

MauriceFlex 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 MauriceFlex cIEF Fractionation Application Kit.

Note: Be sure to prepare batch reagents with crimp top glass vials (PN 110-0018 for the Fluorescence Calibration Standard and PN 110-0019 for all other reagents).

Put Your Reagent and Samples in MauriceFlex

1. Open MauriceFlex's door by touching the metal plate at the top of the door.
2. Install the fractionation adapter in the instrument.
3. Place your prepped batch reagents in Row R and Column K.
 - **R1:** 0.5% Methyl Cellulose
 - **R2:** Fluorescence Calibration Standard
 - **R3:** Water
 - **R4:** Water
 - **R5:** Water
 - **R6:** Air (empty vial)
 - **K2:** Catholyte
 - **K5:** Water

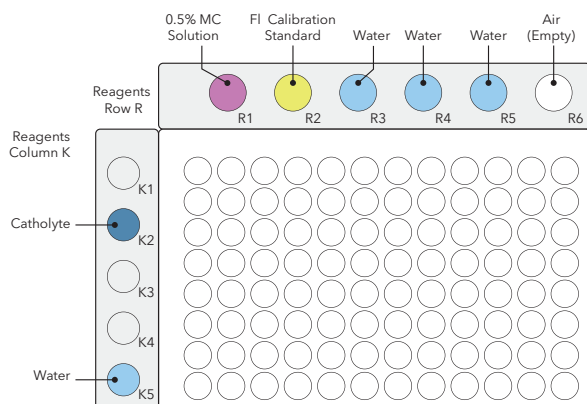


FIGURE 1. Sample and reagent vial placement.

4. Put your prepared sample in the 96-well plate and place the plate on the sample platform.

Install Your Cartridge

1. Take the cIEF Fractionation 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 a failed injection.

Note: Avoid touching the capillary while holding the cartridge.

2. Place the cartridge on a flat surface with its electrolyte tank facing up.
3. Remove the red stopper from the tank.
4. Add 2 mL of analyte to the tank.

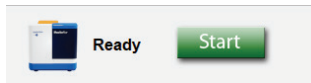
Note: Make sure you don't overfill the tank.



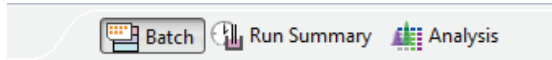
5. Seal the tank with the rubber stopper. If excess liquid comes out of the tank, make sure to wipe it with a lint-free laboratory wipe.
6. Open MauriceFlex's door.
7. Slide the cartridge into the cartridge slot. The lights around the cartridge will turn from **orange** to **blue**, and you'll feel a locking mechanism engage when the cartridge is correctly installed.

Create and Start Your Batch

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



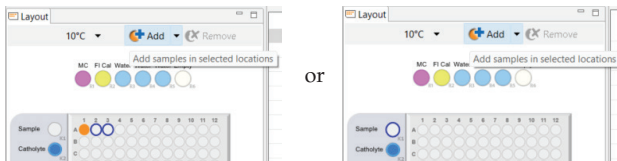
2. Click the **Batch** tab.



3. Set up your batch:

In the File menu, click **New Batch**. Select **MauriceFlex cIEF**.

Adding a sample: Highlight the sample location in the **Layout** window and click **Add**. A sample can be added to any well in the 96-well plate or in vial location K1.



Note: You can perform four injections per MauriceFlex cIEF batch.

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
NIST	10.0 min 500 Volts, 10.0 min ...	1 Exposure	20	7.05, 10.17		

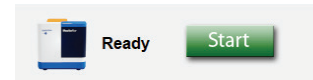
Adding sequential replicate injections: Highlight the injection in the **Injections** window and click **Replicate**.

Injection Name	Sample ID	Location	Method	Notes
1 Sample 01_01	Sample 01	A1	Method1	
2 Sample 01_02	Sample 01	A1	Method1	
3 Sample 02_03	Sample 02	A2	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**.

Injection Name	Sample ID	Location	Method	Notes
1 Sample 01_01	Sample 01	A1	Method1	
2 Sample 02_02	Sample 02	A2	Method1	
3 Sample 03_03	Sample 03	A3	Method1	
4 Sample 02_04	Sample 02	A2	Method1	

4. Click **Start**.



At the End of the Batch

1. Open MauriceFlex's door.
2. Remove your samples.
3. Clean-up and store your cartridge using the instructions in the Maurice User Guide or the MauriceFlex cIEF Fractionation Application Kit.

Note: A cIEF Fractionation cartridge supports a maximum of 15 injections.

biotechne®

Contact Us



Trademarks and registered trademarks are the property of their respective owners.