noted in parenthesis.


Injection Flags

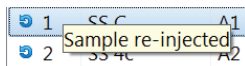
If Compass for iCE detects a potential injection issue, a flag icon will display next to the injection row in the Injections pane.



Past cartridge injection limit notification - This means the injection is over the guaranteed number of injections for the cartridge. Roll your mouse over the icon to display details.

Injections		
	Sample ID	Locat...
3	0.5% Tween	A3
4	TBST	A4
5	SB	A1
6	Past cartridge injection limit	
7	0.5% Tween	A3

-  **Reinjection notification** - This means an abnormal current profile is detected, so the separation was stopped and the sample was reinjected. The second injection always runs to completion even if the current drops again. Roll your mouse over the icon to display details.



Run Status Information

The Status pane shows info specific to each run file:

- Run file name and path (directory location)
- Batch name and type
- Instrument and serial number
- Type of sample tray used
- Run start/complete date and time
- Type of cartridge
- Cartridge serial number
- Cartridge batch injection limit, injections/batches remaining and expiration date
- Recorded run errors or warning. See “Viewing Run Errors or Warnings” on page 424 for more information.

Status	History
run	2021-12-15_11-35-24_Maurice CE-SDS TURBO_96_Inj
path	C:\Users\User\Documents\Compass for iCE\Runs
batch	Maurice CE-SDS TURBO_8_Inj
batch type	CE-SDS
instrument	Maurice : Maurice kf1147 - kf1147
samples	96-well plate
started	Wed 11:35 AM Dec 15, 2021 PST
completed	Thu 1:39 AM Dec 16, 2021 PST
cartridge	Turbo CE-SDS™
serial number	8211213001
injections per batch	96
injections remaining	(4 guaranteed)
batches remaining	20
expires	Dec 2022

run	2021-07-28_14-35-03_2021_0503_Maurice CE-SDS PLUS
path	C:\Users\User\Documents\Compass for iCE\Runs
batch	2021_0503_Maurice CE-SDS PLUS
batch type	CE-SDS
instrument	Maurice : Maurice kf1018 - kf1018
samples	48 vials
started	Wed 2:33 PM Jul 28, 2021 PDT
completed	Wed 5:13 PM Jul 28, 2021 PDT
cartridge	CE-SDS PLUS
serial number	6210115849
injections per batch	48
injections remaining	358 (0 guaranteed)
batches remaining	19
expires	Jan 2022

run	2022-12-20_13-47-10_Maurice FLEX cIEF batch 20221218
path	
batch	Maurice FLEX cIEF batch 20221218
batch type	cIEF Fractionation
instrument	MauriceFlex : Maurice Flex xf1109 - xf1109
samples	96-well plate
started	Tue 1:47 PM Dec 20, 2022 PST
cartridge	MauriceFlex
serial number	9900000100
injections per batch	15
injections remaining	10
batches remaining	5
expires	Jan 2025

run	2022-11-08_11-56-11_Maurice Flex Fractions
path	C:\Users\User\Documents\Compass for iCE\Runs
batch	Maurice Flex Fractions
batch type	cIEF Fractionation
instrument	MauriceFlex : Maurice Flex kf1884 - kf1884
samples	96-well plate
started	Tue 11:56 AM Nov 8, 2022 PST
completed	Tue 1:55 PM Nov 8, 2022 PST
cartridge	MauriceFlex
serial number	5221012106
injections per batch	1
injections remaining	11
batches remaining	10
expires	Oct 2023

run	KF0002_cIEF1160607664_2016-08-09_12-01-50_mAB high low
path	C:\Users\User\Documents\Compass for iCE\Runs
batch	mAB high low_OBM Validation
batch type	cIEF
instrument	Maurice : Maurice kf0002 - kf0002
samples	96-well plate
started	Tue 2:02 PM Aug 9, 2016 CDT
completed	Wed 1:48 AM Aug 10, 2016 CDT
cartridge	cIEF
serial number	1160607664
injections per batch	100
injections remaining	60 (60 guaranteed)
batches remaining	9
expires	Jun 2017

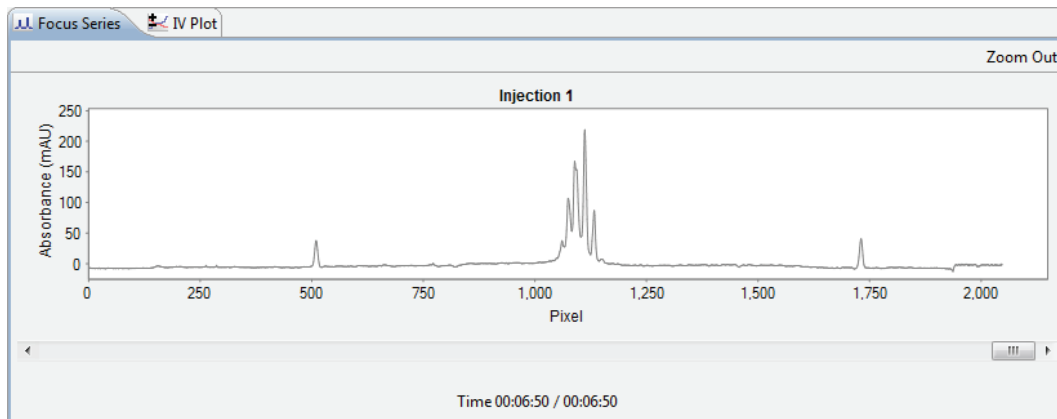
- **To go to the run file directory location** - Double click the path hyperlink, or right-click and select **Open Directory**.
- **To copy the path** - Right-click on the path hyperlink and click **Copy Path**. The path can then be copied into documents. The path can also be copied into the Windows Explorer address bar to launch Compass for iCE and open the run file automatically.

Viewing the Injection Focus Series (cIEF, MauriceFlex cIEF, MauriceFlex Fractionation Only)

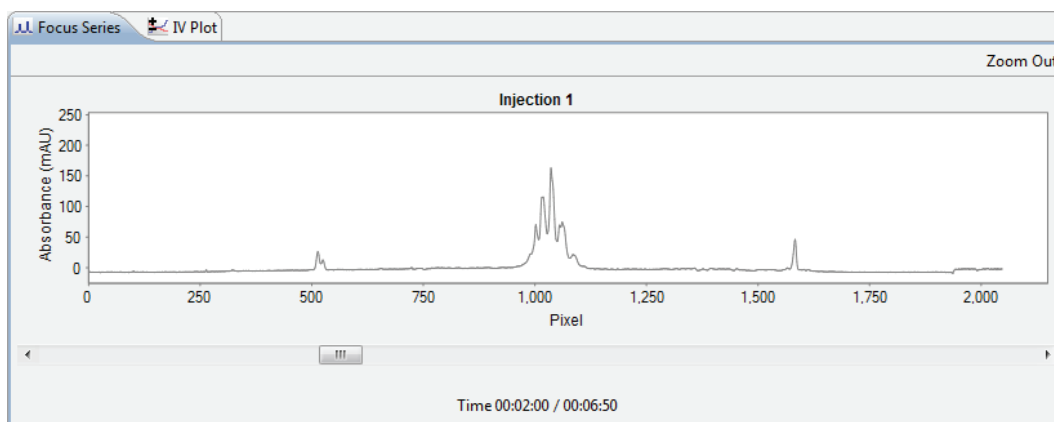
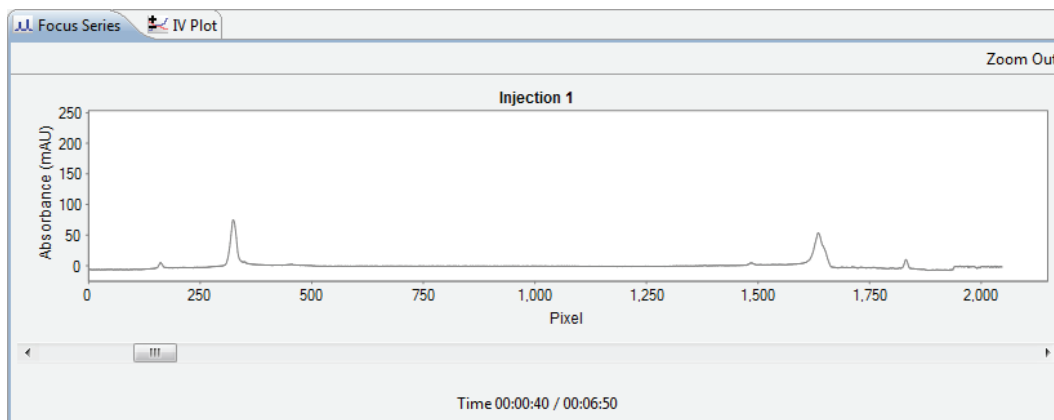
You can view your proteins focusing along the pH gradient in the capillary for each injection in the Focus Series pane.

NOTE: The Focus Series y-axis displays in absorbance for a standard cIEF assay and fluorescence for a MauriceFlex cIEF and MauriceFlex Fractionation run.

1. Select an injection in the Injections pane.
2. Click the Focus Series pane. It'll display the final focusing plot:



3. To view the focusing as it happened, drag the slider bar under the plot to the left or right. To view it frame by frame, click the left/right arrows.



- **To zoom in on an area of the plot** - Hold the mouse button down and draw a box around the area with the mouse.
- **To zoom out** - Click **Zoom Out** in the upper right corner of the pane.

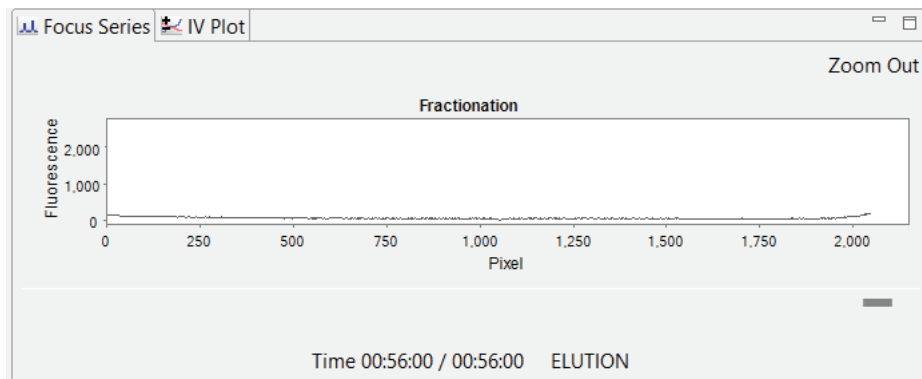
NOTE: Focus Series data for a run in progress won't be available until the injection is executing. Once it starts, the plot displays data up to the current point in time.

Viewing the Fractionation Focus Series (MauriceFlex Fractionation Only)

You can view your separated protein peaks mobilizing out of the capillary in the Focus Series pane.

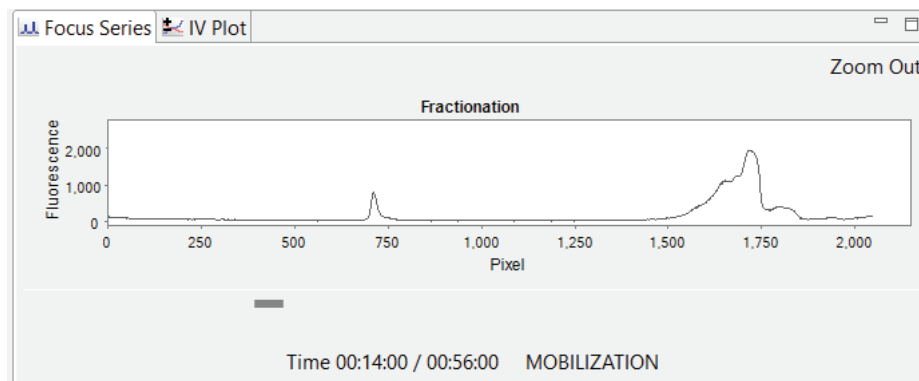
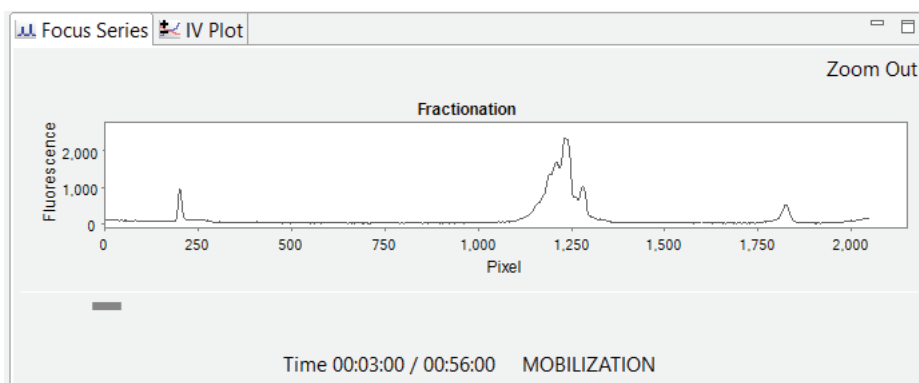
1. Select the Fractionation step in the Method column in the Injections pane.

- Click the Focus Series pane. It will display the final focusing plot:



NOTE: The fractionation status is updated next to Focus Series timestamp.

- To view separated protein peak mobilization as it happened, drag the slider bar under the plot to the left or right. To view it frame by frame, click the left/right arrows.



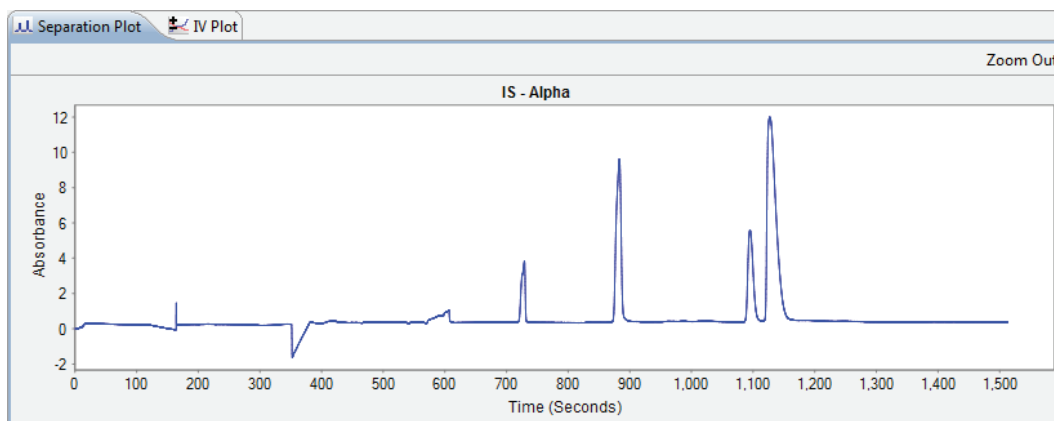
- **To zoom in on an area of the plot** - Hold the mouse button down and draw a box around the area with the mouse.
- **To zoom out** - Click **Zoom Out** in the upper right corner of the pane.

NOTE: Focus Series data for the fractionation step progress won't be available until the IEF separation is complete and mobilization/fractionation is executing. Once it starts, the plot displays data up to the current point in time.

Viewing the Separation Plot (CE-SDS PLUS and Turbo CE-SDS Only)

You can view your protein separation in the capillary for each injection in the Separation Plot pane.

1. Select an injection in the Injections pane.
2. Click the Separation Plot pane. It'll display a plot of the raw separation data.



- **To zoom in on an area of the plot** - Hold the mouse button down and draw a box around the area with the mouse.
- **To zoom out** - Click **Zoom Out** in the upper right corner of the pane.

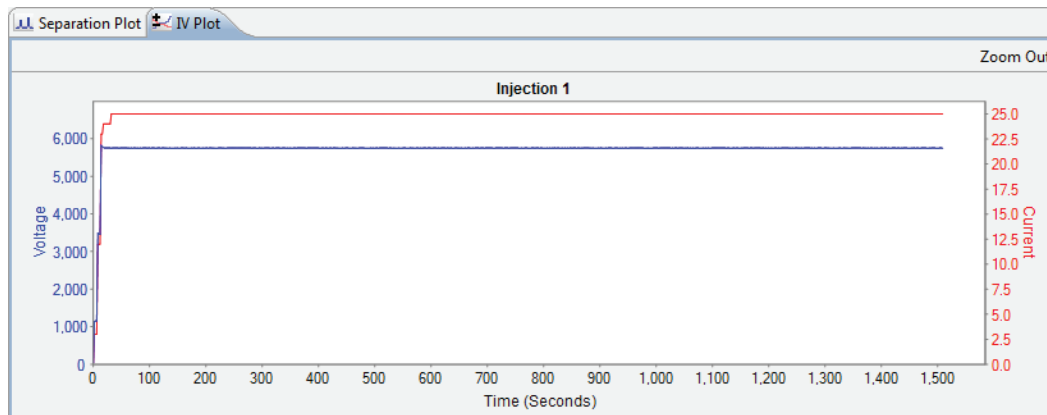
NOTE: Separation data for a run in progress won't be available until the injection is executing. Once it starts, the plot displays data up to the current point in time.

Current and Voltage Plots

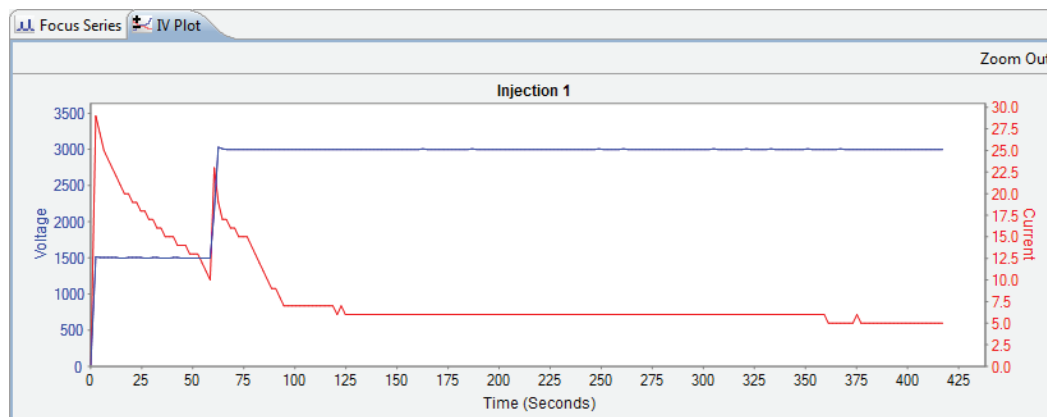
To view plots of the total current and voltage measured during an injection:

1. Select an injection in the Injections pane.

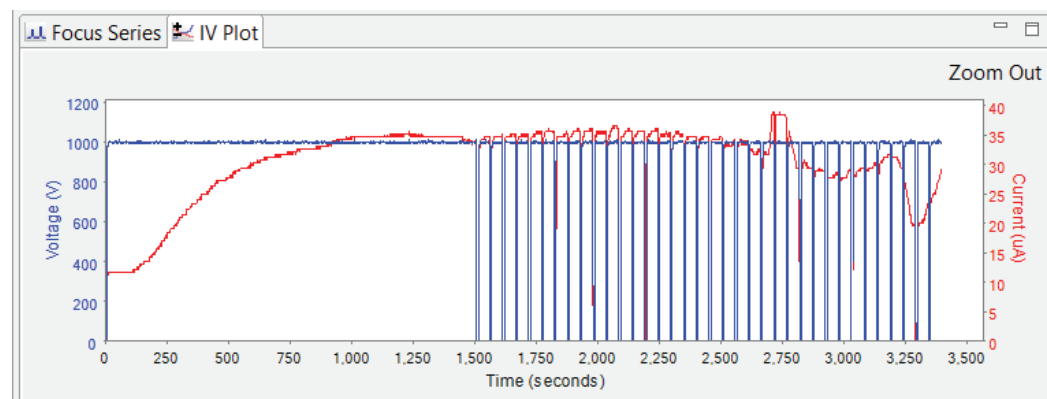
2. Click the IV Plot pane.



CE-SDS Application IV Plot



cIEF Application IV Plot



Fractionation Application Mobilization/Fractionation IV Plot

The blue Y-axis and plot shows the run voltage in volts (V), and the red Y-axis and plot shows the run current in microamps (μA). The X-axis displays time in seconds.

- **To zoom in on an area of the plot** - Hold the mouse button down and draw a box around the area with the mouse.
- **To zoom out** - Click **Zoom Out** in the upper right corner of the pane.

NOTE: IV Plots for a run in progress won't be available until the injection is executing. Once it starts, the plot displays in real time.

Run History

The History pane shows the run file event history, starting with the date and time the run was started through the most current analysis event. Clicking on a row in the table displays the full event details in the box under the table.

Date	User Name	Message	Comment
12/06/2015 3:13 PM		Started run: 2015-12-06_15-13-01_Maurice cIEF_...	
12/08/2015 10:43 AM		Saved analysis changes	
12/08/2015 3:21 PM		Saved analysis changes	
12/09/2015 12:31 PM		Saved analysis changes	
12/14/2015 2:23 PM		Saved analysis changes	
01/05/2016 4:39 PM		Saved analysis changes	

Time	12/06/2015 3:13 PM	User	
Message	Started run: 2015-12-06_15-13-01_Maurice cIEF_Mab11_TechRep Assay: Maurice cIEF.batch		
Comment			

[illegible]

- **Date:** Date and time of the run event.
- **User Name:** User that initiated the event. User names only display if you're using the Access Control feature to help satisfy the 21CFR Part 11 data security requirements.

- **Message:** Description of the event that took place.
- **Comment:** Comments entered by the user when the batch was saved.

Viewing Multiple Events

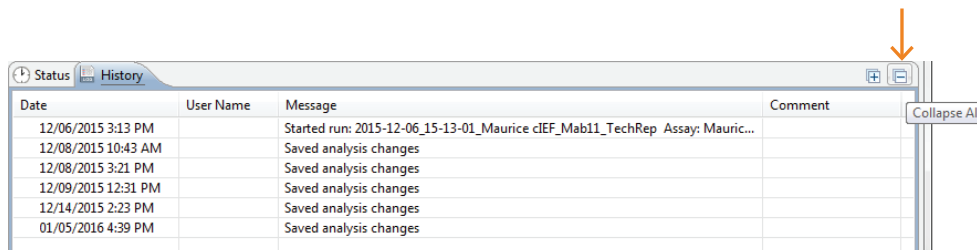
Items in the table with multiple analysis events have an arrow next to the date and time. You can view or hide these details by toggling the arrow:

Status History			
Date	User Name	Message	Comment
12/06/2015 3:13 PM		Started run: 2015-12-06_15-13-01_Maurice cIEF_Mab11_TechRep Assay: Mauric...	
12/08/2015 10:43 AM		Saved analysis changes	
		Changed 1 from "Exposure 1 1 seconds" to "Exposure 3 20 seconds"	
		Changed select: type from absorption to fluorescence	
		Changed Peak Fit Analysis Settings Peak Fit: Range Maximum from 11.0 to 10.1	
		Changed Peak Fit Analysis Settings Peak Fit: Range Minimum from 2.0 to 4.0	
12/08/2015 3:21 PM		Saved analysis changes	
12/09/2015 12:31 PM		Saved analysis changes	
12/14/2015 2:23 PM		Saved analysis changes	
01/05/2016 4:39 PM		Saved analysis changes	
Time 12/08/2015 10:43 AM User			
Message Saved analysis changes			
Comment			

- To view details for all items with multiple analysis events in the run, click the **Expand All** button.

Status History			
Date	User Name	Message	Comment
12/06/2015 3:13 PM		Started run: 2015-12-06_15-13-01_Maurice cIEF_Mab11_TechRep Assay: Mauric...	
12/08/2015 10:43 AM		Saved analysis changes	
		Changed 1 from "Exposure 1 1 seconds" to "Exposure 3 20 seconds"	
		Changed select: type from absorption to fluorescence	
		Changed Peak Fit Analysis Settings Peak Fit: Range Maximum from 11.0 to 10.1	
		Changed Peak Fit Analysis Settings Peak Fit: Range Minimum from 2.0 to 4.0	
12/08/2015 3:21 PM		Saved analysis changes	
		Added Peak Names Apply Settings "apply Peak Names 1 to all"	
		Added Peak Names Group Peak Names 1	
		Control Area: 10000.0	
		Control Reference Capillary: mAb11 Sample 1	
		Protein name: Peak1 pl: 6 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak2 pl: 6 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak3 pl: 6 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak4 pl: 6 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak5 pl: 6 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak6 pl: 6 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak7 pl: 7 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak8 pl: 7 Color: 32512 Range: 0.1 Control: false Show: true	
		Protein name: Peak9 pl: 7 Color: 32512 Range: 0.1 Control: false Show: true	
		Changed Peak Fit Analysis Settings Peak Fit: Range Maximum from 10.1 to 7.5	
		Changed Peak Fit Analysis Settings Peak Fit: Range Minimum from 4.0 to 6.2	

- To hide all items with multiple analysis events, click the **Collapse All** button.



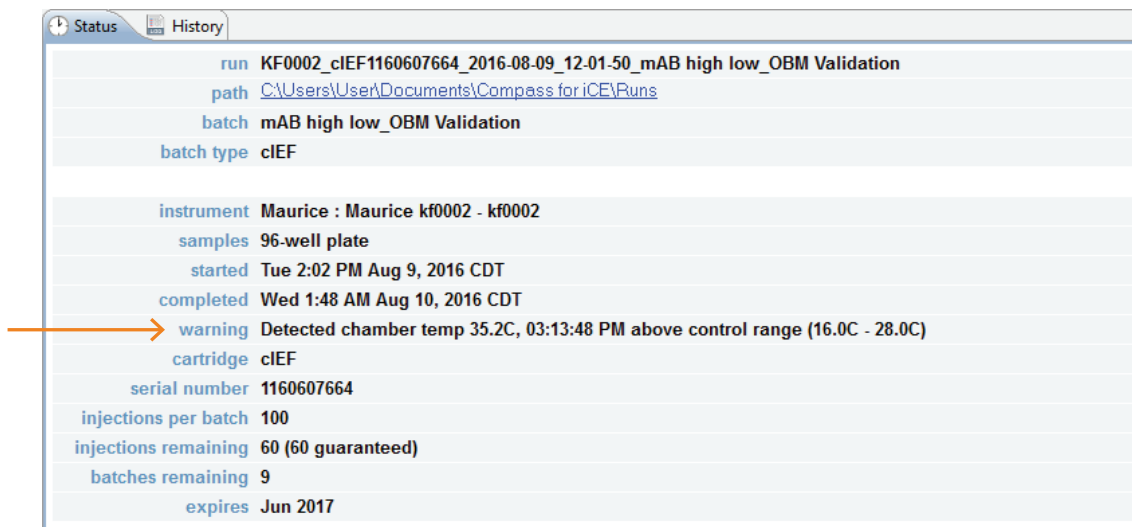
Copying History Info

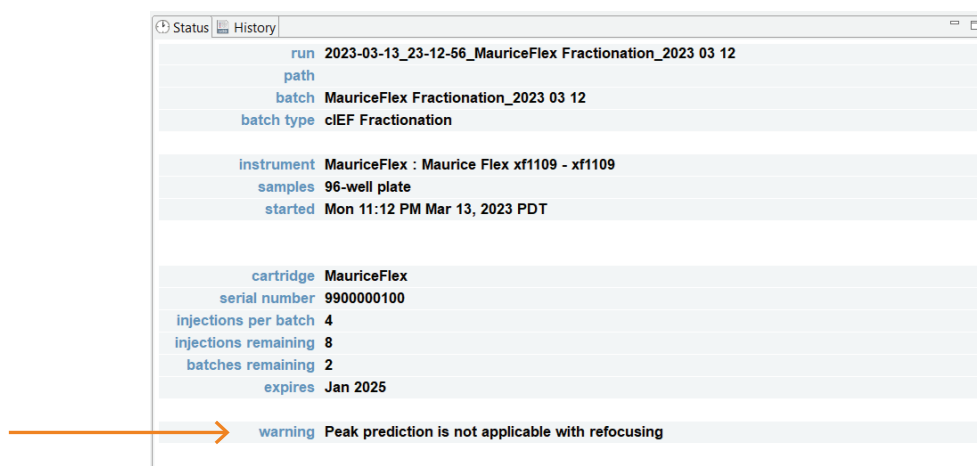
You can copy the information in the History pane to use in other documents:

1. Click the **History** pane to make sure it's active.
2. Click **Edit** in the main menu and select **Copy**.
3. Open a document and click **Paste**.

Viewing Run Errors or Warnings

If an error or warning is detected during the run it will display in the Status pane. You may need to scroll down to view errors or warnings specific to the cartridge.





Injection Reports

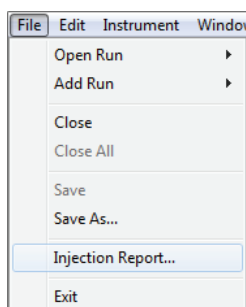
You can export PDF files of the raw and analyzed data, IV plot, peaks table, sample and system info for individual or all injections in a run file. You can also export the run history with all analysis events.

NOTES:

You can have Compass for iCE generate reports for you automatically at the end of each run. See "Setting up Automatic Injection Reports" on page 761 for more information.

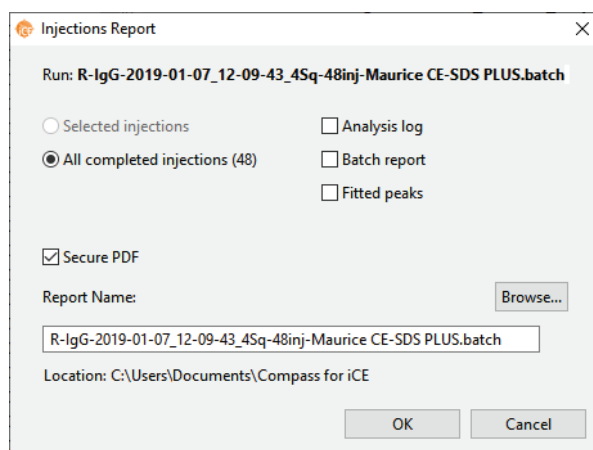
Any time a new report is run, it's saved in a separate run folder. Existing reports aren't overwritten.

1. Click **File > Open Run** and select a run file.
2. If you want reports for all injections, skip to the next step. Otherwise, select the injection in the Injection pane that you want a report for.
3. Select **File** from the main menu in either screen and click **Injection Report**.

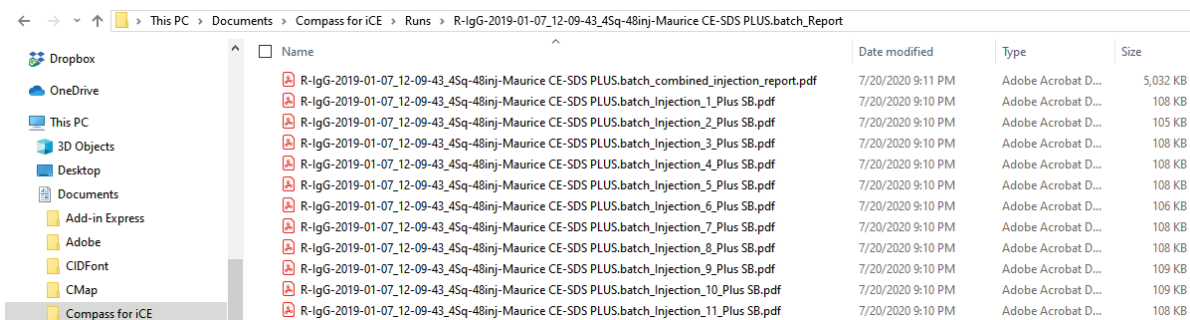


4. In the Injection Reports window:

- Choose either **Selected injections** or **All injections**.
- Select **Analysis log** if you want a run history report with all analysis events.
- Select **Batch Report** if you want to include the sample and method details for each injection in the batch.
- Select **Fitted peaks** if you want to show peak fitting in the electropherograms.
- Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don't want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.
- The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.
- Click **OK**.



- Individual injection report PDF(s) and a combined injection report PDF are exported to the Runs folder in the Compass for iCE directory. They'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Example Analysis and Injection Report: CE-SDS

Run File R-IgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS

Analysis Log

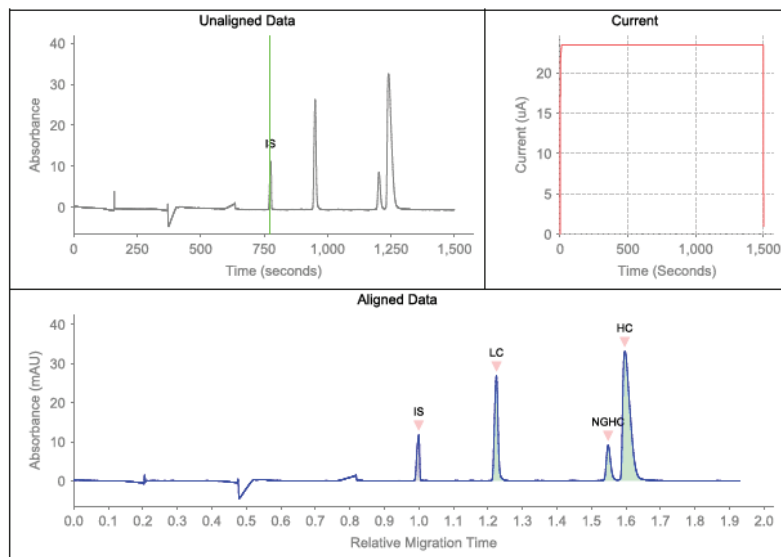
Date	User Name	Message	Comment
2019-01-10 17:25:27		Started run: R-IgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS Batch: 4Sqx12inj-RlgG-Maurice CE-SDS.batch	3180910338
2019-01-14 10:39:07		Saved analysis and methods changes from Compass for iCE v2.1.0-1219	
		Added Peak Names Apply Settings "apply IgG to all"	
		Added Peak Names Group IgG	
		Protein name: LC RMT: 1.21 Color: 32512 Range: 0.1	
		Protein name: NGHC RMT: 1.54 Color: 32512 Range: 0.05	
		Protein name: HC RMT: 1.58 Color: 32512 Range: 0.1	

Created By: Jacquelyn Sat 4:47 PM Feb 2, 2019 CST Software Version: 2.1.0, Build ID: 0130
 C:\Users\Jacquelyn\Documents\Clients\ProteinSimple\Maurice\User Guide\Rev 12 edits\Data from Andrea - no edits\R-IgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS.mbz
 Computer: DESKTOP-C7FPQ08

Page 1 of 2



Uncontrolled Injection 1: R IgG1



Peaks

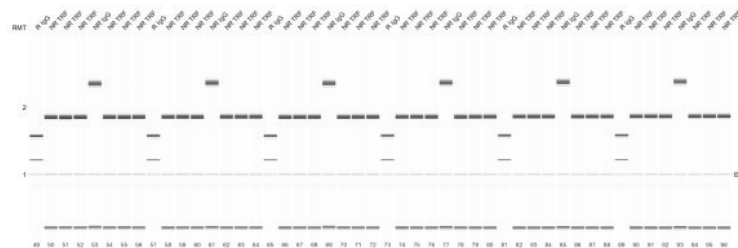
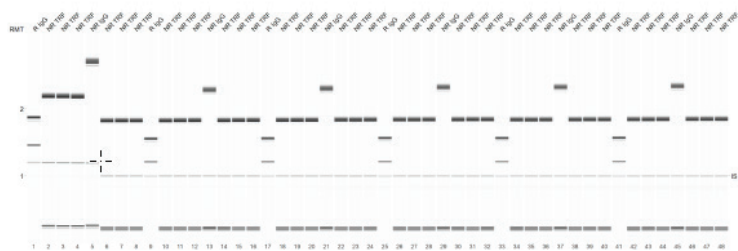
Peak	Name	Time	RMT	Height	Raw Area	Area	%Total	%Area	Width	S/N	Baseline	Resolution
1	IS	776.7	1.000	11.7	164.6	212.4			7.0	658.6	-0.6	
2	LC	951.3	1.225	26.8	498.5	524.1	30.5	30.5	8.8	1507.3	-0.6	15.02
3	NGHC	1202.7	1.549	9.1	205.6	170.8	9.9	9.9	10.2	510.5	-0.6	17.58
4	HC	1240.2	1.597	33.1	1273.4	1021.8	59.5	59.5	17.9	1859.6	-0.6	1.98

Created By: Jacquelyn Bat 4:47 PM Feb 2, 2019 CST
 Software Version: 2.1.0, Build ID: 0130
 C:\Users\Jacquelyn\Documents\Clients\ProteinSimple\Maurice\User Guide\Rev 12 edits\Data from Andrea - no edits\R-Igo-2019-01-10_16-25-12_40qr12m-R-Igo-Maurice CE-606.mtx
 Computer: DESKTOP-C7FPQGB



Page 1 of 2

Lane Image



Created By: Andrea, Mon 8:29 PM Apr 25, 2022 PDT (SECURED)
 C:\Users\Andrea\Documents\Compass for ICE\Runs\Maurice_TURBO_CE-100_Red-NonRed.mlx
 Computer: DESKTOP-IPM7G05 Software Version: Compass for ICE 3.0.0, Build ID: 0218
 Page 1 of 199



Uncontrolled Injection 1: R IgG1

Sample Information

Injection Name	R IgG1
Sample ID	SB
Location	Plate Well B2
Batch Name	4Sgx12inj-RlgG-Maurice CE-SDS
Run Started	Thu 5:25 PM Jan 10, 2019 CST
Run Completed	Fri 8:07 PM Jan 11, 2019 CST
Date Acquired	Thu 6:08 PM Jan 10, 2019 CST
Run Error	None
Reinjection	No

Injection Conditions

Guaranteed Injection	Yes
Focus Period 1	5750V for 25.0 min
Sample Load	20 sec 4600 Volts
Tray Temperature	9.1°C

Maurice Settings

Model	Maurice OBM
Instrument S/N	kf1077
Software Version	2.1.0, Build ID: 0130
Firmware Version	3.1.2019.01.04.20.16.20.5eb707f
Tray Type	96-well plate
Cartridge Type	CE-SDS
Cartridge S/N	3180910338
Cartridge Expiration	Sep 2019
Injections Remaining	152 (52 guaranteed)
Batches Remaining	24

Created By: jacquelyn Sat 4:47 PM Feb 2, 2019 CST
 C:\Users\jacquelyn\Documents\Orients\ProteinSimple\Maurice\User Guide\Rev 12 edit\Data from Andrea - no edit\RIgG-2019-01-10_16-25-12_48lg12inj-RlgG-Maurice CE-SDS.mtz
 Computer: DESKTOP-C7FPQGB

Software Version: 2.1.0, Build ID: 0130

Page 2 of 2



Example Analysis and Injection Report: cIEF

Run File DemoData_Maurice cIEF

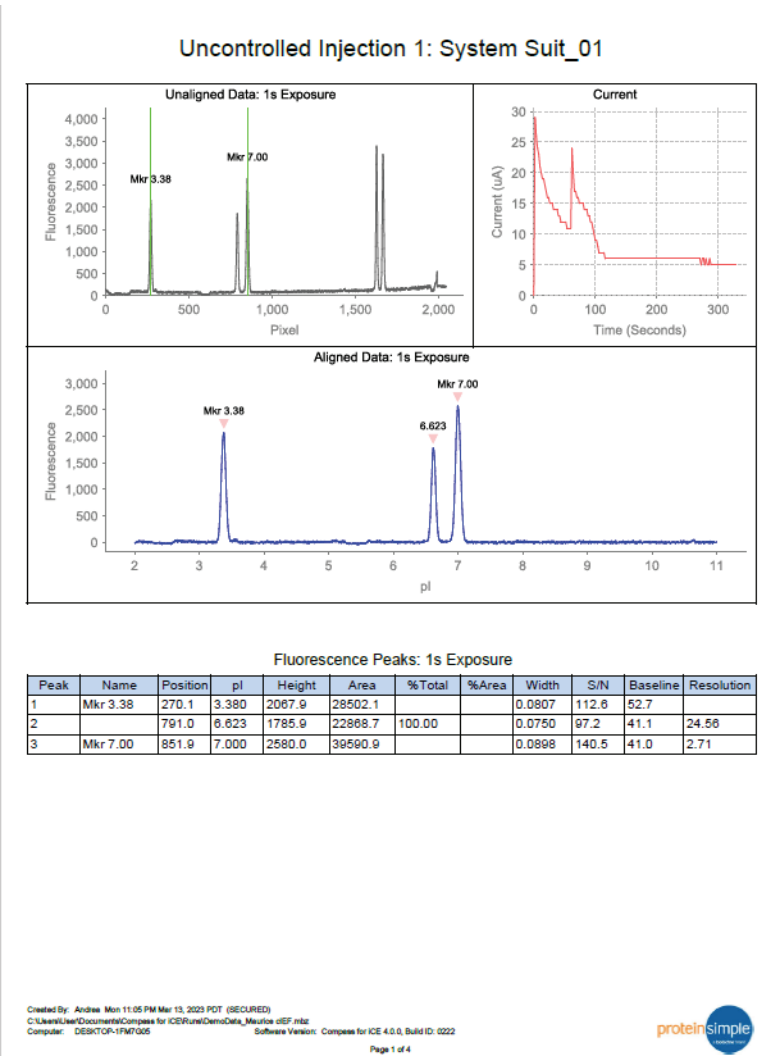
Analysis Log

Date	User Name	Message	Comment
2015-10-27 12:11:27		Started run: 2015-10-27_12-11-12_Maurice cIEF Batch: Maurice cIEF.batch	
2015-10-27 14:46:00		Saved analysis changes	
		Changed MethodName Method Exposure from "Exposure 1 Absorbance 0.005 seconds" to "Exposure 1 Fluorescence 1 seconds"	
		Changed MethodName Method Exposure from "Exposure 1 Absorbance 0.005 seconds" to "Exposure 5 Fluorescence 20 seconds"	
		Added Standards Apply Override "apply Standards 3 to Method1"	
		Added Standards 3	
		pI Marker pI: 3.38 Position: 250	
		pI Marker pI: 7 Position: 800	
2015-10-27 14:54:00		Save run file	

Created By: Andrea Mon 11:05 PM Mar 13, 2023 PDT (SECURED)
 C:\Users\User\Documents\Compass for ICE\Runs\DemoData_Maurice cIEF.mtz
 Computer: DESKTOP-IPMT092 Software Version: Compass for ICE 4.0.0, Build ID: 0222

Page 1 of 2





Uncontrolled Injection 1: System Suit_01

Sample Information

Injection Name	System Suit_01
Sample ID	System Suit
Location	Plate Well A1
Batch Name	Maurice cIEF
Run Started	Tue 12:11 PM Oct 27, 2015 PDT
Run Completed	Tue 1:17 PM Oct 27, 2015 PDT
Date Acquired	Tue 12:17 PM Oct 27, 2015 PDT
Run Error	None
Reinjection	No

Injection Conditions

Guaranteed Injection	Yes
On-board Mixing	Off
Focus Period 1	1500V for 1.0 min
Focus Period 2	3000V for 4.5 min
Sample Load Duration	90.0 Seconds
pI marker 1	3.38
pI marker 2	7.0
Tray Temperature	Not Available

Maurice Settings

Model	Maurice OBM
Instrument S/N	kf1004
Software Version	Compass for iCE 4.0.0, Build ID: 0222
Firmware Version	2.0.2015.10.20.20.35.52.681a5af
Tray Type	48 vials
Cartridge Type	
Cartridge S/N	
Cartridge Expiration	
Injections Remaining	
Batches Remaining	

Created By: Andrea Mon 11:05 PM Mar 13, 2023 PDT (SECURED)
 C:\Users\User\Documents\Compass for ICE\Run\DemoData_Maurice cIEF.mz
 Computer: DESKTOP-1FM7G05 Software Version: Compass for ICE 4.0.0, Build ID: 0222

Page 2 of 4



Example Analysis and Injection Report: MauriceFlex Fractionation

Run File 2023-03-14_14-03-04_MauriceFlex Fractionation_PS IgG_refocus

Analysis Log

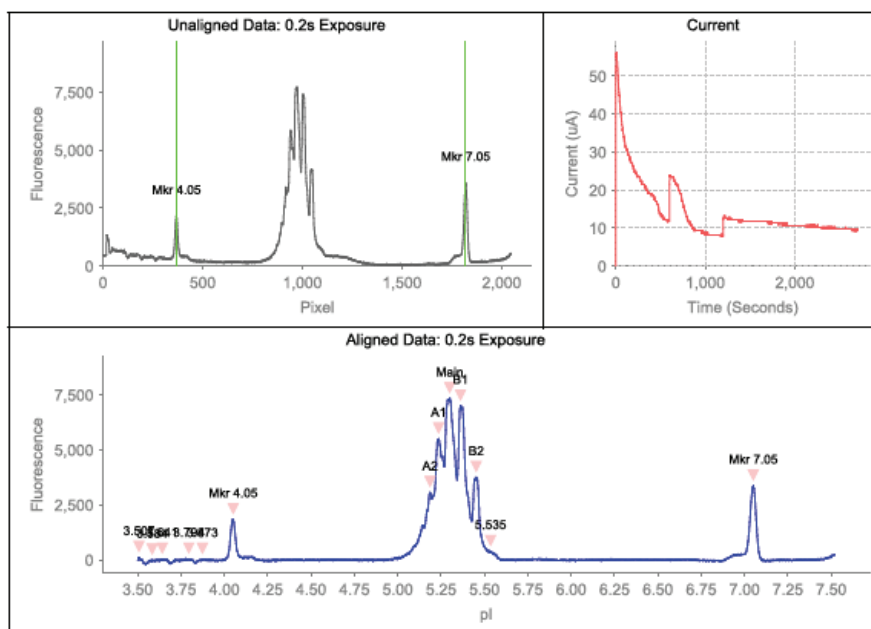
Date	User Name	Message	Comment
2023-03-14 14:08:37		Started run: 2023-03-14_14-03-04_MauriceFlex Fractionation_PS IgG_refocus Batch: MauriceFlex Fractionation(1).batch from Compass for iCE v4.0.0-0222	5230202259
2023-03-15 16:39:29		Saved analysis changes from Compass for iCE v4.0.0-0222	
		Added Peak Names Group Peak Group 1	
		Protein name: Main pl: 5.3 Color: 32512 Range: 0.05	
		Added Peak Names Apply Settings "apply Peak Group 1 to all"	
		Added Protein B1 to Peak Names Group Peak Group 1	
		name: B1	
		pl: 5.363	
		Color: 32512	
		Range: 0.05	
		Added Protein B2 to Peak Names Group Peak Group 1	
		name: B2	
		pl: 5.452	
		Color: 32512	
		Range: 0.05	
		Added Protein A1 to Peak Names Group Peak Group 1	
		name: A1	
		pl: 5.235	
		Color: 32512	
		Range: 0.05	
		Added Protein A2 to Peak Names Group Peak Group 1	
		name: A2	
		pl: 5.187	
		Color: 32512	
		Range: 0.05	

Created By: Andrea Sun 8:11 PM Mar 19, 2023 PDT (SECURED)
 C:\Users\Andrea\Desktop\2023-03-14_14-03-04_MauriceFlex Fractionation_PS IgG_refocus.mtz
 Computer: DESKTOP-1FM7005 Software Version: Compass for iCE 4.0.0, Build ID: 0222

Page 1 of 2



Uncontrolled Injection 1: Sample 01_01



Fluorescence Peaks: 0.2s Exposure

Peak	Name	Position	pI	Height	Area	%Total	%Area	Width	S/N	Baseline	Resolution
1		105.9	3.507	-13.8	0.0			0.0021	-0.5	521.5	
2		143.2	3.584	-68.8	0.0			0.0000	-2.7	474.1	43.83
3		170.7	3.641	36.0	118.3	0.02		0.0070	1.4	412.1	9.62
4		244.7	3.794	38.2	38.2	0.01		0.0000	1.5	332.9	25.92
5		283.1	3.873	-50.4	0.0			0.0041	-2.0	311.1	22.60
6	Mkr 4.05	368.6	4.050	1837.6	31693.9			0.0319	71.3	276.8	5.77
7	A2	918.4	5.187	3068.6	82235.5	11.24	11.4	0.0561	119.0	279.3	15.20
8	A1	941.6	5.235	5506.2	154682.1	21.14	21.4	0.0724	213.6	307.4	0.44
9	Main	973.5	5.300	7339.0	213113.7	29.13	29.5	0.0733	284.7	335.3	0.53
10	B1	1003.5	5.363	6999.9	187342.2	25.61	25.9	0.0526	271.6	363.9	0.58
11	B2	1046.6	5.452	3752.7	85249.6	11.65	11.8	0.0418	145.6	387.0	1.11
12		1087.0	5.535	413.5	8781.7	1.20		0.0478	16.0	396.7	1.10
13	Mkr 7.05	1819.8	7.050	3367.0	71575.3			0.0390	130.6	168.8	20.56

Uncontrolled Injection 1: Sample 01_01

Peak Predictions

Peak	Name	Predicted Wells
1	3.507	G3, G4, G5, G6, G7, G8
2	3.584	G1, G2, G3, G4, G5, G6
3	3.641	F1, G1, G2, G3, G4, G5
4	3.794	F4, F3, F2, F1, G1, G2, G3
5	3.873	F5, F4, F3, F2, F1, G1
6	Mkr 4.05	F8, F7, F6, F5, F4, F3
7	A2	E2, E3, E4, E5, E6, E7
8	A1	E1, E2, E3, E4, E5, E6
9	Main	
10	B1	E1, E2, E3, E4, E5, E6
11	B2	E1, E2, E3, E4, E5
12	5.535	E3, E4, E5, E6, E7, E8
13	Mkr 7.05	

Uncontrolled Injection 1: Sample 01_01

Sample Information

Injection Name	Sample 01_01
Sample ID	Sample 01
Location	Plate Well A3
Batch Name	MauriceFlex Fractionation(1)
Run Started	Tue 2:08 PM Mar 14, 2023 PDT
Run Completed	Tue 4:30 PM Mar 14, 2023 PDT
Date Acquired	Tue 2:22 PM Mar 14, 2023 PDT
Run Error	None
Reinjection	No

Injection Conditions

Guaranteed Injection	Yes
On-board Mixing	Off
Focus Period 1	500V for 10.0 min
Focus Period 2	1000V for 10.0 min
Focus Period 3	1500V for 25.0 min
Detection Exposure	0.2 sec
Detection Interval	5.0 min
Sample Load Duration	20.0 Seconds
pI marker 1	4.05
pI marker 2	7.05
Tray Temperature	10.0°C

Fractionation Conditions

Mobilization	1000 Volts for 45.0 min
Refocus	1500 Volts for 5.0 min
Fractions	1000 Volts for 25.0 sec
Detection Exposure	0.2 sec
Detection Interval	1.0 min

Maurice Settings

Model	MauriceFlex
Instrument S/N	mm0006
Software Version	Compass for iCE 4.0.0, Build ID: 0222
Firmware Version	4.2.2023.02.09.22.55.08.12f36498b
Tray Type	96-well plate
Cartridge Type	MauriceFlex
Cartridge S/N	5230202259

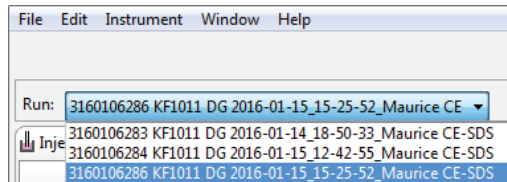
Uncontrolled Injection 1: Sample 01_01

Cartridge Expiration	Feb 2024
Injections Remaining	13 (13 guaranteed)
Batches Remaining	13
Warning	Cartridge post-run cleanup not performed after previous run
Warning	Peak prediction is not applicable with refocusing

Switching Between Open Run Files

If you've got more than one run file open, you can switch between viewing the run information in each.

1. Click the down arrow in the Run box.



2. Select the run you want to view from the drop down list.

Closing Run Files

If you've got more than one run file open, you can close just one file or all the open files at the same time.

- **To close the run file being viewed** - Select **File** from the main menu and click **Close**.
- **To close all open run files** - Select **File** from the main menu and click **Close All**.

Chapter 17:

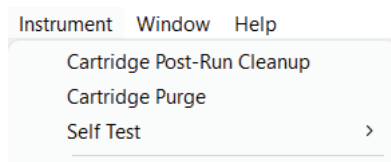
Controlling Maurice, Maurice C., Maurice S. and MauriceFlex

Chapter Overview

- Instrument Control
- Stopping a Run
- Status Modes
- Shutdown
- Instrument Software (Embedded) Updates
- Instrument Self Test
- Viewing and Changing System Properties
- Checking Cartridge Status
- Cartridge Self-Test
- Instrument Reports
- Viewing Log Files
- Sending a Log File to Technical Support

Instrument Control

The Instrument menu lets you to control Maurice, Maurice C., Maurice S. and MauriceFlex



NOTE: Instrument menu options are only active when you've got a computer with Compass for iCE software connected directly to your Maurice system.

Starting a Run

To start your run, click the **Start** button in the Batch screen. You can also start a run by selecting **Instrument** in the main menu and clicking **Start**. For more info on creating and starting batches check out

- Chapter 6: "Running cIEF Applications"
- Chapter 8: "Running MauriceFlex cIEF Applications"
- Chapter 10: "Running MauriceFlex Fractionation Applications"
- Chapter 12: "Running CE-SDS PLUS Applications"
- Chapter 14: "Running Turbo CE-SDS Applications"

Cartridge Purge

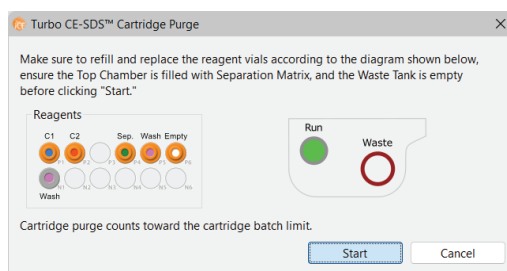
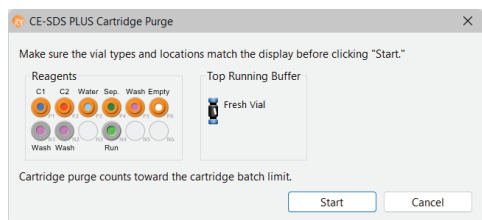
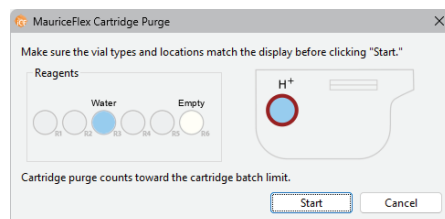
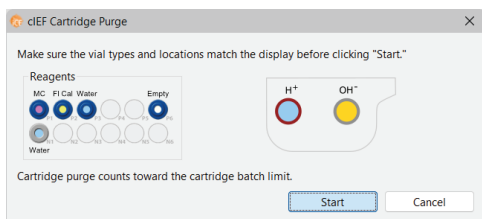
You'll want to run the Cartridge Purge any time you have to stop a run manually or if the run stops because of an error. You'll also want to run the Cartridge Purge if your CE-SDS PLUS or Turbo CE-SDS cartridge hasn't been used in more than 3 months.

Make sure Maurice's door is closed and a cartridge is installed before you run the Cartridge Purge.

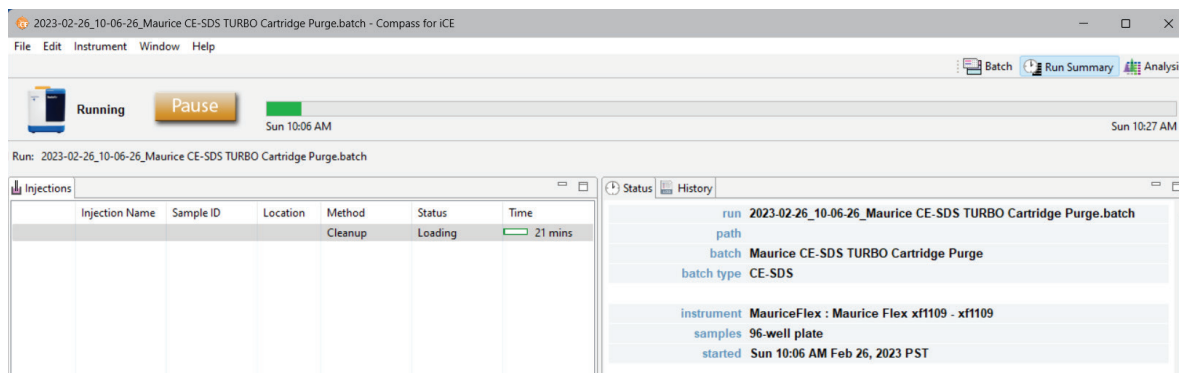
1. Compass for iCE will alert you that a Cartridge Purge is recommended. Click on the brown **Purge** button in the instrument status bar. You can also select **Instrument > Cartridge Purge** in the main menu.

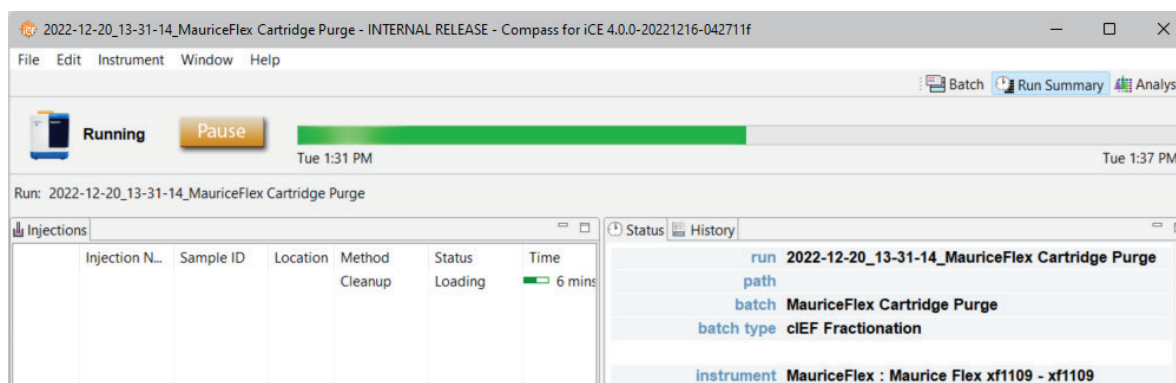
2. Click **Start**.

NOTE: The Cartridge Purge counts toward the batch limit. If there's only one batch left on the cartridge, you'll get a message that it can't be used after purging.



The Cartridge Purge for CE-SDS PLUS or Turbo CE-SDS Cartridges takes about 30 minutes, cIEF Cartridges take a little over 10 minutes and the cIEF Fractionation Cartridge takes about 6 minutes.





3. Once the purge is done, perform a Cartridge Post-Run Cleanup:

- **cIEF Cartridges:** Follow the “Post-batch Procedures” instructions on page 120.
- **cIEF Fractionation Cartridges:** Follow the “Post-batch Procedures” instructions on page 182 for a MauriceFlex cIEF batch and page 248 for a MauriceFlex Fractionation batch.
- **CE-SDS Plus Cartridges:** Follow the “Post-batch Procedures” instructions on page 316.
- **Turbo CE-SDS Cartridges:** Follow the “Post-batch Procedures” instructions on page 385.

4. Once the Cartridge Post-Run Cleanup is completed:

If you’ll be starting a new run right away, prep the cartridge.

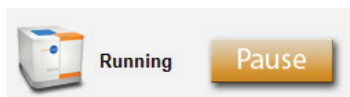
- **cIEF Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 96.
- **cIEF Fractionation Cartridges:** Follow the Step 2: Prep the Cartridge” instructions on page 163 for a MauriceFlex cIEF batch and page 229 for a MauriceFlex Fractionation batch.
- **CE-SDS Plus Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 295.
- **Turbo CE-SDS Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 364.

If you won’t be starting a new run:

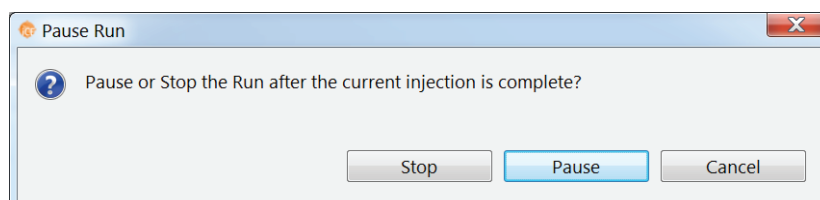
- **cIEF Cartridges:** Place the cartridge and stoppers in its protective packaging and store at room temperature.
- **cIEF Fractionation Cartridges:** Place the cartridge and stopper in its protective packaging and store at room temperature.
- **CE-SDS Plus Cartridges:** Place the cartridge in its protective packaging and store at room temperature.
- **Turbo CE-SDS Cartridges:** Place the cartridge and stoppers in its protective packaging and store at room temperature.

Stopping a Run

1. To stop a run after the current injection, you can either select **Instrument > Stop**, or click the **Pause** button:



2. If you clicked **Pause**, select **Stop** in the pop-up window.



Maurice will stop the run after the current injection is complete and then save the data for the completed instead in the .mbz file. After stopping the run, perform a Cartridge Post-Run Cleanup:

- **cIEF Cartridges:** Follow the “Post-batch Procedures” instructions on page 120.
- **cIEF Fractionation Cartridges:** Follow the “Post-batch Procedures” instructions on page 182 for a MauriceFlex cIEF batch and page 248 for a MauriceFlex Fractionation batch.
- **CE-SDS Plus Cartridges:** Follow the “Post-batch Procedures” instructions on page 316.
- **Turbo CE-SDS Cartridges:** Follow the “Post-batch Procedures” instructions on page 385.

3. Once the post-run cleanup is done:

If you'll be starting a new run right away, prep the cartridge:

- **cIEF Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 96.
- **cIEF Fractionation Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 163 for a MauriceFlex cIEF batch and page 229 for a MauriceFlex Fractionation batch.
- **CE-SDS Plus Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 295.
- **Turbo CE-SDS Cartridges:** Follow the “Step 2: Prep the Cartridge” instructions on page 364.

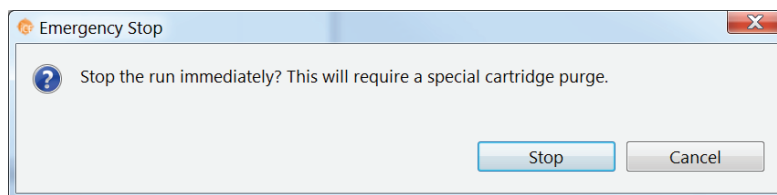
If you won't be starting a new run:

- **cIEF Cartridges:** Place the cartridge and stoppers in its protective packaging and store at room temperature.
- **cIEF Fractionation Cartridges:** Place the cartridge and stopper in its protective packaging and store at room temperature.

- **CE-SDS Plus Cartridges:** Place the cartridge in its protective packaging and store at room temperature.
- **Turbo CE-SDS Cartridges:** Place the cartridge and stoppers in its protective packaging and store at room temperature.

Emergency Stop

1. To stop a run immediately, select **Instrument > Emergency Stop**.
2. Click **Stop** in the pop up window:



Maurice will stop once it's safe to avoid damaging the cartridge. Data for completed injections will be saved in the .mbz file.

3. Perform a Cartridge Purge. Click the brown Purge button in the instrument status bar or select **Instrument > Cartridge Purge**.
4. When the cartridge purge is done, perform a Cartridge Post-Run Cleanup:
 - **cIEF Cartridges:** Follow the "Post-batch Procedures" instructions on page 120.
 - **cIEF Fractionation Cartridges:** Follow the "Post-batch Procedures" instructions on page 182 for a MauriceFlex cIEF batch and page 248 for a MauriceFlex Fractionation batch.
 - **CE-SDS Plus Cartridges:** Follow the "Post-batch Procedures" instructions on page 316.
 - **Turbo CE-SDS Cartridges:** Follow the "Post-batch Procedures" instructions on page 385.
5. Once the Cartridge Post-Run Cleanup is completed:
If you'll be starting a new run right away, prep the cartridge:
 - **cIEF Cartridges:** Follow the "Step 2: Prep the Cartridge" instructions on page 96.
 - **cIEF Fractionation Cartridges:** Follow the Step 2: Prep the Cartridge" instructions on page 163 for a MauriceFlex cIEF batch and page 229 for a MauriceFlex Fractionation batch.
 - **CE-SDS Plus Cartridges:** Follow the "Step 2: Prep the Cartridge" instructions on page 295.
 - **Turbo CE-SDS Cartridges:** Follow the "Step 2: Prep the Cartridge" instructions on page 364.

If you won't be starting a new run:

- **cIEF Cartridges:** Place the cartridge and stoppers in its protective packaging and store at room temperature.
- **cIEF Fractionation Cartridges:** Place the cartridge and stopper in its protective packaging and store at room temperature.
- **CE-SDS Plus Cartridges:** Place the cartridge in its protective packaging and store at room temperature.
- **Turbo CE-SDS Cartridges:** Place the cartridge and stoppers in its protective packaging and store at room temperature.

Status Modes

The instrument status bar displays status, buttons and progress bars depending on what Maurice, Maurice C. or Maurice S. are doing.

- **Ready/Start button** - The instrument is ready and a batch is loaded. Click **Start** to begin a run.
- **Not Ready/Reset button** - The instrument isn't ready and needs to reinitialize. Click **Reset** to start the initialization protocol.
- **Running/Pause button** - The instrument is running. The run name, time it started and when it will be done show in the run progress bar. Click **Pause** to pause or stop the run.
- **Cleaning/Stop button** - The instrument is running a cleaning protocol. The time the cleaning protocol started and when it will be done show in the run progress bar.
- **Error/Reset button** - There's an error. Go to the **Status** pane in the **Run Summary** screen to view details. When you've fixed the source of the error, click **Reset**.

Shutdown

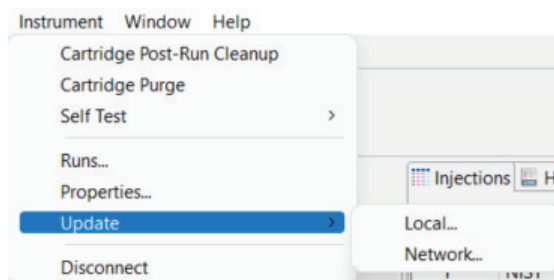
Close Compass for iCE software. Maurice can stay on unless he won't be used for an extended period.

Instrument Software (Embedded) Updates

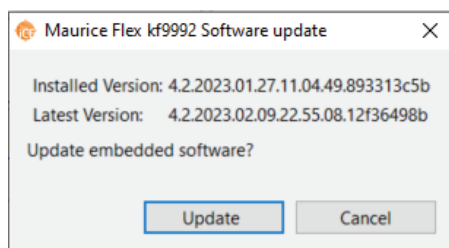
To check for embedded software updates:

If you're on the network:

1. Select **Instrument > Update**, then select **Network**.

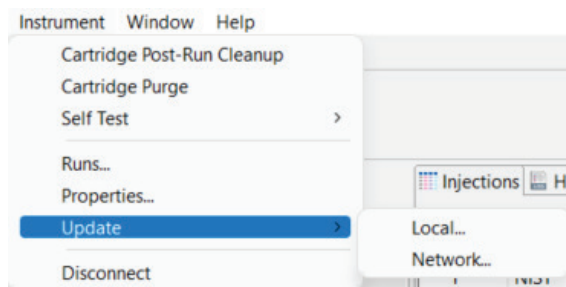


2. The following screen displays. Click **Update**.



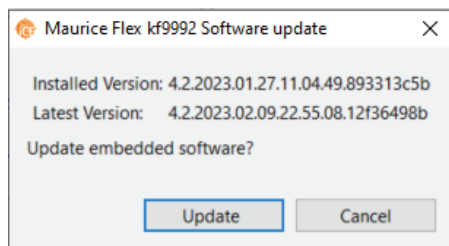
If you're not on the network:

1. Call ProteinSimple Technical Support or your FAS for assistance on getting the latest update.
2. Copy the new embedded software file onto Maurice's computer.
3. Select **Instrument > Update**, then select **Local**.



4. Browse to the location of the embedded software file, select it and click **OK**.

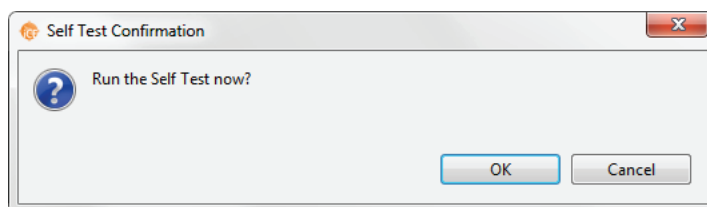
5. The following screen displays. Click **Update**.



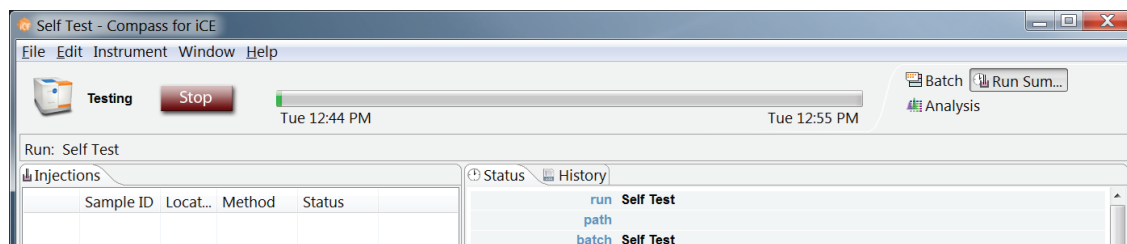
Instrument Self Test

Maurice, Maurice C., Maurice S. and MauriceFlex can run a series of self tests for you to make sure they're operating properly.

1. To start the test, select **Instrument > Self Test > Instrument**.
2. The following screen displays. Click **OK**.



The test takes approximately 11 minutes.



NOTES:

We recommend running the self test before you start a run.

An alert may appear if a cartridge is installed. Remove the cartridge to start the Instrument Self-Test.

Results will automatically display once the test's done. You can also view current or past self test results anytime by selecting **Instrument > Properties** and clicking **Test Log**. See "Self Test Logs" on page 470 for more info.

Viewing and Changing System Properties

Selecting **Instrument > Properties** displays your system properties. They include:

- System Name
- System Location
- Instrument Type
- Serial number
- Instrument software version (firmware)
- Network name and address
- Date and time of the instrument clock
- Adapter block currently in use
- Number of hours on the Deuterium (UV) lamp
- Current sample chiller temperature

Maurice kf0010 - Properties

Name:

Location:

Type: Maurice OBM Network Name: kf0010.local.
 Serial Number: kf0010 Network Address: 10.1.4.118
 Instrument Software: 3.1.2018.11.07.23.43.14.682d616
 Instrument Date and Time: 2018-11-12 10:52:46 -08:00

Adapter Block: plate Sample Chiller: 19.4°C
 D2 Lamp Run Time: 246 hours

Cartridge

Type: CE-SDS PLUS Injections per Batch: 48
 Expires: Nov 2019 Injections Remaining: 452 (52 guaranteed)
 Serial Number: 5181106034 Batches Remaining: 24

Batch	Date	Injections
1	9 Nov 2018	48
2	9 Nov 2018	cleanup

- **To change the system name or location:** Click in the name or location boxes and enter your new info, then click OK.

- **To sync the instrument clock with the computer:** Click **Set to PC time** and then restart Compass for iCE and your Maurice system.

Maurice Flex xf1109 - Properties

Name: Maurice Flex xf1109
Location: 192.168.56.01

Type: MauriceFlex Network Name: 192.168.56.103
Serial Number: xf1109 Network Address: 192.168.56.103
Instrument Software: 4.2.2022.11.03.21.24.04.b15d551f6
Instrument Date and Time: 2023-02-10 21:20:06 -08:00 **Set to PC time**

Restart Compass and MauriceFlex instrument after clicking on "Set to PC time" button.

Adapter Block: vials Sample Chiller: 10.0°C
D2 Lamp Run Time: 85 hours

Cartridge
Type: CE-SDS PLUS Injections per Batch: 48
Expires: Jan 2025 Injections Remaining: 439 (39 guarantee)
Serial Number: 0000000001 Batches Remaining: 20

Batch	Date	Injections
1	1 Sep 2018	48
2	4 Sep 2018	cleanup
3	4 Sep 2018	12
4	5 Sep 2018	cleanup
5	1 Oct 2018	1

Logs Report **OK** Cancel

Checking Cartridge Status

If you've got a cartridge installed in the system, you can see its serial number, the injections and batches it still has available, and a history of batches and injections its run to date. To view this info, select **Instrument > Properties**.

Maurice kf9991 - Properties

Name: Maurice kf9991
Location: 192.168.56.101

Type: Maurice OBM Network Name: 192.168.56.101
Serial Number: kf9991 Network Address: 192.168.56.101
Instrument Software: 4.1.2022.02.17.22.34.36.330f0b19e
Instrument Date and Time: 2022-02-18 16:08:15 -08:00 **Set to PC time**

Adapter Block: vials Sample Chiller: 10.0°C
D2 Lamp Run Time: 162 hours

Cartridge
Type: Turbo CE-SDS™ Injections per Batch: 96
Expires: Dec 2022 Injections Remaining: (50 guaranteed)
Serial Number: 8211213001 Batches Remaining: 2

Batch	Date	Injections
1	22 Dec 2021	cleanup
2	13 Jan 2022	0
3	13 Jan 2022	error
4	14 Jan 2022	0
5	14 Jan 2022	1

Cartridge post-run cleanup recommended or cartridge warranty may be voided

Logs Report **OK** Cancel

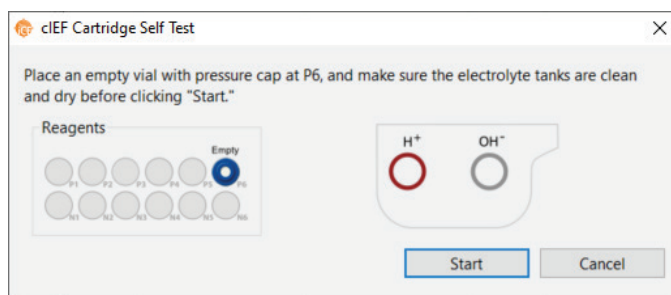
- cIEF, CE-SDS PLUS, and cIEF Fractionation Cartridges will display the number of injections, number of guaranteed injections and number of batches remaining.
- Turbo CE-SDS Cartridges will only display the number of guaranteed injections and batches remaining.
- A warning will appear if a Cartridge Purge or Cartridge Post-Run Cleanup is required.
- The number remaining injections and batches will automatically update as follows:
 - The injections remaining will decrease by one as each injection completes.
 - The batches remaining will decrease by one when a Cartridge Purge is performed on a cIEF, CE-SDS PLUS or Turbo CE-SDS cartridge.
 - The batches and injections remaining will both decrease by one when a Cartridge Purge is performed on a cIEF Fractionation cartridge.
 - The batches and injections remaining will not change when a Cartridge Post-Run Cleanup is performed on all cartridges.

Cartridge Self-Test

You can run a series of self-tests on your cIEF, cIEF Fractionation, CE-SDS PLUS, or Turbo CE-SDS cartridge to make sure they're operating properly.

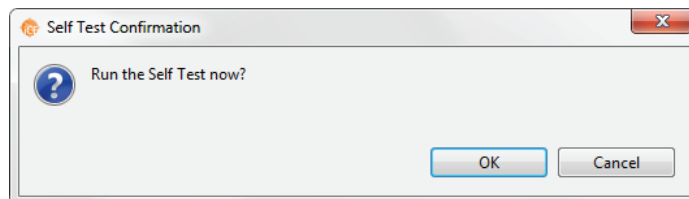
To perform a Cartridge Self-Test on a cIEF cartridge:

1. Confirm the electrolyte tanks are clean and dry.
2. Place an empty vial with a **blue pressure cap** in position P6.



3. Select **Instrument > Self Test > Cartridge**.

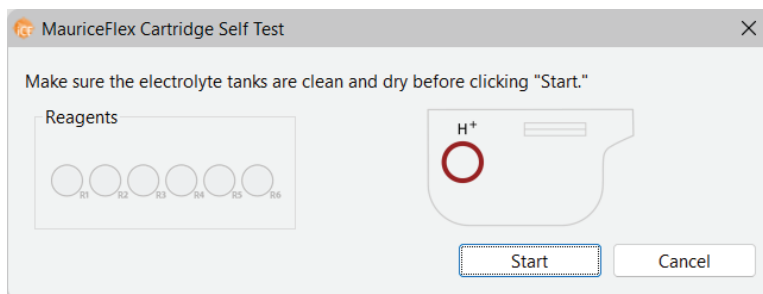
4. The following screen displays. Click OK.



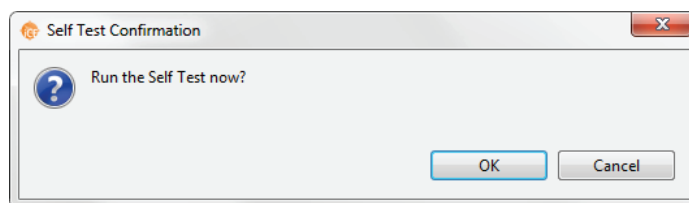
The test takes 3-8 minutes. Results will automatically display once the test's done. You can also view current and past self test results anytime by selecting **Instrument > Properties** and clicking **Test Log**. See "Self Test Logs" on page 470 for more information.

To perform a Cartridge Self-Test on a cIEF Fractionation cartridge:

1. Remove the Fractionation Adapter.
2. Confirm the Anolyte tank is clean and dry.



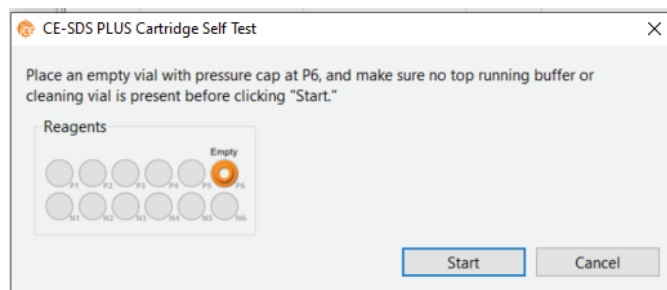
3. Select **Instrument > Self Test > Cartridge**.
4. The following screen displays. Click OK.



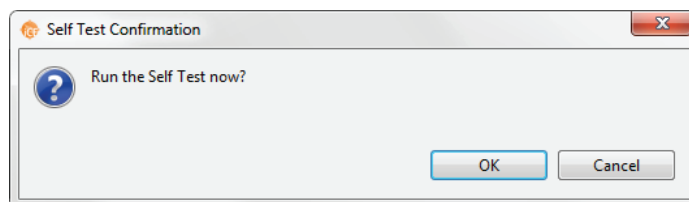
5. The test takes 3-8 minutes. Results will automatically display once the test's done. You can also view current and past self test results anytime by selecting **Instrument > Properties** and clicking **Test Log**. See "Self Test Logs" on page 470 for more information.

To perform a Cartridge Self-Test on a CE-SDS PLUS cartridge:

1. Remove the top Running Buffer and Cleaning vial.
2. Place an empty vial with an **orange pressure cap** in position P6.



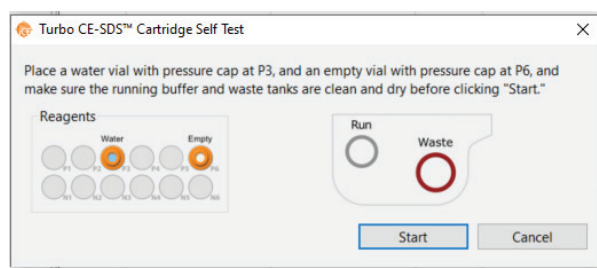
3. Select **Instrument > Self Test > Cartridge**.
4. The following screen displays. Click **OK**.



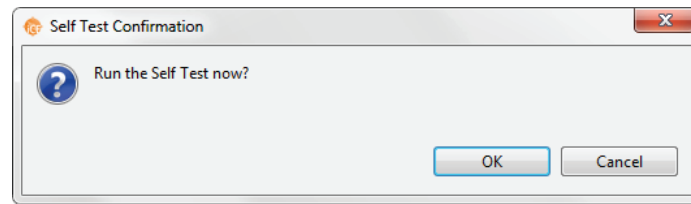
The test takes 3-8 minutes. Results will automatically display once the test's done. You can also view current and past self test results anytime by selecting **Instrument > Properties** and clicking **Test Log**. See "Self Test Logs" on page 470 for more information.

To perform a Cartridge Self-Test on a Turbo CE-SDS cartridge:

1. Confirm the Running Buffer and Waste Tanks are clean and dry.
2. Place an vial of water with an **orange pressure cap** in position P3 and an empty vial with an **orange pressure cap** in position P6.



3. Select **Instrument > Self Test > Cartridge**.
4. The following screen displays. Click **OK**.

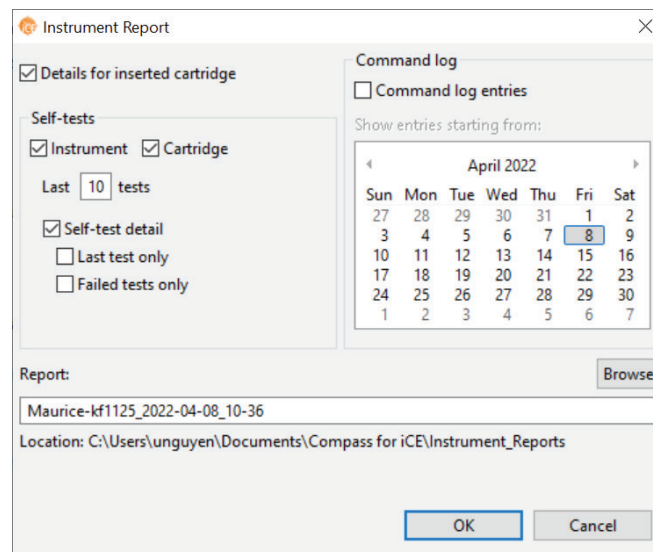


The test takes 3-8 minutes. Results will automatically display once the test's done. You can also view current and past self test results anytime by selecting **Instrument > Properties** and clicking **Test Log**. See "Self Test Logs" on page 470 for more information.

Instrument Reports

When Compass for iCE is connected to a Maurice system, you can generate an instrument report that includes instrument and cartridge info, self test records and details and command logs.

1. To run the report select **Instrument > Properties**.
2. Click **Report**. The Instrument Report window displays:



3. Select the information you want included in the report. You can change the location where the report saves by clicking **Browse**.
4. Click **OK**. The Instrument Report PDF is exported to the folder specified. Once the report is done, the folder opens for you automatically.

NOTE: Report PDFs generated by Compass for iCE are secured. They can be viewed and printed but can't be modified or renamed.

Example Instrument Report

Maurice kf1026

Instrument

Name	Maurice kf1026
Location	
Type	Maurice
Serial Number	kf1026
Instrument Software	4.2.2023.01.13.02.16.24.0f5300503
Adapter Block	plate
D2 Lamp Run Time (hours)	1201
Sample Chiller (°C)	25.3

Cartridge

Type	oIEF
Expires	Mar 2024
Serial Number	000000001
Injections per Batch	100
Injections Remaining	200
Injections Guaranteed	100
Batches Remaining	20

Instrument Self Test Summary - Last 10 Tests

Test Date	Status	Instrument Software	Username
2022-12-13 11:42	FAILED	4.2.2022.12.02.18.45.42.23a67ef7c	
2022-11-17 10:00	PASSED	4.1.2022.06.29.21.58.35.b9d03b73d	
2022-05-23 11:37	STOPPED	4.1.2022.05.11.05.58.39.c54f50194	
2021-03-17 20:54	PASSED	4.0.2020.07.11.04.16.13.ded2bc3	
2021-03-17 18:26	PASSED	4.0.2020.07.11.04.16.13.ded2bc3	
2021-02-08 18:07	PASSED	4.0.2020.07.11.04.16.13.ded2bc3	
2021-02-08 16:59	FAILED	4.0.2020.07.11.04.16.13.ded2bc3	
2021-02-04 17:05	PASSED	4.0.2020.07.11.04.16.13.ded2bc3	
2020-06-14 22:42	PASSED	3.1.2019.02.11.18.58.45.34dd4a5	
2020-05-25 18:21	PASSED	3.1.2019.02.11.18.58.45.34dd4a5	

Instrument Self Test Details

Self Test Date: 2022-12-13 11:42

Instrument Software: 4.2.2022.12.02.18.45.42.23a67ef7c

Username:

Name	Start	Result	Failure Reason
Vacuum Vent	11:42:49	PASSED	
Vacuum Leak	11:41:24	PASSED	
Unregulated Vac level	11:40:59	PASSED	
Temp Sensors	11:40:58	PASSED	
Pressure Vent	11:40:48	PASSED	
Pressure Leak	11:39:31	PASSED	
Low Vacuum Function	11:39:11	PASSED	
High Voltage	11:39:00	PASSED	
Temp Sensor Variance	11:38:58	PASSED	
Tray Move	11:37:07	PASSED	
Tray Jog	11:37:03	PASSED	
Tray Home	11:37:01	PASSED	
Tray Encoders	11:36:58	PASSED	
NFC Service	11:36:55	PASSED	
Disk Storage	11:36:54	PASSED	

Maurice kf1026

Self Test Date: 2022-11-17 10:00

Instrument Software: 4.1.2022.06.29.21.58.35.b9d03b73d

Username:

Name	Start	Result	Failure Reason
Sample Cooler	00:59:53	PASSED	
Sample Coolant Pump	00:59:52	PASSED	
Chamber Heater	00:59:15	PASSED	
Chamber Temperature Spec	00:59:14	PASSED	
Ambient Temperature Spec	00:59:13	PASSED	
Point Detector Light	00:58:32	PASSED	
Point Detector Dark	00:58:07	PASSED	
D2 Lamp On	00:57:36	PASSED	
Image Quality	00:57:12	PASSED	
Filter Wheel Move	00:56:44	PASSED	
Filter Wheel Home	00:56:42	PASSED	
Dark Masters	00:56:41	PASSED	
Camera	00:56:39	PASSED	
Vacuum Vent	00:56:34	PASSED	
Vacuum Leak	00:55:08	PASSED	
Unregulated Vac level	00:54:43	PASSED	
Temp Sensors	00:54:42	PASSED	
Pressure Vent	00:54:33	PASSED	
Pressure Leak	00:53:15	PASSED	
Low Vacuum Function	00:52:55	PASSED	
High Voltage	00:52:44	PASSED	
Temp Sensor Variance	00:52:43	PASSED	
Tray Move	00:51:12	PASSED	
Tray Jog	00:51:08	PASSED	
Tray Home	00:51:06	PASSED	
Tray Encoders	00:51:03	PASSED	
NFC Service	00:51:00	PASSED	

Maurice kf1026

Disk Storage	09:50:59	PASSED	
--------------	----------	--------	--

Self Test Date: 2022-05-23 11:37

Instrument Software: 4.1.2022.05.11.05.58.39.c54f50194

Username:

Name	Start	Result	Failure Reason
Tray Move	11:35:59	PASSED	
Tray Jog	11:35:55	PASSED	
Tray Home	11:35:53	PASSED	
Tray Encoders	11:35:50	PASSED	
NFC Service	11:35:47	PASSED	
Disk Storage	11:35:46	PASSED	

Self Test Date: 2021-03-17 20:54

Instrument Software: 4.0.2020.07.11.04.16.13.ded2bc3

Username:

Name	Start	Result	Failure Reason
Sample Cooler	12:54:03	PASSED	
Sample Coolant Pump	12:54:02	PASSED	
Chamber Heater	12:53:01	PASSED	
Chamber Temperature Spec	12:53:00	PASSED	
Ambient Temperature Spec	12:52:59	PASSED	
Point Detector Light	12:52:18	PASSED	
Point Detector Dark	12:51:53	PASSED	
D2 Lamp On	12:51:22	PASSED	
Camera UV Illumination	12:51:15	PASSED	
Image Quality	12:50:52	PASSED	
Filter Wheel Move	12:50:24	PASSED	
Filter Wheel Home	12:50:23	PASSED	
Dark Masters	12:50:22	PASSED	
Camera	12:50:19	PASSED	
Vacuum Vent	12:50:14	PASSED	

Maurice kf1026

Vacuum Leak	12:48:48	PASSED	
Unregulated Vac level	12:48:23	PASSED	
Temp Sensors	12:48:22	PASSED	
Pressure Vent	12:48:13	PASSED	
Pressure Leak	12:46:51	PASSED	
Low Vacuum Function	12:46:30	PASSED	
High Voltage	12:46:19	PASSED	
Temp Sensor Variance	12:46:18	PASSED	
Tray Move	12:44:47	PASSED	
Tray Jog	12:44:43	PASSED	
Tray Home	12:44:41	PASSED	
Tray Encoders	12:44:38	PASSED	
NFC Service	12:44:35	PASSED	
Disk Storage	12:44:34	PASSED	

Self Test Date: 2021-03-17 18:26

Instrument Software: 4.0.2020.07.11.04.16.13.ded2bc3

Username:

Name	Start	Result	Failure Reason
Sample Cooler	10:25:28	PASSED	
Sample Coolant Pump	10:25:27	PASSED	
Chamber Heater	10:24:28	PASSED	
Chamber Temperature Spec	10:24:27	PASSED	
Ambient Temperature Spec	10:24:26	PASSED	
Point Detector Light	10:23:45	PASSED	
Point Detector Dark	10:23:20	PASSED	
D2 Lamp On	10:22:50	PASSED	
Camera UV Illumination	10:22:39	PASSED	
Image Quality	10:22:17	PASSED	
Filter Wheel Move	10:21:48	PASSED	
Filter Wheel Home	10:21:47	PASSED	
Dark Masters	10:21:46	PASSED	

Maurice kf1026

Camera	10:21:44	PASSED	
Vacuum Vent	10:21:38	PASSED	
Vacuum Leak	10:20:13	PASSED	
Unregulated Vac level	10:19:48	PASSED	
Temp Sensors	10:19:47	PASSED	
Pressure Vent	10:19:37	PASSED	
Pressure Leak	10:18:17	PASSED	
Low Vacuum Function	10:17:56	PASSED	
High Voltage	10:17:45	PASSED	
Temp Sensor Variance	10:17:44	PASSED	
Tray Move	10:16:14	PASSED	
Tray Jog	10:16:10	PASSED	
Tray Home	10:16:08	PASSED	
Tray Encoders	10:16:05	PASSED	
NFC Service	10:16:02	PASSED	
Disk Storage	10:16:01	PASSED	

Self Test Date: 2021-02-08 18:07

Instrument Software: 4.0.2020.07.11.04.16.13.ded2bc3

Username:

Name	Start	Result	Failure Reason
Sample Cooler	11:06:32	PASSED	
Sample Coolant Pump	11:06:31	PASSED	
Chamber Heater	11:05:32	PASSED	
Chamber Temperature Spec	11:05:31	PASSED	
Ambient Temperature Spec	11:05:30	PASSED	
Point Detector Light	11:04:49	PASSED	
Point Detector Dark	11:04:24	PASSED	
D2 Lamp On	11:03:53	PASSED	
Camera UV Illumination	11:03:43	PASSED	
Image Quality	11:03:20	PASSED	
Filter Wheel Move	11:02:52	PASSED	

Maurice kf1026

Filter Wheel Home	11:02:51	PASSED	
Dark Masters	11:02:50	PASSED	
Camera	11:02:48	PASSED	
Vacuum Vent	11:02:43	PASSED	
Vacuum Leak	11:01:17	PASSED	
Unregulated Vac level	11:00:52	PASSED	
Temp Sensors	11:00:51	PASSED	
Pressure Vent	11:00:40	PASSED	
Pressure Leak	10:59:18	PASSED	
Low Vacuum Function	10:58:57	PASSED	
High Voltage	10:58:46	PASSED	
Temp Sensor Variance	10:58:45	PASSED	
Tray Move	10:57:14	PASSED	
Tray Jog	10:57:11	PASSED	
Tray Home	10:57:09	PASSED	
Tray Encoders	10:57:05	PASSED	
NFC Service	10:57:00	PASSED	
Disk Storage	10:56:59	PASSED	

Self Test Date: 2021-02-08 16:59

Instrument Software: 4.0.2020.07.11.04.16.13.ded2bc3

Username:

Name	Start	Result	Failure Reason
Sample Cooler	09:59:09	PASSED	
Sample Coolant Pump	09:59:08	PASSED	
Chamber Heater	09:58:42	PASSED	
Chamber Temperature Spec	09:58:41	FAILED	Detected chamber temp 24.6C, 09:58:41 AM below control range (27.0C - 32.5C)
Ambient Temperature Spec	09:58:40	PASSED	
Point Detector Light	09:57:58	PASSED	
Point Detector Dark	09:57:33	PASSED	
D2 Lamp On	09:57:03	PASSED	
Camera UV Illumination	09:56:55	FAILED	Dark image must be darker than Fluorescence image

Maurice kf1026

Image Quality	09:56:54	FAILED	Camera temperature reads -47.14; must read -49.90 or below to test image quality. Ensure camera has had sufficient time to cool before re-running test.
Filter Wheel Move	09:56:25	PASSED	
Filter Wheel Home	09:56:24	PASSED	
Dark Masters	09:56:23	PASSED	
Camera	09:56:20	PASSED	
Vacuum Vent	09:56:15	PASSED	
Vacuum Leak	09:54:50	PASSED	
Unregulated Vac level	09:54:24	PASSED	
Temp Sensors	09:54:23	PASSED	
Pressure Vent	09:54:14	PASSED	
Pressure Leak	09:52:53	PASSED	
Low Vacuum Function	09:52:32	PASSED	
High Voltage	09:52:21	PASSED	
Temp Sensor Variance	09:52:20	PASSED	
Tray Move	09:50:50	PASSED	
Tray Jog	09:50:46	PASSED	
Tray Home	09:50:44	PASSED	
Tray Encoders	09:50:41	PASSED	
NFC Service	09:50:38	PASSED	
Disk Storage	09:50:37	PASSED	

Self Test Date: 2021-02-04 17:05

Instrument Software: 4.0.2020.07.11.04.16.13.ded2bc3

Username:

Name	Start	Result	Failure Reason
Sample Cooler	10:04:33	PASSED	
Sample Coolant Pump	10:04:32	PASSED	
Chamber Heater	10:03:45	PASSED	
Chamber Temperature Spec	10:03:44	PASSED	
Ambient Temperature Spec	10:03:43	PASSED	
Point Detector Light	10:03:02	PASSED	
Point Detector Dark	10:02:37	PASSED	

Maurice kf1026

D2 Lamp On	10:02:07	PASSED	
Camera UV Illumination	10:01:57	PASSED	
Image Quality	10:01:34	PASSED	
Filter Wheel Move	10:01:06	PASSED	
Filter Wheel Home	10:01:04	PASSED	
Dark Masters	10:01:03	PASSED	
Camera	10:01:01	PASSED	
Vacuum Vent	10:00:56	PASSED	
Vacuum Leak	09:59:30	PASSED	
Unregulated Vac level	09:59:05	PASSED	
Temp Sensors	09:59:04	PASSED	
Pressure Vent	09:58:54	PASSED	
Pressure Leak	09:57:32	PASSED	
Low Vacuum Function	09:57:12	PASSED	
High Voltage	09:57:01	PASSED	
Temp Sensor Variance	09:56:59	PASSED	
Tray Move	09:55:29	PASSED	
Tray Jog	09:55:25	PASSED	
Tray Home	09:55:23	PASSED	
Tray Encoders	09:55:20	PASSED	
NFC Service	09:55:17	PASSED	
Disk Storage	09:55:16	PASSED	

Self Test Date: 2020-06-14 22:42

Instrument Software: 3.1.2019.02.11.18.58.45.34d4a5

Username:

Name	Start	Result	Failure Reason
Sample Cooler	14:41:41	PASSED	
Sample Coolant Pump	14:41:40	PASSED	
Chamber Heater	14:40:59	PASSED	
Chamber Temperature Spec	14:40:58	PASSED	
Ambient Temperature Spec	14:40:57	PASSED	

Maurice kf1026

Point Detector Light	14:40:16	PASSED	
Point Detector Dark	14:39:51	PASSED	
D2 Lamp On	14:39:20	PASSED	
Camera UV Illumination	14:39:10	PASSED	
Image Quality	14:38:47	PASSED	
Filter Wheel Move	14:38:19	PASSED	
Filter Wheel Home	14:38:18	PASSED	
Dark Masters	14:38:17	PASSED	
Camera	14:38:14	PASSED	
Vacuum Vent	14:38:08	PASSED	
Vacuum Leak	14:36:43	PASSED	
Unregulated Vac level	14:36:18	PASSED	
Temp Sensors	14:36:17	PASSED	
Pressure Vent	14:36:07	PASSED	
Pressure Leak	14:34:46	PASSED	
Low Vacuum Function	14:34:26	PASSED	
High Voltage	14:34:15	PASSED	
Temp Sensor Variance	14:34:14	PASSED	
Tray Move	14:32:43	PASSED	
Tray Jog	14:32:40	PASSED	
Tray Home	14:32:38	PASSED	
Tray Encoders	14:32:35	PASSED	
NFC Service	14:32:32	PASSED	
Disk Storage	14:32:31	PASSED	

Self Test Date: 2020-05-25 18:21

Instrument Software: 3.1.2019.02.11.18.58.45.34dd4a5

Username:

Name	Start	Result	Failure Reason
Sample Cooler	10:21:01	PASSED	
Sample Coolant Pump	10:21:00	PASSED	
Chamber Heater	10:20:03	PASSED	
Chamber Temperature Spec	10:20:02	PASSED	

Maurice kf1026

Ambient Temperature Spec	10:20:01	PASSED	
Point Detector Light	10:19:21	PASSED	
Point Detector Dark	10:18:56	PASSED	
D2 Lamp On	10:18:26	PASSED	
Camera UV Illumination	10:18:15	PASSED	
Image Quality	10:17:52	PASSED	
Filter Wheel Move	10:17:24	PASSED	
Filter Wheel Home	10:17:23	PASSED	
Dark Masters	10:17:22	PASSED	
Camera	10:17:21	PASSED	
Vacuum Vent	10:17:15	PASSED	
Vacuum Leak	10:15:50	PASSED	
Unregulated Vac level	10:15:25	PASSED	
Temp Sensors	10:15:24	PASSED	
Pressure Vent	10:15:14	PASSED	
Pressure Leak	10:13:52	PASSED	
Low Vacuum Function	10:13:32	PASSED	
High Voltage	10:13:20	PASSED	
Temp Sensor Variance	10:13:19	PASSED	
Tray Move	10:11:49	PASSED	
Tray Jog	10:11:45	PASSED	
Tray Home	10:11:44	PASSED	
Tray Encoders	10:11:40	PASSED	
NFC Service	10:11:36	PASSED	
Disk Storage	10:11:35	PASSED	

Maurice kf1026**Cartridge Self Test Summary - Last 10 Tests**

Test Date	Status	Cartridge S/N	Instrument Software	Username
2022-11-30 09:51	PASSED	8220511003	4.2.2022.11.23.21.59.20.d530e42a9	
2022-11-29 10:43	PASSED	8220511003	4.2.2022.11.23.21.59.20.d530e42a9	
2022-11-28 10:13	PASSED	8220511003	4.2.2022.11.23.21.59.20.d530e42a9	
2022-11-17 10:43	PASSED	8220615870	4.2.2022.11.07.17.15.11.3643af972	
2022-07-01 13:44	PASSED	8220503001	4.1.2022.06.29.21.58.35.b6d03b73d	admin
2022-06-30 17:47	PASSED	8220503001	4.1.2022.06.29.21.58.35.b6d03b73d	
2022-06-30 17:37	FAILED	8220503001	4.1.2022.06.29.21.58.35.b6d03b73d	
2022-06-30 11:07	PASSED	8220503001	4.1.2022.06.29.21.58.35.b6d03b73d	
2022-06-07 10:50	PASSED	8220503001	4.1.2022.05.28.18.36.41.aa06fd9b0	
2022-05-24 13:06	PASSED	8220503001	4.1.2022.05.11.05.58.39.c54f50194	admin

Cartridge Self Test Details

Self Test Date: 2022-11-30 09:51

Instrument Software: 4.2.2022.11.23.21.59.20.d530e42a9

Username:

Cartridge S/N: 8220511003

Name	Start	Result	Failure Reason
Cartridge Clog	09:44:13	PASSED	

Self Test Date: 2022-11-29 10:43

Instrument Software: 4.2.2022.11.23.21.59.20.d530e42a9

Username:

Cartridge S/N: 8220511003

Name	Start	Result	Failure Reason
Cartridge Clog	10:36:01	PASSED	

Self Test Date: 2022-11-28 10:13

Instrument Software: 4.2.2022.11.23.21.59.20.d530e42a9

Username:

Cartridge S/N: 8220511003

Name	Start	Result	Failure Reason
------	-------	--------	----------------

Maurice kf1026

Cartridge Clog	10:06:17	PASSED	
----------------	----------	--------	--

Self Test Date: 2022-11-17 10:43

Instrument Software: 4.2.2022.11.07.17.15.11.3643af972

Username:

Cartridge S/N: 8220615870

Name	Start	Result	Failure Reason
Cartridge Clog	10:36:03	PASSED	

Self Test Date: 2022-07-01 13:44

Instrument Software: 4.1.2022.06.29.21.58.35.b9d03b73d

Username: admin

Cartridge S/N: 8220503001

Name	Start	Result	Failure Reason
Cartridge Clog	13:37:22	PASSED	

Self Test Date: 2022-06-30 17:47

Instrument Software: 4.1.2022.06.29.21.58.35.b9d03b73d

Username:

Cartridge S/N: 8220503001

Name	Start	Result	Failure Reason
Cartridge Clog	17:40:49	PASSED	

Self Test Date: 2022-06-30 17:37

Instrument Software: 4.1.2022.06.29.21.58.35.b9d03b73d

Username:

Cartridge S/N: 8220503001

Name	Start	Result	Failure Reason
Cartridge Clog	17:30:38	FAILED	Mean reading with water 2267078.1, mean with air 2266837.3, diff 0.01%, minimum 10.0%

Maurice kf1026

Self Test Date: 2022-06-30 11:07

Instrument Software: 4.1.2022.06.29.21.58.35.b9d03b73d

Username:

Cartridge S/N: 8220503001

Name	Start	Result	Failure Reason
Cartridge Clog	11:00:45	PASSED	

Self Test Date: 2022-06-07 10:50

Instrument Software: 4.1.2022.05.26.18.36.41.aa06f6b0

Username:

Cartridge S/N: 8220503001

Name	Start	Result	Failure Reason
Cartridge Clog	10:43:28	PASSED	

Self Test Date: 2022-05-24 13:06

Instrument Software: 4.1.2022.05.11.05.58.39.c54f50194

Username: admin

Cartridge S/N: 8220503001

Name	Start	Result	Failure Reason
Cartridge Clog	12:59:36	PASSED	

Created By: jlevinsky@bruker.com Mon 1:00 PM Mar 13, 2023 PDT
 Maurice kf1026
 Computer: USBKIO-2K792J3-L

Software Version: Compass for ICE 4.0.0, Build ID: 0222

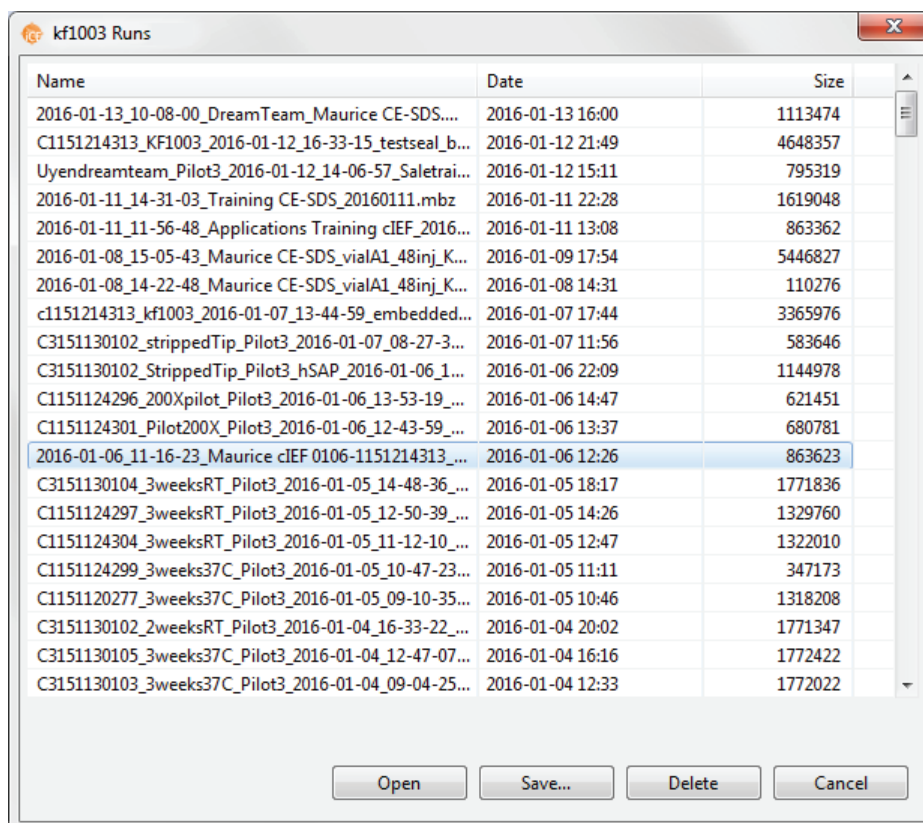


Page 14 of 14

Viewing Log Files

Runs Log

To see a history of all runs your system has performed, select **Instrument > Runs**:

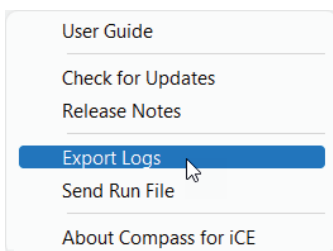


- **To open a run file:** Select a run file from the list and click **Open**.
- **To save a run file:** Select a run file from the list and click **Save**. This lets you save a copy of a completed run or one in progress to either a USB drive or the local computer.
- **To delete a run file:** Select a run file from the list and click **Delete**. The run file will be deleted from the history and from the Run file in the Compass for iCE directory.

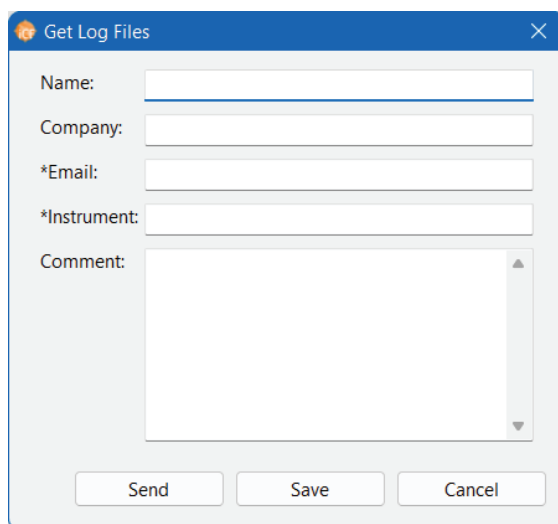
Sending a Log File to Technical Support

If the computer you're using has an internet connection, Compass for iCE can zip and send a log file directly to ProteinSimple Technical Support.

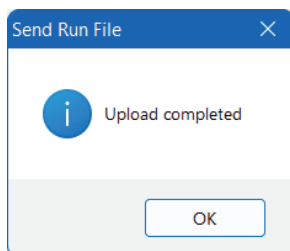
1. Connect to an instrument.
2. Select **Help** in the main menu and click **Export Logs**.



3. A window will display where you can enter information. The email and instrument fields are required.

A screenshot of a dialog box titled 'Get Log Files'. It has a blue header bar with a close button. The dialog contains several input fields: 'Name:', 'Company:', '*Email:', '*Instrument:', and 'Comment:'. The 'Email' and 'Instrument' fields are marked with an asterisk, indicating they are required. At the bottom, there are three buttons: 'Send', 'Save', and 'Cancel'.

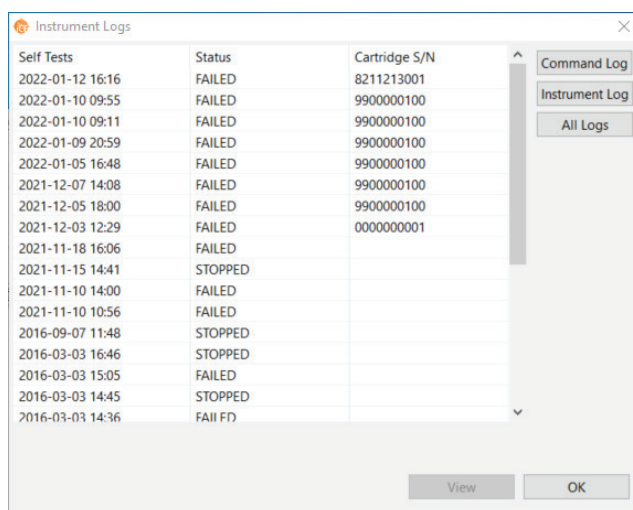
4. Click **Send**. When the file upload is complete, the following message displays. Click **OK**.



5. Click **Save** to save the file directly to the computer. A window will display where you can enter a file name for the zipped log file.

Instrument Log

1. Select **Instrument > Properties**.
2. Click **Logs**. The Instrument Logs window displays:



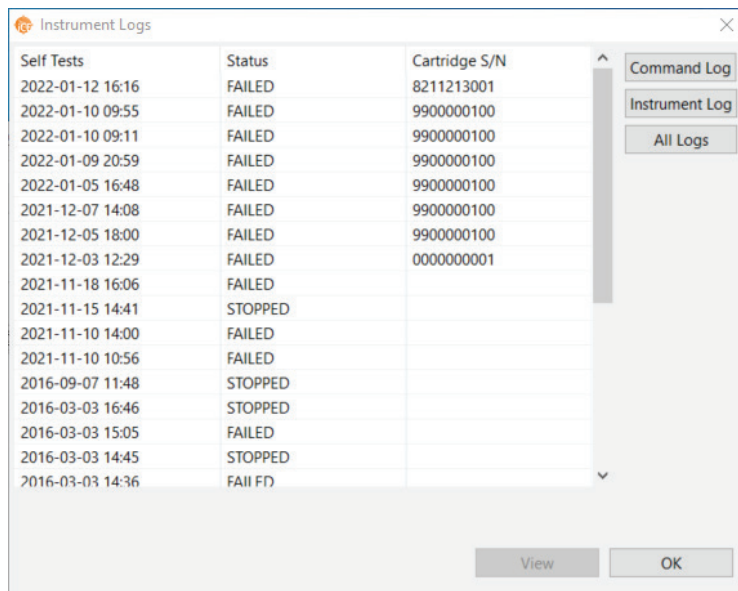
3. Click **Instrument Log**.



4. Click **Save File As** to save a copy of the log file.

Self Test Logs

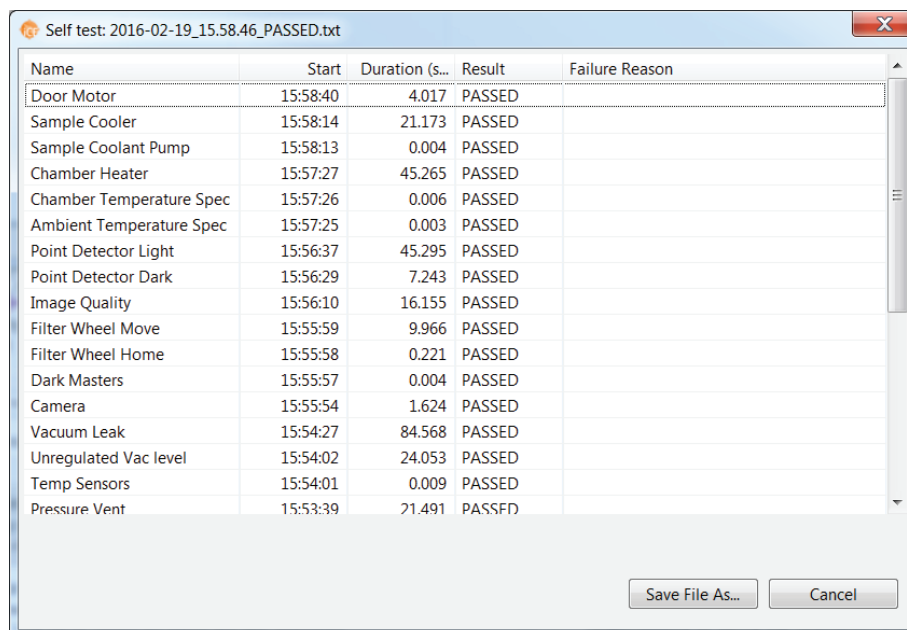
1. Select **Instrument > Properties**.
2. Click **Logs**. A list of dates each self test was run displays:



The screenshot shows the 'Instrument Logs' window. It contains a table with three columns: 'Self Tests', 'Status', and 'Cartridge S/N'. The table lists 20 self-test entries with their respective dates, times, and statuses. To the right of the table are three buttons: 'Command Log', 'Instrument Log', and 'All Logs'. At the bottom right are 'View' and 'OK' buttons.

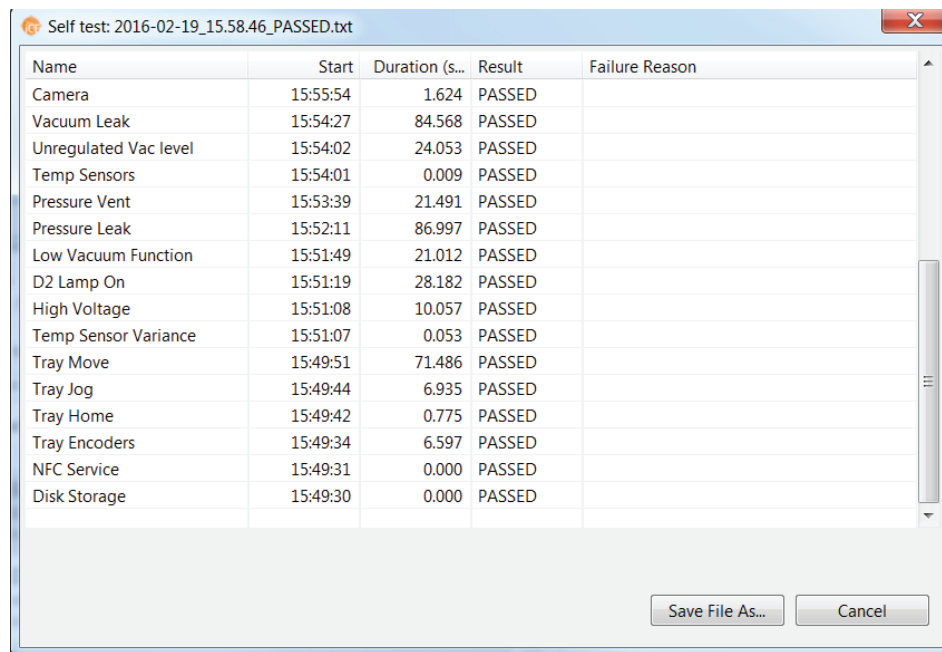
Self Tests	Status	Cartridge S/N
2022-01-12 16:16	FAILED	8211213001
2022-01-10 09:55	FAILED	9900000100
2022-01-10 09:11	FAILED	9900000100
2022-01-09 20:59	FAILED	9900000100
2022-01-05 16:48	FAILED	9900000100
2021-12-07 14:08	FAILED	9900000100
2021-12-05 18:00	FAILED	9900000100
2021-12-03 12:29	FAILED	0000000001
2021-11-18 16:06	FAILED	
2021-11-15 14:41	STOPPED	
2021-11-10 14:00	FAILED	
2021-11-10 10:56	FAILED	
2016-09-07 11:48	STOPPED	
2016-03-03 16:46	STOPPED	
2016-03-03 15:05	FAILED	
2016-03-03 14:45	STOPPED	
2016-03-03 14:36	FAIL FD	

3. Select a test date in the list and click **View** to see the individual test results:



The screenshot shows the 'Self test: 2016-02-19_15.58.46_PASSED.txt' window. It contains a table with five columns: 'Name', 'Start', 'Duration (s...)', 'Result', and 'Failure Reason'. The table lists 20 test items, all of which passed. At the bottom right are 'Save File As...' and 'Cancel' buttons.

