

August 18, 2023

Be happy, extend the life of your cartridge, and save your sample - all for free!

Have you heard about the new cartridge maintenance features included in Compass for iCE version 4.0 software for Maurice? In addition to support for the new Turbo CE-SDS and MauriceFlex cIEF Fractionation cartridges, Compass for iCE v4.0 features new and improved washing and purging steps to reduce maintenance-associated cartridge issues. Users will be notified with various alert messages based on the cartridge cleanup status.

Staying on top of regular and prompted cartridge maintenance has many benefits, such as:

- Reduced sample waste
- Prolonging the life of the cartridge
- Improving the quality and reproducibility of data
- Better instrument uptime by preventing cartridge failures
- Avoiding costly cartridge replacements

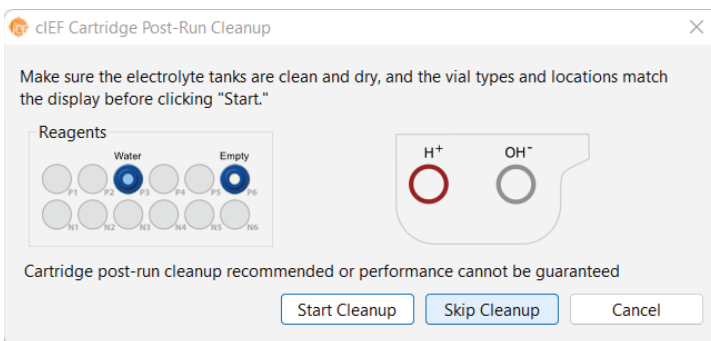
Best of all, **Compass for iCE 4.0 is available free of charge.**

In this software update, a new standalone post-run cleanup method had been added for cIEF cartridges.

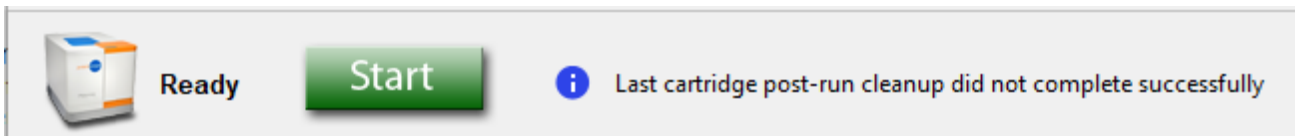
After a batch or Cartridge Purge run, users will now be prompted with a new message recommending a post-run cleanup.



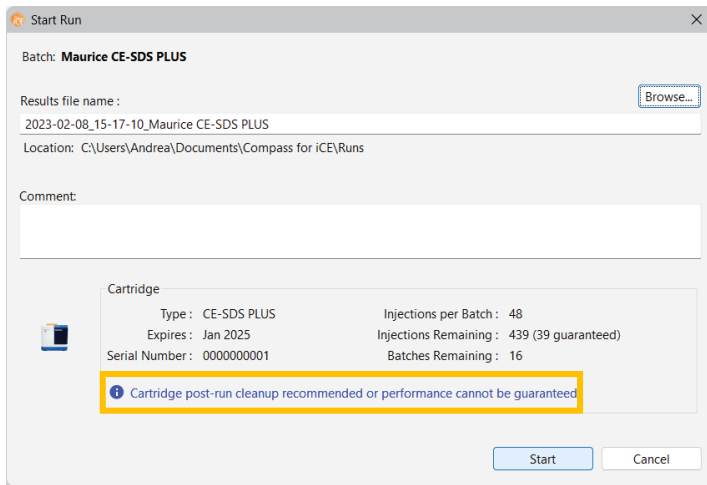
Users have the option to select **Skip Cleanup**; however, this may negatively impact cartridge performance and is not recommended.



If a successful **Cartridge Post-Run Cleanup** is not performed or the clean-up did not run to completion, the following message will be displayed upon insertion of the cartridge into the instrument.

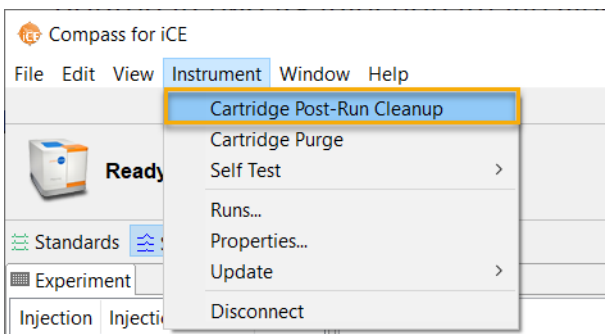


If users select **Start** without performing a cleanup, a secondary message will appear to recommend a cleanup be performed and the decision will be logged in the Injection Report.

