

Maurice cIEF

Quick Reference Guide



PREPARE YOUR REAGENTS AND SAMPLE

1. Prepare your samples and batch reagents using the instructions in the Maurice User Guide or Method Development Kit.

INSTALL YOUR CARTRIDGE

1. Take the cartridge out of its packaging. Save the packaging, you'll need it later.
NOTE: Make sure the tip of the cartridge inlet doesn't come in contact with any surfaces. A damaged inlet may compromise the cartridge and cause failed injection.
2. Place the cartridge on a flat surface with its electrolyte tanks facing up.
3. Remove the red and grey stoppers from each tank.
4. Add 2 mL of catholyte to the OH⁻ tank (white port) and 2 mL of anolyte to the H⁺ tank (red port).
NOTE: Make sure you don't overfill the electrolyte tanks.



5. 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 lint-free laboratory wipe.
6. Open Maurice's door by touching the metal plate on top of the door.
7. Slide the cartridge into the cartridge slot. The lights around the cartridge will turn from orange to blue, and you'll feel a locking mechanism engage when the cartridge is correctly installed.

PUT YOUR REAGENTS AND SAMPLES IN MAURICE

1. Depending on how you prepped your samples, place the metal insert for either the sample vials or a 96-well plate in Maurice.
2. Place your prepped batch reagents in Rows P and N.
 - **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**

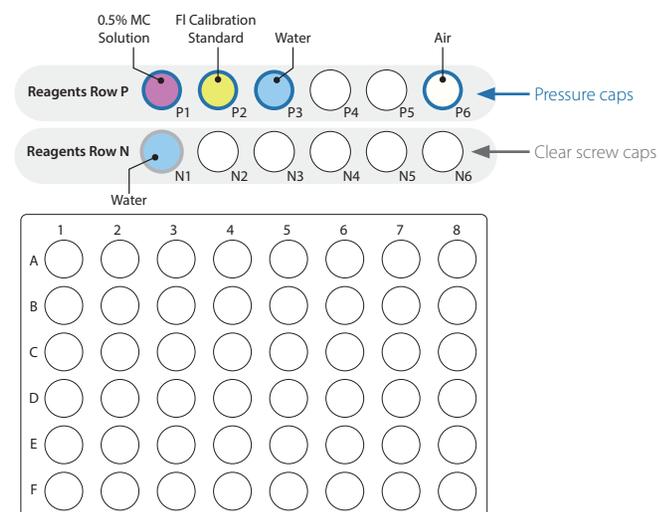


Figure 1. Sample and reagent vial placement when using sample vials.

3. Put your prepared samples into vials or a 96-well plate and place them in the metal insert.

CREATE AND START YOUR BATCH

1. Launch Compass for iCE. The Maurice icon on the top left confirms Maurice is connected to the computer.



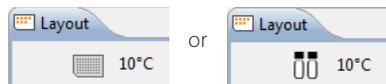
2. Click the **Batch** tab.



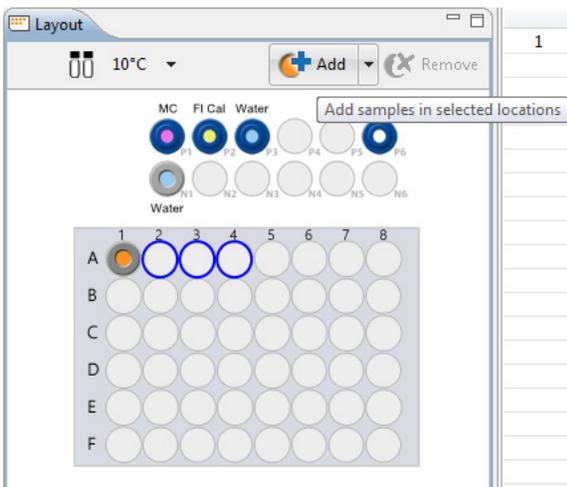
3. Set up your batch:

In the File menu, click **New Batch**. If your Maurice runs both size and charge, select **Maurice cIEF**.

In the **Layout** window, click to toggle the **Plate** icon to the **Vial** icon if you prepped your samples in vials.



Adding a sample: Highlight the sample location in the **Layout** window and click **Add**.



Running multiple methods in the same batch: Click **New** in the **Methods** window, and then enter your method parameters for each method.

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		

Adding sequential replicate injections: Highlight the injection in the **Injections** window and click **Replicate**.

	Sample ID	Location	Method	Notes
1	Sample 1	A1	Method1	
2	Sample 2	A2	Method1	
3	Sample 2	A2	Method1	
4	Sample 3	A3	Method1	

Adding an injection at the end of the batch: Highlight the injection in the **Injections** window and click **Add**. Then use the pulldown menu to assign the method. You can also copy/paste injections or click on the sample position in the **Layout** window and click **Add**.

	Sample ID	Location	Method	
1	Sample 1	A1	Method1	
2	Sample 2	A2	Method1	
3	Sample 3	A3	Method1	
4	Sample 4	A4	Method1	
5	Sample 2	A2	Method1	

4. Click **Start**.



AT THE END OF THE BATCH

As soon as possible:

1. Open Maurice's door.
2. Remove your reagent vials and samples.
3. Remove the cartridge.
4. Clean-up and store your cartridge using the instructions in the Maurice User Guide or cIEF Cartridge Product Insert.

Note: A cartridge is guaranteed for 100 injections and a maximum of 20 batches.