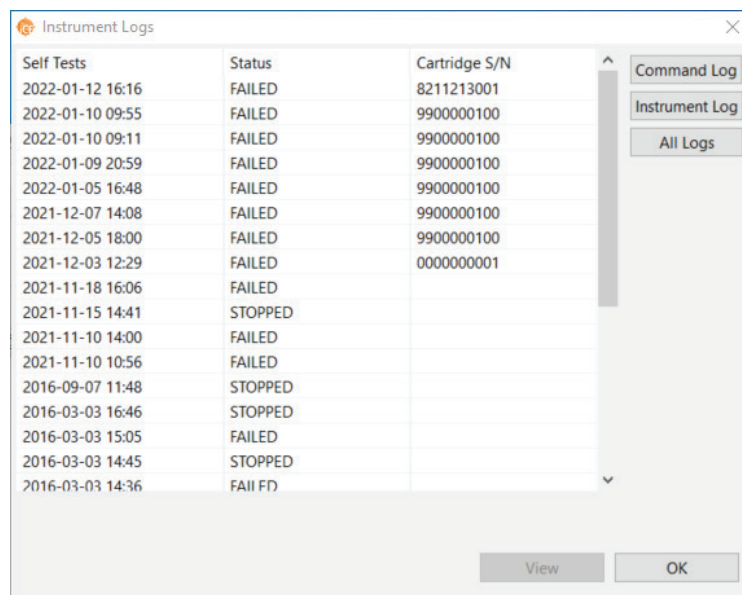
Name	Start	Duration (s...)	Result	Failure Reason
Door Motor	15:58:40	4.017	PASSED	
Sample Cooler	15:58:14	21.173	PASSED	
Sample Coolant Pump	15:58:13	0.004	PASSED	
Chamber Heater	15:57:27	45.265	PASSED	
Chamber Temperature Spec	15:57:26	0.006	PASSED	
Ambient Temperature Spec	15:57:25	0.003	PASSED	
Point Detector Light	15:56:37	45.295	PASSED	
Point Detector Dark	15:56:29	7.243	PASSED	
Image Quality	15:56:10	16.155	PASSED	
Filter Wheel Move	15:55:59	9.966	PASSED	
Filter Wheel Home	15:55:58	0.221	PASSED	
Dark Masters	15:55:57	0.004	PASSED	
Camera	15:55:54	1.624	PASSED	
Vacuum Leak	15:54:27	84.568	PASSED	
Unregulated Vac level	15:54:02	24.053	PASSED	
Temp Sensors	15:54:01	0.009	PASSED	
Pressure Vent	15:53:39	71.491	PASSED	



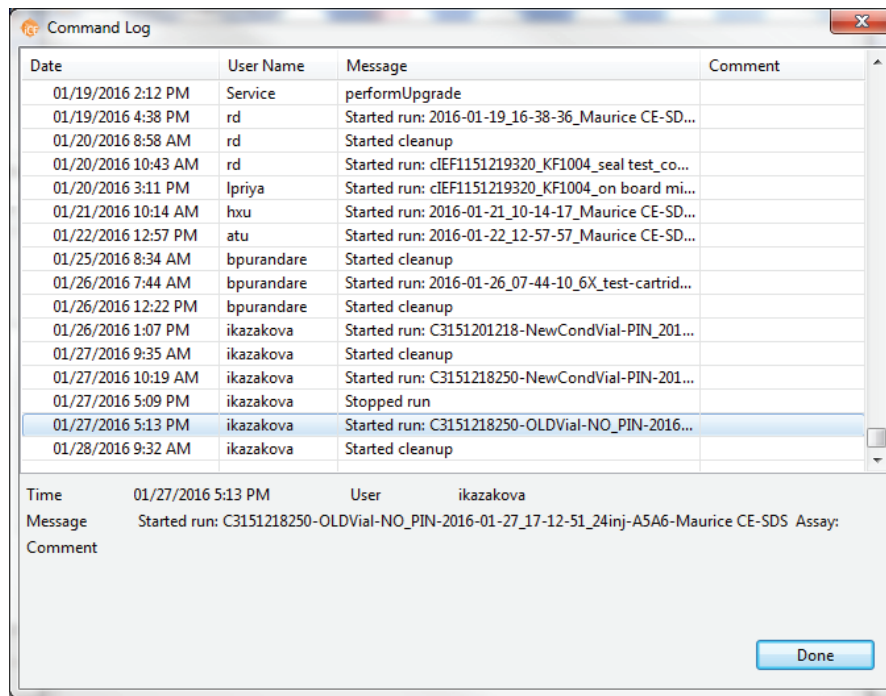
4. Click **Save File As** to save a copy of the test log file.

Command Log

1. Select **Instrument > Properties** to display your system's properties.
2. Click **Logs**. The Instrument Logs window displays:



3. Click **Command Log**. A list of system commands displays:



Date	User Name	Message	Comment
01/19/2016 2:12 PM	Service	performUpgrade	
01/19/2016 4:38 PM	rd	Started run: 2016-01-19_16-38-36_Maurice CE-SD...	
01/20/2016 8:58 AM	rd	Started cleanup	
01/20/2016 10:43 AM	rd	Started run: cIEF1151219320_KF1004_seal test_co...	
01/20/2016 3:11 PM	lpriya	Started run: cIEF1151219320_KF1004_on board mi...	
01/21/2016 10:14 AM	hxu	Started run: 2016-01-21_10-14-17_Maurice CE-SD...	
01/22/2016 12:57 PM	atu	Started run: 2016-01-22_12-57-57_Maurice CE-SD...	
01/25/2016 8:34 AM	bpurandare	Started cleanup	
01/26/2016 7:44 AM	bpurandare	Started run: 2016-01-26_07-44-10_6X_test-cartrid...	
01/26/2016 12:22 PM	bpurandare	Started cleanup	
01/26/2016 1:07 PM	ikazakova	Started run: C3151201218-NewCondVial-PIN_201...	
01/27/2016 9:35 AM	ikazakova	Started cleanup	
01/27/2016 10:19 AM	ikazakova	Started run: C3151218250-NewCondVial-PIN-201...	
01/27/2016 5:09 PM	ikazakova	Stopped run	
01/27/2016 5:13 PM	ikazakova	Started run: C3151218250-OLDVial-NO_PIN-2016...	
01/28/2016 9:32 AM	ikazakova	Started cleanup	

Time	01/27/2016 5:13 PM	User	ikazakova
Message	Started run: C3151218250-OLDVial-NO_PIN-2016-01-27_17-12-51_24inj-ASA6-Maurice CE-SDS Assay;		
Comment			

Done

Chapter 18:

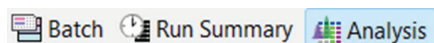
cIEF Data Analysis

Chapter Overview

- Analysis Screen Overview
- Opening Run Files
- How Run Data is Displayed
- Viewing Run Data
- Data Notifications and Warnings
- Checking Your Results
- Naming Peaks
- Group Statistics
- Copying Results Tables and Graphs
- Exporting Run Files
- Changing Sample Protein Identification
- Changing the Electropherogram View
- Closing Run Files
- Analysis Settings Overview
- Detection Settings
- Peak Names Settings
- Peak Fit Analysis Settings
- Manual Peak Integration
- Advanced Analysis Settings
- pI Markers Analysis Settings
- Injection Reports
- Importing and Exporting Analysis Settings

Analysis Screen Overview

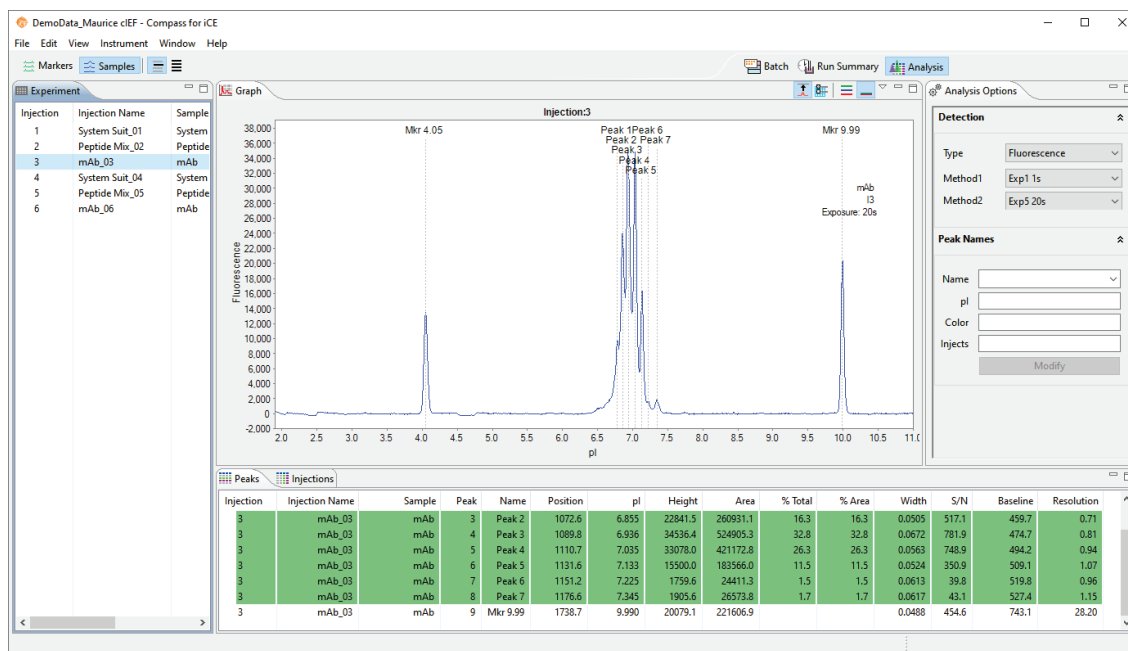
You can use the Analysis screen to view electropherograms and tabulated results for your injections. If any post-run analysis is needed, you can do it here too. To get to this screen, click the **Analysis** screen tab:



Analysis Screen Panes

The Analysis screen has four panes:

- **Experiment** - Lists the injection number, sample IDs, sample locations and methods for each injection in the run and lets you get a quick view of method parameters.
- **Graph** - Displays the electropherograms for sample proteins or pI markers.
- **Peaks** - Shows the tabulated results for sample proteins and pI markers.
- **Injections** - Displays a list of the sample proteins Compass for iCE names automatically using the user-defined peak name analysis parameters.
- **Analysis Options** - Lets you view absorbance or fluorescence data for the run and view, change and add new custom peak name settings.



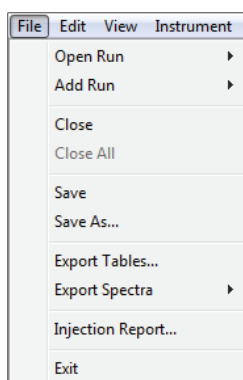
Software Menus Active in the Analysis Screen

These main menu items are active in the Analysis Screen:

- File
- Edit
- View
- Instrument (when Compass for iCE is connected to Maurice, Maurice C. or MauriceFlex.)
- Window
- Help

File Menu

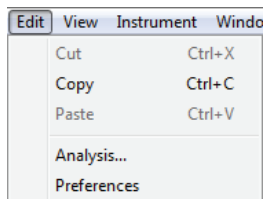
These File menu options are active:



- **Open Run** - Opens a run file.
- **Add Run** - Lets you open and view other run files besides the one that's already open.
- **Close** - Closes the run file currently being viewed.
- **Close All** - Closes all open run files.
- **Save/Save As** - If you made changes in the Analysis screen before you went to the Run Summary screen, this saves your changes to the run file.
- **Export Tables** - Exports the results for all injections in the run in .txt format.
- **Export Spectra** - Exports the raw and analyzed data traces and background for each injection in the run in .txt or .cdf format.
- **Injection Report** - Exports the raw and analyzed data, IV plot, peaks table, sample and system info for individual injections as PDF files. You can also export the run history with all analysis events.
- **Exit** - Closes Compass for iCE.

Edit Menu

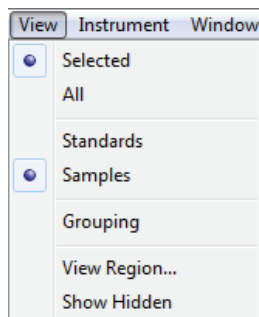
These Edit menu options are active:



- **Copy** - Lets you copy data shown in the graph, lane, peaks or injections pane. See “Copying Results Tables and Graphs” on page 501 for more information.
- **Analysis** - Displays the analysis settings used to analyze the run data and lets you change them as needed. See “Analysis Settings Overview” on page 528 for more information.
- **Preferences** - Lets you set and save your preferences for data export, graph colors, grouped data and Twitter settings. See Chapter 21: “Setting Your Preferences” for more information.

View Menu

These View menu options are active:



- **Selected (Single View)** - Displays the data for only the injections selected.
- **All (Multiple View)** - Displays data for all injections so you can scroll through them.
- **Markers** - Lets you view data just for the pI markers in your injections.
- **Samples** - Lets you view data just for sample proteins in your injections.
- **Grouping** - Displays data for injection groups.
- **View Region** - Lets you change the x-axis range of the data displayed.
- **Show Hidden** - Shows injections that are hidden from the data view.

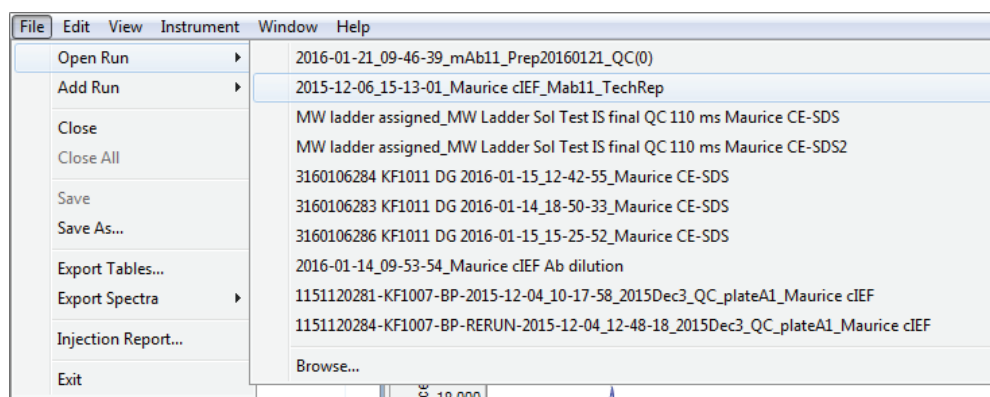
Opening Run Files

You can open one run file or multiple files at the same time to compare information between runs.

NOTE: cIEF run files that use the FL458 nm fluorescence detection filter can only be opened in Compass for iCE 2.1 or later.

Opening One Run File

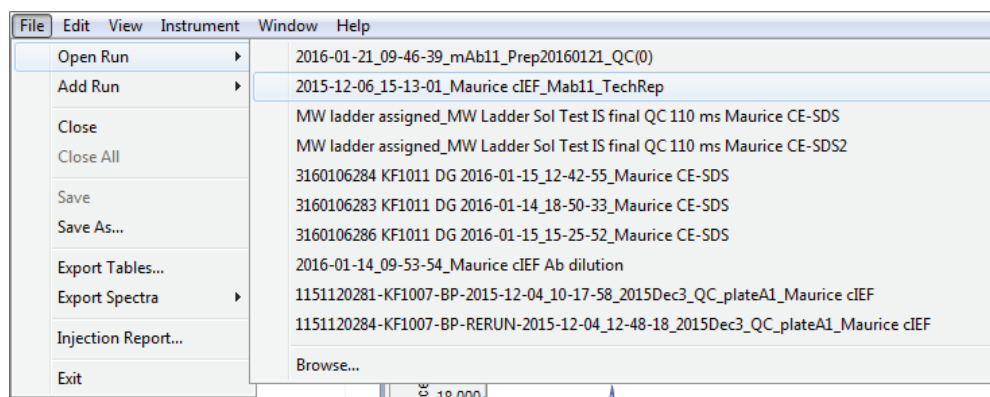
1. Select **File** in the main menu and click **Open Run**.



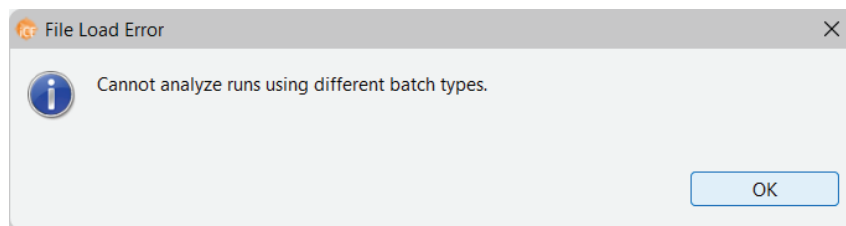
2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.

Opening Multiple Run Files

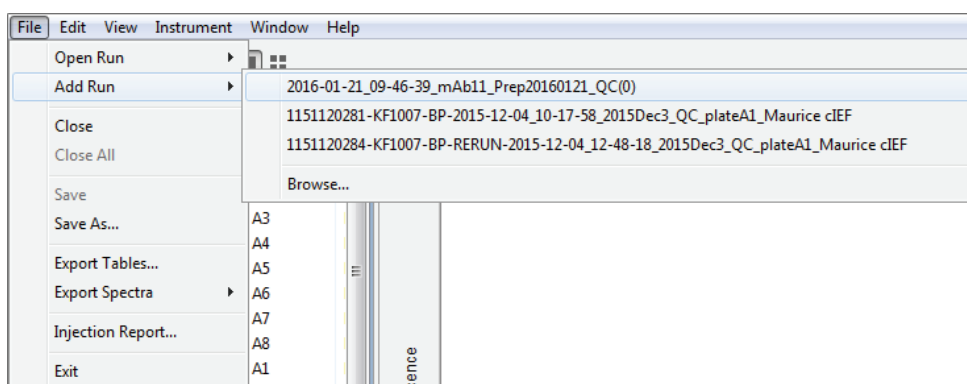
1. To open the first run file, select **File** in the main menu and click **Open Run**.



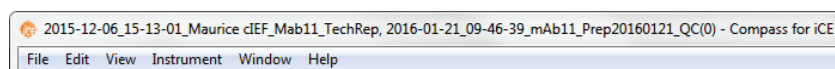
2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file. If the run selected is not a cIEF batch, an alert will appear.



3. To open another run file, select **File** in the main menu and click **Add Run**.



4. A list of cIEF runs will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.
5. When a run is added, its data appends to the open run file and displays as a second set of injections in all screen panes. The second run file name also appears in the title bar:



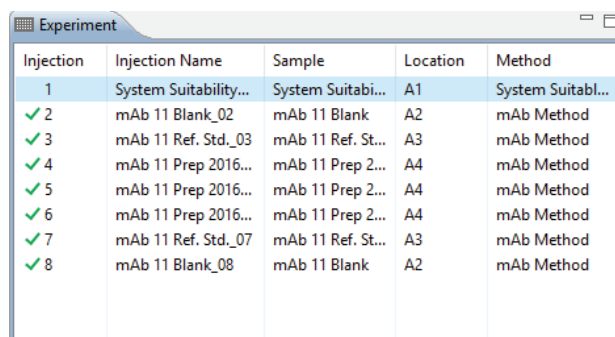
6. Repeat the last two steps to add additional runs.

How Run Data is Displayed

Data in the run file is organized for easy review.

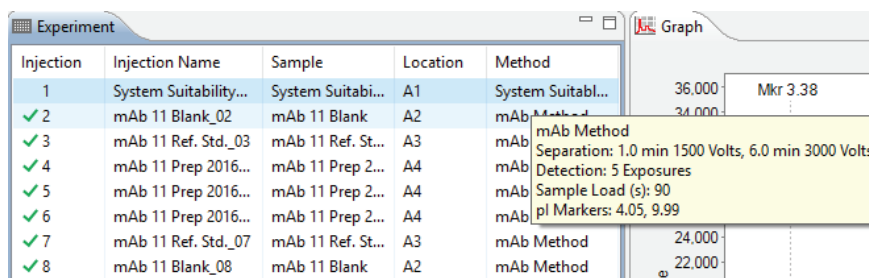
Experiment Pane: Batch Injection Information

The Experiment pane lists all the injections performed in the run, which samples were used for each, the sample location in the 96-well plate or 48-vial tray for cIEF batches and in the 96-well plate for MauriceFlex cIEF batches and the method used.



Injection	Injection Name	Sample	Location	Method
1	System Suitability...	System Suitabi...	A1	System Suitabl...
✓ 2	mAb 11 Blank_02	mAb 11 Blank	A2	mAb Method
✓ 3	mAb 11 Ref. Std._03	mAb 11 Ref. St...	A3	mAb Method
✓ 4	mAb 11 Prep 2016...	mAb 11 Prep 2...	A4	mAb Method
✓ 5	mAb 11 Prep 2016...	mAb 11 Prep 2...	A4	mAb Method
✓ 6	mAb 11 Prep 2016...	mAb 11 Prep 2...	A4	mAb Method
✓ 7	mAb 11 Ref. Std._07	mAb 11 Ref. St...	A3	mAb Method
✓ 8	mAb 11 Blank_08	mAb 11 Blank	A2	mAb Method

- **To view all columns** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.
- **To view method parameters** - Hover the mouse over a method name.



Injection	Injection Name	Sample	Location	Method
1	System Suitability...	System Suitabi...	A1	System Suitabl...
✓ 2	mAb 11 Blank_02	mAb 11 Blank	A2	mAb Method
✓ 3	mAb 11 Ref. Std._03	mAb 11 Ref. St...	A3	mAb Method
✓ 4	mAb 11 Prep 2016...	mAb 11 Prep 2...	A4	mAb Method
✓ 5	mAb 11 Prep 2016...	mAb 11 Prep 2...	A4	mAb Method
✓ 6	mAb 11 Prep 2016...	mAb 11 Prep 2...	A4	mAb Method
✓ 7	mAb 11 Ref. Std._07	mAb 11 Ref. St...	A3	mAb Method
✓ 8	mAb 11 Blank_08	mAb 11 Blank	A2	mAb Method

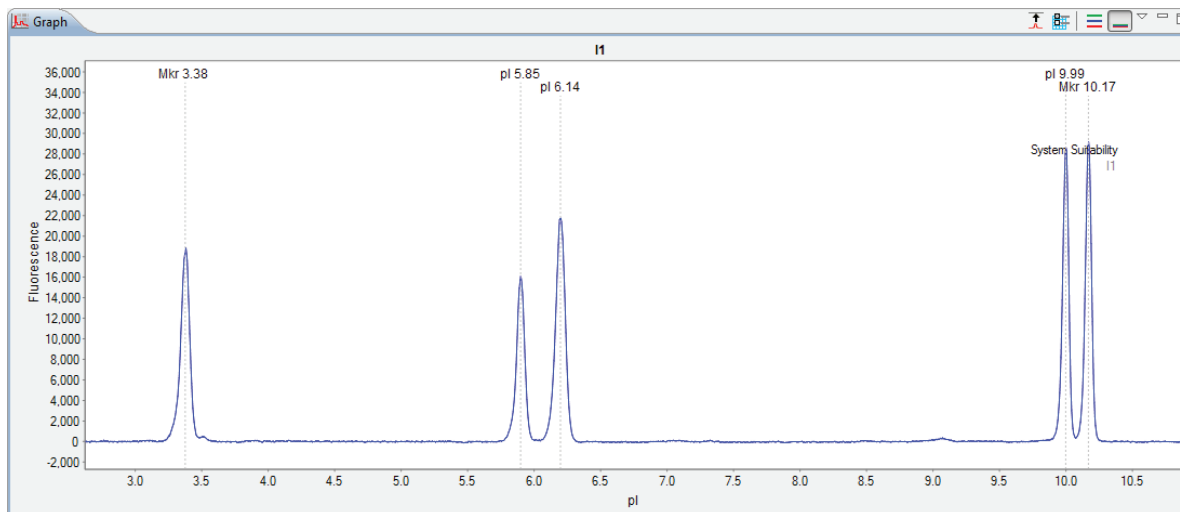
mAb Method
 Separation: 1.0 min 1500 Volts, 6.0 min 3000 Volts
 Detection: 5 Exposures
 Sample Load (s): 90
 pI Markers: 4.05, 9.99

NOTE: Data notification icons will display in the Injection column if Compass for iCE detects a potential analysis issue or data was manually modified by the user. For more information see “Data Notifications and Warnings” on page 411.

Graph Pane: Electropherogram Data

The Graph pane displays the electropherogram(s) for sample proteins or pI markers depending on the view options you’ve selected.

You can get more info on graph view options in “Changing the Electropherogram View” on page 507.



Peaks Pane: Calculated Results

The Peaks pane shows the tabulated results for your sample proteins or pI markers. Each row in the table has the individual results for each peak detected in an injection. Results shown will either be for one injection or multiple injections, samples or pI markers depending on the view options you’re using. Check out “Analysis Options Pane” on page 483 for more info.

Peaks		Injections													
Injection	Injection Name	Sample	Peak	Name	Position	pI	Height	Area	% Total	% Area	Width	S/N	Baseline	Resolution	
3	mAb 11 Ref. St...	mAb 11 Ref. St...	1	Mkr 4.05	542	4.050	8657.5	129440							
3	mAb 11 Ref. St...	mAb 11 Ref. St...	2	Peak1	1082	6.554	548.3	7585	0.5	0.5	0.0621	11.6	84.7	24.28	
3	mAb 11 Ref. St...	mAb 11 Ref. St...	3	Peak2	1107	6.668	1503.9	53212	3.4	3.4	0.1294	31.8	83.1	0.70	
3	mAb 11 Ref. St...	mAb 11 Ref. St...	4	Peak3	1139	6.811	7171.3	111929	7.1	7.1	0.0730	151.7	80.8	0.83	
3	mAb 11 Ref. St...	mAb 11 Ref. St...	5	Peak4	1158	6.881	19395.5	293503	18.6	18.6	0.0745	410.3	79.2	0.56	
3	mAb 11 Ref. St...	mAb 11 Ref. St...	6	Peak5	1177	6.968	28124.1	465924	29.5	29.5	0.0959	595.0	77.5	0.60	
3	mAb 11 Ref. St...	mAb 11 Ref. St...	7	Peak6	1199	7.066	32407.9	438220	27.7	27.7	0.0551	685.6	75.3	0.76	

NOTES:

Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

When the Markers view is selected, the information in the Peaks table includes only injection, sample, peak, position and height. pI markers the software has identified are marked with an **M**.

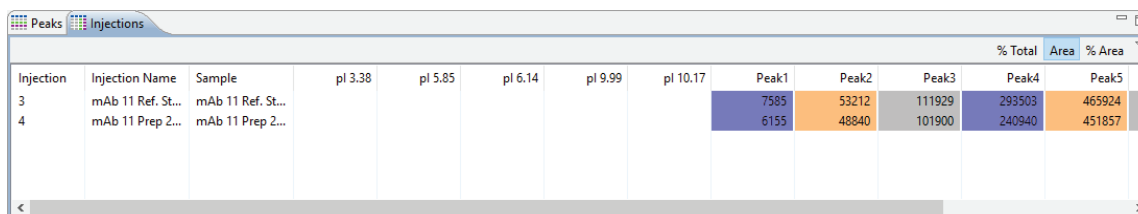
- **To view all rows** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.

The following results and info are listed in the Peaks table:

- **Injection** - Injection number.
- **Injection name** - If injection names were entered in the batch, those names will display here. Otherwise the default 'SampleID_injection number' name will display.
- **Sample** - If sample names were entered in the batch, those names will display here. Otherwise, Sample (default name) will display.
- **Peak** - Peaks are numbered in order of detection.
- **Name** - Displays peaks Compass for iCE named automatically using the user-defined peak name analysis parameters. These cells are blank if the software wasn't able to name the peak or if you didn't enter naming parameters.
- **Position** - Peak location in pixels.
- **pI** - Displays the calculated peak pI based on the migration time of the peak to the pI markers.
- **Height** - The calculated peak height.
- **Area** - Displays the count of the pixel values for dropped line fit and the area of the curve fit for gaussian fit.
- **% Total** - Displays the peak area ratio compared to the sum of all peak areas. This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.
- **% Area** - Displays the calculated percent area for the named peak compared to all named peaks. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100 (shown for named peak sample data only).
- **Width** - Displays the calculated peak width (sample data only).
- **S/N** - Displays the calculated signal to noise. The signal to noise ratio is reported for each peak based on the noise in the injection and the height of the peak. $S/N = 2 \times \text{peak height} / \text{noise}$. For the injection noise calculation, the peak to peak noise is calculated by fitting a quadratic to the test region and taking the difference between the max signal above the fit, and the minimum signal below the fit. A fit is used to mitigate the effect of drift in the signal that would give a falsely high noise value. The software looks for the quietest 50 point region in the entire injection excluding 50 points at each end. The test region is approximately 5 times a typical peak width (FWHM).
- **Baseline** - Displays the raw baseline signal of each peak.
- **Resolution** - Displays resolution of the peak compared to neighboring peaks. Two peaks that are baseline resolved will have a resolution value of 1.5. Smaller values means the peaks are not completely resolved, larger values mean the peaks are fully resolved.

Injections Pane: User-Specified Peak Names

The Injections pane shows tabulated results for sample proteins Compass for iCE labels automatically using user-defined peak name settings. Each row in the table shows the individual results for the named peaks detected in each injection.



Injection	Injection Name	Sample	pI 3.38	pI 5.85	pI 6.14	pI 9.99	pI 10.17	Peak1	Peak2	Peak3	Peak4	Peak5
3	mAb 11 Ref. St...	mAb 11 Ref. St...						7585	53212	111929	293503	465924
4	mAb 11 Prep 2...	mAb 11 Prep 2...						6155	48840	101900	240940	451857

NOTES:

Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

When the Markers view is selected, the information in the Injections table includes only injection, sample and the positions of the pI marker (Mkr) peaks.

- **To view all rows** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.

The following results and info are listed in the Injections table:

- **Injection** - Injection number.
- **Injection name** - If injection names were entered in the batch, those names will display here. Otherwise the default 'SampleID_injection number' name will display.
- **Sample** - If sample names were entered in the batch, those names will display here. Otherwise, Sample (default name) will display.
- **Peak Name Columns** - An individual column per peak name will display for every peak identified by name or as a pI marker peak in the run data. Cells for injections in these columns will be blank if Compass for iCE didn't find peaks automatically using the user-defined peak name analysis and maker parameters (or none were entered).
 - **To view peak area in the peak name columns (default)** - Select **Area** in the upper right corner of the pane. This displays calculated peak area for the individual peak only.
 - **To view % total in the peak name columns** - This displays the calculated percent area for the named peak compared to all named peaks. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100.

NOTE: The sum of the named peak percentages can be less than 100% if some peaks aren't named.

Peaks		Injections										% Total	Area	% Area
Injection	Injection Name	Sample	pI 3.38	pI 5.85	pI 6.14	pI 9.99	pI 10.17	Peak1	Peak2	Peak3	Peak4	Peak5		
3	mAb 11 Ref. St...	mAb 11 Ref. St...						0.5	3.4	7.1	18.6	29.5		
4	mAb 11 Prep 2...	mAb 11 Prep 2...						0.4	3.4	7.2	17.0	31.8		

- **To view % area in the peak name columns** - This displays the peak area ratio compared to the sum of all named peak areas. This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.

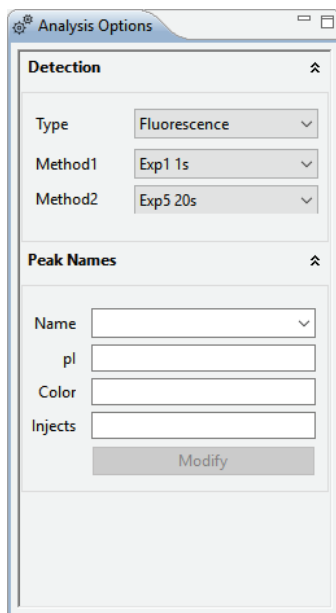
Peaks		Injections										% Total	Area	% Area
Injection	Injection Name	Sample	pI 3.38	pI 5.85	pI 6.14	pI 9.99	pI 10.17	Peak1	Peak2	Peak3	Peak4	Peak5		
3	mAb 11 Ref. St...	mAb 11 Ref. St...						0.5	3.4	7.1	18.6	29.5		
4	mAb 11 Prep 2...	mAb 11 Prep 2...						0.4	3.4	7.2	17.0	31.8		

Analysis Options Pane

The Analysis Options pane gives you a quick way to view data for different exposures and add peak names without having to open and edit the run's analysis settings.

- **Detection** - Lets you choose to view absorbance or fluorescence data for the run and select different fluorescence exposures. For more details see "Switching Between Absorbance and Fluorescence Exposures (For Standard cIEF Runs only)" on page 490.

- **Peak names** - Lets you view, change and add new custom peak name settings for sample proteins. You can apply peak name settings to all, selected, or specified injections. For details on how to use this option, go to “Naming Peaks” on page 493.



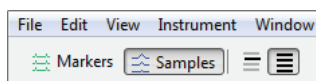
Viewing Run Data

The Analysis screen lets you view data for just one injection, specific injections or all injections in the run. Each run file has data for the sample peaks and the pI markers detected in each injection.

Switching Between Samples and Markers Data Views

Here's how you switch between viewing data for your samples and pI markers:

- **To view sample data** - Click **Samples** in the View bar or select **View** in the main menu and click **Samples**.



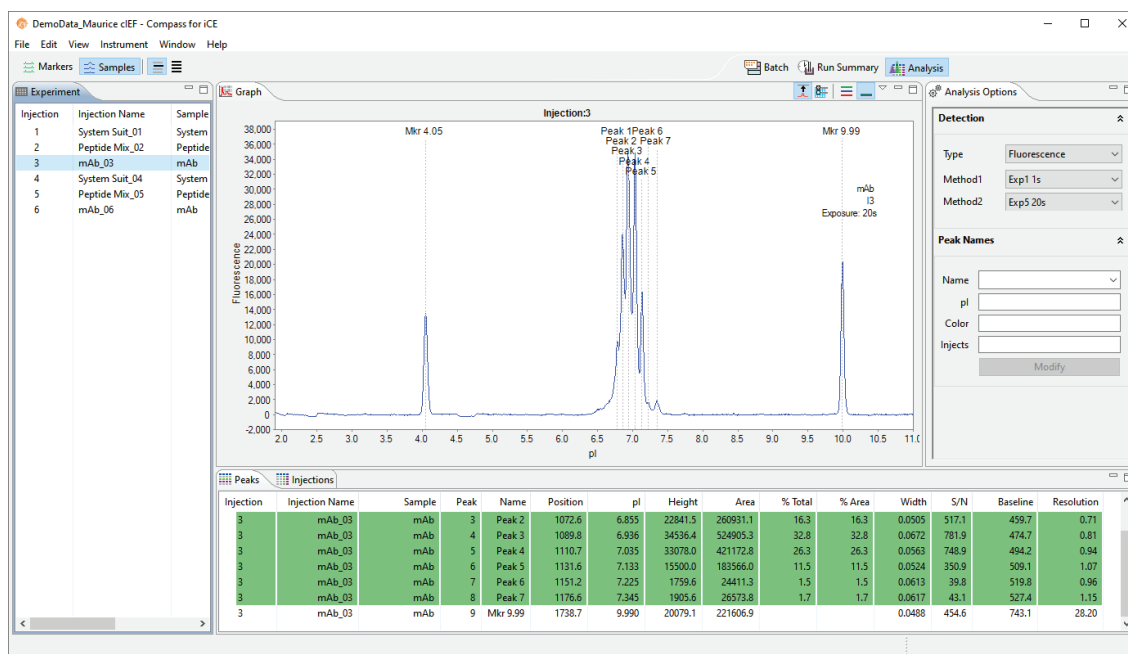
- Data in this view is for sample peaks only.
- The graph displays electropherograms with a y-axis of either Absorbance units (mAU) or Fluorescence units and an x-axis of pI. Go to “Peak Fit Analysis Settings” on page 539 for more info on how to change the detection method to view absorbance, native fluorescence or FL458 fluorescence data.

NOTES:

The FL458 nm fluorescence filter is only available on Maurice, Maurice C. and MauriceFlex systems with the option installed. One MauriceFlex systems, the FL458 nm fluorescence filter is only available for standard cIEF batches.

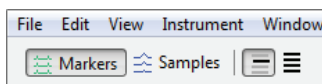
The FL458 nm filter enables detection of fluorescence emission at a longer wavelength to analyze molecules other than proteins, such as small molecule drugs in antibody-drug conjugates (ADCs). When used, ratiometric analysis of fluorescence and absorbance data can be applied to support applications including drug-antibody ratio (DAR) analysis.

- Results for each peak are shown in the Peaks and Injections panes.

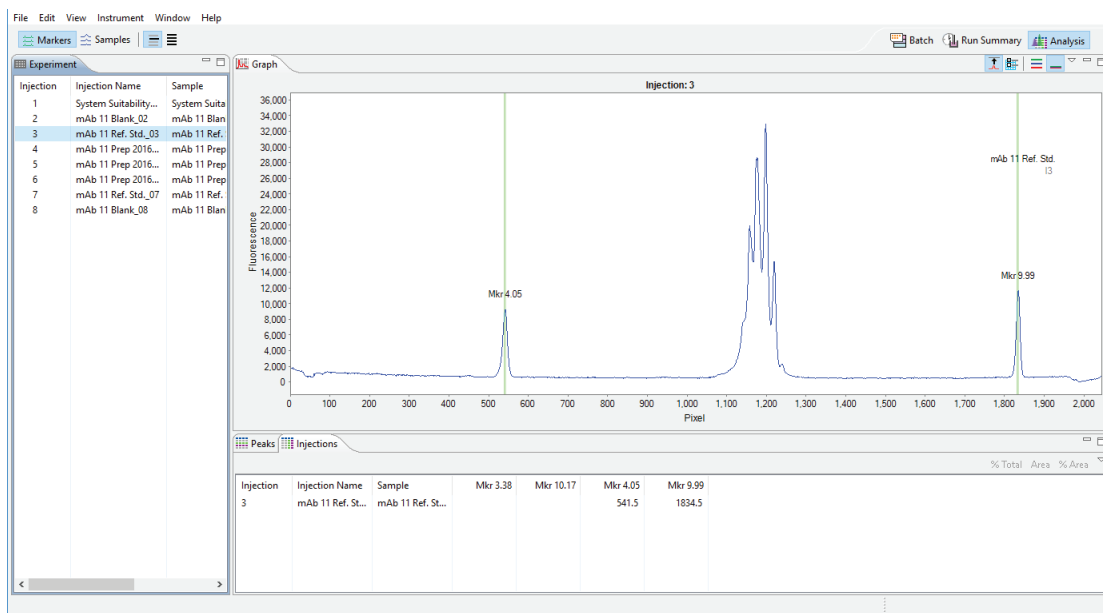


For information on checking and identifying sample peaks, see “Checking Your Data” on page 123 for standard cIEF runs and page 184 for MauriceFlex cIEF runs.

- To view pI marker data - Click **Markers** in the View bar or select **View** in the main menu and click **Markers**.



- Data in this view is for analyzing pI markers only. These are the pI markers you add to your samples during prep.
- The graph displays electropherograms with a y-axis of either Absorbance units (mAU) or Fluorescence units and an x-axis of pixels. Go to “Peak Fit Analysis Settings” on page 539 for more info on how to change the detection method to view either absorbance or native fluorescence data.
- pI markers are identified in the Peaks pane with an **M** and as Mkr in the Injections pane.

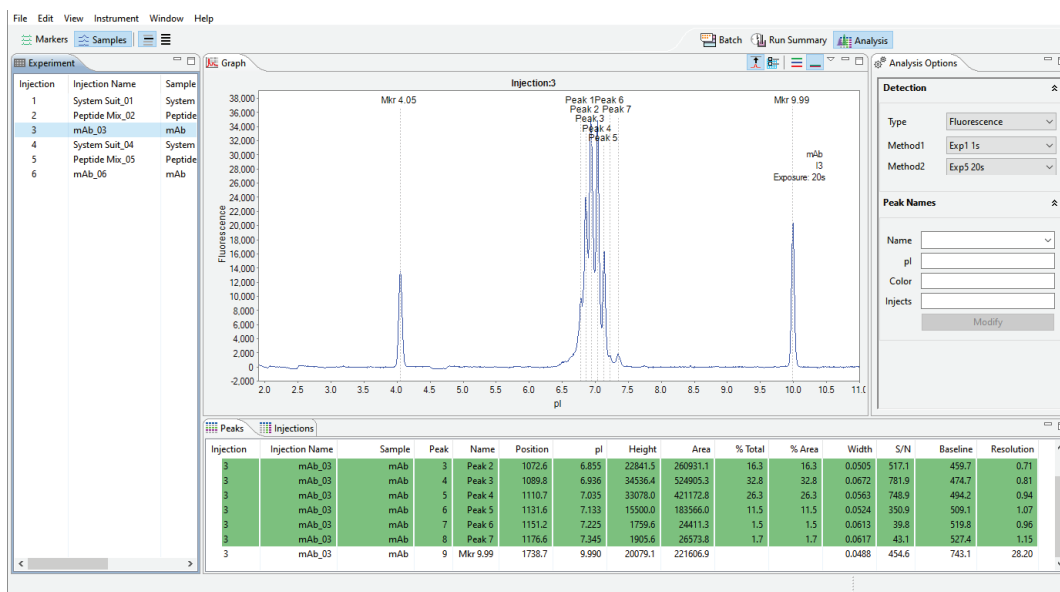


For information on checking and identifying the pI marker peaks, see “Checking Your Data” on page 123 for standard cIEF runs and page 184 for MauriceFlex cIEF runs.

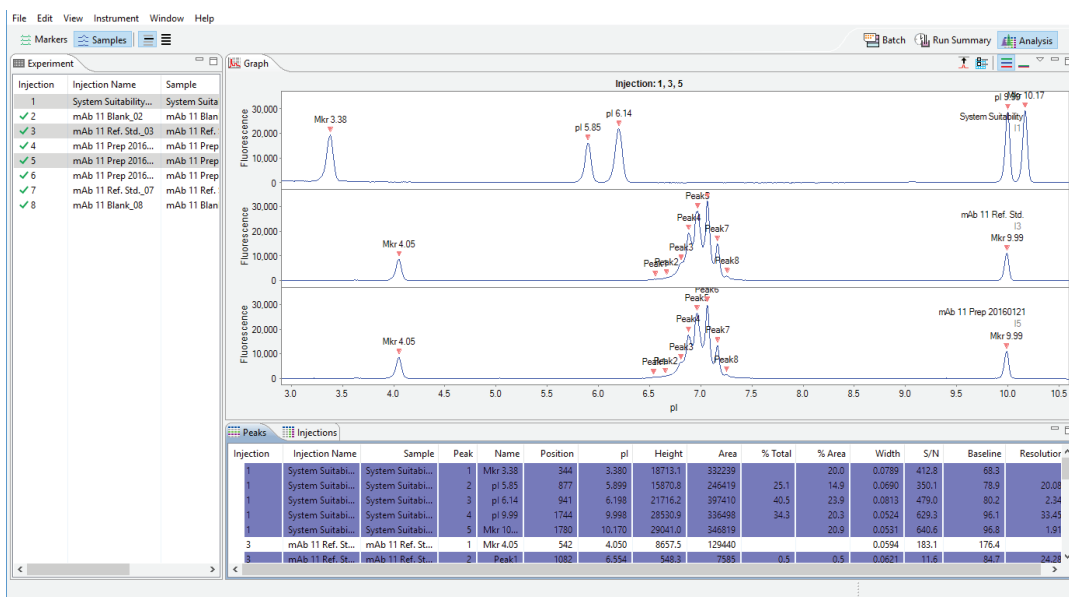
Selecting and Displaying Injection Data

You can view data from one, multiple, or all injections at once.

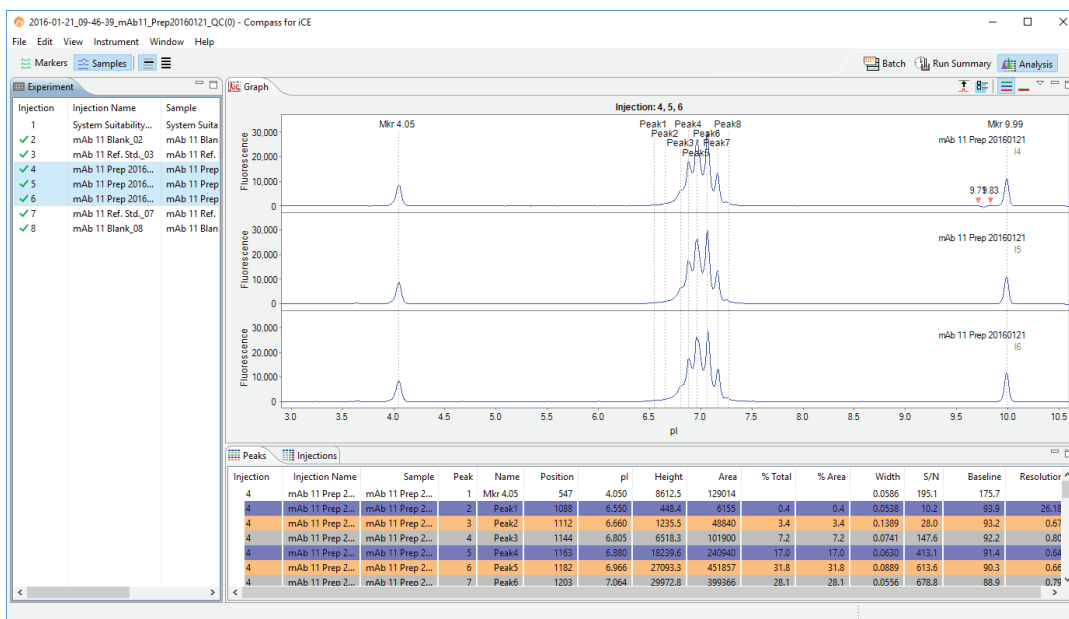
- **To look at data for one injection** - Click an injection row in the Experiment pane. Data for just that injection displays in the graph and tables.



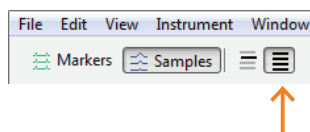
- **To look at data for specific injections** - Hold the **Ctrl** key and select just the injection rows you want to view in the Experiment pane. Data for only the injections selected display in the graph and tables.



- **To look at data for sequential injections** - Select the first injection row in the Experiment pane that you want to view, then hold the **Shift** key and select the last. This selects all rows between the two injections. Data for only the injections selected display in the graph and tables.



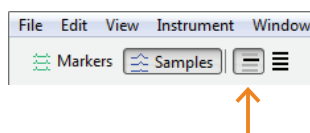
- To look at data for all injections - Just click **View All** in the View bar. Data for all injections displays in the graph and tables.



Switching Between Selected and All Views of Injections

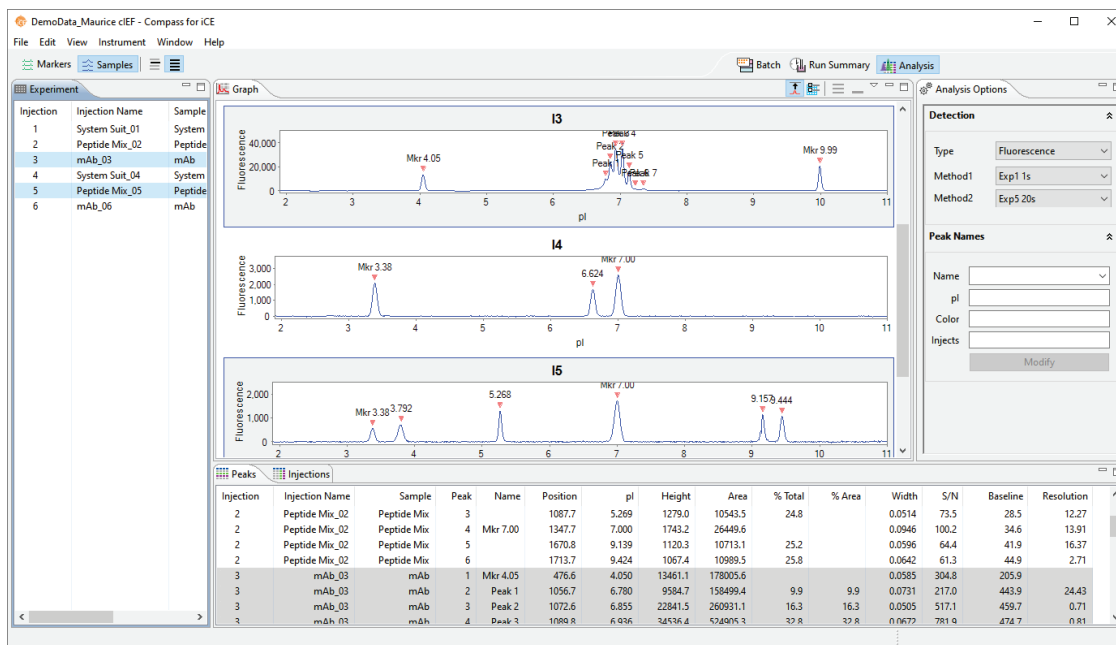
You can switch between displaying run data in a selected, per-injection format or all injections in a multi-injection format.

- To view data per in a per-injection format - Click **Selected View** in the View bar or select **View** in the main menu and click **Selected**.

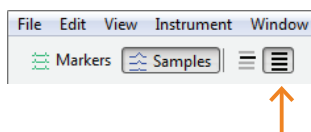


Data for the injection row(s) selected in the Experiment pane:

- Displays with electropherograms either overlaid or stacked in the Graph pane depending on the option you've got chosen.
- Shows only results for the selected row(s) in the Peaks and Injections panes.

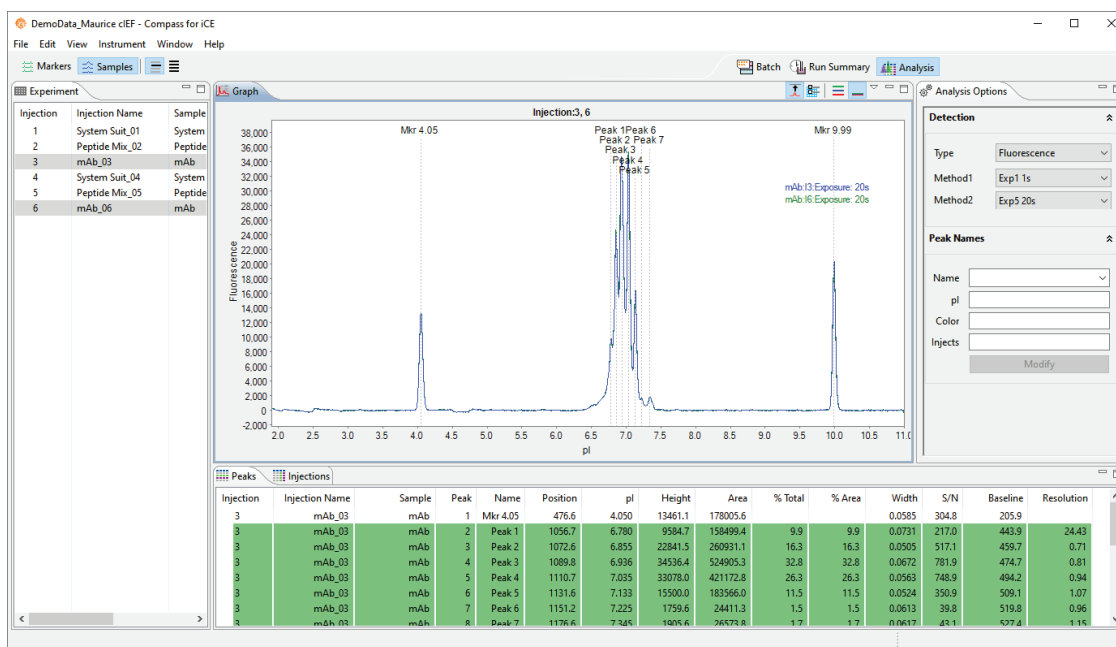


- To view data in a multi-injection format - Click **View All** in the View bar or select **View** in the main menu and click **All**:



Data for the injection row(s) :

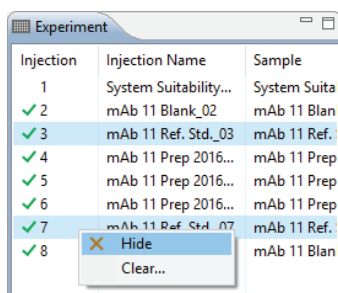
- Displays with the injections in the Graph pane in stacked view.
- Graph pane will scroll to the injection when selected in the Experiment pane.
- Shows the results for injections in the Peaks and Injections panes.



Hiding Injection Data

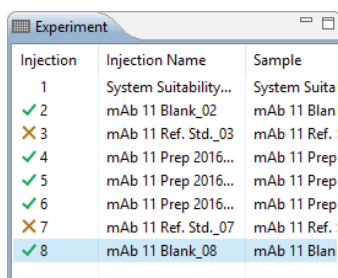
You can hide injection data from the view if needed.

- **To hide injections** - Select the injection rows you want to hide in the Experiment pane, then right click one and select **Hide**.



Data for the injections will be hidden in all data views and results tables.

- **To view hidden injections** - Select **View** in the main menu and click **Show Hidden**. Hidden rows will become visible again in all panes, and are marked with an X in the Experiment pane.



- **To unhide injections** - Select the hidden row(s). Right click on one and click **Unhide**.

Switching Between Absorbance and Fluorescence Exposures (For Standard cIEF Runs only)

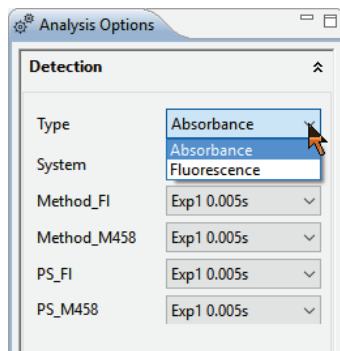
You can choose to display either absorbance or fluorescence data for your standard cIEF run in the Analysis Options pane. Changing detection methods is not an option for MauriceFlex cIEF runs.

NOTES:

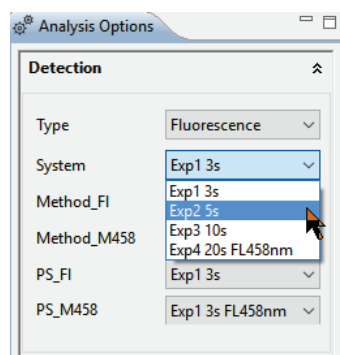
You can also change the detection method and exposure in the analysis settings for the run. See "Detection Settings" on page 529 for more information.

Detection options in the Analysis pane are disabled during a run and when Compass for iCE is reanalyzing data.

- **To change the detection method** - Click the down arrow in the **Type** field and select **Absorbance** or **Fluorescence**.



- **To change the exposure used for the sample data displayed** - Click the down arrow in one of the exposure fields and select an exposure setting.



NOTES:

You'll only be able to choose exposures for the detection method currently selected.

The number of exposures taken and exposure times shown are specified in the method when you set up your batch. They can't be changed after the run has executed.

The Absorbance exposure at 0.005 seconds is an instrument default exposure setting for standard cIEF runs. No other absorbance exposures are available.

The FL458 nm fluorescence filter is only available on Maurice and Maurice C. systems with the option installed. It is only available on MauriceFlex systems with the option installed when running a batch with a cIEF cartridge.

Data Notifications and Warnings

If Compass for iCE detects a potential data issue, a notification or warning icon will display next to the injection row in the Experiment pane.



Manual correction of sample data notification - This means the sample data was manually changed by a user, for example to add or remove a sample peak. Roll your mouse over the icon to display the type of modification that was made.

	System Sensitivity		System Sensitivity
✓ 2	mAb 11 Blank	A2	mAb Method
✓ 3	mAb 11 Ref. Std.	A3	mAb Method
✓ 4	mAb 11 Prep 20160121	A4	mAb Method
✓ 5	Peak Fit Manual rep 20160121	A4	mAb Method



Markers warning - This means one or more of the pI markers may not be identified properly. You can fix this by manually identifying the pI marker using the steps in “Step 2: Check Your pI Markers” on page 124. for standard cIEF runs and page 185 for MauriceFlex cIEF runs. Roll your mouse over the icon to display warning details.

⚠ 10	mAb 25	A3
⚠ 11	mAb 250	A2
⚠ 12	mAb 25	A3
⚠ 13	Markers Error: Not Found	A1
⚠ 14	SS	A1
⚠ 15	SS	A1



Manual correction of markers data notification - This means a user changed the pI marker data manually. Roll your mouse over the icon to display the type of modification that was made.

Injection	Sample	Location	Method
✓ 1	mAb11 Sample 1	A1	Method1
2	mAb11 Sample 2	A2	Method1
3	Markers Manual Sample 3	A3	Method1



Peak fit warning - Means that a peak can't be fit properly. This can sometimes be caused when a broad peak is fitted as multiple narrow peaks. Changing the peak width can help in this case. The warning is also caused by very small peaks around main peaks, or small peaks that are close to the end of the separation range. You can often fix this by removing the peak(s) using the steps in “Step 3: Checking Sample Peaks” on page 126 for standard cIEF runs and page 187 for MauriceFlex cIEF runs. Roll your mouse over the icon to display warning details.

⚠ 6	mAb 25	A3
⚠ 7	mAb 250	A2
⚠ 8	Peak Fit Warning: Too many iterations	
9	mAb 250	A2

Checking Your Results

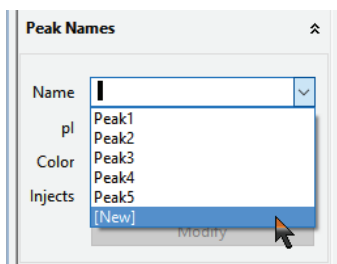
If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues. Compass for iCE detects your sample protein and pI marker peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. Please see the step by step procedure in “Checking Your Data” on page 123 for standard cIEF runs and page 184 for MauriceFlex cIEF runs to do this. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

Naming Peaks

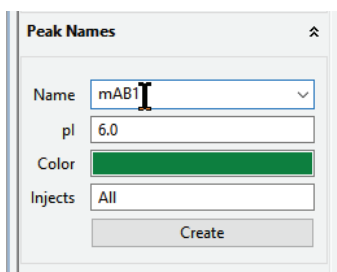
NOTE: Analysis screen options will let you add a new peak name, its pI, identifying color for peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set pI ranges, you can do that in the run analysis settings. For more info go to “Peak Names Settings” on page 532.

Adding New Peak Names in the Analysis Options Pane

1. Click the down arrow in the Name field and select New.



2. Type a name.



- Click in the **pI** field and enter the pI value of your sample protein.

Peak Names

Name: mAB1

pI: 6.78

Color: [Green Bar]

Injects: All

Create

- Click on the **Color** field to display the color selection box.

Peak Names

Name: mAB1

pI: 6.78

Color: [Green Bar]

Injects: All

Create

Color

Basic colors:

Custom colors:

Define Custom Colors >>

OK Cancel

- The color you pick is used to identify the sample protein peak in the Peaks and Injections panes in the Analysis screen. Click a color or define a custom color and click **OK**. The color selection will update in the field:

Peak Names

Name: mAB1

pI: 6.78

Color: [Red Bar]

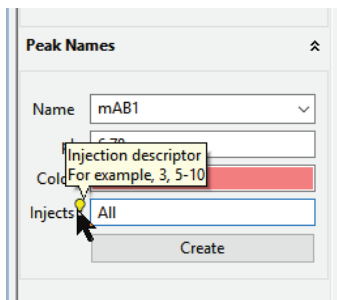
Injects: All

Create

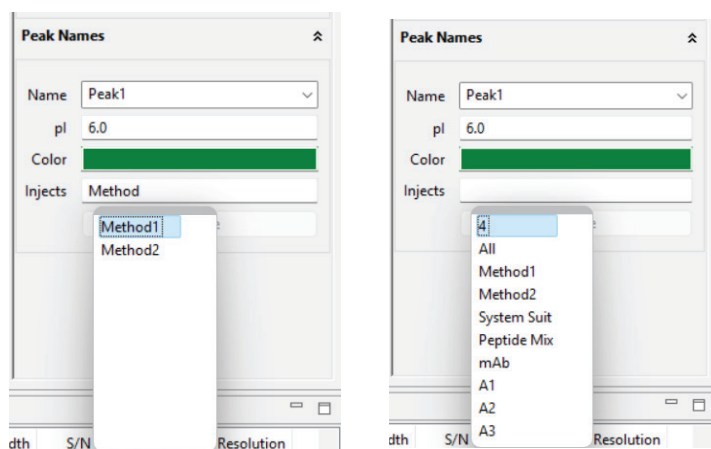
- Click in the **Injects** field:

To apply the new peak name to all injections - leave the default setting of All.

To apply the new peak name to specific injections - enter individual injections separated by commas (for example 2, 5, 8), a range of injections (for example 3-15) or a combination of both (for example 8, 10, 12-15). You can also click the tool tip for more examples.



To apply the new peak name to specific run data - type the first letter of a method, sample name, well or vial location. Compass for iCE will search the run data and display a list of matching options for you to choose from. Alternatively you can highlight the text in the Injections cell, hit **Delete**, then select an option from the drop down list.

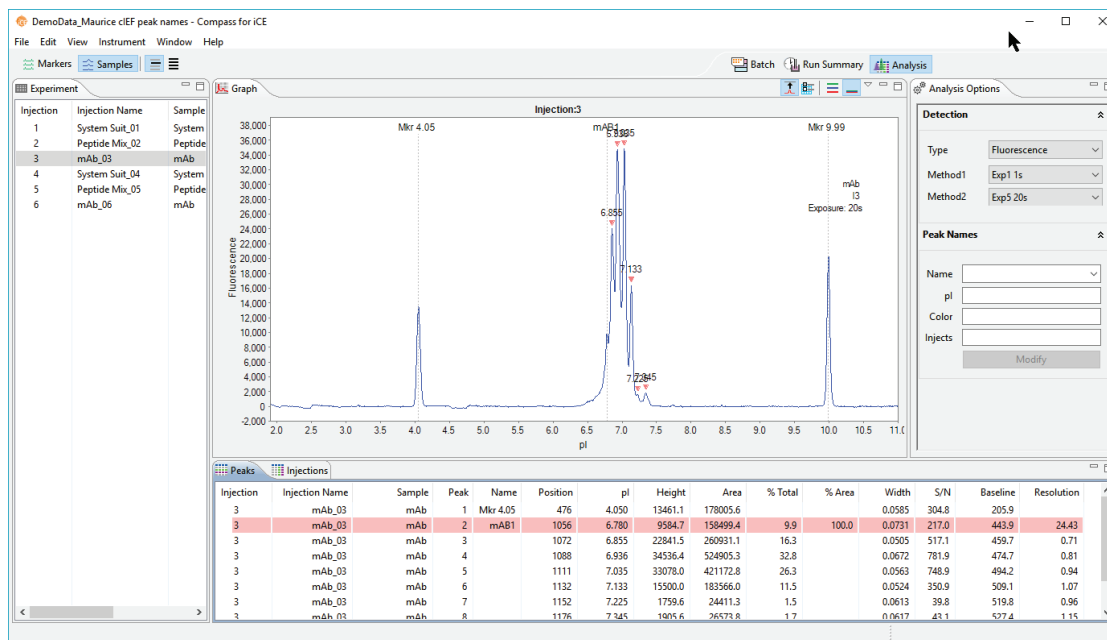


- **Injection** - Applies the peak name to the injection.
- **All** - Applies the peak name to all injections.
- **Methods** - All methods in the run file display in the list. Selecting a method applies the peak name to all injections that used that method.
- **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the peak name to all injections that used that sample name.
- **Wells or vials** - The well or vial numbers currently selected will display in the list. Selecting a well or vial number applies the peak name to all injections that used that well or vial.

7. Click **Create** to save changes.

NOTE: To clear the new peak name information you've entered without saving changes, just delete the peak name in the Name field.

The named peak will be identified with a peak name label in the electropherogram and color-coded in the Peaks and Injections panes:



Adding Peak Names from the Graph or Peaks Table

NOTE: Analysis screen options will let you change an existing peak name, its pI, identifying color in peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set pI ranges, you can do that in the run analysis settings. For more info go to “Manual Peak Integration” on page 547.

1. Right click the peak you want to name in the Graph or Peaks pane.

2. Select Name Peak.

The screenshot shows a chromatogram with a peak labeled 9.424. A context menu is open over the peak, with 'Name Peak' selected. To the right, a 'Peak Names' panel shows fields for Name, pI, and Color. Below the chromatogram is a table with columns: Injection, Injection Name, Sample, Peak, Name, Position, pI, Height, and Resolution.

Injection	Injection Name	Sample	Peak	Name	Position	pI	Height	Resolution
2	Peptide Mix_02	Peptide Mix	1	Mkr 3.38	803	3.380	561.5	
2	Peptide Mix_02	Peptide Mix	2	Peak2	864	3.794	713.6	10
2	Peptide Mix_02	Peptide Mix	3	Peak3	1088	5.269	1279.0	10
2	Peptide Mix_02	Peptide Mix	4	Mkr 7.00	1348	7.000	1743.2	20
2	Peptide Mix_02	Peptide Mix	5	Peak4	1671	9.139	1120.3	10
2	Peptide Mix_02	Peptide Mix	6		1711	9.424	1067.4	10

3. To use an existing peak name - select a name from the list.

To create a new peak name - select New. Type in a name for the peak. Click All to apply to all injections or Selected to apply only to the injections selected.

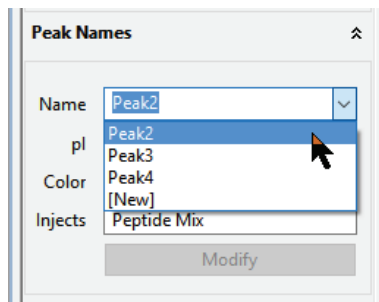
The 'Add New Peak Name' dialog box has a title bar with a close button. It contains a 'Peak Name' text field with 'Peak 5' entered. Below it are two radio buttons: 'All' (selected) and 'Selected'. At the bottom are 'OK' and 'Cancel' buttons.

4. Click OK. If you want to add an identifying color or change the injections the new peak name applies to, you can do that in the Analysis Options pane.

Changing Peak Name Information

NOTE: Analysis screen options will let you change an existing peak name, its pI, identifying color for peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set pI ranges, you can do that in the run analysis settings. For more info go to “Manual Peak Integration” on page 547.

1. In the Analysis Options pane, click the down arrow in the **Name** field and select an existing peak name.



2. Change the name, pI, color and injects as needed then click **Modify**.

Group Statistics

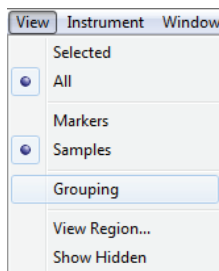
You can use the Grouping view to have Compass for iCE do a statistical analysis of named proteins in your injections (see “Manual Peak Integration” on page 547. for more info on setting named peaks up). Statistics for each protein are also plotted for easy comparison.

Using Groups

1. Groups are automatically created for injections that use the same sample name and method, so to use this feature, you need to make sure you’ve got sample names entered.
 - a. Go to the **Batch** screen.
 - b. Click the **Sample ID** cells in the Injection pane and type a name for any samples you want to calculate statistics for.

Injections History Notes					
	Sample ID	Location	Method	Notes	
1	Sample 1	A1	Method1		
2	Sample 2	A2	Method1		
3	Sample 3	A3	Method1		
4	Sample 4	A4	Method1		

- Go back to the **Analysis** screen. Click **View** in the main menu and select **Grouping**.



NOTE: To turn Grouping off, select **View** in the main menu and deselect **Grouping**.

Viewing Sample Injection Groups

Compass for iCE automatically groups all injections using the same sample name together in the Injection Groups pane.

Injection Groups	
Sample	Method
System Suitability	System Suitability
▶ mAb 11 Blank (2)	mAb Method
▶ mAb 11 Prep 20160121 (3)	mAb Method
▶ mAb 11 Ref. Std. (2)	mAb Method

- To expand a group** - Click the arrow next to a group to see the individual injections in the group and reported data for each

Injection Groups	
Sample	Method
System Suitability	System Suitability
▶ mAb 11 Blank (2)	mAb Method
▲ mAb 11 Prep 20160121 (3)	mAb Method
✓ mAb 11 Prep 20160121	mAb Method
✓ mAb 11 Prep 20160121	mAb Method
✓ mAb 11 Prep 20160121	mAb Method
▶ mAb 11 Ref. Std. (2)	mAb Method

- To expand all groups** - Click **Expand All (+)** in the upper right corner of the pane.
- To collapse all groups** - Click **Collapse All (-)** in the upper right corner of the pane.

Viewing Statistics

Peak and Method Groups

The Peak Groups pane reports statistics for each named protein in every group. Each group shows the statistics for named proteins which includes average area, standard deviation, %CV and SEM (standard error measurement). The number in parenthesis after the sample name is the number of injections in the group.

Peak Groups				Method Groups				Group Plot							
Sample	Method	Injection	Name	Area	Std.Dev.	% CV	SEM					% Total	Area	% Area	
> Peptide Mix (2)	Method1		Peak2	10292	42.81	0.4	30.27								
> Peptide Mix (2)	Method1		Peak3	10496	67.27	0.6	47.56								
> System Suit (2)	Method1		Peak1	22483	544.8	2.4	385.2								

- **To display results using area** - Click **Area** in the upper right corner of the pane.
- **To display results using % total** - Click **% Total** in the upper right corner of the pane to display the calculated percent area for the named peak compared to the total area measured in the injection. This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.
- **To display results using % area** - Click **% Area** in the upper right corner of the pane to display the calculated percent area for the named peak compared to all named peaks. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100 (shown for named peak sample data only).
- **To expand a group** - Click the arrow next to a group to see the individual injections in the group and reported data for each

Peak Groups				Method Groups				Group Plot							
Sample	Method	Injection	Name	Area	Std.Dev.	% CV	SEM					% Total	Area	% Area	
Peptide Mix (2)	Method1		Peak2	10292	42.81	0.4	30.27								
Peptide Mix	Method1	2	Peak2	10322											
Peptide Mix	Method1	5	Peak2	10261											
> Peptide Mix (2)	Method1		Peak3	10496	67.27	0.6	47.56								
> System Suit (2)	Method1		Peak1	22483	544.8	2.4	385.2								

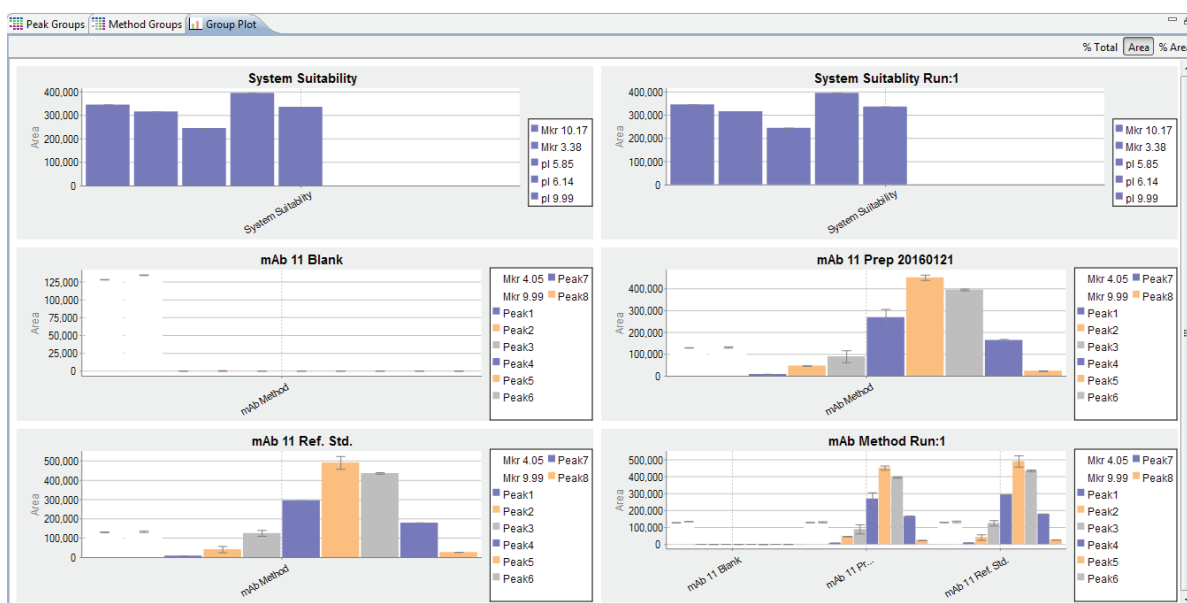
- **To expand all groups** - Click **Expand All (+)** in the upper right corner of the pane.
- **To collapse all groups** - Click **Collapse All (-)** in the upper right corner of the pane.

The Method Groups pane pivots the Peak Groups pane results to show statistics for named protein peaks in individual columns.

Sample	Method	Injection	Peak1:Area	Std.Dev.	%CV	SEM	Peak2:Area	Std.Dev.	%CV	SEM	Peak3:Area	Std.Dev.	%CV	SEM
> Peptide Mix (2)	Method1		0	0.0000	0.0	0.0000	10292	42.81	0.4	30.27	10496	67.27	0.6	47.56
> System Suit (2)	Method1		22483	544.8	2.4	385.2	0	0.0000	0.0	0.0000	0	0.0000	0.0	0.0000
> mAb (2)	Method2		0	0.0000	0.0	0.0000	0	0.0000	0.0	0.0000	0	0.0000	0.0	0.0000

Group Plots

The mean values for named peaks using the same method in each injection group are plotted in bar graphs with error bars showing the standard deviation in the Group Plots pane. You'll also get plots that compare samples using the same method in the run.



Hiding or Removing Injections in Group Analysis

Hidden injections are not included in injection groups. But, hiding injections gives you an easy way to reject individual injections from the statistical analysis. See Hiding Injection Data for details on how to do this.

Copying Results Tables and Graphs

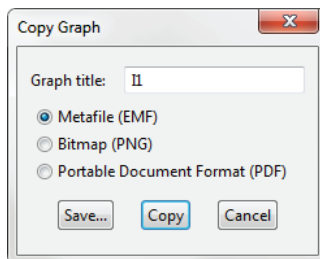
You can copy and paste data and results tables into other documents, or save the electropherogram as a graphic file.

Copying Results Tables

1. Click in the Peaks or Injections pane.
2. Select one or multiple rows.
3. Select **Edit** in the main menu and click **Copy**, or right click on row(s) you selected and click **Copy**.
4. Open a document (Microsoft® Word®, Excel®, PowerPoint®, etc.). Right click in the document and select **Paste**. Data for the rows selected will be pasted into the document.

Copying the Graph

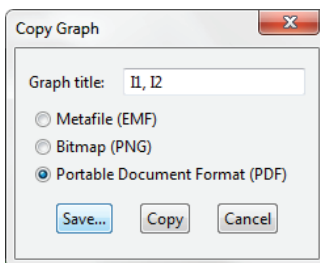
1. Select the Graph pane.
2. Select **Edit** in the main menu and click **Copy**, or right click in the Graph pane and select **Copy**.
3. Select an image option (EMF, PNG or PDF) in the pop-up window, then click **Copy**.



4. Open a document (Microsoft® Word®, Excel®, PowerPoint®, etc.). Right click in the document and select **Paste**. A graphic of the copied electropherogram will be pasted into the document.

Saving the Graph as an Image File

1. Select the Graph pane.
2. Select **Edit** in the main menu and click **Copy**, or right click in the Graph pane and select **Copy**.
3. Select an image option (EMF, PNG or PDF) in the pop-up window, then click **Save**.



4. Select a directory to save the file to, enter a file name, then click **OK**.

Exporting Run Files

Results tables and raw plot data can be exported for use in other applications. Data is organized into these folders when exported:

- Absorbance data: <run file name>Export_ABS
- Fluorescence data:
 - <run file name>Export_FL when the same exposure is selected for all methods
 - <run file name>Export_FL_FL458nm when the FL458 nm filter is selected for all methods
 - <run file name>Export_FL_Various when both the native fluorescence and FL458 nm filter are selected

Exporting Results Tables

To export the information in the Peaks and Injections tables:

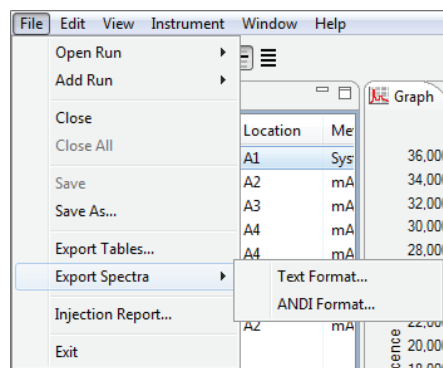
1. Click **File** in the main menu and click **Export Tables**.
2. Select a directory to save the files to and click **OK**. Data will be exported in .txt format.

NOTE: To exclude export of standards (pI markers) data or export results table data in .csv format, see “Setting Data Export Options” on page 760.

Exporting Raw Sample Electropherogram Data

To export raw sample plot and background data:

1. Click **File** in the main menu and click **Export Spectra**.



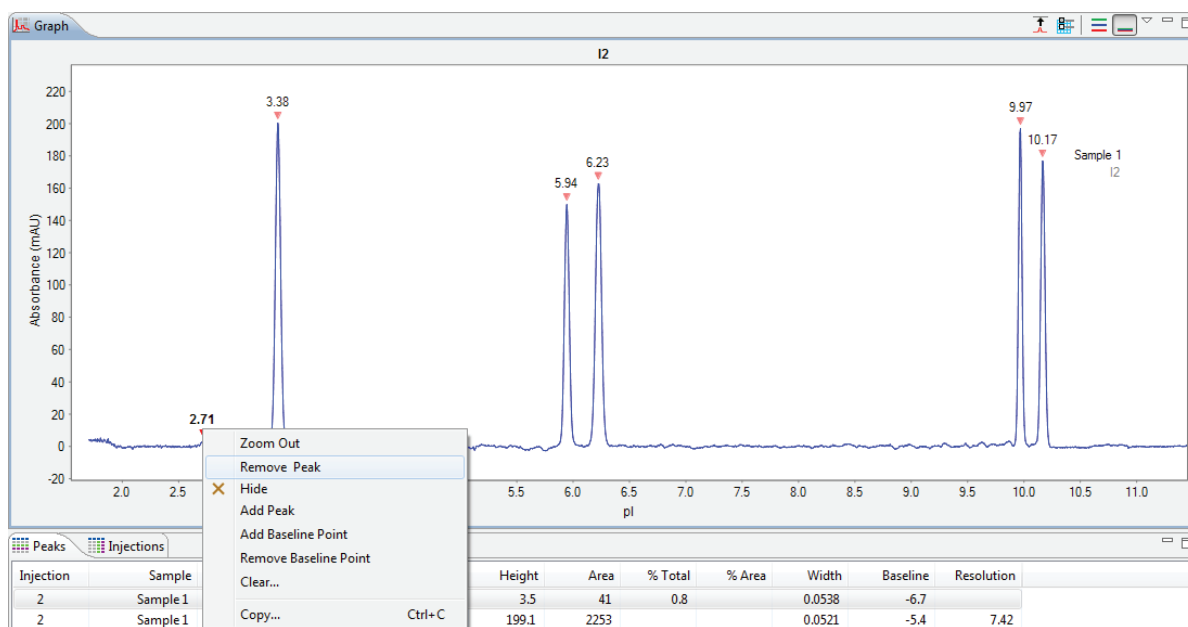
- To export data in .txt format - Select **Text Format**. Data will be exported in one file for all injections.
 - To export data in .cdf format - Select **ANDI Format**. Data will be exported in one file per injection.
2. Select a directory to save the files to and click **OK**. Data will be exported in the selected format.

Changing Sample Protein Identification

Compass for iCE lets you customize what sample proteins are reported in the results tables by making manual adjustments in the electropherogram or Peaks table.

Adding or Removing Sample Data

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that has the injection you want to correct, then click the **Graph** tab.
 - **To remove a peak from the data** - Right click the peak in the electropherogram or Peaks table and select **Remove peak**. The software will no longer identify it as a sample peak in the electropherogram, and the peak data will be removed in the results tables.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment			
Injection	Sample	Location	Method
✓ 1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1

- **To add an unidentified peak to the data** - Right click the peak in the electropherogram or peaks table and select **Add Peak**. The software will calculate and display the results for the peak in the results tables and identify the peak in the electropherogram.

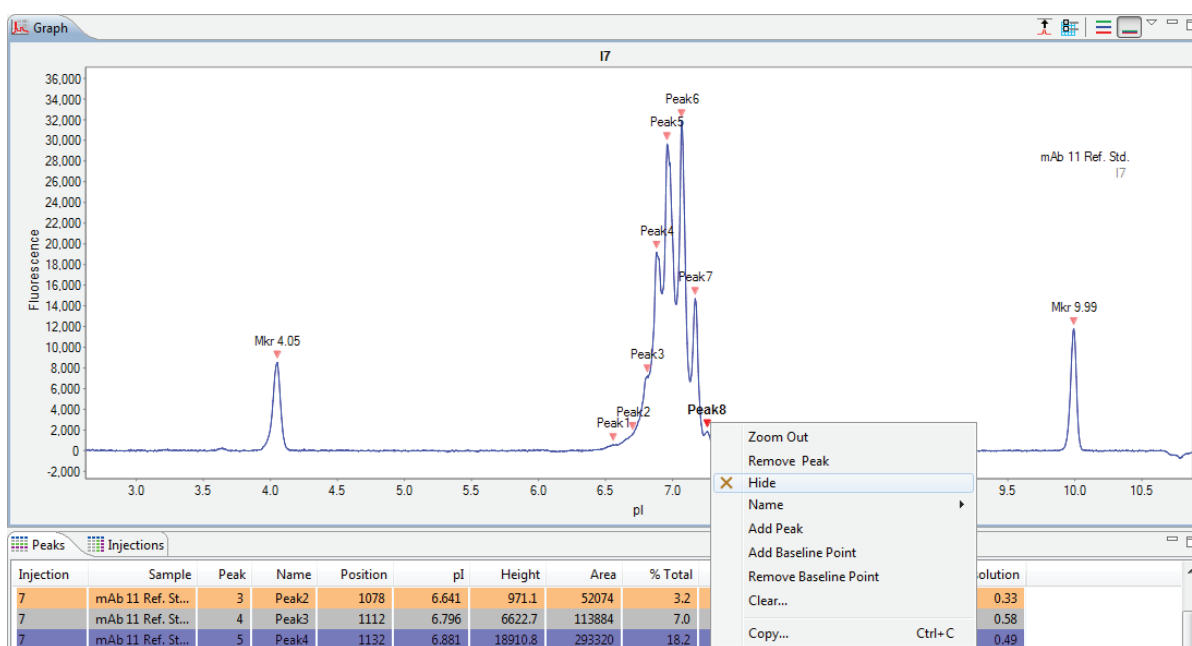
A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear** for the current injection or **Clear All** for all injections in the batch.

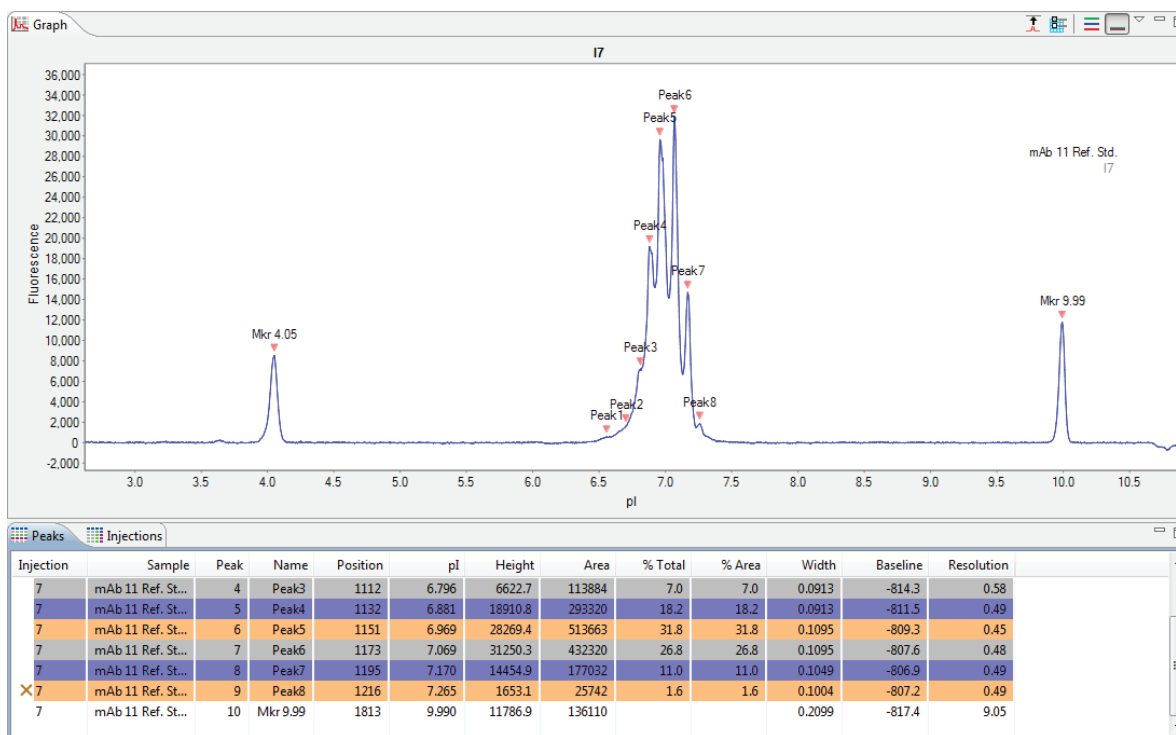
Hiding Sample Data

You can hide the results for a sample protein in the results tables without completely removing it from the reported results.

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that contains the injection you want to correct, then click the **Graph** tab.
4. Right click the peak in the electropherogram or Peaks table and select **Hide**. Compass for iCE will hide the peak data in the results tables.



5. To view hidden peak data, click **View** in the main menu and click **Show Hidden**. Hidden peak data will display in the results table and be marked with an X.



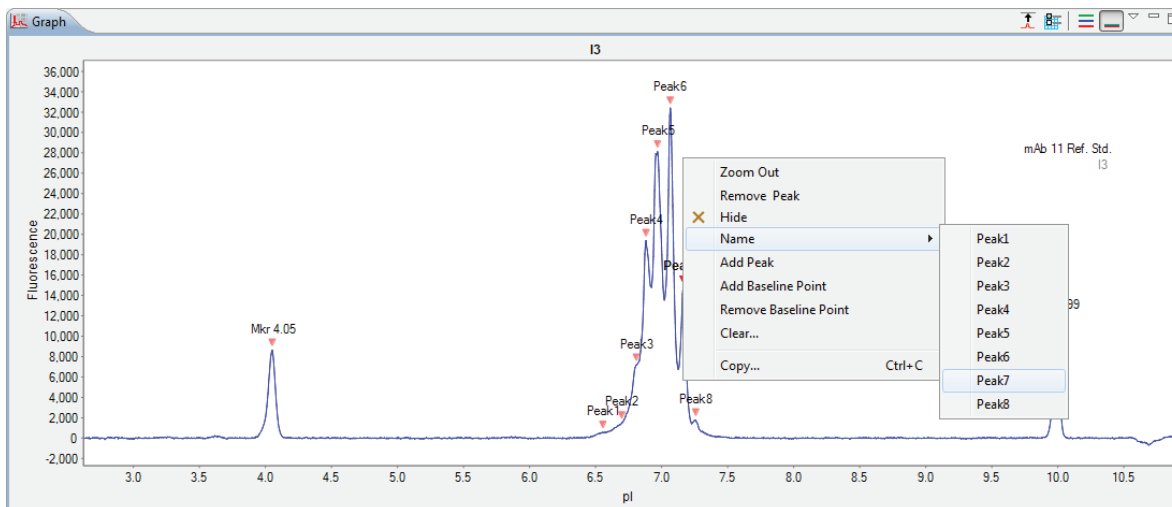
6. To unhide a peak, right click on the peak in the electropherogram or peaks table and select **Unhide**.

Changing Peak Names for Sample Data

If Compass for iCE did not automatically name a sample protein peak, you can do it manually.

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that has the sample you want to correct, then click the **Graph** pane.

- Right click the peak in the electropherogram or Peaks table and click **Name**, then select a name from the list. Compass for iCE will change the peak name in the electropherogram and results tables, and adjust peak names for other sample proteins accordingly.







NOTE: For details on how to specify peak name settings, see “Manual Peak Integration” on page 547.

Changing the Electropherogram View

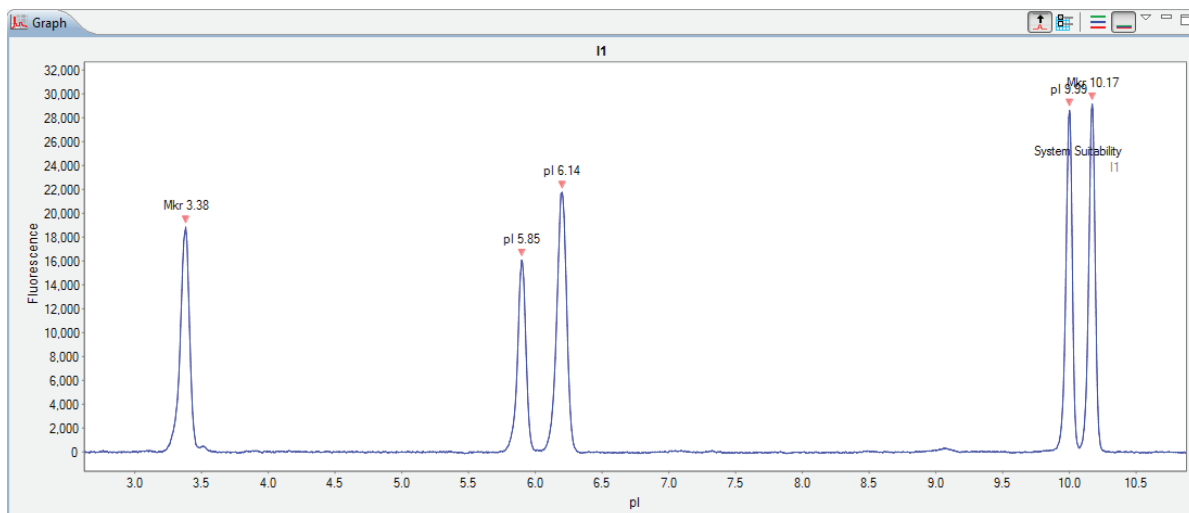
Options in the Graph pane let you zoom and rescale electropherograms, overlay or stack plots and change the peak and plot info displayed.

The Graph pane toolbar has these options:

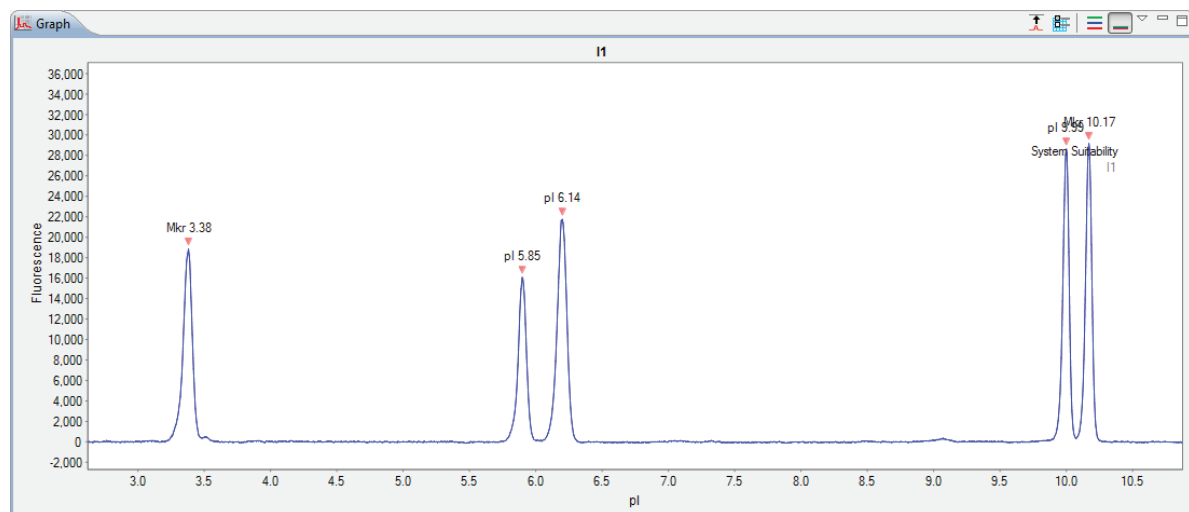
-  Auto Scale
-  Graph Options
-  Stack the Plots
-  Overlay the Plots

Autoscaling the Electropherogram

Click the **Auto Scale** button to scale the y-axis to the largest peak in the electropherogram.

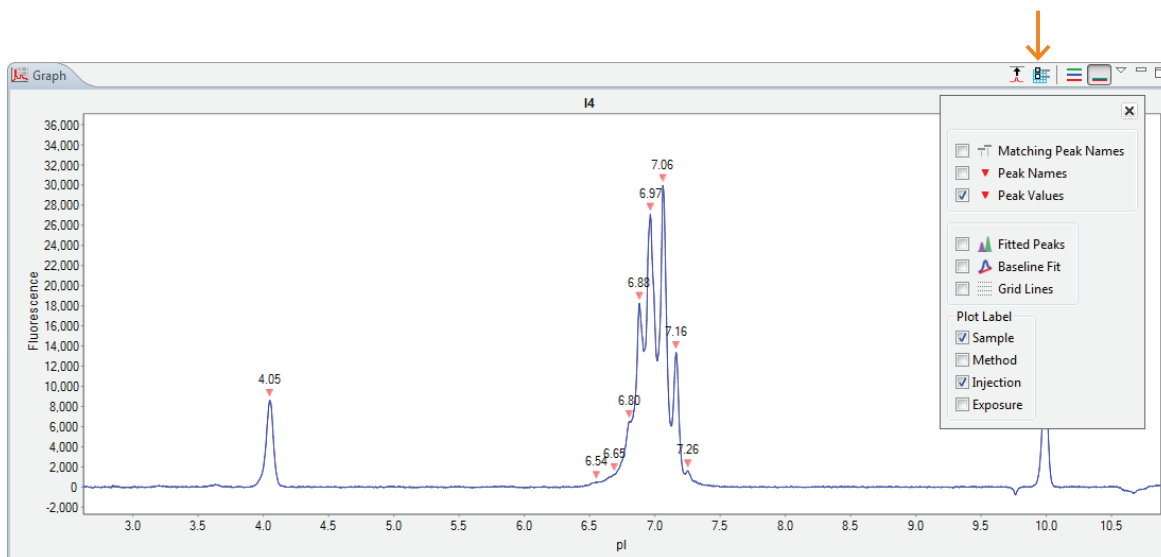


Click the **Auto Scale** button again to return to default scaling.



Customizing the Data Display

You can customize electropherogram peak labels, plot labels and display options. To do this, just select the **Graph Options** button.

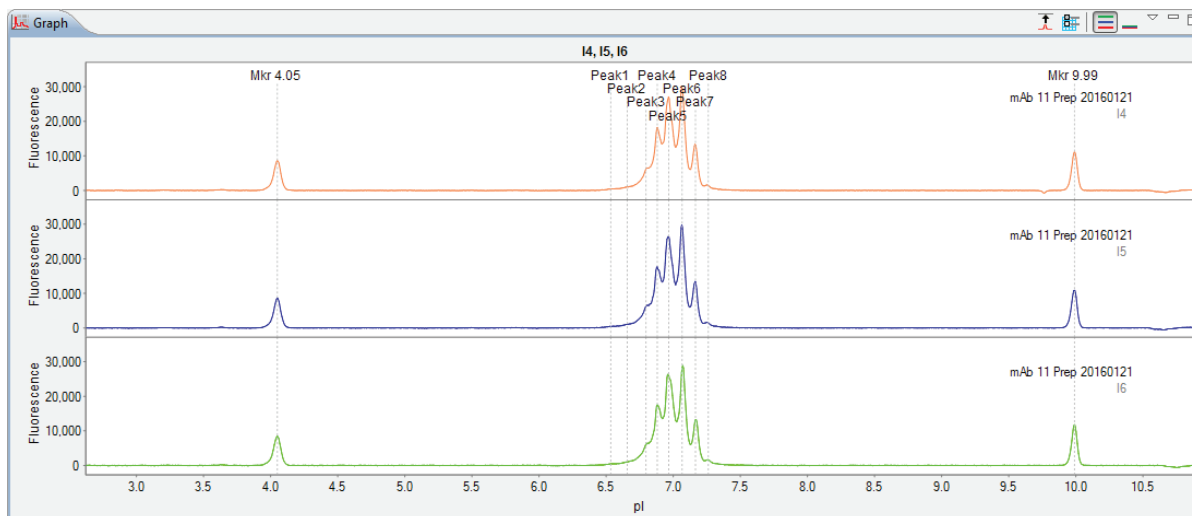


Peak Labels

You can customize the labels used to identify peaks in the electropherogram with these options:

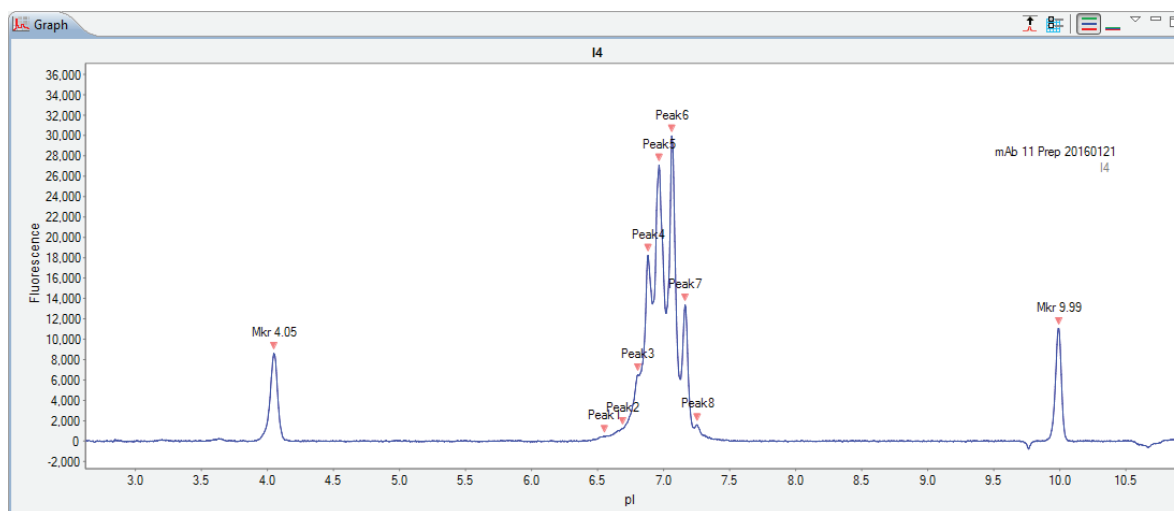
<input type="checkbox"/>	<input type="checkbox"/>	Matching Peak Names
<input type="checkbox"/>	<input checked="" type="checkbox"/>	Peak Names
<input type="checkbox"/>	<input checked="" type="checkbox"/>	Peak Values

- **Matching Peak Names** - Checking this box will draw vertical lines through each named peak. Using this option with Stack the Plots or Overlay the Plots features is helpful for visually comparing your named peaks across multiple injections.



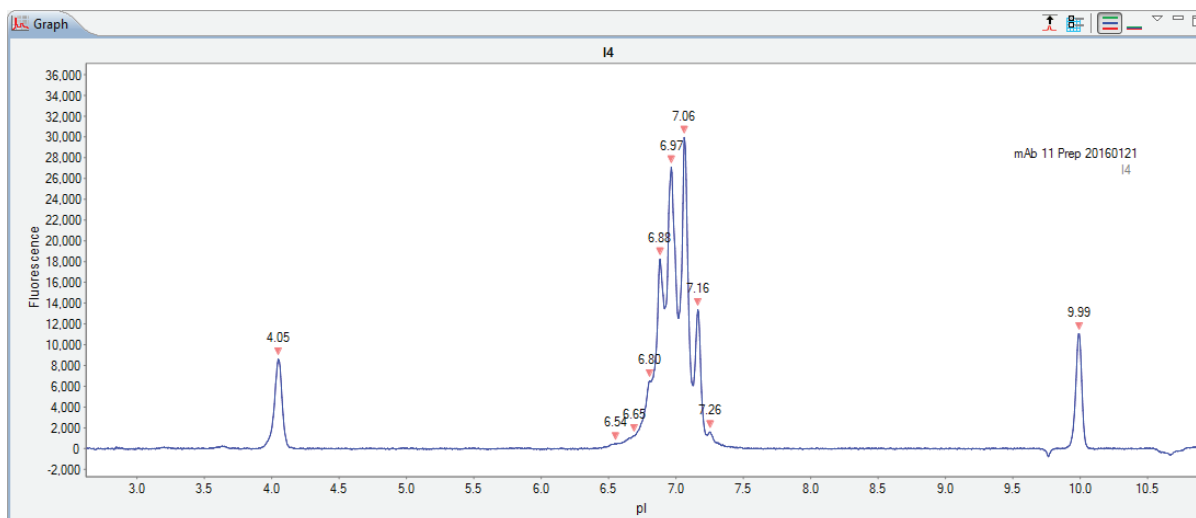
- **Peak Names** - Checking this box displays peak name labels on all named peaks in the electropherogram.

NOTE: If more than one peak label option is selected, peak name labels will always be used for named peaks.



- **Peak Values** - Checking this box will display the pI values on all peaks in the electropherogram.

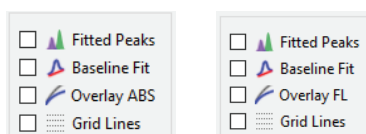
NOTE: If more than one peak label option is selected, peak name labels will always be used for named peaks.



Baseline and Grid Options

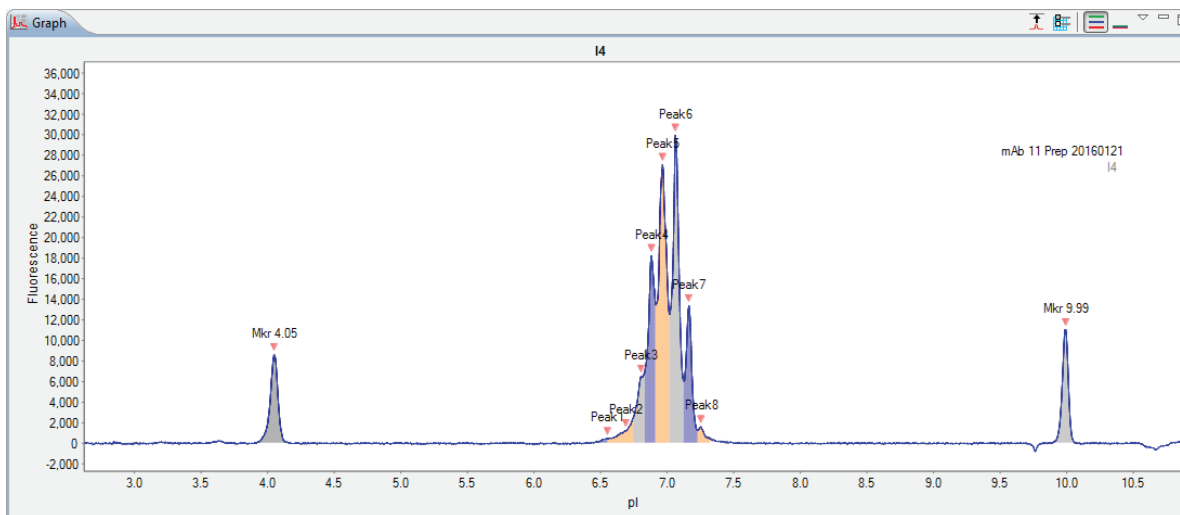
You can view the calculated baseline fit, peak integration, show grid lines and overlay fluorescence and absorbance electropherograms with these options.

NOTE: See “Overlaying Fluorescence and Absorbance Electropherograms” on page 437 for information on the detection mode data overlay.



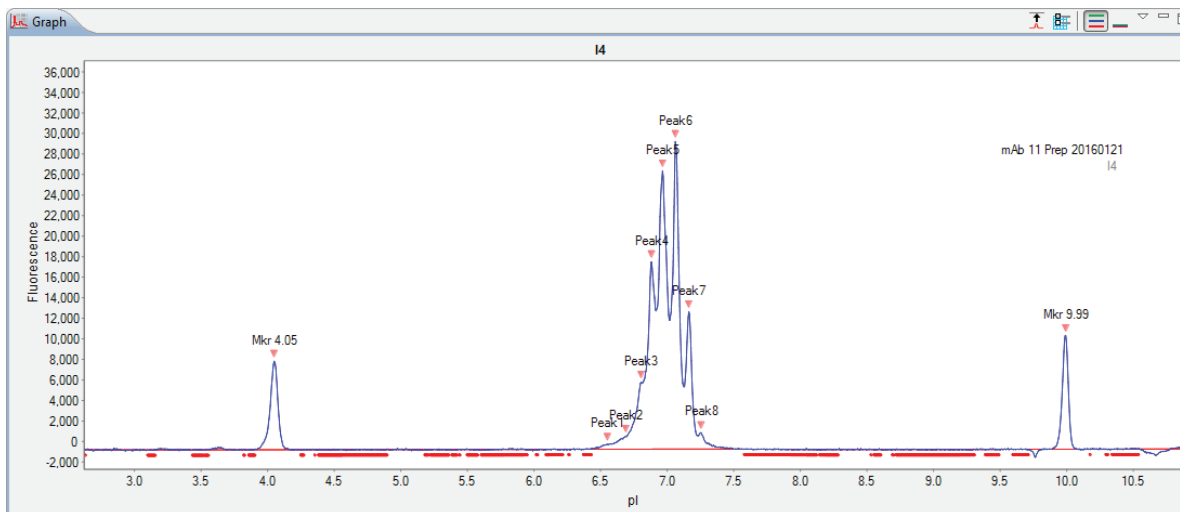
- **Fitted peaks** - Checking this box displays how the peaks were fit by the software. For cIEF runs, the software uses Gaussian Fit by default. For MauriceFlex cIEF runs, the software uses Dropped Line as the default.

NOTE: This option is only available for sample data.

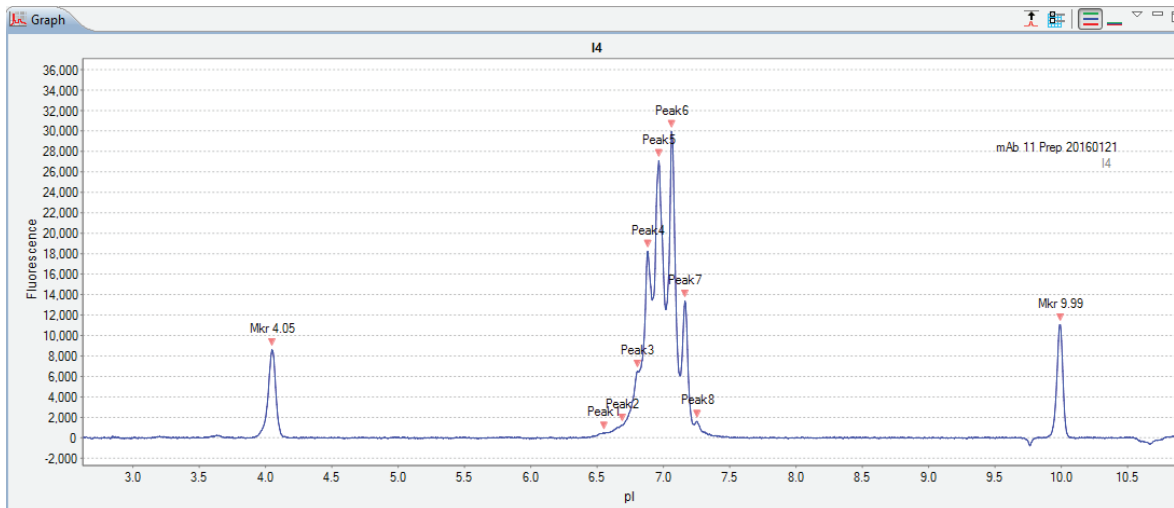


- **Baseline Fit** - Checking this box displays the calculated baseline for the peaks. Baseline points will also display for regions of the electropherogram considered to be at baseline.

NOTE: This option is only available for sample data.



- **Grid Lines** - Checking this box adds grid lines in the graph.



Plot Labels

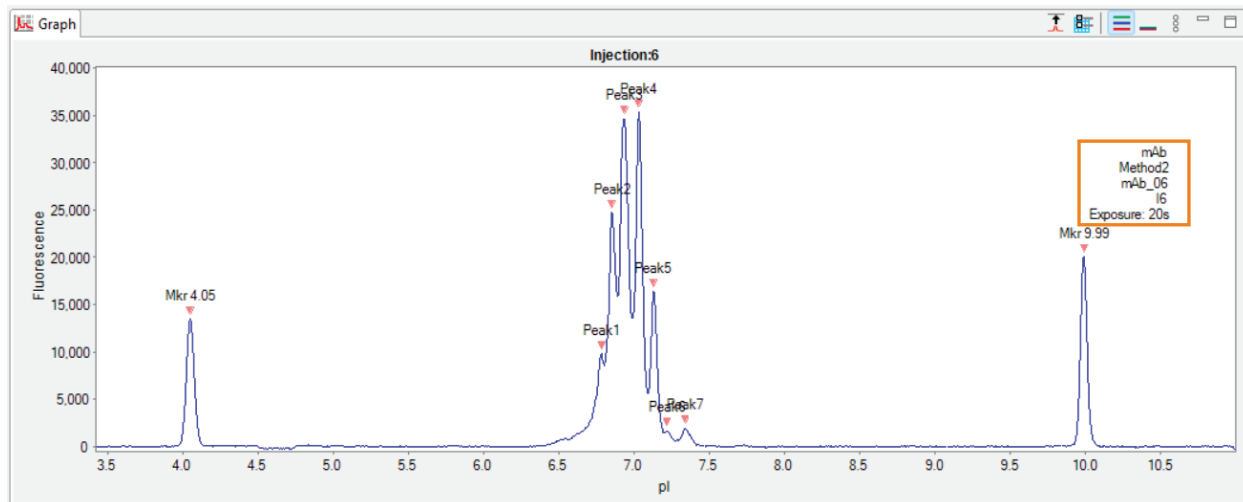
You can customize the plot labels displayed on the electropherogram with these options.

Plot Label	
<input checked="" type="checkbox"/>	Sample
<input type="checkbox"/>	Method
<input checked="" type="checkbox"/>	Injection
<input checked="" type="checkbox"/>	Exposure
<input type="checkbox"/>	Injection Name

Plot labels are shown in the upper right side of the graph.

- **Sample** - Checking this box displays the sample name used for the injection. If sample names were entered with the batch, those names will display here. If not, Sample (default name) displays.
- **Method** - Checking this box displays the method used for the injection.
- **Exposure** - Checking this box display the exposure time(s) used for the data.
- **Injection** - Checking this box displays the injection number. For example, I4 for injection 4 in the run.
- **Injection Name** - Checking this box displays the injection name used for the injection. If injection names were entered with the batch, those names will display here. If not, the default name displays.

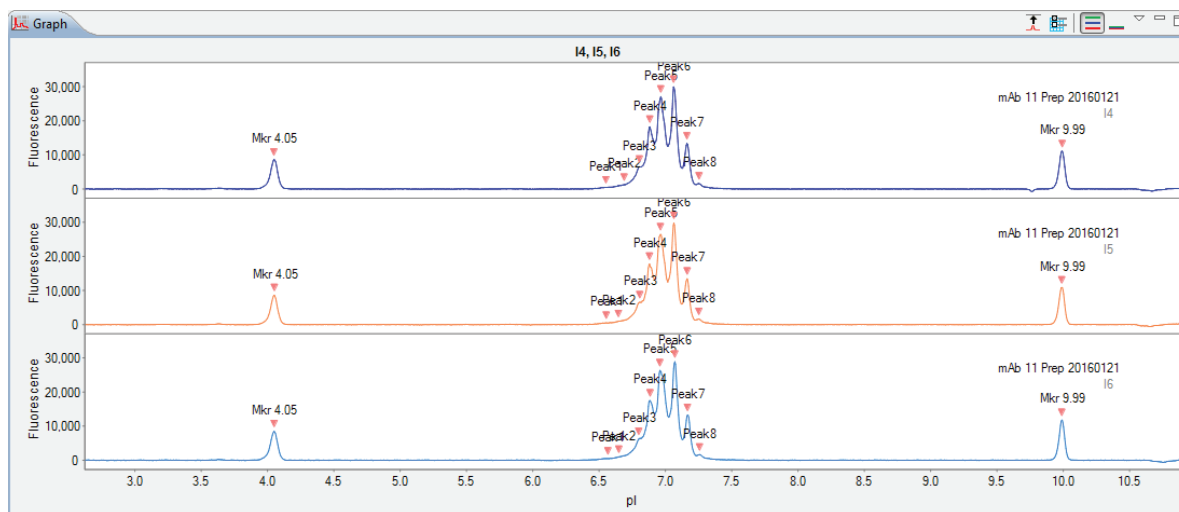
Here's an example of an electropherogram with all plot labels selected:



Stacking Multiple Electropherograms

You can stack electropherograms for multiple injections vertically in the Graph pane for comparison.

1. Click **Single View**.
2. Select multiple injection rows in the Experiment pane.
3. Click the **Stack the Plots** button. The individual electropherograms for each injection you selected will stack in the Graph pane.

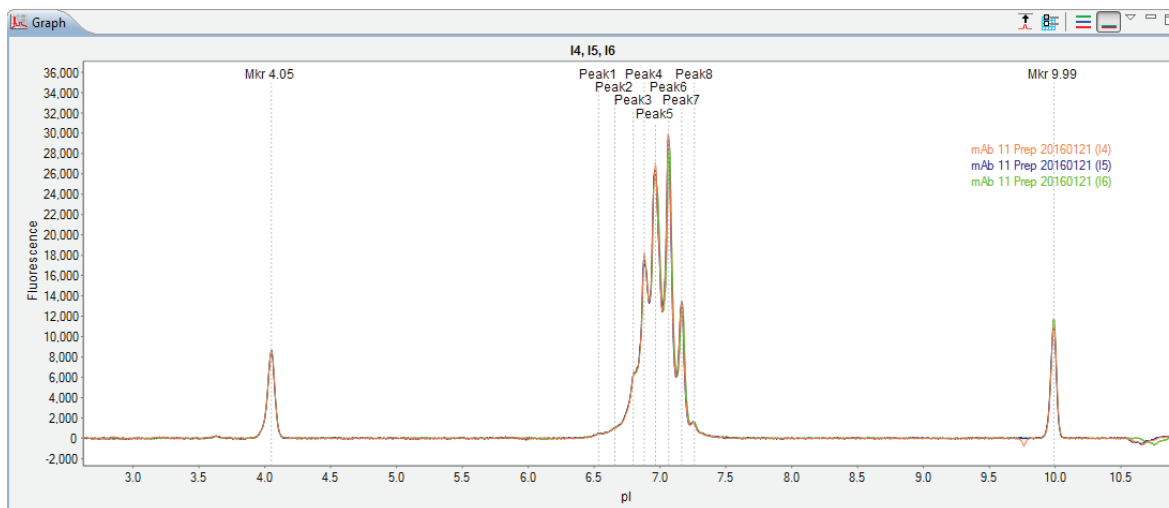


You can also customize the colors used for the stacked plot display. To do that go to “Setting up Automatic Injection Reports” on page 761.

Overlaying Multiple Electropherograms

You can overlay electropherograms for multiple injections on top of each other for comparison in the Graph pane.

1. Click **Single View**.
2. Select multiple injection rows in the Experiment pane.
3. Click the **Overlay the Plots** button. The individual electropherograms for each injection you selected will overlay in the Graph pane.



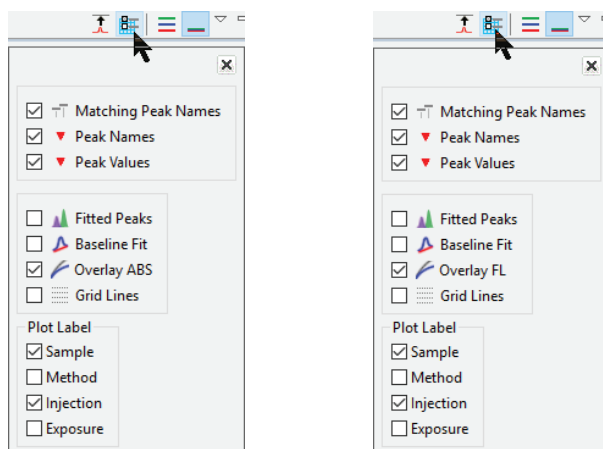
You can also customize the colors used for the overlay plot display. To do that go to “Setting up Automatic Injection Reports” on page 761.

Overlaying Fluorescence and Absorbance Electropherograms (standard cIEF runs only)

You can overlay standard Maurice cIEF fluorescence and absorbance electropherograms for the selected injection in the Graph pane. Due to the different detection methods you’ll see slight differences between the two electropherograms when they’re overlaid.