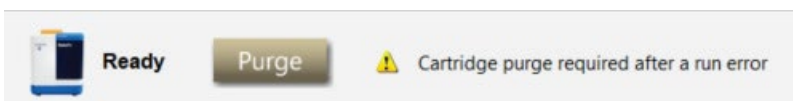


Model	Maurice OBM
Instrument S/N	kf1147
Software Version	Compass for iCE 3.0.0, Build ID: 0124
Firmware Version	4.1.2021.12.09.23.23.49.55b8e524f
Tray Type	96-well plate
Cartridge Type	Turbo CE-SDS™
Cartridge S/N	8211213001
Cartridge Expiration	Dec 2022
Injections Remaining	(4 guaranteed)
Batches Remaining	20
Warning	Cartridge post-run cleanup not performed after previous run.

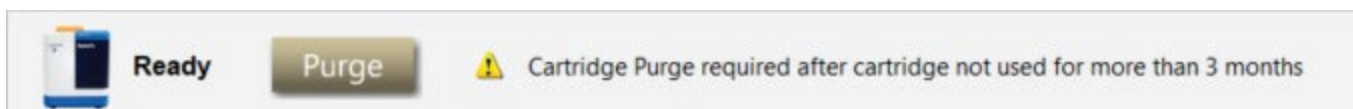
The above-mentioned post-run cleanup methods can be accessed from the **Instrument** drop-down menu bar and by selecting **Cartridge Post-Run Cleanup**.



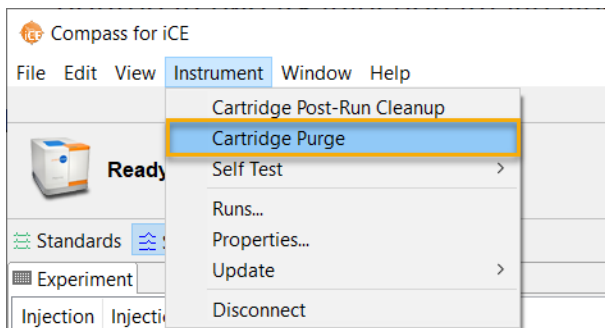
If an error occurred during the run, the following message will be displayed:



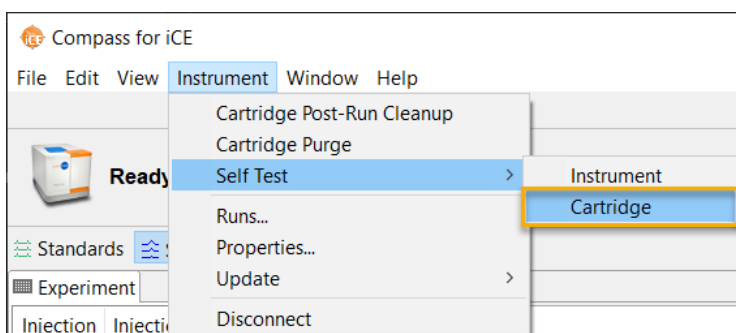
Lastly, if a cartridge has not been used for more than 3 months, a message will be displayed, recommending a cartridge purge be completed to ensure no clogging has occurred while the cartridge was in storage.



Cartridge purge methods can be accessed from the Instrument drop-down menu bar and by selecting the Cartridge Purge option.



Finally, as implemented in Compass for iCE v3.0, users are able to perform a Cartridge Self-Test to ensure the cartridge is functioning properly. You can view current and past self-test results anytime by selecting **Instrument > Properties** and clicking **Test Log**.



Following these new cIEF cartridge cleanup procedures will reduce cartridge failures and downtime. For more information on how to upgrade to Compass for iCE v4.0, please contact your local sales representative, or visit us at www.bio-techne.com.

For any additional questions or concerns, please contact ProteinSimple Technical Support at (US) (888) 607-9692 option 3, or support@proteinsimple.com. For customers in Europe, please contact Technical Support at +44 1235 529449, or Instrument.Support.EMEA@bio-techne.com.

Sincerely,

Yasef Khan

Product Manager Specialist, Biologics Commercial
ProteinSimple, a Bio-Techne brand

WHERE SCIENCE INTERSECTS INNOVATION™

3001 Orchard Parkway, San Jose, CA 95134 | TEL: 1.888.607.9692 | 1.408.510.5500 | proteinsimple.com