

Maurice icIEF 400 Cartridge Handling Guide



Introduction

Handling, preparation, and cleanup of the icIEF 400 cartridge is very similar to that of the cIEF cartridge. Some additional recommendations are presented below to ensure successful usage of the icIEF 400 cartridges.

Prepare your cartridge

Unlike the cIEF cartridge, the capillary in the icIEF 400 cartridge is exposed to air, within the cartridge's optical window.



- Before every run, use filtered, compressed air to clean the cartridge's optical window to reduce particle collection.
- NOTE:** The recommended air output at the nozzle should be around 1–1.5 SCFM with a minimum distance of at least 1 inch away from the cartridge's optical window.

- Apply the air at a side angle towards the top and bottom sides of the cartridge.



- Do not apply the air directly from above of the cartridge



- Do not touch the cartridge's optical window or use Kimwipes, water, etc., to clean it.
- Ensure that objects such as rubber stoppers don't fall onto the exposed capillary.

When pipetting liquid into and out of the cartridge:

- Place the cartridge on a flat surface with the electrolyte tanks facing up to ensure liquid does not drip onto the cartridge's optical window.



- Avoid placing the pipette or aspirator over the cartridge's optical window to avoid dripping onto the exposed capillary.



If you're using Compass for iCE v3.0.0 or higher, go to the next step. If you're using an earlier version of Compass for iCE, skip to step 9.

6. Put the stoppers back on the tanks and install the cartridge in Maurice.
7. In the Compass main menu, select **Instrument** and click **Cartridge Post-Run Cleanup**. It'll only take 7 minutes.
8. Remove the cartridge and reagent vials from Maurice.
9. Leave the stoppers off to allow the tanks to air dry.
10. Store the icIEF 400 cartridge and stoppers back in its original packaging and store at room temperature.

Other design differences

The icIEF 400 cartridge waste tank sponge is not visible compared to the sponge in the cIEF cartridge.



cIEF Cartridge

icIEF 400 Cartridge

At the end of the batch

1. Place the cartridge on a flat surface with the electrolyte tanks facing up.
2. Remove the stoppers on both electrolyte tanks.
3. Using a pipette or low vacuum, aspirate the solution from each tank.
4. Fill each tank with 2.0 mL of DI water, then aspirate it out. Repeat this 3 times.
5. Aspirate all the remaining liquid and make sure the tanks are dry.

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners. PLO-0019, Rev A