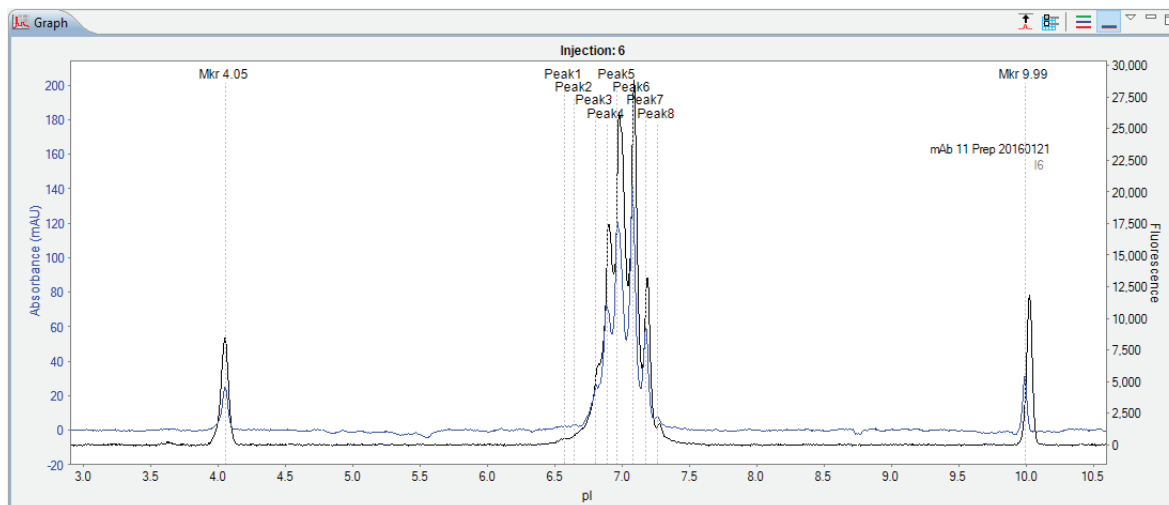
1. Click **Single View**.
2. Select an injection row in the Experiment pane.

3. Click **Graph Options** and select either **Overlay Abs** or **Overlay FL**. The option available depends on what detection mode is currently displaying.



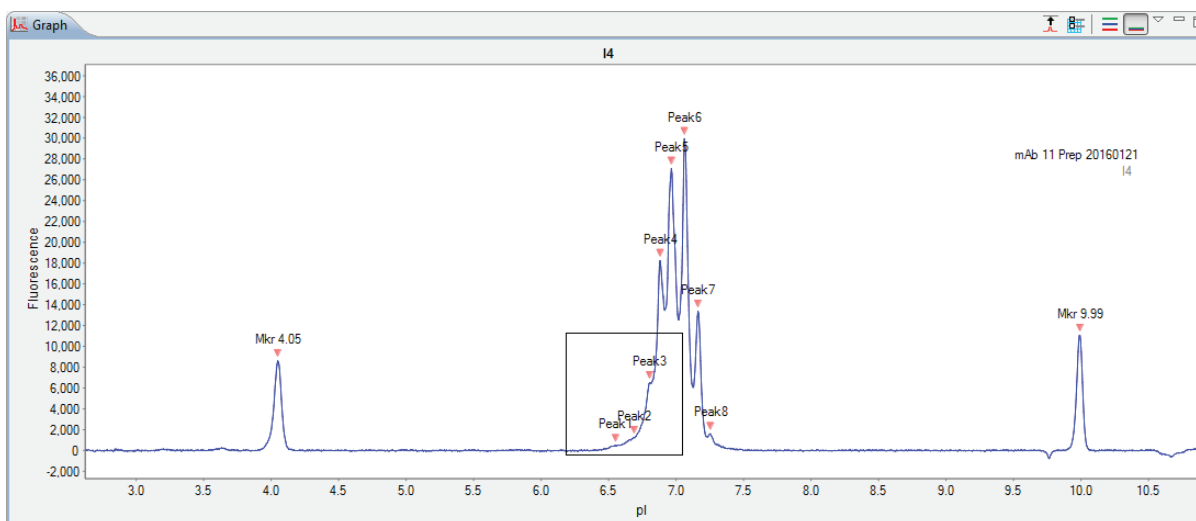
4. The electropherograms for each detection mode will overlay in the Graph pane, and the other detection mode y axis will display on the right.

NOTE: It's helpful to turn on Auto Scale when you're comparing the traces.

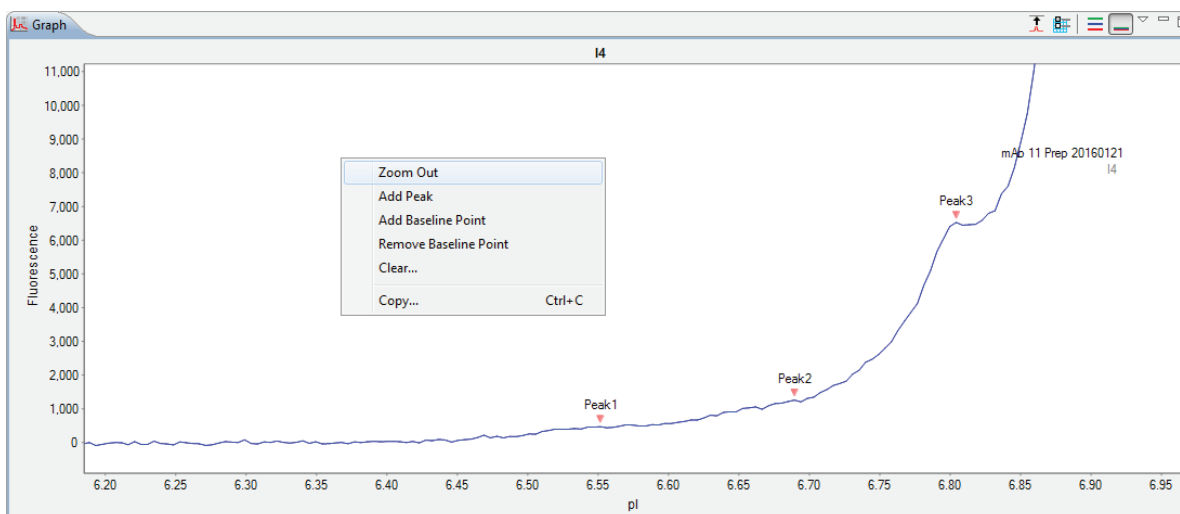


Zooming

To zoom in on a specific area of the electropherogram, hold the mouse button down and draw a box around the area with your mouse:

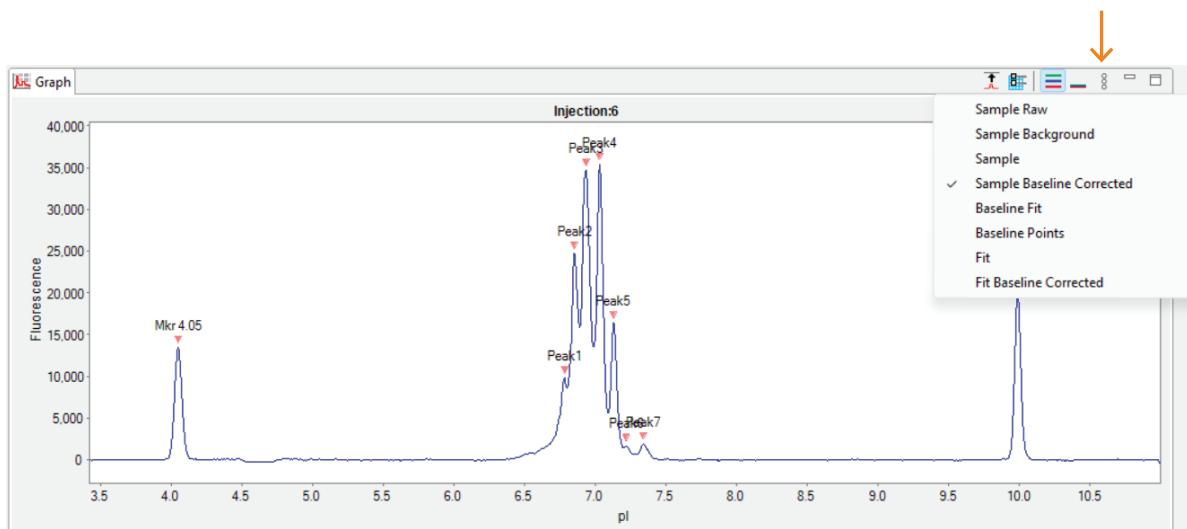


To return to default scaling, right click in the electropherogram and click **Zoom Out**.



Selecting Data Viewing Options

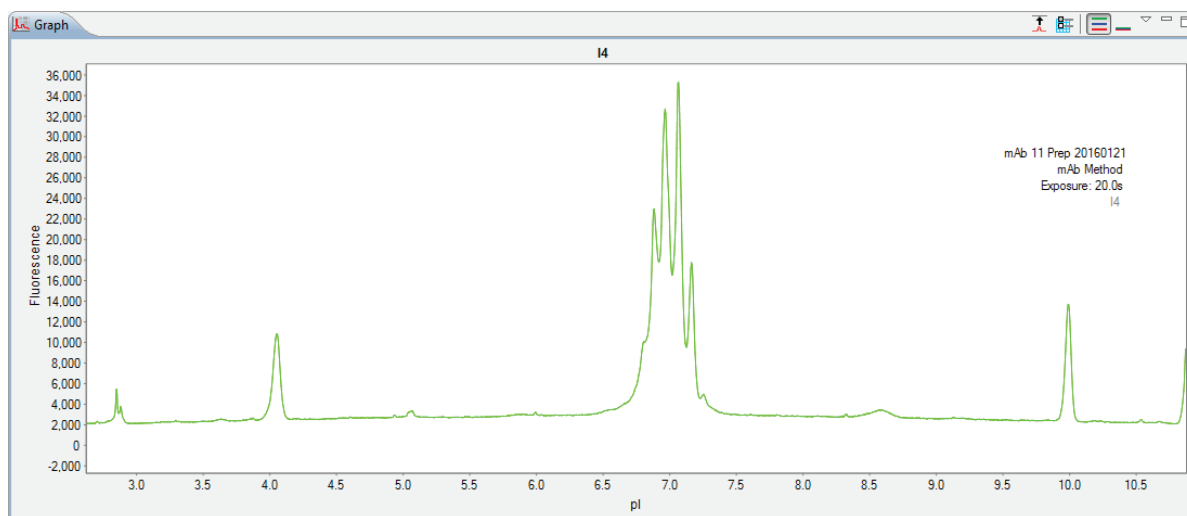
The graph view menu gives you multiple options for changing what type of electropherogram data is displayed. Just click **View Menu** in the graph pane toolbar to view the menu:



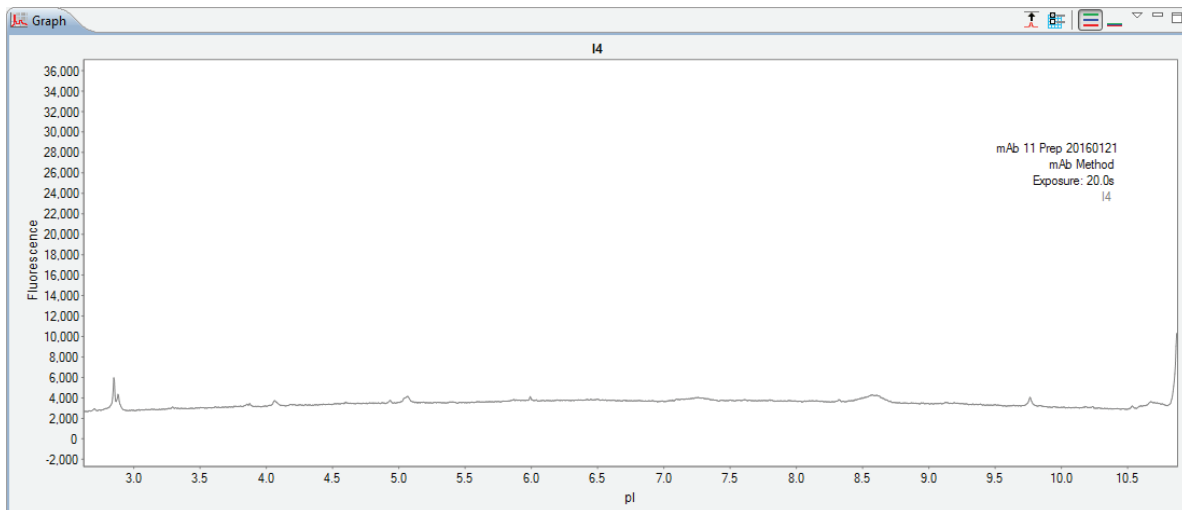
A check mark next to the menu option indicates it's currently selected, and you can select multiple options at once.

NOTE: Unless noted otherwise, graph view menu options are available for sample data only.

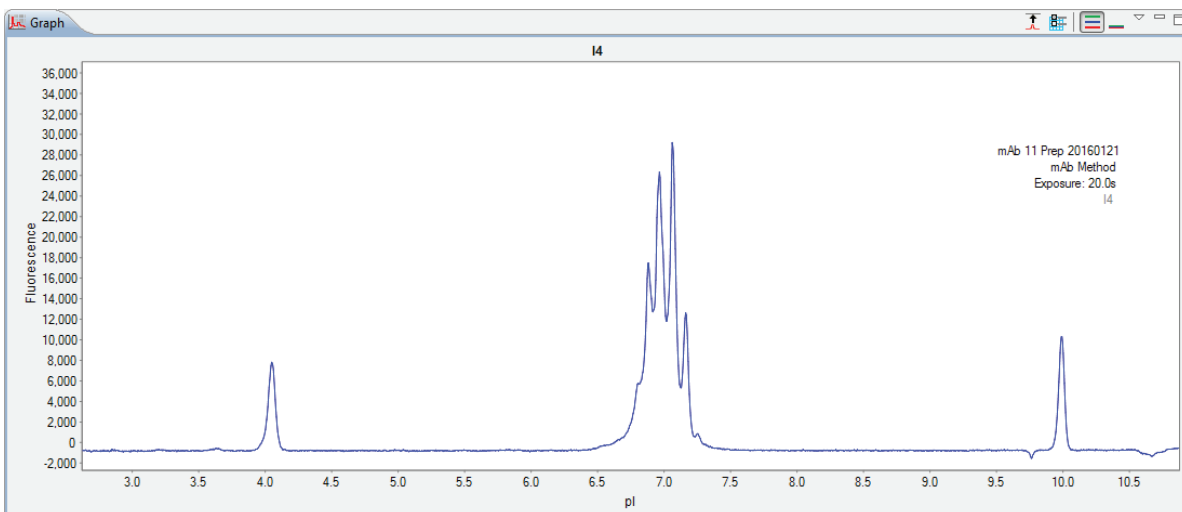
- **Sample Raw** - Clicking this option displays the basic detector values used to calculate peak absorbance or peak fluorescence depending on which mode is being viewed.



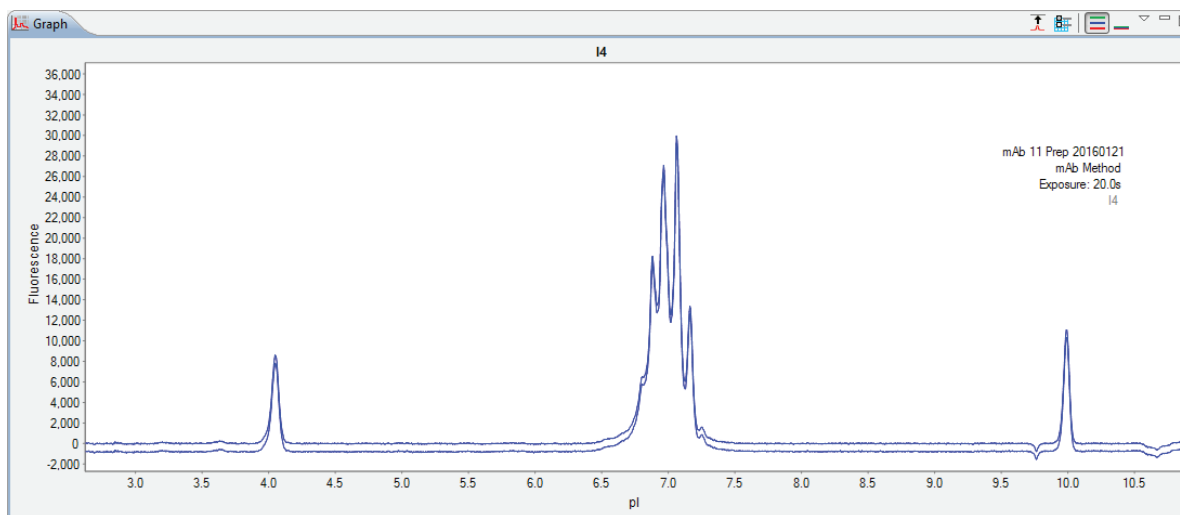
- **Sample Background** - Clicking this option displays the basic detector values used to calculate baseline absorbance or peak fluorescence depending on which mode is being viewed.



- **Sample** - Clicking this option displays raw, uncorrected sample data.

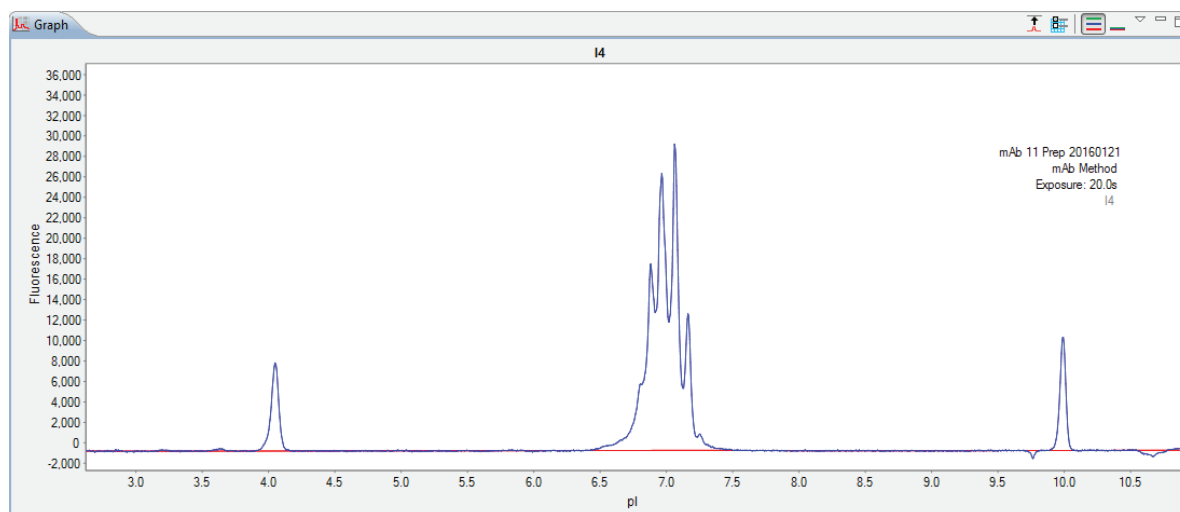


- **Sample Baseline Corrected** - Clicking this option displays sample data with the baseline subtracted (zeroed). This is the default view. In this next example, both Sample and Sample Baseline Corrected are selected.



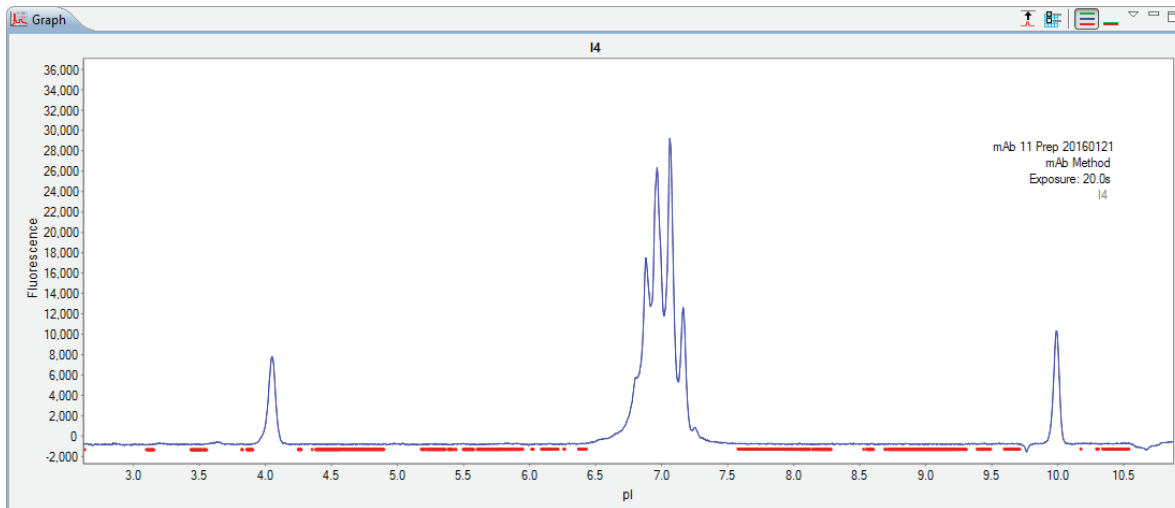
- **Baseline Fit** - Clicking this option displays the calculated baseline for the raw sample data. In this next example, both Baseline Fit and Sample are selected.

NOTE: This option is selected automatically when Baseline Fit is selected in graph options.

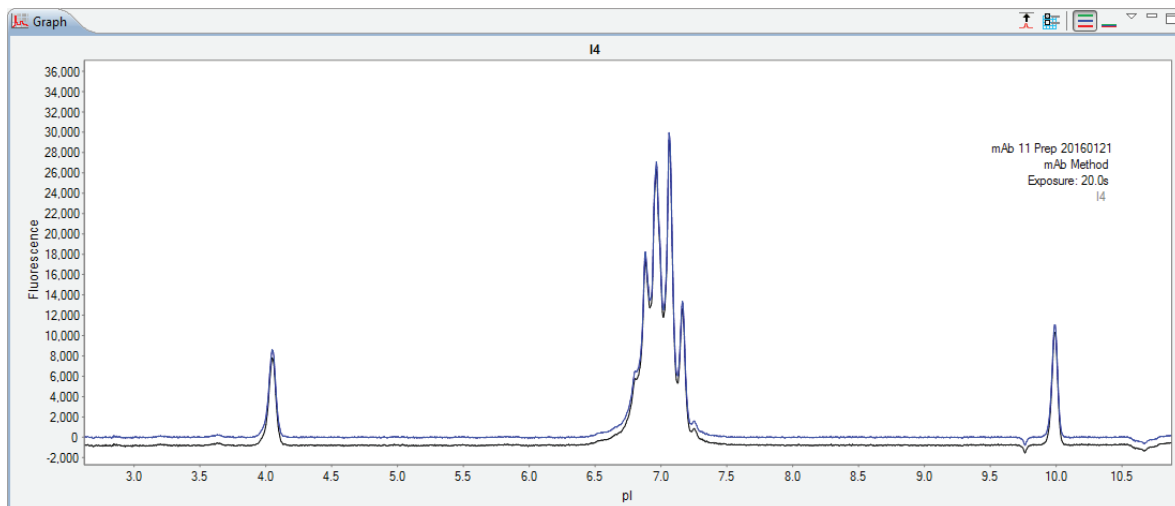


- **Baseline Points** - Clicking this option displays regions of the electropherogram considered to be at baseline. In this example, both Baseline Points and Sample are selected.

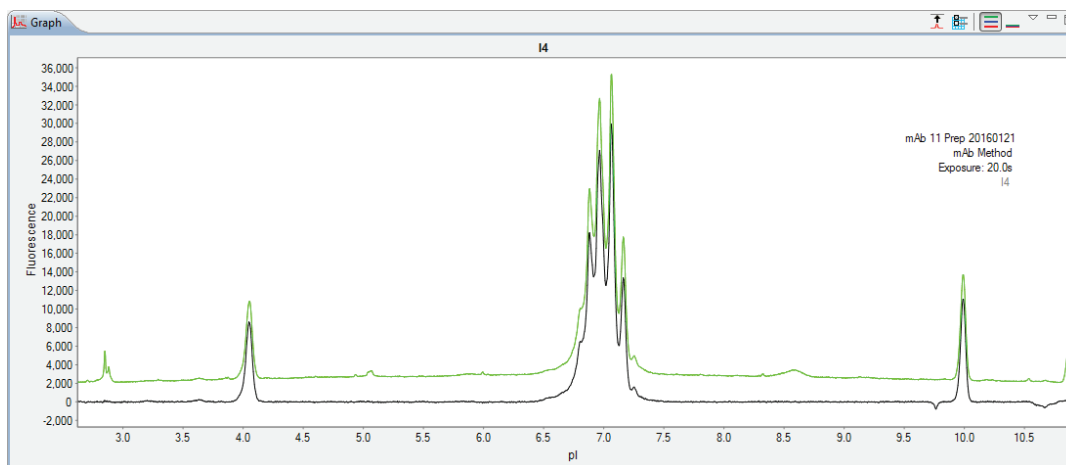
NOTE: This option is selected automatically when Baseline Fit is selected in graph options.



- **Fit** - Clicking this option displays the bounding envelope of the fitted peaks as calculated by the software for the raw sample data. In this example, both Fit and Sample Baseline Corrected are selected.



- **Fit Baseline Corrected** - Clicking this option displays the fitted peaks as calculated by the software for the sample baseline corrected data. In this example, both Fit Baseline Corrected and Sample Raw are selected, the fit plot is on the bottom.

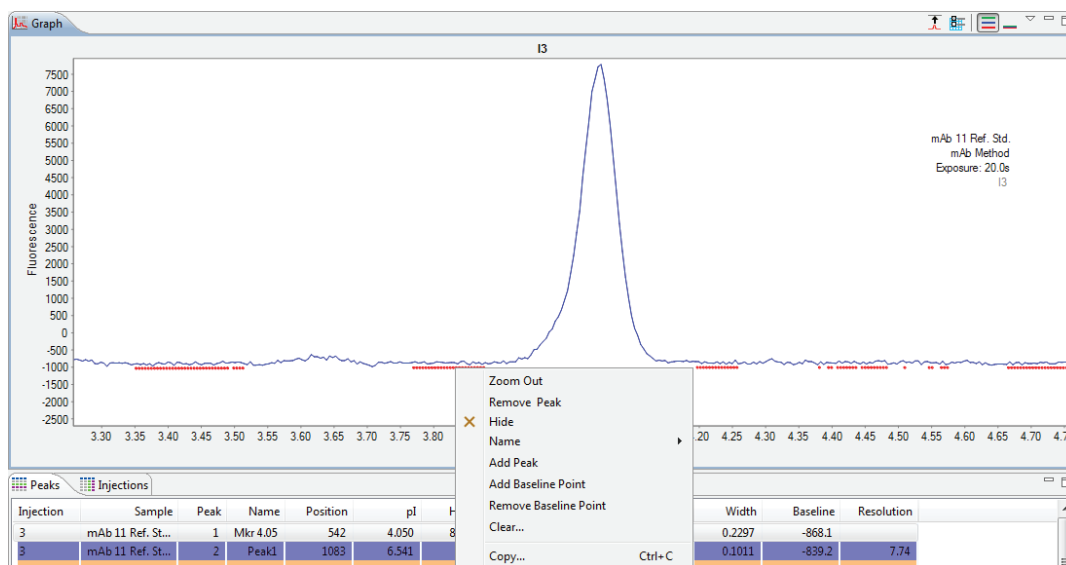


Adjusting the Baseline

Spline Baseline

Points in the baseline can be added or removed as needed.

1. Click the **Graph Options** button in the graph pane toolbar and check **Baseline Points**. This will display baseline points for the raw sample data.
2. Use the mouse to draw a box around the area you want to correct. This will zoom in on the area.
3. Right click a baseline point and select **Add Baseline Point** or **Remove Baseline Point**.



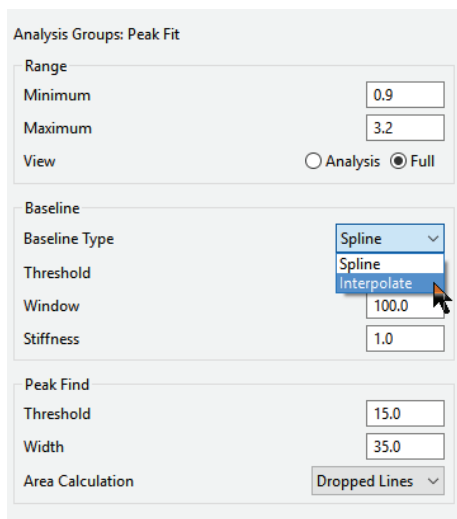
NOTE: To clear the manual addition or removal of baseline points and go back to the original view of the data, right click in the electropherogram and click **Clear All**.

Interpolated Baseline

Compass for iCE lets you interpolate baselines from the base of each peak for named peaks that use Dropped Lines integration.

1. If you haven't already, name your peaks of interest.
2. Select **Edit > Analysis**, and click **Peak Fit** in the left sidebar.
3. Choose **Interpolate** as the Baseline Type if it isn't already selected.

NOTE: Baseline Type selection is only available when Dropped Lines is used for the Area Calculation.



Analysis Groups: Peak Fit

Range

Minimum 0.9

Maximum 3.2

View ☐ Analysis ☒ Full

Baseline

Baseline Type Spline ▼

Threshold 100.0

Window 1.0

Peak Find

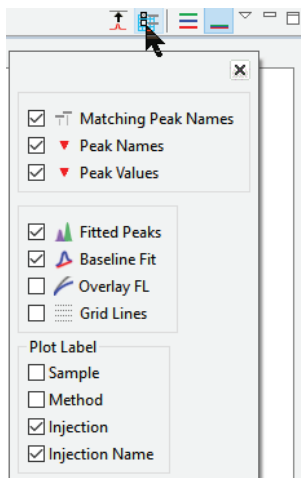
Threshold 15.0

Width 35.0

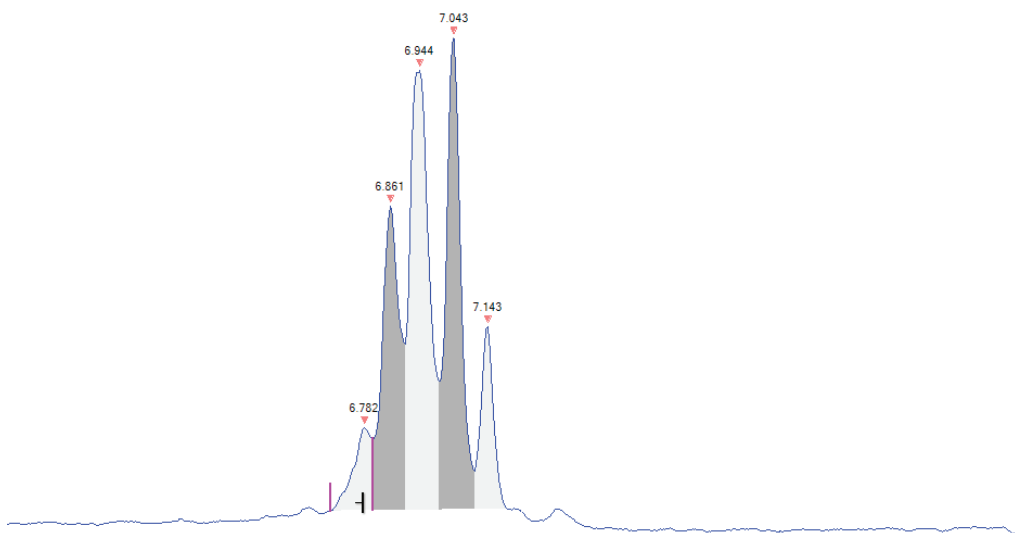
Area Calculation Dropped Lines ▼

4. Click **OK**.

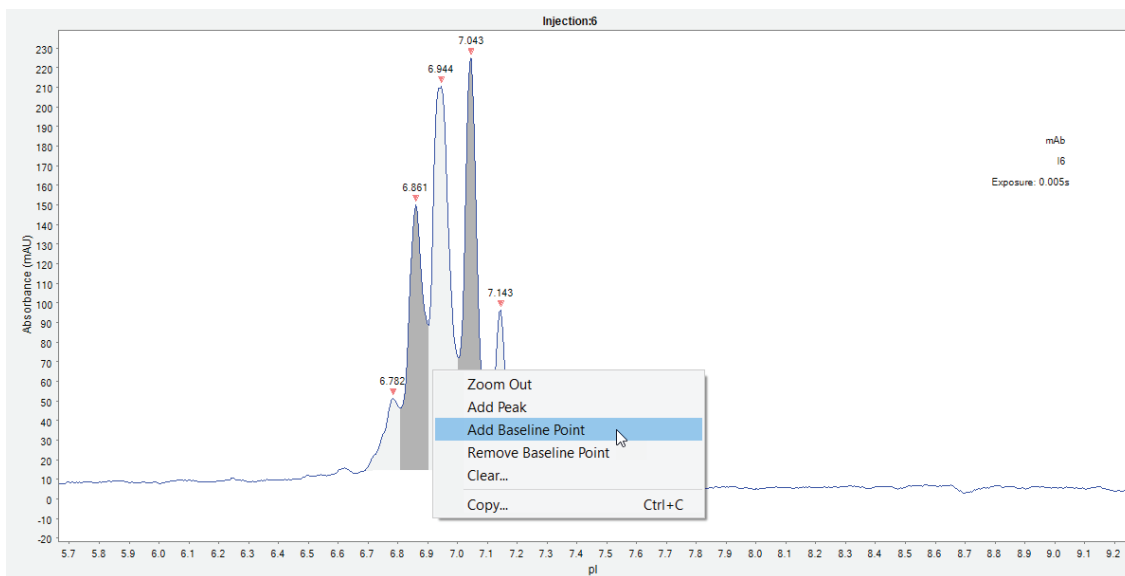
- In the Analysis window Graph Pane, click **Graph Options** and select **Fitted Peaks** and **Baseline Fit**.



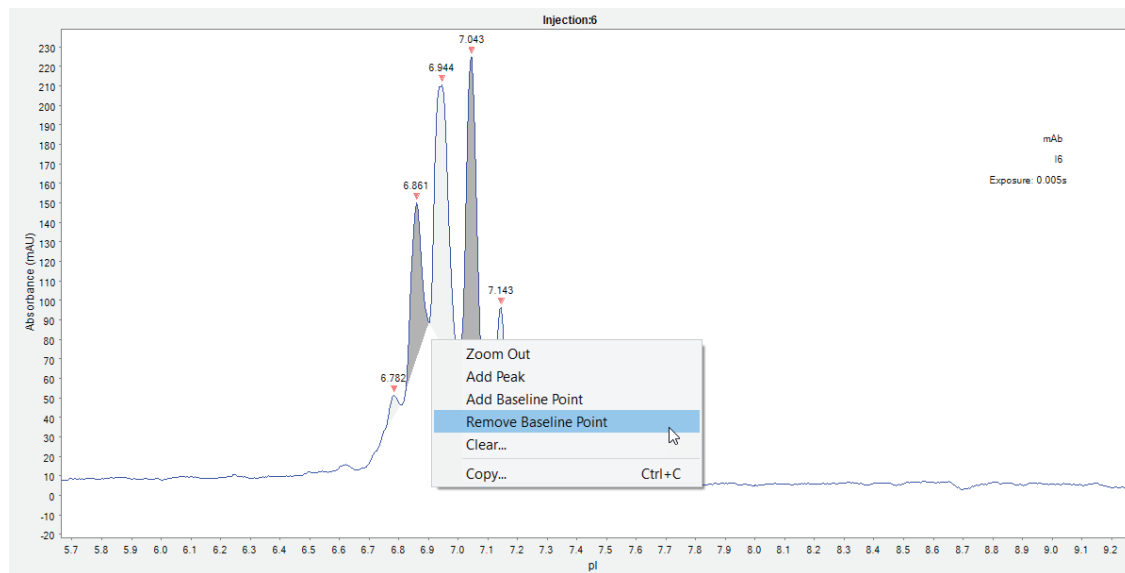
- Select an injection in the Experiment pane.
- Hover the mouse over the peak you want to adjust the baseline on. Two magenta bars will display that show the start and end points of the peak's integration.
- To change the baseline, hover the mouse over one of the magenta bars. Then click the mouse and drag the bar to move it. The cursor changes to indicate which direction you can move the integration point. See "Manual Peak Integration" on page 547 for more info on integrating peaks.



9. Baseline points can also be added between two adjacent peaks for a better fit. Hover the mouse over the peak until the magenta bars appear, then right click and select **Add Baseline Point**.

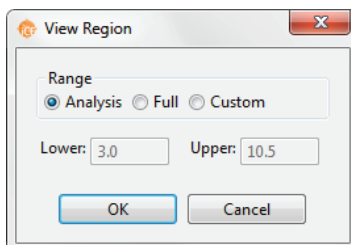


10. To remove the baseline point, hover the mouse over the peak until the magenta bars appear, then right click and select **Remove Baseline Point**.

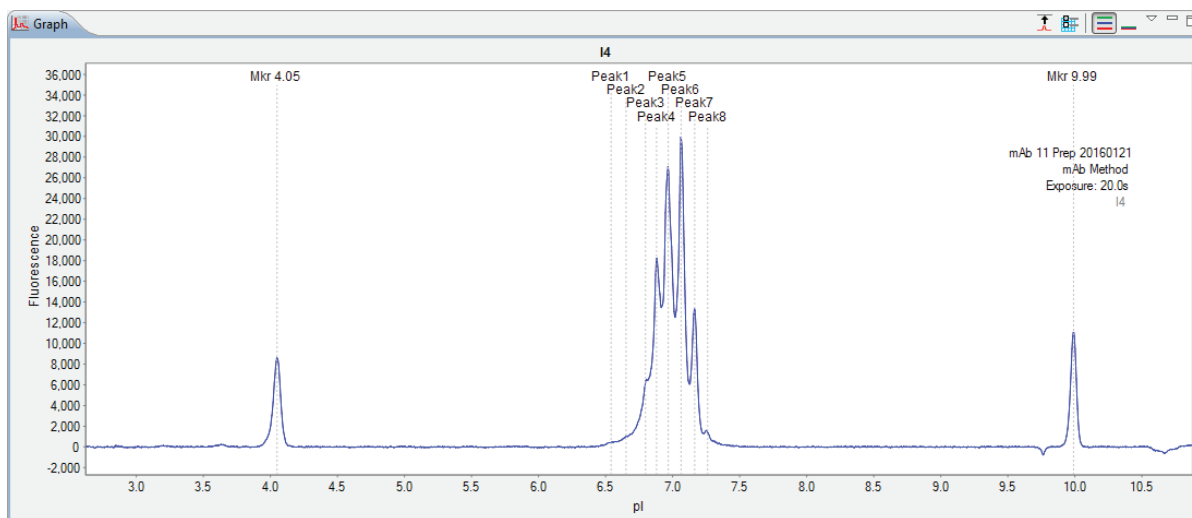


Selecting the Graph X-axis Range

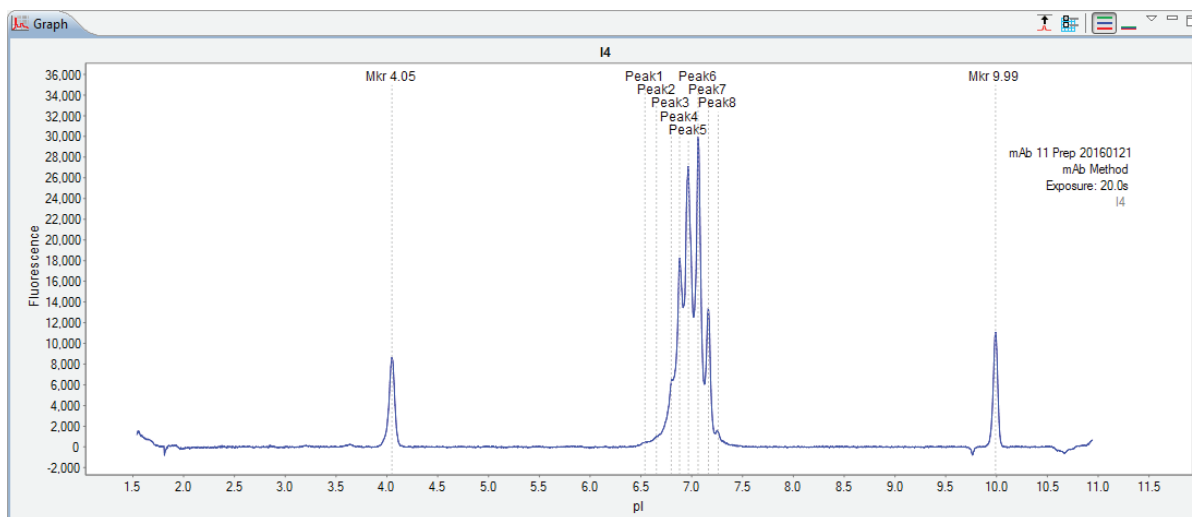
The pI range used for the x-axis can be changed. Just select **View** in the main menu and click **View Region**.



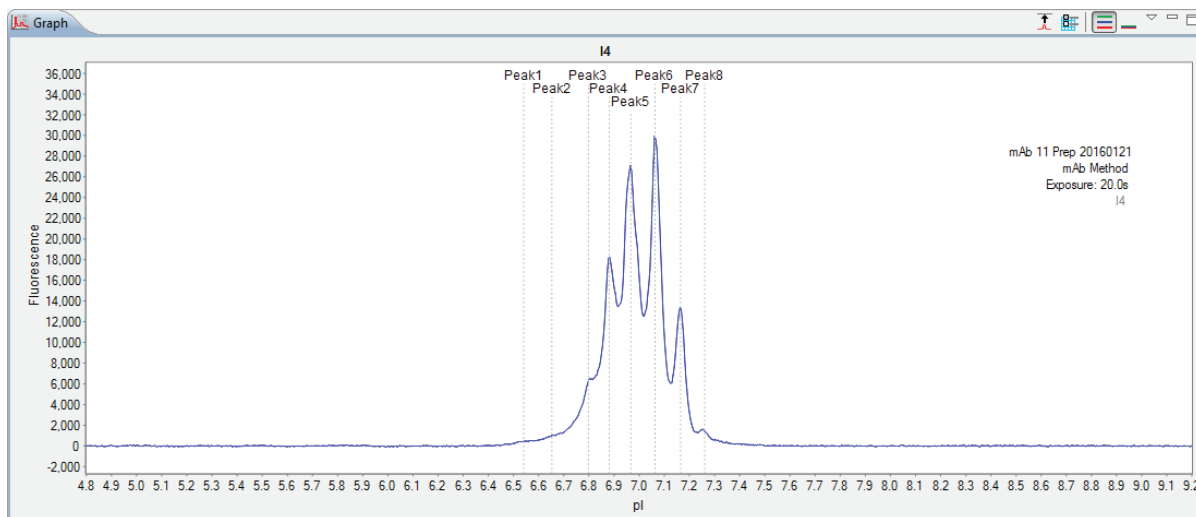
- **Analysis** sets the x-axis range of the electropherogram to what is selected in the Peak Fit range settings. To view or change these analysis settings, go to **Edit > Analysis** and click **Peak Fit** in the left sidebar. In this example, the lower and upper range settings are 3.0 and 10.5.



- **Full** displays the entire separation in the electropherogram. This is the default setting. In this example the lower and upper range settings are 1.5 and 11.4.



- **Custom** lets you manually enter the lower and upper range settings to display in the electropherogram. In this example the lower and upper range settings are 5.0 and 9.0.



NOTE: You can change the default x-axis range that Compass for iCE uses. Go to “Advanced Analysis Settings” on page 551 for more info.

Closing Run Files

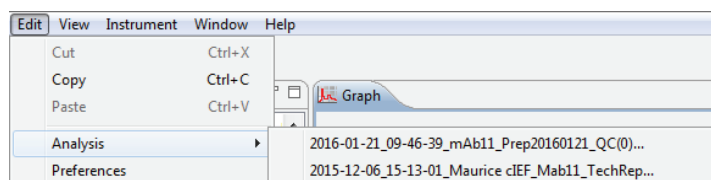
If more than one run file is open, you can close just one file or all the open files at the same time.

- **To close one run file** - In the Experiment pane, click on one of the sample rows in the file. Then click **File** from the main menu and click **Close**.
- **To close all open run files** - Select **File** from the main menu and click **Close All**.

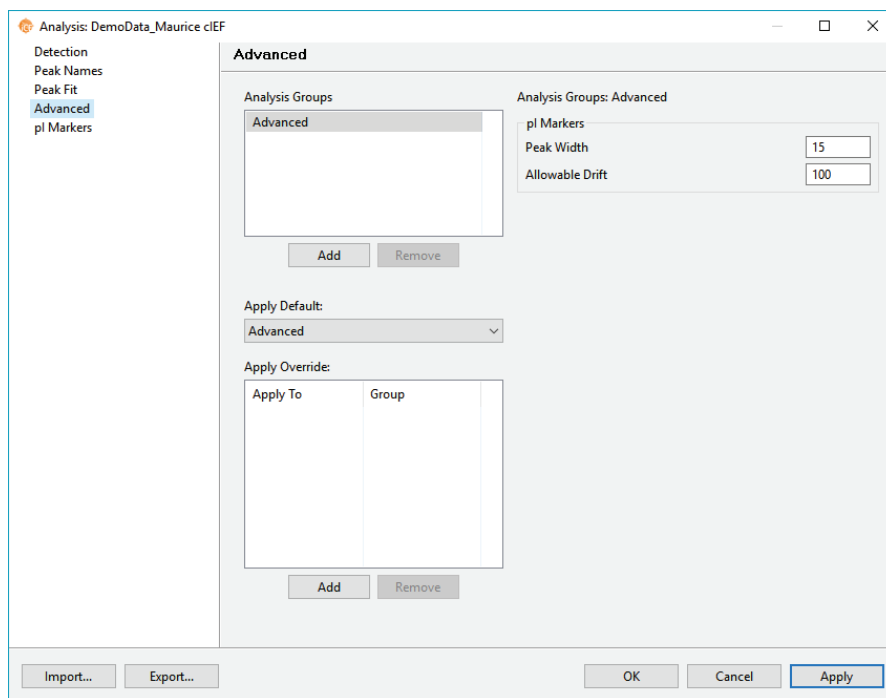
Analysis Settings Overview

Compass for iCE has many analysis features and settings that you can change to enhance your run data.

Select **Edit** in the main menu and click **Analysis**. If more than one run file is open, select the run file you want to view settings for from the list:



This opens the Analysis window:



To move between pages in the window, click on an option in the left sidebar.

- **Detection (Standard Maurice cIEF runs only)** - Lets you choose to view absorbance or native fluorescence data for the run and choose data at different fluorescence exposures.
- **Peak Names** - Lets you enter custom naming settings for sample proteins and have Compass for iCE automatically label the peaks in the run data.
- **Peak Fit** - Lets you customize peak fit settings for sample data.
- **Advanced** - Lets you customize analysis settings for the pI markers.
- **pI Markers** - Lets you customize the pI markers and positions Compass for iCE identifies for each method in your run.

On all pages in the Analysis window:

- Click **Import** to import an analysis settings file. Go to “Importing Analysis Settings” on page 567 to learn how to do this.
- Click **Export** to export the current analysis settings file. Go to “Exporting Analysis Settings” on page 568 to learn how to do this.
- Click **Apply** to apply changes to the run file and update results in real time.
- Click **OK** to save changes to the run file and exit.
- Click **Cancel** to exit without saving changes.

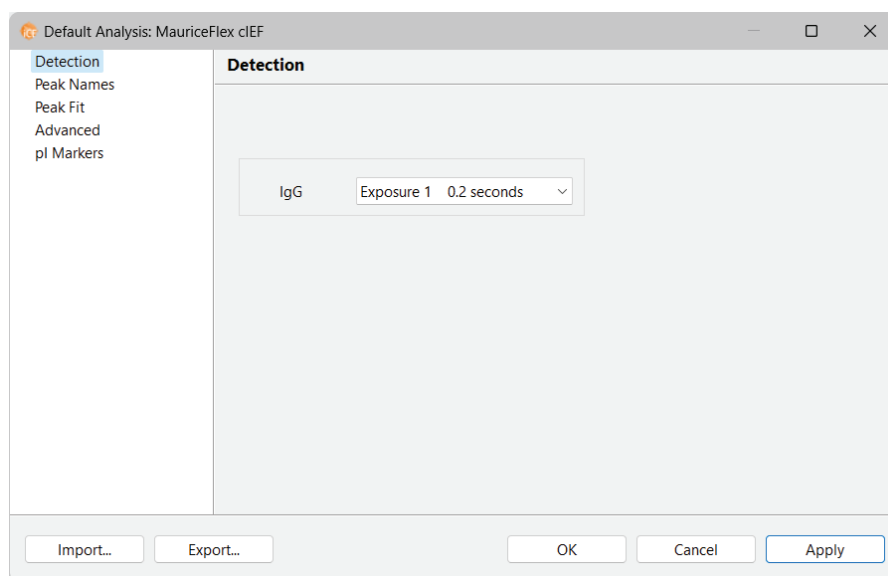
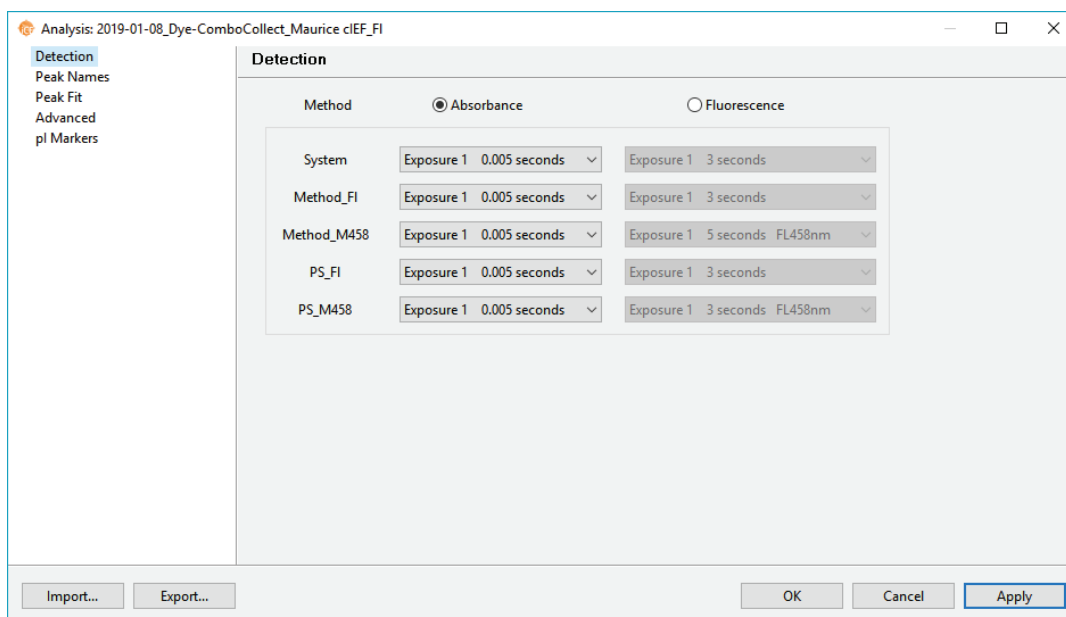
Detection Settings

This page lets you see the absorbance and native fluorescence exposures taken during the run, and select different exposures for data viewing in the Analysis screen. Select **Edit** in the main menu and click **Analysis**, then click **Detection** in the left sidebar.

NOTES:

The FL458 nm fluorescence filter is only available on Maurice, Maurice C. and MauriceFlex systems with the option installed. On MauriceFlex, the filter can only be used with standard Maurice cIEF runs.

Absorbance detection is not available for MauriceFlex cIEF runs.



Changing the Detection Method (Standard Maurice cIEF runs only)

You can choose to display either absorbance or fluorescence data for your run in the Analysis screen.

1. Select **Edit > Analysis**, and select **Detection** in the left sidebar.

2. Select either the Absorbance or Fluorescence radio button.

Detection

Method ☒ Absorbance ☐ Fluorescence

Method	Absorbance	Fluorescence
System	Exposure 1 0.005 seconds	Exposure 1 3 seconds
Method_FI	Exposure 1 0.005 seconds	Exposure 1 3 seconds
Method_M458	Exposure 1 0.005 seconds	Exposure 1 5 seconds FL458nm
PS_FI	Exposure 1 0.005 seconds	Exposure 1 3 seconds
PS_M458	Exposure 1 0.005 seconds	Exposure 1 3 seconds FL458nm

Detection

Method ☐ Absorbance ☒ Fluorescence

Method	Absorbance	Fluorescence
System	Exposure 1 0.005 seconds	Exposure 1 3 seconds
Method_FI	Exposure 1 0.005 seconds	Exposure 1 3 seconds
Method_M458	Exposure 1 0.005 seconds	Exposure 1 5 seconds FL458nm
PS_FI	Exposure 1 0.005 seconds	Exposure 1 3 seconds
PS_M458	Exposure 1 0.005 seconds	Exposure 1 3 seconds FL458nm

Changing the Detection Exposure

You can change the exposure used for the sample data displayed in the Analysis screen.

NOTES:

You'll only be able to choose exposures for the detection method currently selected.

The number of exposures taken and exposure times shown are specified in the method when you set up your batch. They can't be changed after the run has executed.

The Absorbance exposure at 0.005 seconds is an instrument default exposure setting for Maurice cIEF runs. No other absorbance exposures are available.

1. Select **Edit > Analysis**, and select **Detection** in the left sidebar.

- Click the arrow in the exposure button you want to change and select an exposure setting:

The screenshot shows the 'Detection' settings window. It has two tabs: 'Absorbance' and 'Fluorescence'. The 'Fluorescence' tab is selected. Below the tabs, there are several rows of settings for different methods. Each row has a dropdown menu for 'Exposure' and a dropdown menu for 'Time'. The 'Exposure' dropdown is open, showing a list of options: 'Exposure 1 3 seconds', 'Exposure 2 5 seconds', 'Exposure 3 10 seconds', and 'Exposure 4 20 seconds'. The 'Time' dropdown is also open, showing a list of options: 'Exposure 1 3 seconds', 'Exposure 2 5 seconds', 'Exposure 3 10 seconds', and 'Exposure 4 20 seconds'.

- Click OK to save changes. Sample data for the exposure selected will display in the Analysis screen.

Peak Names Settings

This page lets you view and change custom naming settings for sample proteins. Select **Edit** in the main menu and click **Analysis**, then click **Peak Names** in the left sidebar.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.

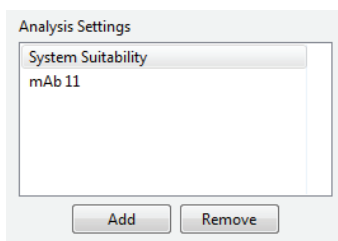
The screenshot shows the 'Peak Names' settings window. It has a sidebar on the left with a list of settings: 'Detection', 'Peak Names', 'Peak Fit', 'Advanced', and 'pI Markers'. The 'Peak Names' setting is selected. The main area of the window is divided into two sections: 'Analysis Groups' and 'Apply Settings'. The 'Analysis Groups' section has a table with columns 'Name', 'pI', 'Color', and 'Range'. Below the table are 'Add' and 'Remove' buttons. The 'Apply Settings' section has a table with columns 'Apply To' and 'Group'. Below the table are 'Add' and 'Remove' buttons. At the bottom of the window are 'Import...', 'Export...', 'OK', 'Cancel', and 'Apply' buttons.

Peak Names Analysis Settings Groups

Peak name settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTE: Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 567.

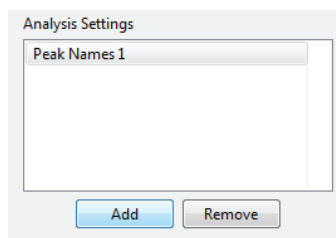
Peak name groups are displayed in the analysis settings box:



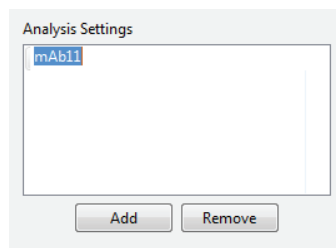
There aren't any Compass for iCE default settings groups, but you can make changes to groups you've created and create new groups. To view settings for a group, click on the group name in the analysis settings box.

Creating a Peak Names Group

1. Select **Edit > Analysis**, and select **Peak Names** in the left sidebar.
2. Click **Add** under the analysis settings box.



3. Enter a new name for the group.




4. Click in the first cell in the **Name** column in the analysis settings peak table and enter a sample protein name.

[illegible]

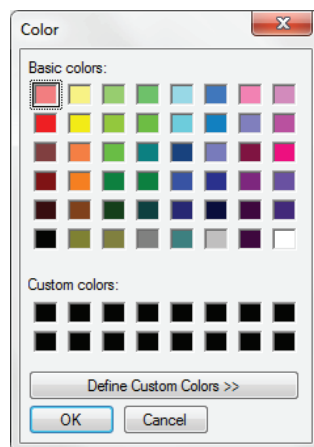
5. Click in the first cell in the **pI** column and enter the expected pI for the sample protein.

Name	pI	Color	Range
Peak1	5.55		0.05

6. Click in the first cell in the **Color** column, then click the button.


Name	pI	Color	Range
Peak1	6.55	 (0.1 ...)	0.05

The color selection box displays:




7. The color you pick is used to identify the sample protein peak in the Peaks and Injections panes in the Analysis screen. Click a color or define a custom color and click OK. The color selection will update in the table:

Analysis Settings: mAb11

Name	pI	Color	Range
Peak1	6.55		0.05








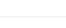
8. Click in the first cell in the **Range** column.

Analysis Settings: mAb11

Name	pI	Color	Range
Peak1	6.55		0.1

9. Enter a \pm range for the pI entered. Compass for iCE will automatically name peaks found within this pI range. For example, if the pI entered is 2 and a 0.1 range is used, all peaks with pIs between 1.9 and 2.1 will be identified with this peak name and color.
10. To add another sample protein, click Add under the peak table. Repeat the previous steps for other sample proteins. In this example, eight proteins were entered:

Analysis Settings: mAb 11

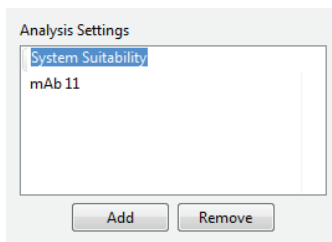
Name	pI	Color	Range
Peak1	6.55		0.1
Peak2	6.65		0.1
Peak3	6.8		0.1
Peak4	6.9		0.1
Peak5	7		0.1
Peak6	7.1		0.1
Peak7	7.2		0.1
Peak8	7.3		0.1

To remove a sample protein, select its row and click **Remove**.

11. Click **OK** to save changes.

Modifying a Peak Names Group

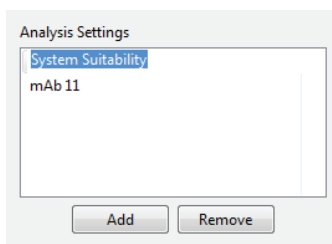
1. Select **Edit > Analysis**, then click **Peak Names** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



3. Change the information in the analysis settings peak table as described in “Creating a Peak Names Group” on page 533.
4. Click **OK** to save changes.

Deleting a Peak Names Group

1. Select **Edit > Analysis**, then click **Peak Names** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.

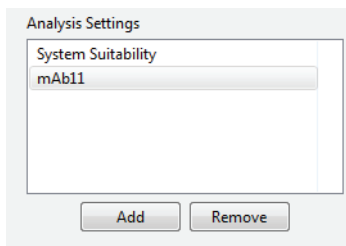


3. Click **OK** to save changes.

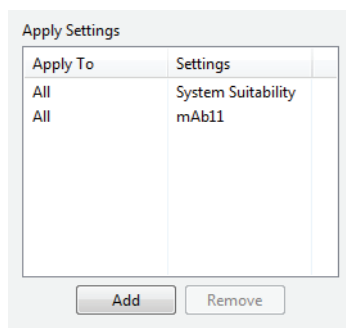
Applying Peak Names Groups to Run Data

1. Select **Edit > Analysis**, then click **Peak Names** in the options list.

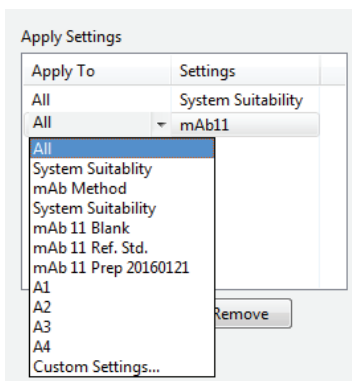
- Click on the group in the analysis settings box you want to apply to specific run data.



- Application of peak names groups to specific run data is done in the apply settings box. A default data set automatically gets created whenever you create a new group and it's applied to all injections in the run. You can either modify the default group or click Add under the box to create a new one.

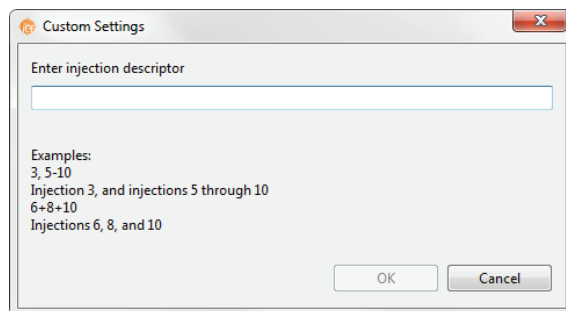


- Click the cell in the **Apply To** column, then click the down arrow.

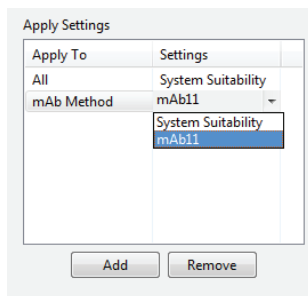


- Select an option from the drop down list. This applies the peak names group selected to specific run data as follows:
 - All** - Selecting this applies peak names group settings to all injections.
 - Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.

- **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.
- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:

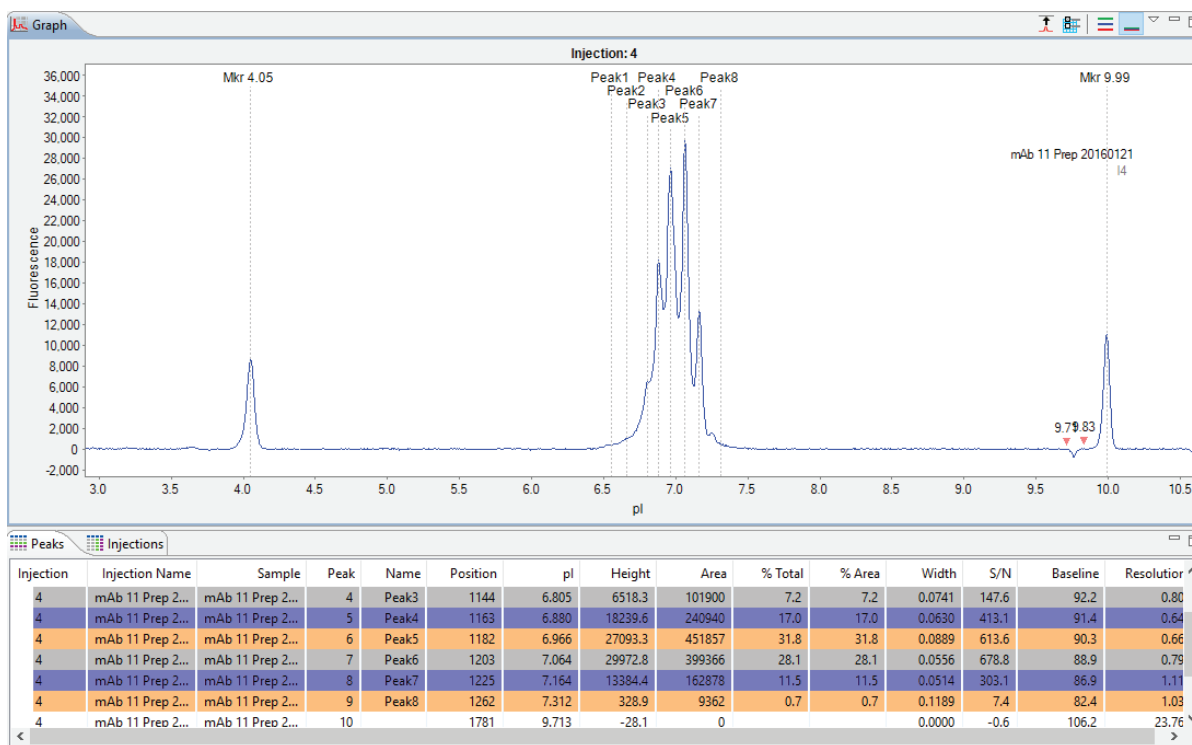


6. If you need to change the peak names group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.

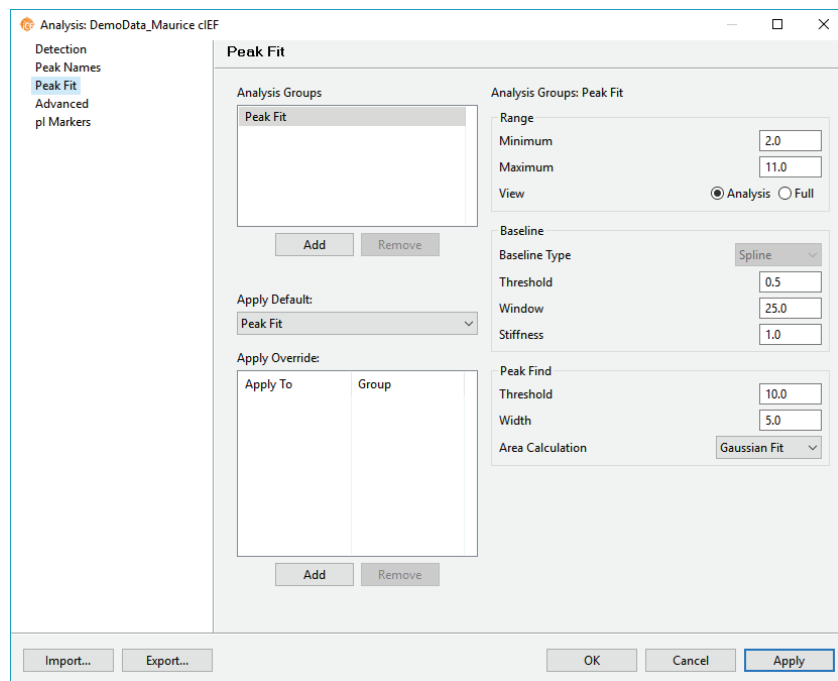
9. Click **OK** to save changes. Named peaks will be identified with a peak name label in the electropherogram and color-coded in the Peaks and Injections panes:



Peak Fit Analysis Settings

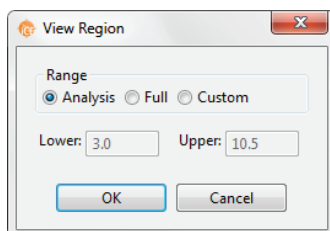
This page lets you view and change peak fit settings for sample data. Select **Edit** in the main menu and click **Analysis**, then click **Peak Fit** in the left sidebar:

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



Range Settings

- **Minimum** - The pI value below which peaks won't be identified. This value is also used as the default lower pI range for data displayed in the electropherogram.
- **Maximum** - The pI value above which peaks won't be identified. This value is also used as the default upper pI range for data displayed in the electropherogram.
- **View** - Sets the default range to either Full or Analysis for the electropherogram x-axis range in the View Region window (select **View** in the main menu and click **View Region**).



- **Analysis** sets the x-axis range of the electropherogram to the Peak Fit range minimum and maximum settings in the electropherogram. This is the default setting.

- **Full** displays the entire separation range of the run data in the electropherogram.

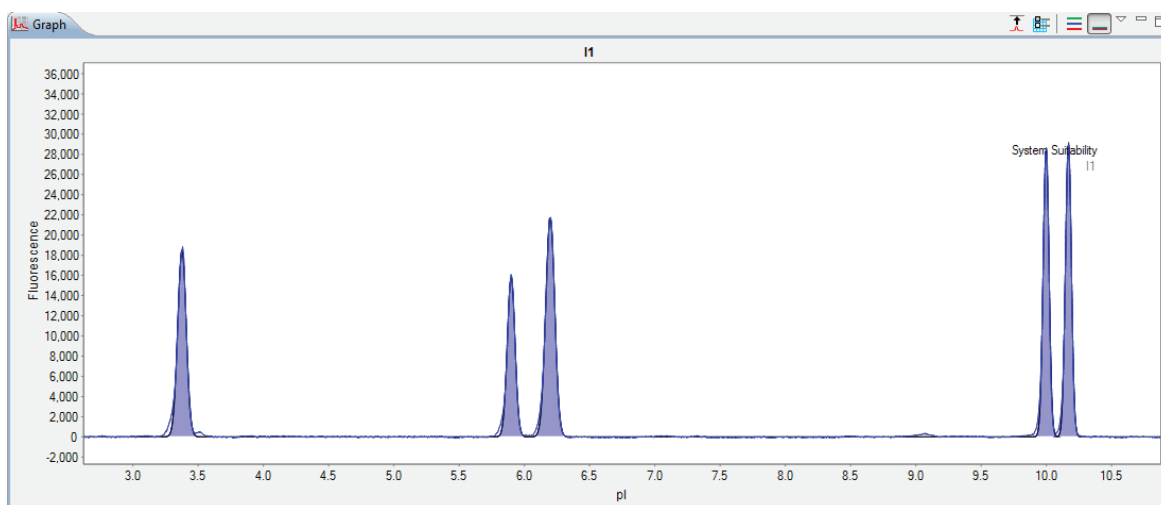
Baseline Settings

These settings apply to spline baselines only.

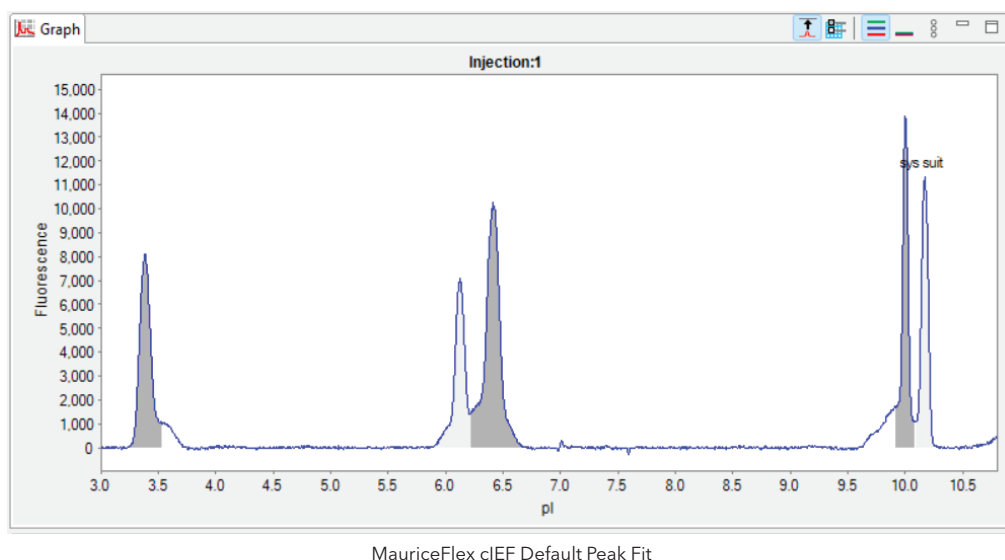
- **Threshold** - The variance, or roughness, in a baseline data segment below which a point is called part of the baseline.
- **Window** - How long baseline data segments are expected to be in pixels. Shorter segments let the baseline follow plateau sections of the signal.
- **Stiffness** - The amount the baseline is allowed to vary from a straight line. Settings between 0.1 and 1.0 make the baseline fit closer to a straight line. Settings from 1.0 to 10.0 will make the baseline fit follow the data more closely.

Peak Find Settings

- **Threshold** - The minimum signal to noise ratio required for a peak to be identified. A setting of 1.0 will detect many peaks, a setting of 10.0 will detect fewer peaks.
- **Width** - The approximate peak width (at full width half max) in pixels used to detect peaks. The minimum value for this setting is 3.0. Larger widths help eliminate the detection of shoulder and noise peaks.
- **Area Calculation** - Two fits are used, either Gaussian Fit or Dropped Lines. These settings can be changed before or after the run is finished.
- For Standard cIEF applications, peak area is calculated using Gaussian Fit by default. For MauriceFlex cIEF applications, peak area is calculated using Dropped Line by default.



Standard Maurice cIEF Default Peak Fit



- The Dropped Line method type of area calculation is also often called the perpendicular drop method. This is the preferred method when peaks overlap or are close to each other. It draws two vertical lines from the left and right bounds of the peak down to the x-axis and then measures the total area bounded by the signal curve, the x-axis ($y=0$ line), and the two vertical lines.

Peak Fit Analysis Settings Groups

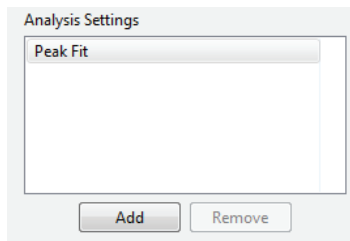
Peak fit settings are saved as a group, and you can create multiple settings groups. Specific group settings can then be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values for peak fit analysis settings. These settings are included in the default Peak Fit group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 567.

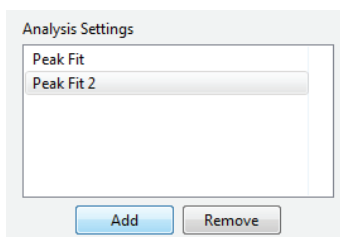
Peak fit groups are displayed in the analysis settings box:



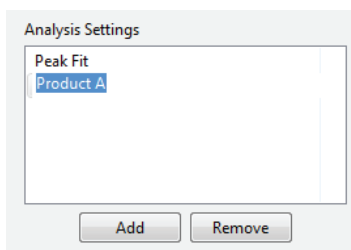
The Peak Fit group shown contains the Compass for iCE default analysis settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Peak Fit Group

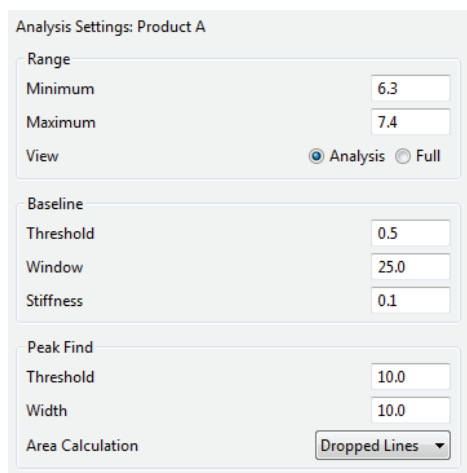
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



3. Click on the new group and enter a new name.

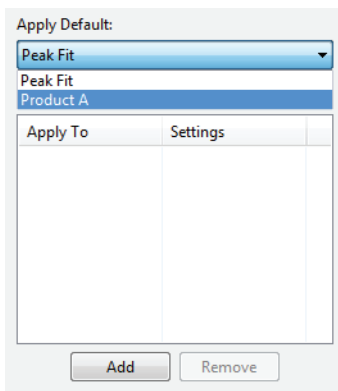


4. Change the settings in the range, baseline or peak find boxes as needed.

The screenshot shows a detailed configuration window titled "Analysis Settings: Product A". It is organized into several sections:

- Range:** Includes input fields for "Minimum" (6.3) and "Maximum" (7.4).
- View:** Includes radio buttons for "Analysis" (selected) and "Full".
- Baseline:** Includes input fields for "Threshold" (0.5), "Window" (25.0), and "Stiffness" (0.1).
- Peak Find:** Includes input fields for "Threshold" (10.0) and "Width" (10.0).
- Area Calculation:** Includes a dropdown menu currently set to "Dropped Lines".

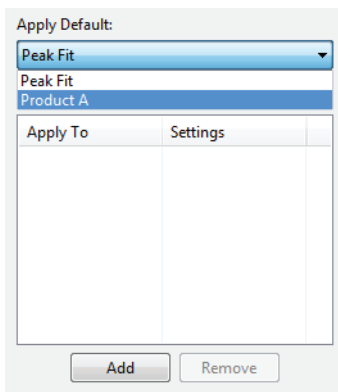
5. To use the new group as the default peak fit settings for the run file data, click the arrow in the drop down list next to Apply Default, then click the new group from the list. Peak fit settings in the new group will then be applied to the run data.



6. Click OK to save changes.

Changing the Default Peak Fit Group

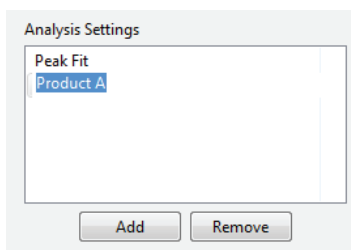
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then click a new default group from the list.



3. Click OK to save changes. Peak fit settings in the group selected will be applied to the run data.

Modifying a Peak Fit Group

1. Select **Edit > Analysis**, and click **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



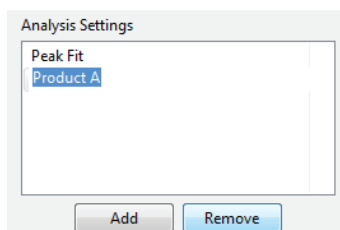
3. Change the settings in the range, baseline or peak find boxes as needed.

The image shows a detailed "Analysis Settings: Product A" dialog box. It is organized into three main sections. The "Range" section includes "Minimum" (6.3) and "Maximum" (7.4) input fields, and a "View" section with "Analysis" (selected) and "Full" radio buttons. The "Baseline" section includes "Threshold" (0.5), "Window" (25.0), and "Stiffness" (0.1) input fields. The "Peak Find" section includes "Threshold" (10.0) and "Width" (10.0) input fields, and an "Area Calculation" dropdown menu currently set to "Dropped Lines".

4. Click **OK** to save changes. The new peak fit settings will be applied to the run data.

Deleting a Peak Fit Group

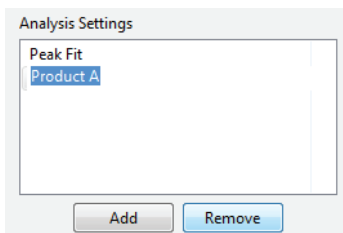
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



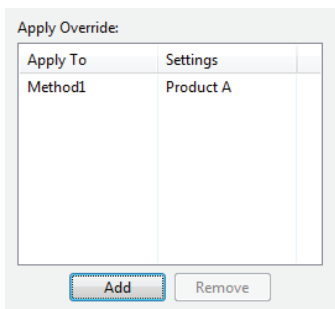
3. Click **OK** to save changes.

Applying Peak Fit Groups to Specific Run Data

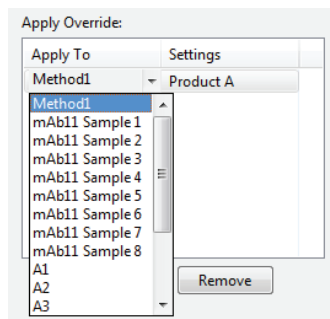
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to apply to specific run data.



3. Application of analysis groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

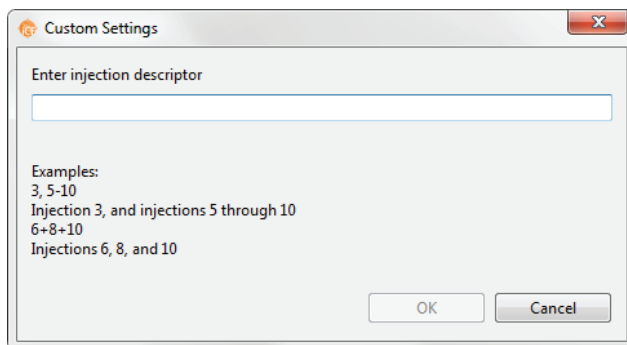


4. Click the cell in the **Apply To** column, then click the down arrow.

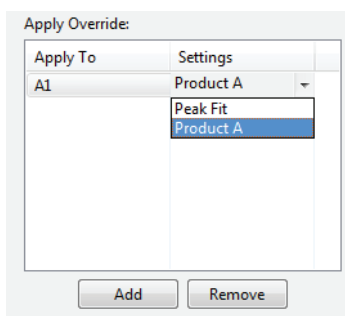


5. Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
 - **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.

- **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.
- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



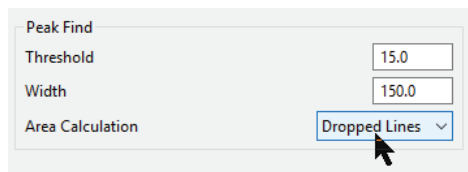
7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click **OK** to save changes.

Manual Peak Integration

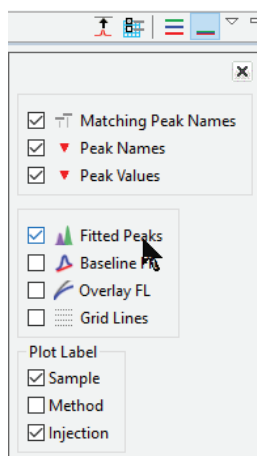
Compass for iCE lets you manually integrate peaks in individual electropherograms.

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.

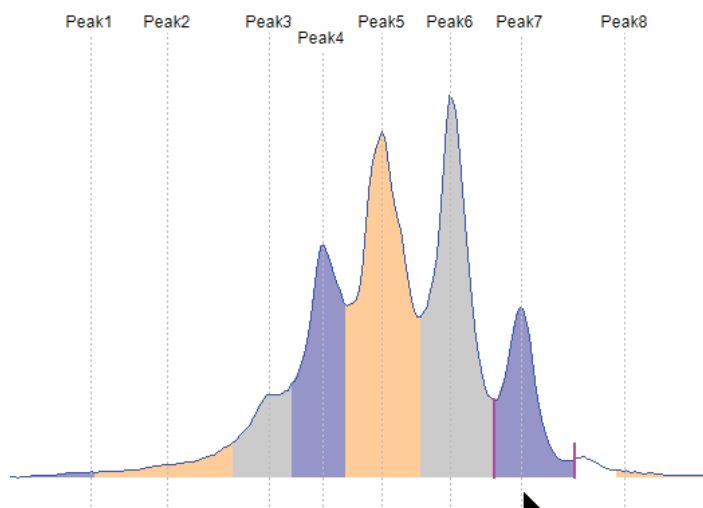
2. Select Dropped lines as the area calculation if it isn't already selected.



3. Select Fitted Peaks in the Graph Options.

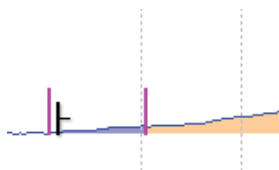



4. Select an injection in the Experiment pane.
5. Hover the mouse over the peak you want to adjust. Two magenta bars will display that indicate the start and end points of its integration.

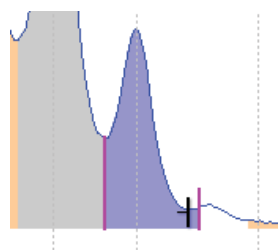


6. To change the integration, hover the mouse over either of the magenta bars. Then click the mouse and drag the bar to move it.

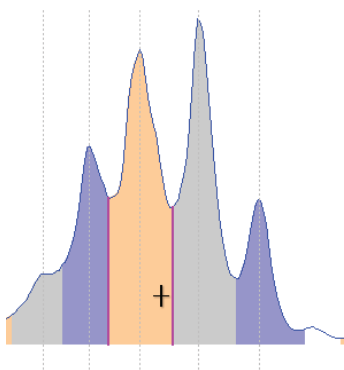
- If the cursor changes to  this is the peak start for the peak on the right.



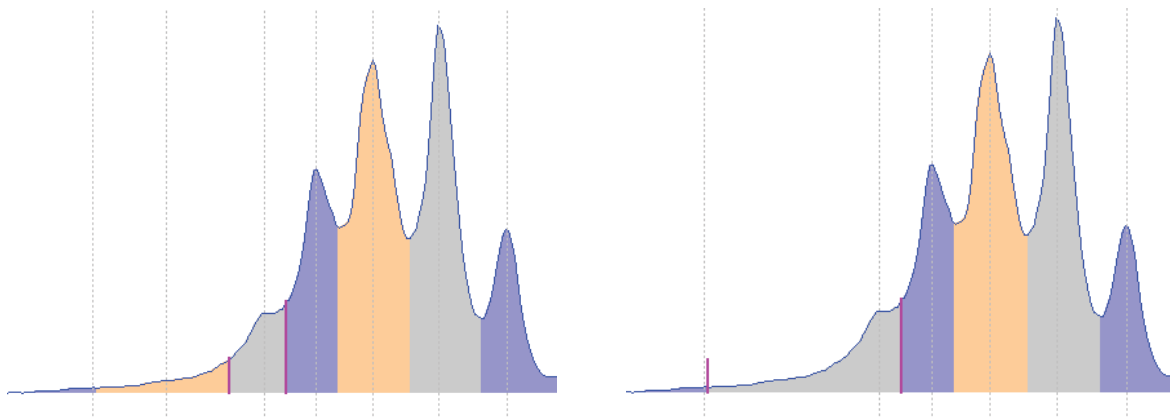
- If the cursor changes to  this is the peak end for the peak on the left.



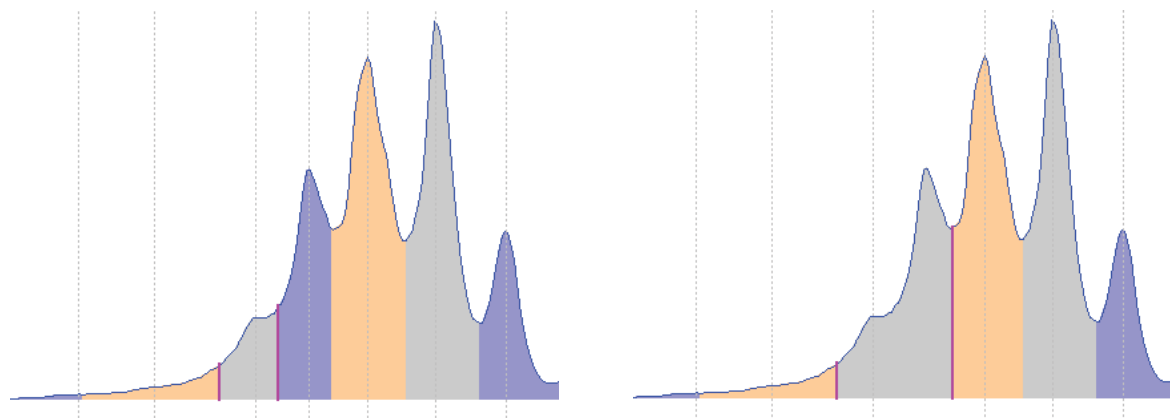
- If the cursor changes to  this is a joint boundary for the peaks on the left and right.


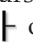
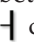


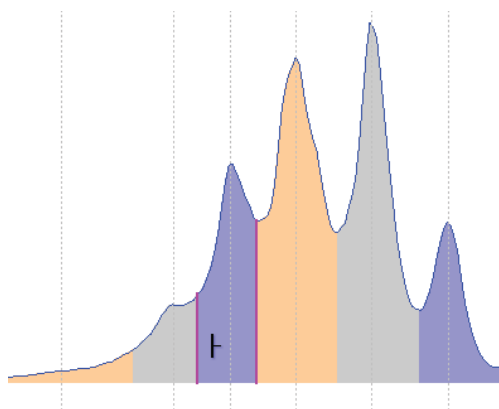
In the example below, we moved the start and end points of the peak to include more area under the peak:



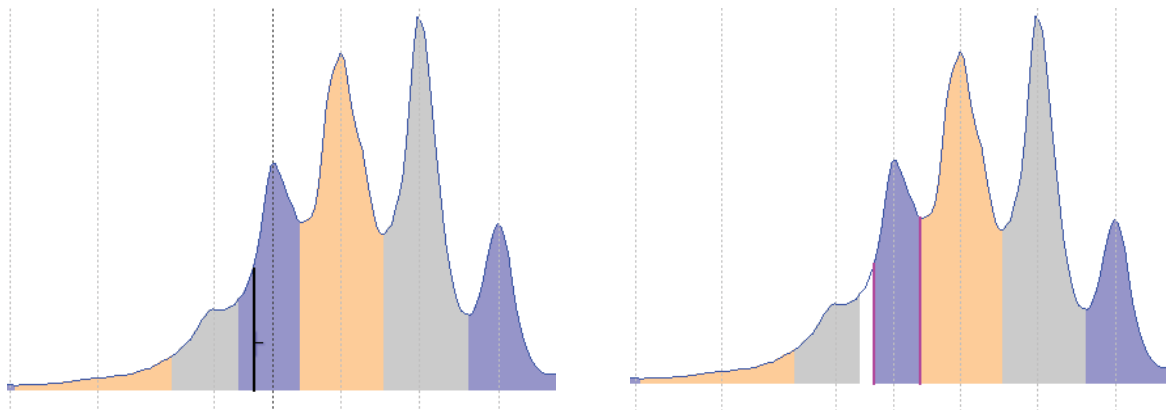
You can also combine peaks by dragging the magenta lines left or right:



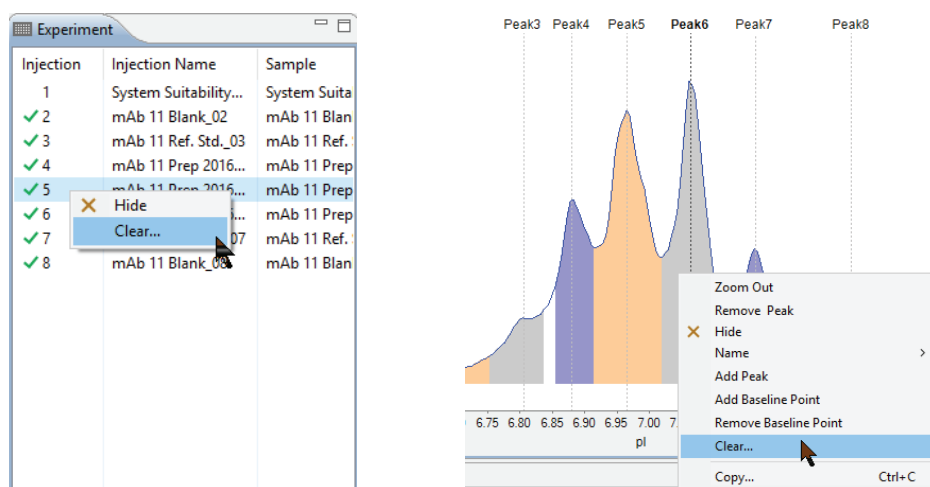
You can also separate areas between peaks. Whenever you have a  cursor between two peaks that aren't baseline resolved, move the mouse slightly to the right or left until you get the  or  cursor.



Click and hold the mouse. A black drop line will display indicating where the current peak start or end point is. While holding the mouse, drag the cursor to the right or left to move the start or end point to where you want it, then release the mouse.



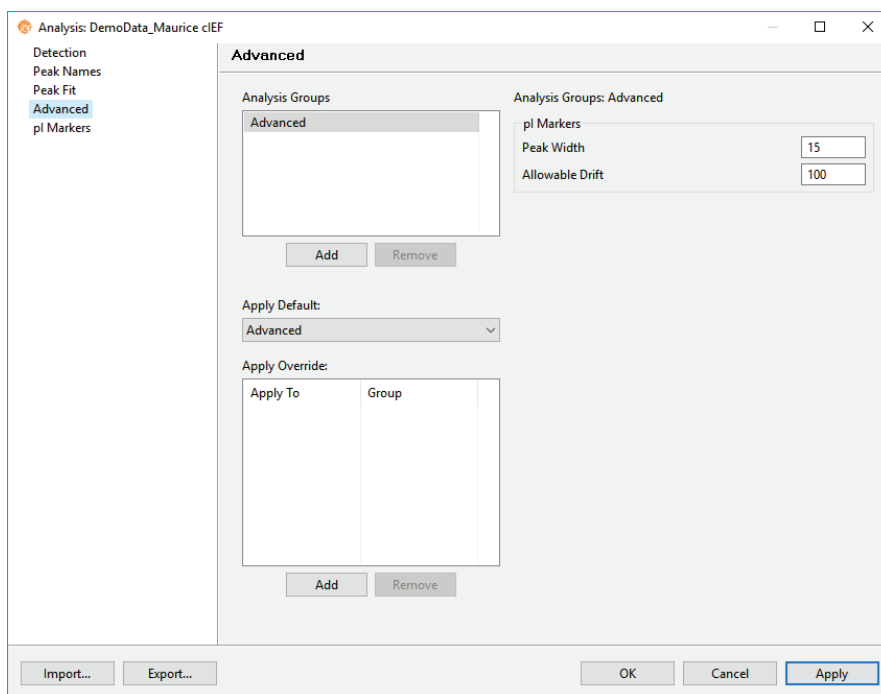
7. To remove the manual integration settings, right click on the graph or on the injection in the Experiment pane and select **Clear**.



Advanced Analysis Settings

This page lets you view and change analysis settings for the pI marker data. Select **Edit** in the main menu and click **Analysis**, then click **Advanced** in the left sidebar:

NOTE: Settings can be changed in batches before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



pI Markers Settings

- **Peak Width** - The approximate width (at full width half max) used to filter out absorbance and fluorescence artifacts which improves recognition of pI markers.
- **Allowable Drift** - The distance the pI marker(s) are expected to move compared to the position entered on the pI Markers page. This setting helps with recognition of the pI marker.

Advanced Analysis Settings Groups

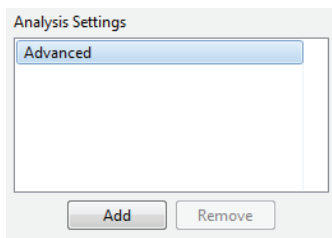
Advanced analysis settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values for advanced analysis settings. These settings are included in the default Advanced group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. See “Importing and Exporting Analysis Settings” on page 567 for more info.

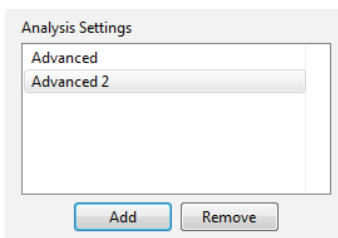
Analysis groups are displayed in the analysis settings box:



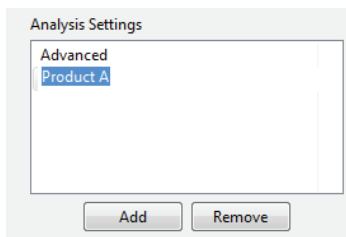
The Advanced group shown contains the Compass for iCE default analysis settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Analysis Group

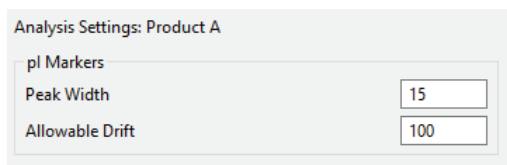
1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



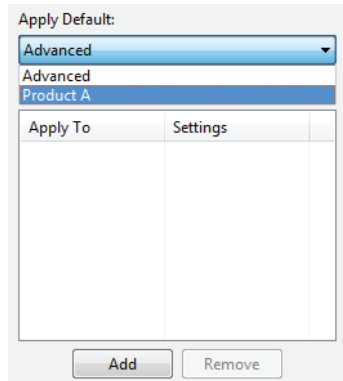
3. Click on the new group and enter a new name.



4. Change the settings in the Markers box as needed.



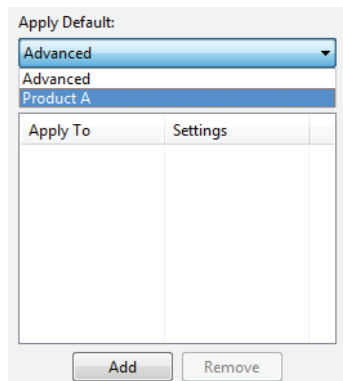
5. To use the new group as the default analysis settings for the run data, click the arrow in the drop down list next to Apply Default, then click the new group from the list. Analysis settings in the new group will then be applied to the run data.



6. Click OK to save changes.

Changing the Default Analysis Group

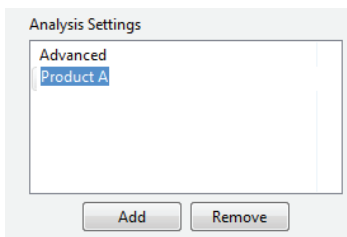
1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then click a new default group from the list.



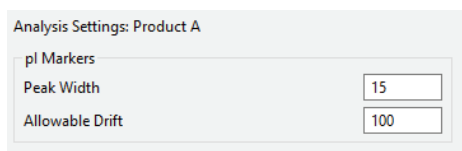
3. Click **OK** to save changes. Analysis settings in the group selected will be applied to the run data.

Modifying an Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



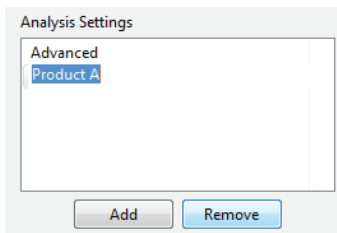
3. Change the settings in the Markers box as needed.



4. Click **OK** to save changes. The new analysis settings will be applied to the run data.

Deleting an Analysis Group

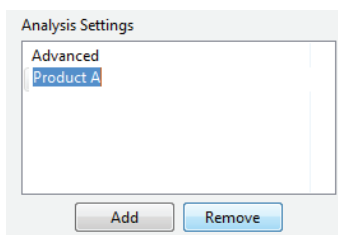
1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



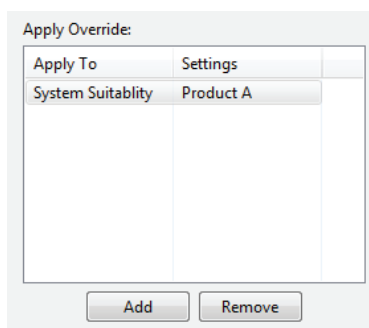
3. Click **OK** to save changes.

Applying Analysis Groups to Specific Run Data

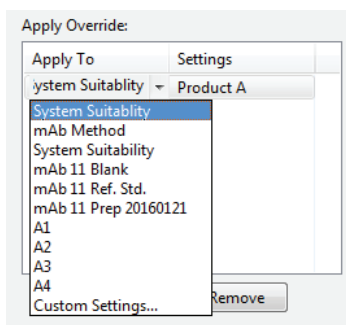
1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to apply to specific run data.



3. Application of analysis groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

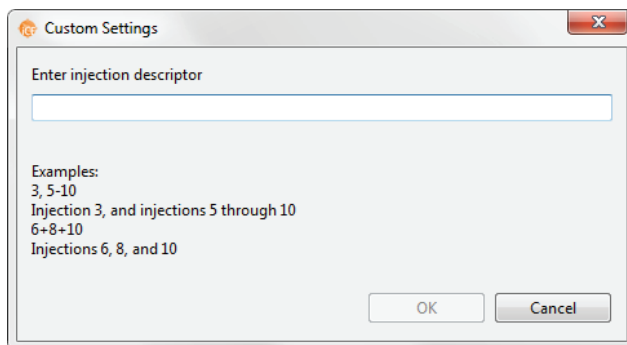


4. Click the cell in the **Apply To** column, then click the down arrow.

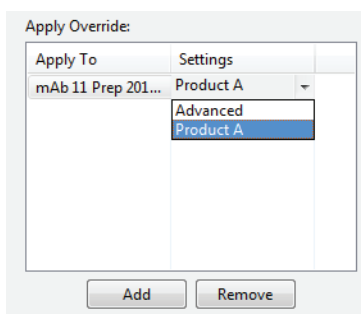


5. Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
 - **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.

- **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.
- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.

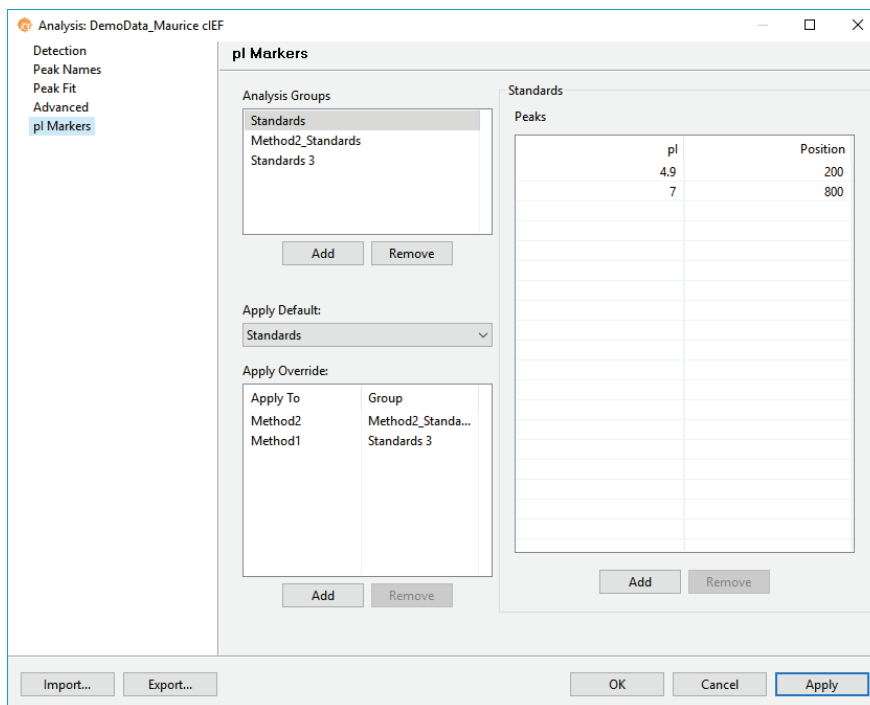


7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click **OK** to save changes.

pI Markers Analysis Settings

This page lets you define the pI and position of the pI Markers you're using in your samples. Select **Edit** in the main menu and click **Analysis**, then click **pI Markers** in the left sidebar.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



pI Markers Analysis Settings Groups

pI marker settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

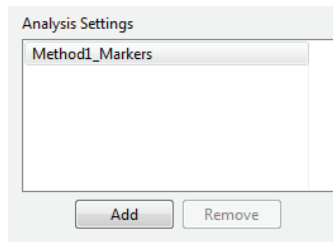
NOTES:

We recommend using the Compass for iCE default values. These settings are included in the default Markers group.

When you edit the pI markers in the method for a batch, Compass for iCE automatically creates a Markers group in the pI Markers Analysis settings for you.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 567.

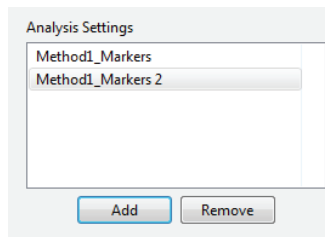
Markers groups are displayed in the analysis settings box:



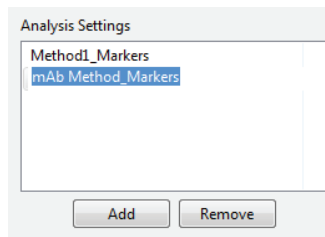
The Markers group shown uses the Compass for iCE default settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Markers Group

1. Select **Edit > Analysis**, and select **pI Markers** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:

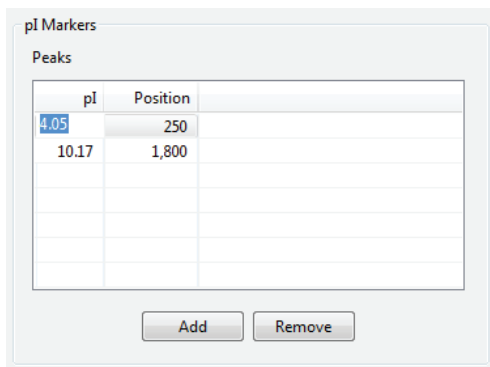


3. Click on the new group and enter a new name.



4. The default Maurice and MauriceFlex cIEF pI marker pI and position values are already populated in the pI Marker Peaks table. If you'd like to use these values, skip to the next step. If you're using different markers, here's how to change the values:

- a. Click in the first cell in the pI column in the table and enter the pI for the marker.

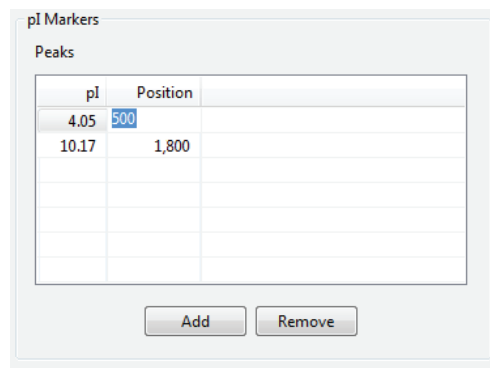


The screenshot shows a window titled "pI Markers" with a sub-header "Peaks". Below it is a table with two columns: "pI" and "Position". The first row has the values "4.05" and "250". The second row has the values "10.17" and "1,800". The "pI" cell in the first row is highlighted with a blue selection box. Below the table are two buttons: "Add" and "Remove".

pI	Position
4.05	250
10.17	1,800

Add Remove

- b. Click in the first cell in the Position column and enter a value for the marker.



The screenshot shows the same "pI Markers" window. In this step, the "Position" cell in the first row (containing "500") is highlighted with a blue selection box. The "pI" cell remains "4.05". The "Add" and "Remove" buttons are still at the bottom.

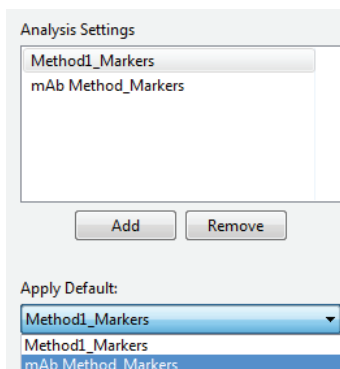
pI	Position
4.05	500
10.17	1,800

Add Remove

NOTE: pI marker peak positions are relative to each other. Only the difference in position is used to help identify them. When entering pI marker peak information for the first time, review the marker data in the Analysis screen to find the correct peak positions.

- c. Repeat the steps above for the remaining markers in the table.
- **To add another marker** - Click **Add** under the table, then change the information in the new row.
 - **To remove a marker** - Select its row and click **Remove**.

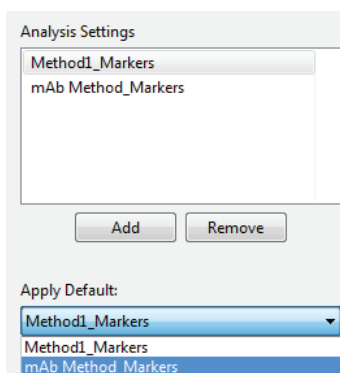
5. To use the new group as the default settings for the run, click the arrow in the drop down list next to Apply Default, then click the new group in the list. The settings in the new group will then be applied to the run data.



6. Click OK to save changes.

Changing the Default Markers Group

1. Select **Edit > Analysis**, and click **pI Markers** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then select a new default group from the list.

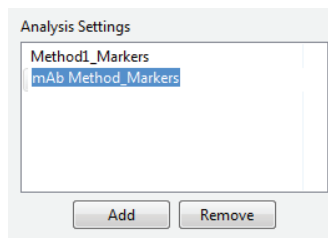


3. Click OK to save changes. Analysis settings in the group selected will be applied to the run data.

Modifying a Markers Group

1. Select **Edit > Analysis**, and click **Markers** in the left sidebar.

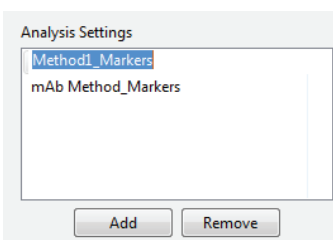
2. Click on the group in the analysis settings box you want to modify.



3. Change the marker info as needed as in Creating a New Markers Group.
4. Click OK to save changes. The new analysis settings will be applied to the run data.

Deleting a Markers Group

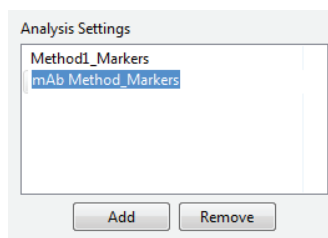
1. Select **Edit > Analysis**, and click **pI Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



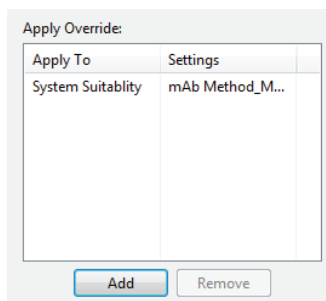
3. Click OK to save changes.

Applying Markers Groups to Specific Run Data

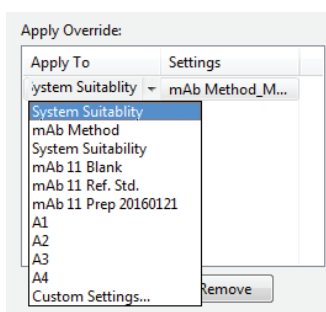
1. Select **Edit > Analysis**, and select **pI Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to apply to specific run data.



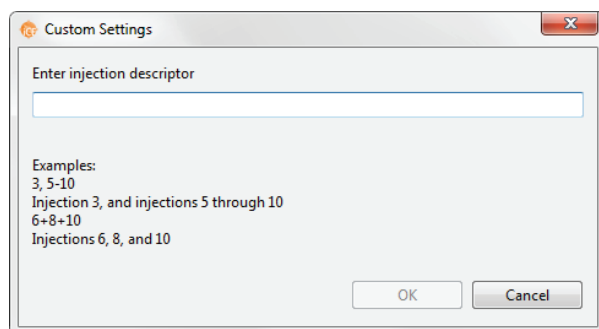
3. Application of markers groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.



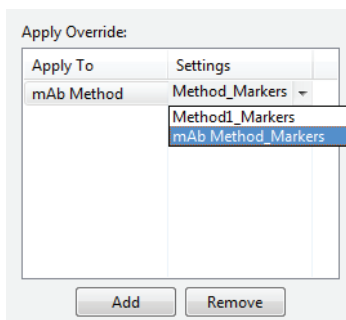
4. Click the cell in the **Apply To** column, then click the down arrow.



5. Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
- **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.
 - **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.
 - **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



- If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



- Repeat the previous steps to apply other groups to specific run data.
- To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
- Click **OK** to save changes.

Injection Reports

You can export PDF files of the raw and analyzed data, IV plot, peaks table, sample and system info for individual or all injections in a run file. You can also export the run history with all analysis events.

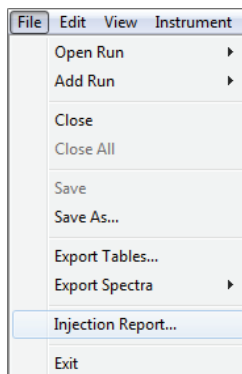
NOTES:

You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it’s saved in a separate run folder. Existing reports aren’t overwritten.

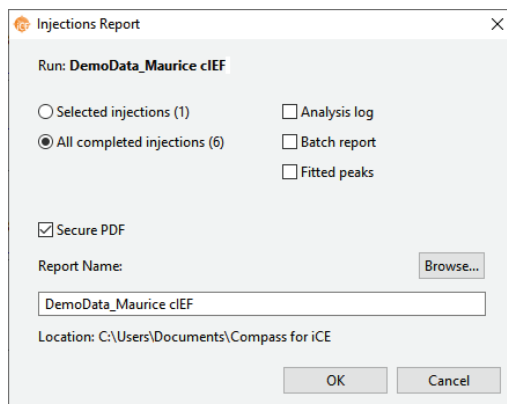
- Click **File > Open Run** and select a run file.
- If you want reports for all injections, skip to the next step. If you only want reports for certain injections, in the Experiment pane:
 - To select sequential injections:** Select the first injection, then hold the **Shift** key and select the last injection you want a report for. This selects all rows between the two injections.
 - To select specific injections:** Hold the **Ctrl** key and select just the injections you want reports for.

3. Select **File** from the main menu in either screen and click **Injection Report**.

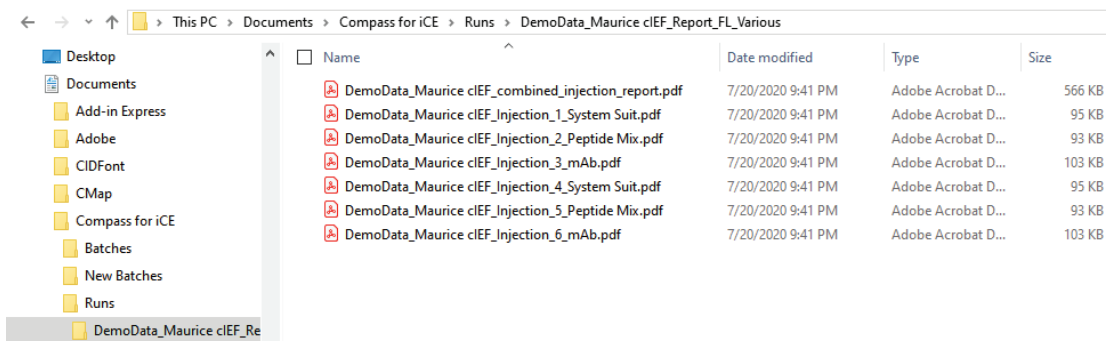


4. In the Injection Reports window:

- a. Choose either **Selected injections** or **All injections**.
- b. Select **Analysis log** if you want a run history report with all analysis events.
- c. Select **Batch Report** if you want to include the sample and method details for each injection in the batch.
- d. Select **Fitted peaks** if you want to show peak fitting in the electropherograms.
- e. Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don't want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.
- f. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.
- g. Click **OK**.

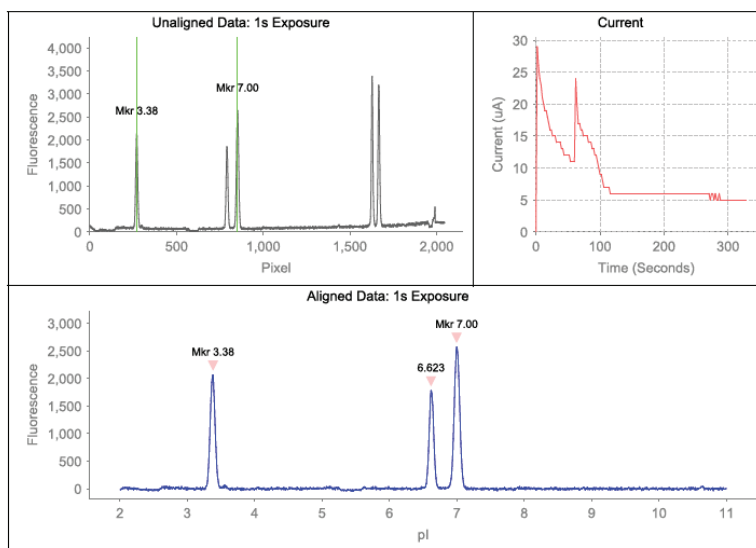


5. The individual injection report PDF(s) and a combined injection report PDF are exported to the Runs folder in the Compass for iCE directory. They'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Example Analysis and Injection Report

Uncontrolled Injection 1: System Suit_01



Fluorescence Peaks: 1s Exposure

Peak	Name	Position	pI	Height	Area	%Total	%Area	Width	S/N	Baseline	Resolution
1	Mkr 3.38	270.1	3.380	2067.9	28502.1			0.0807	112.6	52.7	
2		791.0	6.623	1785.9	22868.7	100.00		0.0750	97.2	41.1	24.56
3	Mkr 7.00	851.9	7.000	2580.0	39590.9			0.0898	140.5	41.0	2.71

Uncontrolled Injection 1: System Suit_01

Sample Information

Injection Name	System Suit_01
Sample ID	System Suit
Location	Plate Well A1
Batch Name	Maurice cIEF
Run Started	Tue 12:11 PM Oct 27, 2015 PDT
Run Completed	Tue 1:17 PM Oct 27, 2015 PDT
Date Acquired	Tue 12:17 PM Oct 27, 2015 PDT
Run Error	None
Reinjection	No

Injection Conditions

Guaranteed Injection	Yes
On-board Mixing	Off
Focus Period 1	1500V for 1.0 min
Focus Period 2	3000V for 4.5 min
Sample Load Duration	90.0 Seconds
pI marker 1	3.38
pI marker 2	7.0
Tray Temperature	Not Available

Maurice Settings

Model	Maurice OBM
Instrument S/N	kf1004
Software Version	Compass for iCE 4.0.0, Build ID: 0222
Firmware Version	2.0.2015.10.20.20.35.52.681a5af
Tray Type	48 vials
Cartridge Type	
Cartridge S/N	
Cartridge Expiration	
Injections Remaining	
Batches Remaining	

Importing and Exporting Analysis Settings

The analysis settings in a run file can be exported as a separate file. This allows the same analysis settings to be imported into other batches or run files at a later time, rather than you having to re-enter them manually.

Importing Analysis Settings

NOTE: Importing an analysis settings file populates the settings in all analysis pages.

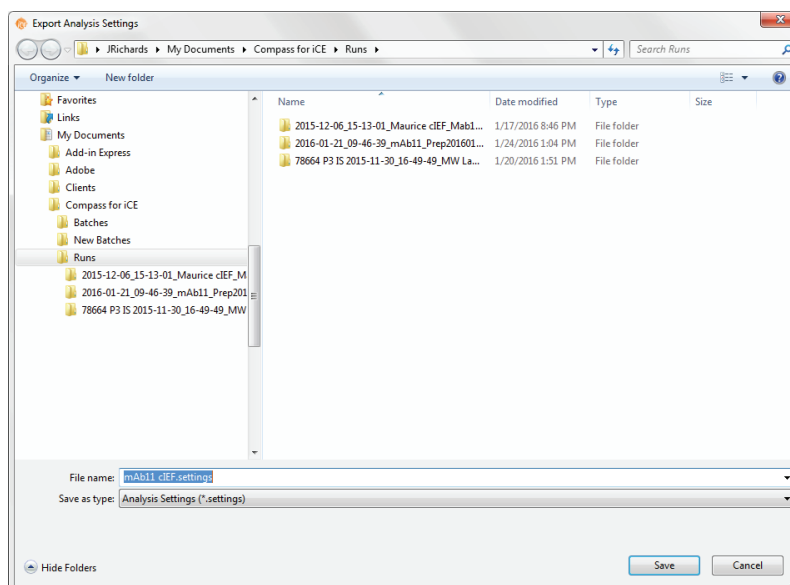
1. Open the run file or batch you want to import analysis settings to.
2. Select **Edit** in the main menu and click **Default Analysis** (Batch screen) or **Analysis** (Analysis screen).
3. Click **Import** on any page.

4. Select a settings file (*.settings) and click OK. The imported settings will display in all analysis pages.

Exporting Analysis Settings

NOTE: Exporting an analysis settings file exports the settings in all analysis pages.

1. Open the run file or batch you want to export analysis settings from.
2. Select **Edit** in the main menu and click **Default Analysis** (Batch screen) or **Analysis** (Analysis screen).
3. Click **Export** on any page. The following window displays:



4. The default directory is Compass for iCE/Runs. Change the directory if needed.
5. Enter a file name and click **Save**. The settings will be saved as a *.settings file.

Chapter 19:

MauriceFlex Fractionation Data Analysis

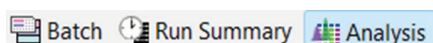
Chapter Overview

- Analysis Screen Overview
- Opening Run Files
- How Run Data is Displayed
- Viewing Run Data
- Data Notifications and Warnings
- Checking Your Results
- Naming Peaks
- Copying Results Tables and Graphs
- Exporting Run Files
- Changing Sample Protein Identification
- Adding or Removing Fractions Data
- Changing the Electropherogram View
- Closing Run Files
- Analysis Settings Overview
- Detection Settings
- Peak Names Settings
- Peak Fit Analysis Settings
- Manual Peak Integration
- Advanced Analysis Settings
- pI Markers Analysis Settings
- Injection Reports
- Importing and Exporting Analysis Settings

NOTE: To learn more how to analyze data from a cIEF run using the fractions from a MauriceFlex Fractionation batch, check out the MauriceFlex cIEF Fractionation Method Development Guide.

Analysis Screen Overview

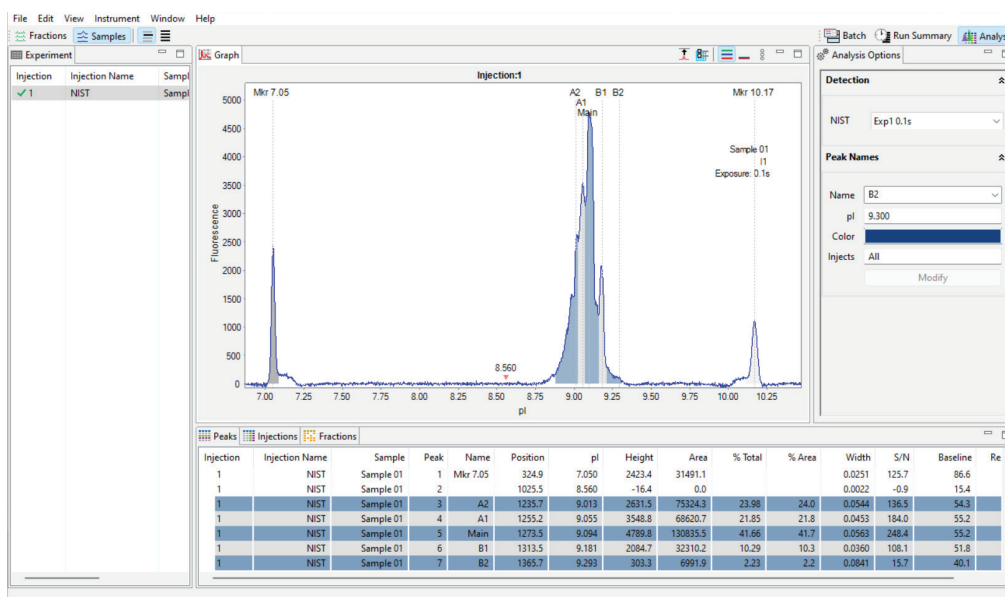
You can use the Analysis screen to view electropherograms and tabulated results for your injections. If any post-run analysis is needed, you can do it here too. To get to this screen, click the **Analysis** screen tab:



Analysis Screen Panes

The Analysis screen has four panes:

- **Experiment** - Lists the injection number, sample IDs, sample locations and methods for each injection in the run and lets you get a quick view of method parameters.
- **Graph** - Displays the electropherograms for sample proteins or pI markers.
- **Peaks** - Shows the tabulated results for sample proteins and pI markers.
- **Injections** - Displays a list of the sample proteins Compass for iCE names automatically using the user-defined peak name analysis parameters.
- **Fractions** - Displays the fractions that each sample peak was collected
- **Analysis Options** - Lets you view absorbance or fluorescence data for the run and view, change and add new custom peak name settings.



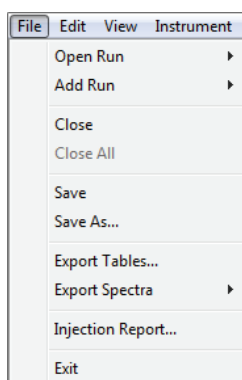
Software Menus Active in the Analysis Screen

These main menu items are active in the Analysis Screen:

- File
- Edit
- View
- Instrument (when Compass for iCE is connected to MauriceFlex.)
- Window
- Help

File Menu

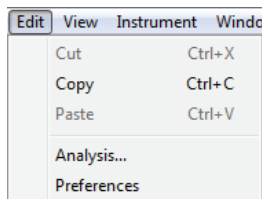
These File menu options are active:



- **Open Run** - Opens a run file.
- **Add Run** - Lets you open and view other run files besides the one that's already open.
- **Close** - Closes the run file currently being viewed.
- **Close All** - Closes all open run files.
- **Save/Save As** - If you made changes in the Analysis screen before you went to the Run Summary screen, this saves your changes to the run file.
- **Export Tables** - Exports the results for all injections in the run in .txt format.
- **Export Spectra** - Exports the raw and analyzed data traces and background for each injection in the run in .txt or .cdf format.
- **Injection Report** - Exports the raw and analyzed data, IV plot, peaks table, sample and system info for individual injections as PDF files. You can also export the run history with all analysis events.
- **Exit** - Closes Compass for iCE.

Edit Menu

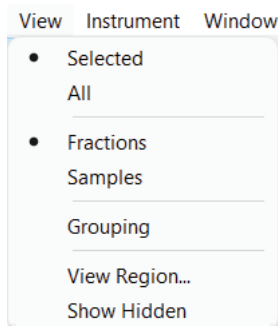
These Edit menu options are active:



- **Copy** - Lets you copy data shown in the graph, lane, peaks or injections panes. See “Copying Results Tables and Graphs” on page 594 for more information.
- **Analysis** - Displays the analysis settings used to analyze the run data and lets you change them as needed. See “Analysis Settings Overview” on page 620 for more information.
- **Preferences** - Lets you set and save your preferences for data export, graph colors, grouped data and Twitter settings. See Chapter 21: “Setting Your Preferences” for more information.

View Menu

These View menu options are active:



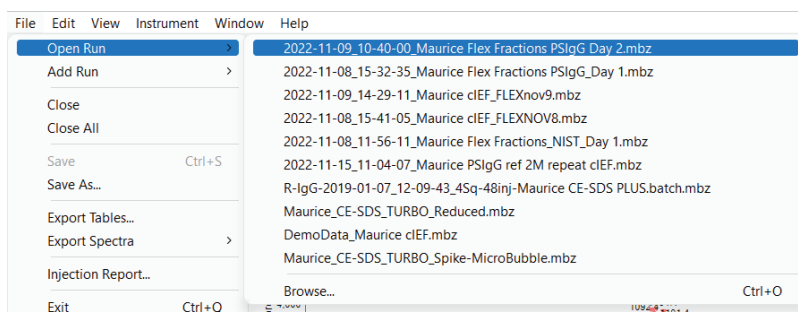
- **Selected (Single View)** - Displays the data for only the injections selected.
- **All (Multiple View)** - Displays data for all injections so you can scroll through them.
- **Fractions** - Lets you view and set pI markers for the sample injection and mobilization electropherograms where you can view the protein mobilizing out of the capillary.
- **Samples** - Lets you view and name peaks for the focused sample injection.
- **Grouping** - Displays data for injection groups.
- **View Region** - Lets you change the x-axis range of the data displayed.
- **Show Hidden** - Shows injections that are hidden from the data view.

Opening Run Files

You can open one run file or multiple files at the same time to compare information between runs.

Opening One Run File

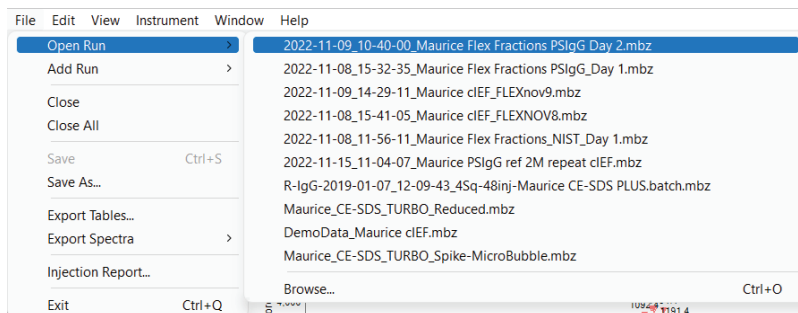
1. Select **File** in the main menu and click **Open Run**.



2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.

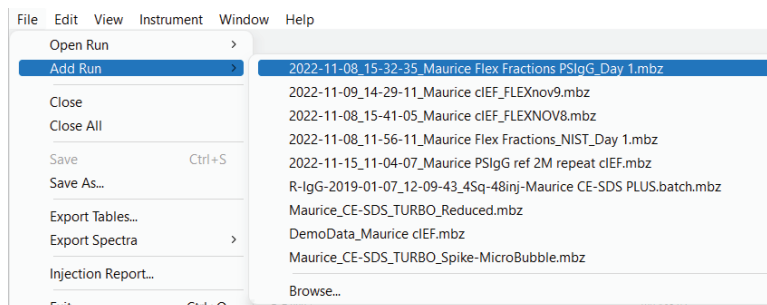
Opening Multiple Run Files

1. To open the first run file, select **File** in the main menu and click **Open Run**.

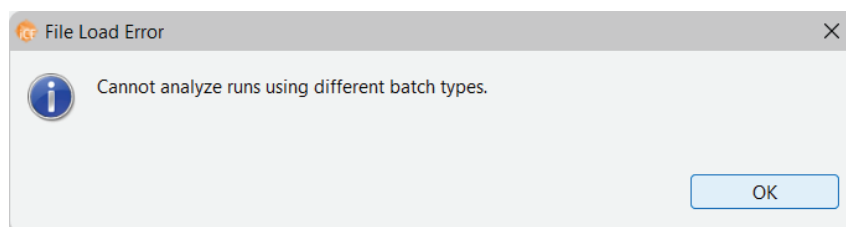


2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.

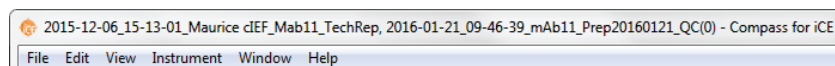
3. To open another run file, select **File** in the main menu and click **Add Run**.



4. A list of MauriceFlex Fractionation runs will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file. If the run selected is not a MauriceFlex Fractionation run, an alert will appear.



5. When a run is added, its data appends to the open run file and displays as a second set of injections in all screen panes. The second run file name also appears in the title bar:



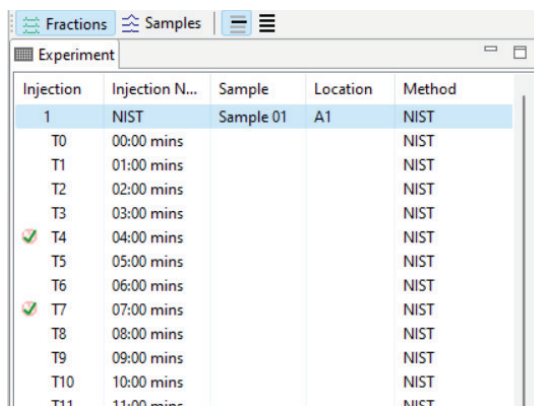
6. Repeat the last two steps to add additional runs.

How Run Data is Displayed

Data in the run file is organized for easy review.

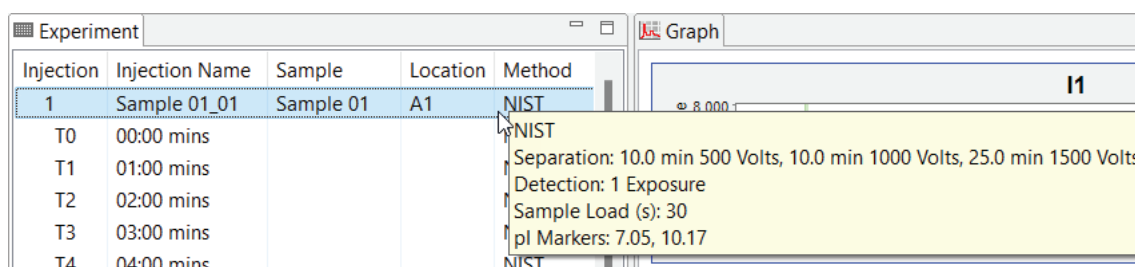
Experiment Pane: Batch Injection Information

The Experiment pane lists the injection performed followed by the fractions collected in the run. Injection information includes the sample used, the sample location and the method used when the Samples view is selected. Fractionation information includes an image number collected during the mobilization series and the time the mobilization electropherogram image was taken and can be viewed when the Fractions view is selected.



Injection	Injection N...	Sample	Location	Method
1	NIST	Sample 01	A1	NIST
T0	00:00 mins			NIST
T1	01:00 mins			NIST
T2	02:00 mins			NIST
T3	03:00 mins			NIST
✓ T4	04:00 mins			NIST
T5	05:00 mins			NIST
✓ T6	06:00 mins			NIST
✓ T7	07:00 mins			NIST
T8	08:00 mins			NIST
T9	09:00 mins			NIST
T10	10:00 mins			NIST
T11	11:00 mins			NIST

- **To view all columns** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.
- **To view method parameters** - Hover the mouse over a method name.



Injection	Injection Name	Sample	Location	Method
1	Sample 01_01	Sample 01	A1	NIST
T0	00:00 mins			
T1	01:00 mins			
T2	02:00 mins			
T3	03:00 mins			
T4	04:00 mins			

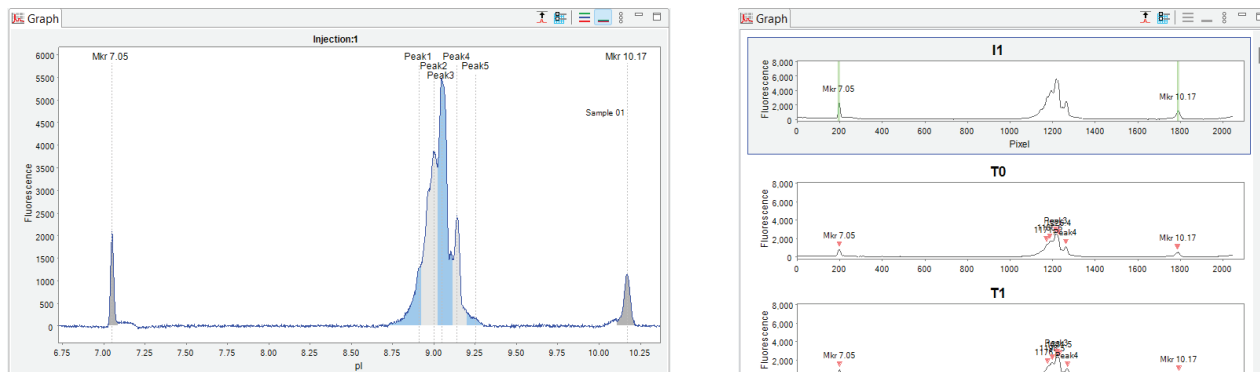
NIST
 Separation: 10.0 min 500 Volts, 10.0 min 1000 Volts, 25.0 min 1500 Volts
 Detection: 1 Exposure
 Sample Load (s): 30
 pl Markers: 7.05, 10.17

NOTE: Data notification icons will display in the Injection column if Compass for iCE detects a potential analysis issue or data was manually modified by the user. For more information see “Data Notifications and Warnings” on page 411.

Graph Pane: Electropherogram Data

The Graph pane displays sample and mobilization electropherograms. When the Fractions view is selected, the sample injection appears first (I1) followed by mobilization electropherograms (T0, T1, T2, etc). The electropherogram viewed depends on the view options you’ve selected.

You can get more info on graph view options in “Changing the Electropherogram View” on page 601.



Peaks Pane: Calculated Results

The Peaks pane shows the tabulated results for your sample proteins or pI markers. Each row in the table has the individual results for each peak detected in an injection. Results shown will either be for one injection or multiple injections, samples or pI markers depending on the view options you’re using. Check out “Analysis Options Pane” on page 580 for more info.

Peaks														
Injection	Injection N...	Sample	Peak	Name	Position	pI	Height	Area	% Total	% Area	Width	S/N	Baseline	Resoluti...
1	Sample 01_...	Sample 01	1	Mkr 7...	196.8	7.050	2089.8	24085.7			0.0211	96.9	153.2	
1	Sample 01_...	Sample 01	2	Peak1	1146.5	8.910	1371.7	43234.2	10.19	10.2	0.0446	63.6	60.4	33.32
1	Sample 01_...	Sample 01	3	Peak2	1193.3	9.002	3882.5	14723...	34.71	34.7	0.0777	180.1	58.6	0.88
1	Sample 01_...	Sample 01	4	Peak3	1216.5	9.047	5489.5	17228...	40.61	40.6	0.0603	254.6	56.6	0.39
1	Sample 01_...	Sample 01	5	Peak4	1263.5	9.139	2435.5	53548.1	12.62	12.6	0.0423	113.0	51.5	1.06
1	Sample 01_...	Sample 01	6	Peak5	1322.8	9.255	297.1	7898.0	1.86	1.9	0.0577	13.8	45.0	1.37

NOTES:

Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

When the Fractions view is selected, the information in the Peaks table includes only injection, injection name, sample, peak, name, position and height. pI markers the software has identified are marked with an **M**.

- **To view all rows** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.

The following results and info are listed in the Peaks table:

- **Injection** - Injection number.
- **Injection name** - If injection names were entered in the batch, those names will display here. Otherwise the default 'SampleID_injection number' name will display.
- **Sample** - If sample names were entered in the batch, those names will display here. Otherwise, Sample (default name) will display.
- **Peak** - Peaks are numbered in order of detection.
- **Name** - Displays peaks Compass for iCE named automatically using the user-defined peak name analysis parameters. These cells are blank if the software wasn't able to name the peak or if you didn't enter naming parameters.
- **Position** - Peak location in pixels.
- **pI** - Displays the calculated peak pI based on the position of the peak relative to the pI markers.
- **Height** - The calculated peak height.
- **Area** - Displays the count of the pixel values for dropped line fit and the area of the curve fit for gaussian fit.
- **% Total** - Displays the peak area ratio compared to the sum of all peak areas. This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.
- **% Area** - Displays the calculated percent area for the named peak compared to all named peaks. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100 (shown for named peak sample data only).
- **Width** - Displays the calculated peak width (sample data only).
- **S/N** - Displays the calculated signal to noise. The signal to noise ratio is reported for each peak based on the noise in the injection and the height of the peak. $S/N = 2 \times \text{peak height} / \text{noise}$. For the injection noise calculation, the peak to peak noise is calculated by fitting a quadratic to the test region and taking the difference between the max signal above the fit, and the minimum signal below the fit. A fit is used to mitigate the effect of drift in the signal that would give a falsely high noise value. The software looks for the quietest 50 point region in the entire injection excluding 50 points at each end. The test region is approximately 5 times a typical peak width (FWHM).
- **Baseline** - Displays the raw baseline signal of each peak.
- **Resolution** - Displays resolution of the peak compared to neighboring peaks. Two peaks that are baseline resolved will have a resolution value of 1.5. Smaller values means the peaks are not completely resolved, larger values mean the peaks are fully resolved.

Injections Pane: User-Specified Peak Names

The Injections pane shows tabulated results for sample proteins Compass for iCE labels automatically using user-defined peak name settings. Each row in the table shows the individual results for the named peaks detected in each injection.

Injection	Injection N...	Sample	Peak1	Peak2	Peak3	Peak4	Peak5
1	Sample 01...	Sample 01	43234.2	147238.8	172281.2	53548.1	7898.0

NOTES:

Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

When the Fractions view is selected, the information in the Injections table includes only injection, sample and the positions of the pI marker (Mkr) peaks.

- **To view all rows** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.

The following results and info are listed in the Injections table:

- **Injection** - Injection number.
- **Injection name** - If injection names were entered in the batch, those names will display here. Otherwise the default 'SampleID_injection number' name will display.
- **Sample** - If sample names were entered in the batch, those names will display here. Otherwise, Sample (default name) will display.
- **Peak Name Columns** - An individual column per peak name will display for every peak identified by name or as a pI marker peak in the run data. Cells for injections in these columns will be blank if Compass for iCE didn't find peaks automatically using the user-defined peak name analysis and maker parameters (or none were entered).
 - **To view peak area in the peak name columns (default)** - Select **Area** in the upper right corner of the pane. This displays calculated peak area for the individual peak only.
 - **To view % total in the peak name columns** - This displays the peak area ratio compared to the sum of all peak areas. This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.

NOTE: The sum of the named peak percentages can be less than 100% if some peaks aren't named.

Peaks Injections Fractions								% Total	Area	% Area
Injection	Injection N...	Sample	Peak1	Peak2	Peak3	Peak4	Peak5			
1	Sample 01_...	Sample 01	10.2	34.7	40.6	12.6	1.9			

- **To view % area in the peak name columns** - This displays the peak area ratio compared to the sum of all named peak areas. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100.

Peaks Injections Fractions								% Total	Area	% Area
Injection	Injection N...	Sample	Peak1	Peak2	Peak3	Peak4	Peak5			
1	Sample 01_...	Sample 01	10.2	34.7	40.6	12.6	1.9			

Fractions Pane: Fraction Peak Predictor

The Fractions Pane appears when the Samples view is selected. It predicts where each sample peak was collected in the 96-well plate. Compass for iCE uses a peak prediction algorithm to model the mobility of sample peaks in the capillary. Results will either be shown for pI markers or named peaks, depending on the view options you're using. Check out the "Analysis Options Pane" on page 580 for more info.

Peaks Injections Fractions										
Injection	Injection Name	Location	Mkr 7.05	8.560	A2	A1	Main	B1	B2	Mkr 10.17
1	NIST	B7							B7	
1	NIST	B8							B8	
1	NIST	B9							B9	
1	NIST	B10					B10	B10	B10	
1	NIST	B11			B11	B11	B11	B11	B11	
1	NIST	B12			B12	B12	B12	B12		
1	NIST	C12			C12	C12	C12			
1	NIST	C11			C11	C11				
1	NIST	C10		C10	C10					
1	NIST	C9								
1	NIST	C8								
1	NIST	C7								
1	NIST	C6								
1	NIST	C5								

NOTES:

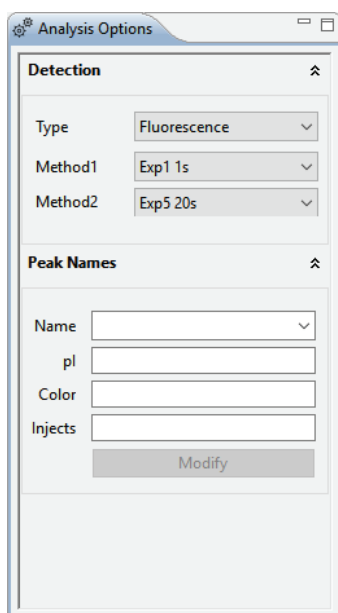
Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

Peak pI values will be used if you have not named peaks yet.

Analysis Options Pane

The Analysis Options pane gives you a quick way to view data for different exposures and add peak names without having to open and edit the run's analysis settings.

- **Peak names** - Lets you view, change and add new custom peak name settings for sample proteins. You can apply peak name settings to all, selected, or specified injections. For details on how to use this option, go to "Naming Peaks" on page 589.



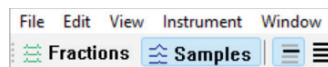
Viewing Run Data

The Analysis screen lets you view one electropherogram, specific electropherograms or all electropherograms in the run. Each run file has data for the sample peaks and pI markers detected. Switching Between Samples and Markers Data Views

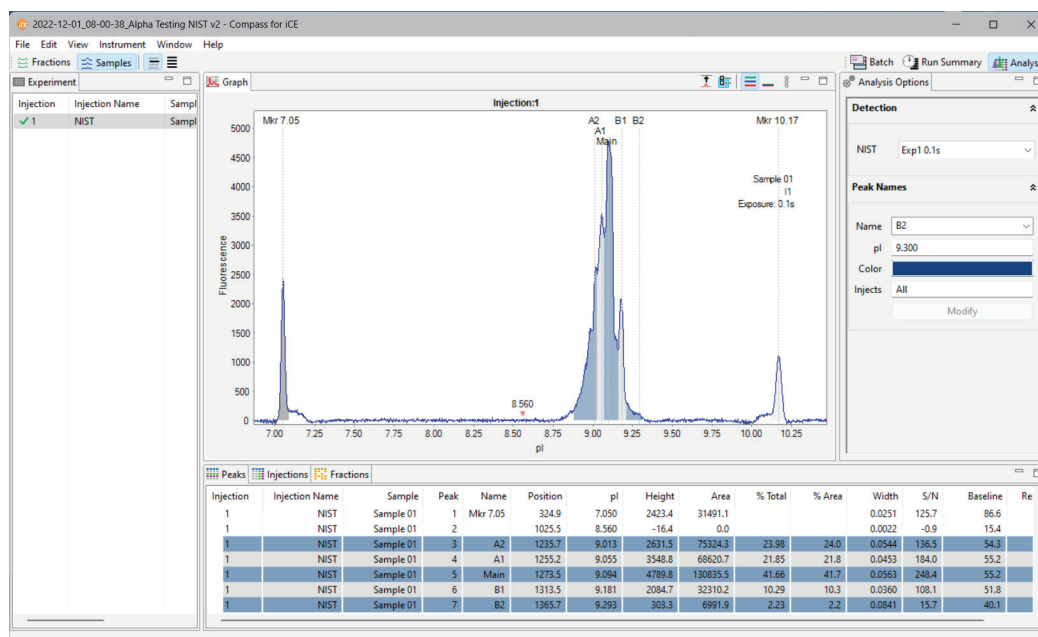
Switching Between Samples and Fractions Data Views

Here's how you switch between viewing data for your sample and pI markers in your final focusing image and viewing mobilization electropherograms:

- To view sample data - Click **Samples** in the View bar or select **View** in the main menu and click **Samples**.

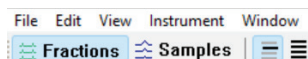


- Data in this view is for the final focusing sample image only.
- The graph displays electropherograms with a y-axis of Fluorescence units and a x-axis of pI.
- Results for each peak are shown in the Peaks and Injections panes. The prediction of which fraction each sample peak was collected in are shown in the Fractions pane.



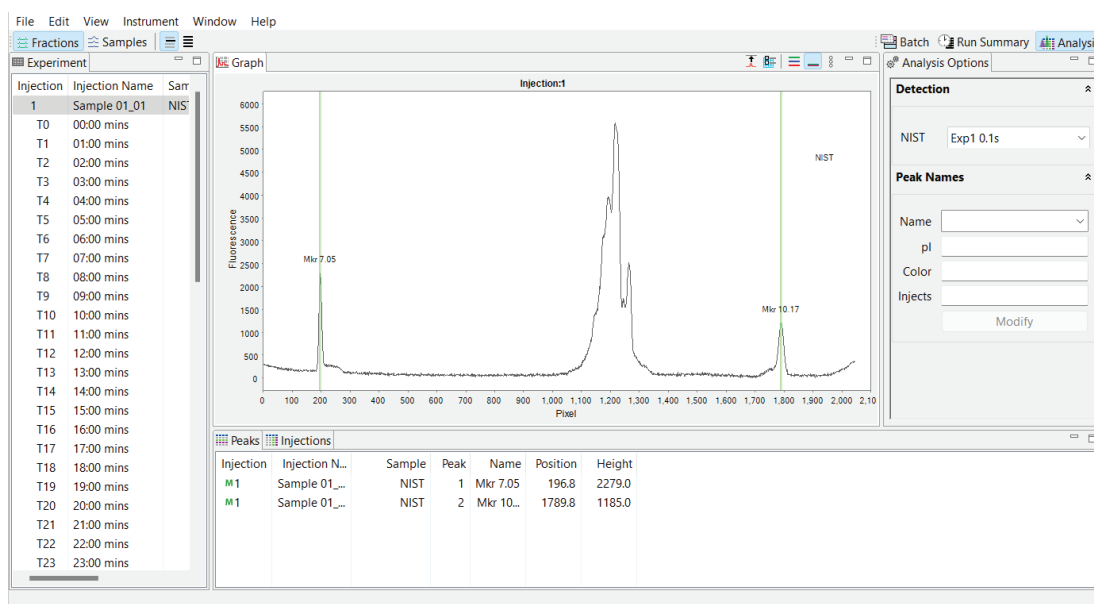
For information on checking and identifying sample peaks, see “Checking Your Data” on page 250.

- To view pI marker data for your sample (I1) - Click **Fractions** in the View bar or select **View** in the main menu and click **Fractions**.



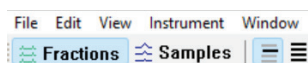
Data in this view is for analyzing pI markers for the final focusing sample image. These are the pI markers you add to your sample during prep. pI markers cannot be adjusted in Samples view. pI markers for mobilization electropherograms cannot be adjusted.

- The graph displays electropherograms with a y-axis of Fluorescence units and an x-axis of pixels.
- pI markers are identified in the Peaks pane with an **M** and as Mkr in the Injections pane.



For information on checking and identifying the pI marker peaks, see “Checking Your Data” on page 250.

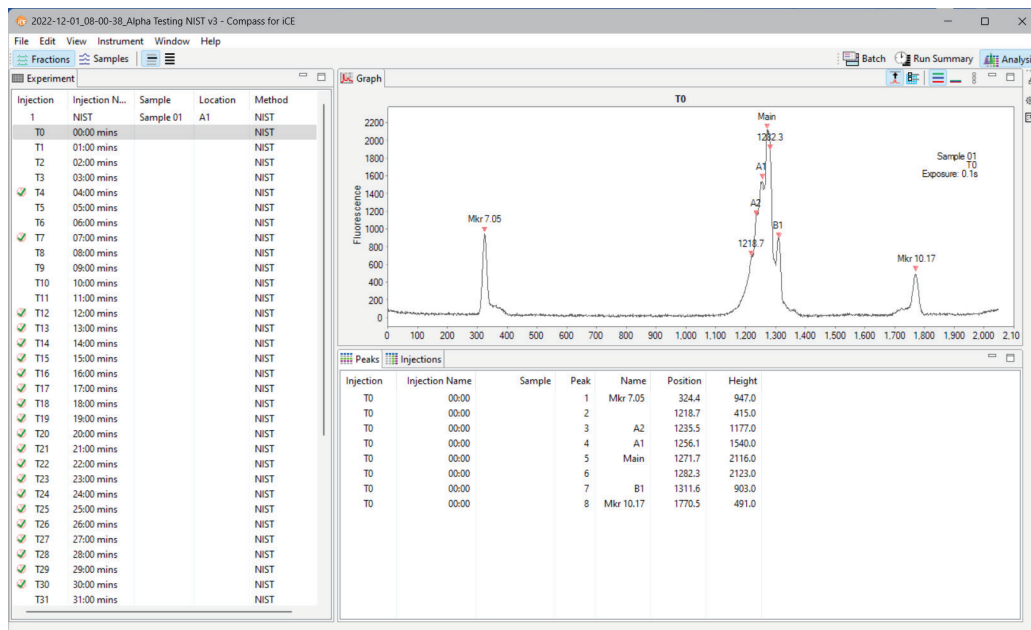
- **To view Fractions data (T0, T1, T2, etc):** Click **Fractions** in the View bar or select **View** in the main menu and click **Fractions**.



Data in the view is for confirming peaks have been correctly identified in the mobilization electropherograms.

- The graph displays electropherograms with a y-axis of Fluorescence units and an x-axis of pixels.
- Incorrectly identified peaks can be removed and missing peaks can be added.
- Peaks will be identified by pixel position by default.
- Compass for iCE will automatically name peaks in the Fractions view using Sample peak names defined by the user in the Samples view. Peak names can be manually adjusted in the Fractions view.

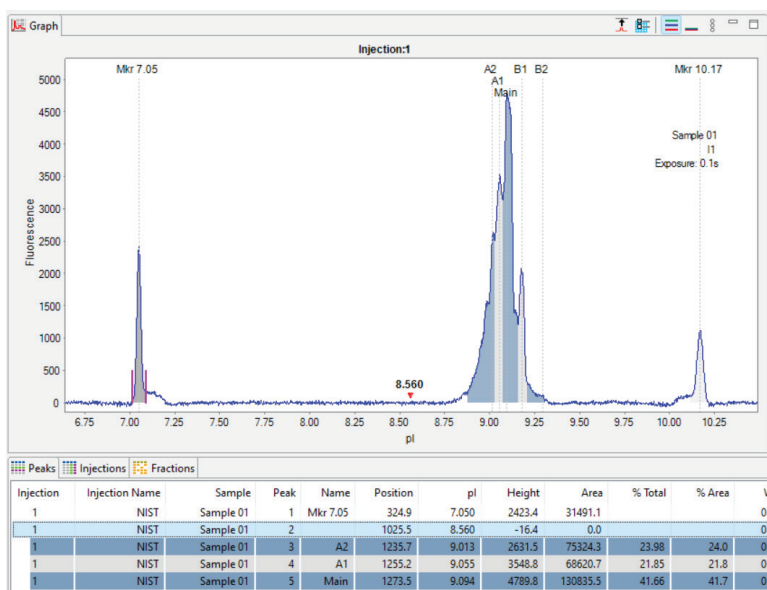
- Information for each peak are shown in the Peaks and Injections panes.

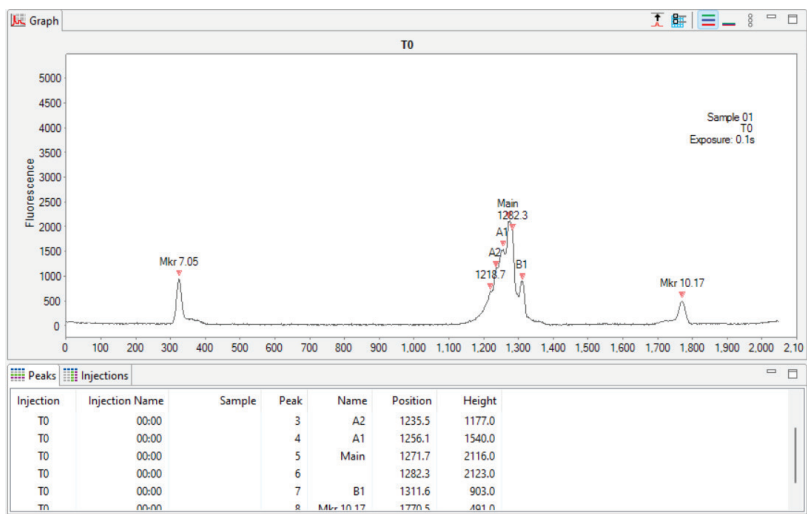


Selecting and Displaying Injection Data

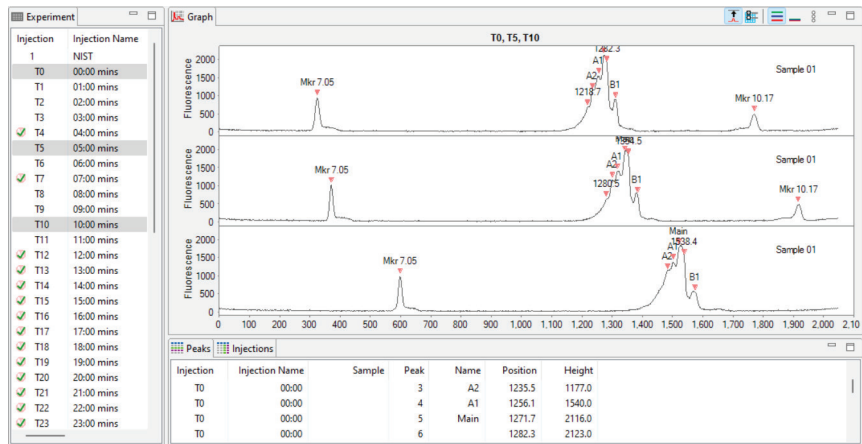
You can view the sample injection and name peaks in the Samples View. You can view the pI markers, injection and mobilization data for the sample and collected fractions in the Fractions View.

- To look at data for one electropherogram** - Select the injection or mobilization electropherogram in the Experiment pane. Data for just that electropherogram displays in the graph and tables.

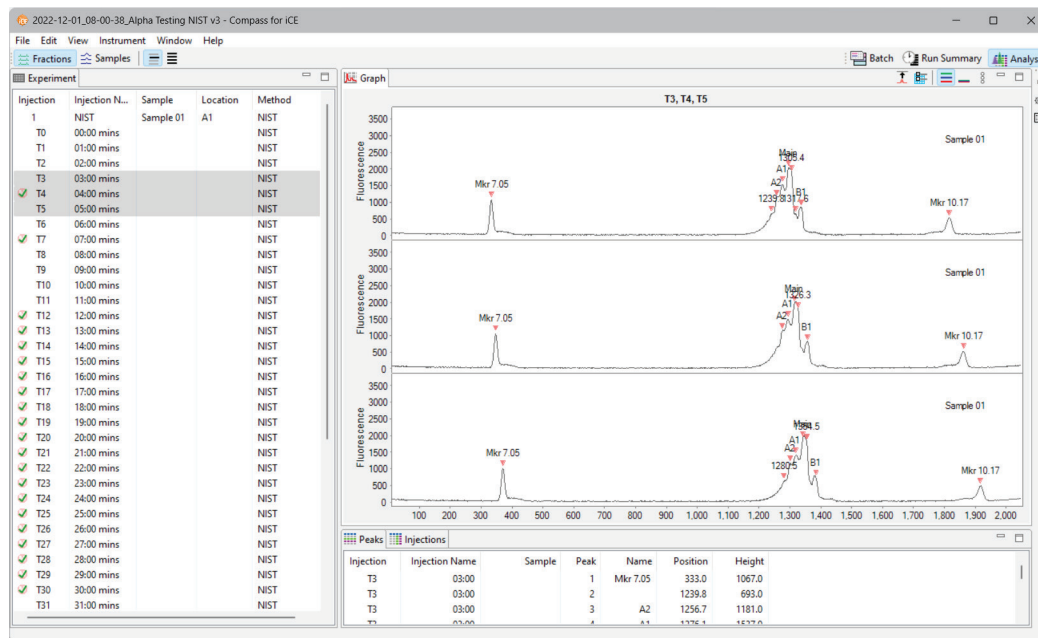




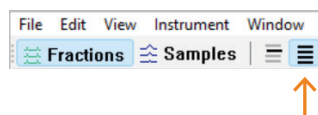
- **To look at data for specific mobilization electropherograms** - In the Fractions view, hold the Ctrl key and select just the rows you want to view in the Experiment pane. Data for only the rows selected display in the graph and tables.



- To look at data for sequential mobilization electropherograms - In the Fractions view, select the first row in the Experiment pane that you want to view, then hold the **Shift** key and select the last. This selects all rows between the two rows. Data for only the rows selected display in the graph and tables.



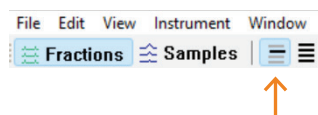
- To look at data for all mobilization electropherograms - Just click **View All** in the View bar. Data for all rows displays in the graph and tables.



Switching Between Selected and All Electropherogram Views (Fractions View)

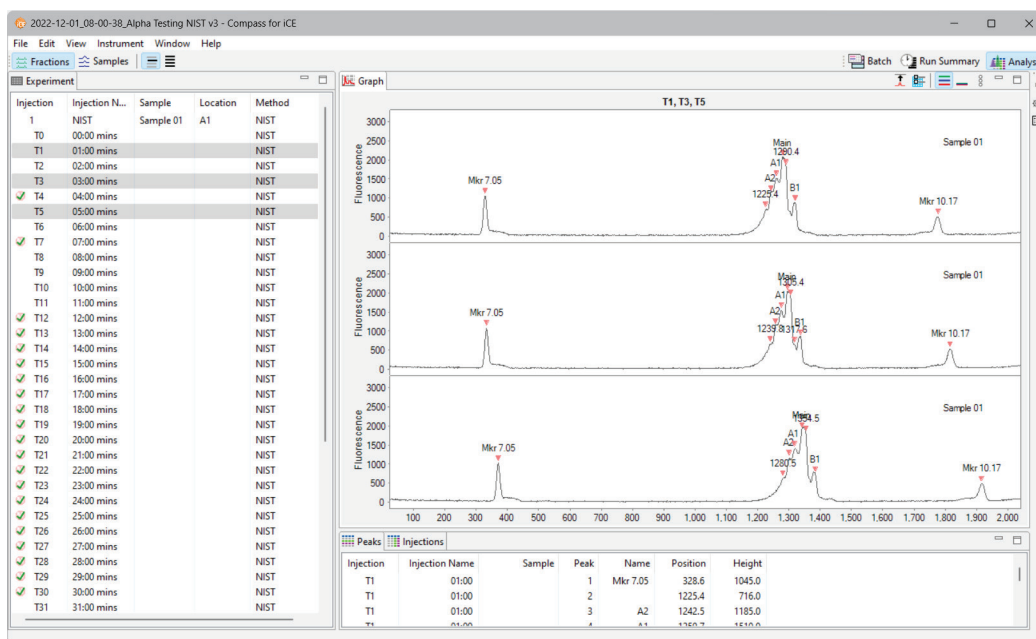
You can switch between displaying run data in a single electropherogram format or a multi-electropherogram format.

- To view data in a selected electropherogram format - Click **Selected View** in the View bar or select **View** in the main menu and click **Selected**.

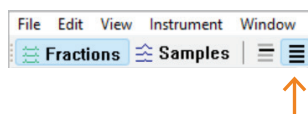


Data for the electropherograms selected in the Experiment pane:

- Displays with electropherograms either overlaid or stacked in the Graph pane depending on the option you've got chosen.
- Shows only results for the selected row(s) in the Peaks and Injections panes.

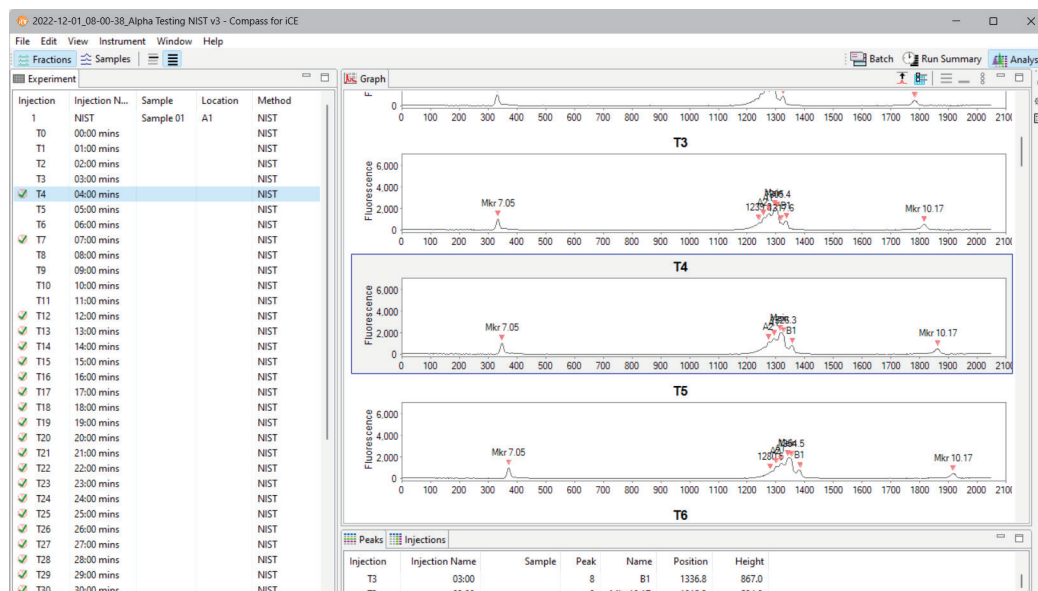


- **To view all electropherograms in a multi-electropherogram format** - Click **View All** in the View bar or select **View** in the main menu and click **All**:



Data for the electropherogram:

- Displays with the electropherograms in the Graph pane in stacked view.
- Graph pane will scroll to the electropherogram when selected in the Experiment pane.
- Shows the results for electropherograms in the Peaks and Injections panes.




Data Notifications and Warnings

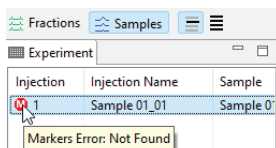
If Compass for iCE detects a potential data issue, a notification or warning icon will display next to the injection row in the Experiment pane.



Manual correction of sample data notification - This means the sample data was manually changed by a user, for example to add or remove a sample peak. Roll your mouse over the icon to display the type of modification that was made.

Experiment		
Injection	Injection Name	Sample
✓ 1	Sample 01_01	Sample 01
T0	00:00 mins	
T1	01:00 mins	

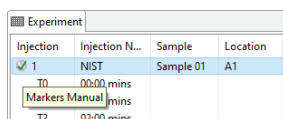
-  **Markers warning** - This means one or more of the pI markers may not be identified properly. You can fix this by manually identifying the pI marker using the steps in “Step 1: Check the pI Markers for the Injected Sample” on page 251. Roll your mouse over the icon to display warning details.



Injection	Injection Name	Sample
1	Sample 01_01	Sample 01


Markers Error: Not Found

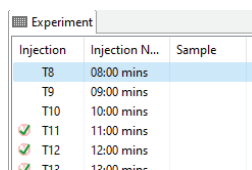
-  **Manual correction of markers data notification** - This means a user changed the pI marker data manually. Roll your mouse over the icon to display the type of modification that was made.




Injection	Injection N...	Sample	Location
1	NIST	Sample 01	A1

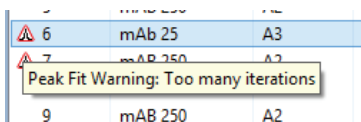
Markers Manual

-  **Manual correction of peak notification** - This means a user added or removed a peak in a mobilization electropherogram.



Injection	Injection N...	Sample	L
T8	08:00 mins		
T9	09:00 mins		
T10	10:00 mins		
T11	11:00 mins		
T12	12:00 mins		

-  **Peak fit warning** - Means that a peak can't be fit properly. This can sometimes be caused when a broad peak is fitted as multiple narrow peaks. Changing the peak width can help in this case. The warning is also caused by very small peaks around main peaks, or small peaks that are close to the end of the separation range. You can often fix this by removing the peak(s) using the steps in “Step 2: Checking Sample Peaks” on page 254. Roll your mouse over the icon to display warning details.



Injection	Injection N...	Sample
6	mAb 25	A3
7	mAb 250	A2
9	mAb 250	A2

Peak Fit Warning: Too many iterations

Checking Your Results

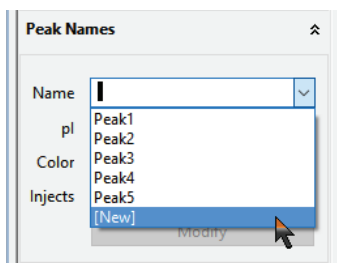
If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues. Compass for iCE detects your sample protein and pI marker peaks and reports results automatically. But, we always recommend you review your data using the steps in this section as a good general practice to make sure your results are accurate. Please see the step by step procedure in “Checking Your Data” on page 250 to do this. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

Naming Peaks

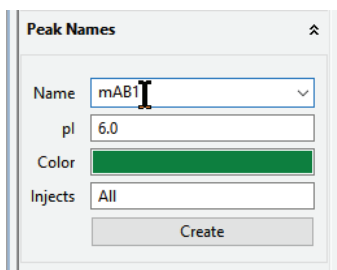
NOTE: Analysis screen options will let you add a new peak name, its pI, identifying color for peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set pI ranges, you can do that in the run analysis settings. For more info go to “Peak Names Settings” on page 622.

Adding New Peak Names in the Analysis Options Pane

1. Click the down arrow in the Name field and select New.



2. Type a name.



- Click in the **pI** field and enter the pI value of your sample protein.

Peak Names

Name: mAB1

pI: 6.78

Color: [Green Bar]

Injects: All

Create

- Click on the **Color** field to display the color selection box.

Peak Names

Name: mAB1

pI: 6.78

Color: [Green Bar]

Injects: All

Create

Color

Basic colors:

Custom colors:

Define Custom Colors >>

OK Cancel

- The color you pick is used to identify the sample protein peak in the Peaks and Injections panes in the Analysis screen. Click a color or define a custom color and click **OK**. The color selection will update in the field:

Peak Names

Name: mAB1

pI: 6.78

Color: [Red Bar]

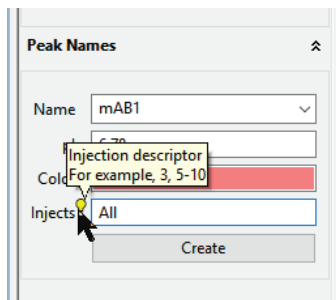
Injects: All

Create

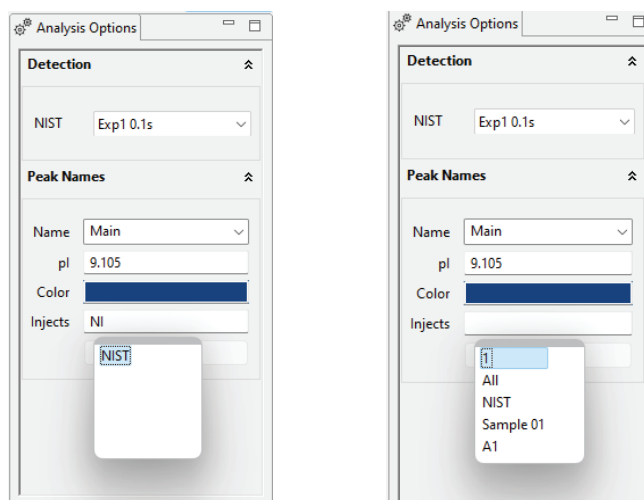
- Click in the **Injects** field:

To apply the new peak name to all injections - leave the default setting of All.

To apply the new peak name to specific injections - enter individual injections separated by commas (for example 2, 5, 8), a range of injections (for example 3-15) or a combination of both (for example 8, 10, 12-15). You can also click the tool tip for more examples.



To apply the new peak name to specific run data - type the first letter of a method, sample name, well or vial location. Compass for iCE will search the run data and display a list of matching options for you to choose from. Alternatively you can highlight the text in the Injections cell, hit **Delete**, then select an option from the drop down list.

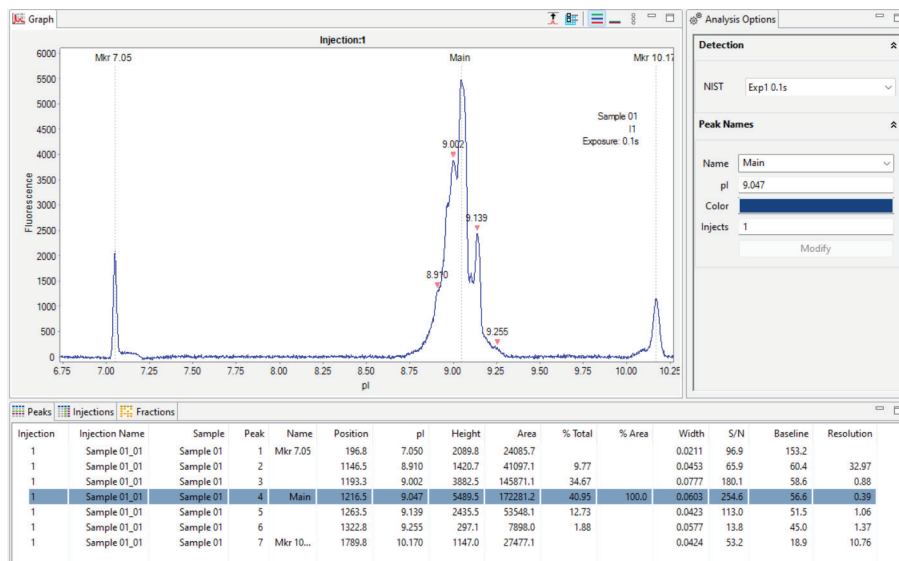


- **Injection** - Applies the peak name to the injection number
- **All** - Applies the peak name to all injections.
- **Methods** - All methods in the run file display in the list. Selecting a method applies the peak name to all injections that used that method.
- **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the peak name to all injections that used that sample name.
- **Wells or vials** - The well or vial numbers currently selected will display in the list. Selecting a well or vial number applies the peak name to all injections that used that well or vial.

7. Click **Create** to save changes.

NOTE: To clear the new peak name information you've entered without saving changes, just delete the peak name in the Name field.

The named peak will be identified with a peak name label in the electropherogram and color-coded in the Peaks and Injections panes:

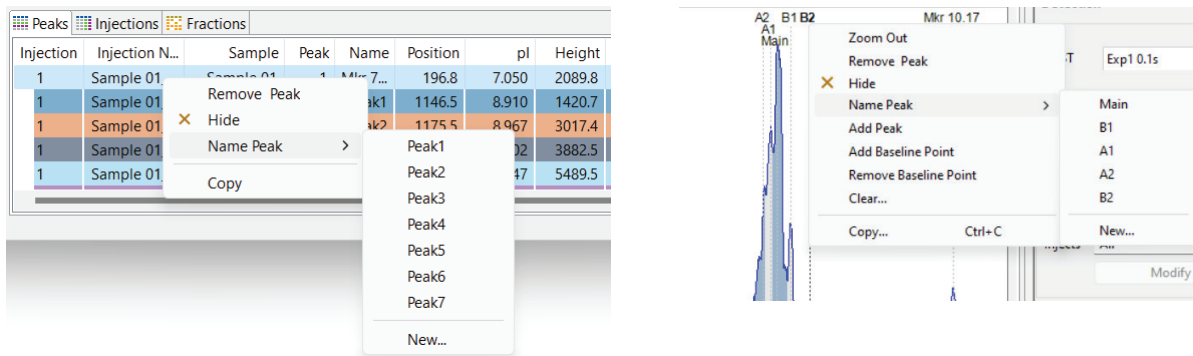


Adding Peak Names from the Graph or Peaks Table

NOTE: Analysis screen options will let you change an existing peak name, its pI, identifying color in peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set pI ranges, you can do that in the run analysis settings. For more info go to "Manual Peak Integration" on page 637.

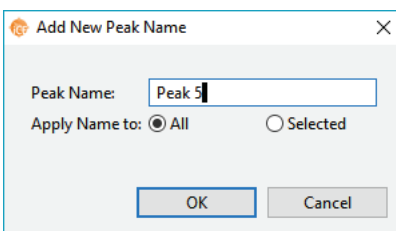
1. Right click the peak you want to name in the Graph or Peaks pane.

2. Select Name Peak.



3. To use an existing peak name - select a name from the list.

To create a new peak name - select **New**. Type in a name for the peak. Click **All** to apply to all injections or **Selected** to apply only to the injections selected.

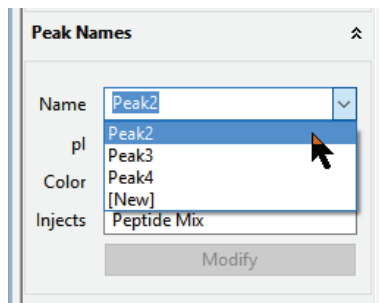


4. Click OK. If you want to add an identifying color or change the injections the new peak name applies to, you can do that in the Analysis Options pane.

Changing Peak Name Information

NOTE: Analysis screen options will let you change an existing peak name, its pI, identifying color for peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set pI ranges, you can do that in the run analysis settings. For more info go to “Manual Peak Integration” on page 637.

1. In the Analysis Options pane, click the down arrow in the Name field and select an existing peak name.



2. Change the name, pI, color and injects as needed then click **Modify**.

Copying Results Tables and Graphs

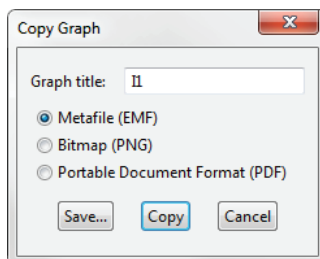
You can copy and paste data and results tables into other documents, or save the electropherogram as a graphic file.

Copying Results Tables

1. Click in the Peaks or Injections pane.
2. Select one or multiple rows.
3. Select **Edit** in the main menu and click **Copy**, or right click on row(s) you selected and click **Copy**.
4. Open a document (Microsoft® Word®, Excel®, PowerPoint®, etc.). Right click in the document and select **Paste**. Data for the rows selected will be pasted into the document.

Copying the Graph

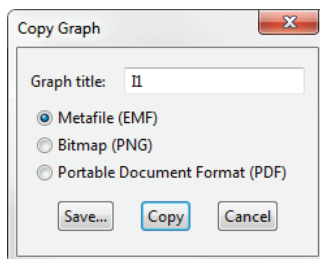
1. Select the Graph pane.
2. Select **Edit** in the main menu and click **Copy**, or right click in the Graph pane and select **Copy**.
3. Select an image option (EMF, PNG or PDF) in the pop-up window, then click **Copy**.



4. Open a document (Microsoft® Word®, Excel®, PowerPoint®, etc.). Right click in the document and select **Paste**. A graphic of the copied electropherogram will be pasted into the document.

Saving the Graph as an Image File

1. Select the Graph pane.
2. Select **Edit** in the main menu and click **Copy**, or right click in the Graph pane and select **Copy**.
3. Select an image option (EMF, PNG or PDF) in the pop-up window, then click **Save**.



4. Select a directory to save the file to, enter a file name, then click **OK**.

Exporting Run Files

Results tables and raw plot data can be exported for use in other applications. Data is organized into these folders when exported:

- <run file name>Export_FL when the same exposure is selected for all methods

Exporting Results Tables

To export the information in the Peaks and Injections tables:

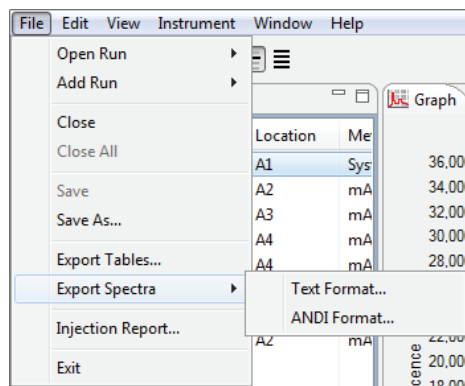
1. Click **File** in the main menu and click **Export Tables**.
2. Select a directory to save the files to and click **OK**. Data will be exported in .txt format.

NOTE: To exclude export of standards (pI markers) data or export results table data in .csv format, see “Setting Data Export Options” on page 760.

Exporting Raw Sample Electropherogram Data

To export raw sample plot and background data:

1. Click **File** in the main menu and click **Export Spectra**.



- **To export data in .txt format** - Select **Text Format**. Data will be exported in one file for the sample injection and all fractions.
- **To export data in .cdf format** - Select **ANDI Format**. Data will be exported in one file per injection/fraction.

2. Select a directory to save the files to and click **OK**. Data will be exported in the selected format.

Changing Sample Protein Identification

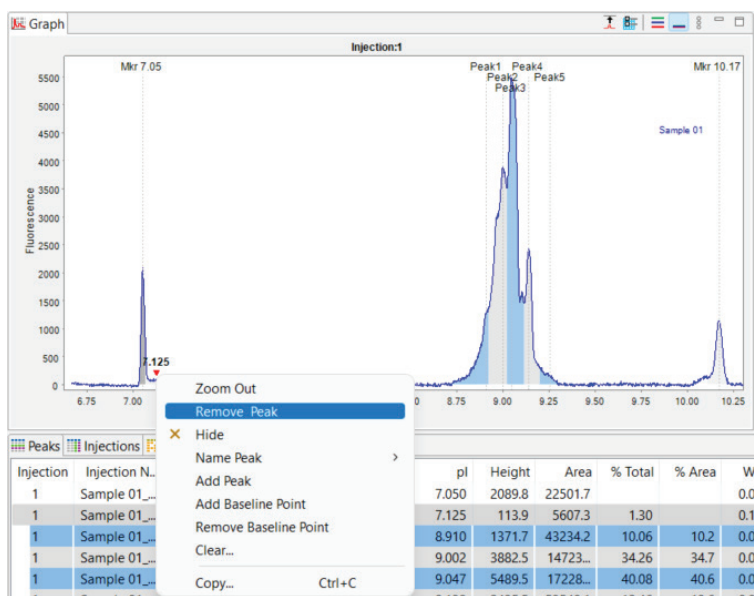
Compass for iCE lets you customize what sample proteins are reported in the results tables by making manual adjustments in the electropherogram or Peaks table.

Adding or Removing Sample Data

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.

3. Click on the row in the experiment pane that has the injection you want to correct, then click the **Graph** tab.

- **To remove a peak from the data** - Right click the peak in the electropherogram or Peaks table and select **Remove peak**. The software will no longer identify it as a sample peak in the electropherogram, and the peak data will be removed in the results tables.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

Experiment		
Injection	Injection Name	Samp
✓ 1	Sample 01_01	Samp

- **To add an unidentified peak to the data** - Right click the peak in the electropherogram or peaks table and select **Add Peak**. The software will calculate and display the results for the peak in the results tables and identify the peak in the electropherogram.

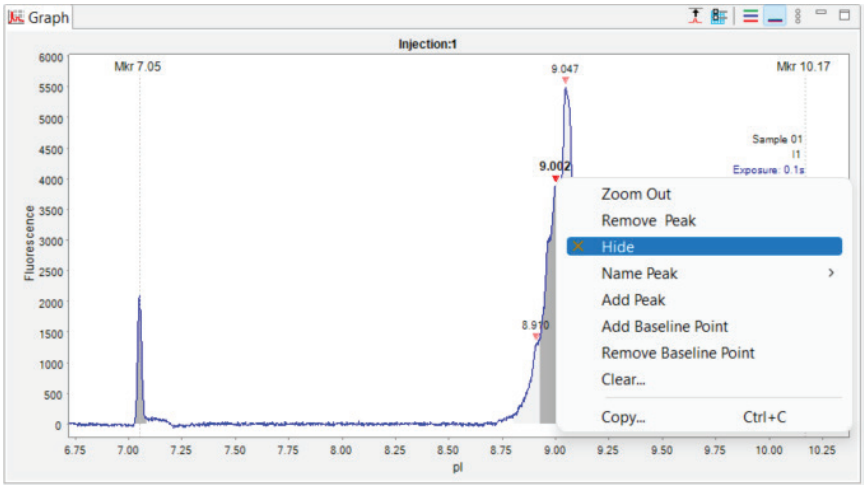
A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear** for the current injection or **Clear All** for all injections in the batch.

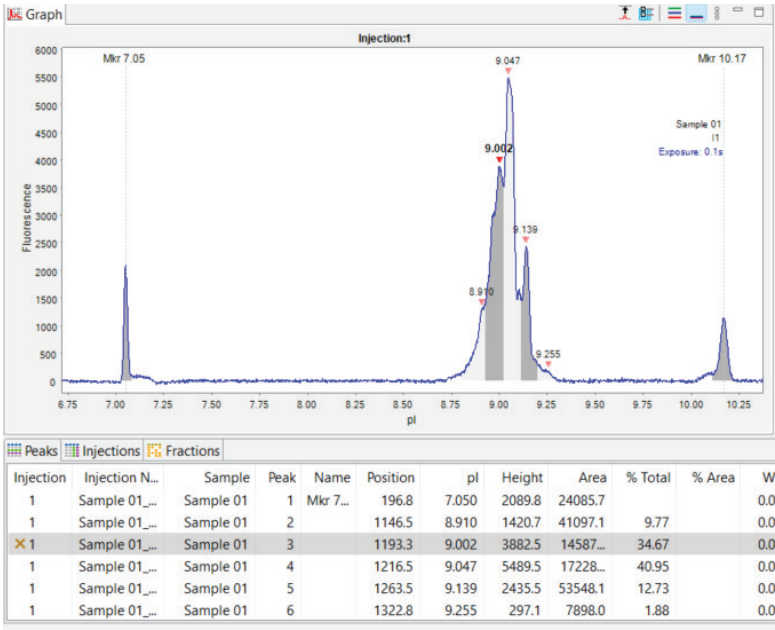
Hiding Sample Data

You can hide the results for a sample protein in the results tables without completely removing it from the reported results.

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that contains the injection you want to correct, then click the **Graph** tab.
4. Right click the peak in the electropherogram or Peaks table and select **Hide**. Compass for iCE will hide the peak data in the results tables.



5. To view hidden peak data, click **View** in the main menu and click **Show Hidden**. Hidden peak data will display in the results table and be marked with an X.

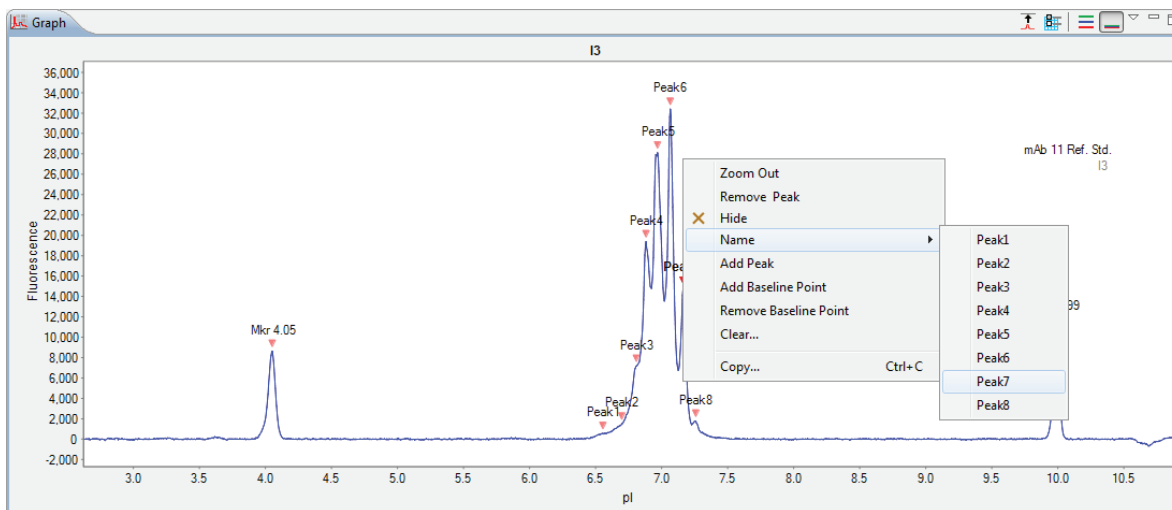


6. To unhide a peak, right click on the peak in the electropherogram or peaks table and select **Unhide**.

Changing Peak Names for Sample Data

If Compass for iCE did not automatically name a sample protein peak, you can do it manually.

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that has the sample you want to correct, then click the **Graph** pane.
4. Right click the peak in the electropherogram or Peaks table and click **Name**, then select a name from the list. Compass for iCE will change the peak name in the electropherogram and results tables, and adjust peak names for other sample proteins accordingly.



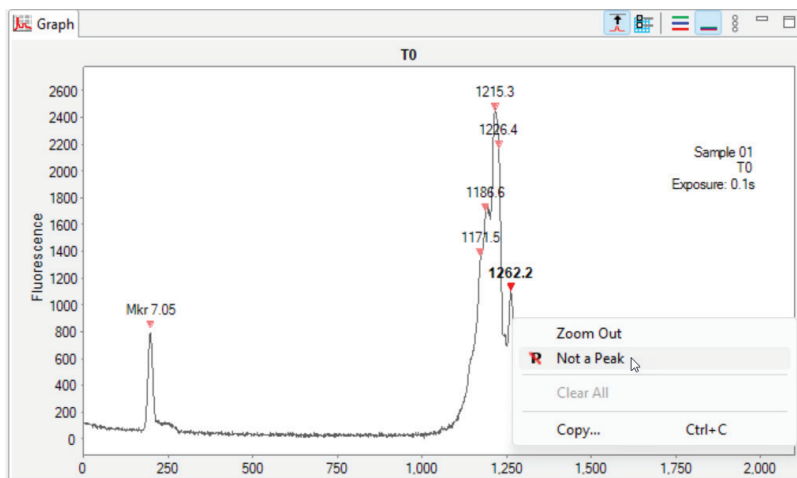
NOTE: For details on how to specify peak name settings, see “Manual Peak Integration” on page 637.

Adding or Removing Fractions Data

1. Click **Fractions** in the View bar.
2. Click **Single View** in the View bar.

3. Click the Graph tab.

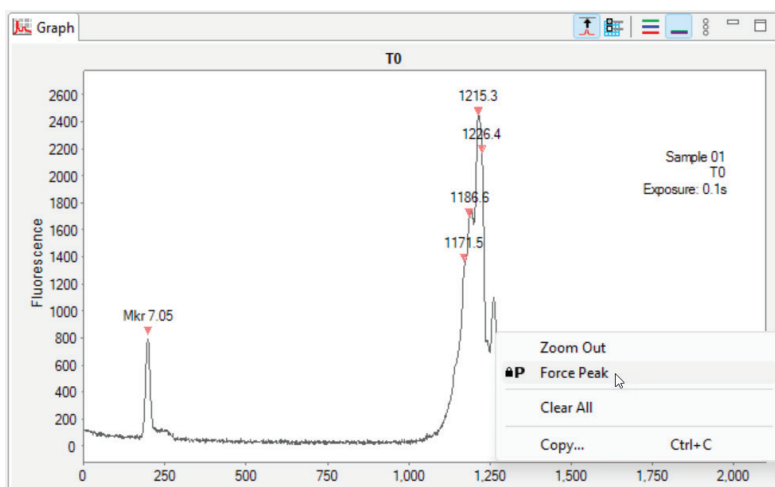
- **To remove a peak** - Right click the peak in the electropherogram or Peaks table and select **Not a Peak**. The software will not longer identify it as a sample peak in the electropherogram, and the peak data will be removed from the results table.



A checkmark will appear next to electropherogram in the Experiment pane to indicate a manual correction was made.

Experiment	
Injection	Injection Name
1	Sample 01_01
✓ T0	00:00 mins
T1	01:00 mins
T2	02:00 mins

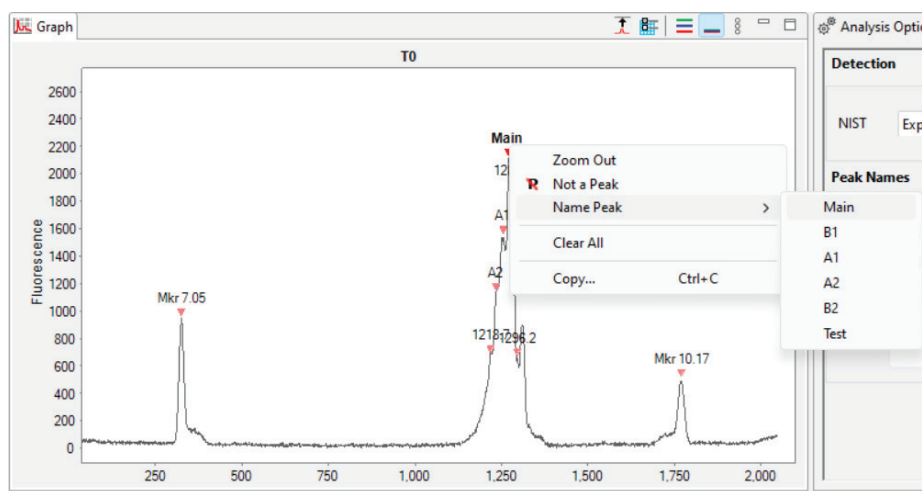
- **To add an unidentified peak**: Right click the peak in the electropherogram and select **Force Peak**. The software will calculate and display the results for the peak in the results table and identify the peak in the electropherogram.



A checkmark will appear next to the electropherogram in the Experiment pane to indicate a manual correct was made.

NOTE: To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear** for the current injection or **Clear All** for all injections in the batch.

- **To name a peak:** Right click the peak in the electropherogram or peaks table and select **Name Peak**. Select the name to assign to the peak. The software will update the peak name in the results table.







IMPORTANT: The sample peaks in the final focusing sample injection image must first be named in the Samples view before peaks can be named in the Fractions view.

Changing the Electropherogram View

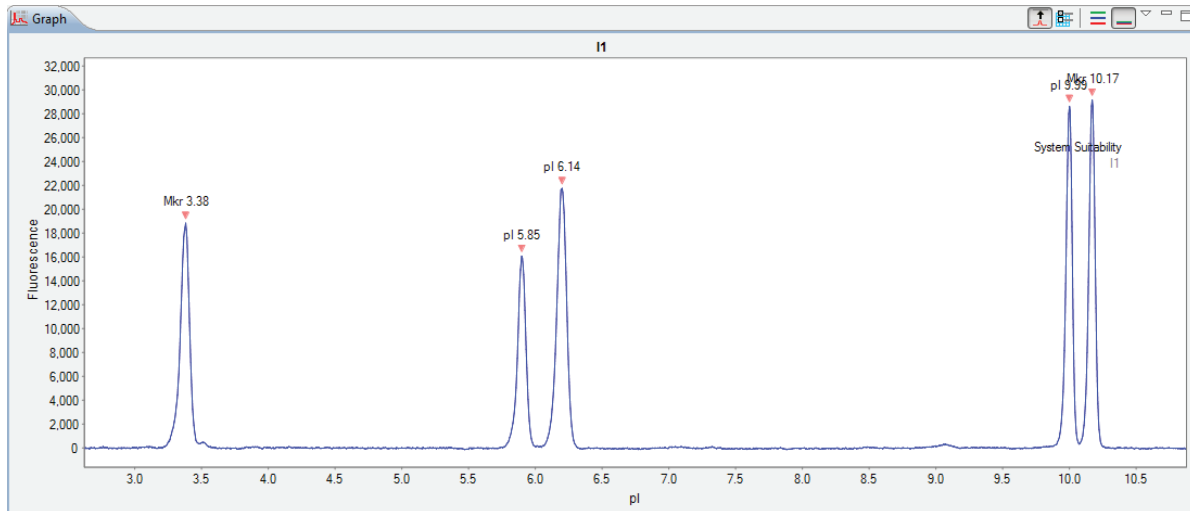
Options in the Graph pane let you zoom and rescale electropherograms, overlay or stack plots and change the peak and plot info displayed.

The Graph pane toolbar has these options:

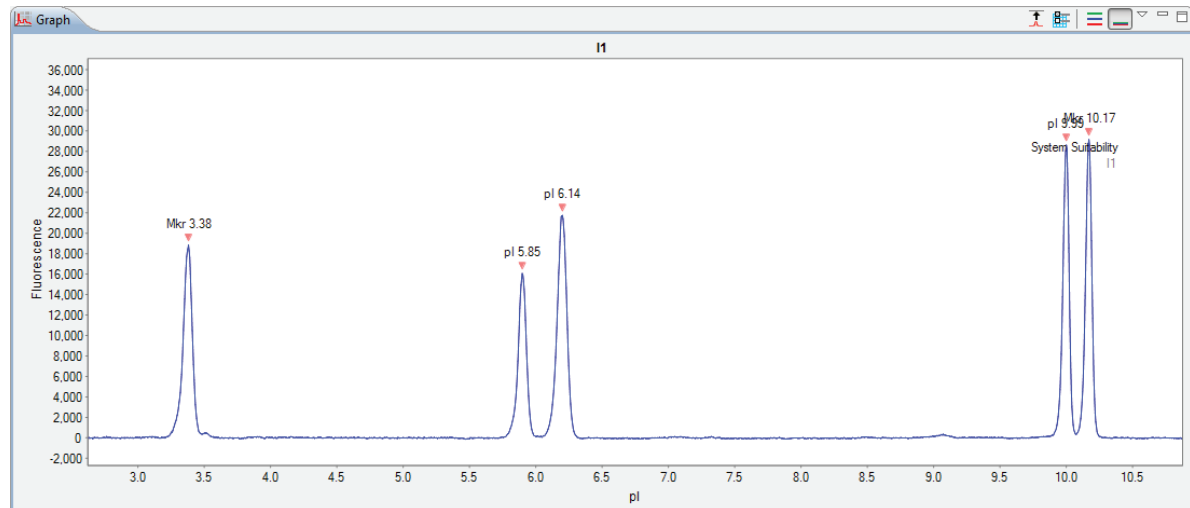
-  Auto Scale
-  Graph Options
-  Stack the Plots
-  Overlay the Plots

Autoscaling the Electropherogram

Click the **Auto Scale** button to scale the y-axis to the largest peak in the electropherogram.

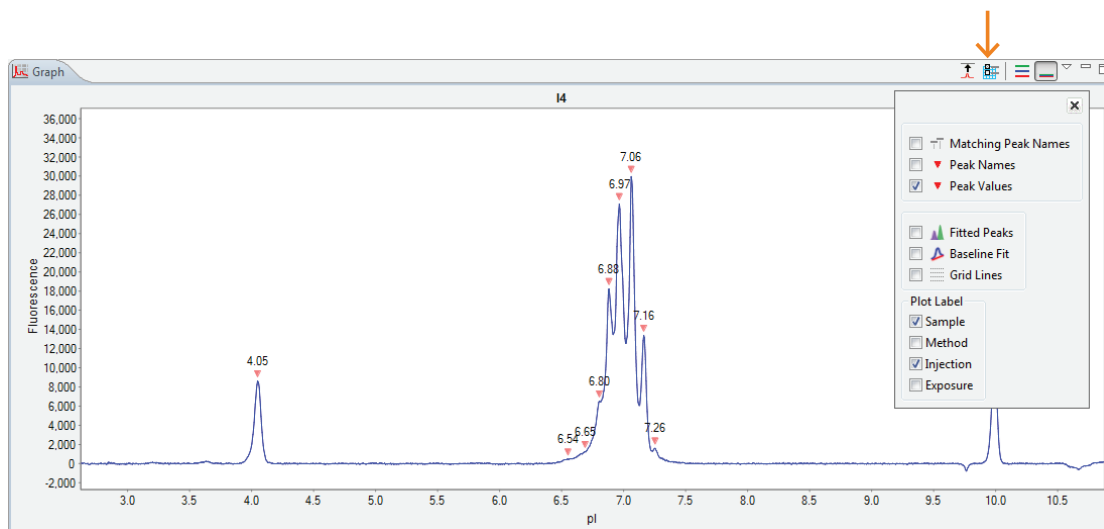


Click the **Auto Scale** button again to return to default scaling.



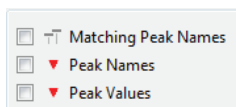
Customizing the Data Display

You can customize electropherogram peak labels, plot labels and display options. To do this, just select the **Graph Options** button.

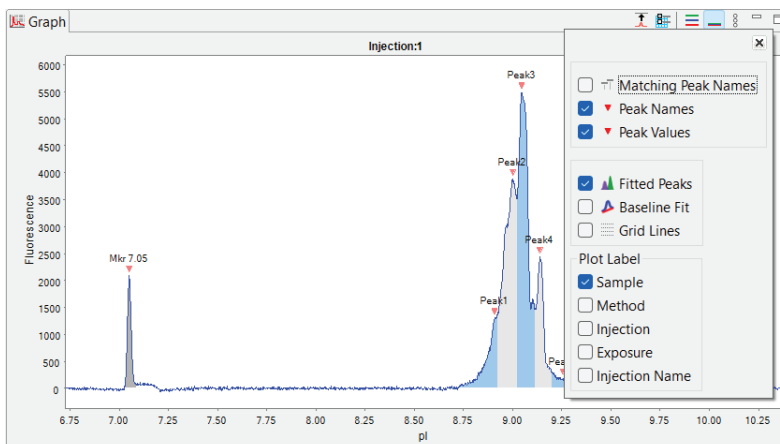


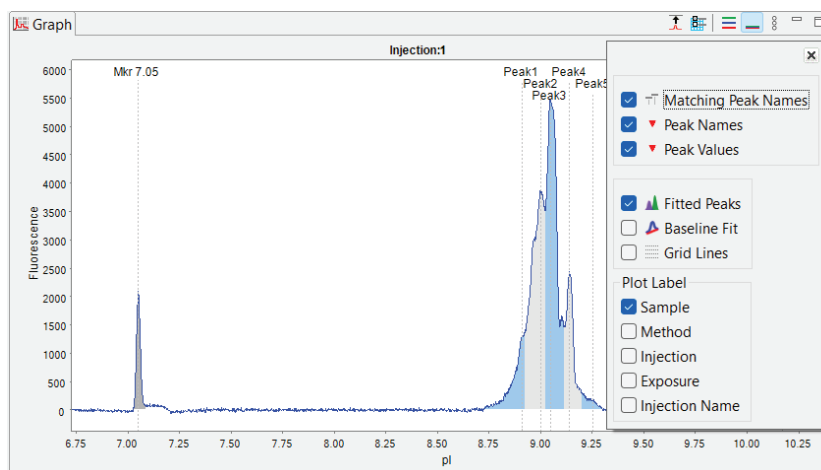
Peak Labels

You can customize the labels used to identify peaks in the electropherogram with these options:



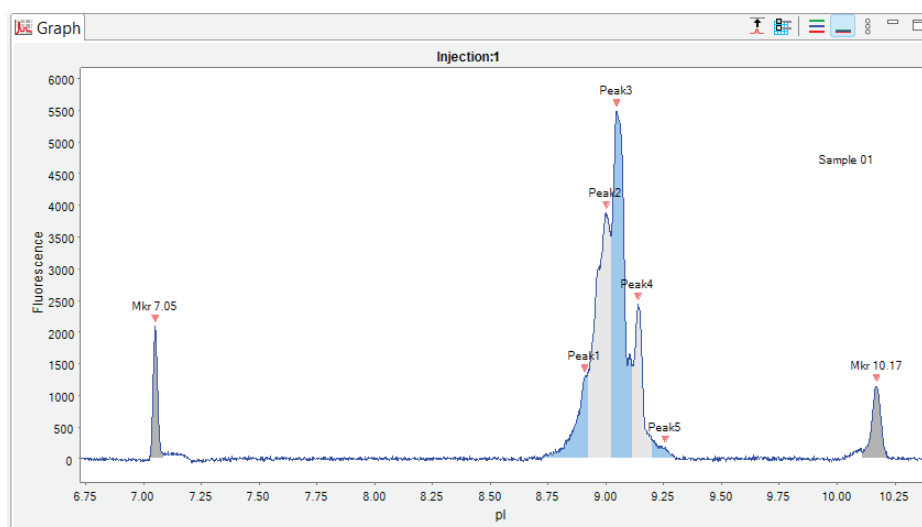
- **Matching Peak Names** - Checking this box will draw vertical lines through each named peak. Using this option with Stack the Plots or Overlay the Plots features is helpful for visually comparing your named peaks across multiple injections.





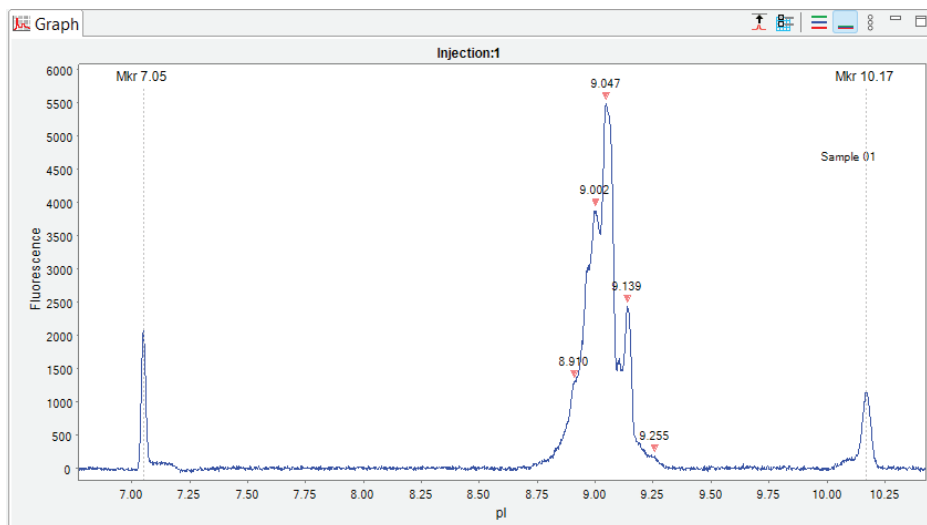
- **Peak Names** - Checking this box displays peak name labels on all named peaks in the electropherogram.

NOTE: If more than one peak label option is selected, peak name labels will always be used for named peaks.



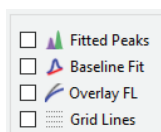
- **Peak Values** - Checking this box will display the pI values on all peaks in the electropherogram.

NOTE: If more than one peak label option is selected, peak name labels will always be used for named peaks.



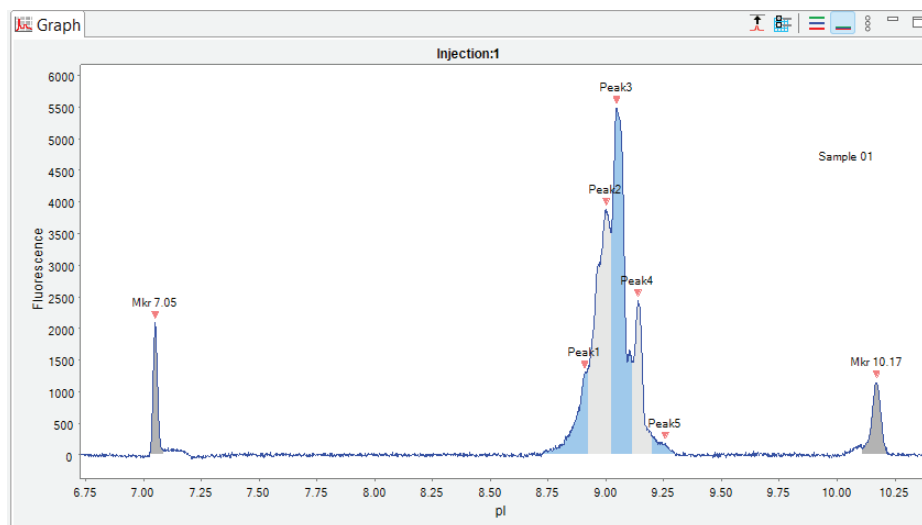
Baseline and Grid Options

You can view the calculated baseline fit, peak integration, show grid lines and overlay fluorescence electropherograms with these options.



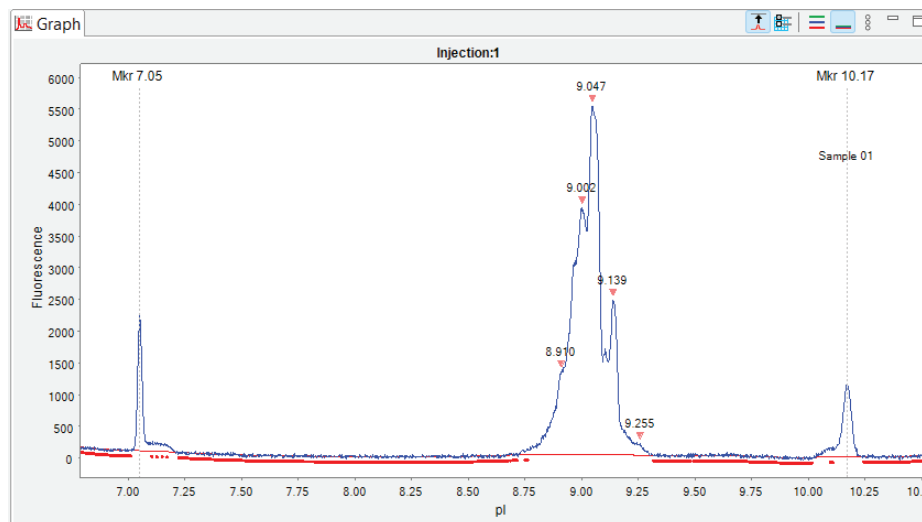
- **Fitted peaks** - Checking this box displays how the peaks were fit by the software. For MauriceFlex Fractionation runs, the software uses Dropped Line as the default.

NOTE: This option is only available for sample data.

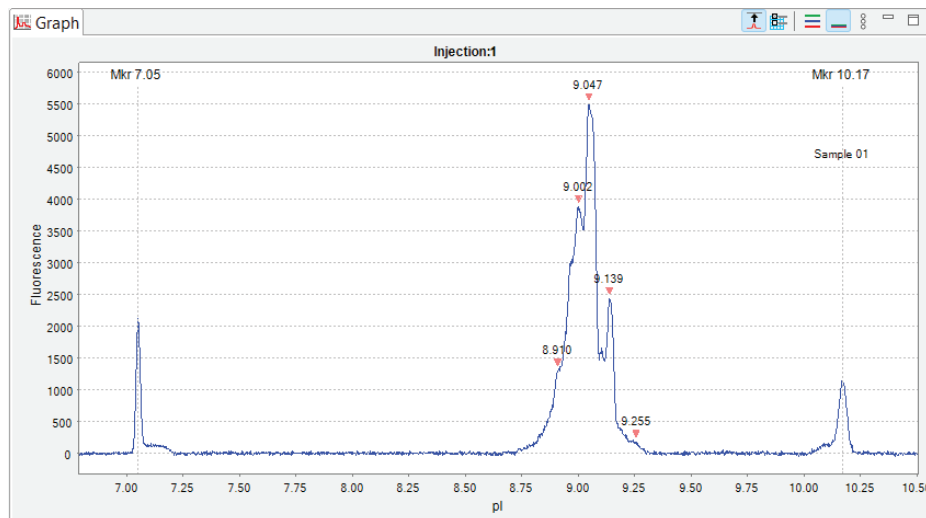


- **Baseline Fit** - Checking this box displays the calculated baseline for the peaks. Baseline points will also display for regions of the electropherogram considered to be at baseline.

NOTE: This option is only available for sample data.



- **Grid Lines** - Checking this box adds grid lines in the graph.



Plot Labels

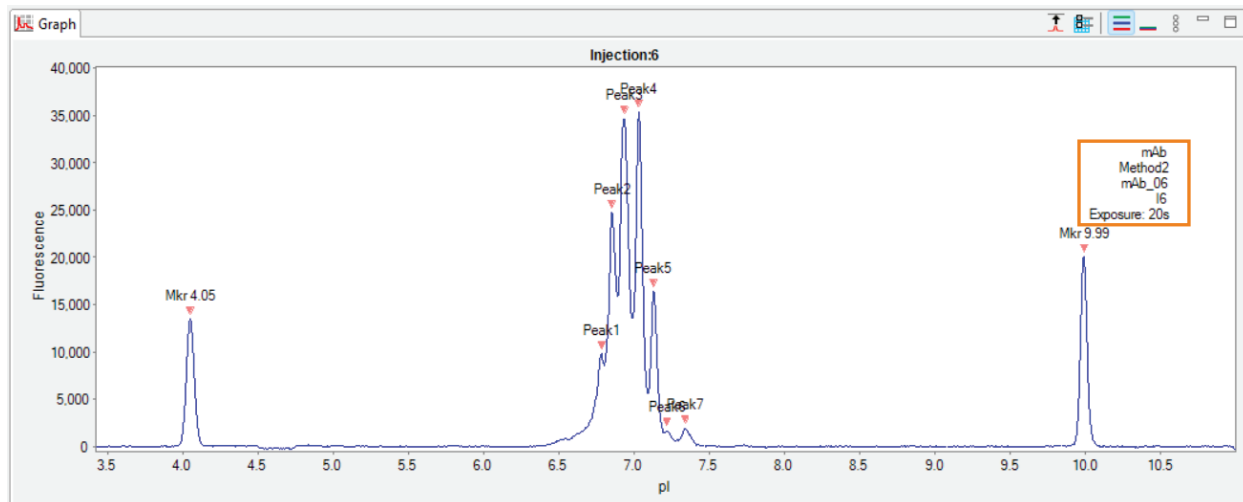
You can customize the plot labels displayed on the electropherogram with these options.

Plot Label	
<input checked="" type="checkbox"/>	Sample
<input type="checkbox"/>	Method
<input checked="" type="checkbox"/>	Injection
<input checked="" type="checkbox"/>	Exposure
<input type="checkbox"/>	Injection Name

Plot labels are shown in the upper right side of the graph.

- **Sample** - Checking this box displays the sample name used for the injection. If sample names were entered with the batch, those names will display here. If not, Sample (default name) displays.
- **Method** - Checking this box displays the method used for the injection.
- **Injection** - Checking this box displays the injection number. For example, I1 for injection 1 in the run.
- **Exposure** - Checking this box display the exposure time(s) used for the data.
- **Injection Name** - Checking this box displays the name used for the injection.
- **Injection Name** - Checking this box displays the injection name used for the injection.

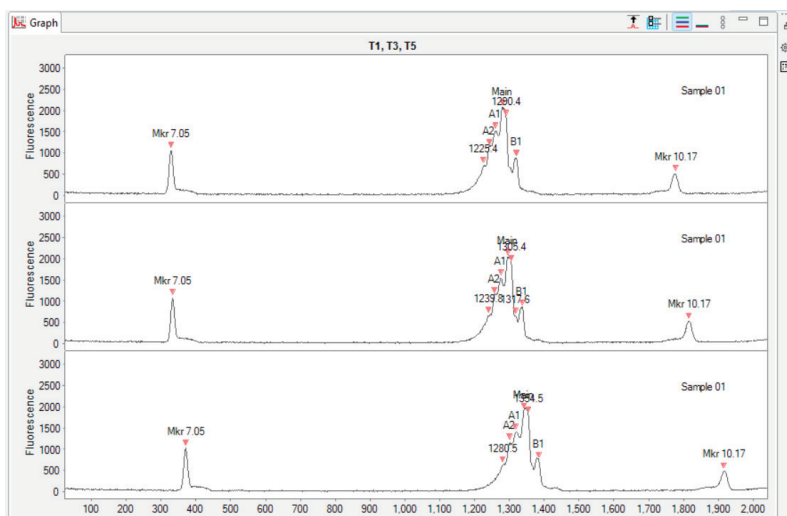
Here's an example of an electropherogram with all plot labels selected:



Stacking Multiple Electropherograms

You can stack electropherograms in the Fractions View for multiple injections or mobilization electropherograms vertically in the Graph pane for comparison. Click **Single View**.

4. Select multiple rows in the Experiment pane.
5. Click the **Stack the Plots** button. The individual electropherograms for each row you selected will stack in the Graph pane.

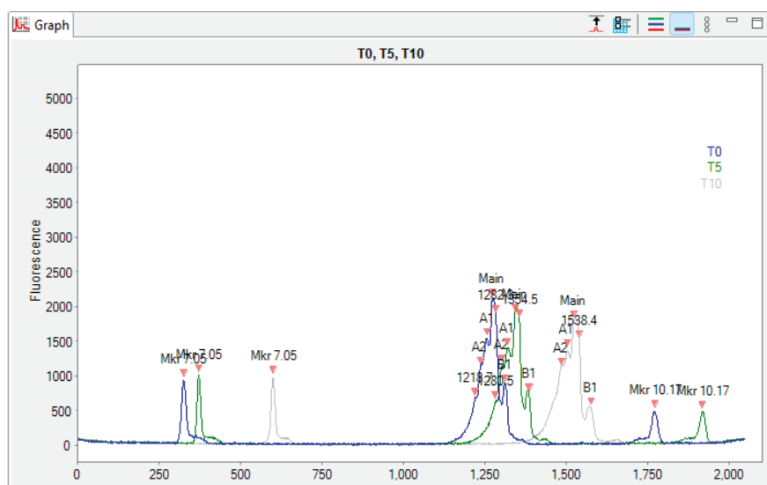


You can also customize the colors used for the stacked plot display. To do that go to "Setting up Automatic Injection Reports" on page 761.

Overlaying Multiple Electropherograms

You can overlay electropherograms in the Fractions View for multiple $\nu\phi\epsilon\chi\tau\iota\omicron\nu\sigma$ $\sigma\rho$ $\mu\omicron\beta\iota\lambda\iota\zeta\alpha\tau\iota\omicron\nu$ electropherograms on top of each other for comparison in the Graph pane.

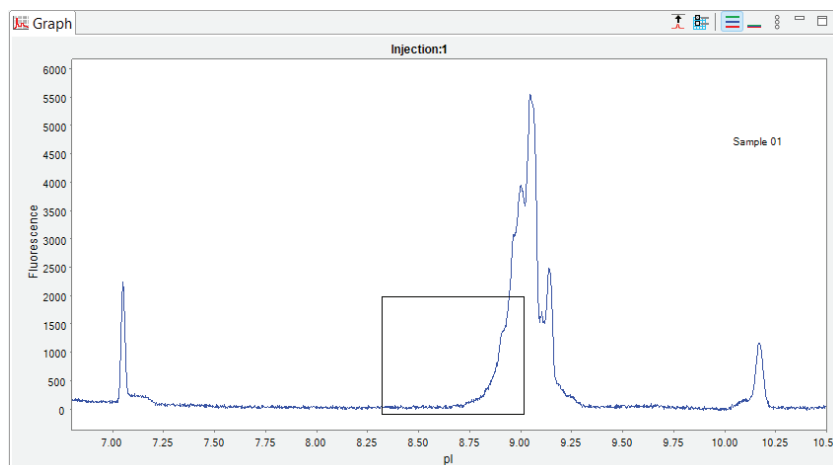
1. Click **Single View**.
2. Select multiple rows in the Experiment pane.
3. Click the **Overlay the Plots** button. The individual electropherograms for each row you selected will overlay in the Graph pane.



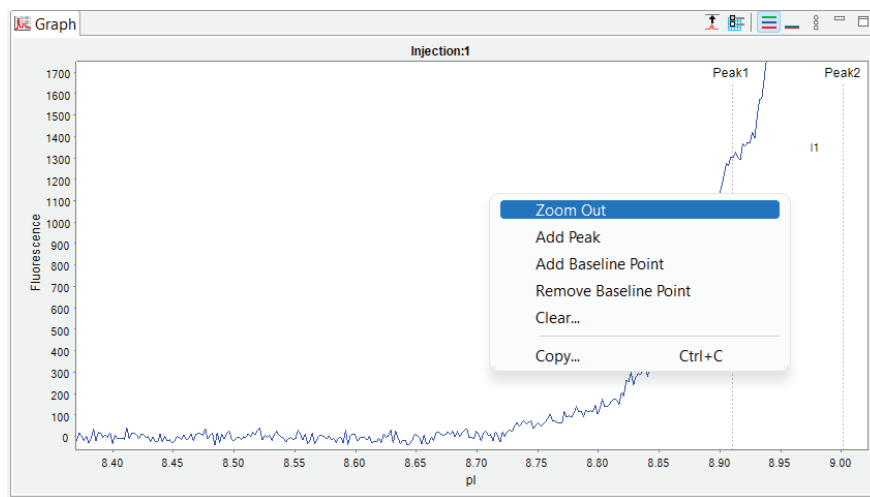
You can also customize the colors used for the overlay plot display. To do that go to “Setting up Automatic Injection Reports” on page 761.

Zooming

To zoom in on a specific area of the electropherogram, hold the mouse button down and draw a box around the area with your mouse:

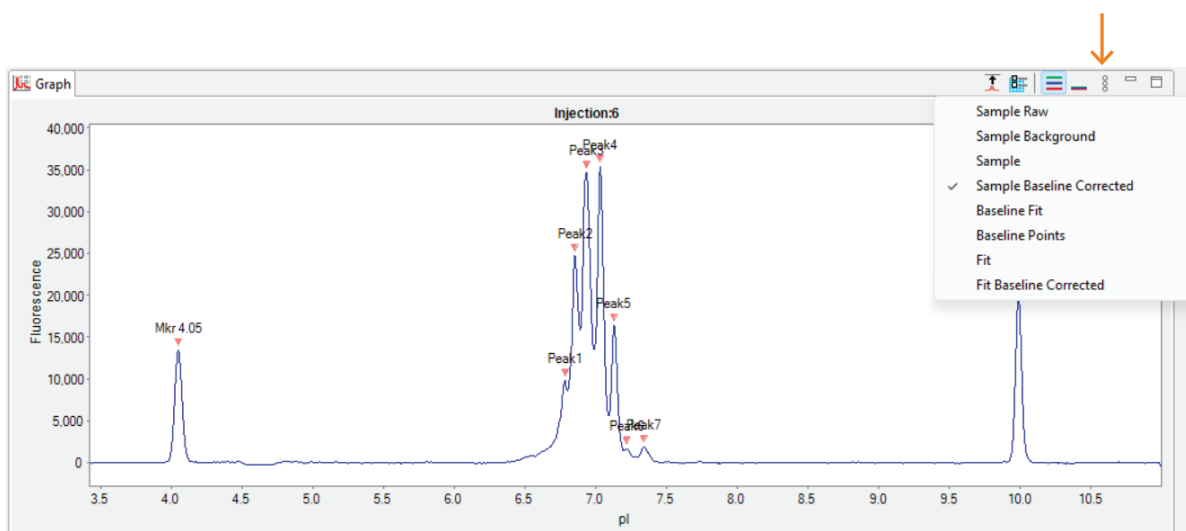


To return to default scaling, right click in the electropherogram and click **Zoom Out**.



Selecting Data Viewing Options

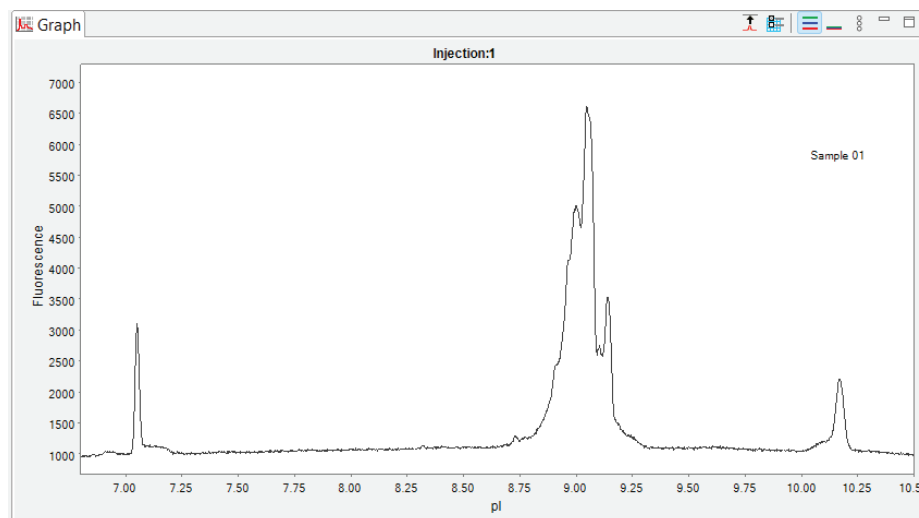
The graph view menu gives you multiple options for changing what type of electropherogram data is displayed. Just click **View More** in the graph pane toolbar to view the menu:



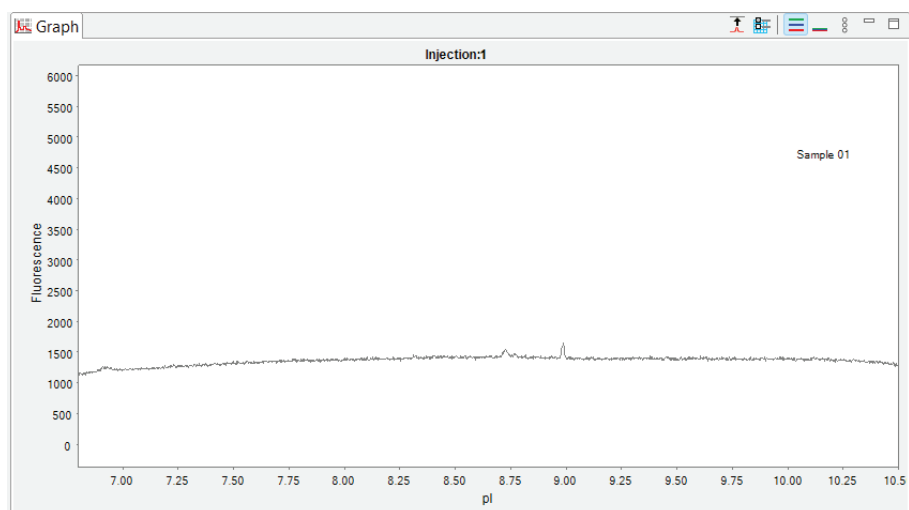
A check mark next to the menu option indicates it's currently selected, and you can select multiple options at once.

NOTE: Unless noted otherwise, graph view menu options are available for sample data only.

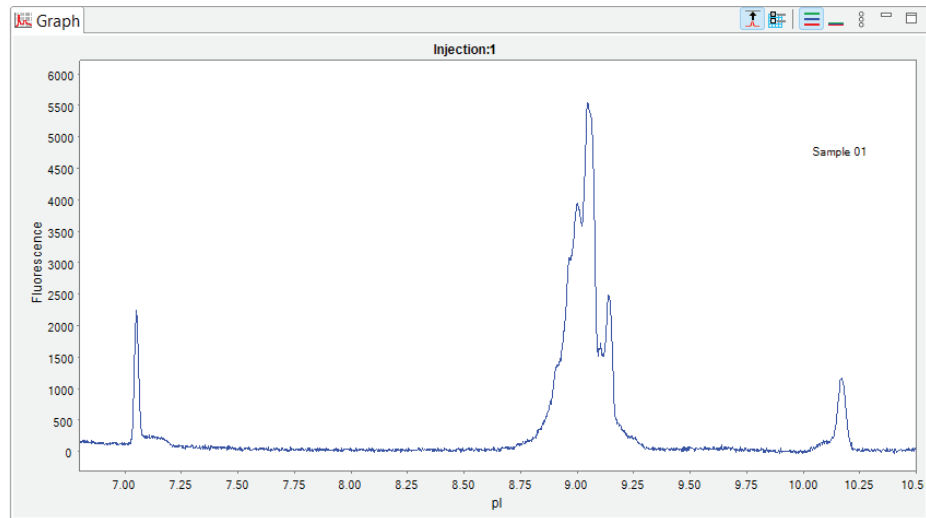
- **Sample Raw** - Clicking this option displays the basic detector values used to calculate peak fluorescence.



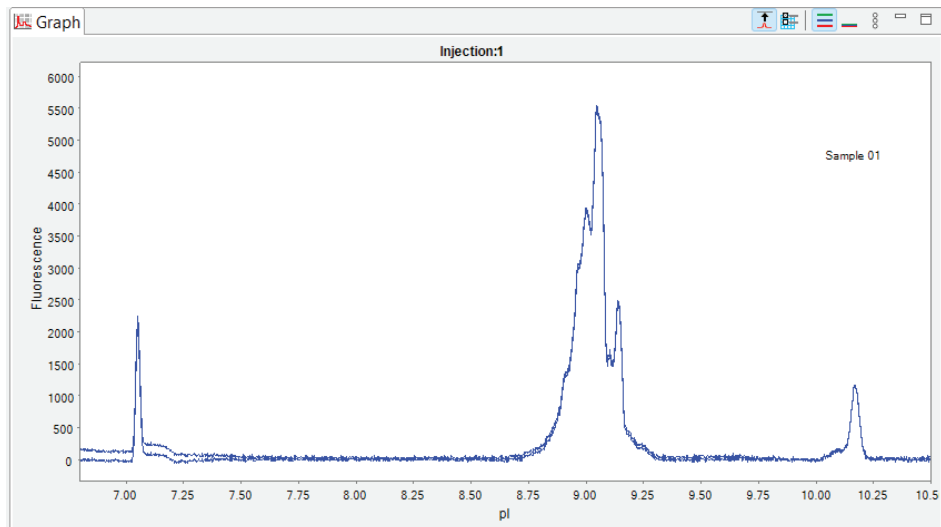
- **Sample Background** - Clicking this option displays the basic detector values used to calculate baseline fluorescence.



- **Sample** - Clicking this option displays raw, uncorrected sample data.

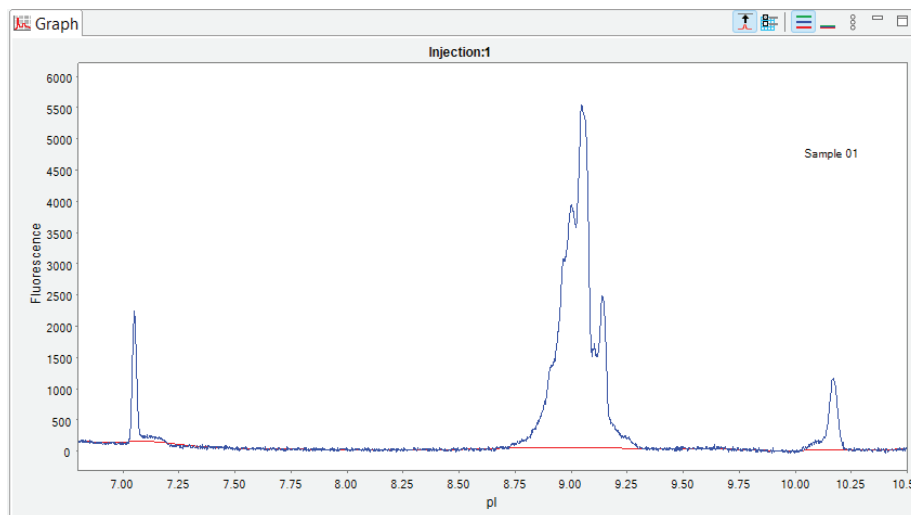


- **Sample Baseline Corrected** - Clicking this option displays sample data with the baseline subtracted (zeroed). This is the default view. In this next example, both Sample and Sample Baseline Corrected are selected.



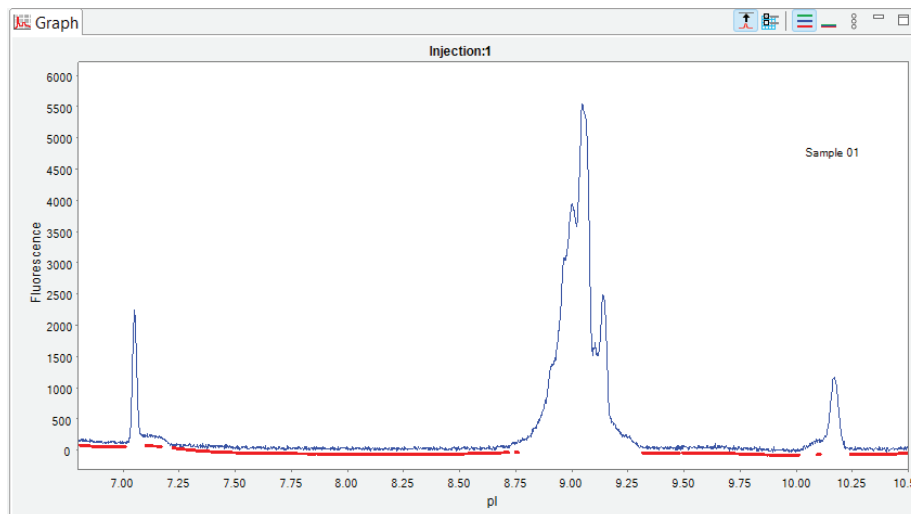
- **Baseline Fit** - Clicking this option displays the calculated baseline for the raw sample data. In this next example, both Baseline Fit and Sample are selected.

NOTE: This option is selected automatically when Baseline Fit is selected in graph options.

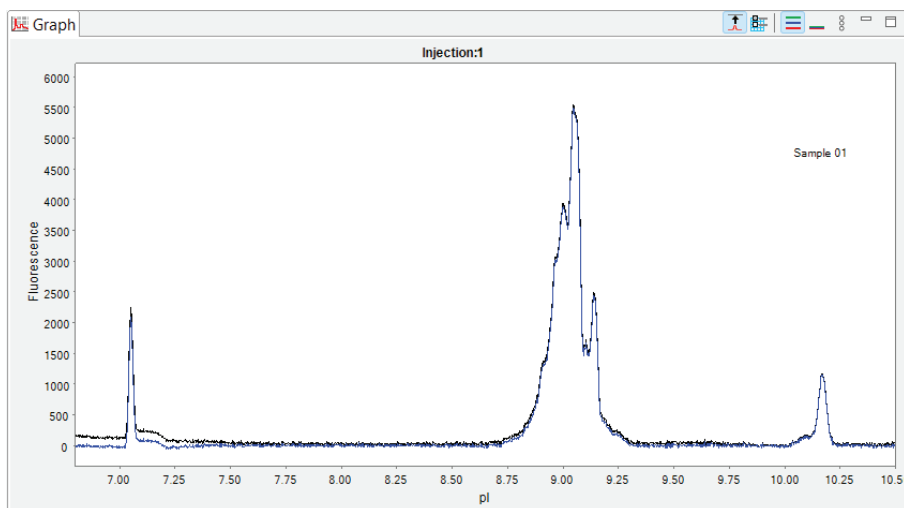


- **Baseline Points** - Clicking this option displays regions of the electropherogram considered to be at baseline. In this example, both Baseline Points and Sample are selected.

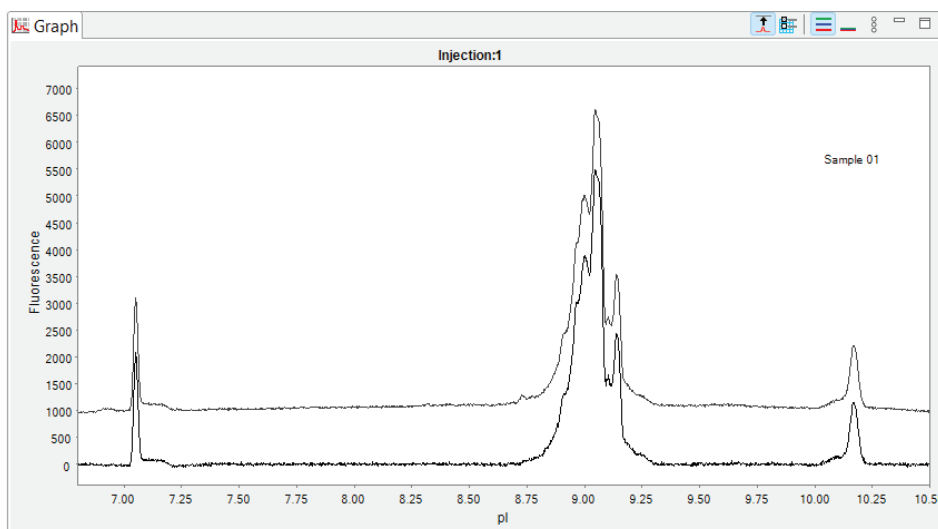
NOTE: This option is selected automatically when Baseline Fit is selected in graph options.



- **Fit** - Clicking this option displays the bounding envelope of the fitted peaks as calculated by the software for the raw sample data. In this example, both Fit and Sample Baseline Corrected are selected.



- **Fit Baseline Corrected** - Clicking this option displays the fitted peaks as calculated by the software for the sample baseline corrected data. In this example, both Fit Baseline Corrected and Sample Raw are selected, the fit plot is on the bottom.

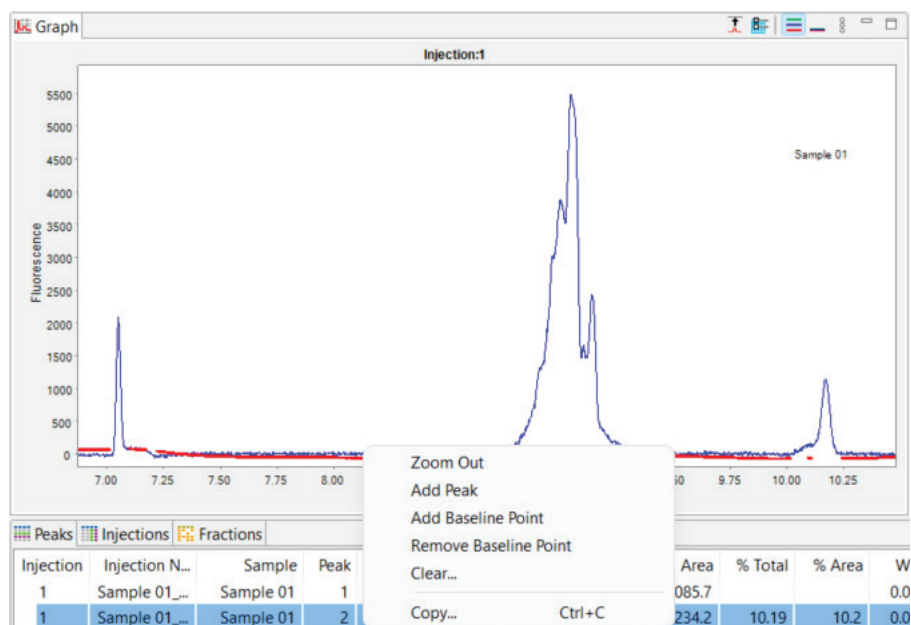


Adjusting the Baseline

Spline Baseline

Points in the baseline can be added or removed as needed.

1. Click the **Graph Options** button in the graph pane toolbar and check **Baseline Points**. This will display baseline points for the raw sample data.
2. Use the mouse to draw a box around the area you want to correct. This will zoom in on the area.
3. Right click a baseline point and select **Add Baseline Point** or **Remove Baseline Point**.



NOTE: To clear the manual addition or removal of baseline points and go back to the original view of the data, right click in the electropherogram and click **Clear All**.

Interpolated Baseline

Compass for iCE lets you interpolate baselines from the base of each peak for named peaks that use Dropped Lines integration.

1. If you haven't already, name your peaks of interest.
2. Select **Edit > Analysis**, and click **Peak Fit** in the left sidebar.

3. Choose **Interpolate** as the Baseline Type if it isn't already selected.

NOTE: Baseline Type selection is only available when Dropped Lines is used for the Area Calculation.

Analysis Groups: Peak Fit

Range

Minimum 3.5

Maximum 10.5

View ☒ Analysis ☐ Full

Baseline

Baseline Type **Interpolate** ▼

Threshold Spline

Window 25.0

Stiffness 1.0

Peak Find

Threshold 10.0

Width 15.0

Area Calculation Dropped Lines ▼

4. Click **OK**.
5. In the Analysis window Graph Pane, click **Graph Options** and select **Fitted Peaks** and **Baseline Fit**.

Graph Options

☒ Matching Peak Names

☒ Peak Names

☒ Peak Values

☒ Fitted Peaks

☒ Baseline Fit

☐ Grid Lines

Plot Label

☒ Sample

☐ Method

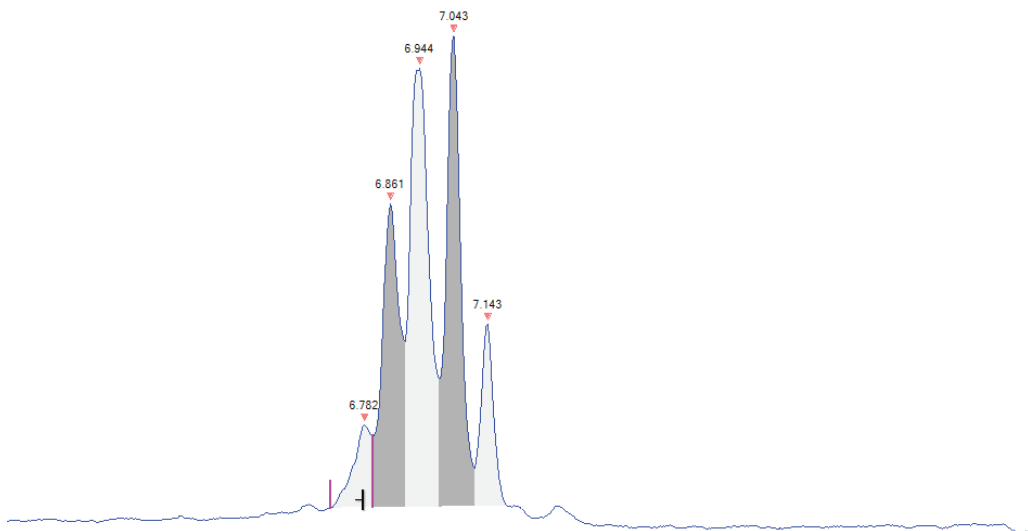
☒ Injection

☒ Exposure

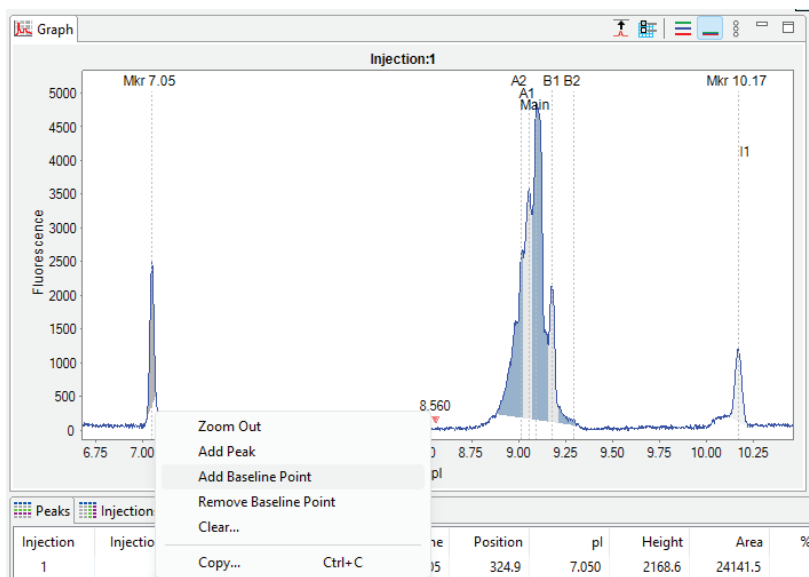
☐ Injection Name

6. Select an injection in the Experiment pane.
7. Hover the mouse over the peak you want to adjust the baseline on. Two magenta bars will display that show the start and end points of the peak's integration.

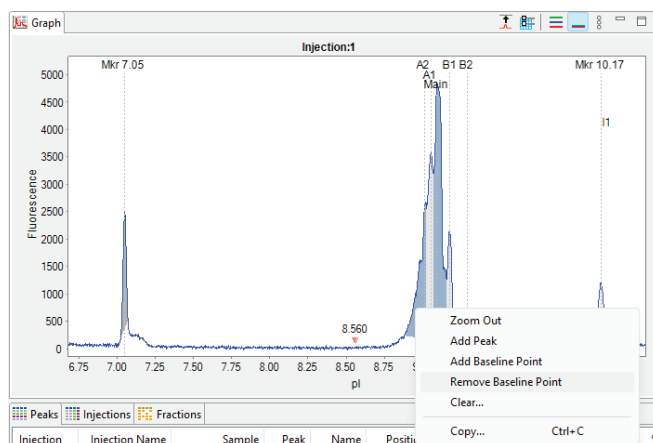
8. To change the baseline, hover the mouse over one of the magenta bars. Then click the mouse and drag the bar to move it. The cursor changes to indicate which direction you can move the integration point. See “Manual Peak Integration” on page 637 for more info on integrating peaks.



9. Baseline points can also be added between two adjacent peaks for a better fit. Hover the mouse over the peak until the magenta bars appear, then right click and select **Add Baseline Point**.

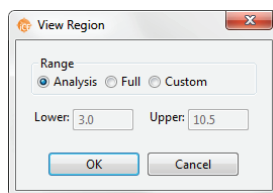


10. To remove the baseline point, hover the mouse over the peak until the magenta bars appear, then right click and select **Remove Baseline Point**.

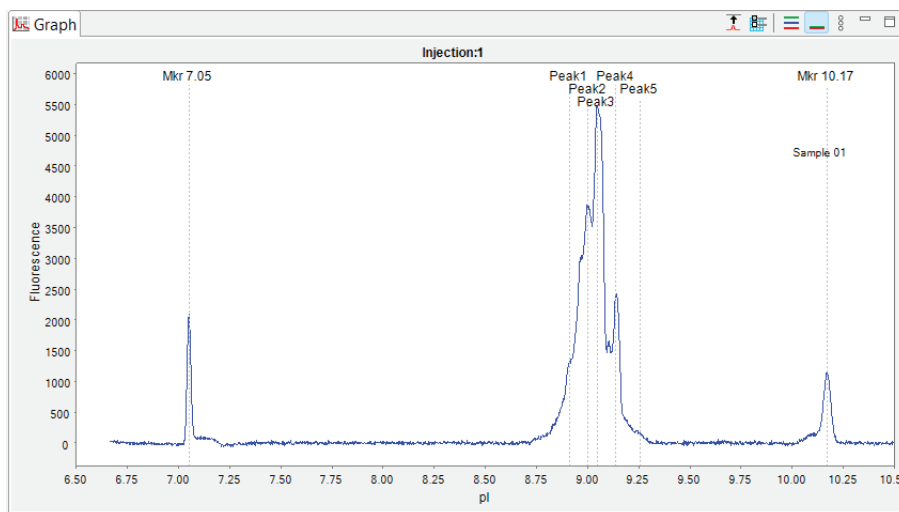


Selecting the Graph X-axis Range

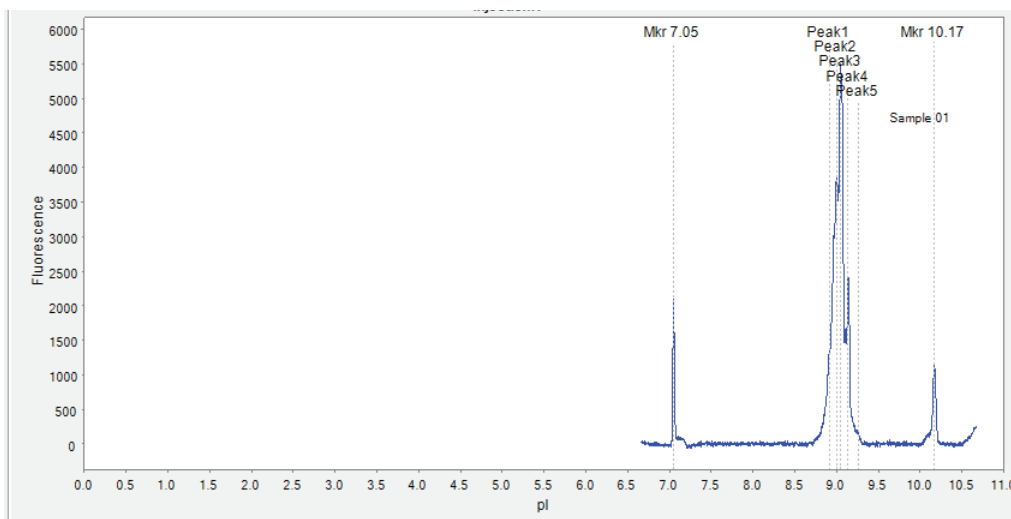
The pI range used for the x-axis can be changed. Just select **View** in the main menu and click **View Region**.



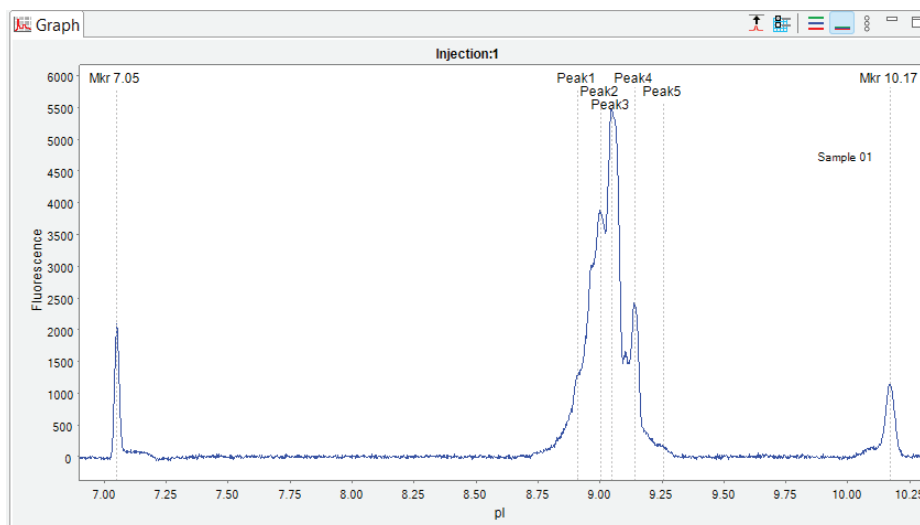
- **Analysis** sets the x-axis range of the electropherogram to what is selected in the Peak Fit range settings. To view or change these analysis settings, go to **Edit > Analysis** and click **Peak Fit** in the left sidebar. In this example, the lower and upper range settings are 6.5 and 10.5.



- **Full** displays the entire separation in the electropherogram. This is the default setting. In this example the lower and upper range settings are 0 and 11.



- **Custom** lets you manually enter the lower and upper range settings to display in the electropherogram. In this example the lower and upper range settings are 6.8 and 10.3.



NOTE: You can change the default x-axis range that Compass for iCE uses. Go to “Advanced Analysis Settings” on page 642 for more info.

Closing Run Files

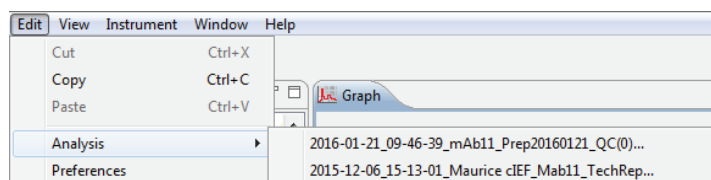
If more than one run file is open, you can close just one file or all the open files at the same time.

- **To close one run file** - In the Experiment pane, click on one of the sample rows in the file. Then click **File** from the main menu and click **Close**.
- **To close all open run files** - Select **File** from the main menu and click **Close All**.

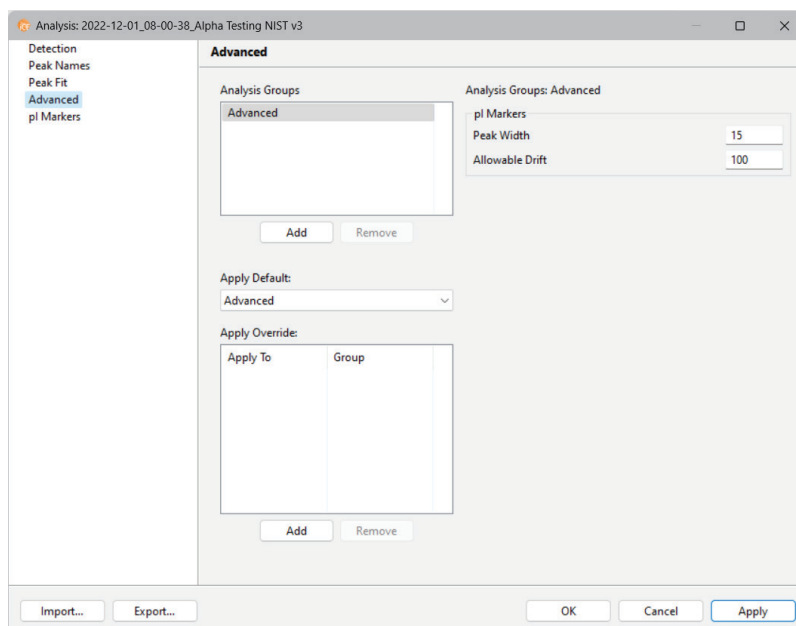
Analysis Settings Overview

Compass for iCE has many analysis features and settings that you can change to enhance your run data.

Select **Edit** in the main menu and click **Analysis**. If more than one run file is open, select the run file you want to view settings for from the list:



This opens the Analysis window:



To move between pages in the window, click on an option in the left sidebar.

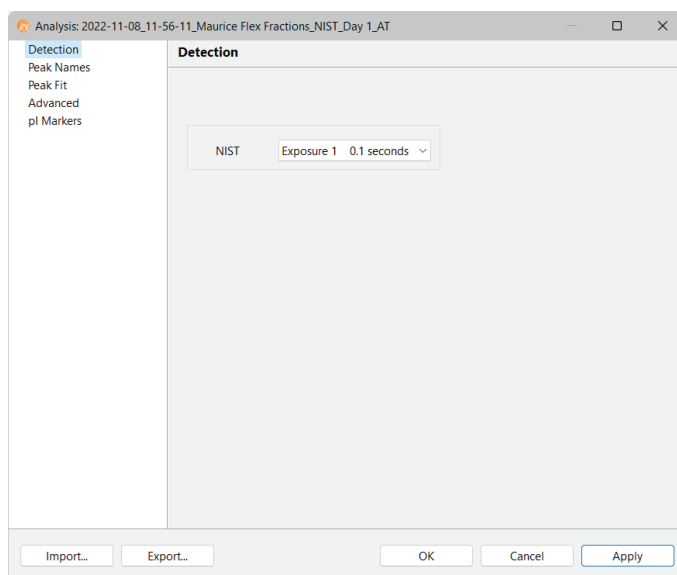
- **Detection** - Lets you view data at different fluorescence exposures.
- **Peak Names** - Lets you enter custom naming settings for sample proteins and have Compass for iCE automatically label the peaks in the run data.
- **Peak Fit** - Lets you customize peak fit settings for sample data.
- **Advanced** - Lets you customize analysis settings for the pI markers.
- **pI Markers** - Lets you customize the pI markers and positions Compass for iCE identifies for each method in your run.

On all pages in the Analysis window:

- Click **Import** to import an analysis settings file. Go to “Importing Analysis Settings” on page 657 to learn how to do this.
- Click **Export** to export the current analysis settings file. Go to “Exporting Analysis Settings” on page 657 to learn how to do this.
- Click **Apply** to apply changes to the run file and update results in real time.
- Click **OK** to save changes to the run file and exit.
- Click **Cancel** to exit without saving changes.

Detection Settings

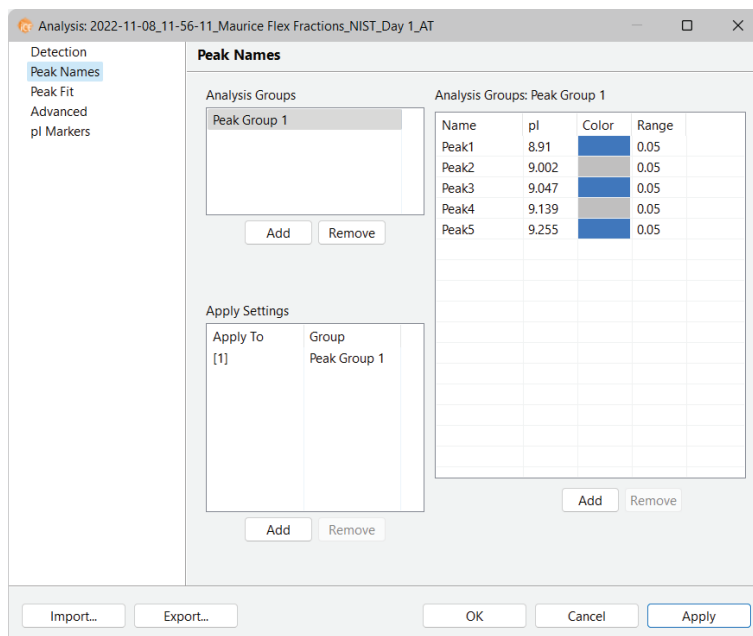
This page lets you see the native fluorescence exposures taken during the run, and select different exposures for data viewing in the Analysis screen. Select **Edit** in the main menu and click **Analysis**, then click **Detection** in the left sidebar.



Peak Names Settings

This page lets you view and change custom naming settings for sample proteins. Select **Edit** in the main menu and click **Analysis**, then click **Peak Names** in the left sidebar.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.

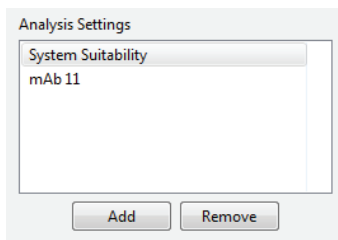


Peak Names Analysis Settings Groups

Peak name settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTE: Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 657.

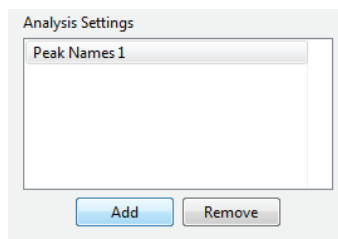
Peak name groups are displayed in the analysis settings box:



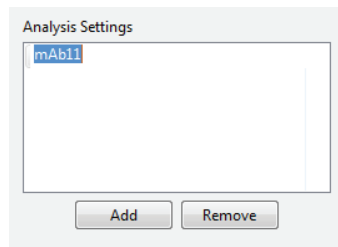
There aren't any Compass for iCE default settings groups, but you can make changes to groups you've created and create new groups. To view settings for a group, click on the group name in the analysis settings box.

Creating a Peak Names Group

1. Select **Edit > Analysis**, and select **Peak Names** in the left sidebar.
2. Click **Add** under the analysis settings box.



3. Enter a new name for the group.




4. Click in the first cell in the **Name** column in the analysis settings peak table and enter a sample protein name.

[illegible]

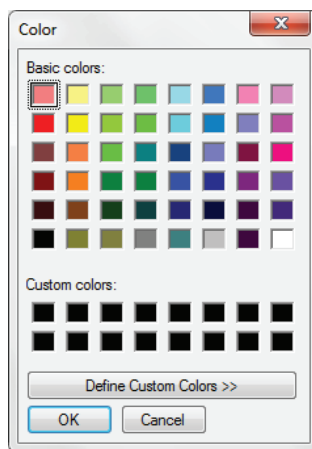
5. Click in the first cell in the **pI** column and enter the expected pI for the sample protein.

Name	pI	Color	Range
Peak1	5.55		0.05

6. Click in the first cell in the **Color** column, then click the button.


Name	pI	Color	Range
Peak1	6.55	 (0,1 ...)	0.05

The color selection box displays:




7. The color you pick is used to identify the sample protein peak in the Peaks and Injections panes in the Analysis screen. Click a color or define a custom color and click OK. The color selection will update in the table:

Analysis Settings: mAb11

Name	pI	Color	Range
Peak1	6.55		0.05








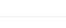
8. Click in the first cell in the **Range** column.

Analysis Settings: mAb11

Name	pI	Color	Range
Peak1	6.55		0.1

9. Enter a \pm range for the pI entered. Compass for iCE will automatically name peaks found within this pI range. For example, if the pI entered is 2 and a 0.1 range is used, all peaks with pIs between 1.9 and 2.1 will be identified with this peak name and color.
10. To add another sample protein, click Add under the peak table. Repeat the previous steps for other sample proteins. In this example, eight proteins were entered:

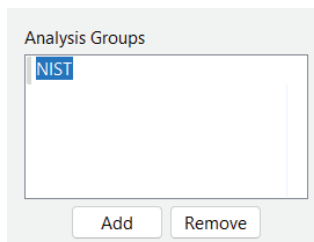
Analysis Settings: mAb 11

Name	pI	Color	Range
Peak1	6.55		0.1
Peak2	6.65		0.1
Peak3	6.8		0.1
Peak4	6.9		0.1
Peak5	7		0.1
Peak6	7.1		0.1
Peak7	7.2		0.1
Peak8	7.3		0.1

11. To remove a sample protein, select its row and click **Remove**.
12. Click **OK** to save changes.

Modifying a Peak Names Group

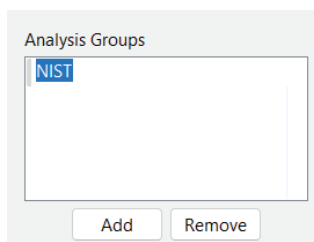
1. Select **Edit > Analysis**, then click **Peak Names** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



3. Change the information in the analysis settings peak table as described in “Creating a Peak Names Group” on page 623.
4. Click **OK** to save changes.

Deleting a Peak Names Group

1. Select **Edit > Analysis**, then click **Peak Names** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.

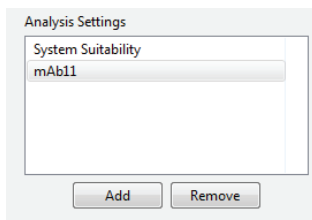


3. Click **OK** to save changes.

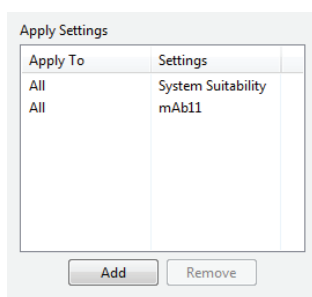
Applying Peak Names Groups to Run Data

1. Select **Edit > Analysis**, then click **Peak Names** in the options list.

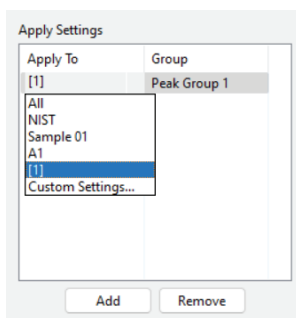
- Click on the group in the analysis settings box you want to apply to specific run data.



- Application of peak names groups to specific run data is done in the apply settings box. A default data set automatically gets created whenever you create a new group and it's applied to all injections in the run. You can either modify the default group or click Add under the box to create a new one.

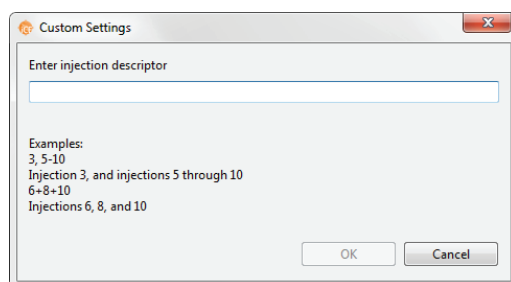


- Click the cell in the Apply To column, then click the down arrow.

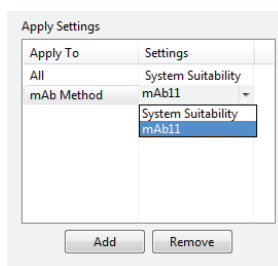


- Select an option from the drop down list. This applies the peak names group selected to specific run data as follows:
 - All** - Selecting this applies peak names group settings to all injections.
 - Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.
 - Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.

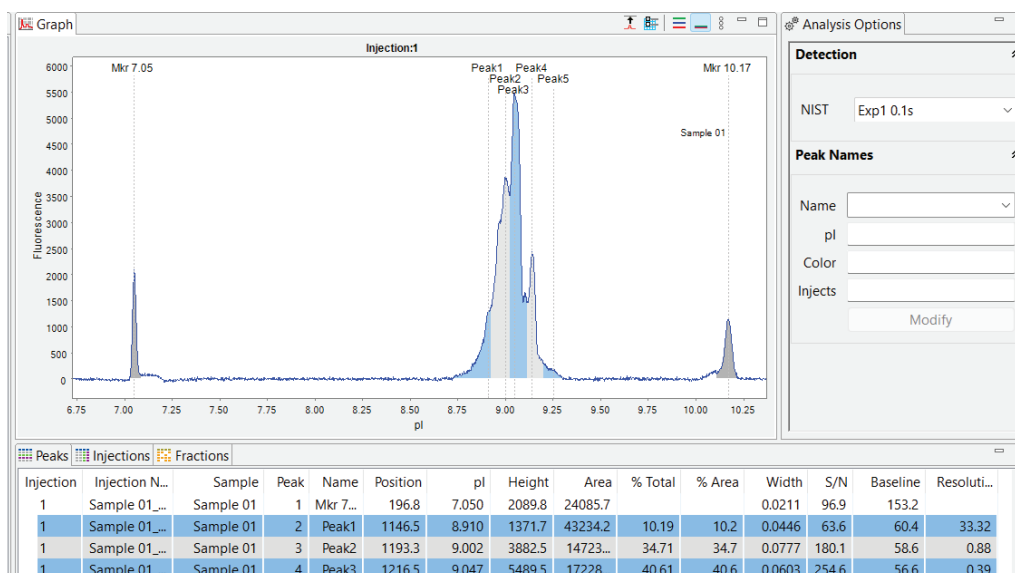
- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the peak names group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click **OK** to save changes. Named peaks will be identified with a peak name label in the electropherogram and color-coded in the Peaks and Injections panes:



Peak Fit Analysis Settings

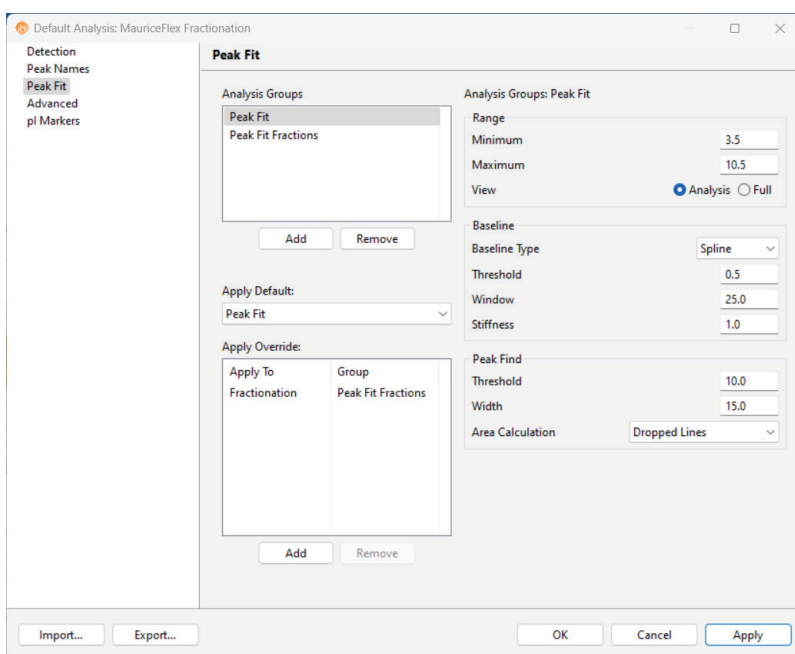
This page lets you view and change peak fit settings for sample data. Select **Edit** in the main menu and click **Analysis**, then click **Peak Fit** in the left sidebar:

IMPORTANT:

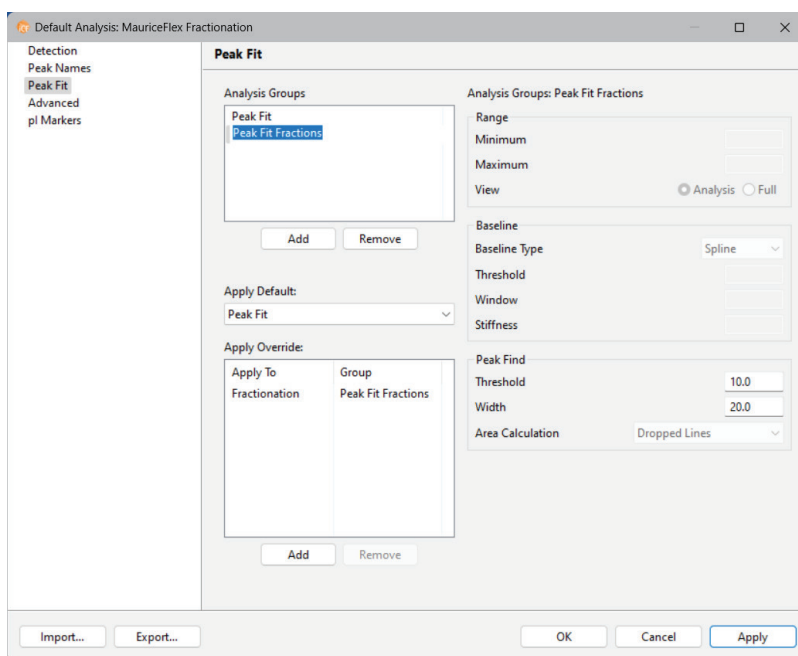
There are two default Peak Fit analysis settings for MauriceFlex Fractionation runs. One set is applied to the final sample injection focusing image and the other set is applied to Fractions data to account for peak shape changes during mobilization.

Only Peak Find settings (threshold and width) can be adjusted when defining analysis settings for Fractions.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



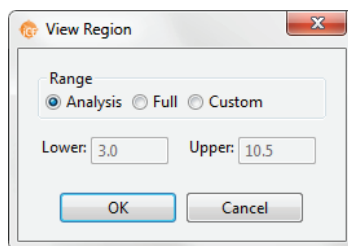
Default Peak Fit settings for sample injection data.



Default Peak Fit settings for Fractions data.

Range Settings

- **Minimum** - The pI value below which peaks won't be identified. This value is also used as the default lower pI range for data displayed in the electropherogram.
- **Maximum** - The pI value above which peaks won't be identified. This value is also used as the default upper pI range for data displayed in the electropherogram.
- **View** - Sets the default range to either Full or Analysis for the electropherogram x-axis range in the View Region window (select **View** in the main menu and click **View Region**).



- **Analysis** sets the x-axis range of the electropherogram to the Peak Fit range minimum and maximum settings in the electropherogram. This is the default setting.
- **Full** displays the entire separation range of the run data in the electropherogram.

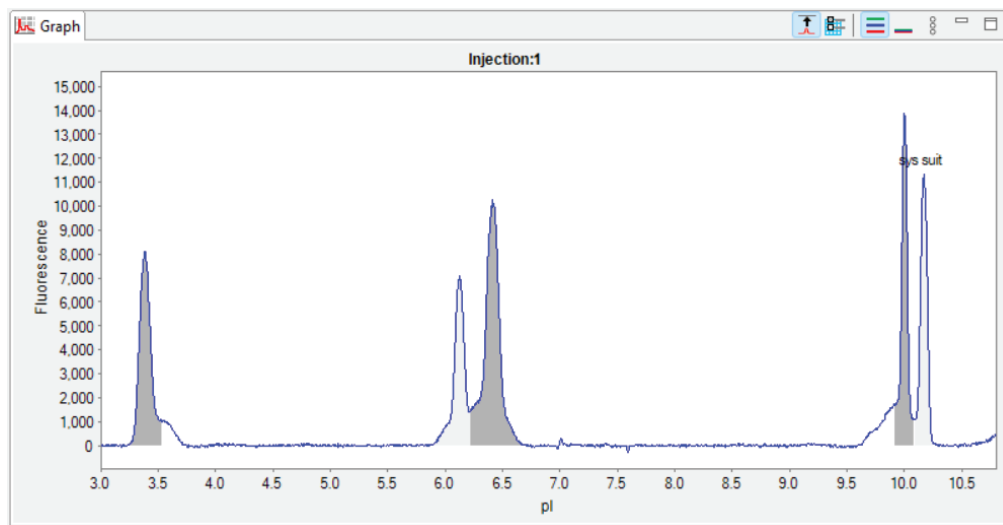
Baseline Settings

These settings apply to spline baselines only.

- **Threshold** - The variance, or roughness, in a baseline data segment below which a point is called part of the baseline.
- **Window** - How long baseline data segments are expected to be in pixels. Shorter segments let the baseline follow plateau sections of the signal.
- **Stiffness** - The amount the baseline is allowed to vary from a straight line. Settings between 0.1 and 1.0 make the baseline fit closer to a straight line. Settings from 1.0 to 10.0 will make the baseline fit follow the data more closely.

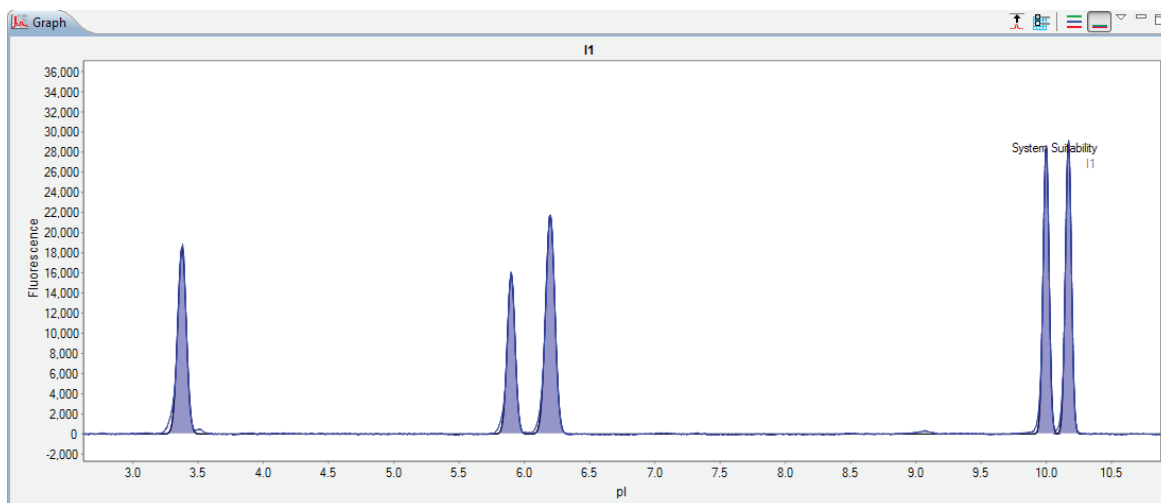
Peak Find Settings

- **Threshold** - The minimum signal to noise ratio required for a peak to be identified. A setting of 1.0 will detect many peaks, a setting of 10.0 will detect fewer peaks.
- **Width** - The approximate peak width (at full width half max) in pixels used to detect peaks. The minimum value for this setting is 3.0. Larger widths help eliminate the detection of shoulder and noise peaks.
- **Area Calculation** - Two fits are used, either Gaussian Fit or Dropped Lines. These settings can be changed before or after the run is finished.
 - For MauriceFlex Fractionation applications, peak area is calculated using Dropped Line by default.



- The Dropped Line method type of area calculation is also often called the perpendicular drop method. This is the preferred method when peaks overlap or are close to each other. It draws two vertical lines from the left and right bounds of the peak down to the x-axis and then measures the total area bounded by the signal curve, the x-axis ($y=0$ line), and the two vertical lines.

- The next view is of the similar data using the Gaussiann peak fit method instead.



Peak Fit Analysis Settings Groups

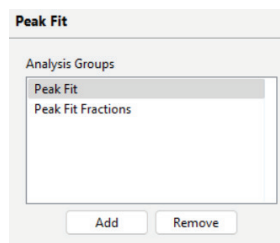
Peak fit settings are saved as a group, and you can create multiple settings groups. Specific group settings can then be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values for peak fit analysis settings. These settings are included in the default Peak Fit group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 657.

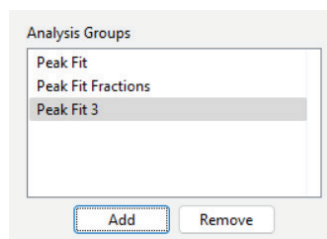
Peak fit groups are displayed in the analysis settings box:



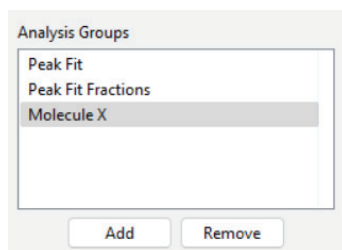
The Peak Fit group shown contains the Compass for iCE default analysis settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Peak Fit Group

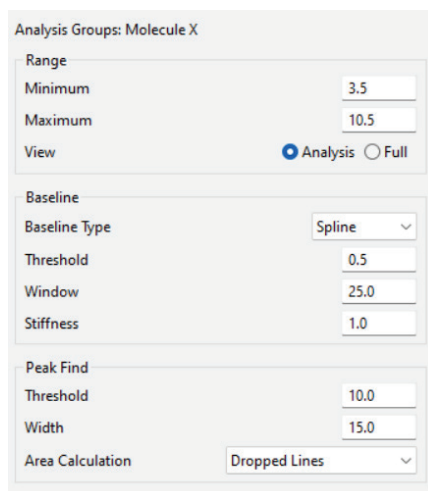
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



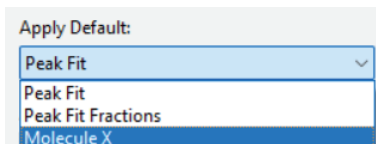
3. Click on the new group and enter a new name.



4. Change the settings in the range, baseline or peak find boxes as needed.

The image shows a settings dialog box titled "Analysis Groups: Molecule X". It has three main sections: "Range", "Baseline", and "Peak Find".
- The "Range" section has "Minimum" set to 3.5 and "Maximum" set to 10.5. There is a "View" section with radio buttons for "Analysis" (selected) and "Full".
- The "Baseline" section has "Baseline Type" set to "Spline", "Threshold" set to 0.5, "Window" set to 25.0, and "Stiffness" set to 1.0.
- The "Peak Find" section has "Threshold" set to 10.0, "Width" set to 15.0, and "Area Calculation" set to "Dropped Lines".

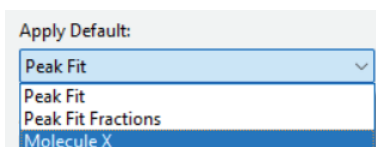
5. To use the new group as the default peak fit settings for the run file data, click the arrow in the drop down list next to Apply Default, then click the new group from the list. Peak fit settings in the new group will then be applied to the run data.



6. Click OK to save changes.

Changing the Default Peak Fit Group

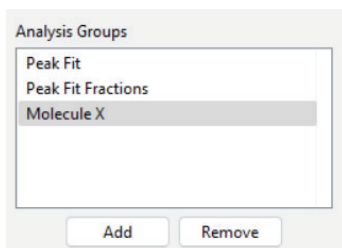
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then click a new default group from the list.



3. Click OK to save changes. Peak fit settings in the group selected will be applied to the run data.

Modifying a Peak Fit Group

1. Select **Edit > Analysis**, and click **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



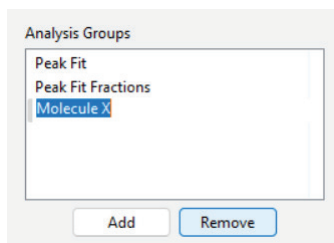
3. Change the settings in the range, baseline or peak find boxes as needed.

The screenshot shows a settings dialog box titled "Analysis Groups: Molecule X". It is divided into three sections: Range, Baseline, and Peak Find. The Range section has input fields for Minimum (3.5) and Maximum (10.5), and radio buttons for View (Analysis is selected, Full is unselected). The Baseline section has a dropdown for Baseline Type (Splines), and input fields for Threshold (0.5), Window (25.0), and Stiffness (1.0). The Peak Find section has input fields for Threshold (10.0) and Width (15.0), and a dropdown for Area Calculation (Dropped Lines).

4. Click OK to save changes. The new peak fit settings will be applied to the run data.

Deleting a Peak Fit Group

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.

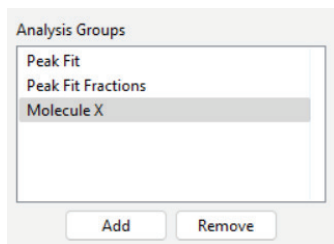


3. Click OK to save changes.

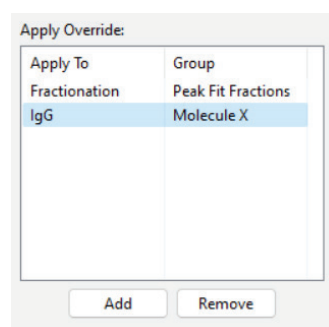
Applying Peak Fit Groups to Specific Run Data

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.

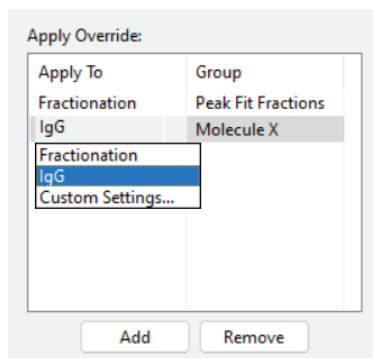
- Click on the group in the analysis settings box you want to apply to specific run data.



- Application of analysis groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.



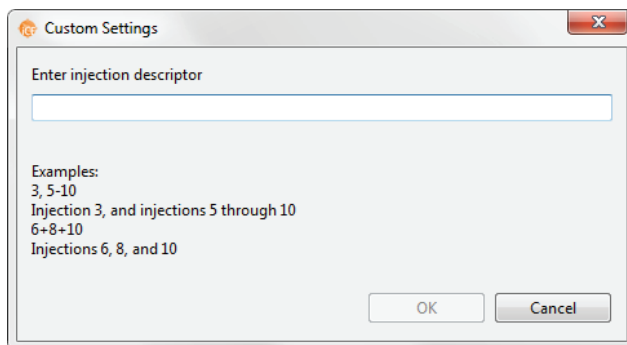
- Click the cell in the **Apply To** column, then click the down arrow.



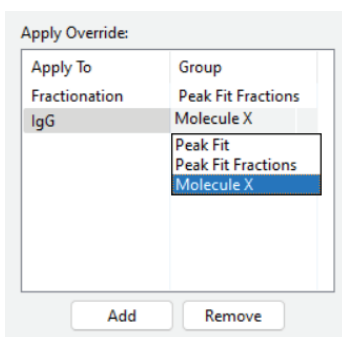
- Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
 - **Fractionation** - All mobilization electropherograms in the run file. Selecting Fractionation applies group settings to all Fractions data.

NOTE: Only the Peak Find Threshold and Width can be adjusted when an Analysis Group is applied to a Fractionation settings group.

- **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
- **Custom settings** - Lets you choose specific electropherograms to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



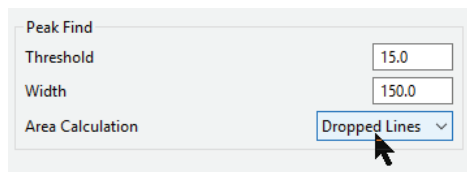
7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click OK to save changes.

Manual Peak Integration

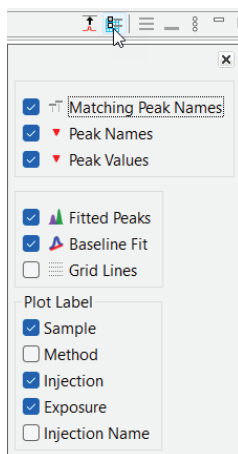
Compass for iCE lets you manually integrate peaks in individual electropherograms when in the Samples view.

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.

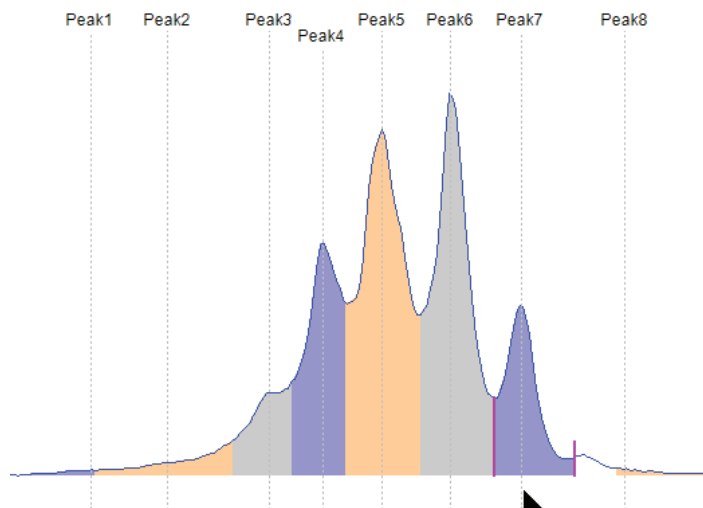
2. Select Dropped lines as the area calculation if it isn't already selected.



3. Select Fitted Peaks in the Graph Options.

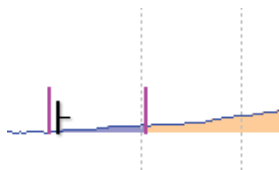



4. Select an injection in the Experiment pane.
5. Hover the mouse over the peak you want to adjust. Two magenta bars will display that indicate the start and end points of its integration.

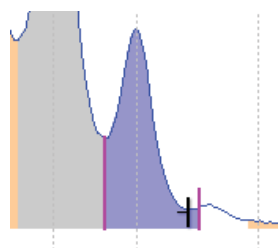


6. To change the integration, hover the mouse over either of the magenta bars. Then click the mouse and drag the bar to move it.

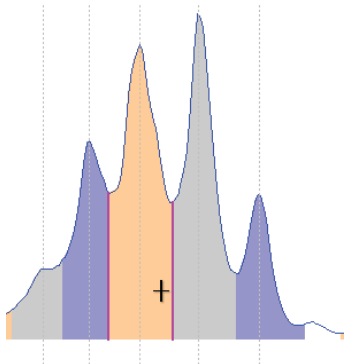
- If the cursor changes to  this is the peak start for the peak on the right.



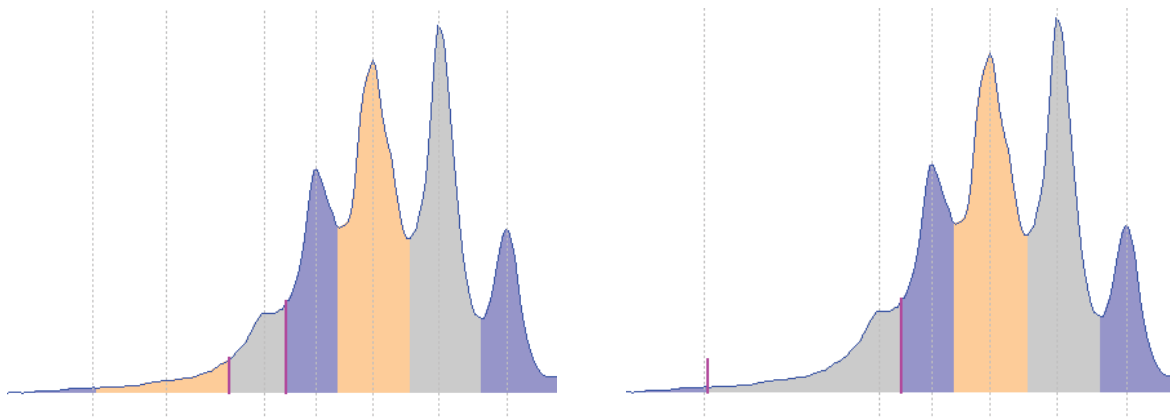
- If the cursor changes to  this is the peak end for the peak on the left.



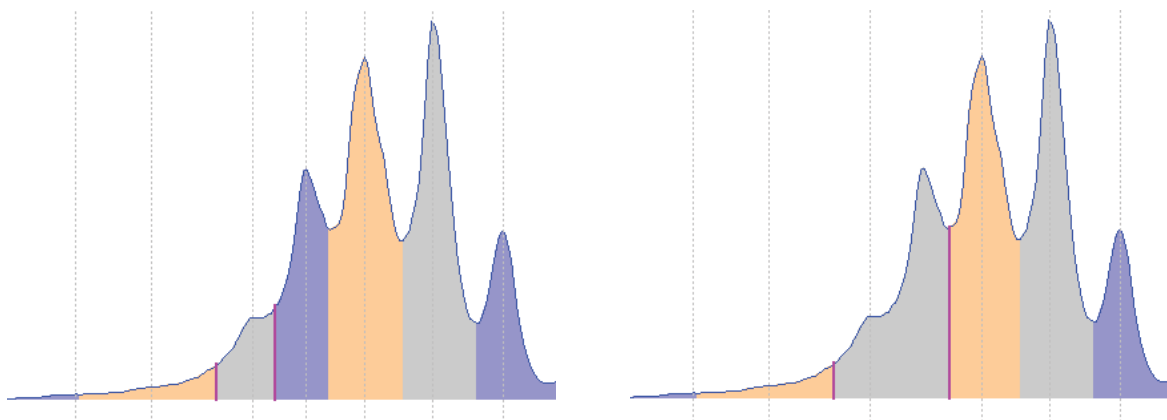
- If the cursor changes to  this is a joint boundary for the peaks on the left and right.


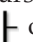



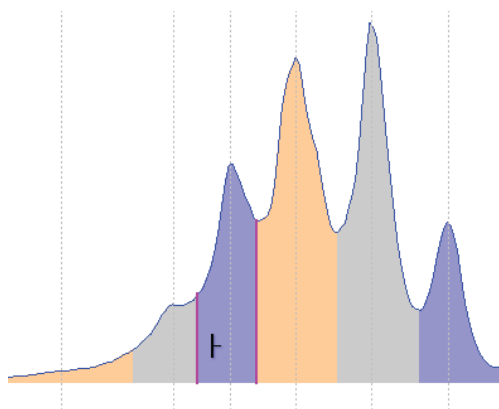
In the example below, we moved the start and end points of the peak to include more area under the peak:



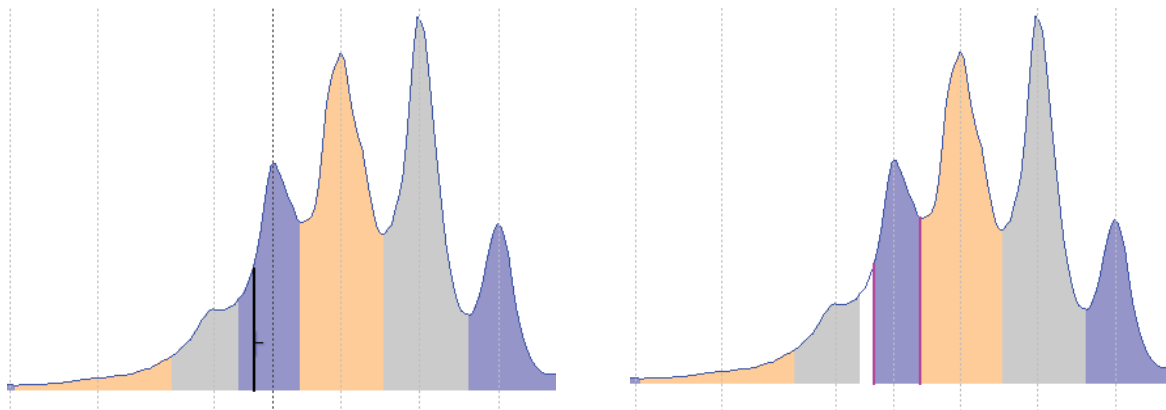
You can also combine peaks by dragging the magenta lines left or right:



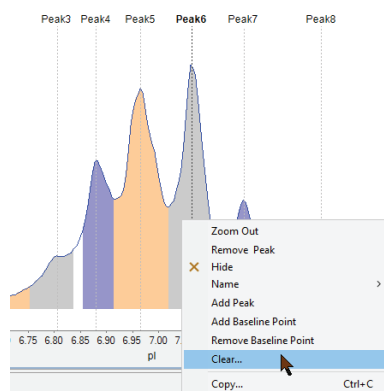
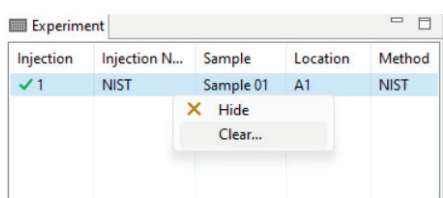
You can also separate areas between peaks. Whenever you have a  cursor between two peaks that aren't baseline resolved, move the mouse slightly to the right or left until you get the  or  cursor.



Click and hold the mouse. A black drop line will display indicating where the current peak start or end point is. While holding the mouse, drag the cursor to the right or left to move the start or end point to where you want it, then release the mouse.



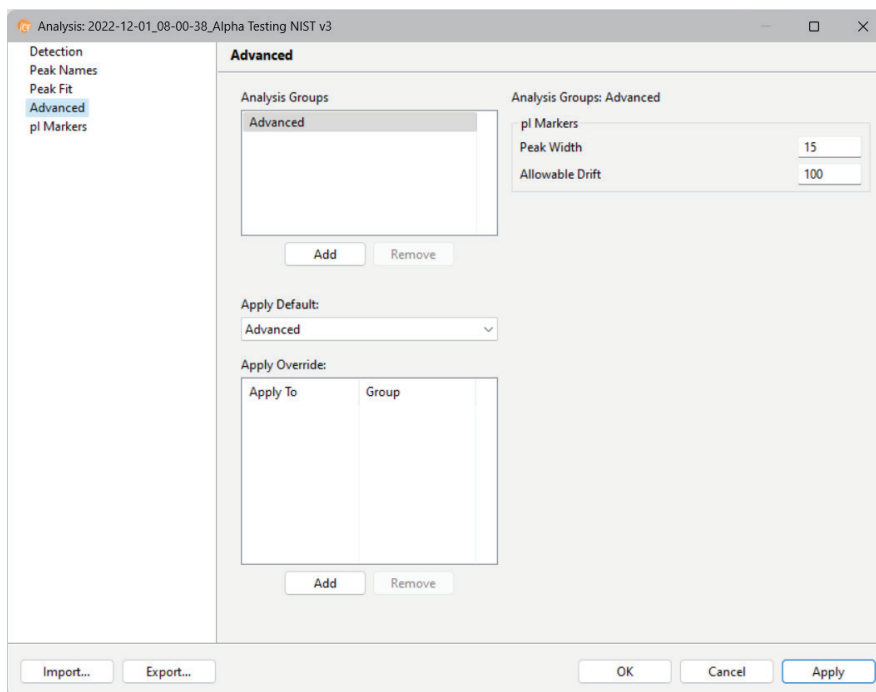
7. To remove the manual integration settings, right click on the graph or on the injection in the Experiment pane and select **Clear**.



Advanced Analysis Settings

This page lets you view and change analysis settings for the pI marker data. Select **Edit** in the main menu and click **Analysis**, then click **Advanced** in the left sidebar:

NOTE: Settings can be changed in batches before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



pI Markers Settings

- **Peak Width** - The approximate width (at full width half max) used to filter out fluorescence artifacts which improves recognition of pI markers.
- **Allowable Drift** - The distance the pI marker(s) are expected to move compared to the position entered on the pI Markers page. This setting helps with recognition of the pI marker.

Advanced Analysis Settings Groups

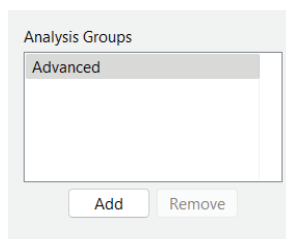
Advanced analysis settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values for advanced analysis settings. These settings are included in the default Advanced group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. See “Importing and Exporting Analysis Settings” on page 657 for more info.

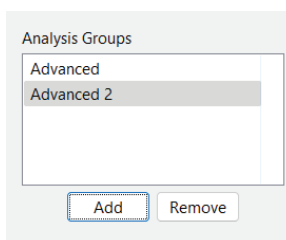
Analysis groups are displayed in the analysis settings box:



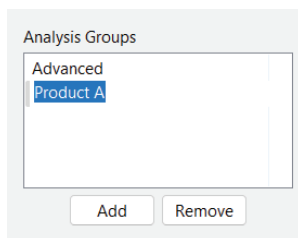
The Advanced group shown contains the Compass for iCE default analysis settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



3. Click on the new group and enter a new name.



4. Change the settings in the Markers box as needed.

Analysis Settings: Product A

pl Markers

Peak Width

Allowable Drift

5. To use the new group as the default analysis settings for the run data, click the arrow in the drop down list next to Apply Default, then click the new group from the list. Analysis settings in the new group will then be applied to the run data.

Apply Default:

▼

Advanced

Product A

Apply To	Settings

6. Click OK to save changes.

Changing the Default Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then click a new default group from the list.

Apply Default:

▼

Advanced

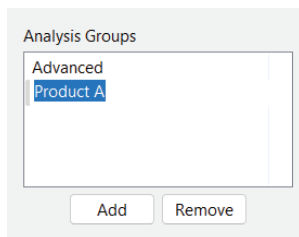
Product A

Apply To	Settings

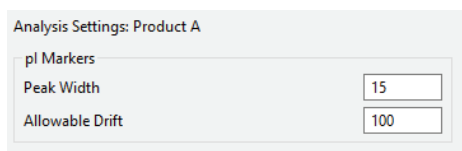
3. Click **OK** to save changes. Analysis settings in the group selected will be applied to the run data.

Modifying an Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



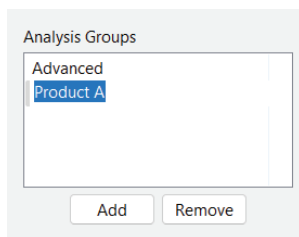
3. Change the settings in the Markers box as needed.



4. Click **OK** to save changes. The new analysis settings will be applied to the run data.

Deleting an Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.

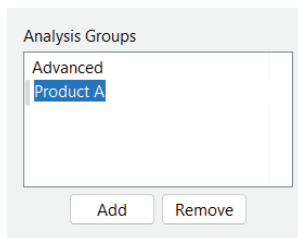


3. Click **OK** to save changes.

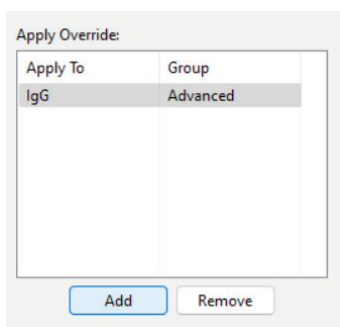
Applying Analysis Groups to Specific Run Data

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.

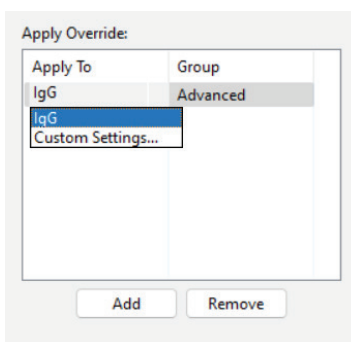
- Click on the group in the analysis settings box you want to apply to specific run data.



- Application of analysis groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

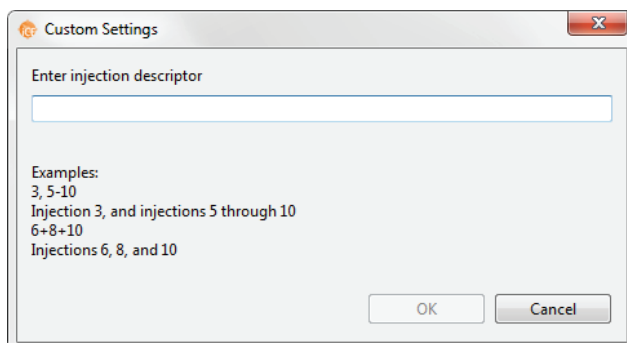


- Click the cell in the **Apply To** column, then click the down arrow.

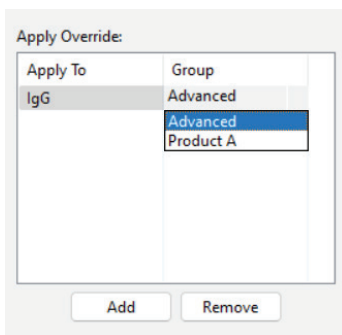


- Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
 - Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.

- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.

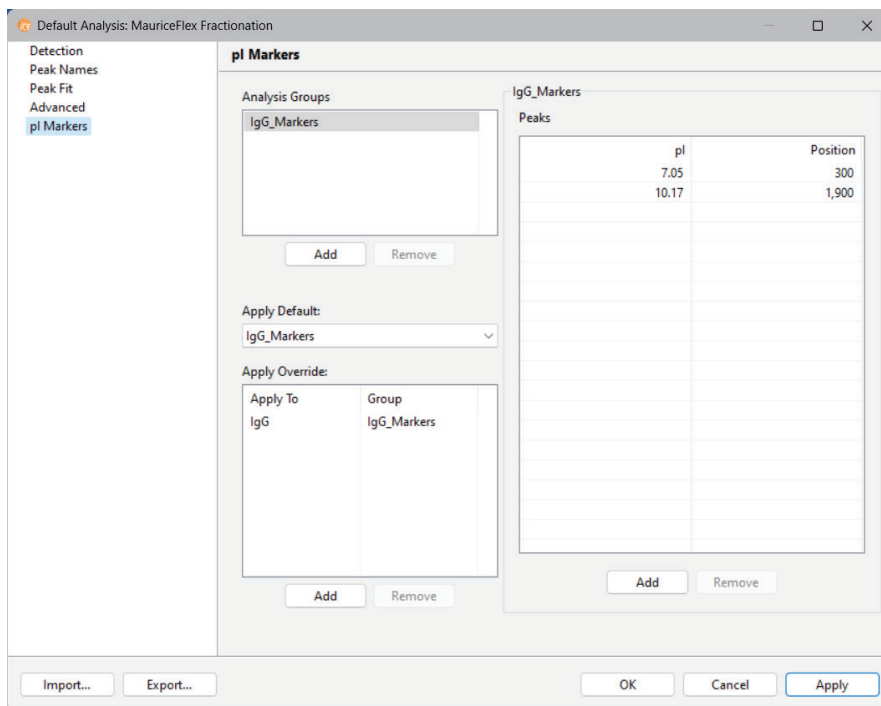


7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click **OK** to save changes.

pI Markers Analysis Settings

This page lets you define the pI and position of the pI Markers you're using in your samples. Select **Edit** in the main menu and click **Analysis**, then click **pI Markers** in the left sidebar.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



pI Markers Analysis Settings Groups

pI marker settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

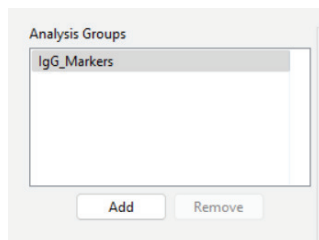
NOTES:

We recommend using the Compass for iCE default values. These settings are included in the default Markers group.

When you edit the pI markers in the method for a batch, Compass for iCE automatically creates a Markers group in the pI Markers Analysis settings for you.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see "Importing and Exporting Analysis Settings" on page 657.

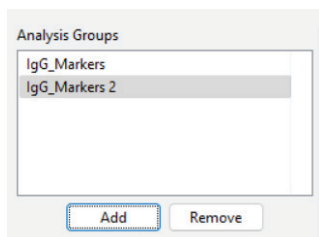
Markers groups are displayed in the analysis settings box:



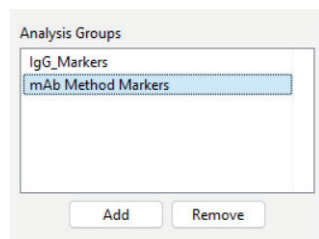
The Markers group shown uses the Compass for iCE default settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Markers Group

1. Select **Edit > Analysis**, and select **pI Markers** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



3. Click on the new group and enter a new name.



4. The default Maurice cIEF pI marker pI and position values are already populated in the pI Marker Peaks table. If you'd like to use these values, skip to the next step. If you're using different markers, here's how to change the values:

- a. Click in the first cell in the pI column in the table and enter the pI for the marker.

pI Markers

Peaks

pI	Position
4.05	250
10.17	1,800

Add Remove

- b. Click in the first cell in the Position column and enter a value for the marker.

pI Markers

Peaks

pI	Position
4.05	500
10.17	1,800

Add Remove

NOTE: pI marker peak positions are relative to each other. Only the difference in position is used to help identify them. When entering pI marker peak information for the first time, review the marker data in the Analysis screen to find the correct peak positions.

- c. Repeat the steps above for the remaining markers in the table.
- **To add another marker** - Click **Add** under the table, then change the information in the new row.
 - **To remove a marker** - Select its row and click **Remove**.
5. To use the new group as the default settings for the run, click the arrow in the drop down list next to Apply Default, then click the new group in the list. The settings in the new group will then be applied to the run data.

Apply Default:

mAb Method Markers

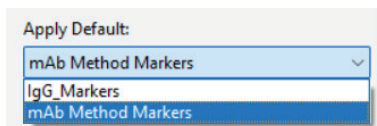
IgG_Markers

mAb Method Markers

6. Click **OK** to save changes.

Changing the Default Markers Group

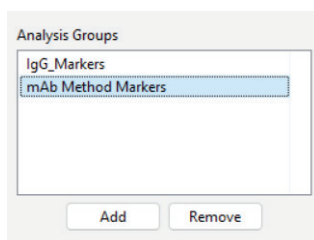
1. Select **Edit > Analysis**, and click **pI Markers** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then select a new default group from the list.



3. Click **OK** to save changes. Analysis settings in the group selected will be applied to the run data.

Modifying a Markers Group

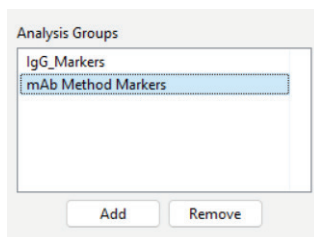
1. Select **Edit > Analysis**, and click **Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



3. Change the marker info as needed as in **Creating a New Markers Group**.
4. Click **OK** to save changes. The new analysis settings will be applied to the run data.

Deleting a Markers Group

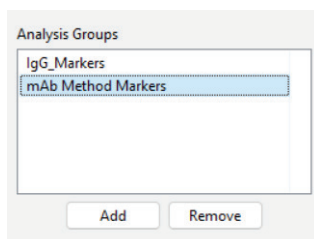
1. Select **Edit > Analysis**, and click **pI Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



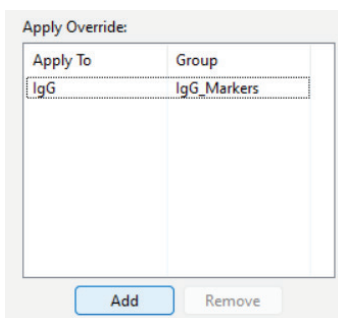
3. Click **OK** to save changes.

Applying Markers Groups to Specific Run Data

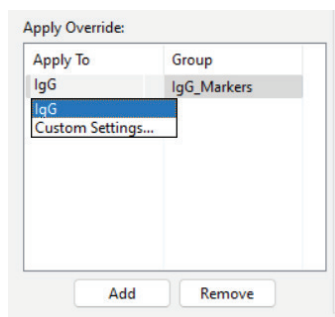
1. Select **Edit > Analysis**, and select **pI Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to apply to specific run data.



3. Application of markers groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

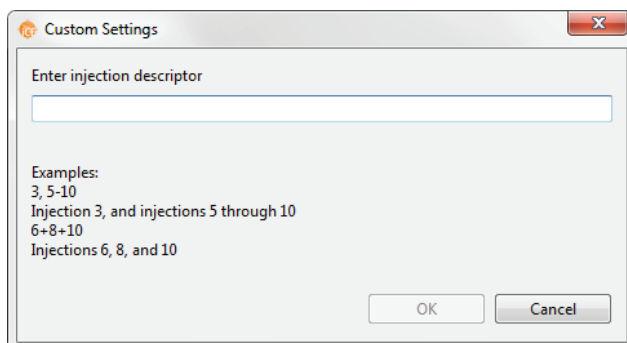


4. Click the cell in the **Apply To** column, then click the down arrow.

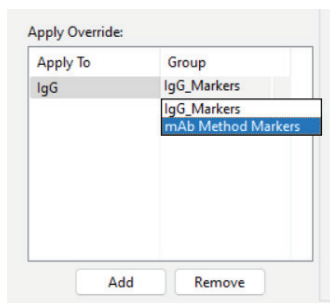


5. Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
 - **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.

- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click OK to save changes.

Injection Reports

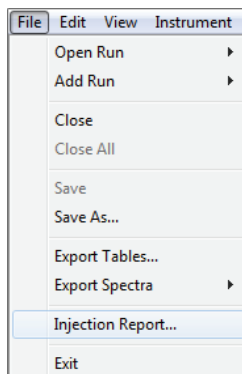
You can export PDF files of the raw and analyzed data, IV plot, peaks table, sample and system info for individual or all injections and mobilization electropherograms in a run file. You can also export the run history with all analysis events.

NOTES:

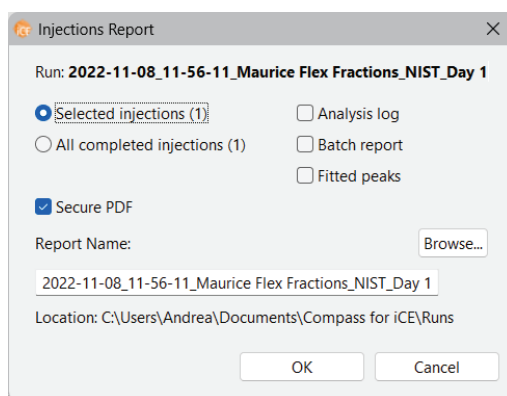
You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it’s saved in a separate run folder. Existing reports aren’t overwritten.

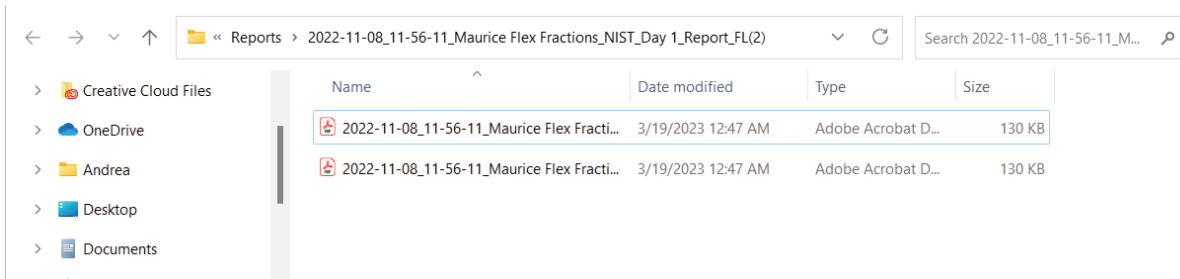
1. Select **File** from the main menu in either screen and click **Injection Report**.



2. In the Injection Reports window:
 - a. Choose either **Selected injections** or **All injections**. Either option will give you the same report since there is only one sample injection in a MauriceFlex Fractionation run.
 - b. Select **Analysis log** if you want a run history report with all analysis events.
 - c. Select **Batch Report** if you want to include the sample and method details for each injection in the batch.
 - d. Select **Fitted peaks** if you want to show peak fitting in the electropherograms.
 - e. Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don't want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.
 - f. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.
 - g. Click **OK**.

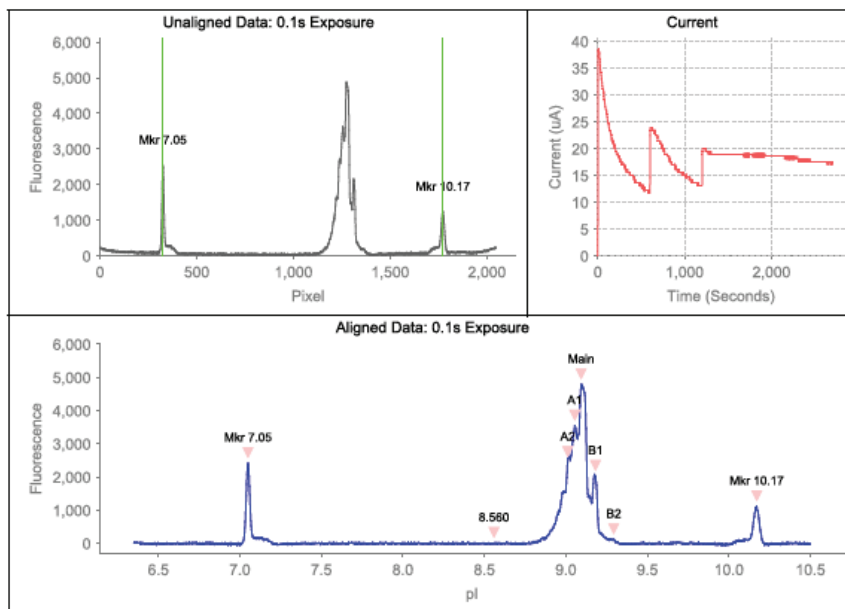


- The individual injection report PDF(s) and a combined injection report PDF are exported to the Runs folder in the Compass for iCE directory. They'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Example Analysis and Injection Report

Uncontrolled Injection 1: NIST



Fluorescence Peaks: 0.1s Exposure

Peak	Name	Position	pI	Height	Area	%Total	%Area	Width	S/N	Baseline	Resolution
1	Mkr 7.05	324.9	7.050	2423.4	31491.1			0.0251	125.7	86.6	
2		1025.5	8.560	-16.4	0.0			0.0022	-0.9	15.4	65.29
3	A2	1235.7	9.013	2631.5	75324.3	24.29	24.3	0.0544	136.5	54.3	9.44
4	A1	1255.2	9.055	3548.8	68620.7	22.13	22.1	0.0453	184.0	55.2	0.49
5	Main	1273.5	9.094	4789.8	121793.8	39.28	39.3	0.0563	248.4	55.2	0.46
6	B1	1313.5	9.181	2084.7	37366.5	12.05	12.0	0.0425	108.1	51.8	1.03
7	B2	1365.7	9.293	303.3	6991.9	2.25	2.3	0.0841	15.7	40.1	1.05
8	Mkr 10.17	1772.5	10.170	1119.3	22427.8			0.0398	58.0	102.3	8.34

Uncontrolled Injection 1: NIST

Peak Predictions

Peak	Name	Predicted Wells
1	Mkr 7.05	
2	8.560	C10, C9, C8, C7, C6
3	A2	B11, B12, C12, C11, C10
4	A1	B11, B12, C12, C11
5	Main	B10, B11, B12, C12
6	B1	B9, B10, B11, B12
7	B2	B7, B8, B9, B10, B11
8	Mkr 10.17	B1, B2

Sample Information

Injection Name	NIST
Sample ID	Sample 01
Location	Plate Well A1
Batch Name	Alpha Testing NIST
Run Started	Thu 8:01 AM Dec 1, 2022 PST
Run Completed	Thu 9:50 AM Dec 1, 2022 PST
Date Acquired	Thu 8:13 AM Dec 1, 2022 PST
Run Error	None
Reinjection	No

Injection Conditions

Guaranteed Injection	Yes
On-board Mixing	Off
Focus Period 1	500V for 10.0 min
Focus Period 2	1000V for 10.0 min
Focus Period 3	1500V for 25.0 min
Detection Exposure	0.1 sec
Detection Interval	5.0 min
Sample Load Duration	30.0 Seconds
pI marker 1	7.05
pI marker 2	10.17
Tray Temperature	11.0°C

Fractionation Conditions

Mobilization	1000 Volts for 25.0 min
Refocus	1500 Volts for 0.0 min
Fractions	1000 Volts for 45.0 sec
Detection Exposure	0.1 sec
Detection Interval	1.0 min

Maurice Settings

Model	MauriceFlex
Instrument S/N	mm0008
Software Version	Compass for iCE 4.0.0, Build ID: 0222
Firmware Version	4.2.2022.11.03.21.24.04.b15d551f8
Tray Type	96-well plate
Cartridge Type	MauriceFlex
Cartridge S/N	5221128022
Cartridge Expiration	Nov 2023
Injections Remaining	16 (16 guaranteed)
Batches Remaining	16

Importing and Exporting Analysis Settings

The analysis settings in a run file can be exported as a separate file. This allows the same analysis settings to be imported into other batches or run files at a later time, rather than you having to re-enter them manually.

Importing Analysis Settings

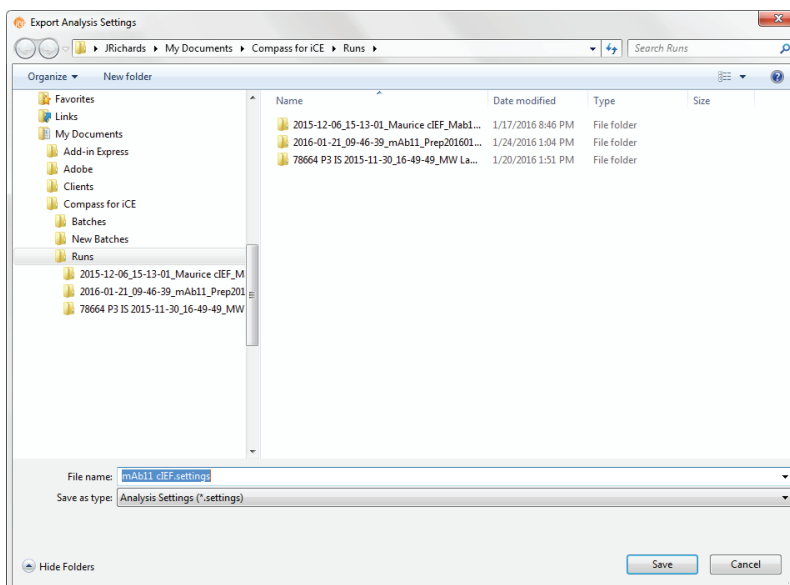
NOTE: Importing an analysis settings file populates the settings in all analysis pages.

1. Open the run file or batch you want to import analysis settings to.
2. Select **Edit** in the main menu and click **Default Analysis** (Batch screen) or **Analysis** (Analysis screen).
3. Click **Import** on any page.
4. Select a settings file (*.settings) and click **OK**. The imported settings will display in all analysis pages.

Exporting Analysis Settings

NOTE: Exporting an analysis settings file exports the settings in all analysis pages.

1. Open the run file or batch you want to export analysis settings from.
2. Select **Edit** in the main menu and click **Default Analysis** (Batch screen) or **Analysis** (Analysis screen).
3. Click **Export** on any page. The following window displays:



4. The default directory is Compass for iCE/Runs. Change the directory if needed.
5. Enter a file name and click **Save**. The settings will be saved as a *.settings file.

Chapter 20:

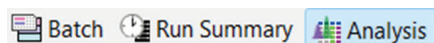
CE-SDS Data Analysis

Chapter Overview

- Analysis Screen Overview
- Opening Run Files
- How Run Data is Displayed
- Viewing Run Data
- Data Notifications and Warnings
- Checking Your Results
- Naming Peaks
- Group Statistics
- Copying Data Views and Results Tables
- Exporting Run Files
- Changing Sample Protein Identification
- Changing the Electropherogram View
- Changing the Virtual Gel View
- Closing Run Files
- Analysis Settings Overview
- Markers Analysis Settings
- Peak Names Settings
- Peak Fit Analysis Settings
- Manual Peak Integration
- Advanced Analysis Settings
- Signal Processing Settings
- Injection Reports
- Importing and Exporting Analysis Settings

Analysis Screen Overview

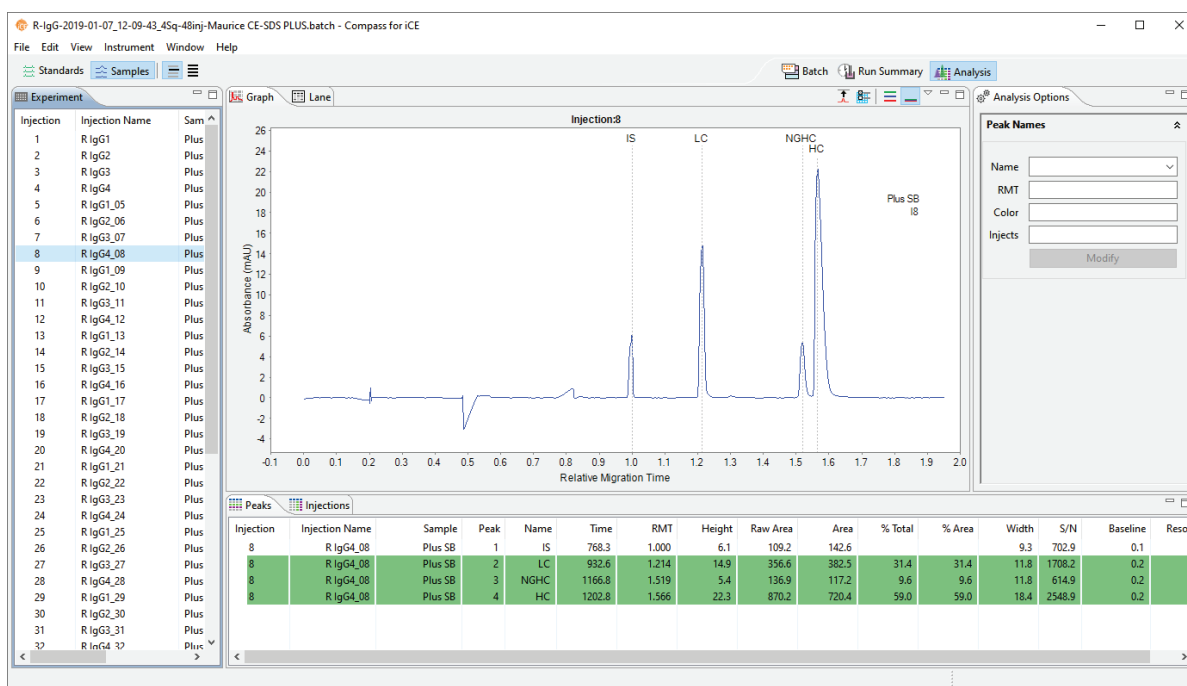
You can use the Analysis screen to view electropherograms, lane view data and tabulated results for your injections. If any post-run analysis is needed, you can do it here too. To get to this screen, click the **Analysis** screen tab:



Analysis Screen Panes

The Analysis screen has four panes:

- **Experiment** - Lists the injection number, sample IDs, sample locations and methods for each injection in the run and lets you get a quick view of method parameters.
- **Graph** - Displays the electropherograms for sample proteins or standards.
- **Lane** - Displays data for sample proteins as bands in individual lanes. This virtual gel-like image is similar to traditional gel results.
- **Peaks** - Shows the tabulated results for sample proteins, internal standards and CE-SDS MW Markers.
- **Injections** - Displays a list of the sample proteins Compass for iCE names automatically using the user-defined peak name analysis parameters.
- **Analysis Options** - Lets you view, change and add new custom peak name settings.



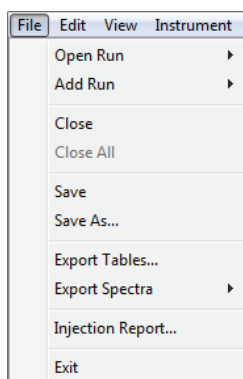
Software Menus Active in the Analysis Screen

These main menu items are active in the Analysis Screen:

- File
- Edit
- View
- Instrument (when Compass for iCE is connected to Maurice, Maurice C., Maurice S. or MauriceFlex)
- Window
- Help

File Menu

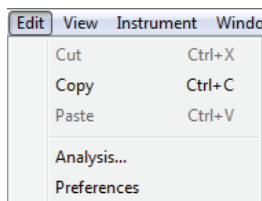
These File menu options are active:



- **Open Run** - Opens a run file.
- **Add Run** - Lets you open and view other run files besides the one that's already open.
- **Close** - Closes the run file currently being viewed.
- **Close All** - Closes all open run files.
- **Save/Save As** - If you made changes in the Analysis screen before you went to the Run Summary screen, this saves your changes to the run file.
- **Export Tables** - Exports the results for all injections in the run in .txt format.
- **Export Spectra** - Exports the raw and analyzed data traces and background for each injection in the run in .txt or .cdf format.
- **Injection Report** - Exports the raw and analyzed data, IV plot, peaks table, sample and system info for individual injections as PDF files. You can also export the run history with all analysis events.
- **Exit** - Closes Compass for iCE.

Edit Menu

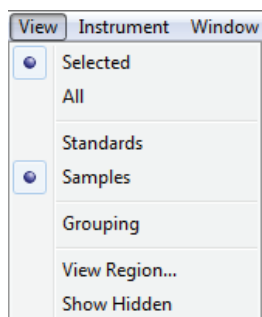
These Edit menu options are active:



- **Copy** - Lets you copy data shown in the graph, lane, peaks or injections panes. See “Copying Data Views and Results Tables” on page 688 for more information.
- **Analysis** - Displays the analysis settings used to analyze the run data and lets you change them as needed. See “Analysis Settings Overview” on page 713 for more information.
- **Preferences** - Lets you set and save your preferences for data export, graph colors, grouped data and Twitter settings. See “Setting Your Preferences” on page 758 for more information.

View Menu

These View menu options are active:



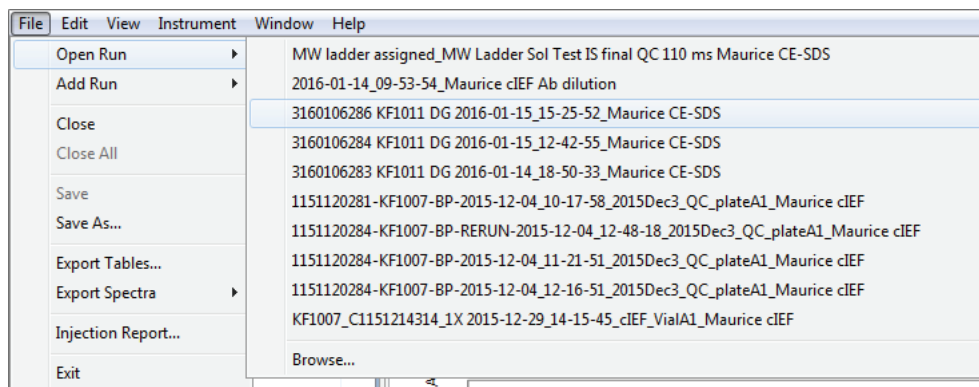
- **Selected** - Displays the data for only the injections selected.
- **All** - Displays data for all injections so you can scroll through them.
- **Standards** - Lets you view data just for the internal standards in your injections.
- **Samples** - Lets you view data just for sample proteins in your injections.
- **Grouping** - Displays data for injection groups.
- **View Region** - Lets you change the x-axis range of the data displayed.
- **Show Hidden** - Shows injections that are hidden from the data view.

Opening Run Files

You can open one run file or multiple files at the same time to compare information between runs.

Opening One Run File

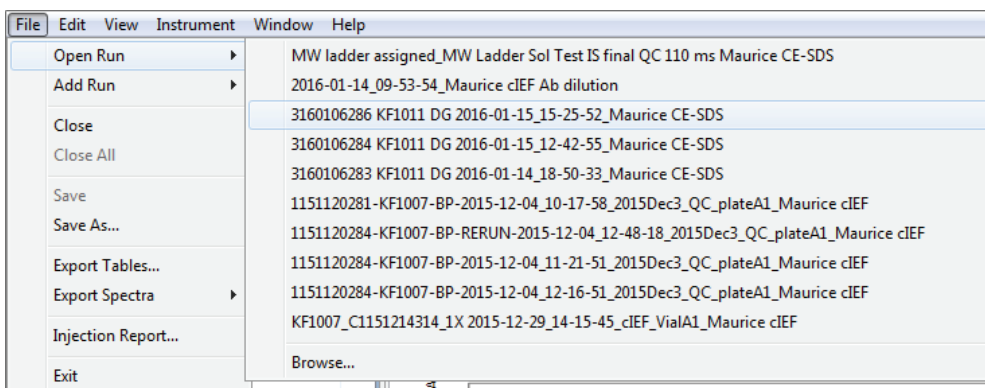
1. Select **File** in the main menu and click **Open Run**.



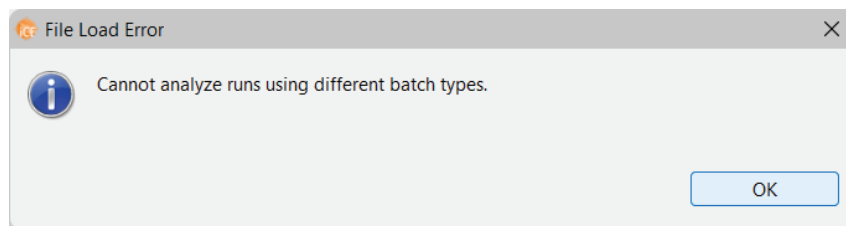
2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.

Opening Multiple Run Files

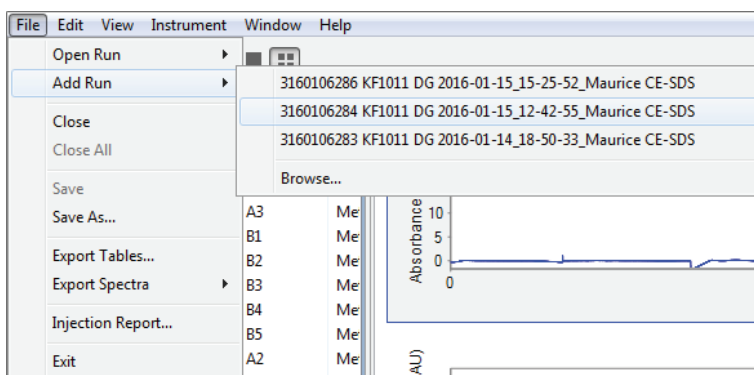
1. To open the first run file, select **File** in the main menu and click **Open Run**.



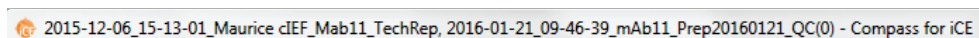
2. A list of the last 10 runs opened will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file. If the run selected is not a CE-SDS PLUS or Turbo CE-SDS batch, an alert will appear.



3. To open another run file, select **File** in the main menu and click **Add Run**.



4. A list of CE-SDS runs will display. Select one of these runs or click **Browse** to open the Runs folder and select a different file.
5. When a run is added, its data appends to the open run file and displays as a second set of injections in all screen panes. The second run file name also appears in the title bar:



6. Repeat the last two steps to add additional runs.

How Run Data is Displayed

Data in the run file is organized for easy review.

Experiment Pane: Batch Injection Information

The Experiment pane lists all the injections performed in the run, which samples were used for each, the sample location in the 96-well plate or 48-vial tray and the method used.

Injection	Injection Name	Sample	Location	Method
1	IgG System Contr...	IgG System Co...	A1	Method1
2	Control Ladder_02	Control Ladder	A2	Method2
3	Test Ladder_03	Test Ladder	A3	Method2
4	IS - Alpha_04	IS - Alpha	B1	Method1
5	IS - Frozen P3_05	IS - Frozen P3	B2	Method1
6	IS - T1 P3_06	IS - T1 P3	B3	Method1
7	IS - T2 P3_07	IS - T2 P3	B4	Method1
8	IS - T3 P3_08	IS - T3 P3	B5	Method1
9	Control Ladder_09	Control Ladder	A2	Method2

- **To view all columns** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.
- **To view method parameters** - Hover the mouse over a method name.

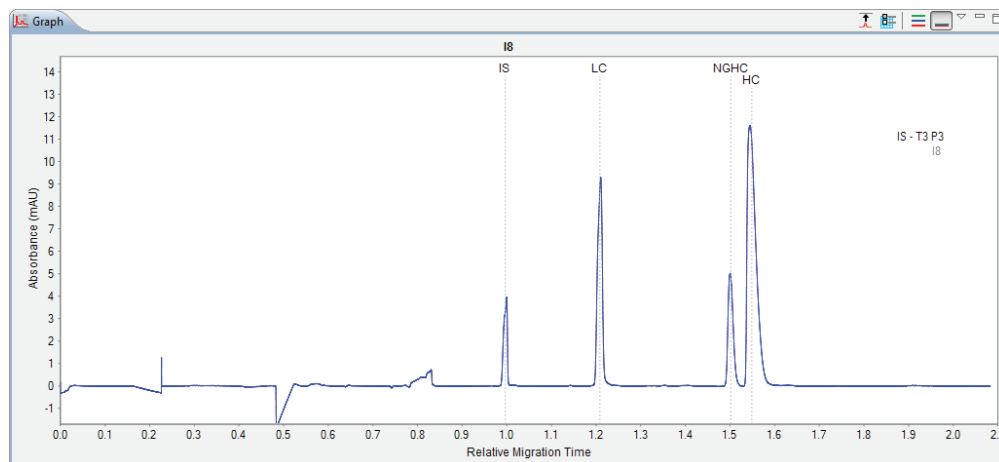
Injection	Injection Name	Sample	Location	Method
1	IgG System Contr...	IgG System Co...	A1	Method1
2	Control Ladder_02	Control Ladder	A2	Method2
3	Test Ladder_03	Test Ladder	A3	Method2
4	IS - Alpha_04	IS - Alpha	B1	Method1
5	IS - Frozen P3_05	IS - Frozen P3	B2	Method1

NOTE: Data notification icons will display in the Injection column if Compass for iCE detects a potential analysis issue or data was manually modified by the user. For more information see “Data Notifications and Warnings” on page 678.

Graph Pane: Electropherogram Data

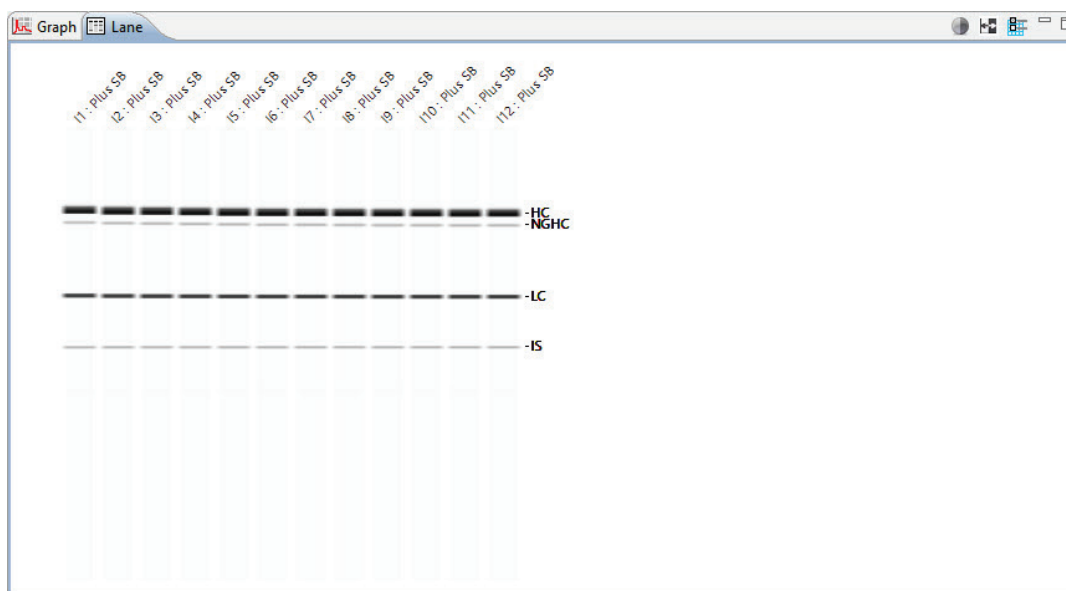
The Graph pane displays the electropherogram(s) for sample proteins or internal standards depending on the view options you’ve selected.

You can get more info on graph view options in “Changing the Electropherogram View” on page 694.



Lane Pane: Virtual Gel-Like Image Data

Click the **Lane** tab to view data for sample proteins as bands in individual lanes. This virtual gel-like image is similar to traditional gel results.



To view information for a band, roll the mouse over a band until the info box appears.

Lane data displayed in the virtual gel is automatically aligned by Compass for iCE. You can get more info on virtual gel view options, see “Changing the Virtual Gel View” on page 711.

Peaks Pane: Calculated Results

The Peaks pane shows the tabulated results for your sample proteins or internal standards. Each row in the table has the individual results for each peak detected in an injection. Results shown will either be for one injection or multiple injections, samples or standards depending on the view options you're using. Check out "Analysis Options Pane: Peak Names" on page 671 for more info.

Injection	Injection Name	Sample	Peak	Name	Time	RMT	MW (kDa)	Height	Raw Area	Area	% Total	% Area	Width	S/N
4	IS - Alpha_04	IS - Alpha	1	IS	729.0	1.000	10	3.5	217	298.4	100.0	100.0	7.2	250.3
4	IS - Alpha_04	IS - Alpha	2	LC	883.1	1.211	25	9.2	762	864.5	30.7	30.7	9.1	665.6
4	IS - Alpha_04	IS - Alpha	3	NGHC	1095.0	1.502	55	5.2	459	418.6	14.8	14.8	9.3	372.1
4	IS - Alpha_04	IS - Alpha	4	HC	1126.9	1.546	63	11.6	1739	1536.6	54.5	54.5	15.6	832.7
5	IS - Frozen P3_...	IS - Frozen P3	1	IS	727.3	1.000	10	3.7	230	317.8	100.0	100.0	7.2	233.1
5	IS - Frozen P3_...	IS - Frozen P3	2	LC	880.8	1.211	25	9.4	773	878.3	30.7	30.7	9.1	592.3
5	IS - Frozen P3_...	IS - Frozen P3	3	NGHC	1092.0	1.502	55	5.2	462	422.2	14.8	14.8	9.3	328.5

NOTES:

Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

When the Standards view is selected, the information in the Peaks table includes only injection, sample, peak, time and height. Internal standards the software has identified are marked with an **S**.

- **To view all rows** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.

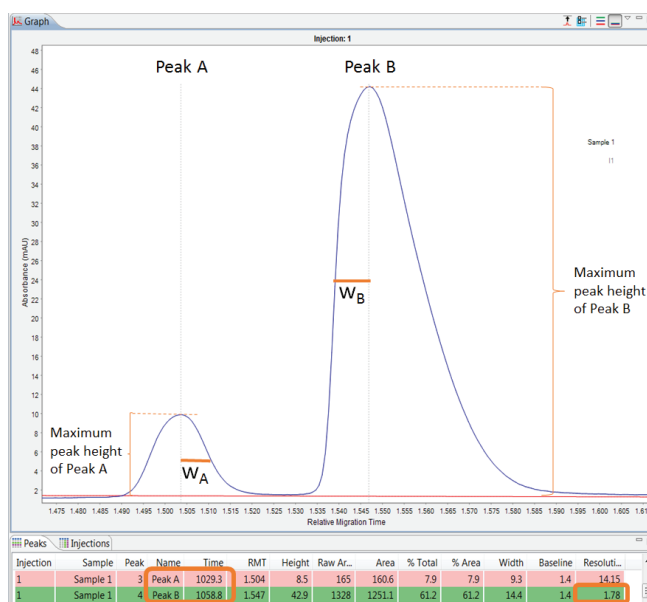
The following results and info are listed in the Peaks table:

- **Injection** - Injection number.
- **Injection name** - If injection names were entered in the batch, those names will display here. Otherwise the default 'SampleID_injection number' name will display.
- **Sample** - If sample names were entered in the batch, those names will display here. Otherwise, Sample (default name) will display.
- **Peak** - Peaks are numbered in order of detection.
- **Name** - Displays peaks Compass for iCE named automatically using the user-defined peak name analysis parameters. These cells are blank if the software wasn't able to name the peak or if you didn't enter naming parameters.
- **Time** - Peak detection time (seconds). This is the elapsed time between the start of the separation and when the peak is detected.
- **RMT** - Relative migration time of the peak to the Internal Standard which has an RMT of 1.0.
- **MW (kDa)** - Displays the relative molecular weight in kDa for sample peaks. MW only displays if you've run the CE-SDS MW Markers as one of the injections in the run and identified that injection in your analysis parameters.

- **Height** - The calculated peak height.
- **Raw Area** - Displays the uncorrected peak area.
- **Area** - Displays the time-corrected peak area. This includes corrections for big and/or slow moving peaks which can be artificially large when uncorrected.
- **% Total** - Displays the peak area ratio compared to the sum of all peak areas (excluding the Internal Standard peak). This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.
- **% Area** - Displays the calculated percent area for the named peak compared to all named peaks. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100 (shown for named peak sample data only).
- **Width** - Displays the calculated peak width (sample data only).
- **S/N** - Displays the calculated signal to noise. The signal to noise ratio is reported for each peak based on the noise in the injection and the height of the peak. $S/N = 2 \times \text{peak height} / \text{noise}$. For the injection noise calculation, the peak to peak noise is calculated by fitting a quadratic to the test region and taking the difference between the max signal above the fit, and the minimum signal below the fit. A fit is used to mitigate the effect of drift in the signal that would give a falsely high noise value. The software looks for the quietest 50 sec region in the first 10 to 600 sec of the injection. The test region is approximately 5 times a typical peak width (FWHM).
- **Baseline** - Displays the raw baseline signal of each peak.
- **Resolution** - Displays resolution of the peak compared to neighboring peaks. Two peaks that are baseline resolved will have a resolution value of 1.5. Smaller values mean the peaks are not baseline resolved. Resolution is calculated using this formula:

$$R = 1.18 * \frac{t_B - t_A}{(2w_B + 2w_A)}$$

- **R** - Resolution between Peak A and Peak B in Peaks Table.
- **w_A** - Right half peak width at half maximum peak height of Peak A in seconds.
- **w_B** - Left half peak width at half maximum peak height of Peak B in seconds.
- **t_A** - Migration time of Peak A in seconds in Peaks Table.
- **t_B** - Migration time of Peak B in seconds in Peaks Table.



Injections Pane: User-Specified Peak Names

The Injections pane shows tabulated results for sample proteins Compass for iCE labels automatically using user-defined peak name settings. Each row in the table shows the individual results for the named peaks detected in each injection.

Injections										
Injection	Injection Name	Sample	IS	HC	NGHC	LC	10kDa	20kDa	33kDa	55
1	IgG System Co...	IgG System Co...	303	1376	381	789				
3	Test Ladder_03	Test Ladder							375	
4	IS - Alpha_04	IS - Alpha	298	1537	419	865				
5	IS - Frozen P3_...	IS - Frozen P3	318	1560	422	878				
6	IS - T1 P3_06	IS - T1 P3	329	1551	415	863				

NOTES:

Peaks that Compass for iCE names automatically with user-defined peak name settings are color-coded.

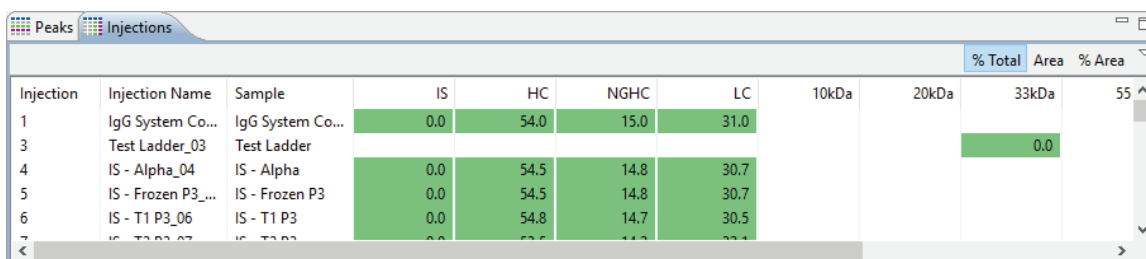
When the Standards view is selected, the information in the Injections table includes only injection, sample and std 1 (the migration time of the standard peak).

- **To view all rows** - Use the scroll bar or click **Maximize** in the upper right corner.
- **To resize columns** - Roll the mouse over a column border until the sizing arrow appears, then click and drag to resize.

The following results and info are listed in the Injections table:

- **Injection** - Injection number.
- **Injection name** - If injection names were entered in the batch, those names will display here. Otherwise the default 'SampleID_injection number' name will display.
- **Sample** - If sample names were entered in the batch, those names will display here. Otherwise, Sample (default name) will display.
- **Peak Name Columns** - An individual column per peak name will display for every peak identified by name or as a MW Marker peak in the run data. Cells for injections in these columns will be blank if Compass for iCE didn't find peaks automatically using the user-defined peak name analysis and maker parameters (or none were entered).
 - **To view peak area in the peak name columns (default)** - Select **Area** in the upper right corner of the pane. This displays calculated peak area for the individual peak only.
 - **To view % total in the peak name columns** - This displays the peak area ratio compared to the sum of all peak areas (excluding the Internal Standard peak). This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.

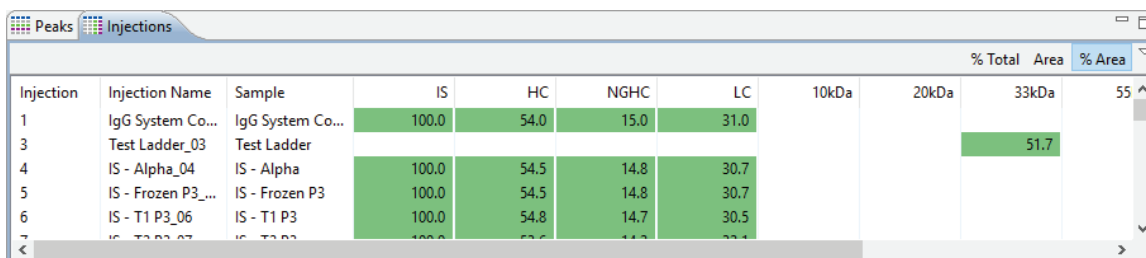
NOTE: The sum of the named peak percentages can be less than 100% if some peaks aren't named.



The screenshot shows the 'Injections' tab in a software interface. The table displays data for several injections. The 'Area' column is selected, showing the calculated peak area for each peak. The 'IS' (Internal Standard) column is highlighted in green for all rows, indicating it is the reference peak.

Injection	Injection Name	Sample	IS	HC	NGHC	LC	10kDa	20kDa	33kDa	55
1	IgG System Co...	IgG System Co...	0.0	54.0	15.0	31.0				
3	Test Ladder_03	Test Ladder							0.0	
4	IS - Alpha_04	IS - Alpha	0.0	54.5	14.8	30.7				
5	IS - Frozen P3_...	IS - Frozen P3	0.0	54.5	14.8	30.7				
6	IS - T1 P3_06	IS - T1 P3	0.0	54.8	14.7	30.5				

- **To view % area in the peak name columns** - This displays the peak area ratio compared to the sum of all named peak areas. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100.

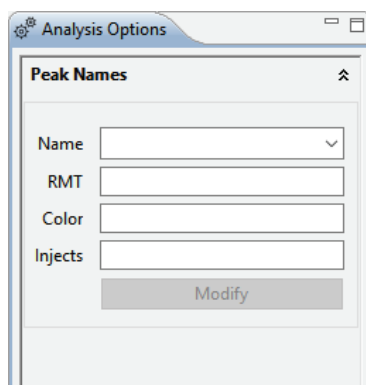


The screenshot shows the 'Injections' tab in a software interface. The table displays data for several injections. The '% Area' column is selected, showing the percentage of the total area for each peak. The 'IS' (Internal Standard) column is highlighted in green for all rows, indicating it is the reference peak.

Injection	Injection Name	Sample	IS	HC	NGHC	LC	10kDa	20kDa	33kDa	55
1	IgG System Co...	IgG System Co...	100.0	54.0	15.0	31.0				
3	Test Ladder_03	Test Ladder							51.7	
4	IS - Alpha_04	IS - Alpha	100.0	54.5	14.8	30.7				
5	IS - Frozen P3_...	IS - Frozen P3	100.0	54.5	14.8	30.7				
6	IS - T1 P3_06	IS - T1 P3	100.0	54.8	14.7	30.5				

Analysis Options Pane: Peak Names

The Analysis Options pane gives you a quick way to view, change and add new custom peak name settings for sample proteins without having to open and edit the run's analysis settings. You can apply peak name settings to all, selected, or specified injections. For details on how to use this option, go to “Naming Peaks” on page 679.



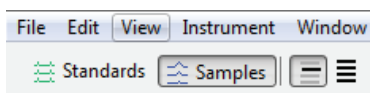
Viewing Run Data

The Analysis screen lets you view data for just one injection, specific injections or all injections in the run. Each run file has data for the sample peaks and the Internal Standard detected in each injection.

Switching Between Samples and Standards Data Views

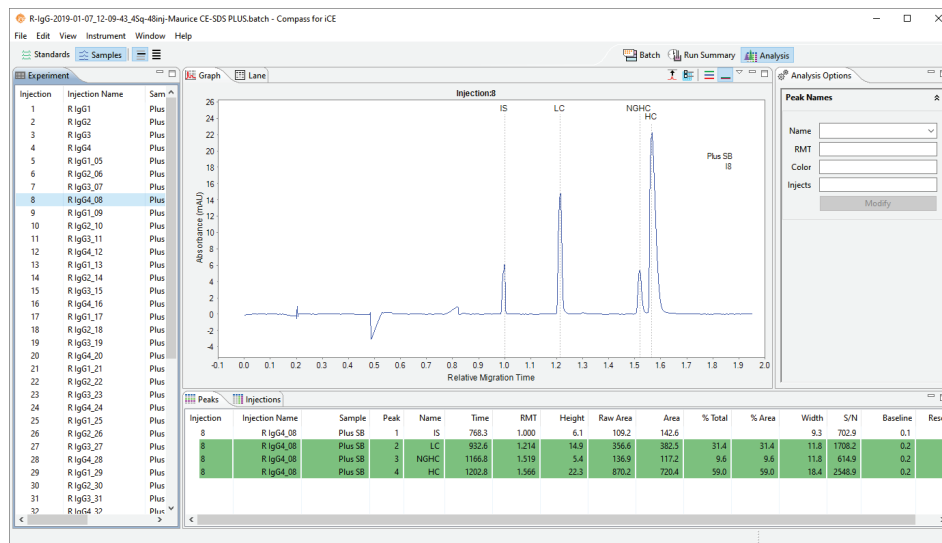
Here's how you switch between viewing data for your samples and standards:

- To view sample data - Click **Samples** in the View bar or select **View** in the main menu and click **Samples**.



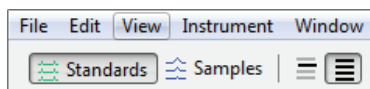
- Data in this view is for sample proteins only.
- The graph displays electropherograms with a y-axis of Absorbance units (mAU) and an x-axis of RMT (relative migration time).
- Lane view data displays sample proteins only.

- Results for each protein are shown in the Peaks and Injections panes.

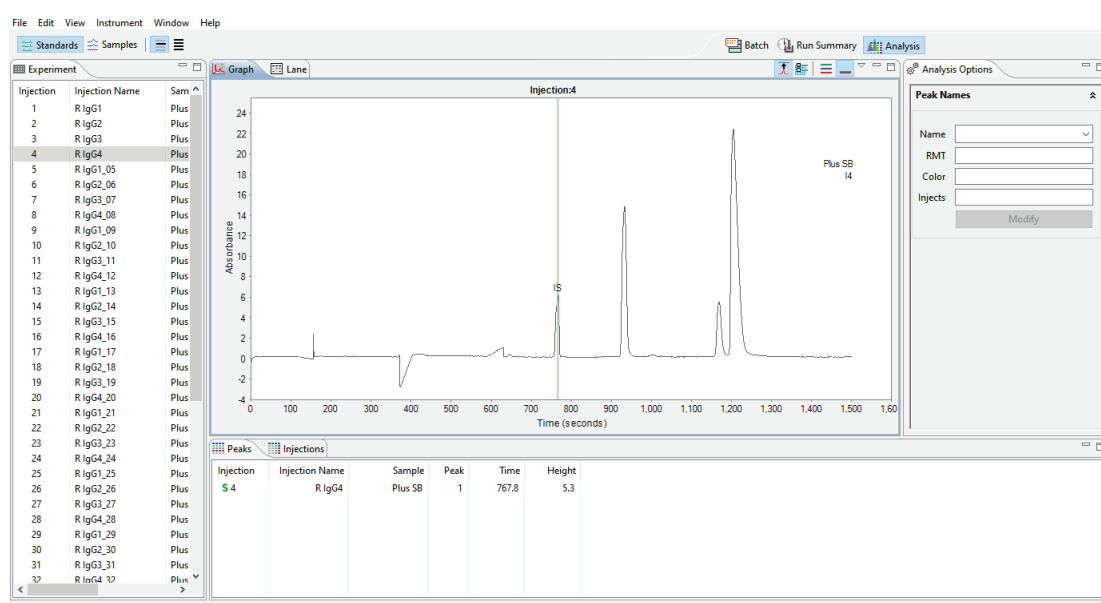


For information on checking and identifying sample peaks, see “Checking Your Data” on page 319 for CE-SDS PLUS runs and page 388 for Turbo CE-SDS runs.

- To view Internal Standard data** - Click **Standards** in the View bar or select **View** in the main menu and click **Standards**.



- Data in this view is for analyzing standards only. This is the Internal Standard you add to your samples during prep.
- The graph displays electropherograms with a y-axis of Absorbance units (mAU) and an x-axis of time in seconds.
- Lane view data displays standards only.
- The Internal Standard is identified in the Peaks pane with an **S** and as IS in the Injections pane.

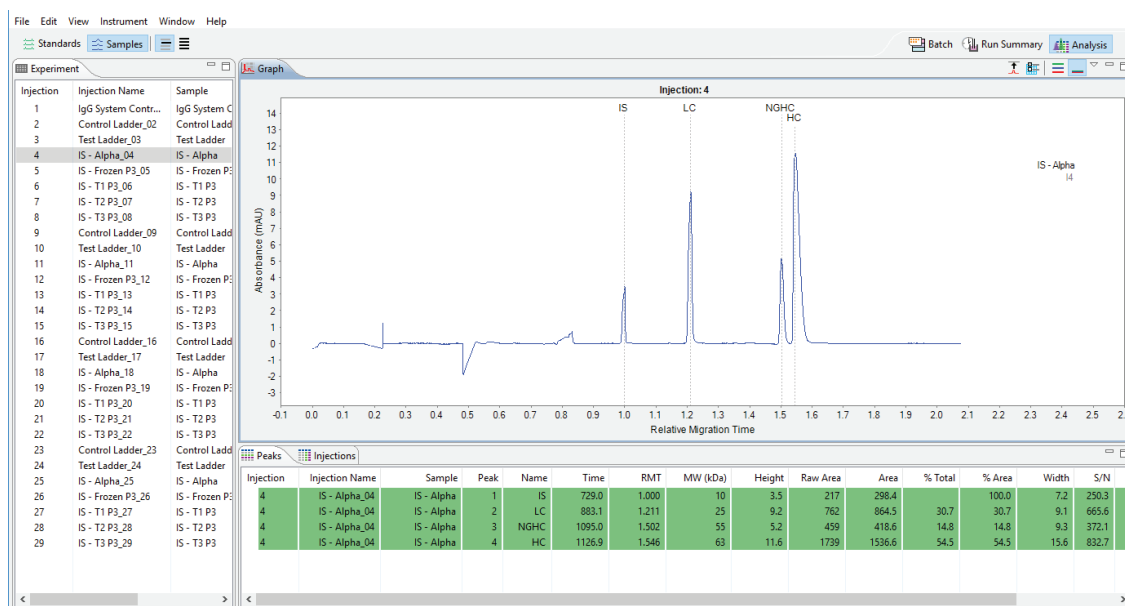


For information on checking and identifying the Internal Standard peak, see “Checking Your Data” on page 319 for CE-SDS PLUS runs and page 388 for Turbo CE-SDS runs.

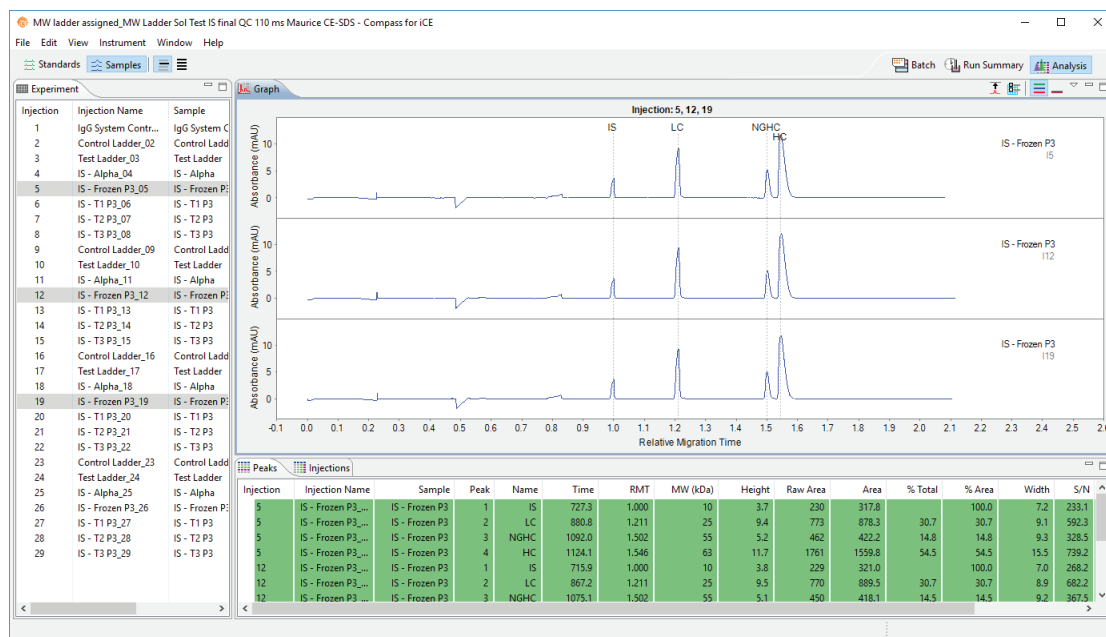
Selecting and Displaying Injection Data

You can view data from one, multiple, or all injections at once.

- **To look at data for one injection** - Click an injection row in the Experiment pane. Data for just that injection displays in the data views and tables.



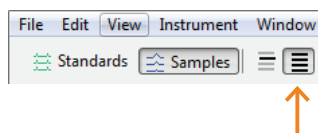
- **To look at data for specific injections** - Hold the **Ctrl** key and select just the injection rows you want to view in the Experiment pane. Data for only the injections selected display in the data views and tables.



- **To look at data for sequential injections** - Select the first injection row in the Experiment pane that you want to view, then hold the **Shift** key and select the last. This selects all rows between the two injections. Data for only the injections selected display in the data views and tables.



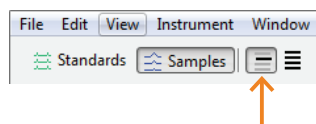
- To look at data for all injections - Just click **View All** in the View bar. Data for all injections displays in the graph and tables.



Switching Between Single and Multiple Views of Injections

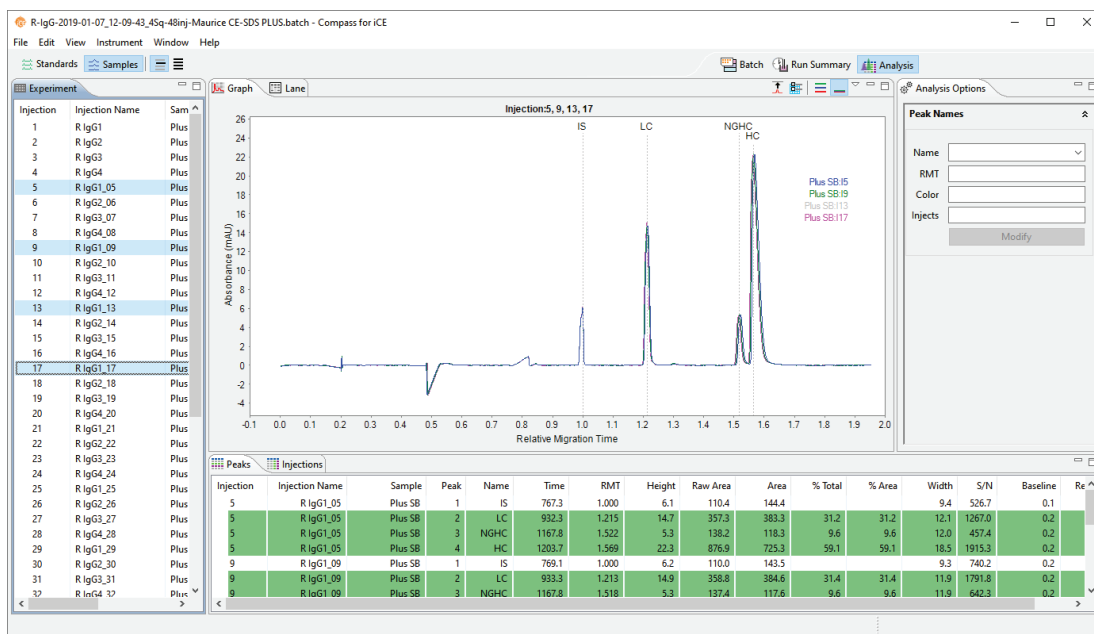
You can switch between displaying run data in a single, per-injection format or a multi-injection format.

- To view data in a per-injection format - Click **View Selected** in the View bar or select **View** in the main menu and click **Selected**.

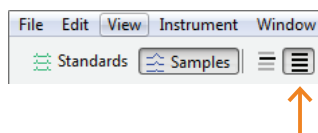


Data for the injection row(s) selected in the Experiment pane:

- Displays with electropherograms either overlaid or stacked in the Graph pane depending on the option you've got chosen.
- Lanes for only the selected row(s) are displayed in the lane pane.
- Shows only results for the selected row(s) in the Peaks and Injections panes.

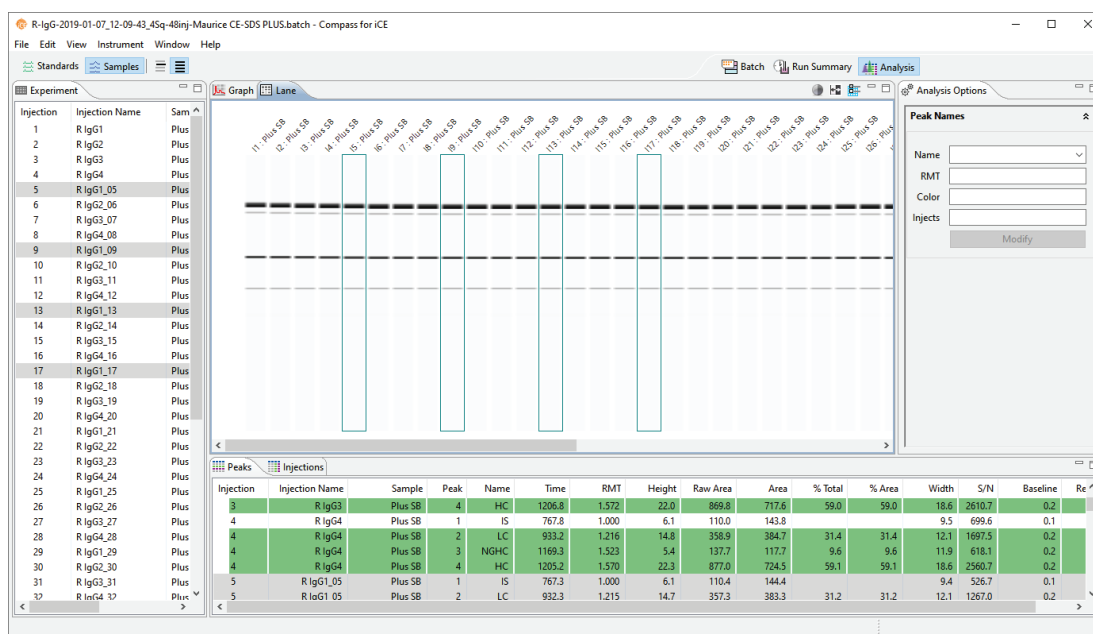


- To view data in a multi-injection format - Click **View All** in the View bar or select **View** in the main menu and click **View All**:



Data for the injection row(s) selected in the Experiment pane:

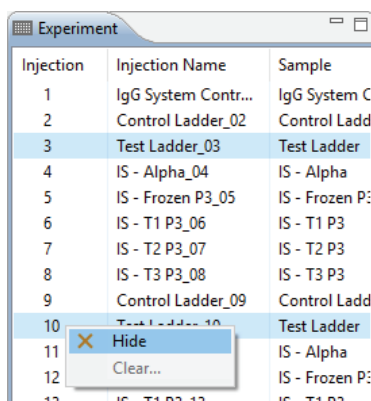
- Displays with the electropherograms of the selected injections highlighted in the Graph pane.
- Displays all lanes in the lane pane, and lanes corresponding to the selected row(s) are highlighted.
- Shows the results for the selected injections highlighted in the Peaks and Injections panes.



Hiding Injection Data

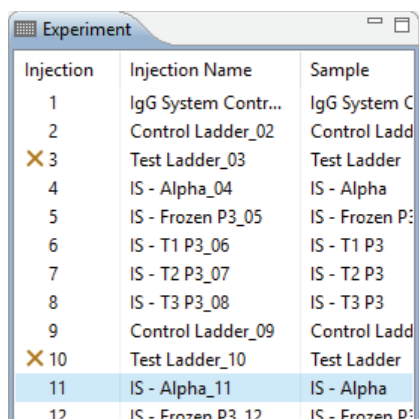
You can hide injection data from the view if needed.

- **To hide injections** - Select the injection rows you want to hide in the Experiment pane, then right click one and select **Hide**.



Data for the injections will be hidden in all data views and results tables.

- **To view hidden injections** - Select **View** in the main menu and click **Show Hidden**. Hidden rows will become visible again in all panes, and are marked with an X in the Experiment pane.



- **To unhide injections** - Select the hidden row(s). Right click on one and click **Unhide**.

Data Notifications and Warnings

If Compass for iCE detects a potential data issue, a notification or warning icon will display next to the injection row in the Experiment pane.



Manual correction of sample data notification - This means the sample data was manually changed by a user, for example to add or remove a sample peak. Roll your mouse over the icon to display the type of modification that was made.

3	Test Ladder	A3	Me
✓ 4	IS - Alpha	B1	Me
5	IS - Frozen P3	B2	Me
6	Peak Fit Manual	B3	Me



Standards warning - This means the Internal Standard may not be identified properly. You can fix this by manually identifying the standard using the steps in “Step 1: Check Your Internal Standard” on page 319 for a CE-SDS PLUS run or page 388 for a Turbo CE-SDS run. Roll your mouse over the icon to display warning details.

11	IgG2B-R	B1	Method1
12	hSAP IgG-R	C1	Method1
Ⓢ 13	IgG2B-NR	D1	Method2
Ⓢ 14	hSAP-IgG-NR	E1	Method2
15	Standards Warning: Low Confidence		Method1



Manual correction of standards data notification - This means a user changed the standards data manually. Roll your mouse over the icon to display the type of modification that was made.

8	IS - T3 P3	B5	Me
9	Control Ladder	A2	Me
✓ 10	Test Ladder	A3	Me
11	IS - Alpha	B1	Me
12	IS - Frozen P3	Standards Manual	Me
13	IS - T1 P3	B3	Me



Peak fit warning - Means that a peak can't be fit properly. This can sometimes be caused when a broad peak is fitted as multiple narrow peaks. Changing the peak width can help in this case. The warning is also caused by very small peaks around main peaks, or small peaks that are close to the end of the separation range. You can often fix this by removing the peak(s) using the steps in “Step 3: Checking Sample Peaks” on page 325 for CE-SDS PLUS runs or page 394 for Turbo CE-SDS runs. Roll your mouse over the icon to display warning details.

6	mAb 25	A3
7	mAb 250	A2
Peak Fit Warning: Too many iterations		
9	mAb 250	A2
10	mAb 25	A3

Checking Your Results

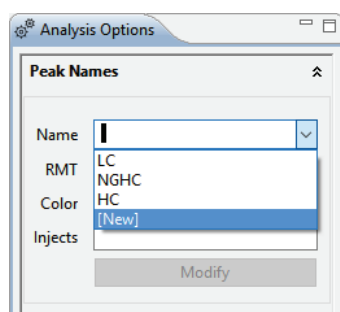
Compass for iCE detects your sample protein, CE-SDS MW Markers and Internal Standard peaks and reports results automatically. But, we always recommend you review and check your data as a good general practice to make sure your results are accurate. Please see the step by step procedure in “Checking Your Data” on page 319 for CE-SDS PLUS runs or page 388 for Turbo CE-SDS runs to do this. If you see a data warning in the Experiment pane, these steps will also help you identify and correct any issues.

Naming Peaks

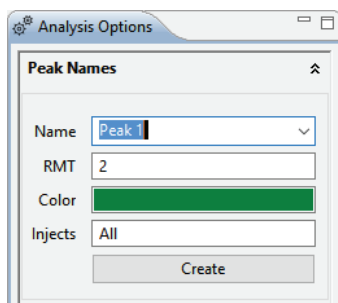
NOTE: Analysis screen options will let you add a new peak name, its RMT, identifying color for peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set RMT ranges, you can do that in the run analysis settings. For more info go to “Peak Names Settings” on page 723.

Adding New Peak Names in the Analysis Options Pane

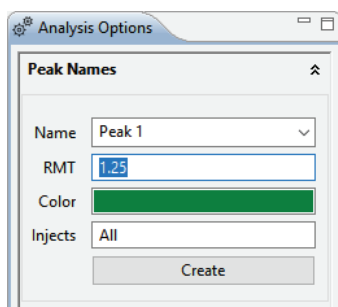
1. Click the down arrow in the Name field and select New.



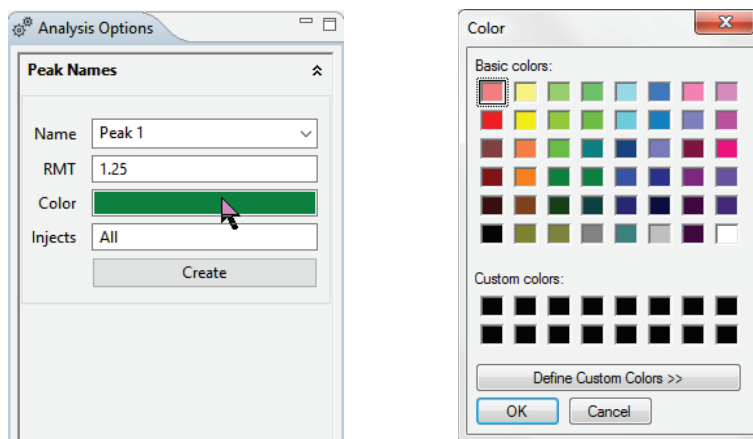
2. Type a name.



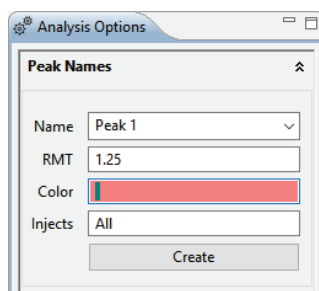
3. Click in the RMT field and enter the relative migration time of your sample protein.



4. Click on the Color field to display the color selection box.



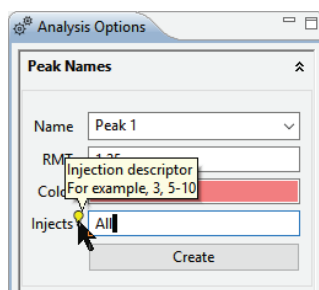
5. The color you pick is used to identify the sample protein peak in the Peaks and Injections panes in the Analysis screen. Click a color or define a custom color and click OK. The color selection will update in the field:



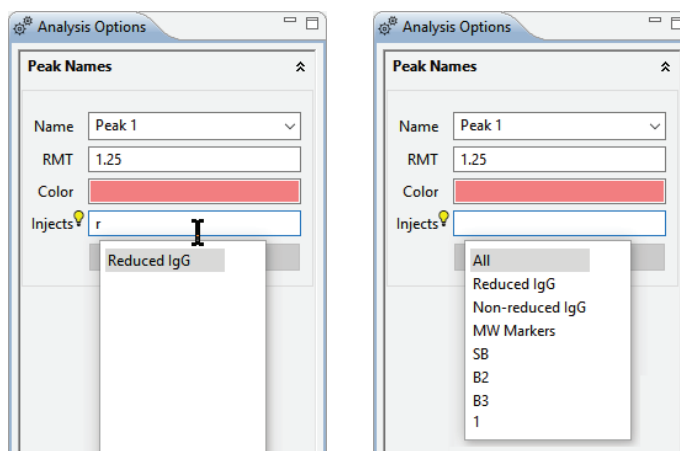
6. Click in the Injects field:

To apply the new peak name to all injections - leave the default setting of All.

To apply the new peak name to specific injections - enter individual injections separated by commas (for example 2, 5, 8), a range of injections (for example 3-15) or a combination of both (for example 8, 10, 12-15). You can also click the tool tip for more examples.



To apply the new peak name to specific run data - type the first letter of a method, sample name, well or vial location. Compass for iCE will search the run data and display a list of matching options for you to choose from. Alternatively you can highlight the text in the Injects cell, hit **Delete**, then select an option from the drop down list.

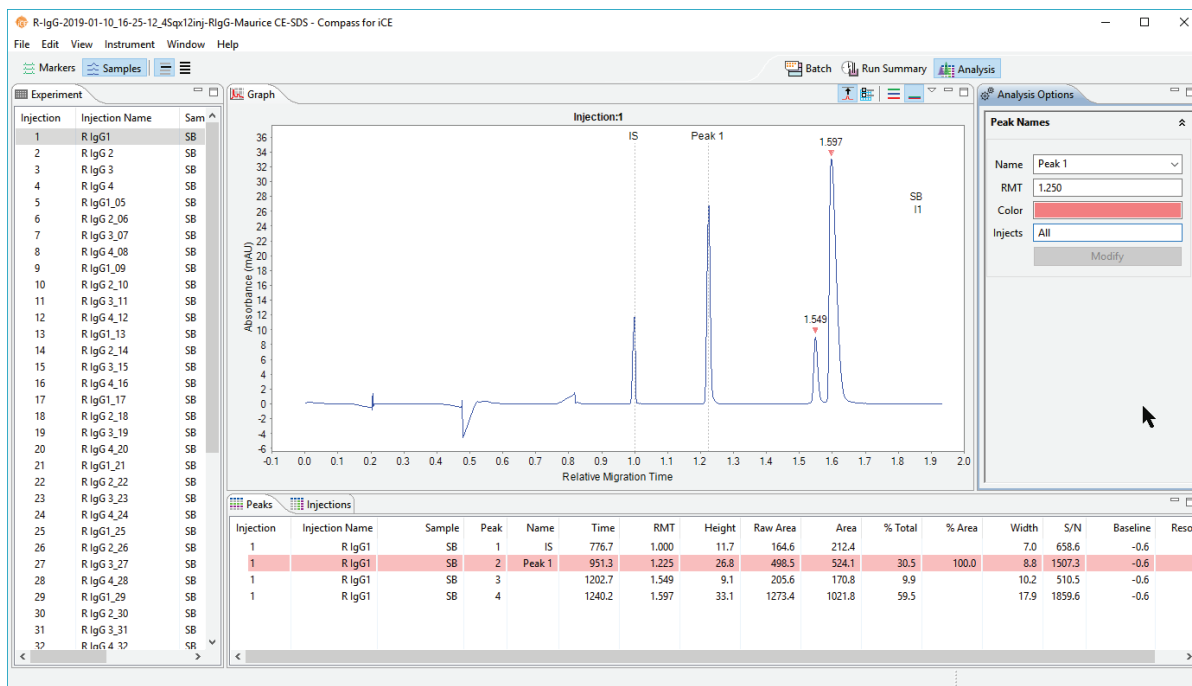


- **All** - Applies the peak name to all injections.
- **Methods** - All methods in the run file display in the list. Selecting a method applies the peak name to all injections that used that method.
- **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the peak name to all injections that used that sample name.
- **Wells or vials** - The well or vial numbers currently selected will display in the list. Selecting a well or vial number applies the peak name to all injections that used that well or vial.

7. Click **Create** to save changes.

NOTE: To clear the new peak name information you've entered without saving changes, just delete the peak name in the Name field.

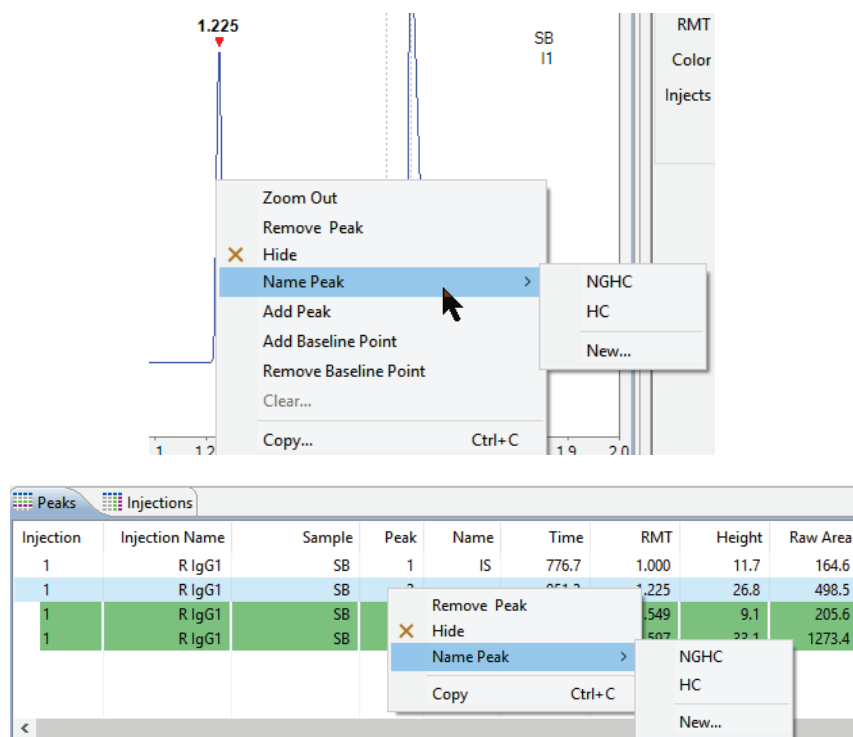
The named peak will be identified with a peak name label in the electropherogram and lane view, and color-coded in the Peaks and Injections panes:



Adding Peak Names from the Graph or Peaks Table

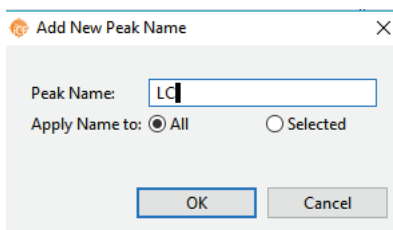
NOTE: Analysis screen options will let you change an existing peak name, its RMT, identifying color in peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set RMT ranges, you can do that in the run analysis settings. For more info go to “Peak Names Settings” on page 723.

1. Right click the peak you want to name in the Graph or Peaks pane.
2. Select Name Peak.



3. To use an existing peak name - select a name from the list.

To create a new peak name - select New. Type in a name for the peak. Click All to apply to all injections or Selected to apply only to the injections selected.

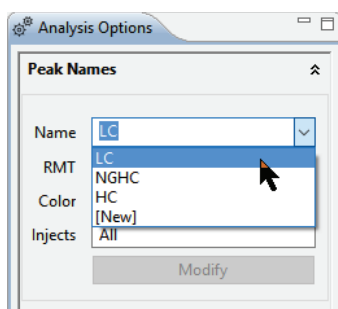


4. Click **OK**. If you want to add an identifying color or change the injections the new peak name applies to, you can do that in the Analysis Options pane.

Changing Peak Name Information

NOTE: Analysis screen options will let you change an existing peak name, its RMT, identifying color for peak tables and what injections it applies to. If you need to remove peak names, add peak name analysis settings groups or set RMT ranges, you can do that in the run analysis settings. For more info go to “Peak Names Settings” on page 723.

1. In the Analysis Options pane, click the down arrow in the **Name** field and select an existing peak name.



2. Change the name, RMT, color and injects as needed then click **Modify**.

Group Statistics

You can use the Grouping view to have Compass for iCE do a statistical analysis of named proteins in your injections (see “Peak Names Settings” on page 723 for more info on setting named peaks up). Statistics for each protein are also plotted for easy comparison.

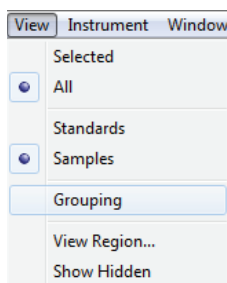
Using Groups

1. Groups are automatically created for injections that use the same sample name and method, so to use this feature, you need to make sure you’ve got sample names entered.
 - a. Go to the **Batch** screen.

- b. Click the **Sample ID** cells in the Injection pane and type a name for any samples you want to calculate statistics for.

Injections						History	Notes	Pause	Stop	Add	Replicate	Remove			
	Injection Name	Sample ID	Location	Method	Notes										
1	Sample 1_01	Product A	A1	Reduced IgG											
2	Sample 02_02	Sample 02	A2	Reduced IgG											
3	Sample 03_03	Sample 03	A3	Reduced IgG											
4	Sample 04_04	Sample 04	A4	Reduced IgG											
5	Sample 05_05	Sample 05	A5	Reduced IgG											
6	Sample 06_06	Sample 06	A6	Reduced IgG											
7	Sample 07_07	Sample 07	A7	Reduced IgG											
8	Sample 08_08	Sample 08	A8	Reduced IgG											
9	Sample 09_09	Sample 09	A9	Reduced IgG											
10	Sample 10_10	Sample 10	A10	Reduced IgG											
11	Sample 11_11	Sample 11	A11	Reduced IgG											
12	Sample 12_12	Sample 12	A12	Reduced IgG											

2. Go back to the **Analysis** screen. Click **View** in the main menu and select **Grouping**.



NOTE: To turn Grouping off, select **View** in the main menu and deselect **Grouping**.

Viewing Sample Injection Groups

Compass for iCE automatically groups all injections using the same sample name together in the Injection Groups pane.

Injection Groups				
Sample		Method		
▶ Control Ladder (3)		Method2		
▶ IS - Alpha (4)		Method1		
▶ IS - Frozen P3 (4)		Method1		
▶ IS - T1 P3 (4)		Method1		
▶ IS - T2 P3 (4)		Method1		
▶ IS - T3 P3 (4)		Method1		
IgG System Control		Method1		
▶ Test Ladder (4)		Method2		

- **To expand a group** - Click the arrow next to a group to see the individual injections in the group and reported data for each

Sample	Method
▶ Control Ladder (3)	Method2
▶ IS - Alpha (4)	Method1
▶ IS - Frozen P3 (4)	Method1
▶ IS - T1 P3 (4)	Method1
▲ IS - T2 P3 (4)	Method1
IS - T2 P3	Method1
IS - T2 P3	Method1
IS - T2 P3	Method1
IS - T2 P3	Method1
▶ IS - T3 P3 (4)	Method1
IgG System Control	Method1
▶ Test Ladder (4)	Method2

- **To expand all groups** - Click Expand All (+) in the upper right corner of the pane.
- **To collapse all groups** - Click Collapse All (-) in the upper right corner of the pane.

Viewing Statistics

Peak and Method Groups

The Peak Groups pane reports statistics for each named protein in every group. Each group shows the statistics for named proteins which includes average area, standard deviation, %CV and SEM (standard error measurement). The number in parenthesis after the sample name is the number of injections in the group.

Sample	Method	Injection	Name	Area	Std.Dev.	% CV	SEM
> SB (48)	Reduced IgG		LC	450	54.73	12.2	7.899
> SB (48)	Reduced IgG		NGHC	181	24.26	13.4	3.502
> SB (48)	Reduced IgG		HC	1134	143.7	12.7	20.74

- **To display results using area** - Click Area in the upper right corner of the pane.
- **To display results using % total** - Click % Total in the upper right corner of the pane to display the calculated percent area for the named peak compared to the total area measured in the injection. This value results from dividing the individual peak area by the sum of all peak areas for the injection and multiplying by 100.
- **To display results using % area** - Click % Area in the upper right corner of the pane to display the calculated percent area for the named peak compared to all named peaks. This value results from dividing the individual peak area by the sum of all named peak areas for the injection and multiplying by 100 (shown for named peak sample data only).

- **To expand a group** - Click the arrow next to a group to see the individual injections in the group and reported data for each

Sample	Method	Injection	Name	Area	Std.Dev.	% CV	SEM
SB (48)	Reduced IgG		LC	450	54.73	12.2	7.899
SB	Reduced IgG	1	LC	499			
SB	Reduced IgG	2	LC	584			
SB	Reduced IgG	3	LC	495			
SB	Reduced IgG	4	LC	487			
SB	Reduced IgG	5	LC	488			
SB	Reduced IgG	6	LC	568			
SB	Reduced IgG	7	LC	472			

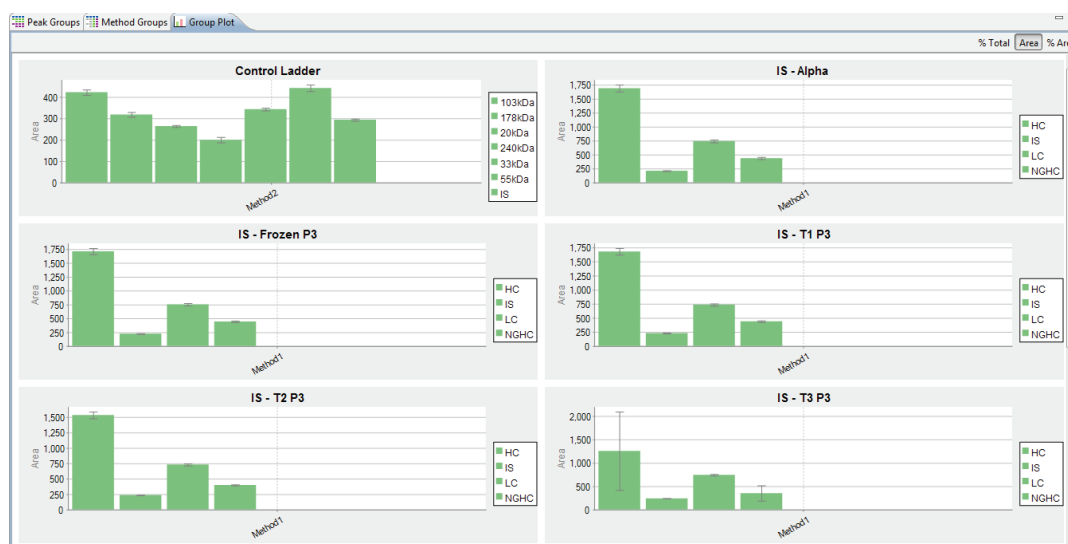
- **To expand all groups** - Click Expand All (+) in the upper right corner of the pane.
- **To collapse all groups** - Click Collapse All (-) in the upper right corner of the pane.

The Method Groups pane pivots the Peak Groups pane results to show statistics for named protein peaks in individual columns.

Sample	Method	Injection	IntSt:Area	Std.Dev.	%CV	SEM	BSA:Area	Std.Dev.	%CV	SEM
> 0 (3)	Reduced IgG		0	0.0000	0.0	0.0000	0	0.05963	87.7	0.03443
> 0 PLUS (3)	IS		0	0.06329	21.6	0.03654	0	0.0000	0.0	0.0000
> 0.13 PLUS (3)	IS		1	0.2110	41.7	0.1218	0	0.0000	0.0	0.0000
> 0.41 (3)	Reduced IgG		0	0.0000	0.0	0.0000	1	0.3290	51.1	0.1899
> 0.41 PLUS (3)	IS		1	0.1279	17.2	0.07382	0	0.0000	0.0	0.0000
> 1.23 (3)	Reduced IgG		0	0.0000	0.0	0.0000	1	0.2445	17.3	0.1412
> 1.23 PLUS (3)	IS		2	0.09553	3.9	0.05515	0	0.0000	0.0	0.0000
> 100 PLUS (3)	IS		148	4.119	2.8	2.378	0	0.0000	0.0	0.0000

Group Plots

The mean values for named peaks using the same method in each injection group are plotted in bar graphs with error bars showing the standard deviation in the Group Plots pane. You'll also get plots that compare samples using the same method in the run.



Hiding or Removing Injections in Group Analysis

Hidden injections are not included in injection groups. But, hiding injections gives you an easy way to reject individual injections from the statistical analysis. See Hiding Injection Data for details on how to do this.

Copying Data Views and Results Tables

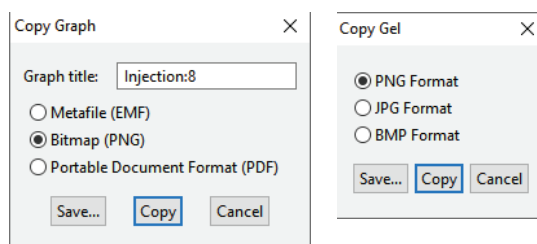
You can copy and paste data and results tables into other documents, or save a data view as a graphic file.

Copying Results Tables

1. Click in the Peaks or Injections pane.
2. Select one or multiple rows.
3. Select **Edit** in the main menu and click **Copy**, or right click on row(s) you selected and click **Copy**.
4. Open a document (Microsoft® Word®, Excel®, PowerPoint®, etc.). Right click in the document and select **Paste**. Data for the rows selected will be pasted into the document.

Copying Data Views

1. Click in the Graph or Lane pane.
2. Select **Edit** in the main menu and click **Copy**, or right click and select **Copy**.
3. Select an image option in the pop-up window, then click **Copy**.

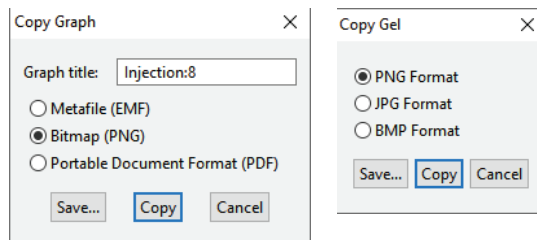


4. Open a document (Microsoft® Word®, Excel®, PowerPoint®, etc.). Right click in the document and select **Paste**. A graphic of the copied electropherogram will be pasted into the document.

Saving the Data Views as an Image File

1. Click in the Graph or Lane pane.
2. Select **Edit** in the main menu and click **Copy**, or right click and select **Copy**.

3. Select an image option in the pop-up window, then click **Save**.



4. Select a directory to save the file to, enter a file name, then click **OK**.

Exporting Run Files

Results tables and raw plot data can be exported for use in other applications.

Exporting Results Tables

To export the information in the Peaks and Injections tables:

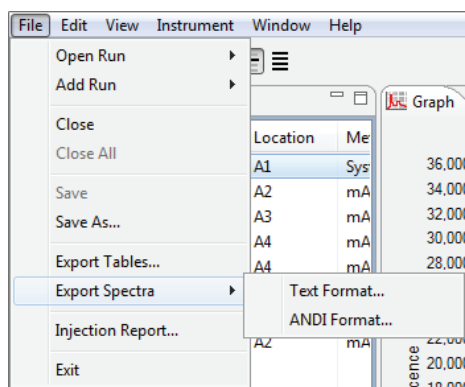
1. Click **File** in the main menu and click **Export Tables**.
2. Select a directory to save the files to and click **OK**. Data will be exported in .txt format.

NOTE: To exclude export of standards data or export results table data in .csv format, see “Setting Data Export Options” on page 760.

Exporting Raw Sample Electropherogram Data

To export raw sample plot and background data:

1. Click **File** in the main menu and click **Export Spectra**.



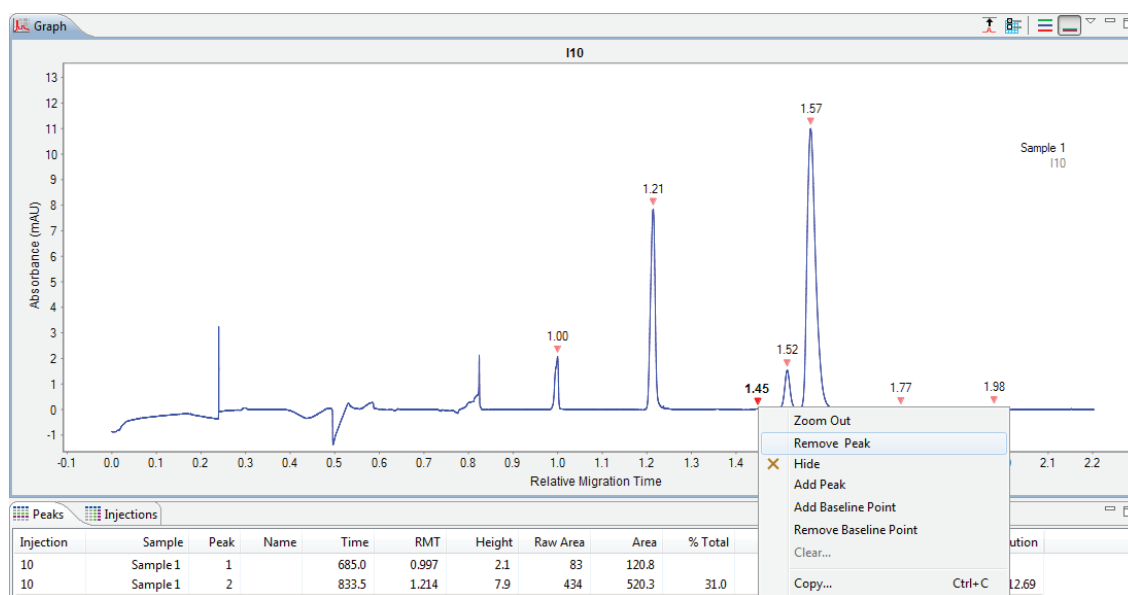
- **To export data in .txt format** - Select **Text Format**. Data will be exported in one file for all injections.
 - **To export data in .cdf format** - Select **ANDI Format**. Data will be exported in one file per injection.
2. Select a directory to save the files to and click **OK**. Data will be exported in the selected format.

Changing Sample Protein Identification

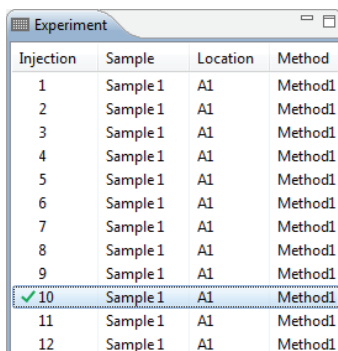
Compass for iCE lets you customize what sample proteins are reported in the results tables by making manual adjustments in the electropherogram or Peaks table.

Adding or Removing Sample Data

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that has the injection you want to correct, then click the **Graph** tab.
 - **To remove a peak from the data** - Right click the peak in the electropherogram or Peaks table and select **Remove peak**. The software will no longer identify it as a sample peak in the electropherogram, and the peak data will be removed in the results tables.



A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.



Injection	Sample	Location	Method
1	Sample 1	A1	Method1
2	Sample 1	A1	Method1
3	Sample 1	A1	Method1
4	Sample 1	A1	Method1
5	Sample 1	A1	Method1
6	Sample 1	A1	Method1
7	Sample 1	A1	Method1
8	Sample 1	A1	Method1
9	Sample 1	A1	Method1
✓ 10	Sample 1	A1	Method1
11	Sample 1	A1	Method1
12	Sample 1	A1	Method1

- **To add an unidentified peak to the data** - Right click the peak in the electropherogram or peaks table and select **Add Peak**. The software will calculate and display the results for the peak in the results tables and identify the peak in the electropherogram.

A check mark will appear next to the injection in the Experiment pane to indicate a manual correction was made.

NOTES:

To remove sample peak assignments that were made manually and go back to the original peak data, right-click the peak in the electropherogram and select **Clear** for the current injection or **Clear All** for all injections in the batch.

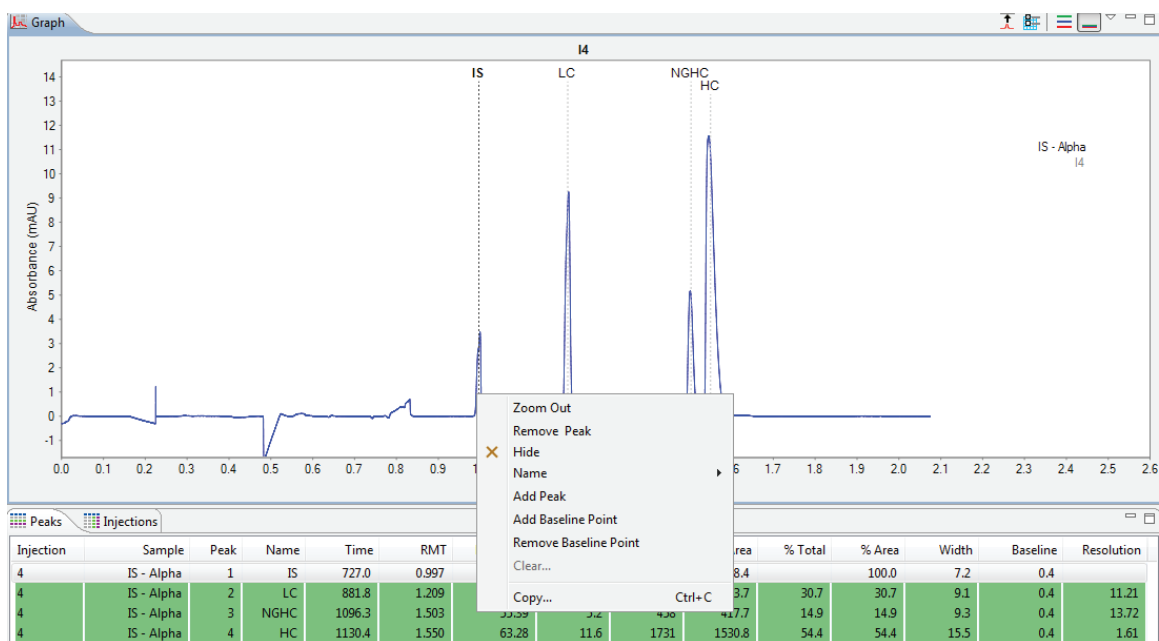
Virtual gel data in the lane pane will also update to reflect changes made in the graph pane.

Hiding Sample Data

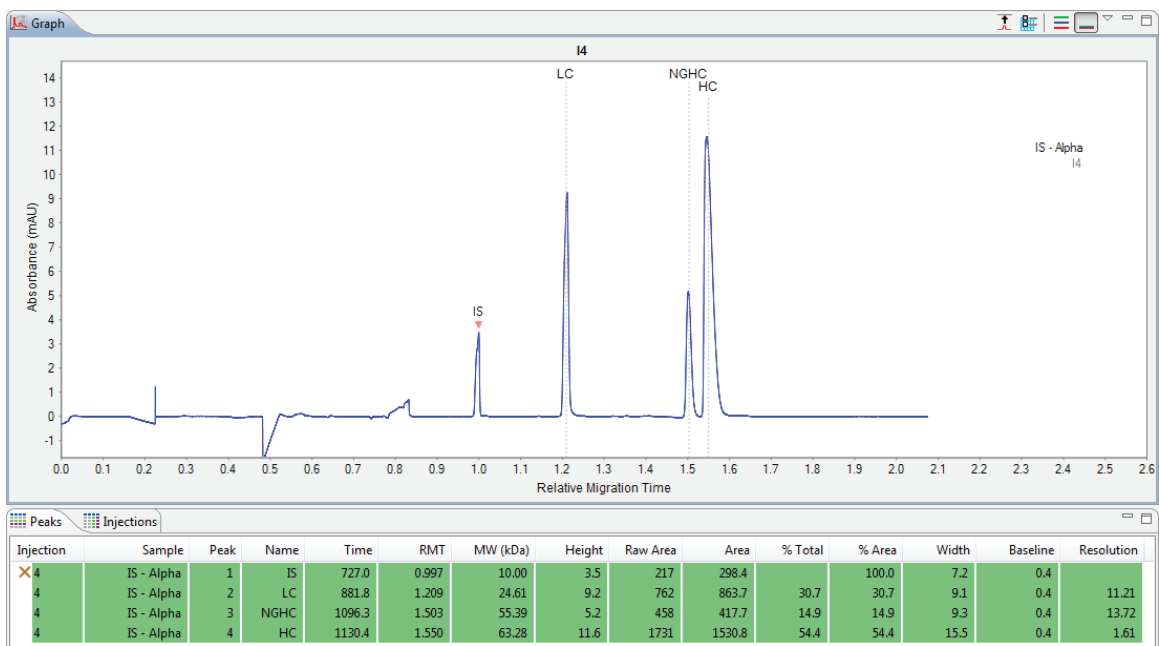
You can hide the results for a sample protein in the results tables without completely removing it from the reported results.

1. Click **Show Samples** in the View bar.
2. Click **Single View** in the View bar.
3. Click on the row in the experiment pane that contains the injection you want to correct, then click the **Graph** tab.

- Right click the peak in the electropherogram or Peaks table and select **Hide**. Compass for iCE will hide the peak data in the results tables.



- To view hidden peak data, click **View** in the main menu and click **Show Hidden**. Hidden peak data will display in the results table and be marked with an X.

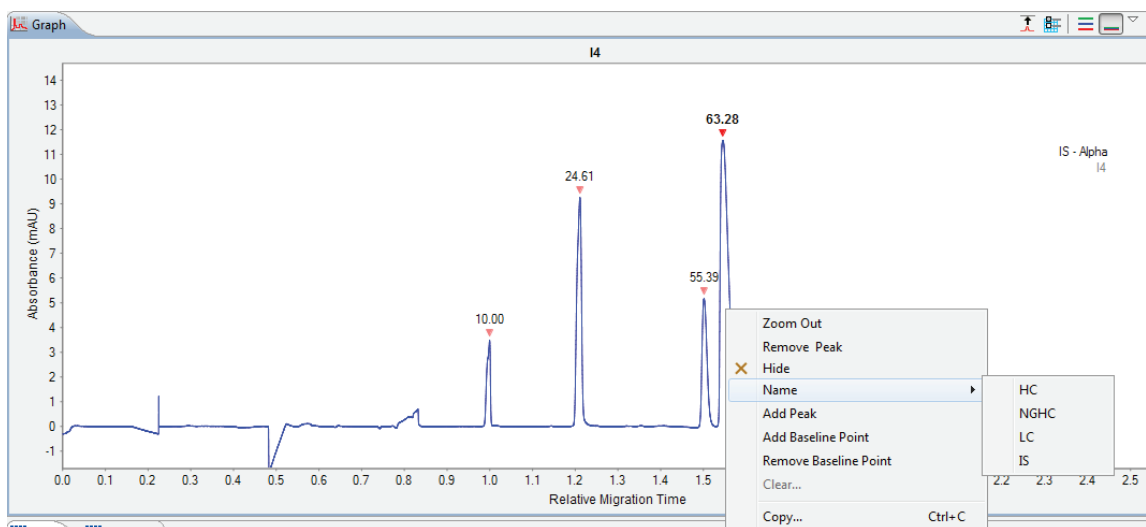


- To unhide a peak, right click on the peak in the electropherogram or peaks table and select **Unhide**.

Changing Peak Names for Sample Data

If Compass for iCE did not automatically name a sample protein peak, you can do it manually.

- Click **Show Samples** in the View bar.
- Click **Single View** in the View bar.
- Click on the row in the experiment pane that has the sample you want to correct, then click the **Graph** pane.
- Right click the peak in the electropherogram or Peaks table and click **Name**, then select a name from the list. Compass for iCE will change the peak name in the electropherogram and results tables, and adjust peak names for other sample proteins accordingly.



NOTES:

For details on how to specify peak name settings, see “Peak Names Settings” on page 723.

Virtual gel data in the lane pane will also update to reflect changes made in the graph pane.

Changing the Electropherogram View

Options in the Graph pane let you zoom and rescale electropherograms, overlay or stack plots and change the peak and plot info displayed.

The Graph pane toolbar has these options:



Auto Scale



Graph Options



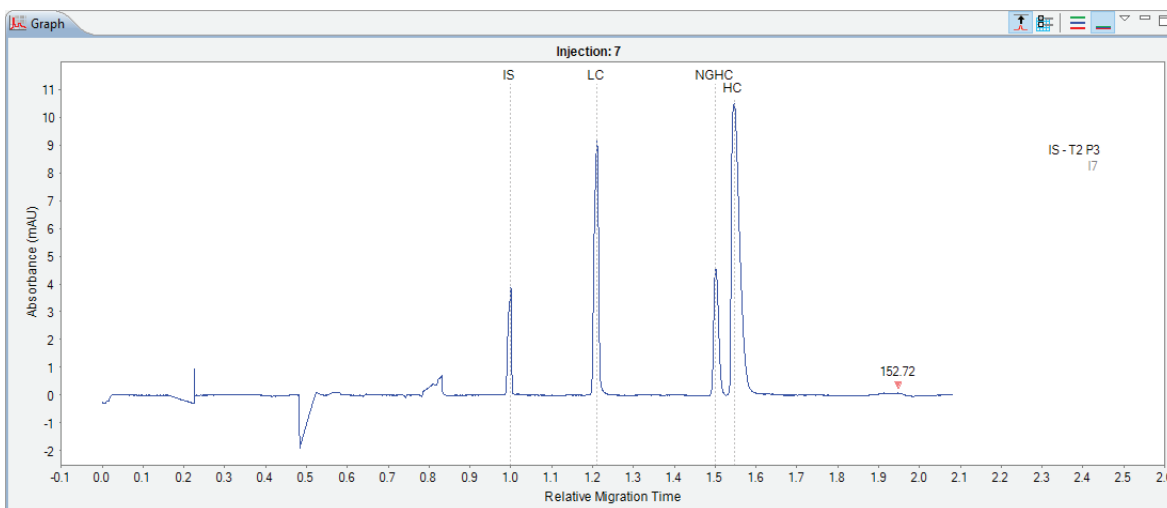
Stack the Plots



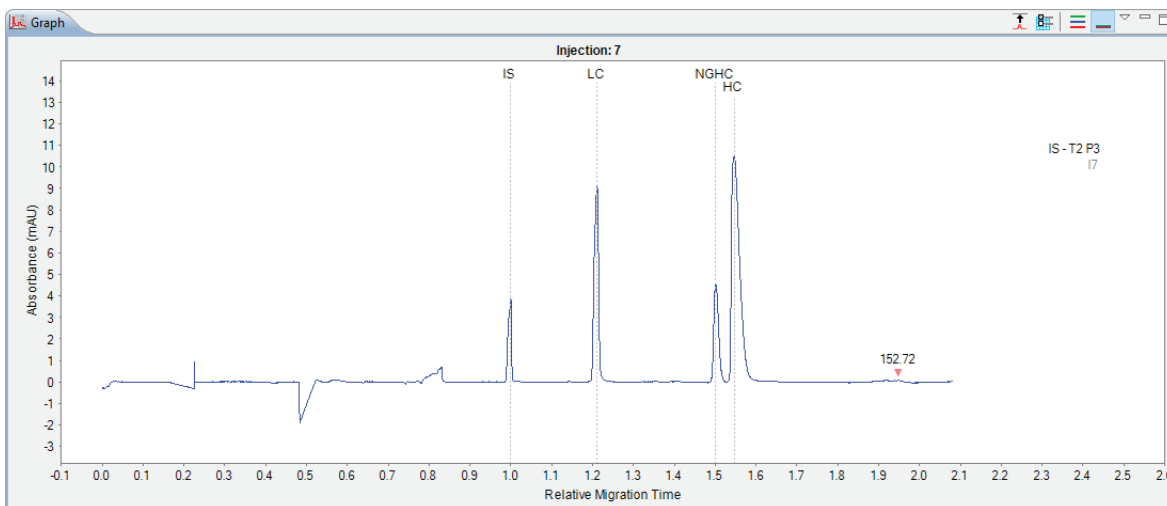
Overlay the Plots

Autoscaling the Electropherogram

Click the **Auto Scale** button to scale the y-axis to the largest peak in the electropherogram.

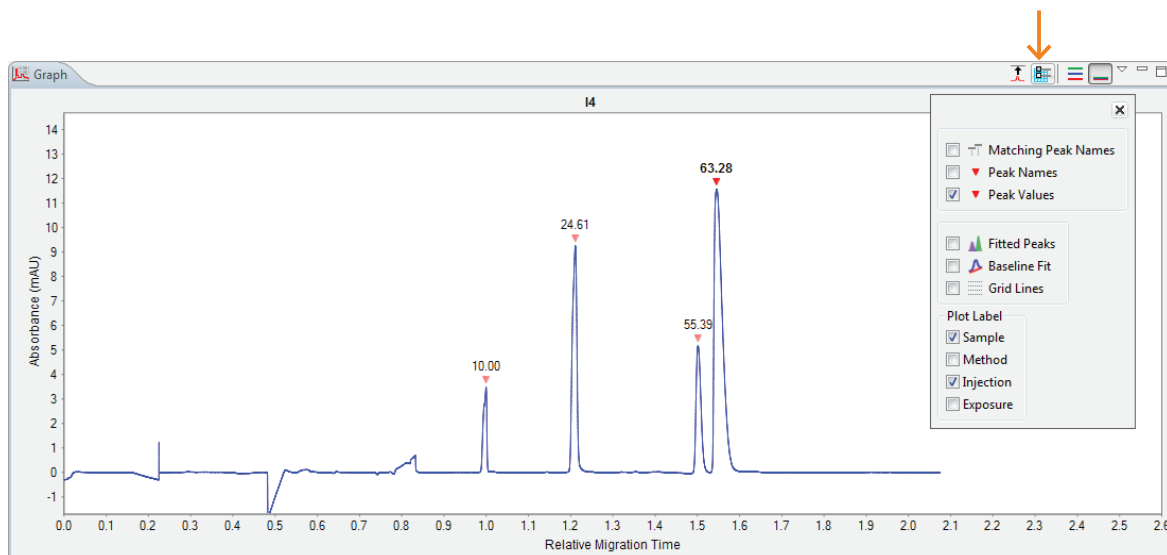


Click the **Auto Scale** button again to return to default scaling.



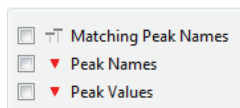
Customizing the Data Display

You can customize electropherogram peak labels, plot labels and display options. To do this, just select the **Graph Options** button.

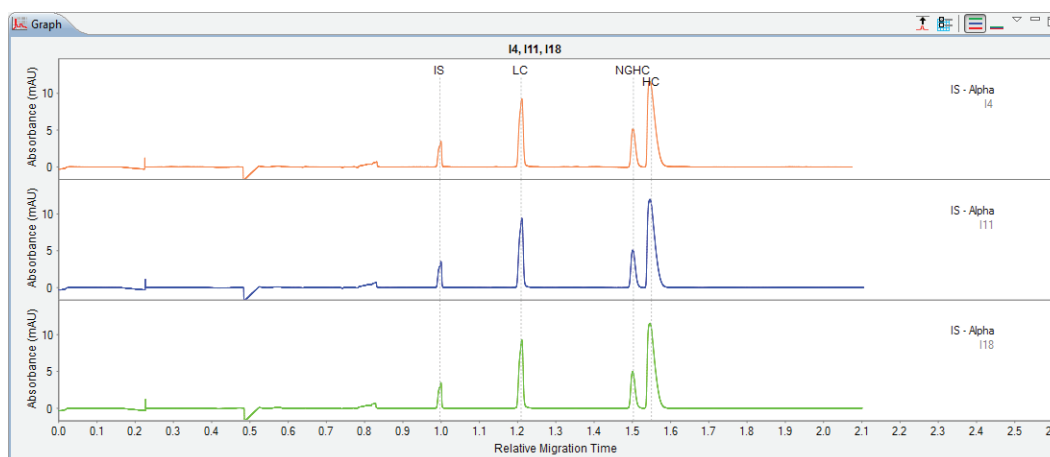


Peak Labels

You can customize the labels used to identify peaks in the electropherogram with these options:

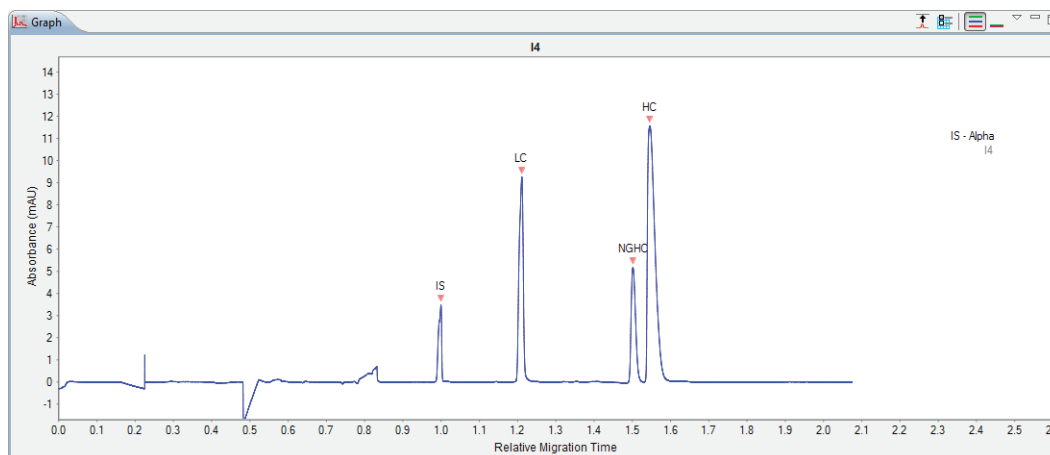


- **Matching Peak Names** - Checking this box will draw vertical lines through each named peak. Using this option with Stack the Plots or Overlay the Plots features is helpful for visually comparing your named peaks across multiple injections.



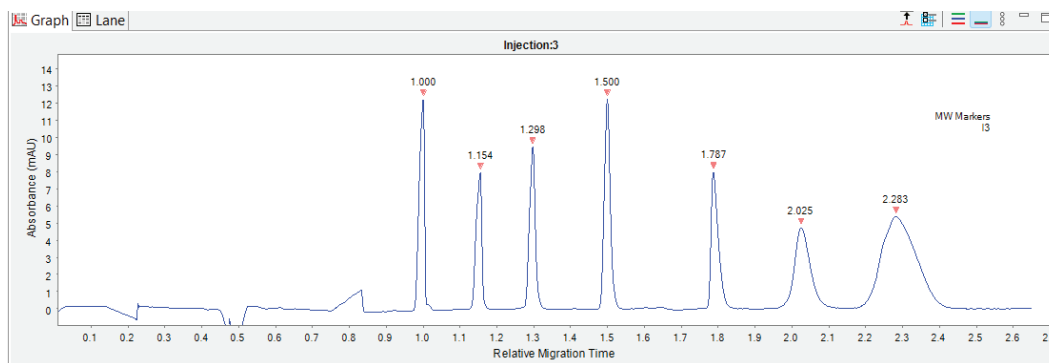
- **Peak Names** - Checking this box displays peak name labels on all named peaks in the electropherogram.

NOTE: If more than one peak label option is selected, peak name labels will always be used for named peaks.



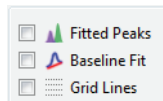
- **Peak Values** - Checking this box will display the molecular weight labels on all peaks in the electropherogram. If a MW Marker wasn't run, RMT values are displayed.

NOTE: If more than one peak label option is selected, peak name labels will always be used for named peaks.



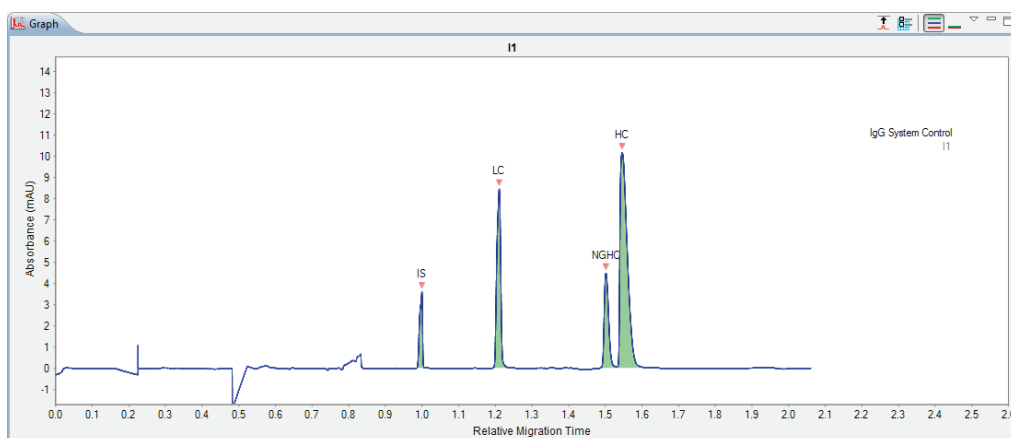
Baseline and Grid Options

You can view the calculated baseline fit, peak integration and show grid lines with these options.



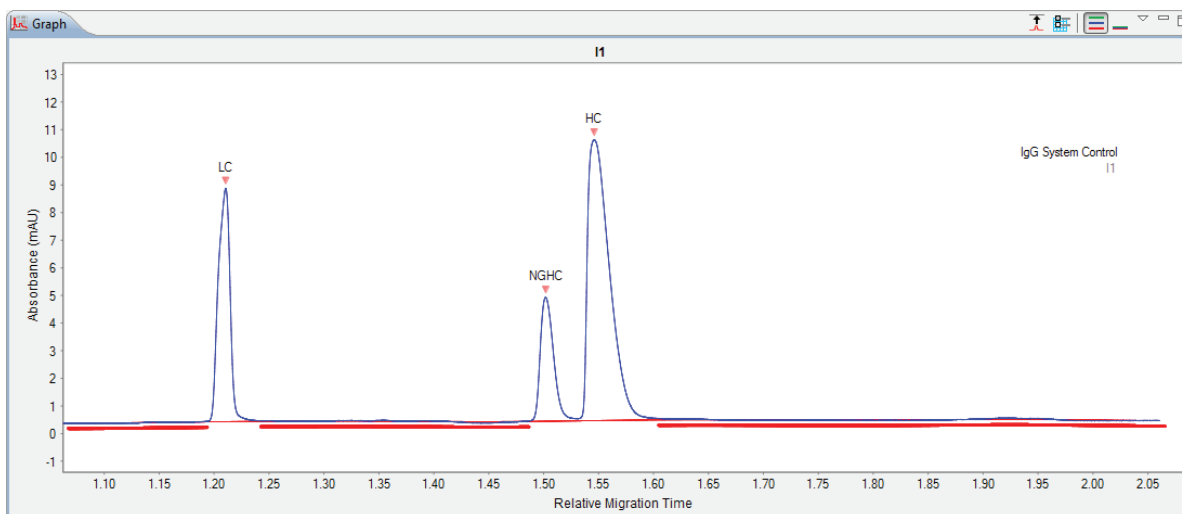
- **Fitted peaks** - Checking this box displays how the peaks were fit by the software. For CE-SDS runs, the software uses Dropped Lines by default.

NOTE: This option is only available for sample data.

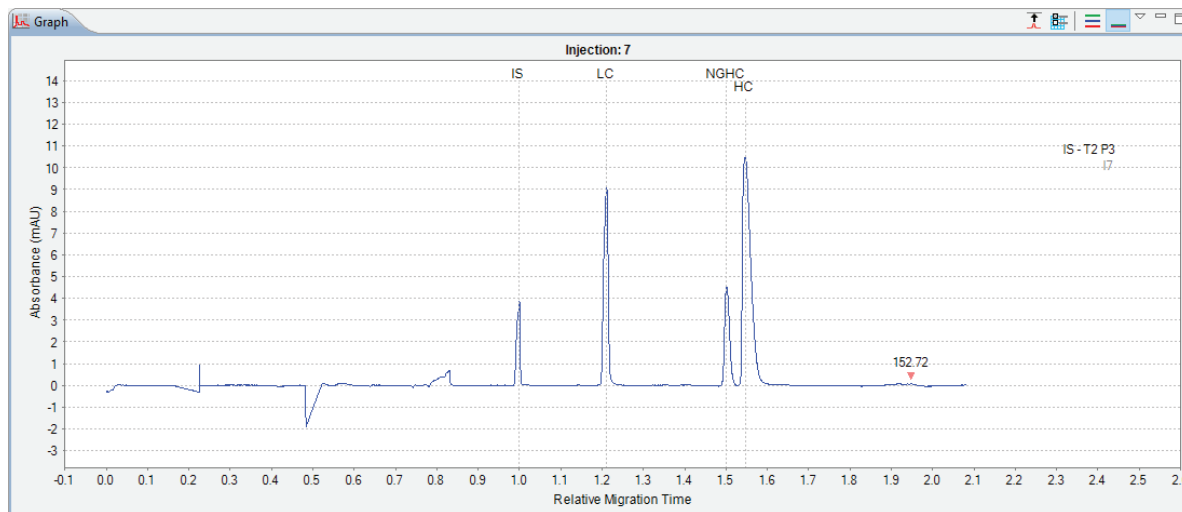


- **Baseline Fit** - Checking this box displays the calculated baseline for the peaks. Depending on your baseline type setting (baseline points or interpolated baseline) baseline points will also display for regions of the electropherogram considered to be at baseline.

NOTE: This option is only available for sample data.



- **Grid Lines** - Checking this box adds grid lines in the graph.



Plot Labels

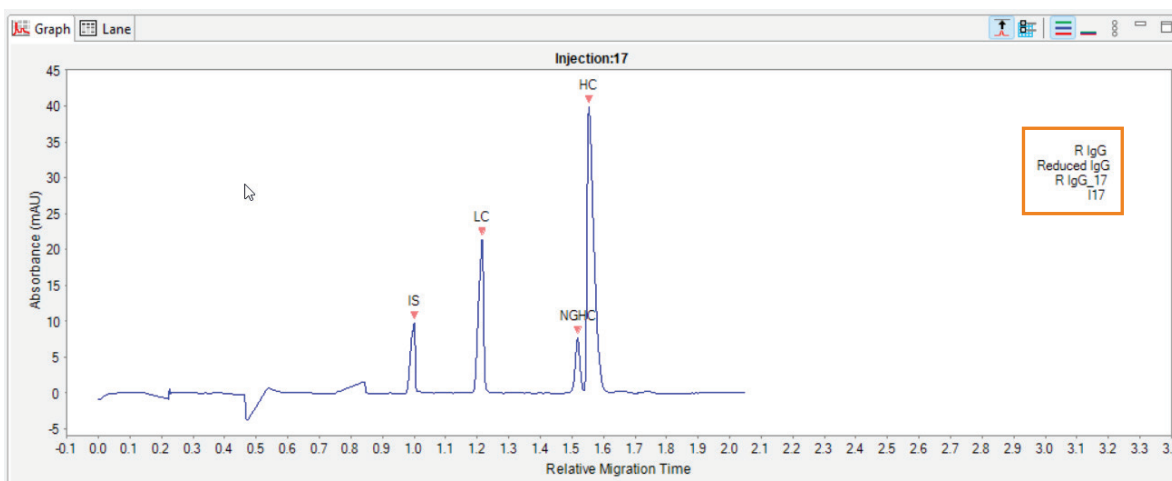
You can customize the plot labels displayed on the electropherogram with these options.

Plot Label	
<input checked="" type="checkbox"/>	Sample
<input type="checkbox"/>	Method
<input checked="" type="checkbox"/>	Injection
<input type="checkbox"/>	Injection Name

Plot labels are shown in the upper right side of the graph.

- **Sample** - Checking this box displays the sample name used for the injection. If sample names were entered with the batch, those names will display here. If not, Sample (default name) displays.
- **Method** - Checking this box displays the method used for the injection.
- **Injection** - Checking this box displays the injection number. For example, I4 for injection 4 in the run.
- **Injection Name** - Checking this box displays the injection name used for the injection. If injection names were entered with the batch, those names will display here. If not, the default name displays.

Here's an example of an electropherogram with all plot labels selected:

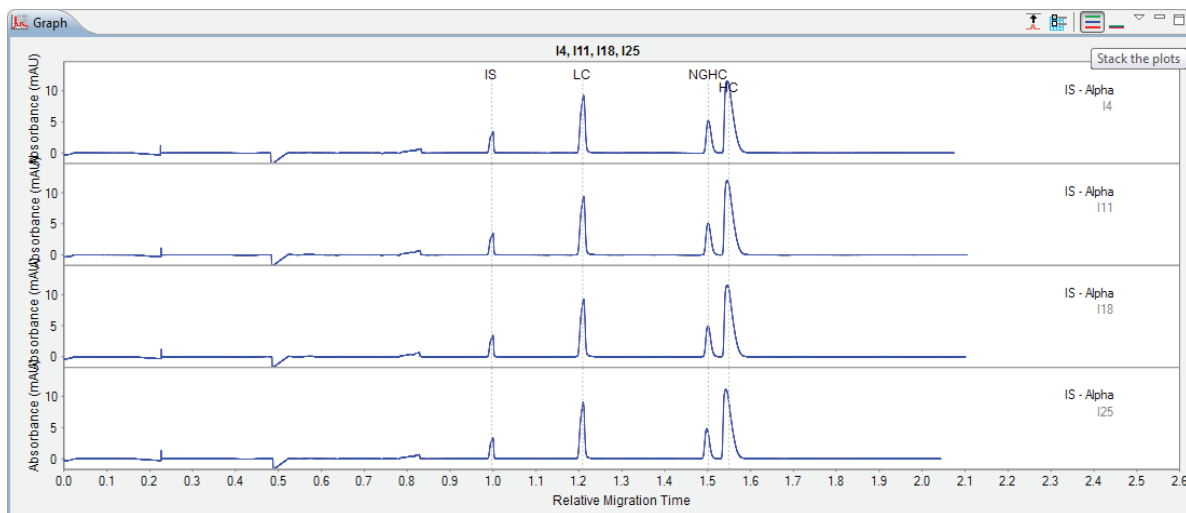


Stacking Multiple Electropherograms

You can stack electropherograms for multiple injections vertically in the Graph pane for comparison.

1. Click **Single View**.
2. Select multiple injection rows in the Experiment pane.

- Click the **Stack the Plots** button. The individual electropherograms for each injection you selected will stack in the Graph pane.

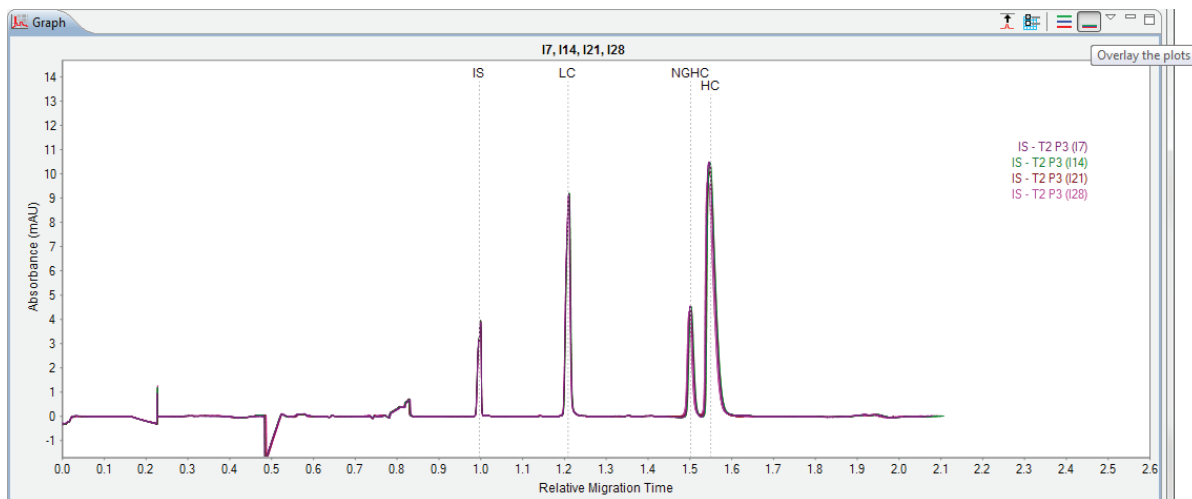


You can also customize the colors used for the stacked plot display. To do that go to “Setting up Automatic Injection Reports” on page 761.

Overlaying Multiple Electropherograms

You can overlay electropherograms for multiple injections on top of each other for comparison in the Graph pane.

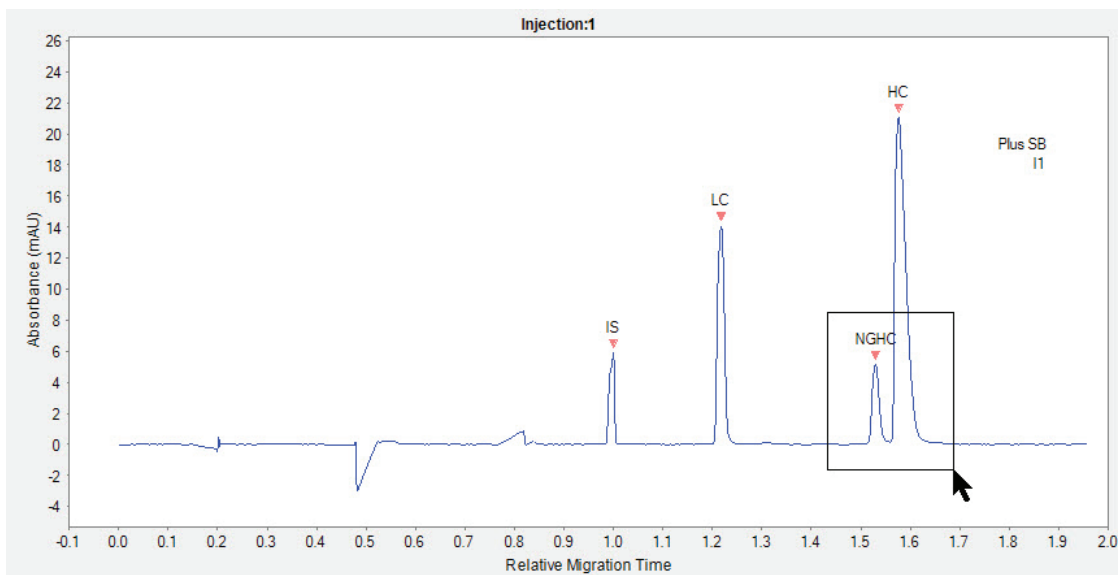
- Click **Single View**.
- Select multiple injection rows in the Experiment pane.
- Click the **Overlay the Plots** button. The individual electropherograms for each injection you selected will overlay in the Graph pane.



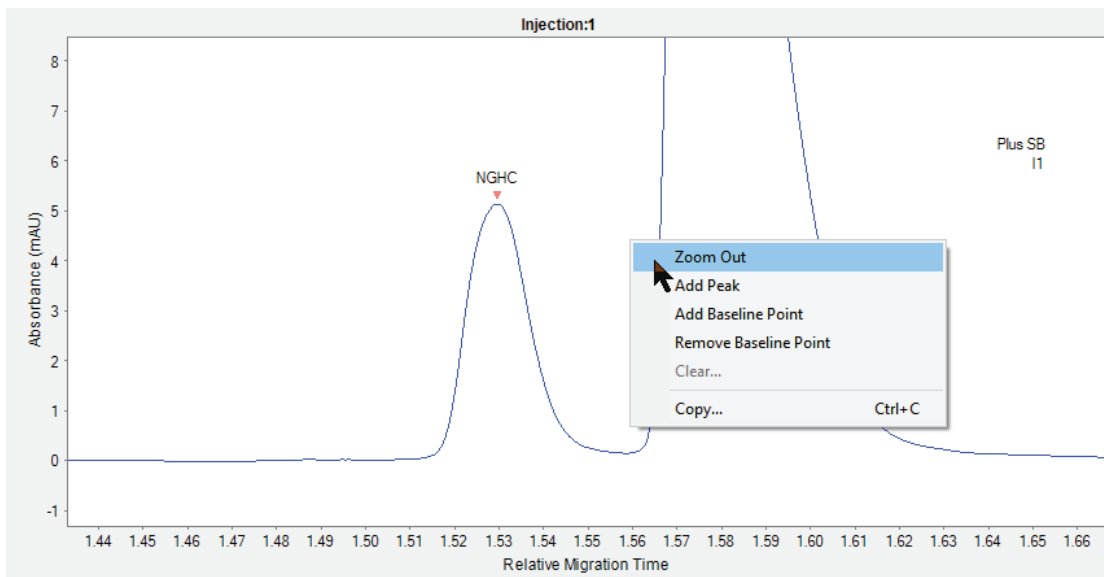
You can also customize the colors used for the overlay plot display. To do that go to “Setting up Automatic Injection Reports” on page 761.

Zooming

To zoom in on a specific area of the electropherogram, hold the mouse button down and draw a box around the area with your mouse:

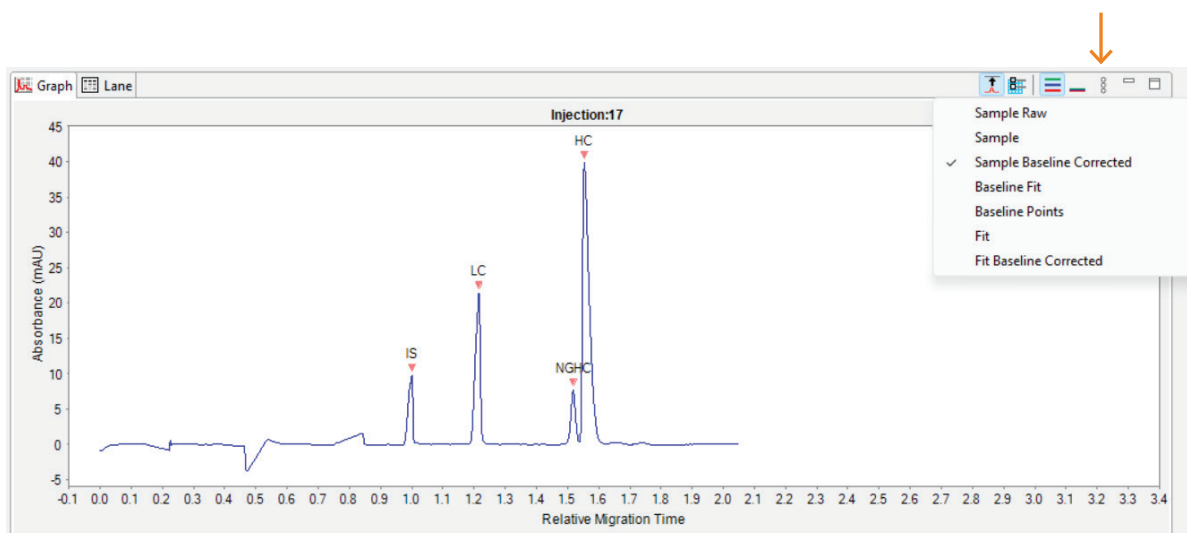


To return to default scaling, right click in the electropherogram and click **Zoom Out**.



Selecting Data Viewing Options

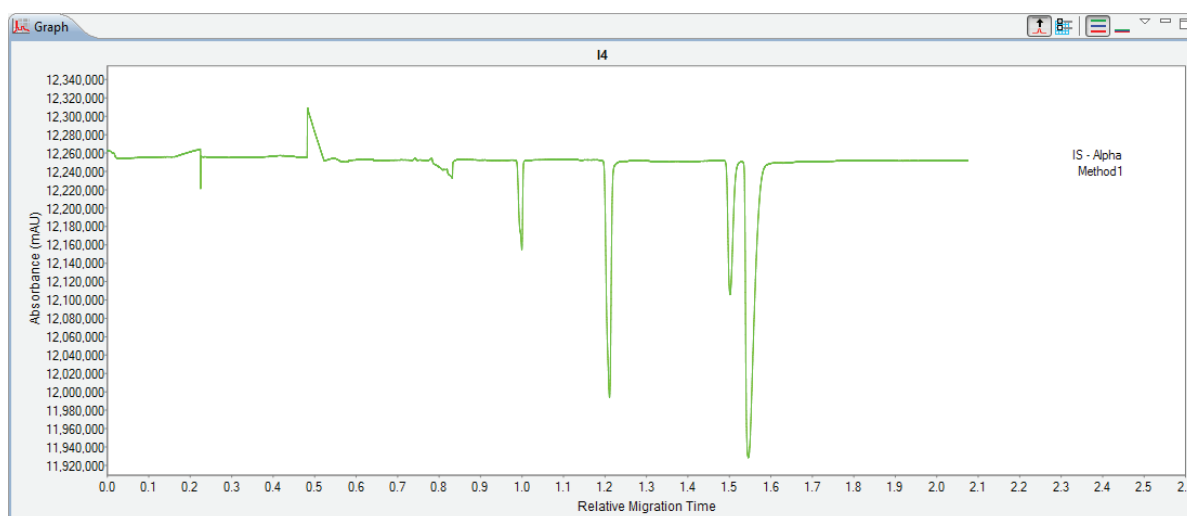
The graph view menu gives you multiple options for changing what type of electropherogram data is displayed. Just click **View More** in the graph pane toolbar to view the menu:



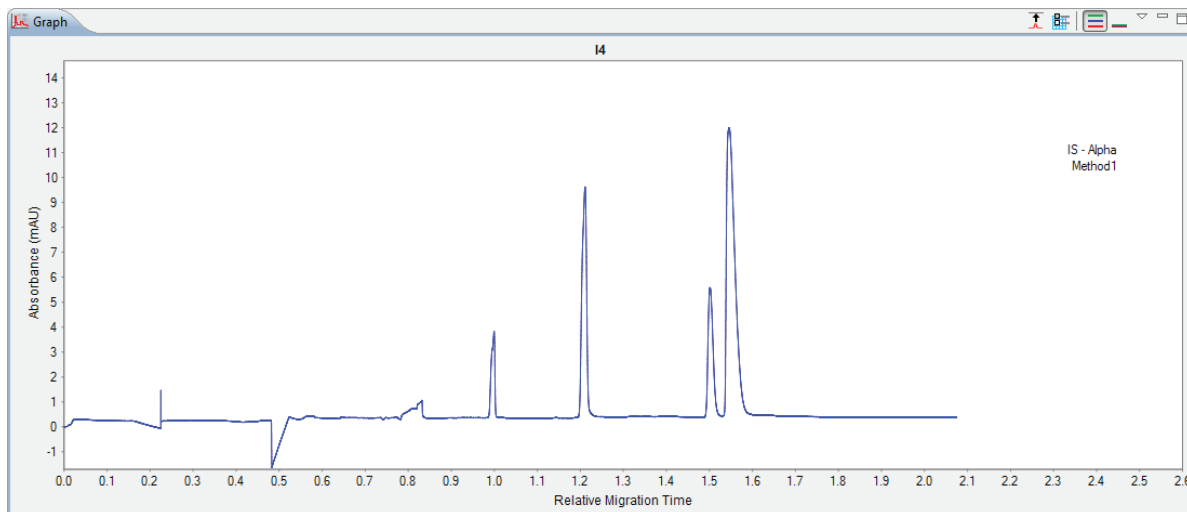
A check mark next to the menu option indicates it's currently selected, and you can select multiple options at once.

NOTE: Unless noted otherwise, graph view menu options are available for sample data only.

- **Sample Raw** - Clicking this option displays the basic detector values used to calculate peak absorbance. To view this you'll need to select **Auto Scale** in the Graph pane tool bar.

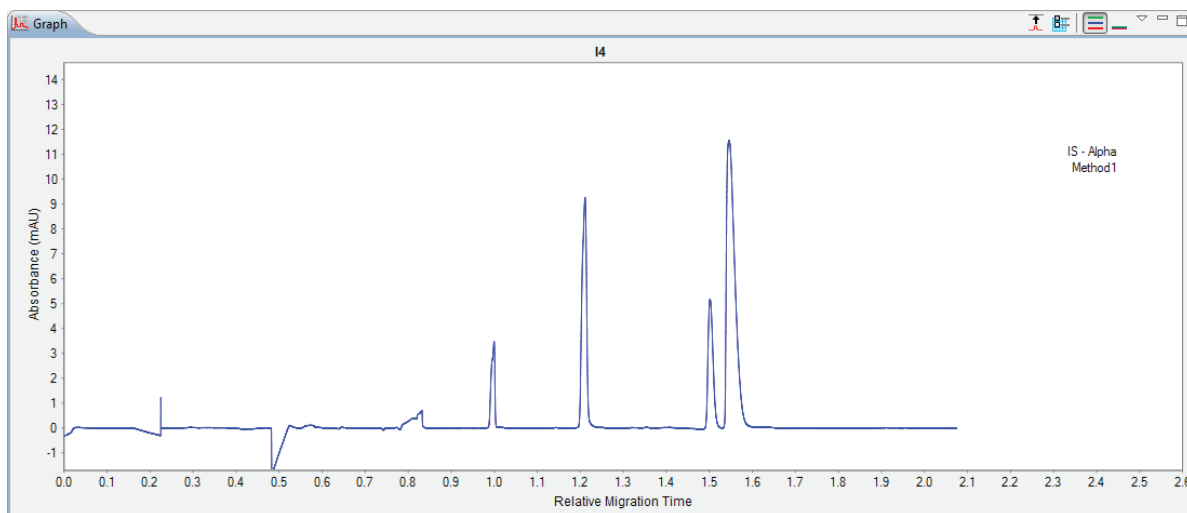


- **Sample** - Clicking this option displays raw, uncorrected sample data.



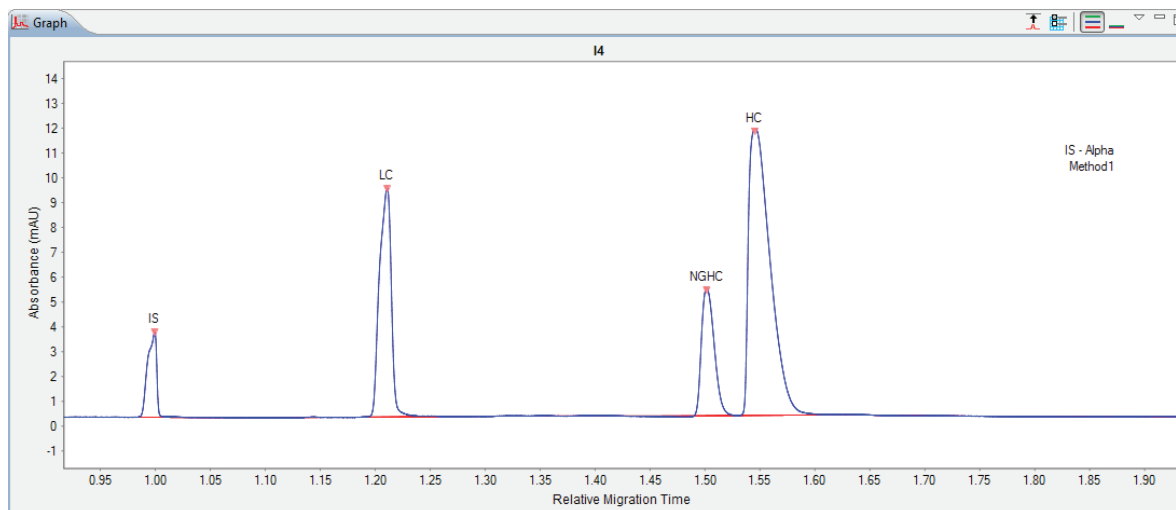
- **Sample Baseline Corrected** - Clicking this option displays sample data with the baseline subtracted (zeroed). This is the default view.

NOTE: The Sample Baseline Corrected data line will not be displayed if Baseline Type is set to Interpolate in the analysis menu option.



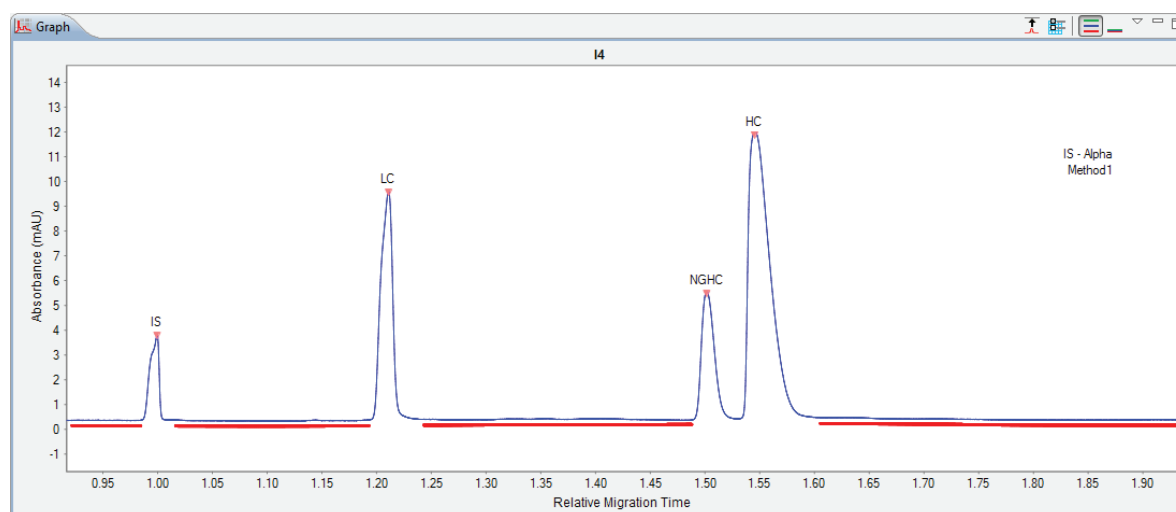
- **Baseline Fit** - Clicking this option displays the calculated baseline for the raw sample data. In this next example, both Baseline Fit and Sample are selected.

NOTE: This option is selected automatically when Baseline Fit is selected in graph options.

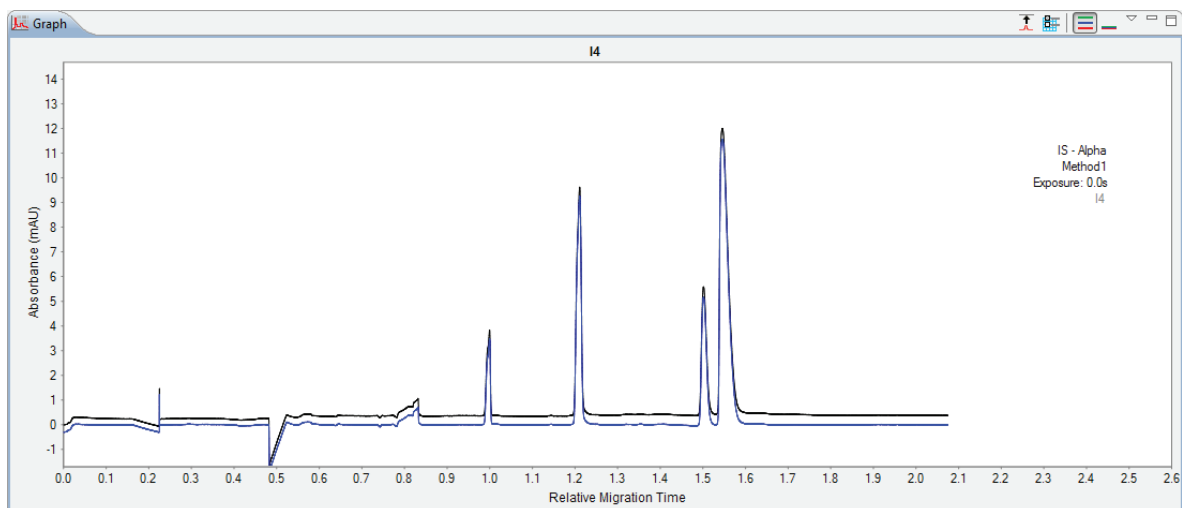


- **Baseline Points** - Clicking this option displays regions of the electropherogram considered to be at baseline. In this example, both Baseline Points and Sample are selected.

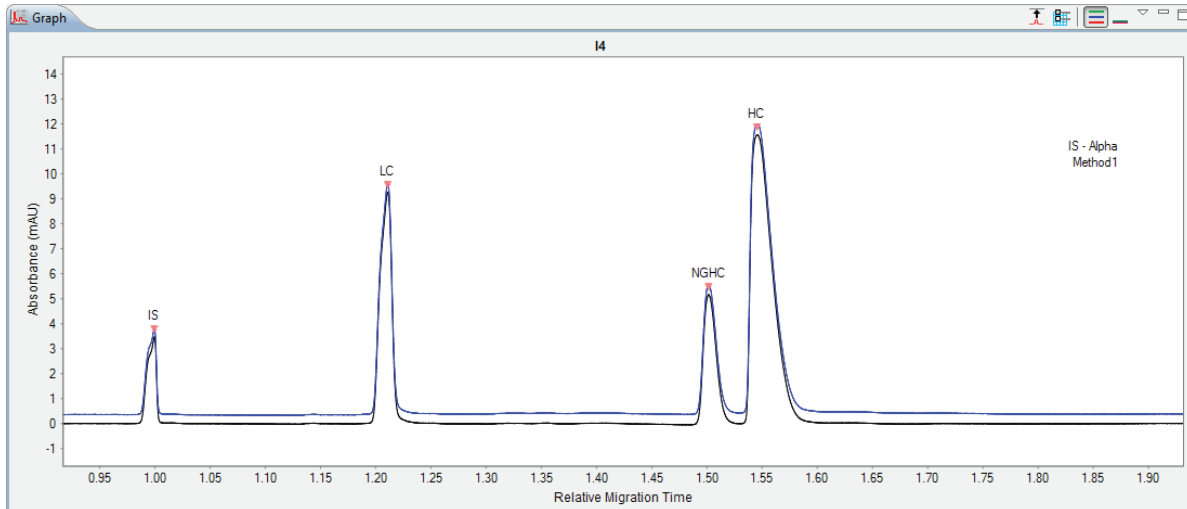
NOTE: This option is selected automatically when Baseline Fit is selected in graph options.



- **Fit** - Clicking this option displays the bounding envelope of the fitted peaks as calculated by the software for the raw sample data. In this example, both Fit and Sample Baseline Corrected are selected.



- **Fit Baseline Corrected** - Clicking this option displays the fitted peaks as calculated by the software for the sample baseline corrected data. In this example, both Fit Baseline Corrected and Sample are selected, the fit plot is on the bottom.

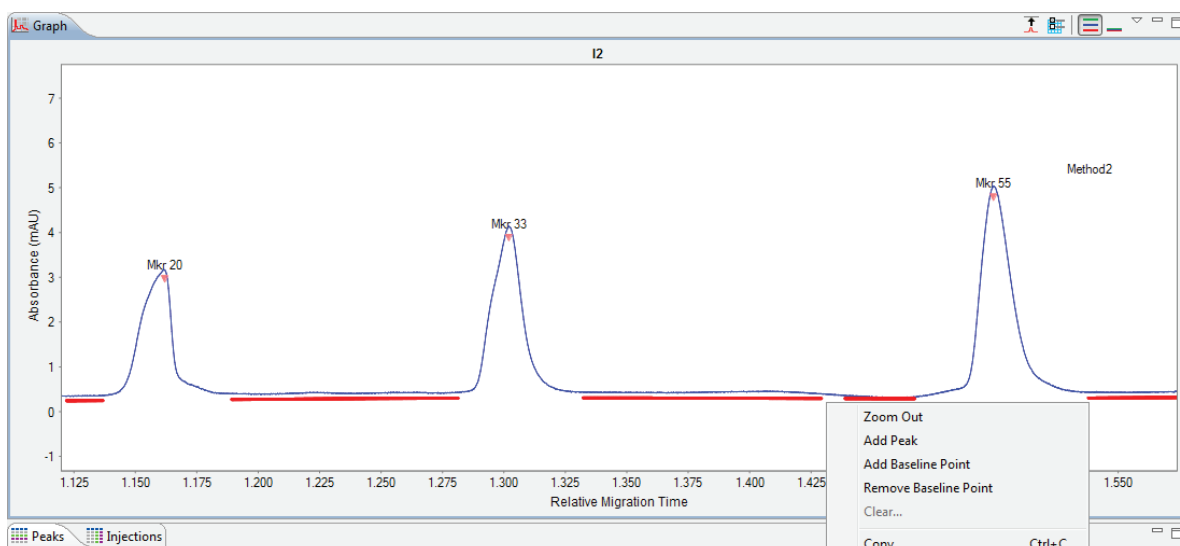


Adjusting the Baseline

Spline Baseline

Points in the baseline can be added or removed as needed.

1. Click the **Graph Options** button in the graph pane toolbar and check **Baseline Points**. This will display baseline points for the raw sample data.
2. Use the mouse to draw a box around the area you want to correct. This will zoom in on the area.
3. Right click a baseline point and select **Add Baseline Point** or **Remove Baseline Point**.



NOTE: To clear the manual addition or removal of baseline points and go back to the original view of the data, right click in the electropherogram and click **Clear All**.

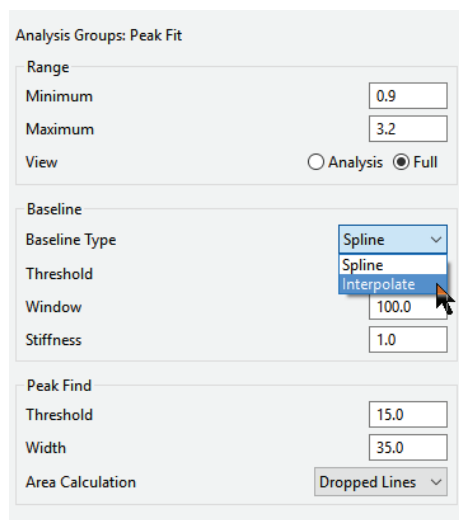
Interpolated Baseline

Compass for iCE lets you interpolate baselines from the base of each peak for named peaks that use Dropped Lines integration.

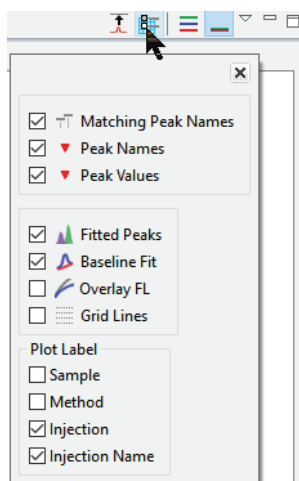
1. If you haven't already, name your peaks of interest.
2. Select **Edit > Analysis**, and click **Peak Fit** in the left sidebar.

- Choose **Interpolate** as the Baseline Type if it isn't already selected.

NOTE: Baseline Type selection is only available when Dropped Lines is used for the Area Calculation.

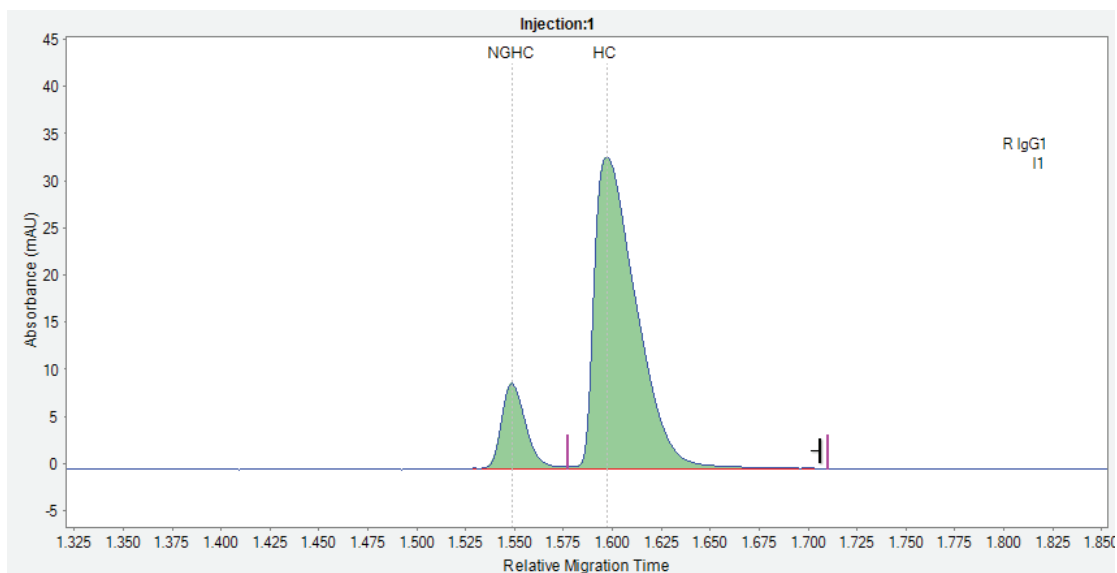


- Click **OK**.
- In the Analysis window Graph Pane, click **Graph Options** and select **Fitted Peaks** and **Baseline Fit**.

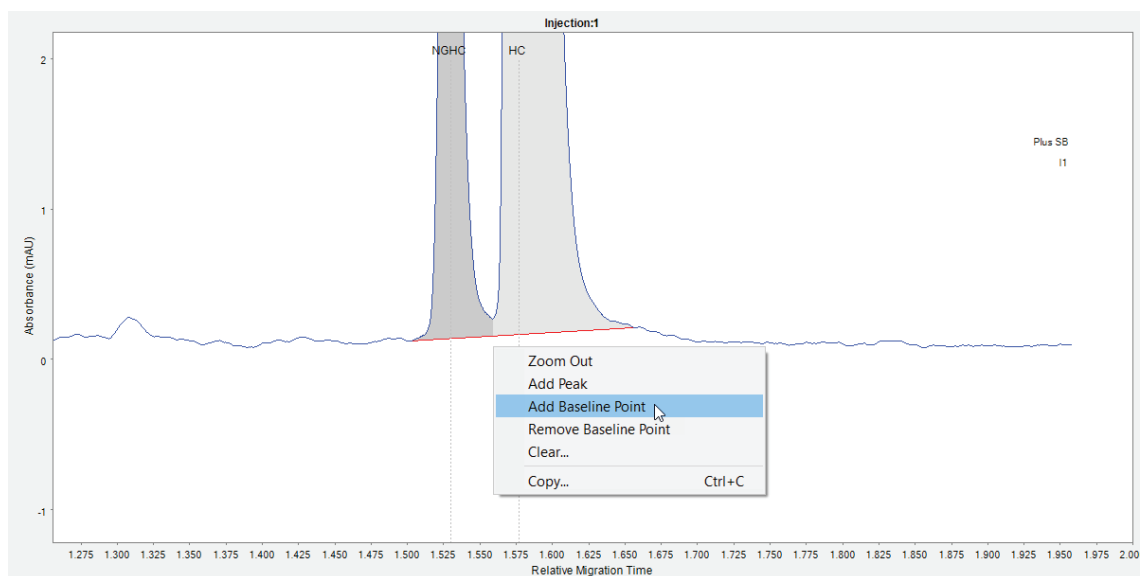


- Select an injection in the Experiment pane.
- Hover the mouse over the peak you want to adjust the baseline on. Two magenta bars will display that show the start and end points of the peak's integration.

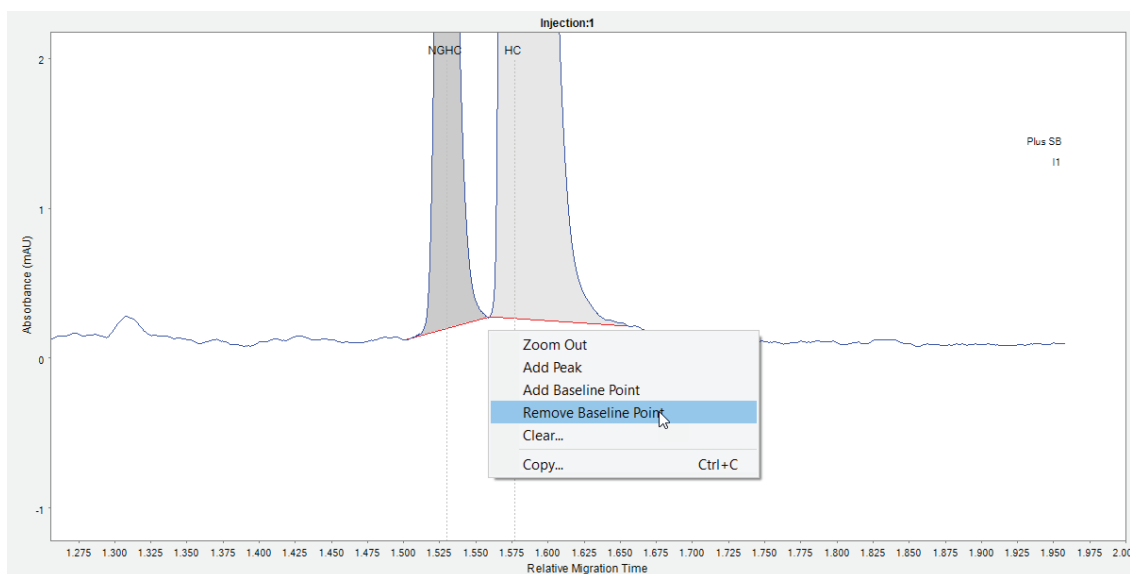
8. To change the baseline, hover the mouse over one of the magenta bars. Then click the mouse and drag the bar to move it. The cursor changes to indicate which direction you can move the integration point. See “Manual Peak Integration” on page 739 for more info on integrating peaks.



9. Baseline points can also be added between two adjacent peaks for a better fit. Hover the mouse over the peak until the magenta bars appear, then right click and select **Add Baseline Point**.

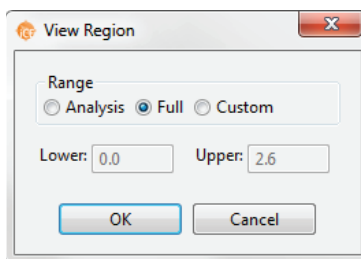


10. To remove the baseline point, hover the mouse over the peak until the magenta bars appear, then right click and select **Remove Baseline Point**.

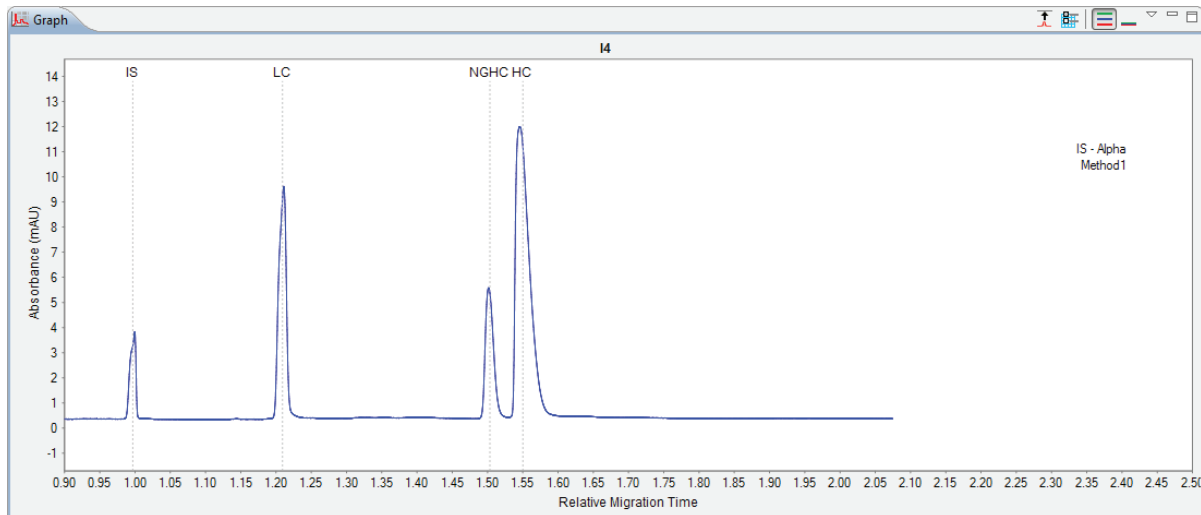


Selecting the Graph X-axis Range

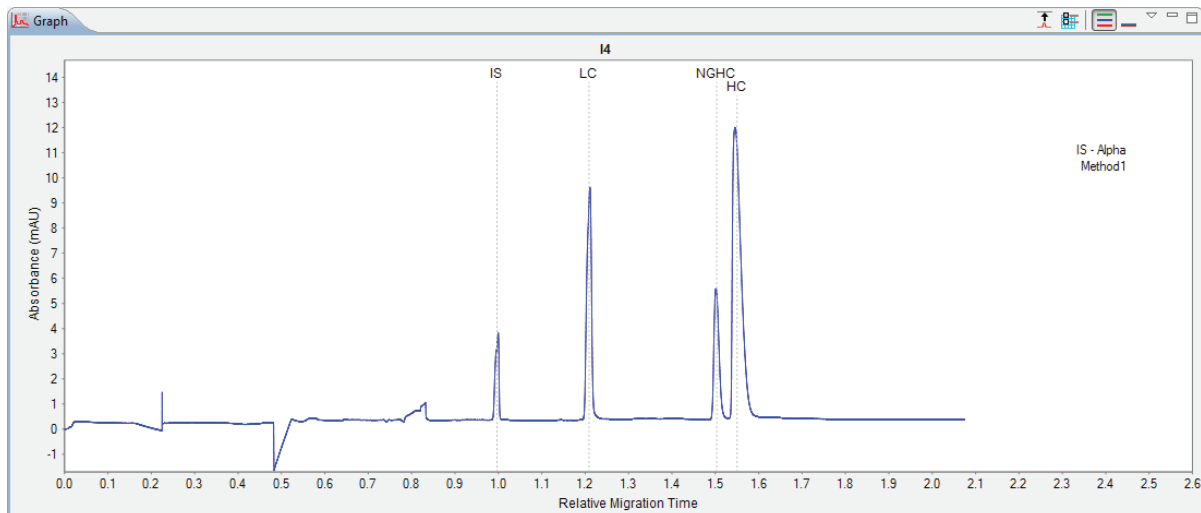
The RMT (relative migration time) range used for the x-axis can be changed. Just select **View** in the main menu and click **View Region**.



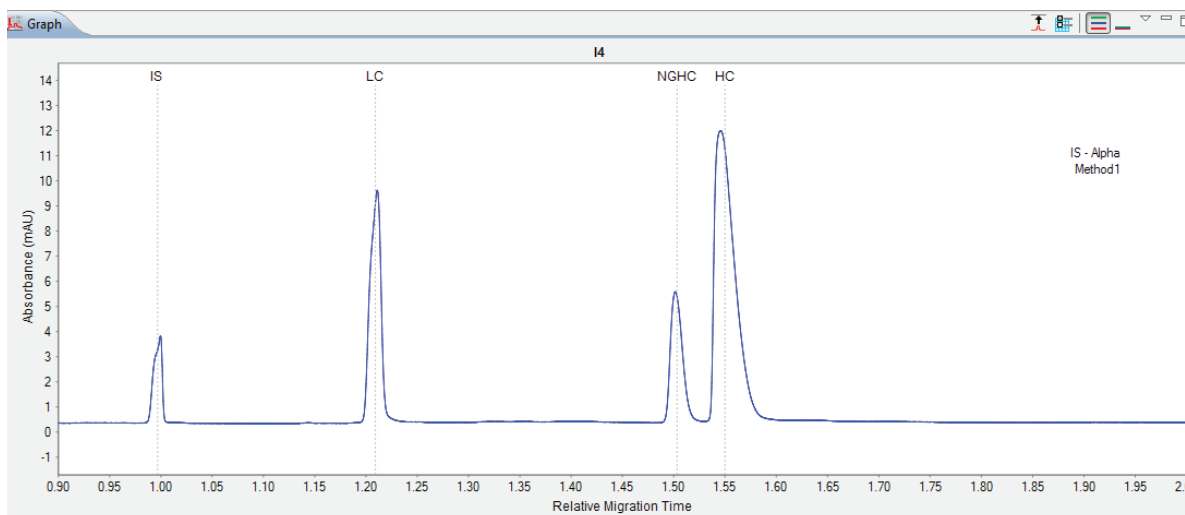
- **Analysis** sets the x-axis range of the electropherogram to what is selected in the Peak Fit range settings. To view or change these analysis settings, go to **Edit > Analysis** and click **Peak Fit** in the left sidebar. In this example, the lower and upper range settings are 0.9 and 2.5.



- **Full** displays the entire separation in the electropherogram. This is the default setting. In this example the lower and upper range settings are 0 and 2.6.



- **Custom** lets you manually enter the lower and upper range settings to display in the electropherogram. In this example the lower and upper range settings are 0.9 and 2.0.



NOTE: You can change the default x-axis range that Compass for iCE uses. Go to “Advanced Analysis Settings” on page 743 for more info.

Changing the Virtual Gel View

Options in the lane pane let you change the contrast or invert the virtual gel or change lane labels.

The Lane pane toolbar has the following options:



Contrast Adjustment



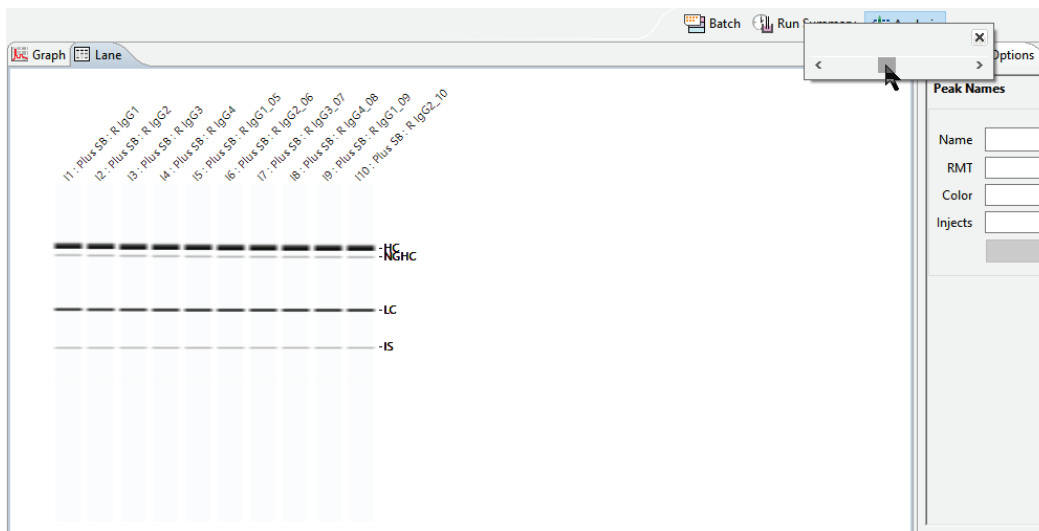
Invert



Lane Options

Adjusting the Contrast

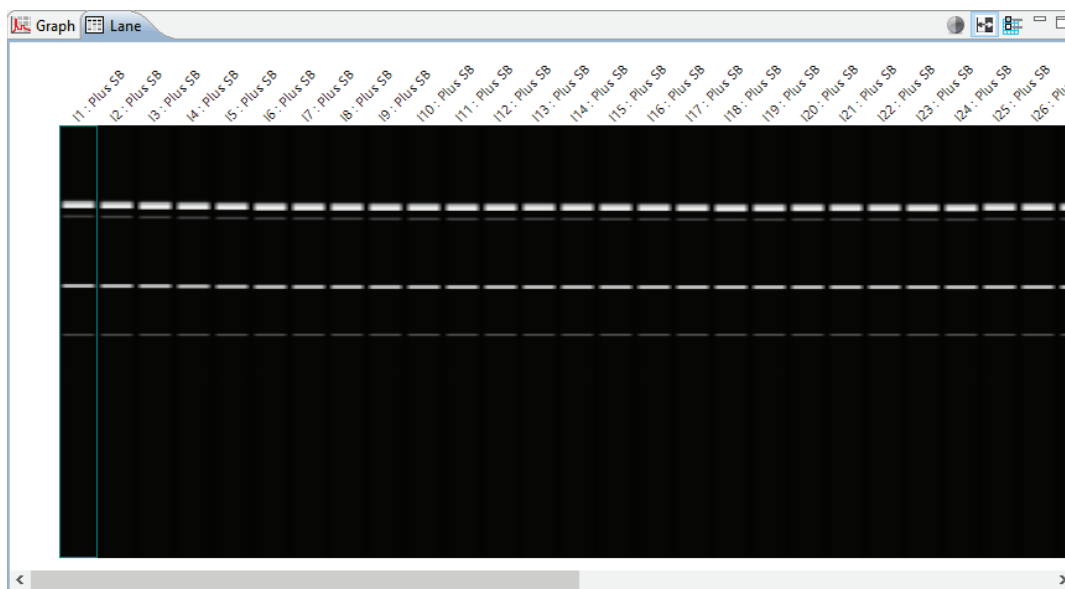
1. Click the **Contrast Adjustment** button. The contrast tool will display:



2. Use the mouse to move the slider left or right until the desired contrast is reached.

Inverting the Virtual Gel

1. Click the **Invert** button. The virtual gel image will invert:

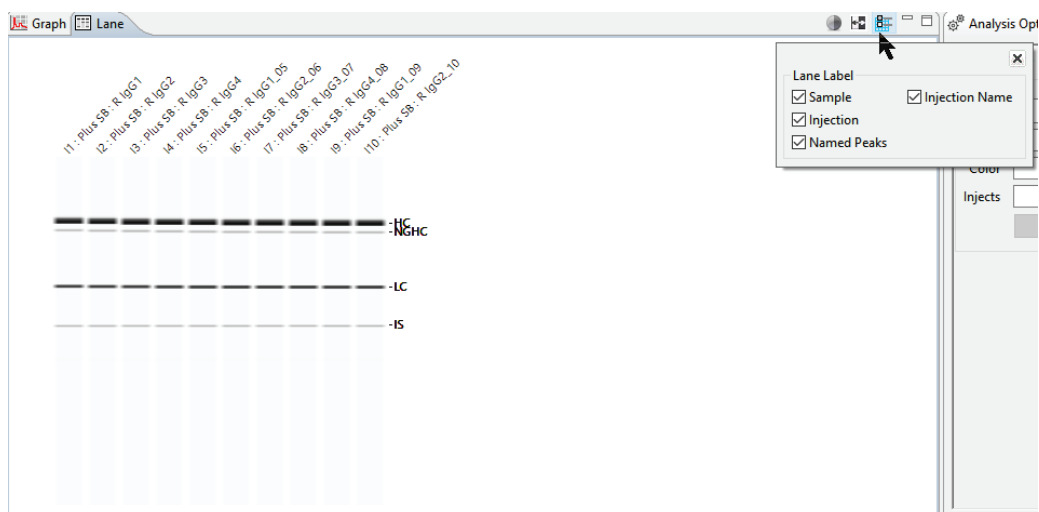


2. Click the **Invert** button again to return to the default view.

Selecting Lane Labels

The labels shown above the lanes in the virtual gel can be customized. To do this:

1. Click the **Lane Options** button. The label box will display:



2. The following lane display options are available:

- **Sample** - Displays the sample IDs entered in the batch above each lane.
- **Injection** - Displays the batch injection number above each lane.
- **Named peaks** - Displays peak names (if entered) next to the individual bands in the lane view.
- **Injection name** - Displays the injection name entered in the batch above each lane.

Closing Run Files

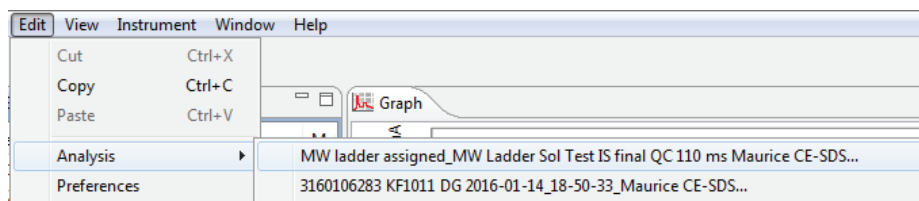
If more than one run file is open, you can close just one file or all the open files at the same time.

- **To close one run file** - In the Experiment pane, click on one of the sample rows in the file. Then click **File** from the main menu and click **Close**.
- **To close all open run files** - Select **File** from the main menu and click **Close All**.

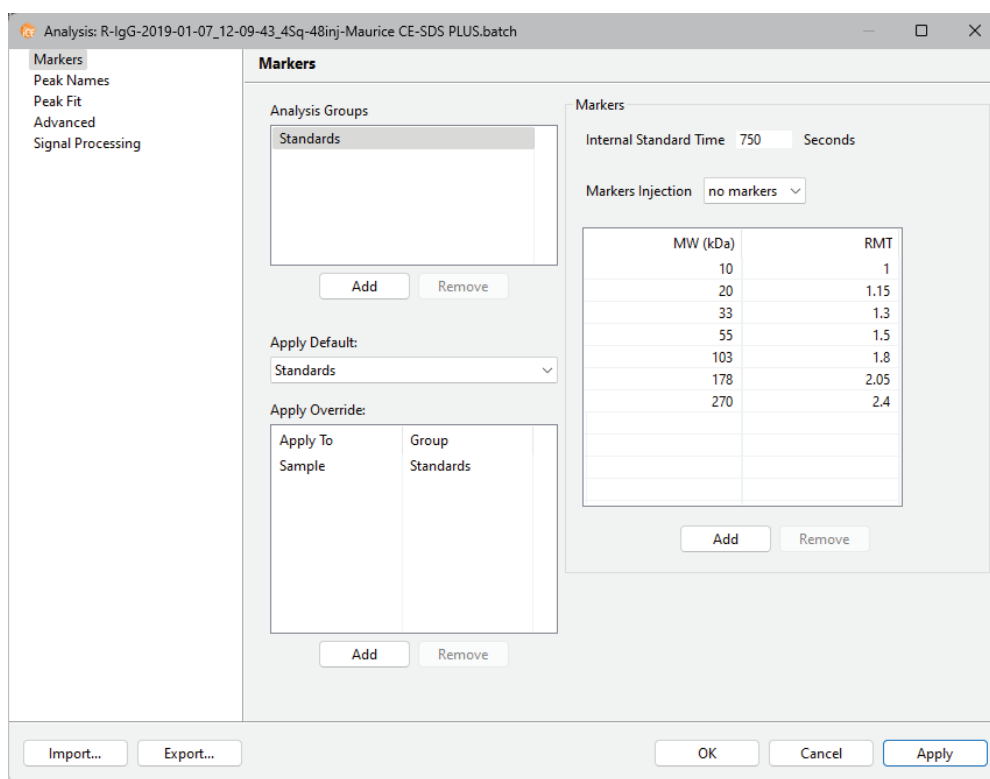
Analysis Settings Overview

Compass for iCE has many analysis features and settings that you can change to enhance your run data.

Select **Edit** in the main menu and click **Analysis**. If more than one run file is open, select the run file you want to view settings for from the list:



This opens the Analysis window:



To move between pages in the window, click on an option in the left sidebar.

- **Markers** - Lets you customize the Internal Standard migration time, and the molecular weight and RMT Compass for iCE uses to identify your CE-SDS MW Markers.
- **Peak Names** - Lets you enter custom naming settings for sample proteins and have Compass for iCE automatically label the peaks in the run data.
- **Peak Fit** - Lets you customize peak fit settings for sample data.
- **Advanced** - Lets you customize analysis settings for the Internal Standard.
- **Signal Processing** - Lets you apply reference signal corrections and baseline smoothing options to your data.

On all pages in the Analysis window:

- Click **Import** to import an analysis settings file. Go to “Importing Analysis Settings” on page 756 to learn how to do this.
- Click **Export** to export the current analysis settings file. Go to “Exporting Analysis Settings” on page 757 to learn how to do this.
- Click **Apply** to apply changes to the run file and update results in real time.
- Click **OK** to save changes to the run file and exit.
- Click **Cancel** to exit without saving changes.

Markers Analysis Settings

This page lets you select the injection for your CE-SDS MW Markers, enter a list of molecular weights and RMTs for each marker peak, and set the expected migration time of the Internal Standard for all injections. Select **Edit** in the main menu and click **Analysis**, then click **Markers** in the left sidebar.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they’re completed. If you make analysis settings changes to an executing run, they won’t be saved to the final run file.

Analysis: R-IgG-2019-01-07_12-09-43_45q-48inj-Maurice CE-SDS PLUS.batch

Markers

Peak Names
Peak Fit
Advanced
Signal Processing

Markers

Analysis Groups

Standards

Add Remove

Apply Default:
Standards

Apply Override:

Apply To	Group
Sample	Standards

Add Remove

Markers

Internal Standard Time 750 Seconds

Markers Injection no markers

MW (kDa)	RMT
10	1
20	1.15
33	1.3
55	1.5
103	1.8
178	2.05
270	2.4

Add Remove

Import... Export... OK Cancel Apply

Markers Settings

- **Internal Standard Time** - The approximate migration time (in seconds) of the Internal Standard. This is applied to all injections.

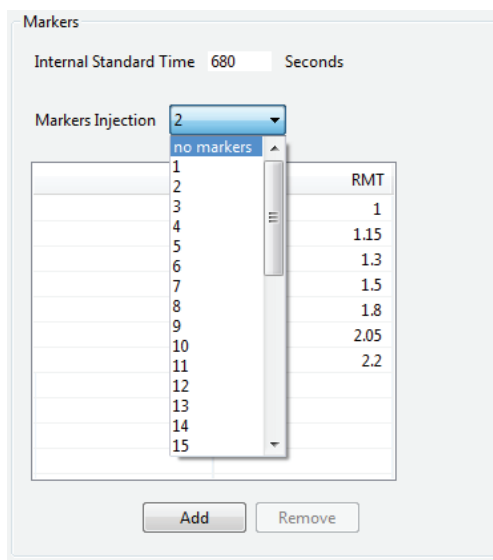
Changing the Injection Used for the CE-SDS MW Markers

You can use known markers to calculate molecular weights of your unknown sample proteins. You can select the injection you ran your CE-SDS MW Markers in, or opt to not use one.

NOTE: When the markers injection is set to no markers, the molecular weight for sample proteins in the run isn't displayed.

To change the markers injection:

1. Select **Edit > Analysis**, and select **Markers** in the left sidebar.
2. Click the arrow in the drop down list next to Markers Injection, then select an **injection number** or **no markers** from the list.



Compass for iCE will use the data in the selected injection to calculate molecular weights for sample proteins in the run data using the information in the table. If no markers is selected, Compass for iCE doesn't display molecular weight for sample proteins.

Standards Analysis Settings Groups

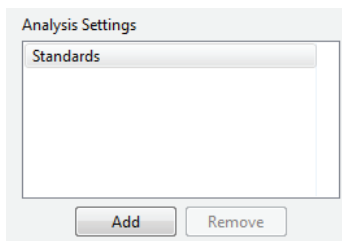
Standards settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values. These settings are included in the default Standards group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 756.

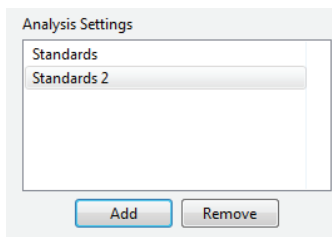
Standards groups are displayed in the analysis settings box:



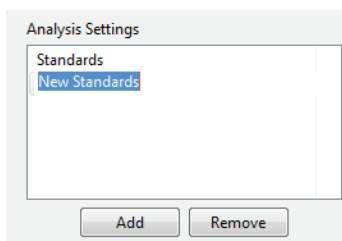
The Standards group shown uses the Compass for iCE default settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Standards Group

1. Select **Edit > Analysis**, and select **Markers** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



3. Click on the new group and enter a new name.



4. Change the Internal Standard time as needed.

Markers

Internal Standard Time 750 Seconds

5. Click the arrow in the drop down list next to Markers Injection, then click an injection number or no markers from the list.

Markers

Internal Standard Time 680 Seconds

Markers Injection 2

no markers

	RMT
1	
2	
3	1
4	1.15
5	1.3
6	1.5
7	1.8
8	2.05
9	2.2
10	
11	
12	
13	
14	
15	

Add Remove

Compass for iCE will use the data in the selected injection to recalculate molecular weights for sample proteins in the run data using the information in the table. If no markers is selected, Compass for iCE doesn't display molecular weight for sample proteins.

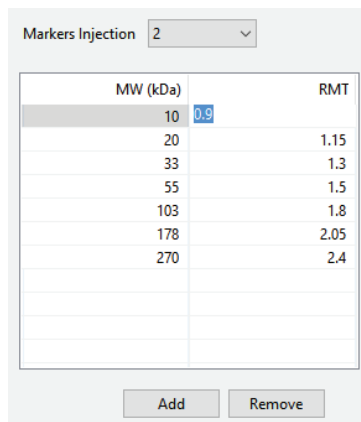
6. If a markers injection was selected, the default Maurice CE-SDS MW Markers molecular weights and relative migration time (RMT) values are already populated in the table. If you'd like to use these values, skip to the next step. If you're using different markers, here's how to change the values:
 - a. Click in the first cell in the MW column in the table and enter the molecular weight (in kDa) for the marker.

Markers Injection 2

MW (kDa)	RMT
15	1
20	1.15
33	1.3
55	1.5
103	1.8
178	2.05
270	2.4

Add Remove

- b. Click in the first cell in the RMT column and enter a value for the marker.



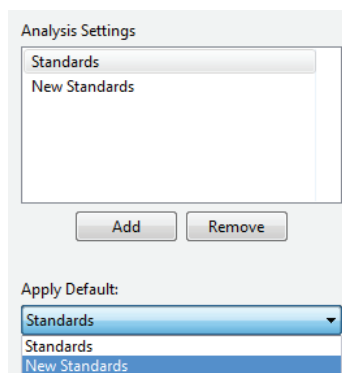
MW (kDa)	RMT
10	0.9
20	1.15
33	1.3
55	1.5
103	1.8
178	2.05
270	2.4

Markers Injection: 2

Add Remove

NOTE: Marker peak positions are relative to each other. Only the difference in RMT is used to help identify them. When entering marker peak information for the first time, review the marker data in the Analysis screen to find the correct peak RMT.

- c. Repeat the steps above for the remaining markers in the table.
- **To add another marker** - Click **Add** under the table, then change the information in the new row.
 - **To remove a marker** - Select its row and click **Remove**.
7. To use the new group as the default settings for the run, click the arrow in the drop down list next to Apply Default, then click the new group in the list. The settings in the new group will then be applied to the run data.



Analysis Settings

Standards

New Standards

Add Remove

Apply Default:

Standards

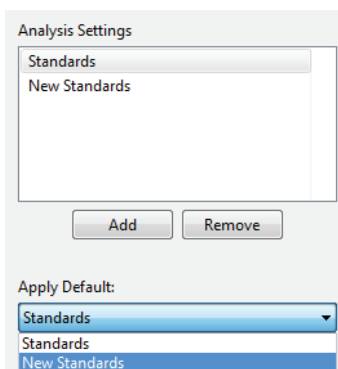
Standards

New Standards

8. Click **OK** to save changes.

Changing the Default Standards Group

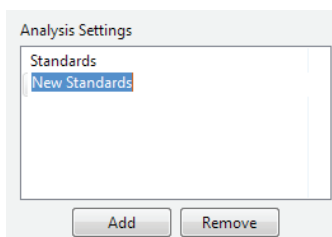
1. Select **Edit > Analysis**, and click **Markers** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then select a new default group from the list.



3. Click **OK** to save changes. Analysis settings in the group selected will be applied to the run data.

Modifying a Standards Group

1. Select **Edit > Analysis**, and click **Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.

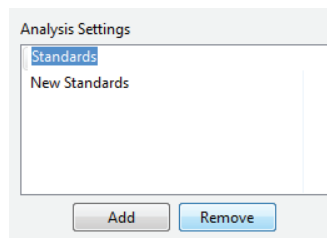


3. Change the marker info as needed as in **Creating a New Standards Group**.
4. Click **OK** to save changes. The new analysis settings will be applied to the run data.

Deleting a Standards Group

1. Select **Edit > Analysis**, and click **Markers** in the left sidebar.

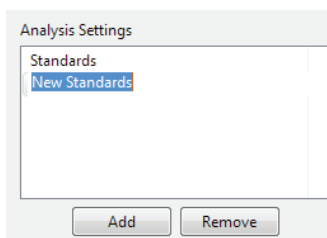
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



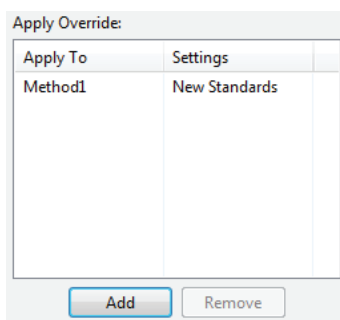
3. Click **OK** to save changes.

Applying Standards Groups to Specific Run Data

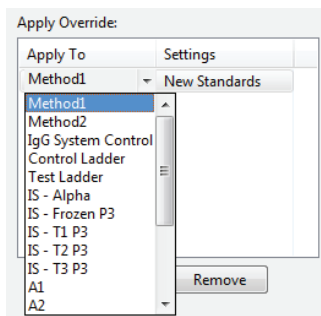
1. Select **Edit > Analysis**, and select **Markers** in the left sidebar.
2. Click on the group in the analysis settings box you want to apply to specific run data.



3. Application of standards groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

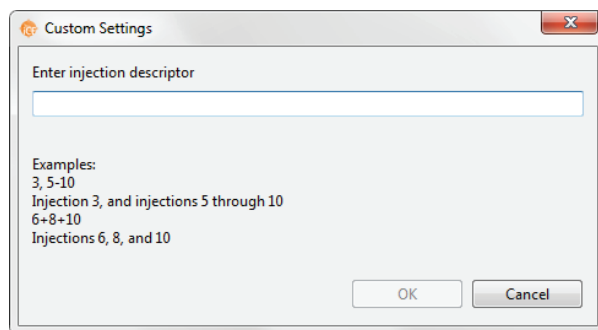


4. Click the cell in the **Apply To** column, then click the down arrow.

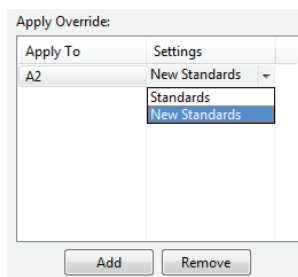


5. Select an option from the drop down list. This applies the settings group selected to specific run data as follows:

- **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
- **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.
- **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.
- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.

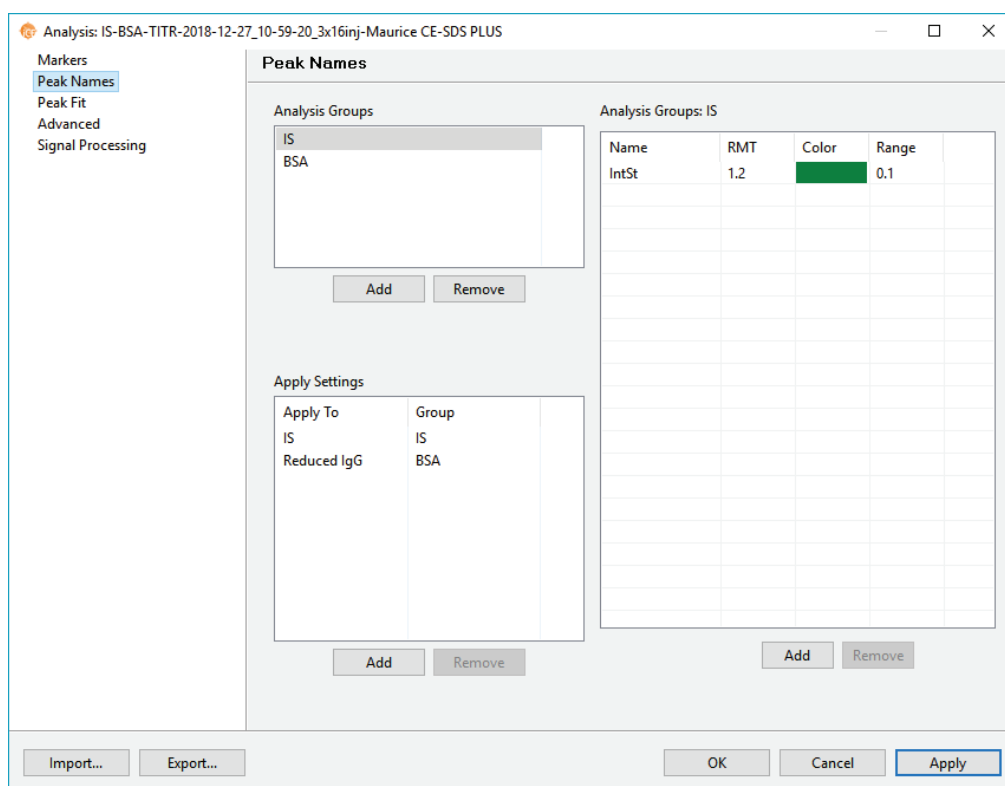


7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click **OK** to save changes.

Peak Names Settings

This page lets you view and change custom naming settings for sample proteins. Select **Edit** in the main menu and click **Analysis**, then click **Peak Names** in the left sidebar.

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.

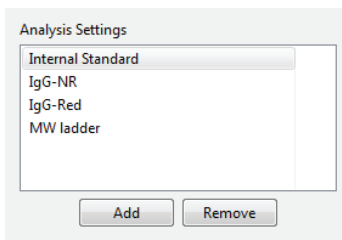


Peak Names Analysis Settings Groups

Peak name settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTE: Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 756.

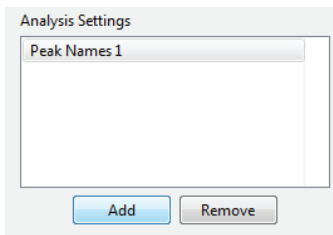
Peak name groups are displayed in the analysis settings box:



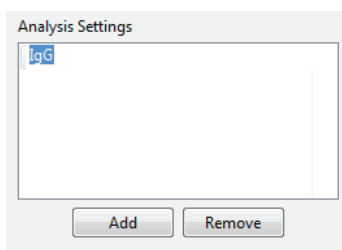
There aren't any Compass for iCE default settings groups, but you can make changes to groups you've created and create new groups. To view settings for a group, click on the group name in the analysis settings box.

Creating a Peak Names Group

1. Select **Edit > Analysis**, and select **Peak Names** in the left sidebar.
2. Click **Add** under the analysis settings box.



3. Enter a new name for the group.



4. Click in the first cell in the **Name** column in the analysis settings peak table and enter a sample protein name.

Analysis Settings: IgG

Name	RMT	Color	Range
HC	2		0.1

Add Remove

5. Click in the first cell in the **RMT** column and enter the relative migration time for the sample protein.

Analysis Settings: IgG

Name	RMT	Color	Range
HC	1.55		0.1

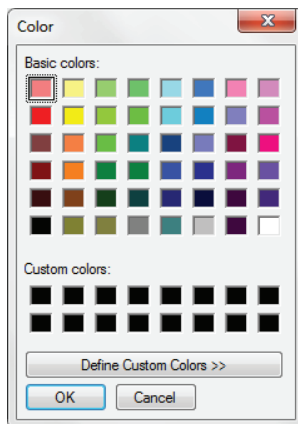
6. Click in the first cell in the **Color** column, then click the button.

Analysis Settings: IgG

Name	RMT	Color	Range
HC	1.55	■ (0.1...)	0.1

Line:

The color selection box displays:



7. The color you pick is used to identify the sample protein peak in the Peaks and Injections panes in the Analysis screen. Click a color or define a custom color and click OK. The color selection will update in the table:

Analysis Settings: IgG				
Name	RMT	Color	Range	
HC	1.55		0.1	

8. Click in the first cell in the **Range** column.

Analysis Groups: IgG				
Name	RMT	Color	Range	
HC	1.55		0.1	

9. Enter a \pm range for the RMT entered. Compass for iCE will automatically name peaks found within this RMT range. For example, if the RMT entered is 2 and a 0.1 range is used, all peaks with RMTs between 1.9 and 2.1 will be identified with this peak name and color.
10. To add another sample protein, click **Add** under the peak table. Repeat the previous steps for other sample proteins. In this example, three proteins were entered:

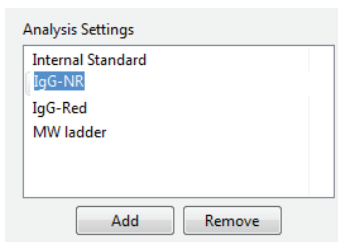
Analysis Groups: IgG				
Name	RMT	Color	Range	
HC	1.55		0.1	
NGHC	1.5		0.1	
LC	1.2		0.1	

To remove a sample protein, select its row and click **Remove**.

11. Click **OK** to save changes.

Modifying a Peak Names Group

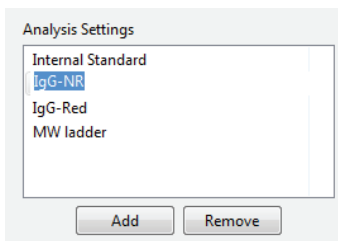
1. Select **Edit > Analysis**, then click **Peak Names** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



3. Change the information in the analysis settings peak table as described in [Creating a Peak Names Group](#).
4. Click **OK** to save changes.

Deleting a Peak Names Group

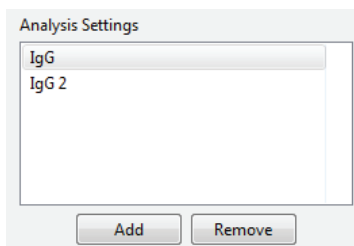
1. Select **Edit > Analysis**, then click **Peak Names** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



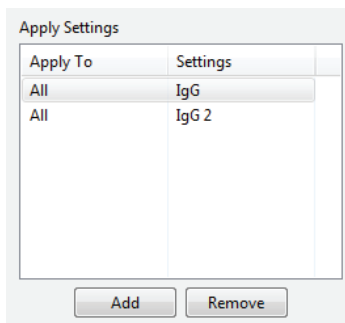
3. Click **OK** to save changes.

Applying Peak Names Groups to Run Data

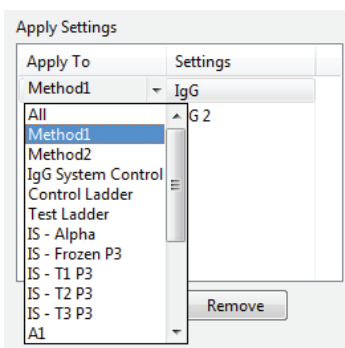
1. Select **Edit > Analysis**, then click **Peak Names** in the options list.
2. Click on the group in the analysis settings box you want to apply to specific run data.



3. Application of peak names groups to specific run data is done in the apply settings box. A default data set automatically gets created whenever you create a new group and it's applied to all injections in the run. You can either modify the default group or click **Add** under the box to create a new one.

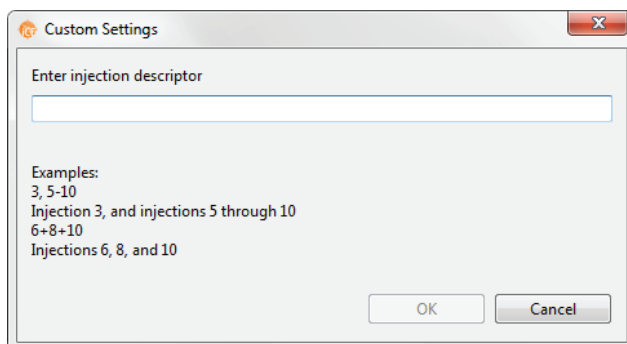


4. Click the cell in the **Apply To** column, then click the down arrow.

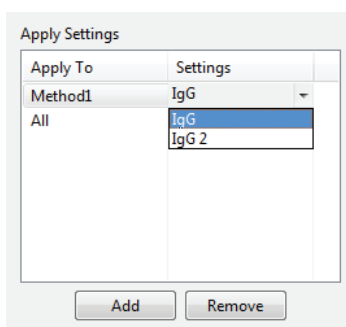


5. Select an option from the drop down list. This applies the peak names group selected to specific run data as follows:
- **All** - Selecting this applies peak names group settings to all injections.
 - **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.
 - **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.

- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:

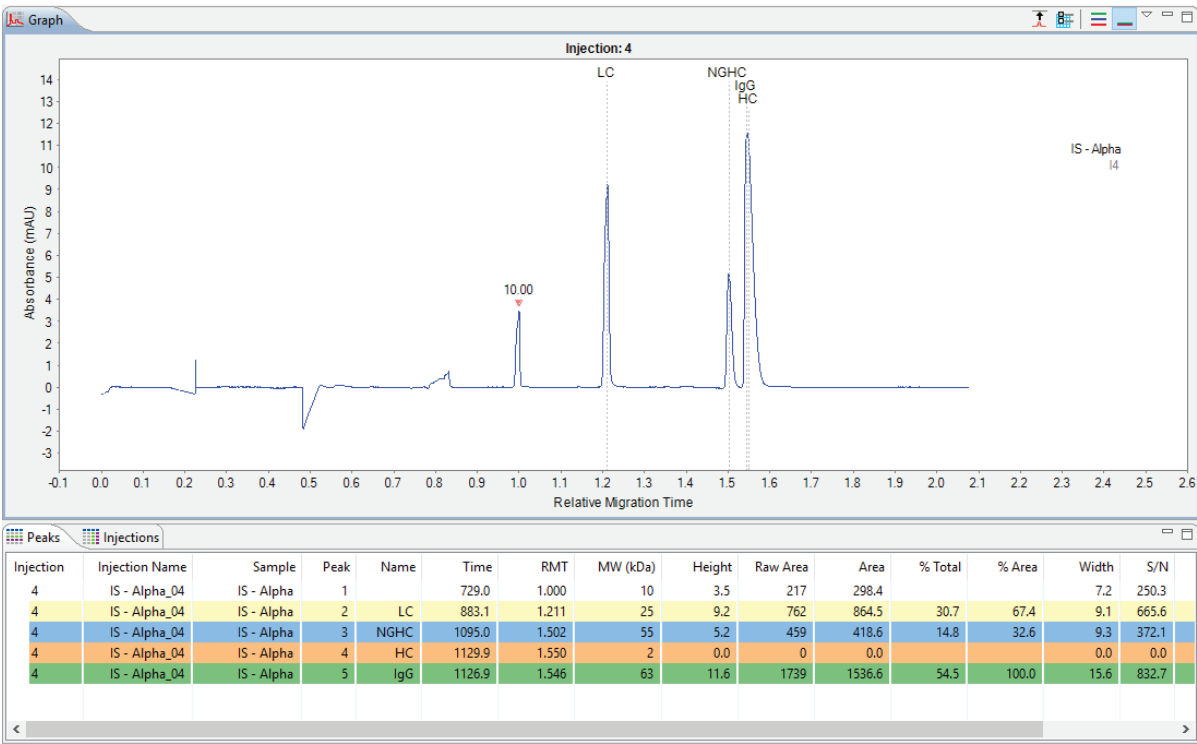


6. If you need to change the peak names group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.

9. Click OK to save changes. Named peaks will be identified with a peak name label in the electropherogram and color-coded in the Peaks and Injections panes:



Peak Fit Analysis Settings

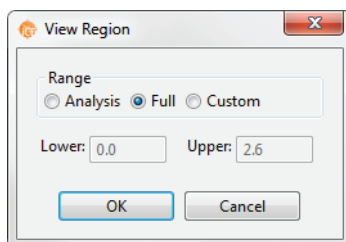
This page lets you view and change peak fit settings for sample data. Select **Edit** in the main menu and click **Analysis**, then click **Peak Fit** in the left sidebar:

NOTE: Settings can be changed in the batch default analysis before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.

Range Settings

- **Minimum** - The RMT value below which peaks won't be identified. This value is also used as the default lower RMT range for data displayed in the electropherogram.
- **Maximum** - The RMT value above which peaks won't be identified. This value is also used as the default upper RMT range for data displayed in the electropherogram.

- **View** - Sets the default range to either Full or Analysis for the electropherogram x-axis range in the View Region window (select **View** in the main menu and click **View Region**).



- **Analysis** sets the x-axis range of the electropherogram to the Peak Fit range minimum and maximum settings in the electropherogram.
- **Full** displays the entire separation range of the run data in the electropherogram. This is the default setting.

Baseline Settings

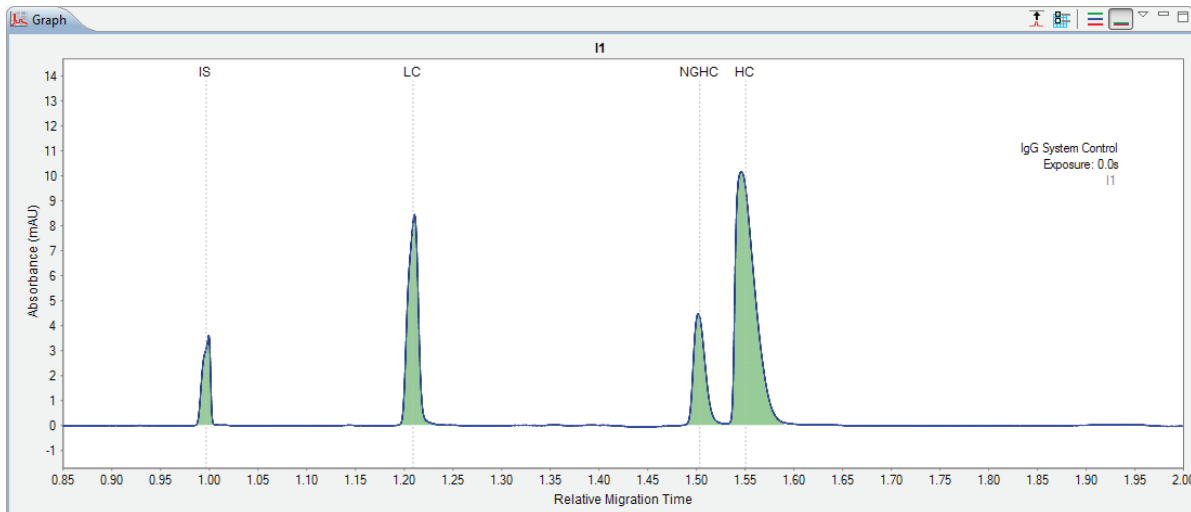
These settings apply to spline baselines only.

- **Threshold** - The variance, or roughness, in a baseline data segment below which a point is called part of the baseline.
- **Window** - How long baseline data segments are expected to be in pixels. Shorter segments let the baseline follow plateau sections of the signal.
- **Stiffness** - The amount the baseline is allowed to vary from a straight line. Settings between 0.1 and 1.0 make the baseline fit closer to a straight line. Settings from 1.0 to 10.0 will make the baseline fit follow the data more closely.

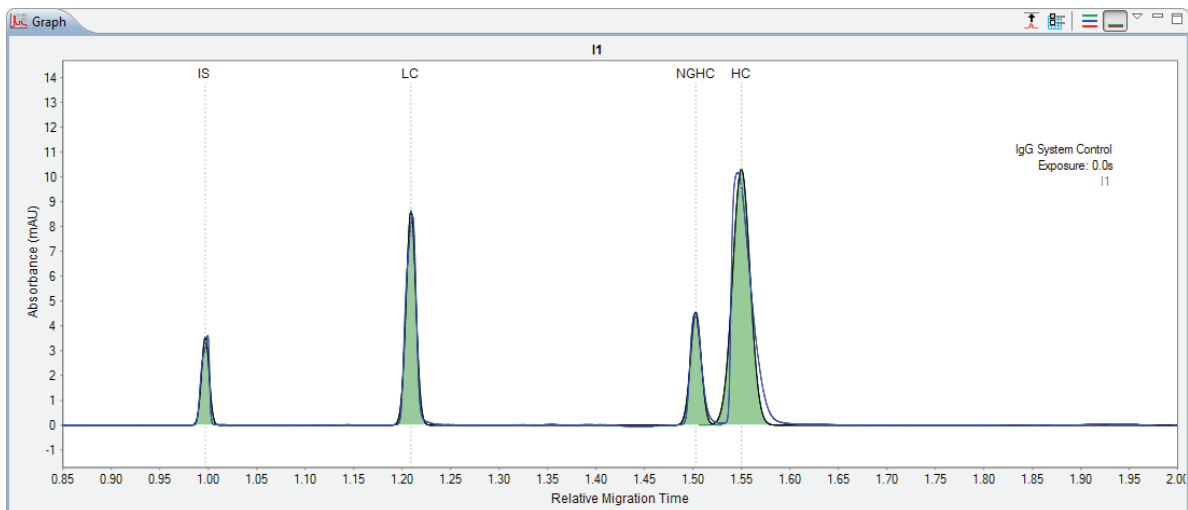
Peak Find Settings

- **Threshold** - The minimum signal to noise ratio required for a peak to be identified. A setting of 1.0 will detect many peaks, a setting of 10.0 will detect fewer peaks.
- **Width** - The approximate peak width (at full width half max) in pixels used to detect peaks. The minimum value for this setting is 3.0. Larger widths help eliminate the detection of shoulder and noise peaks.
- **Area Calculation** - Two fits are used, either Gaussian Fit or Dropped Lines. These settings can be changed before or after the run is finished.

- For CE-SDS applications, peak area is calculated using Dropped Lines by default. This type of area calculation is also often called the perpendicular drop method. This is the preferred method when peaks overlap or are close to each other. It draws two vertical lines from the left and right bounds of the peak down to the x-axis and then measures the total area bounded by the signal curve, the x-axis ($y=0$ line), and the two vertical lines.



- This next view is of the same data using Gaussian fit instead:



Peak Fit Analysis Settings Groups

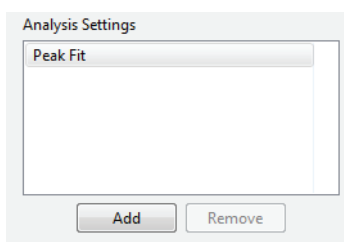
Peak fit settings are saved as a group, and you can create multiple settings groups. Specific group settings can then be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values for peak fit analysis settings. These settings are included in the default Peak Fit group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. For more information see “Importing and Exporting Analysis Settings” on page 756.

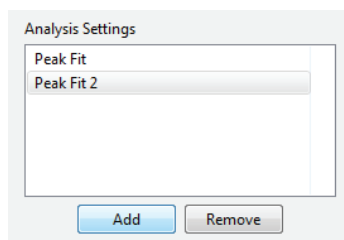
Peak fit groups are displayed in the analysis settings box:



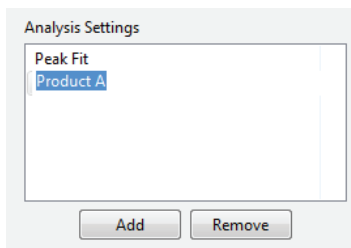
The Peak Fit group shown contains the Compass for iCE default analysis settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Peak Fit Group

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



- Click on the new group and enter a new name.



- Change the settings in the range, baseline or peak find boxes as needed.

Analysis Settings: Product A

Range

Minimum 0.9

Maximum 2.5

View ☐ Analysis ☒ Full

Baseline

Threshold 2.0

Window 100.0

Stiffness 0.7

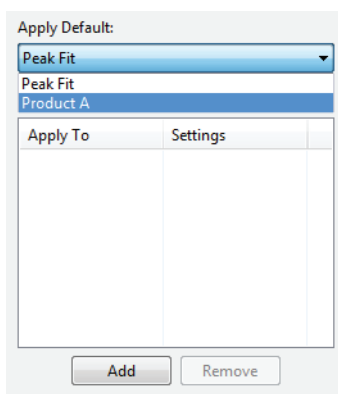
Peak Find

Threshold 15.0

Width 150.0

Area Calculation **Dropped Lines**

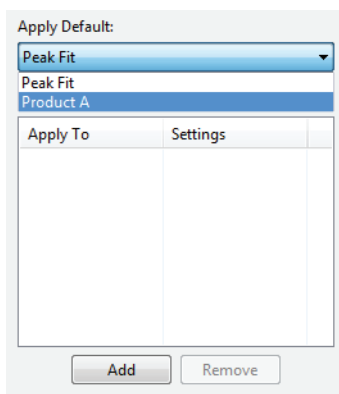
- To use the new group as the default peak fit settings for the run file data, click the arrow in the drop down list next to Apply Default, then click the new group from the list. Peak fit settings in the new group will then be applied to the run data.



- Click **OK** to save changes.

Changing the Default Peak Fit Group

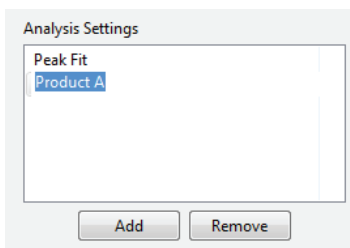
1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click the arrow in the drop down list next to **Apply Default**, then click a new default group from the list.



3. Click **OK** to save changes. Peak fit settings in the group selected will be applied to the run data.

Modifying a Peak Fit Group

1. Select **Edit > Analysis**, and click **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



3. Change the settings in the range, baseline or peak find boxes as needed.

Analysis Settings: Product A

Range

Minimum 0.9

Maximum 2.5

View ☐ Analysis ☒ Full

Baseline

Threshold 2.0

Window 100.0

Stiffness 0.7

Peak Find

Threshold 15.0

Width 150.0

Area Calculation Dropped Lines ▼

4. Click OK to save changes. The new peak fit settings will be applied to the run data.

Deleting a Peak Fit Group

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.

Analysis Settings

Peak Fit

Product A

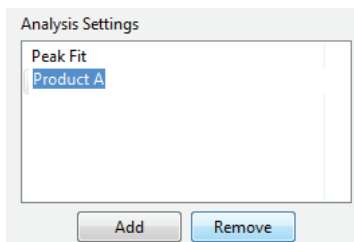
Add Remove

3. Click OK to save changes.

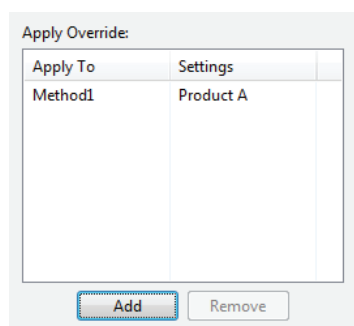
Applying Peak Fit Groups to Specific Run Data

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.

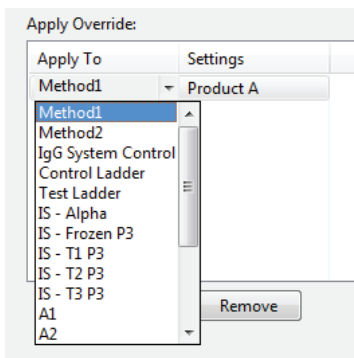
- Click on the group in the analysis settings box you want to apply to specific run data.



- Application of analysis groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

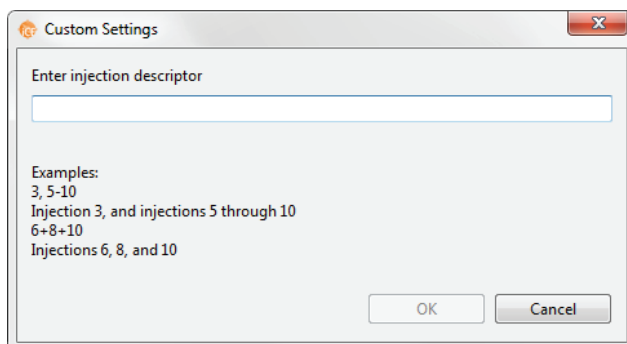


- Click the cell in the **Apply To** column, then click the down arrow.

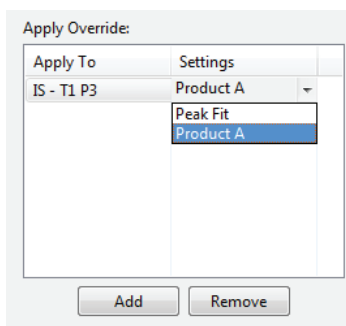


- Select an option from the drop down list. This applies the settings group selected to specific run data as follows:
 - Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
 - Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.
 - Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.

- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.

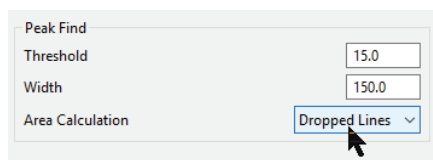


7. Repeat the previous steps to apply other groups to specific run data.
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.
9. Click OK to save changes.

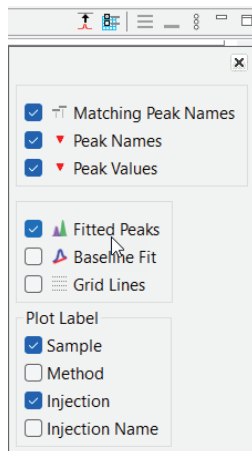
Manual Peak Integration

Compass for iCE lets you manually integrate peaks in individual electropherograms.

1. Select **Edit > Analysis**, and select **Peak Fit** in the left sidebar.
2. Select **Dropped lines** as the area calculation if it isn't already selected.

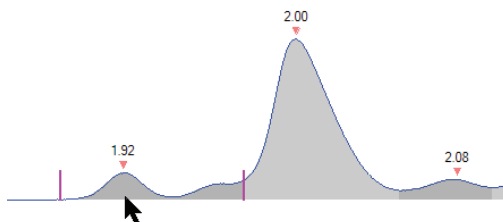


3. Select **Fitted Peaks** in the Graph Options.



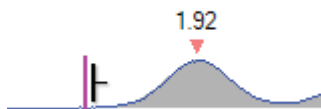
4. Select an injection in the Experiment pane.

5. Hover the mouse over the peak you want to adjust. Two magenta bars will display that indicate the start and end points of its integration.

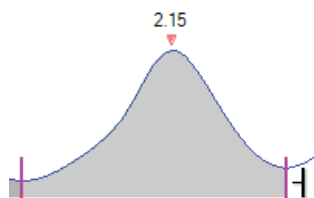


6. To change the integration, hover the mouse over either of the magenta bars. Then click the mouse and drag the bar to move it. The cursor changes to indicate which direction you can move the integration point:

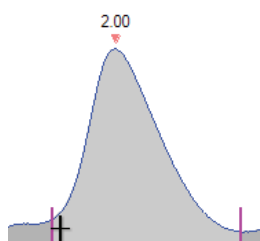
- If the cursor changes to  this is the peak start for the peak on the right.



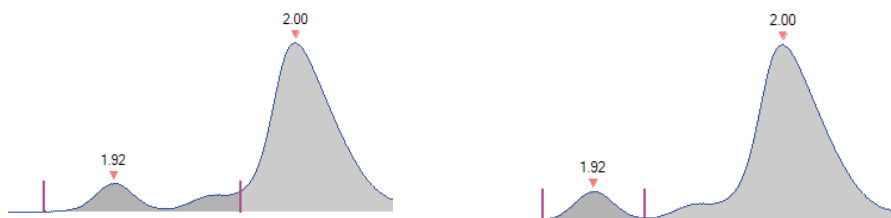
If the cursor changes to  this is the peak end for the peak on the left.



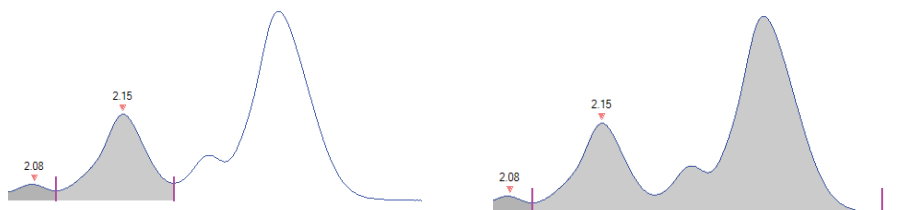
- If the cursor changes to  this is a joint boundary for the peaks on the left and right.






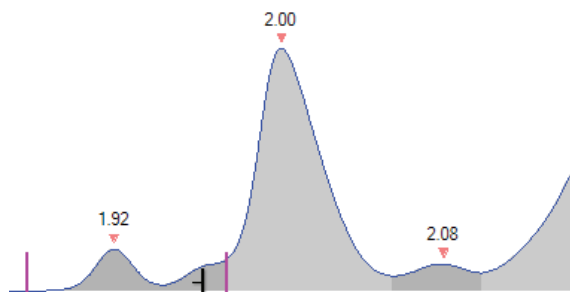
In the example below, we moved the start and end points of the peak at 1.92 RMT to just include the area under the peak:



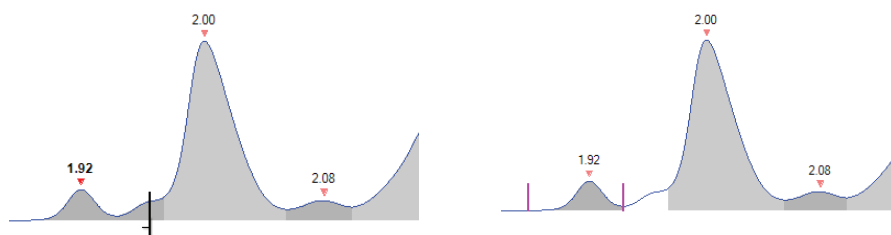
You can also combine peaks:



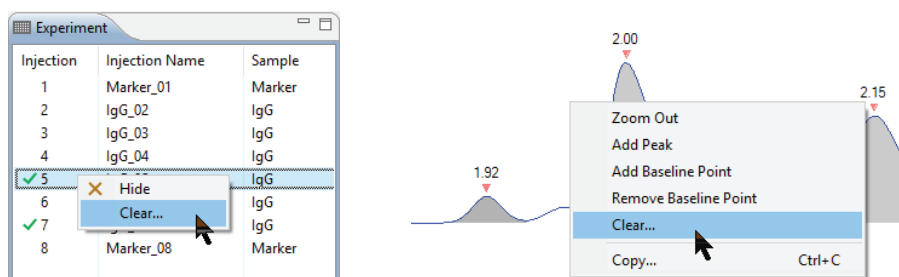
You can also separate areas between peaks. Whenever you have a  cursor between two peaks that aren't baseline resolved, move the mouse slightly to the right or left until you get the  or  cursor.



Click and hold the mouse. A black drop line will display indicating where the current peak start or end point is. While holding the mouse, drag the cursor to the right or left to move the start or end point to where you want it, then release the mouse.



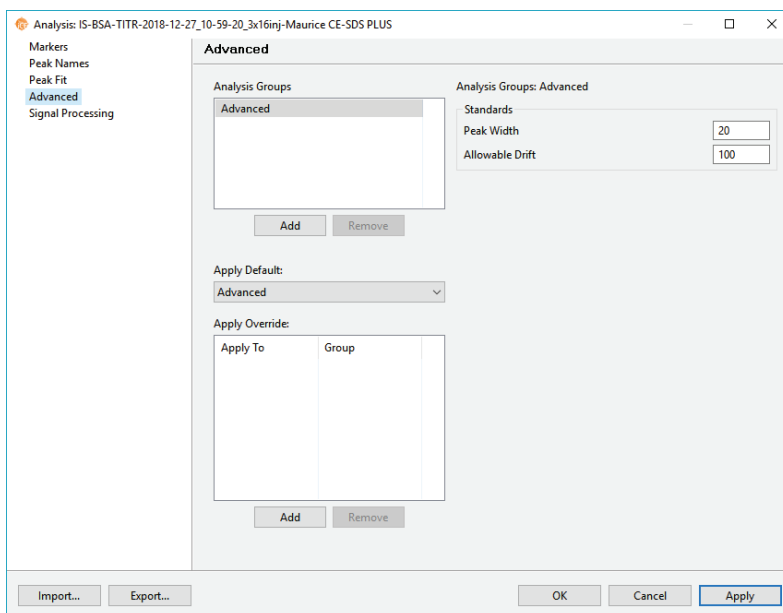
7. To remove the manual integration settings, right click on the graph or on the injection in the Experiment pane and select **Clear**.



Advanced Analysis Settings

This page lets you view and change analysis settings for the Internal Standard data. Select **Edit** in the main menu and click **Analysis**, then click **Advanced** in the left sidebar:

NOTE: Settings can be changed in batches before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



Internal Standard Settings

- **Peak Width** - The approximate width (at full width half max) used to filter out absorbance artifacts which improves recognition of standards.
- **Allowable Drift** - The distance the Internal Standard is expected to move compared to the entered number of seconds on the Markers page. This setting helps with recognition of the Internal Standard.

Advanced Analysis Settings Groups

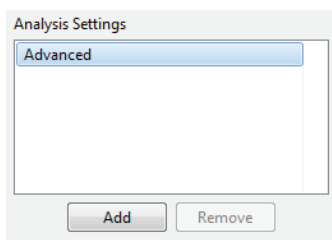
Advanced analysis settings are saved as a group, and you can create multiple settings groups. Specific group settings can be applied to methods, injections, sample names or other attributes in the run data.

NOTES:

We recommend using the Compass for iCE default values for advanced analysis settings. These settings are included in the default Advanced group.

Analysis settings are run-file specific. But, settings can be imported or exported for use with other run files. See “Importing and Exporting Analysis Settings” on page 756.

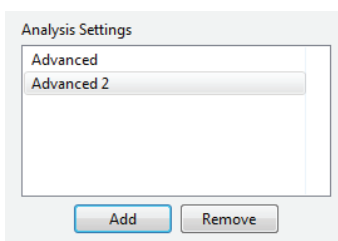
Analysis groups are displayed in the analysis settings box:



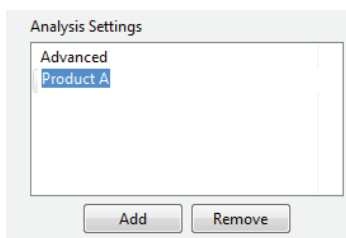
The Advanced group shown contains the Compass for iCE default analysis settings. You can make changes to this group and create new groups. To view settings for a group, click on the group name.

Creating a New Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click **Add** under the analysis settings box. A new group will be created:



3. Click on the new group and enter a new name.



4. Change the settings in the Standards box as needed.

Analysis Settings: Product A

Standards

Peak Width	10
Allowable Drift	50

5. To use the new group as the default analysis settings for the run data, click the arrow in the drop down list next to Apply Default, then click the new group from the list. Analysis settings in the new group will then be applied to the run data.

Apply Default:

Advanced

Advanced

Product A

Apply To	Settings
----------	----------

Add Remove

6. Click OK to save changes.

Changing the Default Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click the arrow in the drop down list next to Apply Default, then click a new default group from the list.

Apply Default:

Advanced

Advanced

Product A

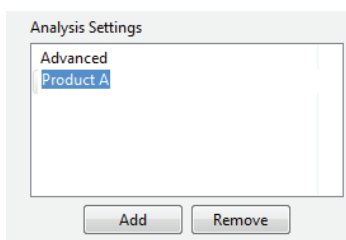
Apply To	Settings
----------	----------

Add Remove

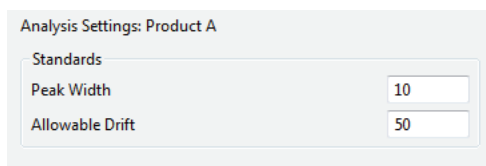
3. Click **OK** to save changes. Analysis settings in the group selected will be applied to the run data.

Modifying an Analysis Group

1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to modify.



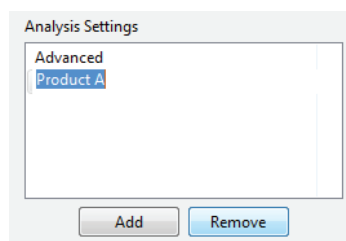
3. Change the settings in the Standards box as needed.



4. Click **OK** to save changes. The new analysis settings will be applied to the run data.

Deleting an Analysis Group

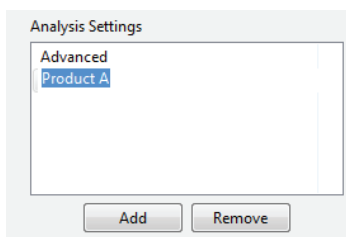
1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to delete and click **Remove**.



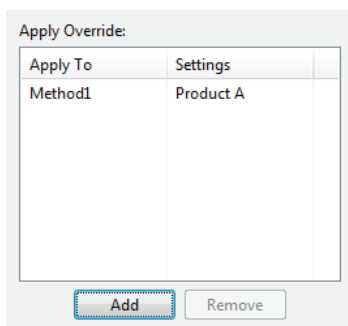
3. Click **OK** to save changes.

Applying Analysis Groups to Specific Run Data

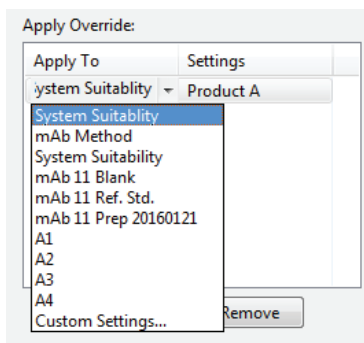
1. Select **Edit > Analysis**, and select **Advanced** in the left sidebar.
2. Click on the group in the analysis settings box you want to apply to specific run data.



3. Application of analysis groups to specific run data is done in the override box. Click **Add** under the override box. A default override data set will be created from sample information found in the run file.

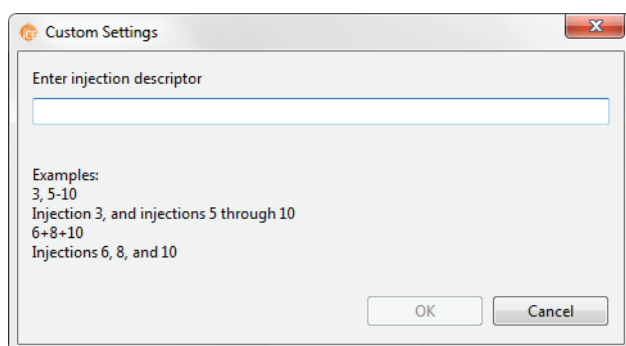


4. Click the cell in the **Apply To** column, then click the down arrow.

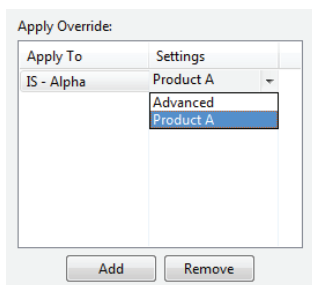


5. Select an option from the drop down list. This applies the settings group selected to specific run data as follows:

- **Methods** - All methods in the run file display in the list. Selecting a method applies the group settings to all injections that used that method.
- **Sample names** - All sample names in the run file display in the list, otherwise the default name of Sample shows. Selecting a sample name applies the group settings to all injections that used that sample name.
- **Wells or vials** - All well or vial numbers used in the run display in the list. Selecting a well/vial number applies the group settings to all injections that used that well/vial.
- **Custom settings** - Lets you choose specific injections to apply the group settings to. When you select this in the list, a pop-up box displays to let you enter a specific injection number or range of injections:



6. If you need to change the analysis group used for a data set, click the cell in the **Settings** column and click the down arrow. Select a group from the drop down list.



7. Repeat the previous steps to apply other groups to specific run data.

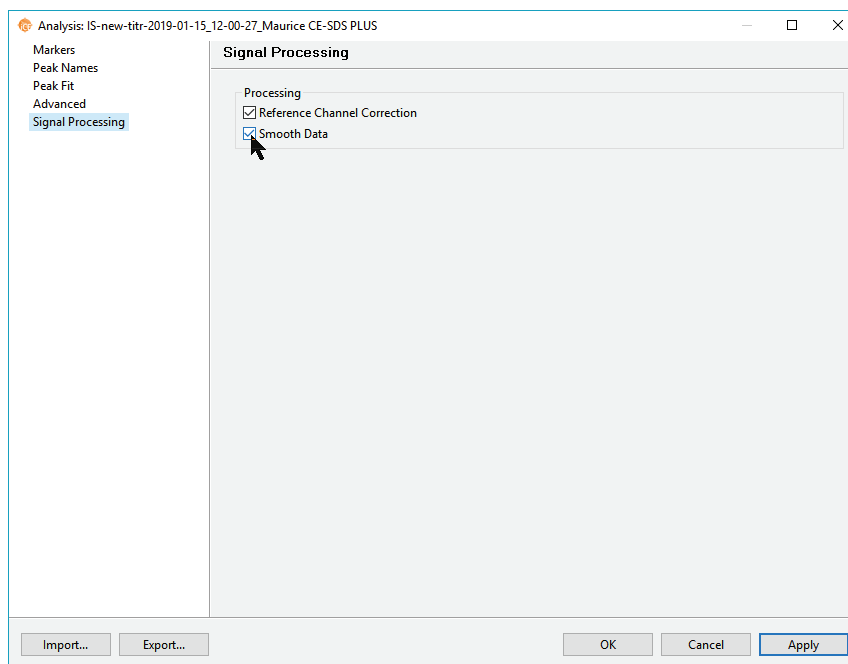
8. To remove a data set, click on its cell in the **Apply To** column, then click **Remove**.

9. Click **OK** to save changes.

Signal Processing Settings

This page lets you apply reference signal corrections and baseline smoothing options to your data. Select **Edit** in the main menu and click **Analysis**, then click **Signal Processing** in the left sidebar:

NOTE: Settings can be changed in batches before you start the run, or in run files once they're completed. If you make analysis settings changes to an executing run, they won't be saved to the final run file.



Reference Channel Correction

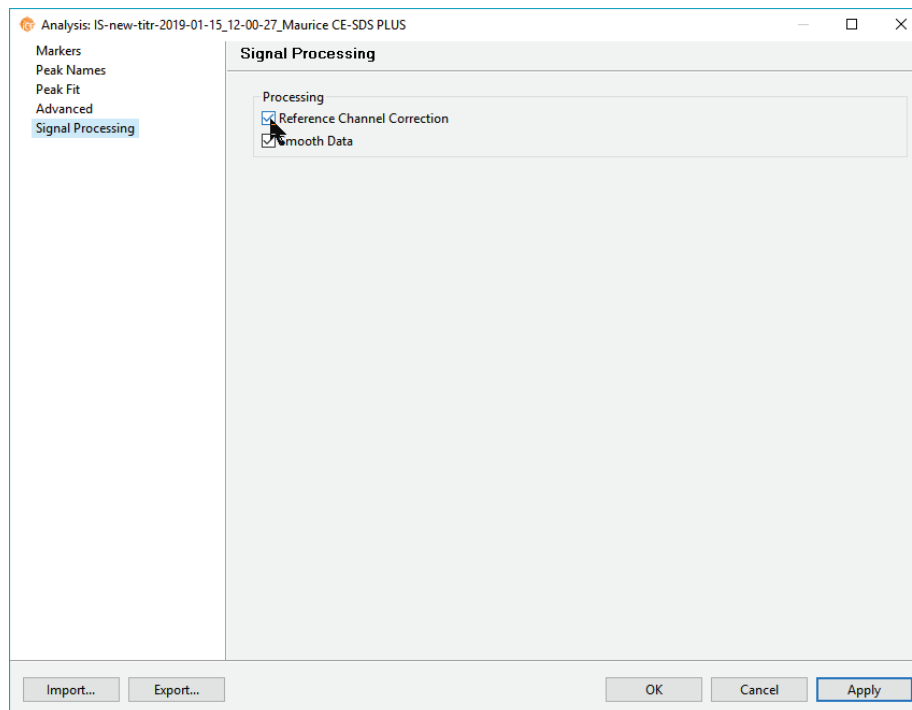
The Reference Channel Correction improves CE-SDS data quality by automatically removing any signal within a certain range that can be attributed to lamp variability. It's on by default for new CE-SDS PLUS and Turbo CE-SDS batches and runs in Compass for iCE 2.1 or higher, and off when opening runs collected in previous versions of the software. It can be turned on when analyzing older runs if preferred.

NOTE: We recommend keeping the Reference Channel Correction on for the highest data quality. For more information, contact Protein Simple Technical Support.

To turn reference channel correction on or off:

1. From the Batch screen, select **Edit > Default Analysis**. From the Analysis screen, select **Edit > Analysis**.
2. Click **Signal Processing** in the left sidebar.

3. Select or deselect the **Reference Channel Correction** checkbox.



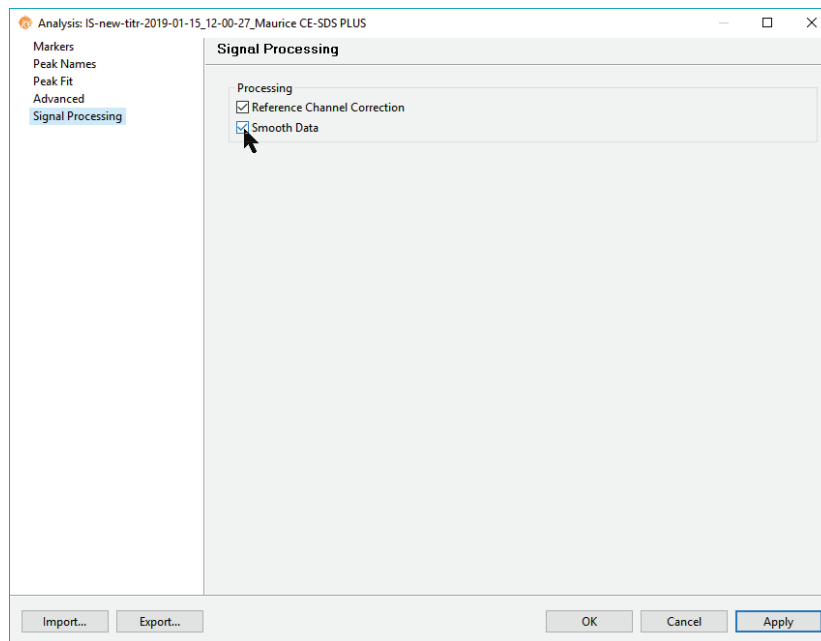
4. Click **Apply**.

Baseline Smoothing

Compass for iCE can apply baseline smoothing to CE-SDS data to decrease noise and improve signal-to-noise ratios. Smoothing is on by default but can be turned off in the Default Analysis Settings before you start a batch or during analysis after a batch is completed.

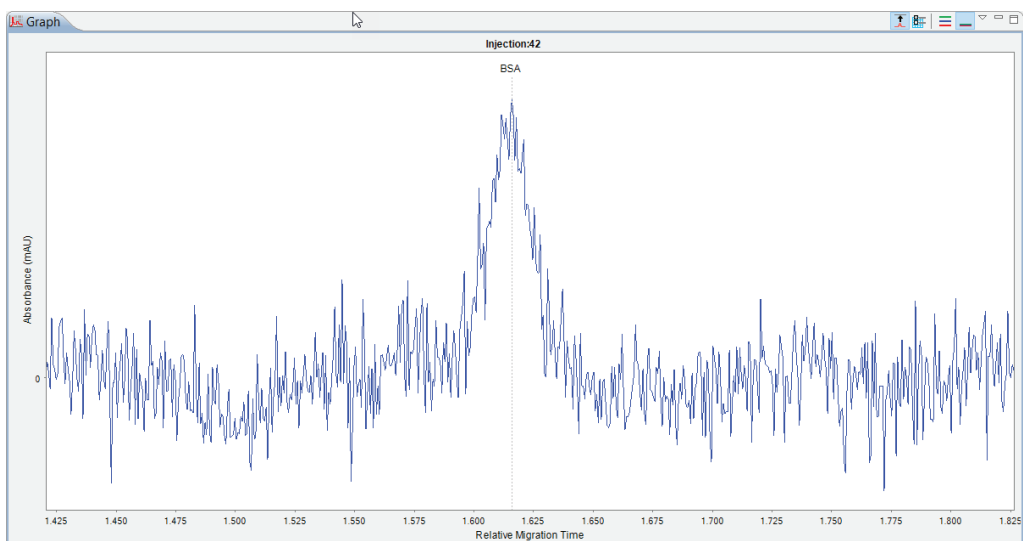
To turn baseline smoothing on or off:

1. From the Batch screen, select **Edit > Default Analysis**. From the Analysis screen, select **Edit > Analysis**.
2. Click **Signal Processing** in the left sidebar.

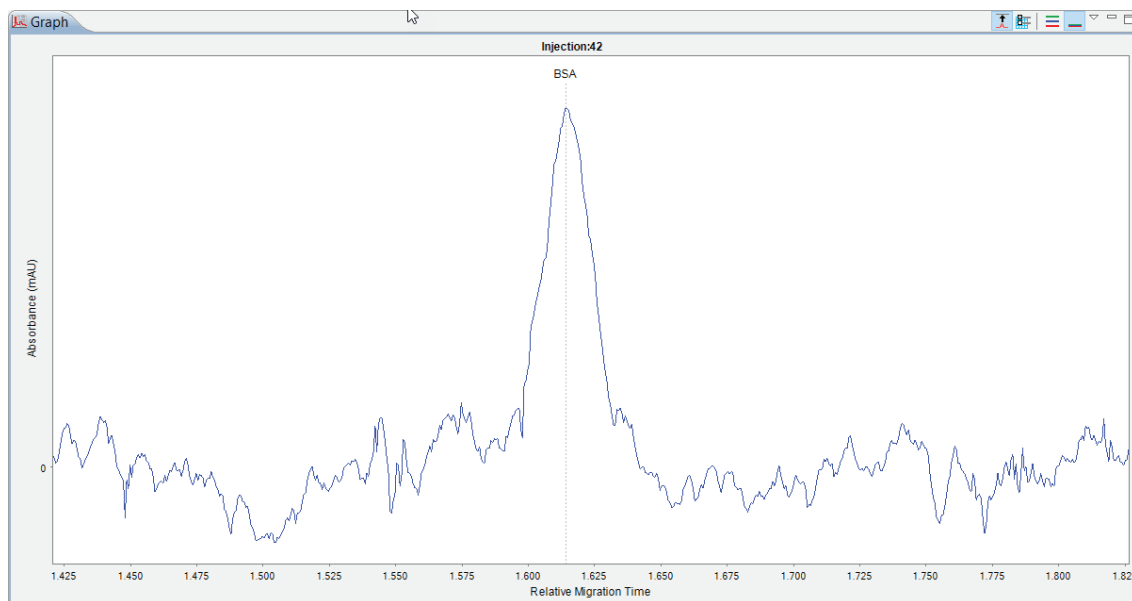
3. Select or deselect the **Smooth Data** checkbox.4. Click **Apply**.

The examples below show data with and without the Smooth Data option:

- Smooth Data off



- Smooth Data on



Injection Reports

You can export PDF files of the raw and analyzed data, IV plot, peaks table, sample and system info for individual or all injections in a run file. You can also export the run history with all analysis events.

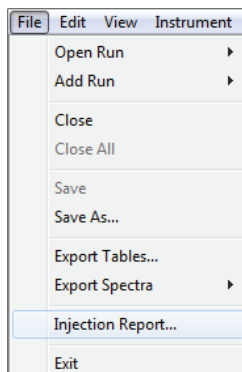
NOTES:

You can have Compass for iCE generate reports for you automatically at the end of each run. See “Setting up Automatic Injection Reports” on page 761 for more information.

Any time a new report is run, it’s saved in a separate run folder. Existing reports aren’t overwritten.

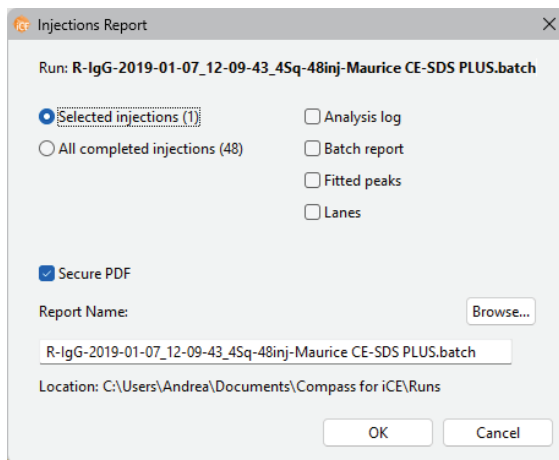
1. Click **File > Open Run** and select a run file.
2. If you want reports for all injections, skip to the next step. If you only want reports for certain injections, in the Experiment pane:
 - **To select sequential injections:** Select the first injection, then hold the **Shift** key and select the last injection you want a report for. This selects all rows between the two injections.
 - **To select specific injections:** Hold the **Ctrl** key and select just the injections you want reports for.

3. Select **File** from the main menu in either screen and click **Injection Report**.

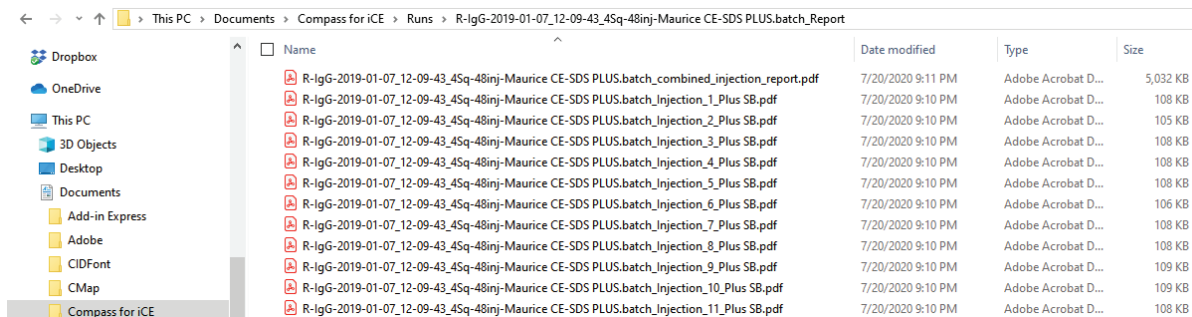


4. In the Injection Reports window:

- a. Choose either **Selected injections** or **All injections**.
- b. Select **Analysis log** if you want a run history report with all analysis events.
- c. Select **Batch Report** if you want to include the sample and method details for each injection in the batch.
- d. Select **Fitted peaks** if you want to show peak fitting in the electropherograms.
- e. Select **Lanes** if you want to show lane view data.
- f. Report PDFs generated by Compass for iCE are secured by default, which means they can be viewed and printed but not modified or renamed. Uncheck **Secure PDF** if you don't want to generate a secure report. To learn more about permissions for secured vs. unsecured reports, see page 800.
- g. The report name defaults to the run file name. If you want to change it, type in the **Report Name** box to make updates.
- h. Click **OK**.



5. The individual injection report PDF(s) and a combined injection report PDF are exported to the Runs folder in the Compass for iCE directory. They'll be in a folder with the report name used in the prior step. When the reports are done, the folder opens for you automatically.



Example Analysis and Injection Report

Run File R-IgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS

Analysis Log

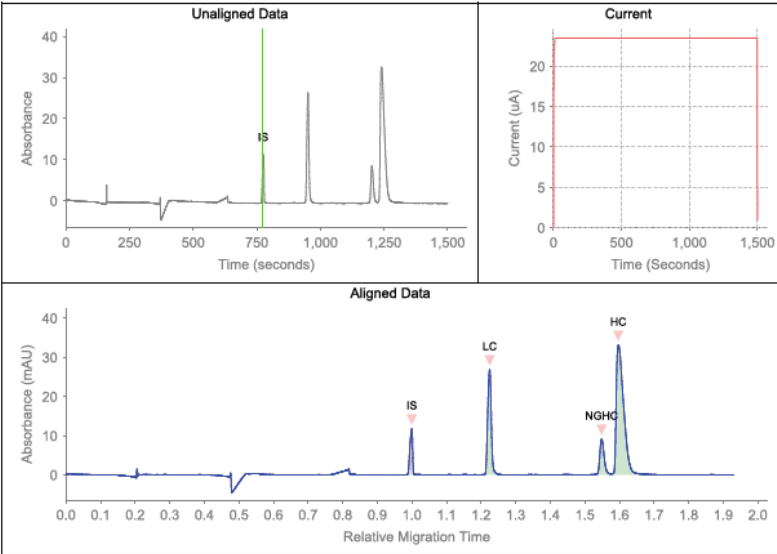
Date	User Name	Message	Comment
2019-01-10 17:25:27		Started run: R-IgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS Batch: 4Sqx12inj-RlgG-Maurice CE-SDS.batch	3180910338
2019-01-14 10:39:07		Saved analysis and methods changes from Compass for iCE v2.1.0-1219	
		Added Peak Names Apply Settings "apply IgG to all"	
		Added Peak Names Group IgG	
		Protein name: LC RMT: 1.21 Color: 32512 Range: 0.1	
		Protein name: NGHRC RMT: 1.54 Color: 32512 Range: 0.05	
		Protein name: HC RMT: 1.58 Color: 32512 Range: 0.1	

Created By: Jacquelyn Sat 4:47 PM Feb 2, 2019 CST Software Version: 2.1.0 Build ID: 0130
 C:\Users\Jacquelyn\Documents\Clients\ProteinSimple\Maurice\User Guide\Rev 12 edits\Data from Andrea - no edits\R-IgG-2019-01-10_16-25-12_4Sqx12inj-RlgG-Maurice CE-SDS.mlx
 Computer: DESKTOP-C7FPOGB

Page 1 of 2



Uncontrolled Injection 1: R IgG1



Peaks											
Peak	Name	Time	RMT	Height	Raw Area	Area	%Total	%Area	Width	S/N	Baseline Resolution
1	IS	776.7	1.000	11.7	164.6	212.4			7.0	658.6	-0.6
2	LC	951.3	1.225	26.8	498.5	524.1	30.5	30.5	8.8	1507.3	-0.6
3	NGHC	1202.7	1.549	9.1	205.6	170.8	9.9	9.9	10.2	510.5	-0.6
4	HC	1240.2	1.597	33.1	1273.4	1021.8	59.5	59.5	17.9	1859.6	-0.6

Uncontrolled Injection 1: R IgG1

Sample Information

Injection Name	R IgG1
Sample ID	SB
Location	Plate Well B2
Batch Name	4Sgx12inj-RlgG-Maurice CE-SDS
Run Started	Thu 5:25 PM Jan 10, 2019 CST
Run Completed	Fri 8:07 PM Jan 11, 2019 CST
Date Acquired	Thu 6:08 PM Jan 10, 2019 CST
Run Error	None
Reinjection	No

Injection Conditions

Guaranteed Injection	Yes
Focus Period 1	5750V for 25.0 min
Sample Load	20 sec 4600 Volts
Tray Temperature	9.1°C

Maurice Settings

Model	Maurice OBM
Instrument S/N	kf1077
Software Version	2.1.0, Build ID: 0130
Firmware Version	3.1.2019.01.04.20.16.20.5eb707f
Tray Type	96-well plate
Cartridge Type	CE-SDS
Cartridge S/N	3180910338
Cartridge Expiration	Sep 2019
Injections Remaining	152 (52 guaranteed)
Batches Remaining	24

Created By: jacquelyn Bat 4:47 PM Feb 2, 2019 CST Software Version: 2.1.0, Build ID: 0130
 C:\Users\jacquelyn\Documents\Clients\ProteinSimple\Maurice\User Guide\Rev 12 edits\Data from Andrea - no edits\RlgG-2019-01-10_16-25-12_4Sgx12inj-RlgG-Maurice CE-SDS.mz
 Computer: DESKTOP-C7FPQGB

Page 2 of 2



Importing and Exporting Analysis Settings

The analysis settings in a run file can be exported as a separate file. This allows the same analysis settings to be imported into other batches or run files at a later time, rather than you having to re-enter them manually.

Importing Analysis Settings

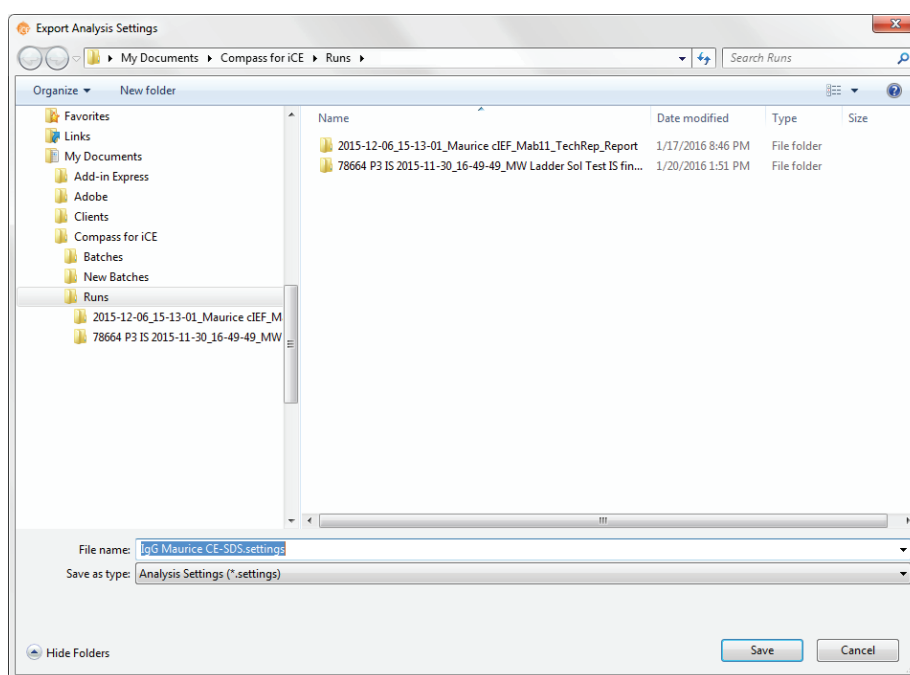
NOTE: Importing an analysis settings file populates the settings in all analysis pages.

1. Open the run file or batch you want to import analysis settings to.
2. Select **Edit** in the main menu and click **Default Analysis** (Batch screen) or **Analysis** (Analysis screen).
3. Click **Import** on any page.
4. Select a settings file (*.settings) and click **OK**. The imported settings will display in all analysis pages.

Exporting Analysis Settings

NOTE: Exporting an analysis settings file exports the settings in all analysis pages.

1. Open the run file or batch you want to export analysis settings from.
2. Select **Edit** in the main menu and click **Default Analysis** (Batch screen) or **Analysis** (Analysis screen).
3. Click **Export** on any page. The following window displays:



4. The default directory is Compass for iCE/Runs. Change the directory if needed.
5. Enter a file name and click **Save**. The settings will be saved as a *.settings file.

Chapter 21:

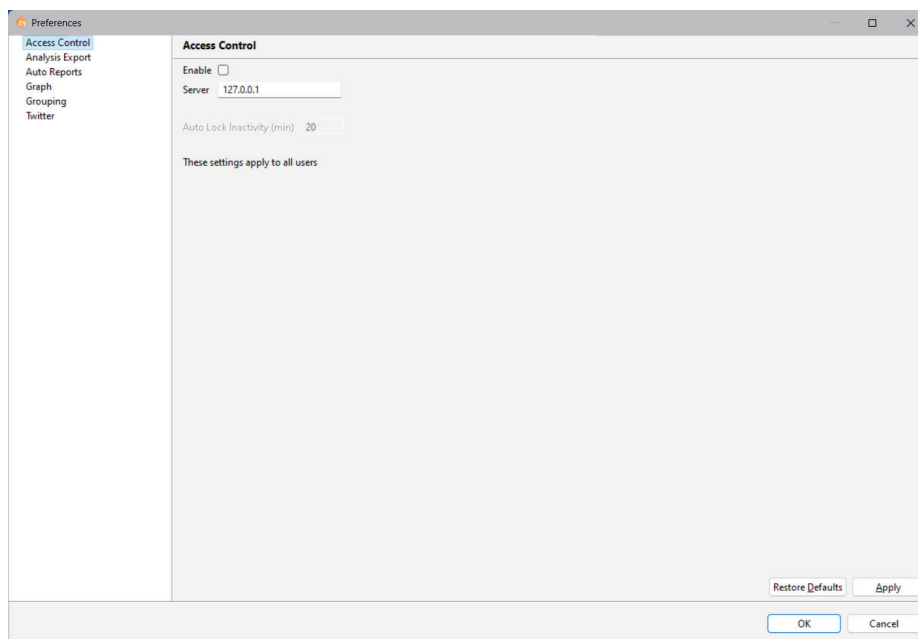
Setting Your Preferences

Chapter Overview

- Customize Your Preferences
- Enabling Access Control
- Setting Data Export Options
- Setting up Automatic Injection Reports
- Selecting Custom Plot Colors for Graph Overlay
- Grouping Options
- Setting Up Maurice Systems to Send Tweets

Customize Your Preferences

You can set and save several custom preferences in Compass for iCE. To view and change these settings, select **Edit** in the main menu and click **Preferences**.



To move between preferences pages, click on an option in the left sidebar. Here's what you can customize:

- **Access Control** - Lets you log on to Compass for iCE through an Authorization Server.
- **Analysis Export** - Lets you customize data export options.
- **Auto Reports** - Lets you set reports to generate automatically at the end of a run.
- **Graph** - Lets you customize graph color displays.
- **Grouping** - Groups samples with the same name together across runs, so you can get statistics for the same sample in multiple runs.
- **Twitter** - Lets you configure Compass for iCE to tweet Maurice, Maurice C., Maurice S. and MauriceFlex. run status.

In all preferences windows:

- Click **Apply** to apply changes to any open run files in Compass for iCE.
- Click **Restore Defaults** to restore the values on the page to default settings.
- Click **OK** to save changes and exit.
- Click **Cancel** to exit without saving changes.

Enabling Access Control

You can use the Access Control feature to help satisfy 21CFR Part 11 data security requirements when using Maurice instruments. Please go to “Enabling Access Control” on page 781 to get more info.

Setting Data Export Options

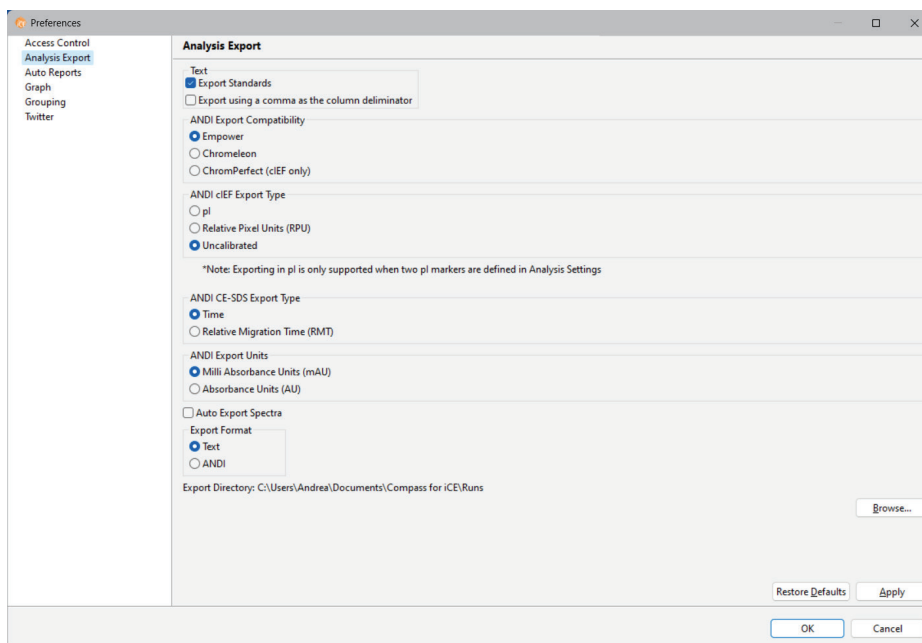
Select **Analysis Export** in the sidebar.

- **Text:**
 - **Export Standards** - This option exports data for the standards in each injection when run data is exported. It's selected by default. If it's not selected, only sample injection data is exported.
 - **Export using a comma as the column delimiter** - This option exports run data with a comma separator in .csv format. When it's not selected, data is exported in .txt format with a tab separator (this is the default setting).
- **ANDI Export Compatibility:**
 - **Empower** - This option exports run data in a format that is compatible for further analysis in Empower and other analysis programs that use standard formats.
 - **Chromeleon** - This option exports run data in a format that is compatible for further analysis in Chromeleon.
 - **ChromPerfect (cIEF only)** - This option exports standard cIEF, MauriceFlex cIEF and MauriceFlex Fractionation run data in a format that is compatible for further analysis in ChromPerfect.
- **ANDI cIEF Export Type:**
 - **pI** - Exports the pI of identified peaks in the cIEF run data.
 - **Relative Pixel Units (RPU)** - Exports the RPU of identified peaks in the cIEF run data.
 - **Uncalibrated** - Exports uncalibrated cIEF run data.

NOTE: ANDI spectra data for standard cIEF, MauriceFlex cIEF and MauriceFlex Fractionation can be exported. For MauriceFlex Fractionation, only sample injection data can be exported. Mobilization electropherograms can not be exported.

- **ANDI CE-SDS Export Type:**
 - **Time** - Exports the peak detection time
 - **Relative Migration Time (RMT)** - Exports the RMT of identified peaks in the CE-SDS run data.
- **ANDI Export Units:**
 - **Milli Absorbance Units** - Exports peak height in milli absorbance units (mAU).
 - **Absorbance Units** - Exports peak height in absorbance units (AU).

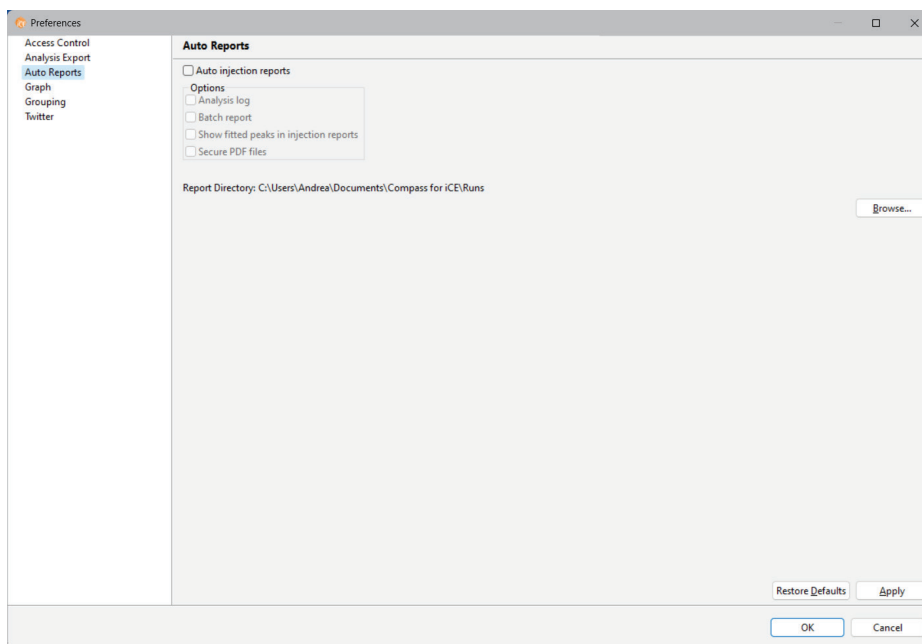
- Auto Export Spectra:
 - When Text or ANDI options are selected, spectra will automatically export at the end of every run. Data is saved in \Documents\Compass for iCE\Runs by default. To save the files in another folder, click Browse and select a different location.
 - **Text** - Exports spectra in text format.
 - **ANDI** - Exports spectra in ANDI format.



Setting up Automatic Injection Reports

1. Select **Auto Reports** in the sidebar.
2. Check the **Auto Injection Reports** checkbox.
3. Check any additional reports you'd like to have automatically generated at the end of each run
4. Check **Secure PDF** to have the software automatically generate secure report PDFs that can be viewed and printed but not modified or renamed.

Reports include all injections in the batch, and are saved in \Documents\Compass for iCE\Runs by default. To save reports in another folder, click Browse and select a different location.

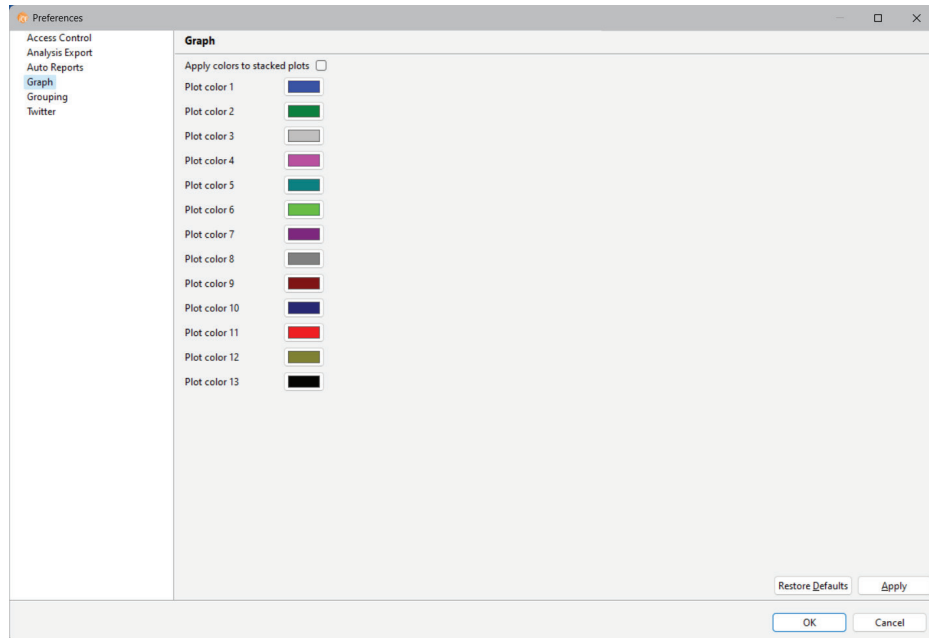


Selecting Custom Plot Colors for Graph Overlay

Select **Graph** in the sidebar.

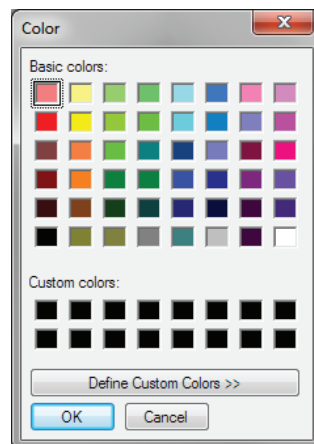
- **Apply colors to stacked plots** - This option applies the color scheme shown to individual plots when Stack the plots is selected in the Analysis screen's Graph pane. When this option isn't selected, all plots use the same color (this is the default setting).

NOTE: If **Apply colors to stacked plots** isn't selected, the colors shown are only applied to plots when Overlay the plots is selected in the Graph pane.



Changing Plot Colors

1. Click the button next to a Plot color number. You'll get a color selection box:



2. Select a color or define a custom color and click **OK**. The color button will update to the new color selected.
3. Repeat the steps above for any other plot colors.

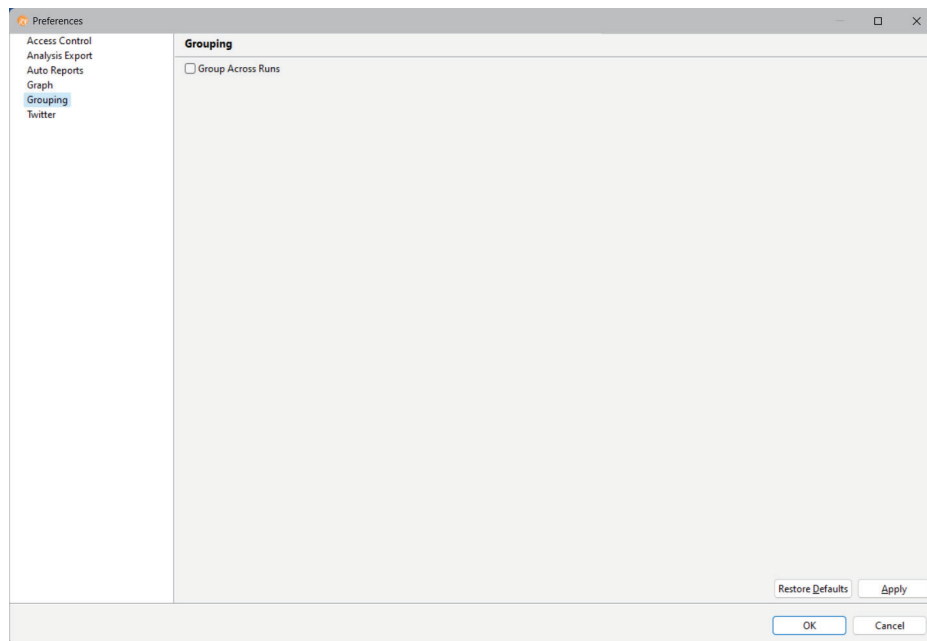
4. Check **Apply Colors to Stacked Plots** if you also want the new color settings to be used for the Stack the plots option in the Graph pane.
5. Click **Apply** to apply the new color settings to the plots currently displayed. This lets you see the changes without having to close the Graph window.
6. Click **OK** to save changes and exit.
7. Select **Overlay the plots** in the Graph pane. The new color scheme will be used.

Grouping Options

Select **Grouping** in the sidebar.

Selecting the **Group Across Runs** box groups samples with the same name together even if they're in different runs, so you can get statistics for the same samples across multiple runs. When the box isn't selected, only samples with the same name within the same run are grouped for statistics (this is the default setting).

NOTE: To activate grouping and get statistics for runs you have open in the Analysis Screen, select **View** in the main menu and click **Grouping**.



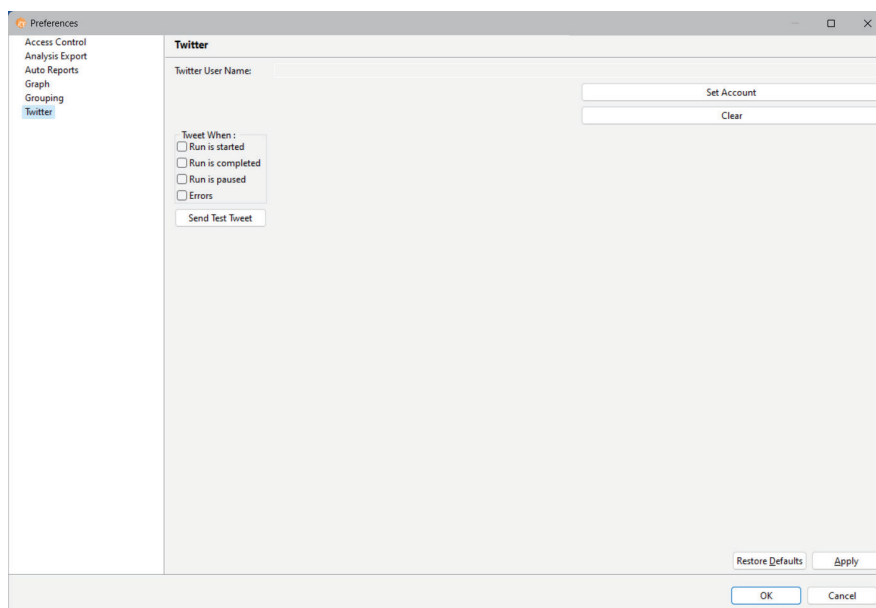
Setting Up Maurice Systems to Send Tweets

Select **Twitter** in the sidebar.

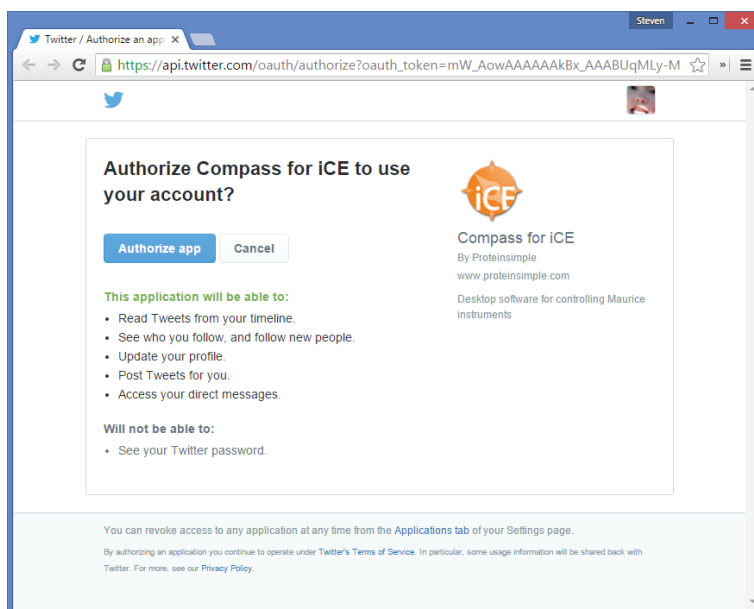
NOTES:

To set your Maurice system up to tweet, the computer you're using needs to be connected to the internet through a network connection or the local lab computer.

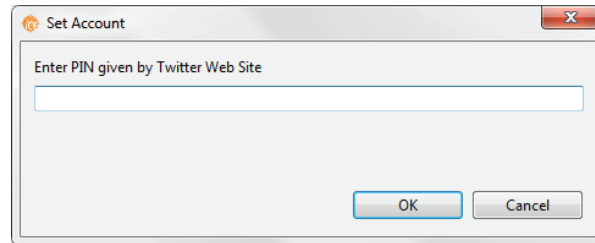
We recommend setting up separate Twitter accounts for each system. This lets multiple people in the lab follow run progress.



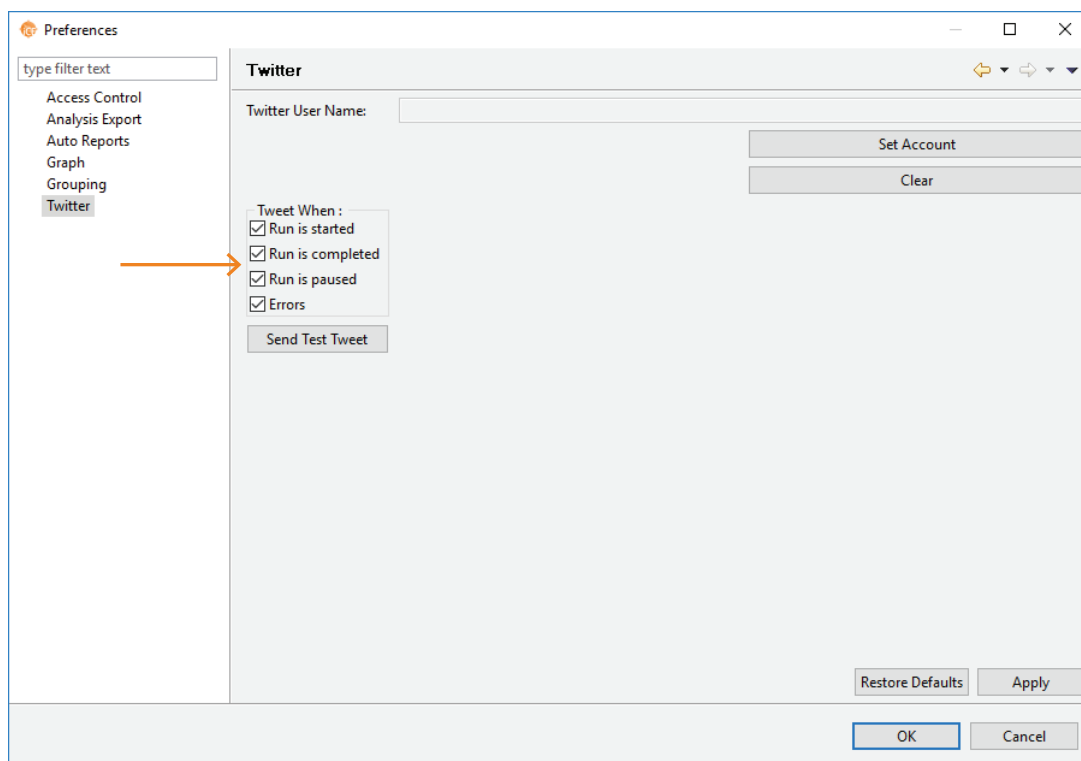
1. Click **Set Account**. A set account window will display in Compass for iCE and a browser window will open:



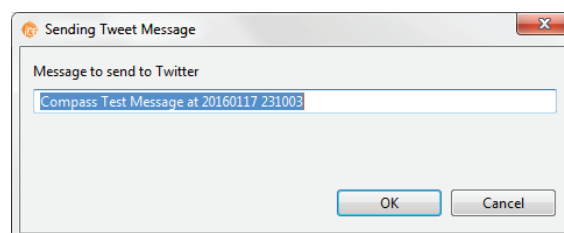
2. Enter a user name or email and password, then click **Authorize app**. A new page will display in the browser with a PIN number.
3. Enter the PIN number in the Compass for iCE set account window and click **OK**:



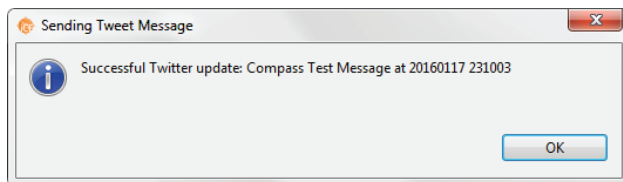
4. The user name will now appear in the Twitter User Name box. Select your Tweet When options and click **Apply**.



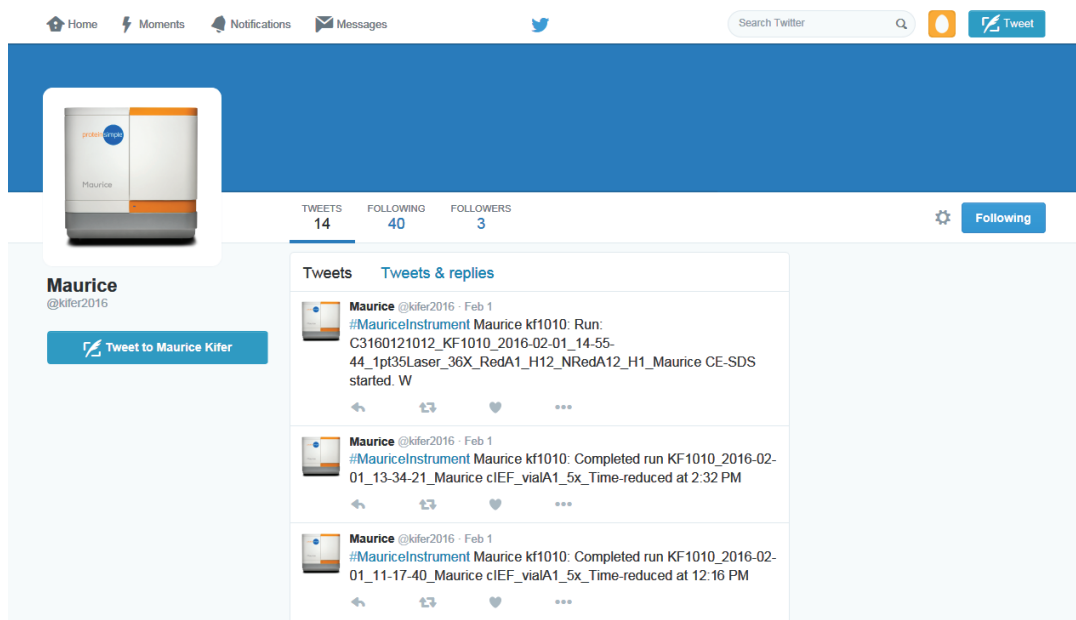
5. To confirm the Twitter account is receiving messages, click **Tweet Message**. Enter a test message and click **OK**.



6. If the test Tweet was successful, you'll get this message:



7. Click OK to save changes and exit. Maurice, Maurice C., Maurice S. and MauriceFlex will automatically tweet as the selected options occur:



Changing the Twitter Account

To change the Twitter account your system uses:

1. Select **Edit > Preferences**, then select **Twitter** in the left sidebar.
2. Click **Clear**.
3. Follow the same steps to set up the account as in “Setting Up Maurice Systems to Send Tweets”.

Chapter 22:

Compass Access Control and 21 CFR Part 11 Compliance

Chapter Overview

- Overview
- Compass Authorization Server for iCE
- Enabling Access Control
- Changing the Software Inactivity Auto Lock
- Logging In to Compass for iCE
- Saving Changes
- Signing Files
- Exporting Uncontrolled Files
- Instrument Command Log
- Run File History
- Troubleshooting Problems and Suggested Solutions



Overview

The Compass Access Control feature can be used to help satisfy the 21CFR Part 11 data security requirements when using Maurice instruments. When Access Control is enabled and the Compass Authorization Server for iCE has been installed:

- Users are required to log in to Compass for iCE when the software is launched
- A history of all actions is maintained
- Data files are signed and encrypted to prevent unauthorized changes (*e.g.*, all files are controlled)
- Each instrument maintains a history of user commands
- Each batch and data file includes a history of signed changes to the file

Compass for iCE can be run with or without Access Control enabled. When Access Control is disabled, no user log in is required and files are not encrypted or signed. The instrument history and file history are still maintained but the entries are not signed.

Compass Authorization Server for iCE

The Compass Authorization Server for iCE controls the log in access to Compass for iCE. In the simplest configuration, the server is run on the same computer as Compass for iCE and only that copy of Compass for iCE is controlled. A single server can also be used to control access to multiple copies of Compass for iCE running on different computers, so long as they have network access to the server. Multiple copies of the server may be run on the same network, and each server will have its own user database.

To enable Compass for iCE to use a particular Authorization Server for iCE, click **Edit**, then **Preferences** and **Access Control** and enter the server IP address using format X.X.X.X.

NOTES:

Always use the default port setting of 8443, this should not be changed.

If the server is installed on the same computer as Compass for iCE (*e.g.*, the local machine), enter localhost:8443 instead of the IP address. Contact your local IT Administrator to assist with installing the Compass Authorization Server for iCE in your preferred format.

Server Administration

The Authorization Server for iCE is configured through a web interface at the IP address of the server on port 8443. To access the Server home page, open any browser and type the IP address on port 8443 in a X.X.X.X:8443 or http://X.X.X.X:8443 format. Use localhost:8443 instead of the IP address if the Server is installed on the local machine.

NOTE: If you have upgraded your system from a previous version of Compass Authorization Service, the localhost login page will have changed from http://localhost:8000/admin/login/?next=/ to http://localhost:8443/admin/login/?next=/

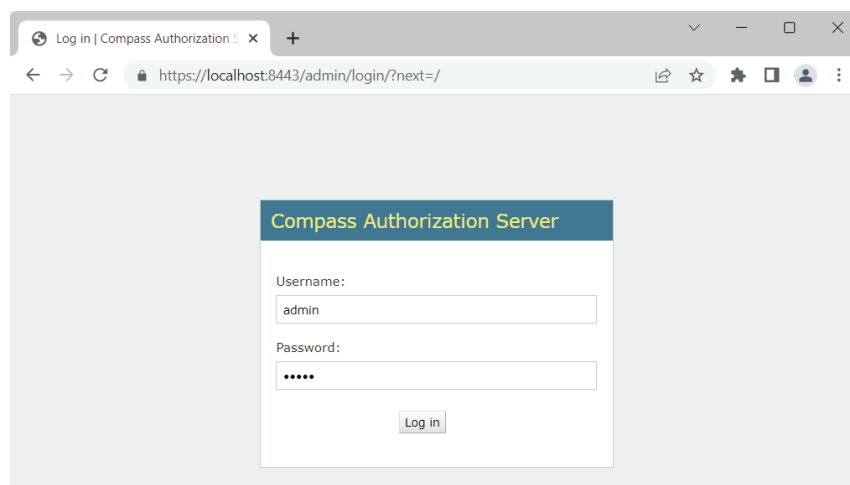
The default server administrator is:

- User: admin
- Password: admin

NOTE: The user account “admin” is not an active Compass for iCE login account. In order to log into Compass for iCE in the 21 CFR Part 11 compliant mode, at least one user account must be created in the Authorization Server for iCE application.

After installing the Compass Authorization Server for iCE, the administrator user name and password can be changed.

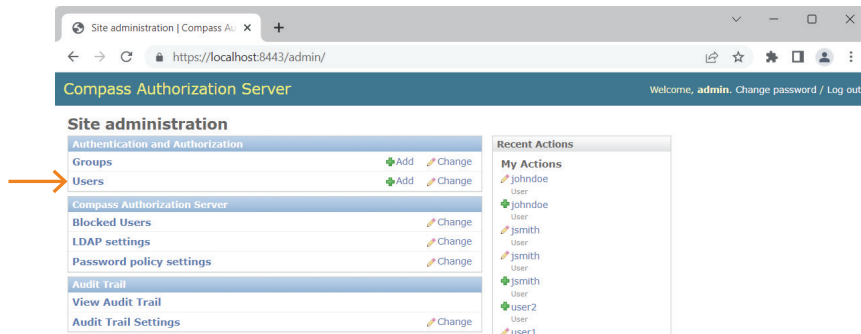
NOTE: Sites are responsible for maintaining their own admin and user account credentials. Record these and keep them safe. ProteinSimple can’t provide access once the admin account has been edited.



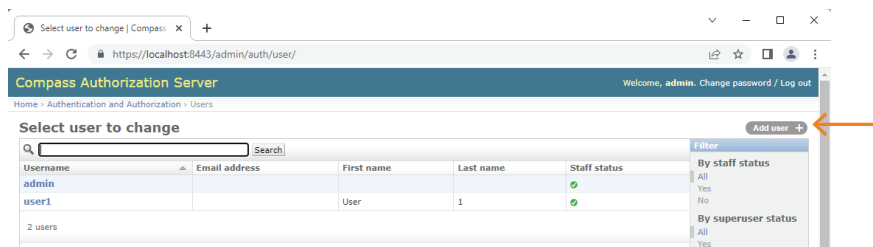
Adding Non-admin Users

Add a user to the server to allow that user to log in to Compass for iCE. To do this:

1. Select **Users** from the Site Administration home page:



2. From the Users page, select **Add User**:



3. Fill in the fields to create a new user:

The screenshot shows the 'Add user' form of the Compass Authorization Server. The form includes fields for Username, First name, Last name, Password, and Password confirmation. There is a checkbox for 'LDAP User' and a checkbox for 'Staff' (which is checked). The 'Staff' checkbox is accompanied by the text 'Staff users may log in and view audit trail'. At the bottom, there are three buttons: 'Save and add another', 'Save and continue editing', and 'Save'.

After adding a new user more information can be added:

The screenshot displays the 'Change user' page in the Compass Authorization Server. The page is titled 'Change user' and includes a 'History' button. The 'Username' field is set to 'user1'. The 'Password' field has a 'Click here to change' link. The 'Personal info' section contains fields for 'First name', 'Last name', and 'Email address'. The 'Permissions' section has three checkboxes: 'Active' (checked), 'Staff status', and 'Superuser status'. The 'Groups' section shows 'Available groups' and 'Chosen groups'. The 'Chosen groups' list contains 'Administrator'.

NOTE: Users are blocked after the number of login failures defined in the Password policy setting.

Permissions

All users can log in to Compass for iCE, but the commands available within Compass for iCE are controlled by Permission settings. Commands a user does not have permission to use will be disabled. After user permissions have been changed on the server the user must close and re-open Compass for iCE to use the new permissions.

Users can belong to groups that have multiple permissions such as Operator or Scientist:

The screenshot displays the Compass Admin interface. At the top, the browser address bar shows the URL `https://localhost:8443/admin/auth/user/5/`. The main content area is divided into two sections: 'Groups' and 'User permissions'.

Groups Section:

- Available groups:** A list with a search filter. The 'Chosen groups' list on the right contains: Administrator, Operator, Reviewer, and Scientist.
- Buttons:** 'Choose all' and 'Remove all'.
- Text:** 'The groups this user belongs to. A user will get all permissions granted to each of their groups. Hold down "Control", or "Command" on a Mac, to select more than one.'

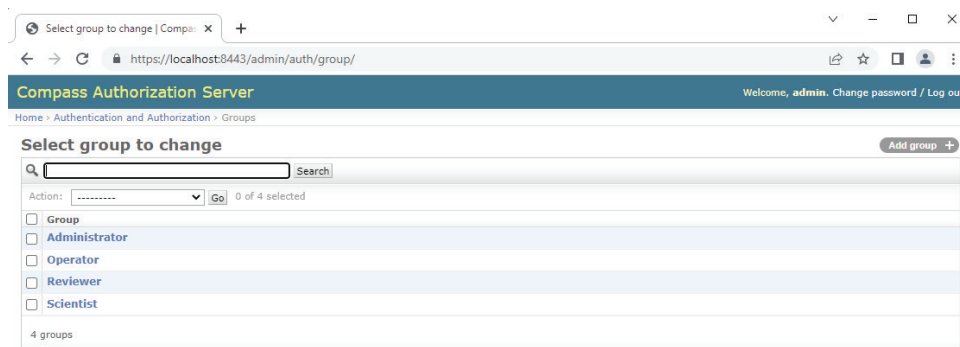
User permissions Section:

- Available user permissions:** A list with a search filter. The 'Chosen user permissions' list on the right is currently empty.
- Buttons:** 'Choose all' and 'Remove all'.

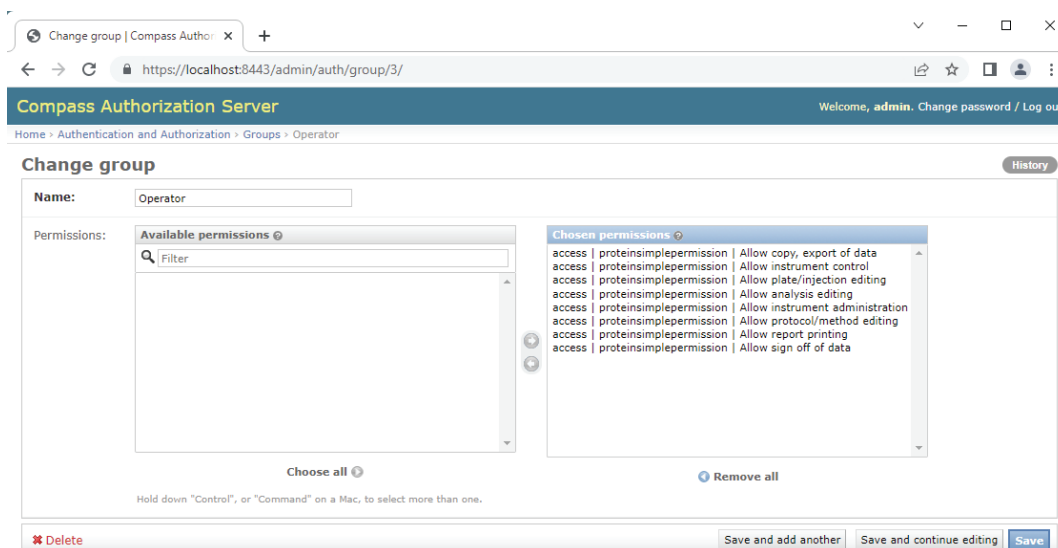
Use the Groups page to change the permissions in a group or create new groups:

A screenshot of a web browser showing the Compass Authorization Server administration page. The address bar shows 'https://localhost:8443/admin/'. The page has a dark blue header with the title 'Compass Authorization Server' and a welcome message 'Welcome, admin. Change password / Log out'. Below the header, there's a section titled 'Site administration' with a sub-section 'Authentication and Authorization'. An orange arrow points to the 'Groups' link in this sub-section. Other links include 'Users', 'Blocked Users', 'LDAP settings', 'Password policy settings', 'Audit Trail', 'View Audit Trail', and 'Audit Trail Settings'. To the right, there's a 'Recent Actions' panel titled 'My Actions' listing several user actions like 'johnndoe User' and 'jsmith User'.

To change permissions for a group click **Change**, then select a group:



Move individual group permissions in or out of the Available Permissions and Chosen Permissions boxes by selecting a permission in either box. Click the **left** or **right** arrow button to move the permission into the other box.



Here are the current user and group permissions:

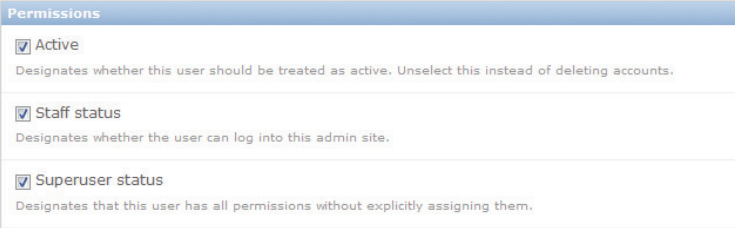
- **Allow analysis editing** - lets users change analysis settings
- **Allow copy, export of data** - lets users copy and export data
- **Allow instrument control** - gives users permission to connect to an instrument
- **Allow plate editing** - lets users update the plate layout and edit injections
- **Allow protocol/method editing** - gives users permission to edit methods
- **Allow report printing** - lets users print the Batch and Injection reports
- **Allow sign off of data** - gives users permission to sign off on changes using e-signatures

- **Allow instrument administration** - lets users update the Maurice embedded software, delete run dates, update the Compass for iCE software and turn Access Control on or off

Adding Admin Users

To create a user with administrator permissions:

1. Follow the steps described in “Adding Non-admin Users” on page 771” to create the admin user.
2. Under permissions, select **Staff status** and **Superuser status**:



Permissions	
<input checked="" type="checkbox"/> Active	Designates whether this user should be treated as active. Unselect this instead of deleting accounts.
<input checked="" type="checkbox"/> Staff status	Designates whether the user can log into this admin site.
<input checked="" type="checkbox"/> Superuser status	Designates that this user has all permissions without explicitly assigning them.

3. Assign the admin user to a group.

NOTE: Selecting Superuser status enables server permissions only. Admin users must be also be assigned to a group to in order to have Compass for iCE permissions.

Resetting User Passwords

NOTE: Users are blocked after the number of login failures defined in the Password policy setting.

To reset a user password:

1. Select **Users** from the Site Administration home page, then select the user to change. The following screen displays:

The screenshot shows a web browser window with the address bar displaying `https://localhost:8443/admin/auth/user/2/`. The page title is "Change user | Compass Authoriz...". The breadcrumb trail is "Home > Authentication and Authorization > Users > user1". The page content includes a "Change user" section with a "History" button. The "Username" field is "user1" with a note: "Required. 30 characters or fewer. Letters, digits and @/./+/-/_ only." The "Password" field has a link "Click here to change". Below this are sections for "Personal info" (First name, Last name, Email address) and "Permissions" (Active, Staff status, Superuser status).

2. Click the text link to access the password change form:

The screenshot shows a web browser window with the address bar displaying `https://localhost:8443/admin/auth/user/2/password/`. The page title is "Change password: user1 | Compass Authoriz...". The breadcrumb trail is "Home > Authentication and Authorization > Users > user1 > Change password". The page content includes a "Change password: user1" section with the prompt "Enter a new password for the user user1." Below this are two input fields: "New password:" and "New password confirmation:". A "Change password" button is located at the bottom right.

3. Enter the new password, then click **Change password**.

Audit Trail

Admin users with Staff Status can view, print and download the Audit Trail. Select **View Audit Trail** from the Site Administration home page to access it.

The image shows two screenshots of the Compass Authorization Server web application. The top screenshot shows the 'Site administration' menu with an orange arrow pointing to the 'View Audit Trail' link. The bottom screenshot shows the 'Audit Trail' page, which includes a table of logged actions.

Audit Trail
(current server version 3.0.0.189)
[Download as PDF](#)

Page: 1

datetime	machine	user	first name	last name	action	description	comment
2022-05-05 14:56 -0700	Compass Authorization Server	—	—	—	install	Version 3.0.0.189	—
2022-05-05 15:47 -0700	Compass Authorization Server	admin	—	—	login	success	—
2022-05-05 15:48 -0700	Compass Authorization Server	admin	—	—	User Created	Tu, Andrea as user1	—
2022-05-05 15:49 -0700	Compass Authorization Server	admin	—	—	User Modified	Removed group Operator from user user1	—
2022-05-05 15:49 -0700	Compass Authorization Server	admin	—	—	User Modified	Added group Administrator to user user1	—
2022-05-05 15:49 -0700	Compass Authorization Server	admin	—	—	User Modified	Added permission Allow analysis editing to user user1	—
2022-05-05 15:49 -0700	Compass Authorization Server	admin	—	—	User Modified	Added permission Allow copy, export of data to user user1	—
2022-05-05 15:49 -0700	Compass Authorization Server	admin	—	—	User Modified	Added permission Allow instrument administration to user user1	—
2022-05-05 15:49 -0700	Compass Authorization Server	admin	—	—	User Modified	Added permission Allow instrument control to user user1	—

Here are the actions currently logged in the Audit Trail:

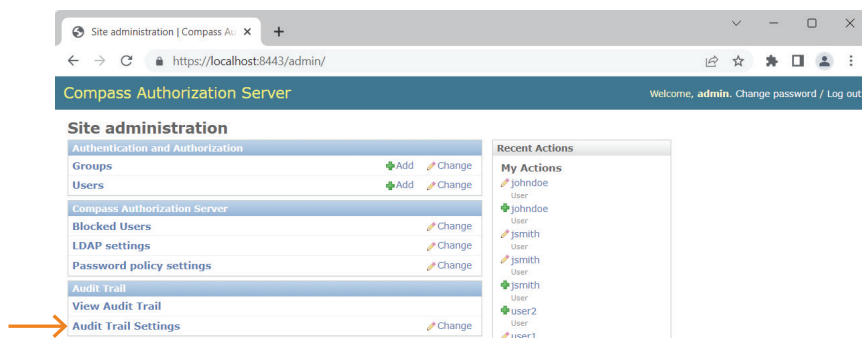
- **Login.** User logged into CAS or Compass.
- **Logout.** User logged out of CAS or Compass.
- **Start run.** User started a run in Compass.
- **Stop run.** User stopped a run in Compass.
- **Delete run.** Run file was deleted.
- **Cartridge Cleanup.** User initiated a cartridge cleanup routine from Compass.

- **Self test.** User initiated a self-test on an instrument.
- **Open runfile.** User opened a run file (*.mbz) in Compass.
- **Save runfile.** User saved a run file (*.mbz) in Compass to the file location shown.
- **Open batchfile.** User opened a batch file (*.batch) in Compass.
- **Save batchfile.** User saved a batch file (*.batch) in Compass to the file location shown.
- **Instrument upgrade.** User initiated an update of instrument software from Compass.
- **Change Control Settings.** User disabled Access Control from Compass.
- **Generate Batch Report.** User generated Batch Report from open batch in Compass.
- **Generate Injection Report.** User generated Injection Report from run in Compass.
- **Export spectra andi format.** User exported the raw and analyzed data traces and background for each injection in the run in .cdf format from Compass.
- **Export spectra text format.** User exported the raw and analyzed data traces and background for each injection in the run in .txt format from Compass.
- **Export tables.** User exported the results for all injections in the run in .txt format from Compass.
- **User created.** A new user was created on CAS.
- **User modified.** A user was modified on CAS.
- **User deleted.** A user was deleted on CAS.
- **Group created.** A group was created on CAS.
- **Group modified.** A group was modified on CAS.
- **Group deleted.** A group was deleted on CAS.
- **Change password.** A user's password was changed.
- **Sign off.** A user signed off on changes to the file location shown utilizing e-signatures in Compass.
- **Edit LDAP settings.** LDAP settings in CAS were modified.
- **Edit password policy.** Password policy settings were modified in CAS.
- **Install.** CAS software was installed.
- **Upgrade.** CAS software was upgraded to new version.
- **Blocked.** A user has been blocked from logging in.

- **Unblock.** A user has been unblocked, allowed to log in.
- **Emergency Stop.** User initiated the Emergency Stop command to instrument from Compass.
- **Pause.** User added a Pause operation to the running batch on the instrument from Compass.
- **Continue.** User initiated the Continue command to the instrument from Compass.
- **Stop events.** User initiated a stop to the instrument from Compass.

Audit Trail Settings

The Secure Audit Trail PDF setting (selected by default), allows users to download audit trail PDFs securely. Secure audit trail PDFs can be viewed and printed, but content cannot be copied or modified.



Password Policy Settings

These settings let administrators set password policies. Select **Password policy settings** from the Site Administration home page to make changes.

Select password policy setting to change

Display name	Value
Number of previous passwords to compare to	3
Minimum number of uppercase characters	1
Minimum amount of symbol characters	0
Minimum amount of number characters	1
Minimum number of lowercase characters	1
Minimum password length	8
Number of login attempts permitted	3
Days password is valid	180

8 password policy settings

Save

LDAP Settings

LDAP settings allow you to connect the Compass Authorization Server for iCE to your own network's domain controller, so users can log on with their existing network password. With LDAP, passwords are not maintained by the Compass Authorization Server for iCE they are administered by the network admin.

First select **LDAP settings** from the Site Administration page and set your LDAP settings.

Change LDAP setting

☐ Enabled

Address:

Domain:

SSL Port:

☒ Sync Information

Save and continue editing Save

Next, add users as described in “Adding Non-admin Users” on page 771 and select the LDAP User checkbox. Passwords aren't required for LDAP users.

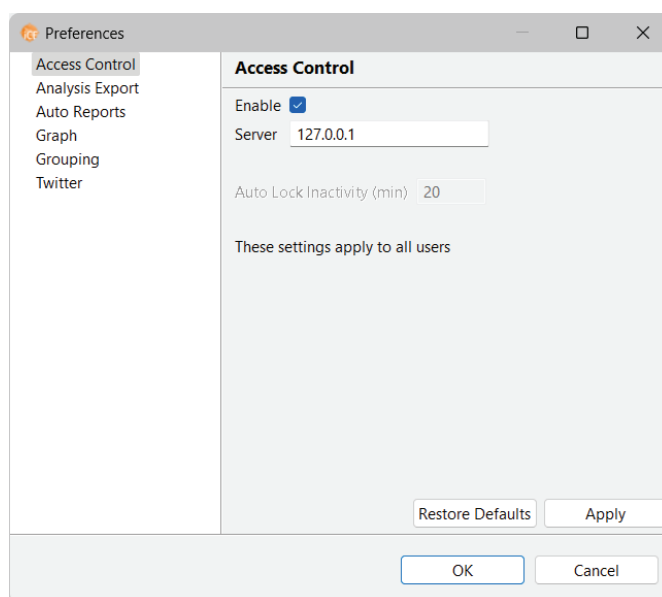
Encryption Details

Compass for iCE uses the SHA1 hash algorithm to generate a 160 bit hash code that is unique for all files. All files saved by Compass for iCE are encrypted with a digital key. This key along with the hash codes guarantees the file history is correct and no other edits were made. All changes saved to a file have the electronic signature of the user who saved the file. The **e-Signature** command allows a user to sign off on a state such as approved or verified.

There is no individual ownership of files, all users who log into Compass for iCE can open any file.

Enabling Access Control

Access Control is enabled in **Preferences**. Select **Edit** in the main menu, click **Preferences**, then select **Access Control**.



To enable Access Control:

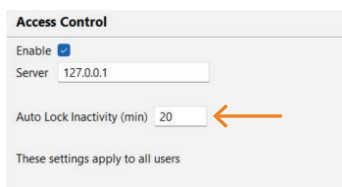
1. Check the **Enable** box.
2. Enter the IP address of the Authorization Server for iCE. Use format X.X.X.X:8443 or localhost:8443 if installing the server on the local machine.
3. Close Compass for iCE. The next time the software is launched, a user log in will be required.

NOTE: Access Control can only be disabled by logging into Compass for iCE and deselecting the **Enable** box in the Access Control page of Preferences.

Changing the Software Inactivity Auto Lock

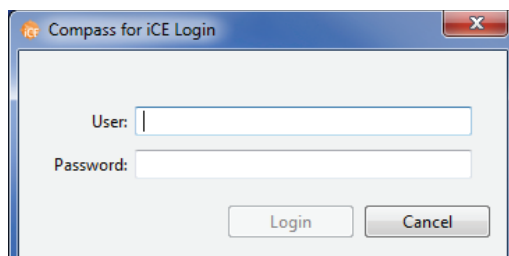
Compass for iCE software automatically locks to prevent user access after a period of inactivity. Once the software is locked, users must login again. Only users with instrument administrator permission can set the auto lock inactivity time.

1. Select **Edit > Preferences > Access Control**.
2. Enter an **Auto Lock Inactivity** time in minutes. The default setting is 20 minutes.



Logging In to Compass for iCE

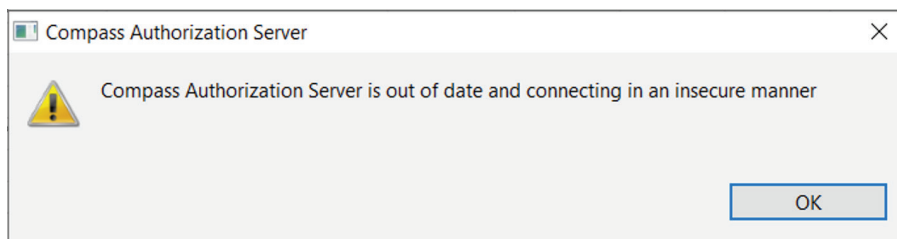
With Access Control enabled, all users must log in to Compass for iCE whenever the software is launched.



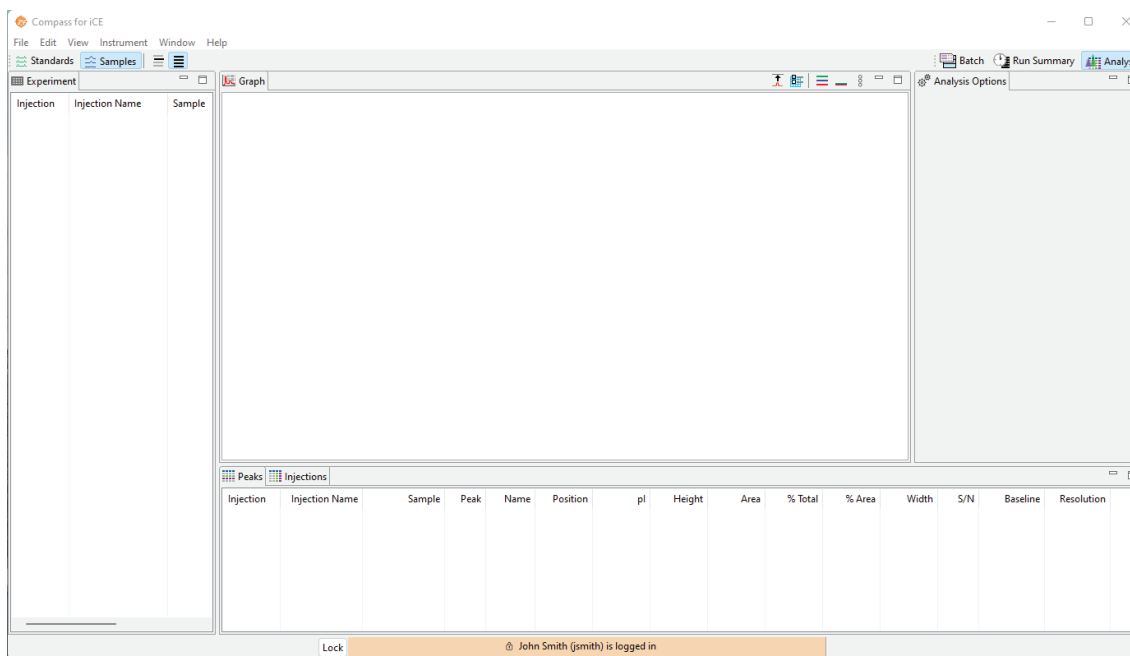
Enter your user name and password previously setup by your Compass for iCE Administrator.

NOTE: Your account will be blocked after a certain number of login failures. If this happens, contact your administrator to unblock the account.

A window will appear if Compass for iCE is connecting to a previous version of the Compass Authorization Server that does not support a secure connection.

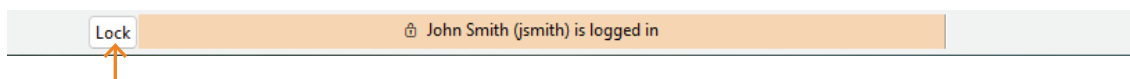


A successful log in will display the Compass for iCE main window with the user information in the lower status bar. The full user name is displayed with the unique user ID in parenthesis:

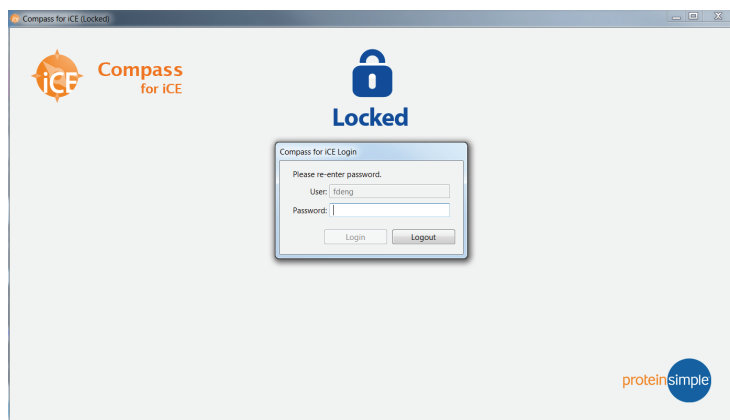


Locking and Unlocking the Application

You can click the **Lock** button to lock Compass for iCE and prevent access by other users. To unlock the application, users must re-enter their password.



If there is no activity in Compass for iCE for 20 minutes, the application automatically locks. Users must re-enter their passwords to perform any controlled actions:

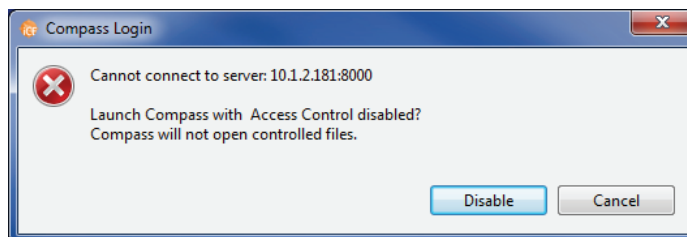


Resolving Log In Issues

Log in failures may occur when:

- The server is temporarily unavailable
- Compass for iCE is using the wrong IP address

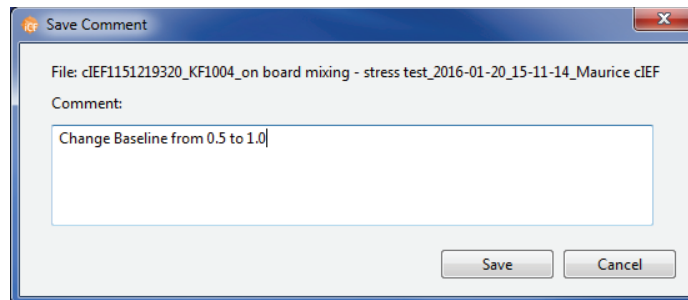
When this happens, the following message displays:



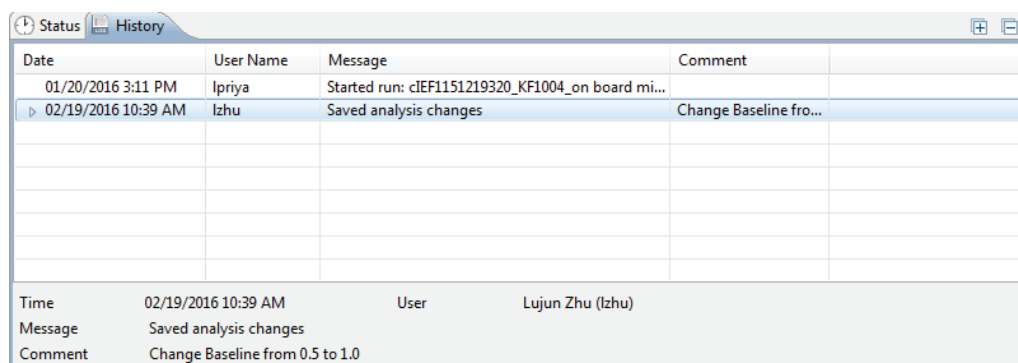
Click **Disable** to restart Compass for iCE with Access Control disabled. Verify or correct the server IP address then close and restart the software to log in with Access Control enabled.

Saving Changes

When **Save** is selected from the **File** menu, a dialog box will display to allow you to enter a comment before saving the signed file:



The comment is added to the signature entry in the file History:

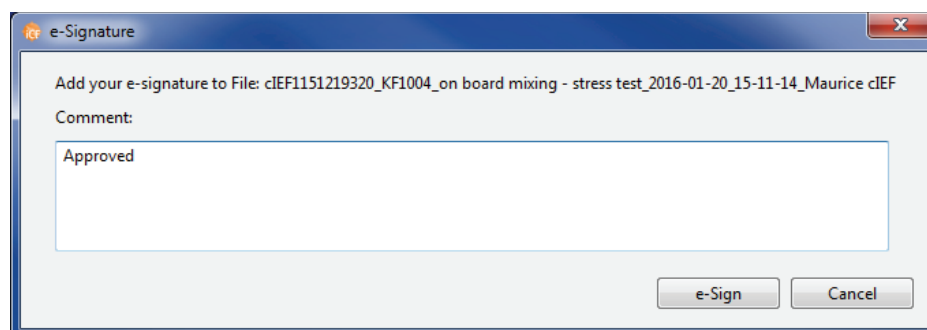
A screenshot of a 'History' window. It has a tabbed interface with 'Status' and 'History' tabs. The 'History' tab is active, showing a table with columns: Date, User Name, Message, and Comment. The table has two rows. The first row is: 01/20/2016 3:11 PM, Ipriya, Started run: cIEF1151219320_KF1004_on board mi..., and an empty comment. The second row is highlighted: 02/19/2016 10:39 AM, Izhu, Saved analysis changes, and Change Baseline fro... Below the table, there is a summary section with labels: Time (02/19/2016 10:39 AM), User (Lujun Zhu (Izhu)), Message (Saved analysis changes), and Comment (Change Baseline from 0.5 to 1.0).

Date	User Name	Message	Comment
01/20/2016 3:11 PM	Ipriya	Started run: cIEF1151219320_KF1004_on board mi...	
02/19/2016 10:39 AM	Izhu	Saved analysis changes	Change Baseline fro...

Time	02/19/2016 10:39 AM	User	Lujun Zhu (Izhu)
Message	Saved analysis changes		
Comment	Change Baseline from 0.5 to 1.0		

Signing Files

Select **e-Signature** from the **File** menu to add an electronic signature to a file.

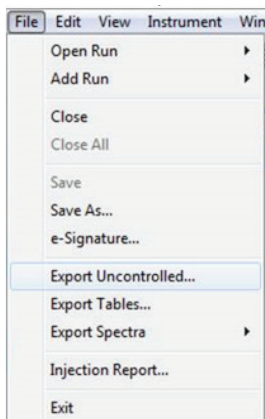


The signed entry will be added to the file History with the meaning of the signature entered in the comment, such as *Approved* or *Verified*.

Status		History		
Date	User Name	Message	Comment	
01/20/2016 3:11 PM	Ipriya	Started run: cIEF1151219320_KF1004_on board mi...		
02/19/2016 10:39 AM	Izhu	Saved analysis changes	Change Baseline fro...	
02/19/2016 10:50 AM	Izhu	e-Signature applied	Approved	
Time	02/19/2016 10:50 AM		User	Lujun Zhu (Izhu)
Message	e-Signature applied			
Comment	Approved			

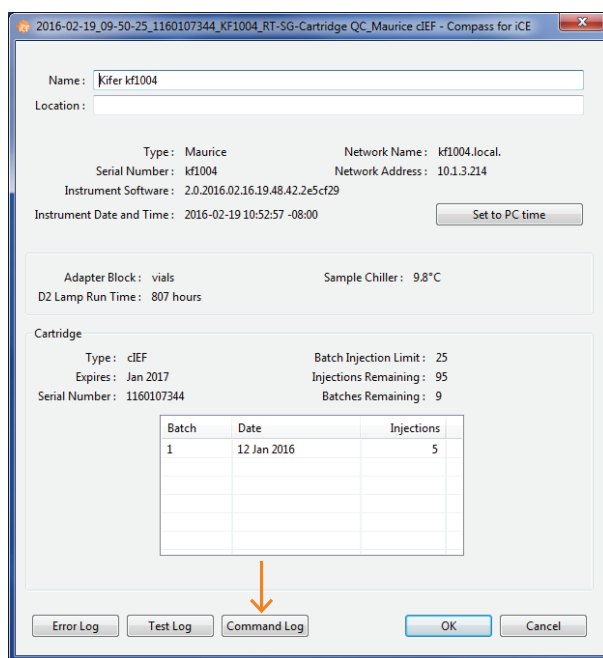
Exporting Uncontrolled Files

You can export an uncontrolled file by selecting **File > Export Uncontrolled....** The export will also be logged in the Audit Trail and the History log for the exported file.

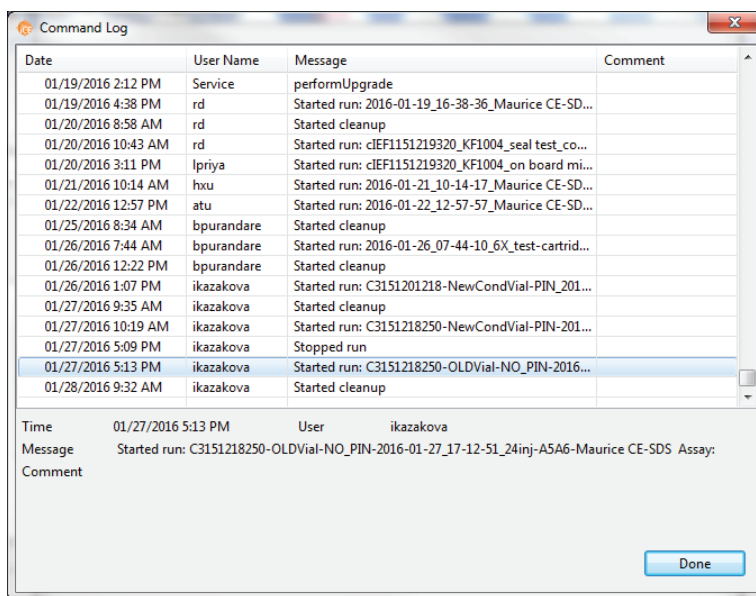


Instrument Command Log

The Instrument Command Log can be viewed at any time by selecting the **Instrument** menu and clicking **Properties**, and then clicking the **Command Log** button:



The Command Log lists all the commands sent to the instrument that were signed by the user who sent the command. If you want to copy the Command Log at any time, right click in the table and select **Copy**, then paste into another document.



Run File History

Select the **Run Summary** screen tab and then the **History** tab to see the file History. To copy the file History, right click in the table and select **Copy**, then paste into another document.

The screenshot displays the 'Run File History' window in the Compass for iCE software. The window has a menu bar (File, Edit, Instrument, Window, Help) and a toolbar with 'Batch', 'Run Summary', and 'Analysis' buttons. The main area is divided into two panes. The left pane, titled 'Injections', contains a table with columns: Sample ID, Location, Method, and Status. It lists 24 injection events, all with a status of 'Completed'. The right pane, titled 'History', contains a table with columns: Date, User Name, Message, and Comment. It shows three events: 'Started run: cIEF1151219320_KF1004_on board mi...' by 'Ipriya' on 01/20/2016, 'Saved analysis changes' by 'Izhu' on 02/19/2016, and 'e-Signature applied' by 'Izhu' on 02/19/2016. Below the 'History' table, there is a section for 'Time', 'Message', and 'Comment' for the selected event, showing '02/19/2016 10:50 AM', 'e-Signature applied', and 'Approved' respectively. At the bottom of the window, a status bar indicates 'Lujun Zhu (Izhu) is logged in.' and a 'Lock' button is present.

Troubleshooting Problems and Suggested Solutions

If any of the following error messages are encountered, follow the recommended steps below to resolve the issue.

- **Unknown user name or password.**
 - Check if the Caps Lock is on, user name and password are case sensitive.
 - Ask a Compass for iCE administrator to confirm your user name. If your password is unknown then the administrator can reset your password (see Resetting User Passwords for more information).
- **Server not available.**
 - From the **Edit** menu, click **Preferences** and then **Access Control** to confirm the server address is set to the correct Authorization Server for iCE address. Compass for iCE must be able to reach the server on the network.
- **Controlled file cannot be opened without log in.** To open a controlled Run file, enable Access Control by clicking **Edit**, then **Preferences** and **Access Control**. Select **Enable**, close Compass for iCE, then re-launch the software with a valid log in.
- **Uncontrolled file cannot be opened when logged in.** To open an uncontrolled Run file, disable Access Control by clicking **Edit**, then **Preferences** and **Access Control**. Deselect **Enable**, close Compass for iCE then re-launch the software.

NOTES:

Only users with Instrument Administrator permission can turn Access Control on or off. This event will also be logged in the Audit Trail.

Uncontrolled files cannot be opened when Compass Access Control is enabled (controlled mode).

- **Command disabled.** Certain commands are only available when a user with the correct permissions is logged in. To change user permissions, use a web browser to log in to the Authorization Server for iCE web interface at the address shown on the **Access Control** page in **Preferences**, such as: 10.1.3.231:8443.
- **Compass for iCE does not prompt for log in.** Compass for iCE will only prompt for a log in on launch when Access Control is enabled in Preferences. Enable Access Control by clicking **Edit**, then **Preferences** and **Access Control**. Select **Enable**, close Compass for iCE, then re-launch the software. You should now be prompted for a log in.

Chapter 23:

Maintenance and Troubleshooting

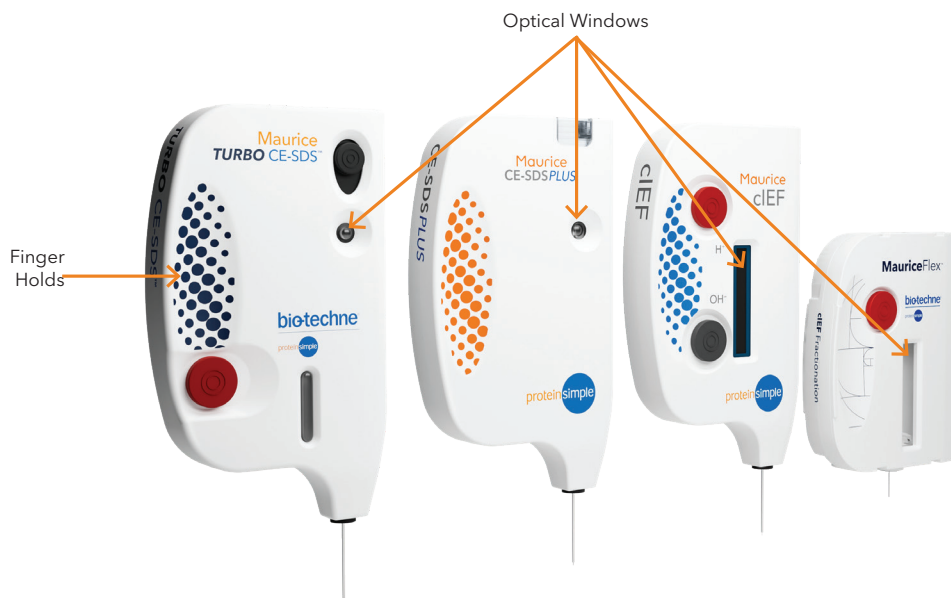
Chapter Overview

- Cartridge Handling and Care
- Installing or Replacing the System Filter
- Maintenance
- Spare Parts
- Software Updates
- Instrument Software (Embedded) Updates
- Frequently Asked Questions: General
- Frequently Asked Questions: cIEF Applications
- Frequently Asked Questions: cIEF Fractionation Applications
- Frequently Asked Questions: CE-SDS Applications
- Best Practices
- Troubleshooting

Cartridge Handling and Care

The cIEF, cIEF Fractionation, CE-SDS PLUS and Turbo CE-SDS Cartridges were developed for use with Maurice systems. Each cartridge is individually tested and shipped with a Certificate of Analysis, are also online at <https://www.bio-techne.com/resources/cofa-finder-tool>. The cartridges are shipped dry and should be stored free of liquid.

- **Cartridges need to be used, cleaned and stored properly to reach their maximum lifetime.** Additional cleaning vials (PN 046-125) and cartridge inserts (PN 046-124) for CE-SDS PLUS cartridges may be purchased separately.
- Store cartridges in their original packaging at room temperature when you receive them.
- Hold cartridges using the blue or orange finger holds on either side of the cartridge.
- Don't touch the recessed optical windows of the cartridge.



- Whenever you handle the cartridge or remove it from its packaging, make sure the cartridge inlet doesn't come in contact with any surfaces. A damaged inlet may compromise the cartridge and cause a failed injection.



- cIEF Cartridges are guaranteed for 100 injections and an absolute maximum of 200 injections.
- CE-SDS PLUS Cartridges are guaranteed for 100 injections and an absolute maximum of 500 injections.

- Turbo CE-SDS Cartridges are guaranteed for 100 injections.
- cIEF Fractionation Cartridges support 15 injections.
- Maurice reads the cartridge's RFID and keeps track of how many injections are left for you automatically.
- **Always clean the cartridge before storing.** See page 120 for cIEF Cartridge cleaning steps, page 182 or page 248 for cIEF Fractionation Cartridge cleaning steps, page 316 for CE-SDS PLUS Cartridge cleaning steps and page 385 for Turbo CE-SDS Cartridge cleaning steps.
- Always store the cartridge in its original packaging at room temperature when not in use.

cIEF Cartridge

A cIEF Cartridge is guaranteed for 100 injections with an absolute maximum of 200 injections and a maximum of 20 batches. Its RFID will keep track of how many are left for you.

- The cartridge is designed for use with common cIEF reagents like methyl cellulose, ampholytes, urea and other protein solubilizing agents, anolyte and catholyte, but over exposure or high concentrations of certain components can harm it. The cartridge is optimized for 0.35% methyl cellulose in the sample mixture.
- Make sure to not get any liquid on the cartridge's optical window.
- **Always clean the electrolyte tank and run the cartridge post-run cleanup at the end of a run.** See page 120 for cIEF Cartridge post-run cleaning steps.

Compatibility with Sample Components

- **Methyl Cellulose (MC):** The sample mix must contain 0.35% methyl cellulose. The cartridge must be flushed with 0.5% methyl cellulose between runs.
- **Solvents:** The cartridge is not compatible with organic solvents. Do not rinse with untested solvents and minimize the amount of solvent in the sample mix.
- **Salt and surfactants:** High current can harm the internal coating in the cartridge capillary. High concentrations of salt and surfactants in the sample mix can generate high currents above 40 microamps. This high current will compress the pH gradient and also damage the cartridge. Please take care to minimize the concentration of salts in the final sample mix to below 15 mM. To keep current at a minimum, we suggested using only non-ionic or zwitterionic surfactants. Don't use aromatic surfactants as they can interfere with sample detection.

Cleaning the Outside of the Cartridge

If you see spikes in your data, the outside of the cartridge should be cleaned with canned air. You'll need to use residue- and moisture-free canned air to prevent fouling of the optical path through the separation capillary.

1. Place the can's nozzle or tube opening 10–12 inches from the cartridge surface. Then depress the aerosol actuator down about halfway so you get a gentle flow of air.
2. Sweep the air stream across the entire length of the optical window.
3. Flip the cartridge over and repeat the prior steps.

4. Flip the cartridge over again and gently clean the top surface one last time before reinstalling in Maurice.

cIEF Fractionation Cartridge

A cIEF Fractionation Cartridge supports up to 15 injections and a maximum of four injections per MauriceFlex cIEF batch and one injection per MauriceFlex Fractionation batch. Its RFID will keep track of how many are left for you.

- The cartridge is designed for use with common cIEF reagents like methyl cellulose, ampholytes, urea and other protein solubilizing agents, anolyte and catholyte, but over exposure or high concentrations of certain components can harm it. The cartridge also requires 0.35% methyl cellulose in the sample mixture. We do not recommend using more than 0.8% methyl cellulose in the sample mixture.
- Make sure to not get any liquid on the cartridge's optical window.
- Always clean the electrolyte tanks and run the cartridge post-run cleanup at the end of a run. See page 182 or page 248 for cIEF Fractionation Cartridge post-run cleanup steps.

Compatibility with Sample Components

- **Methyl Cellulose (MC):** The sample mix must contain 0.35% methyl cellulose. The cartridge must be flushed with 0.5% methyl cellulose between runs.
- **Solvents:** The cartridge is not compatible with organic solvents. Do not rinse with solvents and minimize the amount of solvent in the sample mix.
- **Salt and surfactants:** High current can harm the internal coating in the cartridge capillary. High concentrations of salt and surfactants in the sample mix can generate high currents above 90 microamps. This high current will compress the pH gradient and also damage the cartridge. Please take care to minimize the concentration of salts in the final sample mix to below 10 mM. To keep current at a minimum, we suggested using only non-ionic or zwitterionic surfactants. Don't use aromatic surfactants as they can interfere with sample detection.

Cleaning the Outside of the Cartridge

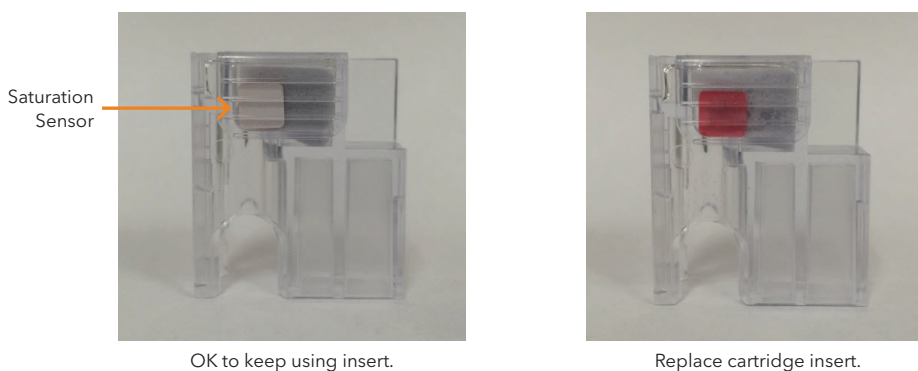
If you see spikes in your data, the outside of the cartridge should be cleaned with canned air. You'll need to use residue- and moisture-free canned air to prevent fouling of the optical path through the separation capillary.

1. Place the can's nozzle or tube opening 10-12 inches from the cartridge surface. Then depress the aerosol actuator down about halfway so you get a gentle flow of air.
2. Sweep the air stream across the entire length of the optical window.
3. Flip the cartridge over and repeat the prior steps.
4. Flip the cartridge over again and gently clean the top surface one last time before reinstalling in MauriceFlex.

CE-SDS PLUS Cartridges

A CE-SDS PLUS Cartridge is guaranteed for 100 injections with an absolute maximum of 500 injections, a maximum of 48 injections per batch and a maximum of 25 batches. Its RFID will keep track of how many are left for you.

- Once you've inserted the Top Running Buffer vial in the cartridge insert, the CE-SDS PLUS Cartridge **must** be kept in an upright position at all times.
- If you see any separation matrix sticking to the cartridge inlet, soak the inlet in DI water for 5 minutes. Then wipe it using a lint-free laboratory wipe that's been moistened with DI water. Other than this, no external cleaning of the cartridge is required.
- Check the saturation sensor on the back of the CE-SDS PLUS Cartridge insert after every run. If it's red, you'll need to use a new cartridge insert for your next batch. If the saturation sensor isn't red, you can keep using the current cartridge insert.



- We highly recommend running a Cartridge Purge if the cartridge has not been used in more than 3 months. See "Cartridge Purge" on page 439 for instructions.
- Always run the post-run cleanup at the end of a run. See page 316 for CE-SDS PLUS Cartridge post-run steps.

Compatibility with Sample Components

- **Salt:** The salt concentration in your sample should be <50 mM. Higher concentrations will adversely affect electrokinetic injections. Dilute your sample with CE-SDS Sample Buffer to reach the recommended salt concentration. If the protein concentration in your samples is low, we recommend desalting the sample.

Turbo CE-SDS Cartridges

A Turbo CE-SDS Cartridge is guaranteed for 100 injections with a maximum of 96 injections per batch and a maximum of 25 batches. Its RFID will keep track of how many are left for you.

- Prepare your samples and reagents before preparing the cartridge. Allowing the Separation Matrix to sit in the cartridge for longer than 15 minutes may result in cartridge clogs.
- If you see any separation matrix sticking to the cartridge inlet, soak the inlet in DI water for 5 minutes. Then wipe it using a lint-free laboratory wipe that's been moistened with DI water. Other than this, no external cleaning of the cartridge is required.
- We highly recommend running a Cartridge Purge if the cartridge has not been used in more than 3 months. See "Cartridge Purge" on page 439 for instructions.

- **Always clean the Running Buffer and Waste Tanks and run the post-run cleanup at the end of a run.** See page 385 for Turbo CE-SDS Cartridge post-run cleanup steps.

Compatibility with Sample Components

- **Salt:** The salt concentration in your sample should be <50 mM. Higher concentrations will adversely affect electrokinetic injections. Dilute your sample with CE-SDS Sample Buffer to reach the recommended salt concentration. If the protein concentration in your samples is low, we recommend desalting the sample.

Installing or Replacing the System Filter

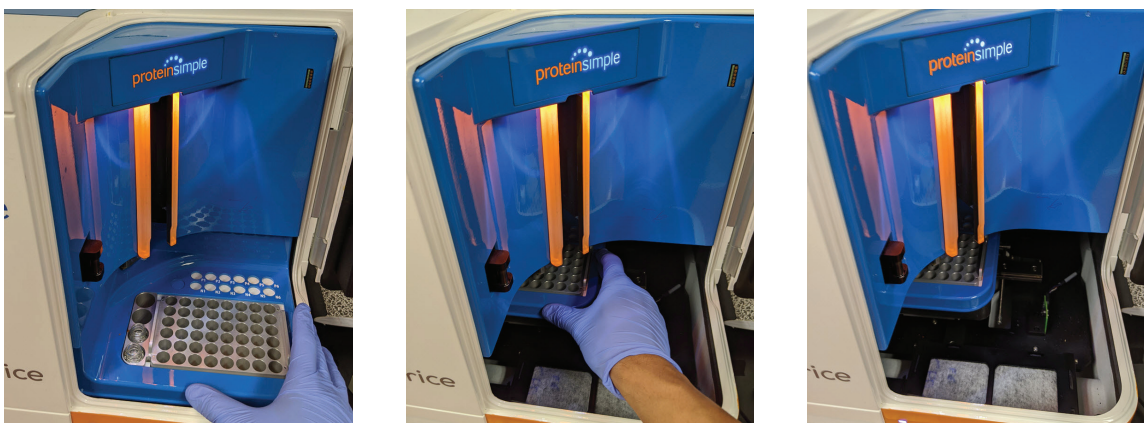
If the optional Maurice System Filter Upgrade is installed on your Maurice, Maurice S. or MauriceFlex system, we recommend changing the filter every 3 months.

1. Make sure Maurice isn't running a batch (light should be steady blue).
2. Open Maurice's door by touching the metal plate on top of the door. If a cartridge is installed, remove it. For more info on removing and storing cartridges, see "Post-batch Procedures" on page 316 for a CE-SDS PLUS Cartridge and "Post-batch Procedures" on page 385 for a Turbo CE-SDS Cartridge.

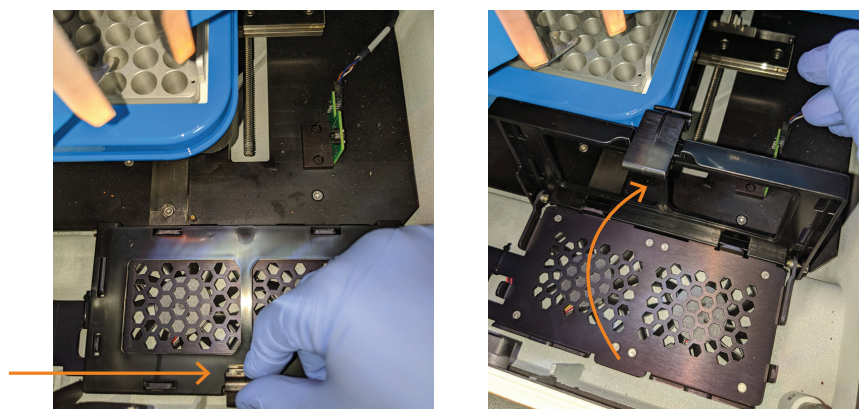


NOTE: The indicator light on Maurice's front panel will blink rapidly as the door disengages.

3. Swing the door open to access the reagents and samples platform.
4. Gently apply pressure to the front right corner of the reagents and samples platform and push the platform to the left and then back until the front corner is below the cartridge slot (**orange** lights). The filter holder will now be accessible.



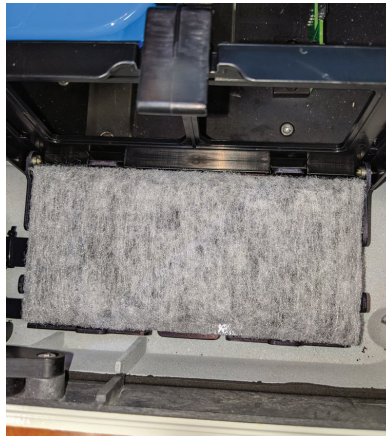
5. Open the filter holder by squeezing the front tab inward (orange arrow) and lifting the hinged holder door up and open. The images below show the filter holder without a filter installed.



6. Remove the old filter (if present) and discard.
7. Remove the new filter from the plastic wrapping.

NOTE: Use scissors to cut open the filter packaging, taking care not to cut the filter itself. Tearing open the package may bend or damage the filter. We recommend handling the filter with gloves to prevent contamination.

8. Place the new filter in the holder, it doesn't matter which side is up or down. It should lie flat and within the raised edges on the sides of the holder.

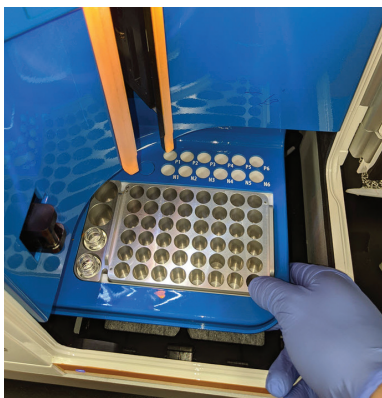


9. Close the holder lid until the latch clicks into place. Gently tug on the latch to make sure the filter holder is fully closed.



10. Gently move the reagents and sample platform back into position by pulling on the corner of the platform until the reagent and sample locations are visible.

NOTE: The exact positioning of the platform isn't critical. Maurice will automatically home the platform to the correct position once the door has been closed.



11. Close Maurice's door. A steady blue light will indicate that the system is ready to use.

Maintenance

Daily

- Empty out any condensation in the sample plate or sample tray inserts. Wipe out the sample block too if needed.
- Dispose of your samples and reagent vials after each run. Compass for iCE will let you know when a cartridge is at the end of its useful life and should be discarded.

Yearly

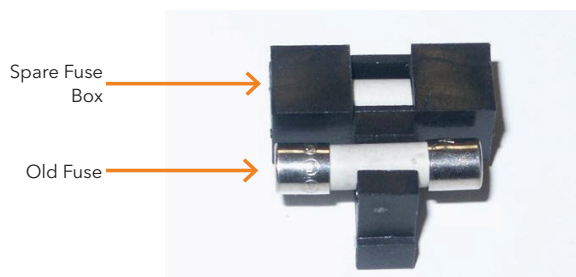
We recommend Maurice systems have annual preventive maintenance performed by an authorized ProteinSimple engineer. Please contact Technical Support to schedule a visit.

Changing the Fuse

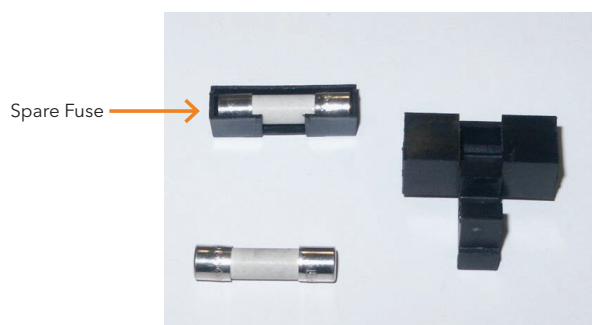
1. Power down Maurice and unplug his power cord.
2. Locate the fuse holder on his rear panel.



3. Use a flat-head screwdriver to gently pry the fuse holder open and remove the fuse holder.



4. Remove the old fuse.
5. There's a spare fuse in the small box. Pull the box out, pull the new fuse out and use it to replace defective fuse.



6. Reinsert the fuse holder.
7. Plug Maurice's power cord back in and turn his power on.

Spare Parts

Maurice, Maurice C., Maurice S. and MauriceFlex don't have any user replaceable parts except for the filter and power fuse. Please contact ProteinSimple Technical Support if they get sick and need repair.

- **Maurice Filter Kit** - PN 046-576, includes 4 filters and 8 plate lids.

!WARNING!

You can't replace or service any parts on Maurice systems except for the filter and power entry fuse.

Software Updates

To check for software updates, first make sure the computer you're using has an active internet connection. Then go to Compass for iCE software, select **Help** in the main menu and click **Check for Updates**. If you don't have internet access, call your FAS for assistance on getting the latest update.

Instrument Software (Embedded) Updates

To check for embedded updates, go to Compass for iCE software, select **Instrument** in the main menu, then **Update** and select **Network**. If you're not on the network, contact Technical Support for assistance on getting the latest update.

Frequently Asked Questions: General

Should I download a secured or non-secured batch or injection report?

See the table below for what you can do with a secured PDF (Compass default) vs. a non-secured PDF.

Parameter	Secured PDF	Non-Secured PDF
Security Method	Random password security	No security
Printing	Allowed	Allowed
Document Assembly	Not allowed	Allowed
Content Copying	Not allowed	Allowed
Content Copying for Accessibility	Not allowed	Allowed
Page Extraction	Not allowed	Allowed
Commenting	Not allowed	Allowed
Filling of Form Fields	Not allowed	Allowed
Signing	Not allowed	Allowed
Creation of Template Pages	Not allowed	Allowed
Encryption Level	128-bit RC4	N/A

Frequently Asked Questions: cIEF Applications

NOTE: Please refer to the Maurice cIEF Method Development Guide for info on initial application conditions and method optimization.

I have a new protein sample to analyze. What starting conditions should I use?

Begin with the following initial sample conditions:

- Carrier ampholytes: pH 3-10 Pharmalytes (4%)
- Additive: 0.35% methyl cellulose
- Sample analyte: 0.1 mg/mL concentration in final solution. The balance of the solution should be HPLC-grade deionized water.
- 10 mM arginine
- 1X pI 4.05 and 9.99 markers

NOTES:

If you want to use pI markers below pI 4.05, we suggest adding iminodiacetic acid 10 mM IDA to the sample solution

ProteinSimple provides a 1% methyl cellulose solution (PN 101876).

Another way to start is to simply use the same sample conditions used if you were successful in running this sample on slab gel IEF. Use the same carrier ampholytes and additives for analysis on Maurice systems.

What carrier ampholytes are commercially available, and which one is best for my sample?

At present, carrier ampholytes are commercially available from four different manufacturers under the following brand names:

- Pharmalytes (Cytiva - formerly GE)
- Servalyts (Serva)

Other carrier ampholytes exist, however, they are all repackaged and resold using one of the products listed above.

Each brand may give slightly different separation resolution due to slight differences in ampholytic compositions. Identification of the optimal carrier ampholytes for a given protein sample is best determined experimentally.

Along with native fluorescence, Maurice systems use UV absorption detection at 280 nm. All carrier ampholytes exhibit some degree of absorption at this wavelength, which causes some baseline noise. Pharmalytes have low and uniform UV absorption and produce no background signal in native fluorescence along the entire pH range. Because of the low background noise of Pharmalytes, these ampholytes are recommended for initial sample conditions.

Does my sample matrix affect my results?

Yes. However, the sample is diluted 20X in carrier ampholytes, methyl cellulose and HPLC-grade deionized water, which minimizes matrix effects. For example, if the concentration of your sample stock solution is 2 mg/ μ L, 10 μ L of the sample can be directly diluted by adding 112 μ L of HPLC-grade deionized water, 8 μ L of pH 3–10 Pharmalytes and 70 μ L of 1% methyl cellulose. The final solution is 200 μ L with a sample concentration of 0.1 mg/ μ L. In this example, the original sample matrix will not affect analysis.

If the original stock sample concentration is >2 mg/mL and contains high salt concentrations, then diluting further and using native fluorescence to boost signal is recommended. If detection in both absorbance and fluorescence is required, desalting may be necessary.

I cannot get reproducible peaks due to sample precipitation, what should I do?

1. Dilute the sample.
2. The native fluorescence detection mode provides higher sensitivity and can be used for low concentration samples.

Doing either or both of the above reduces the potential for aggregation or precipitation. If the issue is still observed, several additives can be tested to increase protein solubility. The following additives have been successfully tested with Maurice systems and should help stabilize proteins during analysis:

- Up to 40% SimpleSol
- Up to 4 M urea
- Up to 20% formamide
- Up to 25% sorbitol
- Up to 25% sucrose
- Up to 25% glycerol
- Denaturing conditions, such as 8 M urea

In rare cases, sample precipitation may be caused by the carrier ampholytes. To avoid this problem, try using a different brand of carrier ampholytes. If additive conditions for stable sample runs have been established for gel IEF, then these additive conditions can often be successfully used for cIEF analysis on Maurice systems.

NOTE: All additives may change the pI value of the protein slightly, especially if the method uses pI markers in the acidic range of the pH gradient.

How do I prepare sample solutions in 8 M Urea?

For a 200 µL final sample solution:

- Weigh 96 mg of urea powder in a 1.5 mL centrifuge tube.
- Add 32 µL HPLC-grade deionized water, 70 µL of 1% methyl cellulose, 8 µL of carrier ampholytes and 10 µL of sample to the urea powder in the centrifuge tube.

This will make a final volume of 200 µL (96 mg urea powder and 120 µL liquid). If more than 10 µL or less than 10 µL of sample is added, the volume of water should be adjusted to ensure a final volume of 120 µL.

NOTE: Urea must always be prepared fresh before use. Do not heat to dissolve urea in solution.

When running samples in 8 M urea, the focusing time should be increased 1-2 minutes compared to normal conditions. This is due to the higher viscosity of the urea-containing solution.

How do I choose a protein solubilizer?

Protein precipitation or aggregation may occur during the isoelectric focusing step as proteins concentrate close to their pIs, leading to non-reproducible peak profiles. One of the strategies to address this issue is the addition of protein solubilizers to the protein-ampholyte mixture. Urea, a chaotropic agent, is a popularly used protein solubilizer that disrupts non-covalent bonds to reduce protein precipitation and aggregation. However, a number of known issues exist for urea, including potential induction of carbamylation, anodic gradient drift causing shift of apparent pI values, and protein

denaturation above 4 M. In addition, urea needs to be prepared fresh daily which creates an additional step in sample analysis.

SimpleSol can effectively solubilize proteins for icIEF but is a significantly more stable agent than urea, eliminating the need for analysts to prepare urea fresh every day. SimpleSol is also stable when pre-mixed with methyl cellulose as opposed to urea, and is compatible with absorbance and native fluorescence detection on Maurice icIEF. In addition, SimpleSol has less of an impact on the acidic portion of the pH gradient formed during icIEF compared to urea, resulting in more stable pI values of protein peaks relative to using urea. Consider SimpleSol if you are concerned about solution stability, pH gradient formation, or ease-of-use for your protein solubilizer.

How can I identify peaks in different runs and different samples?

A reliable way to identify peaks in electropherograms is to use internal pI Markers. First run the sample without internal pI Markers. The pI values of sample peaks can be estimated from their peak positions relative to the full pI range of the carrier ampholytes.

In Compass for iCE graphs, the left side of the electropherogram is the anodic end of the capillary (acidic) and the right side is the cathodic end (basic). For example, if pH 3-10 Pharmalytes are used as the carrier ampholytes, the x-axis of the electropherogram represents pI 3 to pI 10 from left to right. The pI value of a peak at the middle of the trace should be about 6.5.

Two pI Markers are mixed into the sample solution. Ideally, the peaks of the two markers should bracket the sample peaks and the two marker peaks should be as close as possible in order to achieve good precision in peak identification.

The electropherograms of the sample mixed with pI Markers are processed using Compass for iCE for pI determination. The software uses the method settings to automatically identify the pI markers to convert from pixel position in the Markers View to pI in the Samples View.

In this way, the sample peaks are identified by their measured pI values. The precision of peak identification by measuring the pI values using Maurice systems is less than +0.03 pH units.

Since the measured pI value of a protein is affected by many factors such as sample matrix and the type of carrier ampholytes used, to correctly identify peaks in different samples or different runs, all runs should be done under the exact same conditions.

What kind of pI markers can I use?

ProteinSimple recommends using low molecular weight amphoteric compounds with well defined isoelectric points and strong UV absorbance when using Maurice systems. Conversely, we do not recommend using protein pI markers since they often produce multiple isoelectric points and, on occasion, may interact with the sample analyte.

ProteinSimple offers a selection of absorbance- and fluorescence-compatible pI markers at pI 3.38, 4.05, 5.85, 6.14, 7.05, 8.40, 9.50, 9.99 and 10.17.

The distance between the two pI Markers in my sample electropherograms is different from run to run even though I use the same pI Markers and carrier ampholytes. What is the reason for this?

Usually this is caused by different salt concentrations in the sample solutions. Salt can compress the pH gradient created by the carrier ampholytes. So, the higher the salt concentration, the shorter the distance between the two pI Markers.

However, since the whole pH gradient is compressed by the salt, this will not affect peak identification results as long as two pI Markers are used and their peaks bracket the sample peaks.

Can I use narrow pH range carrier ampholytes to improve the resolution for my sample?

Yes. The most efficient way to do this is to use a mixture of narrow pH range carrier ampholytes and wide pH range carrier ampholytes. The proportion of carrier ampholytes can be from 1:1 (narrow range: wide range) up to 5:1 depending on the resolution requirement. Focusing time should be increased with the increasing proportion of the narrow pH range carrier ampholytes, from 6 to 12 minutes.

The measured pI value of my sample peak is slightly different when I use different pI Markers or different carrier ampholytes with the same pI markers. What is the reason for this?

When using different pI Markers, the small difference in the measured pI value is due to the slight non-linearity of the pH gradient established by the carrier ampholytes along the separation capillary. Compass for iCE pI calibration assumes that the pH gradient is perfectly linear between the two pI Markers. In reality, carrier ampholytes are not perfectly linear throughout their pH gradient.

When different carrier ampholytes are used, their pH gradients may be slightly different causing a small difference in measured pI value. This effect is most obvious when using a carrier ampholyte mixture (i.e. narrow and wide pH range carrier ampholytes). Under these conditions, the pH gradient will not be linear at the edges of the overlapping pH regions of the different carrier ampholytes. Changing the ratio of the different carrier ampholytes in the mixture will affect the measured pI values of a protein.

In conclusion, only measured pI values obtained using the same carrier ampholytes and the same pI markers can be compared. Also, as long as the run conditions are the same, the measured pI values can be used to identify protein peaks.

Frequently Asked Questions: cIEF Fractionation Applications

NOTE: Please refer to the MauriceFlex cIEF Fractionation Method Development Guide for info on initial application conditions and method optimization. Refer to the Maurice cIEF Method Development Guide and “Frequently Asked Questions: cIEF Applications” on page 801 for initial sample and method optimization of Maurice cIEF prior to running cIEF fractionation.

How do I optimize a cIEF fractionation batch?

Method development starts with a previously developed Maurice cIEF method, and you'll typically only need to optimize a few parameters from there. For many molecules, adding an additional 20 mM arginine to the final sample with the separation mix used for Maurice cIEF methods provides sufficient fractionation performance and don't need further development.

For more challenging molecules that do not fractionate well and/or have limited solubility refer to the MauriceFlex cIEF Fractionation Method Development Guide. In summary, the first step is to start with a generic method and evaluate the

separation current and focused peak profile. Once you have the desired peak profile, fractionation can be optimized for the desired peak purity and concentration.

What is the recommended sample concentration for cIEF fractionation?

cIEF fractionation typically uses higher sample concentrations than used for Maurice cIEF. ProteinSimple recommends initial sample concentrations of 1 mg/mL. Increasing the sample concentration can improve concentrations of the fractions, but may also result in aggregation. Addition of protein solubilizers, such as 2 M urea, will eliminate sample aggregation most of the time.

How are sample peaks mobilized and collected during cIEF fractionation?

Samples migrate out of the cartridge capillary and into the fraction wells using chemical mobilization with 5 mM ammonium acetate (Mobilizer Solution). During the chemical mobilization, acetate ions from the Mobilizer Solution enter the capillary to displace the hydroxy ions when a voltage is applied. This results in the change of the local pH and charged molecules in the capillary, including protein charge variants, become positively charged. The positively charged protein moves toward the cathodic end of the capillary, eventually into the fraction wells with Mobilizer Solution. Other charged molecules in the sample, like ampholytes, will also be mobilized and collected in fractions.

How do I prepare sample solutions in 4 M urea?

The total sample volume should be 200 µL. You'll need the following amount of components per 160 µL of master mix prepared with 48 mg urea powder:

- 70 µL of 1% Methyl Cellulose
- 8 µL of carrier ampholytes (total)
- 1.5 µL of each pI marker
- 8 µL of 500 mM Arginine
- Add DI water to reach 160 µL total volume

NOTE: Urea must always be prepared fresh before use. Do not heat to dissolve urea in solution.

How do I check the purity and identity of the sample in the fractions?

To verify the presence and identity of charge variants in specific fractions, we recommend performing a Maurice cIEF batch with the fractions and a sample of the unfractionated molecule as a reference. Peaks detected in the fractions can be aligned based on pI with peaks in the unfractionated reference sample to verify identity. Peak purity can be assessed in the fractions by measuring percent peak area for each peak detected in a fraction. We recommend using fluorescence detection in a Maurice cIEF batch for calculating peak purity and estimating concentration of charge variants within a fraction.

Can I use the cIEF Fractionation Cartridge for standard Maurice cIEF analysis/batches?

No, the cIEF Fractionation Cartridge cannot be used with a Maurice cIEF batch. In addition, the cIEF Fractionation Cartridge is not designed for quantitative cIEF analysis of proteins – use the Maurice cIEF Cartridge for this purpose. Focused images on a cIEF Fractionation Cartridge are provided for qualitative purposes only.

Frequently Asked Questions: CE-SDS Applications

NOTE: Please refer to the Maurice CE-SDS Application Guide for info on initial application conditions and method optimization.

What is the molecular weight size range?

The CE-SDS application lets you size proteins between sizing between 10 and 270 kDa.

What wavelength is used for detection?

Proteins are detected via absorbance at 220 nm.

How good is the resolution?

The CE-SDS application was optimized to achieve baseline resolution (>1.5) between the glycosylated and non-glycosylated heavy chain.

Can I run non-reduced samples?

You can run both reduced and non-reduced IgG samples on Maurice. See the instructions in Chapter 12: “Running CE-SDS PLUS Applications” or Chapter 14: “Running Turbo CE-SDS Applications” for sample preparation. Compass for iCE also has default methods for running reduced and non-reduced samples, so you can select the appropriate method for your sample.

Best Practices

Maurice cIEF

Follow the setup and cleanup instructions in the Maurice cIEF Cartridge Product Insert (PN 046-298).

Important Tips:

- Immediately empty and rinse the electrolyte tanks on the cartridges after opening the Maurice door at the end of a run. After clean-up, leave the stoppers off the cartridge for storage. Store the stoppers and the cartridge together in its original packaging.
- Confirm you are using the correct glass reagents for the application you are running. cIEF batch reagents should be prepared using 2 mL glass reagent vials (PN 046-017).
- Use only ProteinSimple-recommended plate seals (<https://biochromato.com/slit-seal/>). Use plate seals for standard Maurice cIEF ONLY, DO NOT use plate seals for MauriceFlex cIEF, MauriceFlex Fractionation or CE-SDS applications.

MauriceFlex cIEF and MauriceFlex Fractionation

Follow the setup and cleanup instructions in the MauriceFlex cIEF Fractionation Cartridge Product Insert (PN PL3-0058).

Important Tips:

- Immediately empty and rinse the electrolyte tank on the cartridge after opening the MauriceFlex door at the end of a run. After clean-up, leave the stopper off the cartridge for storage. Store the stopper and the cartridge together in its original packaging.
- When installing the fractionation adapter, confirm the adapter is seated and sitting flat in MauriceFlex.
- Remove batch reagent vials from the reagent platform prior to placing the fractionation adapter in MauriceFlex if a Maurice cIEF or CE-SDS application was previously performed.
- Confirm you are using the correct glass reagent vials for the application you are running. MauriceFlex cIEF and MauriceFlex Fractionation batch reagents should be prepared using glass reagent vials, 0.3 mL (PN 110-0018) for the Fluorescence Calibration Standard and 2 mL crimp top glass vials (PN 110-0019) for all other reagents.
- DO NOT use plate seals for MauriceFlex cIEF or MauriceFlex Fractionation applications.

Maurice CE-SDS

Follow the setup and cleanup instructions in the Maurice CE-SDS PLUS Cartridge Product Insert (PN PL3-0013) of the Maurice Turbo CE-SDS Product Insert (PN PL3-055) depending on the cartridge you are using.

Important Tips:

- Clean the cartridge according to the procedure outlined in the Product Insert immediately after opening the Maurice door at the end of a run.
- The Top Running Buffer (TRB) vial should be inserted into the CE-SDS PLUS cartridge only once. Do not remove and re-insert the vial.
- Once a TRB vial is inserted into a cartridge, keep the cartridge upright.
- CE-SDS PLUS Cleaning vials may only be used up to 5 times.
- Immediately empty and rinse the Top Chamber on the Turbo CE-SDS cartridge after opening the Maurice door at the end of a run. After clean-up, leave the stopper off the cartridge for storage. Store the stopper and the cartridge together in its original packaging.
- DO NOT use plate seals for CE-SDS applications.

General Tips

- Use only the recommended sample vials, sample plates and reagent vials from ProteinSimple.
- Avoid touching the cartridge capillary tip to any surfaces.
- Store all cartridges in their original packaging when not in use.
- Use only compressed air to clean dust off the Maurice system. Do not clean with alcohol.
- Verify batch reagents are prepared using the appropriate glass vial.
- If a run is stopped midway, always perform a cleanup or purge.
- DO NOT use plate seals for CE-SDS applications.
- For Maurice cIEF, CE-SDS PLUS and Turbo CE-SDS batches only:
 - Verify that pressure caps have O-rings installed prior to use.
 - Verify that vials and caps are sitting flush and level in the sample tray. Check that caps are on correctly and vials are seated correctly.
 - Verify that the correct vial caps are being used in the correct positions according to the Assay map.
- For batches run on a MauriceFlex system:
 - Verify batch reagents are locked in the reagent platform or fractionation adapter before starting a batch.

Troubleshooting

Compass for iCE lets you run a self test on Maurice, Maurice C., Maurice S. and MauriceFlex systems. These diagnostic tests check many internal components and can help you determine if you have an instrument issue or not. Go to “Instrument Self Test” on page 446 for details on how to get started.

For more Maurice and application troubleshooting information, please contact ProteinSimple Technical Support at (888) 607-9692 (option 3), support@proteinsimple.com or visit <https://www.bio-technne.com/support/instrument-support>. You can also contact your local Field Application Specialist for help.

Maurice System Troubleshooting

Problem	Solution
Maurice does not initialize The status light on Maurice's front panel does not turn blue.	Restart the instrument. If Maurice still does not initialize, contact Technical Support at support@proteinsimple.com .
Maurice's door won't open Maurice's door is unresponsive and won't open when you touch the Touch Plate at the top of the door.	Turn off Maurice and engage the emergency door release on the lower, left-hand side of the machine. Remove the cap and press up on the bar inside to pop the door open. Return the cap to its original position and try to power on Maurice. If the door is unresponsive contact Technical Support at support@proteinsimple.com .

cIEF Application Troubleshooting

Problem	Solution
Error message: Calibration standard not detected The Fluorescence Calibration Standard was not stored at the proper temperature, the reagent vial is in the wrong position, or the reagent vial doesn't contain the right volume of solution. The cartridge may be clogged or the tip may be damaged.	<ul style="list-style-type: none"> The Fluorescence Calibration Standard should always be stored at 4 °C. If it's been stored at another temperature, replace the bottle with a new one then prepare a fresh reagent vial with the new solution. Make sure reagent vials are placed in the right positions in the sample and reagent platform. Confirm that there is 500 µL of Fluorescence Calibration Standard in the reagent vial. Ensure the cap on the reagent vial for the Fluorescent Calibration Standard is properly aligned and tightened. Ensure that you are using the newer pressure caps: cIEF blue pressure caps (PN 046-573) or CE-SDS orange pressure caps (PN 046-572). Examine the tip of the cartridge. If damaged, discard the cartridge and replace with a new one. <p>If the cartridge tip is undamaged, soak the tip in warm DI water overnight, with the cartridge positioned vertically (prop against a wall or stable object). Perform a Cartridge Purge after the overnight soaking.</p>

Abnormal Focusing

Problem	Solution
cIEF Cartridge tank level low If the anolyte or catholyte tank fluid level is not high enough to make good contact with the electrodes, the current will drop to < 2 μ A.	2 mL of electrolyte is needed in each tank. Aspirate out all the electrolyte solution and add 2 mL of anolyte and catholyte into their designated tanks.
Electrolyte contamination	Replace the anolyte and catholyte solutions in the electrolyte tanks.
Current keeps increasing beyond 80 μA	<p>Either the electrolytes are in the wrong tanks or the cartridge is defective.</p> <ol style="list-style-type: none"> 1. Immediately stop the system and run the Maurice Cartridge Purge in Compass for iCE. 2. When the purge is done, remove the cartridge, wash out the tanks and transfer new electrolyte solutions into the proper tanks. 3. Rerun the System Suitability test to confirm the internal coating is intact. 4. Replace the cartridge if the System Suitability run fails to meet resolution specifications or the current still remains high at >80 μA.

Artificial Peaks

Problem	Solution
Dust or particulates on the optical window	<p>Remove the cartridge. Use compressed air or nitrogen to gently clean the optical window, then reinstall the cartridge.</p> <hr/> <p>NOTE: Don't wash or submerge the cartridge's optical window in water or solvent.</p> <hr/>
Particles or precipitate in sample	Use an aqueous additive to stabilize the sample solution.
Air bubble in capillary	<ul style="list-style-type: none"> • Always spin samples for 5 minutes at 1000 xg before adding them to sample vials or wells of a 96-well plate. • Dispense the solution at the bottom of the vial/well to avoid trapping any air bubbles.

cIEF Fractionation Application Troubleshooting

Problem	Solution
Error message: Calibration standard not detected The Fluorescence Calibration Standard was not stored at the proper temperature, the reagent vial is in the wrong position, or the reagent vial doesn't contain the right volume of solution. The cartridge may be clogged or the tip may be damaged.	<ul style="list-style-type: none"> The Fluorescence Calibration Standard should always be stored at 4 °C. If it's been stored at another temperature, replace the bottle with a new one then prepare a fresh reagent vial with the new solution. Make sure reagent vials are placed in the right positions in the fractionation adapter. Confirm that there is 350 µL of Fluorescence Calibration Standard in the reagent vial. Confirm the Fluorescence Calibration Standard is prepared in the glass reagent vial with insert, 0.3 mL (PN 110-0018) Examine the tip of the cartridge. If damaged, discard the cartridge and replace with a new one.

Abnormal Focusing

Problem	Solution
Unstable current: cIEF Fractionation Cartridge tank level low. If the anolyte tank fluid level is not high enough to make good contact with the electrodes, the current will drop to < 2 µA.	2 mL of anolyte is needed in the tank. Aspirate out all the anolyte solution and add 2 mL of fresh to the tank.
Unstable current, large peak shift or poor resolution: Electrolyte contamination. If the electrolytes are contaminated or aged, the separation current could be unstable, peak position could shift significantly, or the peak resolution could be poor.	Replace the anolyte and catholyte solutions in the anolyte tank in the cartridge and catholyte vial. Do not reuse electrolytes between batches.
Current keeps increasing beyond 80 µA. Either the electrolytes are in the wrong locations or the cartridge is defective.	<ol style="list-style-type: none"> Immediately stop the system and run the Maurice Cartridge Purge in Compass for iCE. When the purge is done, remove the cartridge, wash out the tank and transfer new anolyte solution into the tank. Add new catholyte solution to a new reagent vial. Inspect the cartridge to ensure there is no anolyte leaking out of the tank. Replace the cartridge if the current still remains high at >80 µA.

Artificial Peaks

Problem	Solution
Dust or particulates on the optical window	Remove the cartridge. Use compressed air or nitrogen to gently clean the optical window, then reinstall the cartridge. NOTE: Don't wash or submerge the cartridge's optical window in water or solvent.
Particles or precipitate in sample	Use an aqueous additive to stabilize the sample solution.
Air bubble in capillary	<ul style="list-style-type: none"> Always spin samples for 5 minutes at 1000 xg before adding them to sample vials or wells of a 96-well plate. Dispense the solution at the bottom of the vial/well to avoid trapping any air bubbles

CE-SDS Application Troubleshooting**Error Message: Detected current below minimum threshold**

Problem	Solution
Run stops before the first injection Capillary is likely clogged.	Run the Cartridge Purge then start the batch again. If the error reoccurs, replace the cartridge with a new one.

Spikes, Poor Resolution

Problem	Solution
Air bubble in capillary	Always spin samples for 10 min at 1000 xg and use fresh reagents to minimize bubble occurrence.

Late Peak Arrival, Poor Resolution

Problem	Solution
Top Running Buffer vial leak	Use a new Top Running Buffer vial.
Insufficient conditioning	Use fresh Conditioning Buffers with each run.
Partial capillary clog	Run the Cartridge Purge.

Current Drifts to 0, No Signal, Saturation Sensor is Red

Problem	Solution
Non-viscous liquid in Separation Buffer vial	Make sure reagent vials are placed in the right positions in the sample and reagent platform.
Top Running Buffer vial is overloaded	Use a new cartridge insert and new Top Running Buffer vial.
Top Running Buffer vial leak	Use a new cartridge insert and new Top Running Buffer vial.

Low Signal

Problem	Solution
Sample composition is affecting the electrokinetic injection	Make sure the salt and protein concentrations of your sample are within the recommended ranges.

Rising Baseline

Problem	Solution
UV lamp is approaching the life time limit	Replace the lamp. Call ProteinSimple Technical Support for assistance.

Chapter 24:

General Information

Chapter Overview

- Compliance
- Safety Guidelines
- Consumables and Reagents
- Customer Service and Technical Support
- Legal Notices

Compliance

Maurice complies with:

- **UL 61010-1:2001:** Safety requirements for electrical equipment for measurement, control, and laboratory use - Part 1: General requirements (US)
- **EN 61010-1:2010:** Safety requirements for electrical equipment for measurement, control, and laboratory use - Part 1: General requirements (EU)
- **CAN/CSA 22.2 No. 61010-1-04:** Safety requirements for electrical equipment for measurement, control, and laboratory use - Part 1: General requirements (CA)
- **EN 61326-1:2013:** Electrical equipment for measurement, control and laboratory use. EMC Requirements. General requirements (EU)



This device complies with Part 15 of the FCC Rules. Operation is subject to the following two conditions:

1. This device may not cause harmful interference.
2. This device must accept any interference received, including interference that may cause undesired operation.

Note: This equipment has been tested and found to comply with the limits for a Class B digital device, pursuant to Part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference in a residential installation. This equipment generates, uses and can radiate radio frequency energy and, if not installed and used in accordance with the instructions, may cause harmful interference to radio communications. However, there is no guarantee that interference will not occur in a particular installation. If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try to correct the interference by one or more of the following measures:

- Reorient or relocate the receiving antenna.
- Increase the separation between the equipment and receiver.
- Connect the equipment into an outlet on a circuit different from that to which the receiver is connected.
- Consult the dealer or an experienced radio/television technician for help.

Modifications: Any modifications made to this device that are not approved by ProteinSimple, Inc. may void the authority granted to the user by the FCC to operate this equipment.

FCC ID: 2AHGG-MAURICE

Safety Guidelines

!WARNING!

If Maurice is not used as specified by ProteinSimple, overall safety will be impaired.

!WARNING!

If Maurice is damaged and doesn't function properly, stop him safely and contact ProteinSimple Technical Support right away.

!WARNING!

You can't replace or service any parts on Maurice except for the power entry fuse.

!WARNING! SHARPS HAZARD

The capillary inlet of the cartridge may present a potential sharps hazard. Dispose of used cartridges according to your organizations health and safety regulations.

CAUTION

Avoid using Maurice ways not specified by ProteinSimple. Although Maurice has been designed to protect you, this protection may not be effective if he isn't used properly.

Chemical Hazards

!WARNING! CHEMICAL HAZARD

Some chemicals used can be potentially hazardous, and can cause injury or illness.

- Read and understand the Product Inserts and Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials.
- Minimize contact with and inhalation of chemicals. Wear appropriate personal protective equipment when handling chemicals (e.g., safety glasses, gloves, or clothing). For additional safety guidelines, consult the SDS.
- Do not leave chemical containers open.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended on the SDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

Chemical Waste Hazards



!WARNING! BIOHAZARD

Cartridges, sample plates and vials should be handled by procedures recommended in the CDC/NIH manual: Biosafety in Microbiological and Biomedical Laboratories (BMBL). The manual is available from the U.S. Government Printing Office or online at <http://www.cdc.gov/biosafety/publications/bmbl5/>.

Depending on the samples used, the cartridges, plates and vials may constitute a chemical or a biohazard. Dispose of the cartridges, plates and vials in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations. Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste content before you store, handle, or dispose of chemical waste.

- Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste container before you store, handle, or dispose of chemical waste.
- Minimize contact with chemical waste. Wear appropriate personal protective equipment when handling chemicals (e.g., safety glasses, gloves, or clothing).
- Use precaution when emptying waste.
- Dispose of waste in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.

Waste Production and Disposal

cIEF Application on Maurice, Maurice C. and MauriceFlex

Maurice produces approximately 0.75 mL of waste per 100 injections run on a single cIEF cartridge and will contain the following:

- Samples
- Methyl cellulose (~0.5%)
- Carrier ampholytes
- Fluorescence calibration standard
- pI markers
- Sample additives

Additionally, the Anolyte and Catholyte used in the cIEF cartridge will also need to be discarded and replaced after each batch and during cartridge cleanup prior to storage.

- Catholyte: 100 mM sodium hydroxide in 0.1% methyl cellulose, 1.5 mL
- Anolyte: 80 mM phosphoric acid in 0.1% methyl cellulose, 1.5 mL

MauriceFlex cIEF and MauriceFlex Fractionation Application on MauriceFlex

MauriceFlex produces approximately 0.35 mL of waste per injection run on a single cIEF Fractionation Cartridge and will contain the following:

- Water
- Samples
- Methyl cellulose (~0.5%)
- Carrier ampholytes
- Fluorescence calibration standard
- pI markers
- Sample additives

Additionally, the Anolyte used in the cIEF Fractionation Cartridge will also need to be discarded and replaced after each batch and during cartridge cleanup prior to storage.

- Anolyte: 80 mM phosphoric acid in 0.1% methyl cellulose, 2.0 mL

CE-SDS Application on Maurice, Maurice S. and MauriceFlex

Maurice produces approximately 0.75 mL of waste per 48 injections that is contained within the Top Running Buffer vial or Waste Tank. It contains the following:

- Samples
- Conditioning Solution 1 and 2
- Separation Matrix
- Running Buffer
- Wash Solution
- Additives such as β -mercaptoethanol or iodoacetamide.

Waste should be disposed of in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.

Safety Data Sheets




Some chemicals used with Maurice may be listed as hazardous. Warnings are displayed on the labels of all chemicals when hazards exist.

SDSs provide users with safety information needed to store, handle, transport and dispose of the chemicals safely. We recommend updating laboratory SDS records periodically.

Safety Data Sheets for ProteinSimple reagents are available online at <https://www.bio-technne.com/> on the reagent product page or by calling (888) 607-9692. Otherwise, call the chemical manufacturer directly or visit their web site.

Instrument Safety Labels

The following safety labels are located on Maurice. Each label will display a safety alert symbol indicating a potential safety hazard.

Symbol	Description
	Risk of Electric Shock.
	Refer to Maurice User Guide before proceeding.
	Danger of hazardous waste. Use caution in these areas. This warning only applies if using hazardous material. Maurice reagents are not considered hazardous waste. If you are using hazardous materials, please contact your field service representative to place labels in the appropriate locations.

Consumables and Reagents

Maurice CE-SDS Consumables, Kits and Reagents

Item	PN	Description
Maurice CE-SDS PLUS Cartridges (2-pack)	PS-MC02-SP	Cartridges to run CE-SDS PLUS batches, 2 cartridges/pk. For use with Maurice, Maurice S. and MauriceFlex systems only.
Maurice Turbo CE-SDS Cartridges (2-pack)	PS-MC02-TS	Cartridges to run Turbo CE-SDS batches, 2 cartridges/pk. For use with Maurice. Maurice S. and MauriceFlex systems only.
Maurice Turbo CE-SDS Cartridges (1-pack)	PS-MC01-TS	Cartridges to run Turbo CE-SDS batches, 1 cartridge/pk. For use with Maurice, Maurice S. and MauriceFlex systems only.
Maurice glass reagent vials, 2 mL	046-017	Screw-top glass vials, 2 mL. 100/pk. Suitable for use with Maurice, Maurice S. and Maurice C. systems only.
Maurice sample vials with integrated inserts	046-083	Vials with integrated inserts for samples, accommodate 200 µL sample volume. 100/pk. Suitable for use with Maurice, Maurice S., Maurice C. and MauriceFlex systems only.
Maurice CE-SDS Orange Pressure Caps	046-572	Pressure screw tops with o-ring for glass reagent vials (PN 046-017), 12/pk. Suitable for use with Maurice and Maurice S. systems only.

Item	PN	Description
Maurice clear screw caps	046-138	Clear screw caps with round opening, 100/pk. Suitable for use with Maurice, Maurice S., Maurice C systems only.
96-well plates	046-021	96-well plate for high-throughput analysis, 10/pk. Suitable for use with Maurice, Maurice S., Maurice C. and MauriceFlex systems only.
Maurice CE-SDS PLUS Application Kit	PS-MAK03-S	The CE-SDS PLUS Application Kit contains all components required to run CE-SDS PLUS batches. The kit includes 2 CE-SDS PLUS Cartridges, Separation Matrix, Running Buffer, CE-SDS PLUS Sample Buffer, Wash Solution, Conditioning Solutions, Internal Standard, Reagent Vials, Pressure Caps and 96-Well Plates. 200 samples/kit. Molecular weight markers can be ordered separately.
Maurice CE-SDS PLUS Reagent Kit	PS-MRK01-S	The CE-SDS PLUS Reagent Kit contains reagents for running CE-SDS PLUS batches. The kit includes Separation Matrix, Running Buffer, CE-SDS PLUS Sample Buffer, Wash Solution, Conditioning Solutions, and Internal Standard. 200 samples/kit. Molecular weight markers can be ordered separately.
Maurice Turbo CE-SDS Application Kit	PS-MAK01-TS	The Turbo CE-SDS Application Kit contains all components required to run a Turbo CE-SDS batch. The kit includes one (1) Turbo CE-SDS Cartridge, Separation Matrix, Running Buffer, CE-SDS 1X Sample Buffer, CE-SDS PLUS 1X Sample Buffer, Wash Solution, Conditioning Solutions, Internal Standard, Reagent Vials, Pressure Caps and 96-Well Plates. 200 samples/kit. Molecular weight markers can be ordered separately.
Maurice Turbo CE-SDS Reagent Kit	PS-MRK01-TS	The Turbo CE-SDS Reagent Kit contains reagents for running Turbo CE-SDS batches. The kit includes Separation Matrix, Running Buffer, CE-SDS Sample Buffer, CE-SDS PLUS Sample Buffer, Wash Solution, Conditioning Solutions, and Internal Standard. 200 samples/kit. Molecular weight markers can be ordered separately.
Maurice CE-SDS IgG Standard	046-039	Lyophilized antibody standard for 8 runs.
Maurice CE-SDS Molecular Weight Markers	046-432	Lyophilized molecular weight markers (10, 20, 33, 55, 103, 178 and 270 kDa) for 8 runs.
Maurice CE-SDS Separation Matrix	046-386	Separation Matrix for CE-SDS application, 15 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP) or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Running Buffer - Top	046-384	Pre-assembled vial insert containing Top Running Buffer for CE-SDS application. For use with the cartridge insert (046-124) of CE-SDS PLUS cartridge (PS-MC02-SP) on Maurice, Maurice S. and MauriceFlex systems only. 12 vials. Top Running Buffer should be stored at 2-8 °C when not in use.
Maurice CE-SDS Running Buffer - Bottom	046-385	Running Buffer for CE-SDS application, 15 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice Turbo CE-SDS Running Buffer - Bottom	046-579	Buffer for sample preparation for CE-SDS application, 25 mL. For use with Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.

Item	PN	Description
Maurice CE-SDS 1X Sample Buffer	046-012	Buffer for sample preparation for CE-SDS application, 25 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP) or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice and Maurice S. systems only.
Maurice CE-SDS PLUS 1X Sample Buffer	046-567	Buffer for sample preparation for CE-SDS application, 25 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP), or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Wash Solution	046-013	Wash Solution for CE-SDS application, 20 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP) or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Conditioning Solution 1	046-014	Conditioning Solution 1 CE-SDS application, 20 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP) or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Conditioning Solution 2	046-015	Conditioning Solution 2 CE-SDS application, 20 mL. For use with CE-SDS PLUS cartridge (PS-MC02-SP) or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Internal Standard	046-144	Internal Standard for addition to each sample for CE-SDS application. 2 vials/pk. For use with CE-SDS PLUS cartridge (PS-MC02-SP) or Turbo CE-SDS cartridge (PS-MC02-TS, PS-MC01-TS) on Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Cartridge inserts	046-124	Cartridge Inserts for holding the Top Running Buffer vial assembly (046-010) in the CE-SDS PLUS cartridge (PS-MC02-SP). 2 inserts/pk. For use with Maurice, Maurice S. and MauriceFlex systems only.
Maurice CE-SDS Cartridge Cleaning Vial	046-125	Clear cleaning vial for use with the Maurice CE-SDS Cartridge insert (046-124). 2 vials/pk.

Maurice cIEF Consumables, Kits and Reagents

Item	PN	Description
Maurice cIEF Cartridges	PS-MC02-C	Cartridges for cIEF application, 2 cartridges/pk. For use with Maurice, Maurice C. and MauriceFlex systems only.
Maurice glass reagent vials, 2 mL	046-017	Screw-top glass vials, 2 mL. 100/pk. Suitable for use with Maurice, Maurice S. and Maurice C. systems only.

Item	PN	Description
Maurice sample vials with integrated inserts	046-083	Vials with integrated inserts for samples, accommodate 200 µL sample volume. 100/pk. Suitable for use with Maurice, Maurice S., Maurice C. and MauriceFlex systems only.
Maurice cIEF Blue Pressure Caps	046-573	Pressure screw tops with o-ring for glass reagent vials (P/N 046-017), 12/pk. Suitable for use with Maurice, Maurice C. and MauriceFlex systems only.
Maurice clear screw caps	046-138	Clear screw caps with round opening, 100/pk. Suitable for use with Maurice, Maurice S., Maurice C systems only.
96-well plates	046-021	96-well plate for high-throughput analysis, 10/pk. Suitable for use with Maurice, Maurice S., Maurice C. and MauriceFlex systems only.
Maurice cIEF Method Development Kit	PS-MDK01-C	The cIEF Method Development Kit provides all the reagents and instructions to help you develop fast and robust cIEF methods on Maurice, Maurice C. and MauriceFlex systems. The kit includes a Method Development Guide as well as a selection of reagents required for cIEF method development on the system. This kit includes a Fluorescence Calibration Standard, System Suitability Kit, Anolyte, Catholyte, Methyl Cellulose, five types of Ampholytes (Pharmalyte pH ranges 3-10, 2.5-5, 5-8 and 8 to 10.5 and Servalyte pH range 2-9), nine pI Markers (3.38, 4.05, 5.85, 6.14, 7.05, 8.40, 9.50, 9.99 and 10.17) and additives (SimpleSol Protein Solubilizer, PN 046-574, lyophilized urea and arginine). 30 samples/kit. The expiration date for this kit is 12 months from date of manufacture. For use with the cIEF cartridge (PS-MC02-C) on Maurice, Maurice C. and MauriceFlex systems only.
Maurice cIEF Chemical Test Kit	046-036	The cIEF Chemical Test Kit is designed to confirm the overall performance of the Maurice, Maurice C. and MauriceFlex. systems with the cIEF cartridge (PS-MC02-C) on 8 separate occasions. The kit contains Anolyte, Catholyte, Methyl Cellulose, Fluorescence Calibration Standard and a System Suitability kit. Each time, the performance can be checked up to 24 hours. The expiration date for this kit is 12 months from date of manufacture. 8 runs/kit. For use with the cIEF cartridge (PS-MC02-C) on Maurice, Maurice C. and MauriceFlex systems only.
Maurice cIEF System Suitability Kit	046-044	The cIEF System Suitability Kit is used to run a control test for cIEF application on Maurice, Maurice C. and MauriceFlex systems on 8 separate occasions. Each time, the performance can be checked up to 24 hours. The kit contains a vial of lyophilized Peptide Panel and a vial of System Suitability Test mix. 8 runs/kit. The expiration date for this kit is 12 months from date of manufacture. For use with the cIEF cartridge (PS-MC02-C) on Maurice, Maurice C. and MauriceFlex systems only.
Maurice cIEF Fluorescence Calibration Standard	046-025	Fluorescence Calibration Standard for cIEF application, 5.5 mL. For use with Maurice, Maurice C. and MauriceFlex systems only.
SimpleSol Protein Solubilizer	046-575	SimpleSol is a protein solubilizer, 200 mL. This reagent is more stable than urea, eliminating the need to prepare urea fresh every day. For use with Maurice, Maurice C., MauriceFlex and iCE3 systems.

Item	PN	Description
0.5% Methyl Cellulose Solution	102505	0.5% Methyl Cellulose Solution, 100 mL. Use this concentration for the wash cycle between injections. The solution is filtered to ensure consistent viscosity to coat the capillary lumen to minimize electroosmotic flow (EOF). Conveniently packaged in 2 x 100 mL bottles. The expiration date for this kit is 12 months from date of manufacture. Suitable for use with Maurice, Maurice C., MauriceFlex, iCE3 and iCE280.
1% Methyl Cellulose Solution	101876	1% Methyl Cellulose Solution, 100 mL. This solution is used to prepare samples for cIEF applications. The expiration date for this kit is 12 months from date of manufacture. Suitable for use with Maurice, Maurice C., MauriceFlex, iCE3 and iCE280.
iCE Electrolyte Kit	102506	These 100 mL Anolyte and Catholyte solutions in 0.1% MC are used to fill the electrolyte tanks on Maurice cIEF cartridge as well as FC and HT cartridges. The labels and bottles are color coded to improve safety. The kit contains 2 x 100 mL bottles. The expiration date for this kit is 12 months from date of manufacture. Suitable for use with Maurice, Maurice C., MauriceFlex, iCE3 and iCE280.
Maurice cIEF pI Marker - 3.38	046-028	Maurice cIEF pI Marker - 3.38, 210 µL, lyophilized.
Maurice cIEF pI Marker - 4.05	046-029	Maurice cIEF pI Marker - 4.05, 210 µL, lyophilized.
Maurice cIEF pI Marker - 5.85	046-030	Maurice cIEF pI Marker - 5.85, 210 µL, lyophilized.
Maurice cIEF pI Marker - 6.14	046-031	Maurice cIEF pI Marker - 6.14, 210 µL, lyophilized.
Maurice cIEF pI Marker - 7.05	046-032	Maurice cIEF pI Marker - 7.05, 210 µL, lyophilized.
Maurice cIEF pI Marker - 8.40	046-033	Maurice cIEF pI Marker - 8.4, 210 µL, lyophilized.
Maurice cIEF pI Marker - 9.50	046-047	Maurice cIEF pI Marker - 9.50, 210 µL, lyophilized.
Maurice cIEF pI Marker - 9.99	046-034	Maurice cIEF pI Marker - 9.99, 210 µL, lyophilized.
Maurice cIEF pI Marker - 10.17	046-035	Maurice cIEF pI Marker - 10.17, 210 µL, lyophilized.
Maurice cIEF electrolyte tank caps	046-123	Electrolyte tank caps, 20 mm. Red is used for the Anolyte tank and the grey cap is used for the Catholyte tank. 5 pairs/pk.
Electrolyte Pipette	101788	Pipettes with soft tips for adding Anolyte and Catholyte into the electrolyte tanks in the cIEF cartridge. 10/pk.

MauriceFlex cIEF Fractionation Consumables, Kits and Reagents

Item	PN	Description
MauriceFlex cIEF Fractionation Cartridge	PS-MC02-F	The MauriceFlex cIEF Fractionation Cartridge is designed for icIEF-based fractionation on the MauriceFlex system. Each pack contains 2 Cartridges, and each Cartridge supports up to 15 fractionation runs. Suitable for use with MauriceFlex systems only. Not for analytical icIEF analysis.
MauriceFlex crimp top glass reagent vials, 2 mL, 100/pkg	110-0019	Glass vial, crimp top, 2mL, 100/pkg. Reagent vials for MauriceFlex cIEF fractionation use. Suitable for use with MauriceFlex systems only.
MauriceFlex glass vials with insert, 0.3 mL, 15/pkg	110-0018	Glass vials, 2mL, with 0.3 mL insert, 15/pkg. For MauriceFlex cIEF fractionation use. Suitable for use with MauriceFlex systems only.
96-well plates	046-021	96-well plate for high-throughput analysis, 10/pk. Suitable for use with Maurice, Maurice S., Maurice C. and MauriceFlex systems only.
MauriceFlex cIEF Fractionation Method Development Kit	PS-MDK01-F	The MauriceFlex cIEF Fractionation Method Development Kit provides all the reagents and instructions to help you develop both analytical icIEF method as well as the icIEF fractionation method. The kit includes a Maurice cIEF Method Development Kit (PS-MDK01-C) and additional reagents and consumables for cIEF fractionation method development. A single MauriceFlex cIEF Fractionation Cartridge is also included to support up to 15 icIEF fractionation runs. The expiration date for this kit is 12 months from date of manufacture. For use with the cIEF cartridge (PS-MC02-C) for analytical icIEF analysis. The kit is to be ordered with a new MauriceFlex system.
MauriceFlex cIEF Fractionation Reagent Kit	PS-MRK01-F	The MauriceFlex cIEF Fractionation Reagent Kit contains Fluorescence Calibration Standard, Anolyte, Catholyte, Methyl Cellulose, arginine and ammonium acetate to support up to 15 cIEF fractionation runs on the MauriceFlex system with a MauriceFlex cIEF Fractionation Cartridge.
MauriceFlex cIEF Fractionation Application Kit	PS-MAK01-F	The MauriceFlex cIEF Fractionation Application Kit contains one MauriceFlex cIEF Fractionation Reagent Kit (PS-MRK01-F), one cIEF Fractionation Cartridge, Reagent Vials, and 96-well Plates to support up to 15 icIEF fractionation runs. For initial icIEF fractionation method development on a new MauriceFlex system, order MauriceFlex cIEF Fractionation Method Development Kit (PS-MDK-01) instead.
2 M Ammonium Acetate, 350 µL	046-580	Ammonium acetate aqueous solution, 2 M, 350 µL. Stock for preparing chemical mobilizer for MauriceFlex cIEF fractionation use.

Customer Service and Technical Support

Need pricing information or want to know who your sales rep is? Our Customer Service team can help.

Email: orders@proteinsimple.com

Telephone: 1-408-510-5500, option 1

Toll-free (US and Canada): 1-888-607-9692, option 1

Fax: 1-408-520-4831

Have product-related questions? Ping our Tech Support group, they'll be happy to help!

Use our online technical support request

Email: support@proteinsimple.com

Telephone: 1-408-510-5500

Toll-free (US and Canada): 1-888-607-9692, option 3

Web

www.bio-techne.com

Address

ProteinSimple
3001 Orchard Parkway
San Jose, CA 95134
USA

Legal Notices

NOTE: Read the Legal Notices carefully before using Maurice.

Maurice Disclaimer of Warranty

EXCEPT AS EXPRESSLY PROVIDED IN ANY ProteinSimple SOFTWARE LICENSE AGREEMENT OR QUOTATION, THE PRODUCTS SOLD AND SERVICES PROVIDED BY ProteinSimple ARE PROVIDED ON AN “AS IS” AND “AS AVAILABLE” BASIS WITHOUT WARRANTY OF ANY KIND. ProteinSimple AND ITS SUPPLIERS DO NOT WARRANT THE SECURITY, PRIVACY, OR ACCURACY OF ANY DATA PROVIDED VIA THE PRODUCTS OR SERVICES, AND YOU AGREE THAT THE USE OF ANY SUCH DATA BY YOU IS AT YOUR SOLE RISK. TO THE MAXIMUM EXTENT ALLOWED UNDER APPLICABLE LAW, ProteinSimple AND ITS SUPPLIERS DISCLAIM ANY AND ALL WARRANTIES, WHETHER EXPRESS, IMPLIED OR STATUTORY, INCLUDING, WITHOUT LIMITATION, ANY WARRANTY OF MERCHANTABILITY, TITLE, NON-INFRINGEMENT, OR FITNESS FOR A PARTICULAR PURPOSE.

Compass Software and Compass Authorization Server for iCE License Agreement

IMPORTANT - PLEASE READ CAREFULLY THE TERMS OF THIS COMPASS SOFTWARE AND AUTHORIZATION SERVER LICENSE AGREEMENT (“AGREEMENT”). BY CLICKING ON THE “I AGREE” BUTTON, (1) YOU ACKNOWLEDGE THAT YOU HAVE READ, UNDERSTAND, AND AGREE TO BE BOUND BY THIS AGREEMENT AND (2) YOU REPRESENT THAT YOU HAVE THE AUTHORITY TO ENTER INTO THIS AGREEMENT, PERSONALLY OR IF YOU HAVE NAMED A COMPANY AS CUSTOMER, ON BEHALF OF THAT COMPANY (YOU OR ANY SUCH COMPANY, THE “CUSTOMER”), AND TO BIND THE CUSTOMER TO THE TERMS OF THIS AGREEMENT. IF YOU DO NOT AGREE TO ALL TERMS AND CONDITIONS OF THIS AGREEMENT, OR IF YOU DO NOT HAVE SUCH AUTHORITY, YOU SHOULD CLICK ON THE “CANCEL” BUTTON TO DISCONTINUE THE DOWNLOAD OF THE LICENSED SOFTWARE.

1. Definitions

- 1.1. **“Authorized Use Parameters”** means the following usage restrictions, which restrict the operation of the Licensed Software to a particular set of conditions: Customer shall (a) limit simultaneous use of the Licensed Software to a maximum of ten (10) Authorized Users; and (b) use the Licensed Software only in connection with the accompanying System purchased by Customer pursuant to the System Quotation and located at the Site.
- 1.2. **“Authorized User”** means one (1) User who initiates the execution of the Licensed Software and/or interacts with or directs the Licensed Software in the performance of its functions. Multiple Authorized Users may work simultaneously with one installation of the Licensed Software, as on a server, or they may each have their own installation on single-user machines, or a mix of these, provided that in all cases the total number of simultaneous Users does not exceed the applicable Authorized Use Parameters.
- 1.3. **“Company”** means ProteinSimple.
- 1.4. **“Documentation”** means Company’s then-current manuals, guides, and on-line help pages, if any, applicable to the Licensed Software and made generally available by Company to its customers.
- 1.5. **“Enterprise”** means those organizations that have Internet addresses located at top level and second-level domain names set forth in the System Quotation.
- 1.6. **“Error”** means a reproducible error in the Licensed Software that prevents such Licensed Software from operating substantially in accordance with its Documentation.
- 1.7. **“Executable Code”** means the fully compiled binary version of Licensed Software that can be executed by a computer and used by an end user without further compilation.
- 1.8. **“Intellectual Property Rights”** means all copyrights, trade secrets, patents, patent applications, moral rights, contract rights, and other proprietary rights, but specifically excluding any trademarks or service marks.

- 1.9. **“Licensed Software”** means the Compass software program in Executable Code form, and any Updates that Company makes available to Customer in accordance with this Agreement.
- 1.10. **“Site”** means the facility or campus set forth in the System Quotation.
- 1.11. **“System”** means the proprietary NP1000, NP100, Simon, Sally, Peggy, Wes, Sally Sue, Peggy Sue and Maurice protein analysis system or any future model or successor thereto that is provided to Customer by Company pursuant to a separate agreement between the parties (the “System Quotation”).
- 1.12. **“Update”** means those releases of the Licensed Software that Company provides to customers to correct Errors, fix bugs, or create minor improvements, incremental features, or enhancements of existing features which Company designates by a change in the number to the right of the first or second decimal point. Updates do not include those releases of the Licensed Software that provide substantial new features or additional functionality which Company designates by a change in the number to the left of the first decimal point.
- 1.13. **“User”** means any individual that has an e-mail address within the Enterprise.

2. License and Restrictions

- 2.1. **License Grant.** Subject to the terms and conditions of this Agreement and the payment of the required fees set forth in the System Quotation, Company grants to Customer a nontransferable, nonexclusive, royalty-free, revocable, worldwide license (without the right to sublicense) to (a) install the Licensed Software on any computer located at any Site; (b) use, execute, and display the Licensed Software, in Executable Code form only; and (c) copy the Licensed Software and Documentation, solely as necessary to support Authorized Users; in each of the foregoing, solely in accordance with the Documentation and the Authorized Use Parameters. Customer agrees that it will comply with the Authorized Use Parameters.
- 2.2. **License Restrictions.** Customer acknowledges that the Licensed Software and its structure and organization constitute valuable trade secrets of Company. Accordingly, the license granted in this Agreement is subject to the following restrictions: Customer and its Authorized Users (a) may not reverse engineer, disassemble, decompile, or otherwise attempt to derive the source code of Licensed Software; (b) may not modify, adapt, alter, translate, or create derivative works from the Licensed Software; (c) may not merge the Licensed Software with other software; (d) may not use the Licensed Software in any service bureau or time-sharing arrangement, license, sell, rent, lease, transfer, assign, distribute, host, outsource, disclose, or otherwise commercially exploit or make the Licensed Software or Documentation available to any third party; (e) shall only make that number of exact copies of the Licensed Software and Documentation as delivered by Company that are necessary to support Customer’s use of the Licensed Software in accordance with this Agreement; (f) shall include any titles, trademarks, and copyright and restricted rights notices that are included on or in the Licensed Software as delivered by Company on and in any copies of the Licensed Software that it makes; and (g) shall ensure that Customer’s use of the Licensed Software does not exceed the scope of the license that Customer has purchased pursuant to this Agreement.
- 2.3. **Open Source Software.** Certain items of independent, third-party code may be included in the Licensed Software that are subject to open source licenses (“Open Source Software”). Such Open Source Software is licensed under the terms of the license that accompanies such Open Source Software. Nothing in this Agreement limits Customer’s rights under, or grants Customer rights that supersede, the terms and conditions of any applicable end user license for such Open Source Software. In particular, nothing in this Agreement restricts Customer’s right to copy, modify, and distribute such Open Source Software that is subject to the terms of such open source licenses.
- 2.4. **Ownership.** Company reserves all rights not expressly granted to Customer in this Agreement. Without limiting the generality of the foregoing, Customer acknowledges and agrees that, except as expressly set forth in this Agreement, Company and its suppliers retain all Intellectual Property Rights, title and interest in and to the Licensed Software and Documentation.

3. Support and Maintenance Services

- 3.1. **Services.** Subject to Customer’s payment of the Services fees, as set forth in the System Quotation, and to the terms and conditions herein, Company will use commercially reasonable efforts to provide to Customer the following support and maintenance services (the “Services”) for the Licensed Software: (a) Company will answer technical questions concerning functions and features of the Licensed Software; (b) Company will provide Error verification, analysis and corrective efforts for the Licensed Software; and (c) Company will provide, without charge, Updates of the software released during the term of this Agreement. Customer will be responsible for providing, in a manner consistent with good industry practice, all Services to Users. Customer acknowledges that Company may not be able to correct all reported Errors. Any Update of the Licensed Software will be deemed part of the Licensed Software and Customer will use such Updates in accordance with the requirements and obligations in this Agreement.

- 3.2. **Service Conditions.** Company's obligation to provide the Services is conditioned on Customer: (a) notifying Company of any Error within a reasonable period of time; (b) providing Company all information relating to the Error; (c) providing access to the Licensed Software and Customer's facility where the Licensed Software is located and informing Company of any potential hazards which may be encountered while servicing the Licensed Software. Customer may contact Company via telephone at 1-888-607-9692 or e-mail at support@proteinsimple.com during the hours of 8 a.m. (Pacific Time) and 5 p.m. (Pacific Time) Monday through Friday, excluding holidays, to report any Error. A list of standard holidays will be provided to Customer upon request. Company shall have the right to determine in its sole discretion what corrective action Company will perform to support the Licensed Software. Company may subcontract the Services to a third party contractor provided that Company will be responsible for the third party contractor's compliance with this Agreement.
- 3.3. **Service Exclusions.** Company will not be obligated to provide the Services if (a) Company determines that an Error is caused by malfunction of any hardware (other than malfunction of the System) or third party software used with the Licensed Software; or (b) Customer has failed to incorporate the latest Update previously released to Customer.

4. Warranty

- 4.1. **Licensed Software Warranty.** Company warrants that the Licensed Software, as properly installed, and under normal use, will perform substantially in accordance with its Documentation during the Warranty Period. The "Warranty Period" for the Licensed Software begins on date Customer downloads the Licensed Software and ends twelve (12) months thereafter.
- 4.2. **Remedy.** If Customer notifies Company in writing during the Warranty Period of an Error, Company will, at its expense and as its sole obligation for any breach of the foregoing warranty, use commercially reasonable efforts to correct the Error or replace the Licensed Software. Any Error correction or replacement of the Licensed Software will not extend the original Warranty Period. The warranty and the remedies provided above will not apply to the Licensed Software if (a) Company determines that an Error is caused by accident, abuse, misuse, negligence, fire, earthquake, flood, other force majeure event, failure of electrical power, the use of unauthorized products, or unauthorized repairs or modifications; (b) Company determines that an Error is caused during or as a result of delivery; (c) a problem arises from or is based on Company's compliance with Customer's specifications; or (d) Company determines that an Error is caused by malfunction of any hardware (other than malfunction of the System) or third party software used with the Licensed Software.
- 4.3. **Disclaimer.** THE WARRANTIES ABOVE ARE EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER EXPRESS, IMPLIED OR STATUTORY, INCLUDING WITHOUT LIMITATION THE IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, TITLE, AND NONINFRINGEMENT.

5. **Limitation of Liability.** NEITHER COMPANY NOR ITS SUPPLIERS SHALL BE RESPONSIBLE OR LIABLE WITH RESPECT TO ANY SUBJECT MATTER OF THIS AGREEMENT OR TERMS OR CONDITIONS RELATED THERETO UNDER ANY CONTRACT, NEGLIGENCE, STRICT LIABILITY OR OTHER THEORY (A) FOR LOSS OR INACCURACY OF DATA, LOSS OF PROFITS OR COST OF PROCUREMENT OF SUBSTITUTE GOODS, SERVICES OR TECHNOLOGY, OR (B) FOR ANY INDIRECT, INCIDENTAL OR CONSEQUENTIAL DAMAGES INCLUDING, BUT NOT LIMITED TO LOSS OF REVENUES AND LOSS OF PROFITS. COMPANY'S AGGREGATE CUMULATIVE LIABILITY HEREUNDER SHALL NOT EXCEED THE GREATER OF FIVE HUNDRED DOLLARS (\$500.00).

6. Term and Termination

- 6.1. **Term of Agreement.** The Agreement is effective on the date Customer downloads the Licensed Software and shall remain in effect until terminated by either party as provided in this section.
- 6.2. **Termination For Material Breach.** Either party may terminate this Agreement upon written notice if the other party materially breaches this Agreement and fails to cure such breach within thirty (30) calendar days following receipt of written notice from the other party specifying the breach in detail. Notwithstanding the foregoing, Company may immediately terminate this Agreement and all licenses granted hereunder if Customer breaches Section 2 (License and Restrictions) hereof or upon termination of the System Quotation. The foregoing rights of termination are in addition to any other rights and remedies provided in this Agreement or by law.
- 6.3. **Effect of Termination.** Upon termination of this Agreement (or termination or expiration of any license granted hereunder), all rights of Customer to use the Licensed Software and Documentation will cease and (a) all license rights granted under this Agreement will immediately terminate and Customer shall promptly stop all use of the Licensed Software and Documentation; (b) all Services will terminate immediately; (c) Customer shall promptly erase all copies of the Licensed Software from Customer's computers, and destroy all copies of the Licensed Software and Documentation on tangible media in Customer's possession or control or return such copies to Company; and (d) upon request by Company, Customer shall certify in writing to Company that it has returned or destroyed such Licensed Software and Documentation. The parties' rights and obligations under Sections 1 (Definitions), 2.4 (Ownership), 4.3 (Disclaimer), 5 (Limitation of Liability), 6 (Term and Termination), and 7 (General) shall survive termination of this Agreement.

7. General

- 7.1. **Assignment.** This Agreement and Customer's rights hereunder may not be assigned to any third party by Customer except with the prior written approval of Company. Any attempted assignment of this Agreement or any rights or obligations hereunder will be null and void.
- 7.2. **Governing Law.** This Agreement is made in, governed by, and shall be construed in accordance with the laws of the State of California, without regard to any conflicts of law principles that would result in application of laws of any other jurisdiction. The United Nations Convention on Contracts for the International Sale of Goods does not apply to this contract. Any legal action or other legal proceeding relating to this contract or the enforcement of any provision of this contract must be brought in any state or federal court located in Santa Clara County, California. Customer and Company expressly and irrevocably consents and submits to the jurisdiction of such courts.
- 7.3. **Injunctive Relief.** Customer acknowledges that the Licensed Software contains valuable trade secrets and proprietary information of Company, that any actual or threatened breach of this Agreement will cause harm to Company for which monetary damages would be an inadequate remedy, and that injunctive relief is an appropriate remedy for such breach.
- 7.4. **Modifications.** Company reserves the right to change the terms and conditions of this Agreement or its policies relating to the Licensed Software at any time. Company will notify Customer of any material changes to this Agreement by sending Customer an e-mail to the last e-mail address Customer provided to Company or by prominently posting notice of the changes on Company's website. Any material changes to this Agreement will be effective upon the earlier of thirty (30) calendar days following Company's dispatch of an e-mail notice to Customer or thirty (30) calendar days following Company's posting of notice of the changes on Company's website. These changes will be effective immediately for new users of our Licensed Software. Please note that at all times Customer is responsible for providing Company with its most current e-mail address. In the event that the last e-mail address that Customer has provided Company is not valid, or for any reason Company is not capable of delivering to Customer the notice described above, Company's dispatch of the e-mail containing such notice will nonetheless constitute effective notice of the changes described in the notice. If Customer does not agree with the changes to this Agreement, Customer must notify Company prior to the effective date of the changes that Customer wishes to terminate its license to the Licensed Software. Continued use of the Licensed Software, following notice of such changes, shall indicate Customer's acknowledgement of such changes and agreement to be bound by the terms and conditions of such changes.
- 7.5. **Severability.** In the event any provision of this Agreement is held to be invalid or unenforceable, the remaining provisions of this Agreement will remain in full force.
- 7.6. **Waiver.** The waiver by either party of any default or breach of this Agreement shall not constitute a waiver of any other or subsequent default or breach.
- 7.7. **Export.** Customer agrees not to export, reexport, or transfer, directly or indirectly, any U.S. technical data acquired from Company, or any products utilizing such data, in violation of the United States export laws or regulations.
- 7.8. **Force Majeure.** Company shall not be liable, directly or indirectly, for any delay or failure in performance of any obligation under this Agreement, including any delivery obligation, where such delay or failure arises or results from a cause beyond Company's reasonable control, or beyond the reasonable control of Company's suppliers or contractors, including, but not limited to strike, boycott or other labor disputes, embargo, governmental regulation, inability or delay in obtaining materials, acts of God, war, earthquake, fire, or flood. In the event of such force majeure, the time for delivery or other performance will be extended for a period equal to the duration of the delay caused thereby, provided that Company notifies Customer of the nature and duration of such force majeure event.
- 7.9. **Entire Agreement; Notice.** This Agreement constitutes the complete agreement between the parties and supersedes all prior or contemporaneous agreements or representations, written or oral, concerning the subject matter of this Agreement. Except as otherwise expressly provided in this Agreement, any modifications of this Agreement must be in writing and agreed to by both parties. Company may provide any notice to Customer by e-mail. Customer may provide notice to Company by sending an e-mail to info@proteinsimple.com or a letter by United States mail to ProteinSimple, 3040 Oakmead Village Drive, Santa Clara, CA 95051, or to such other address as Company may specify in writing by posting the new address on the Company website.
- 7.10. **Relationship of the Parties.** The parties are acting hereunder as independent contractors and not as partners, agents, fiduciaries, or joint venturers. Neither party has the power or authority represent, act for, bind, or otherwise create or assume any obligation on behalf of the other party.

Where Science Intersects Innovation™

Bio-Techne® | R&D Systems™ Novus Biologicals™ Tocris Bioscience™ ProteinSimple™ ACD™ ExosomeDx™ Asuragen®



Contact Us

Global info@bio-techne.com bio-techne.com/find-us/distributors

North America TEL 800 343 7475

Europe | Middle East | Africa TEL +44 (0)1235 529449

China info.cn@bio-techne.com TEL +86 (21) 52380373

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners.

bio-techne®