

MauriceFlex Fractionation Quick Reference Guide

Prepare Your Reagents and Sample

1. Prepare your sample 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 Sample 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
- **K3:** Mobilizer Solution (5 mM ammonium acetate)
- **K4:** Mobilizer Solution (5 mM ammonium acetate)
- **K5:** Water

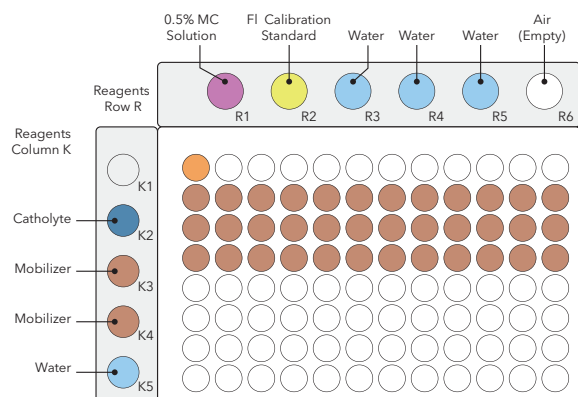


FIGURE 1. Sample and reagent vial placement.

4. Put your prepared Mobilizer Solution and 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.

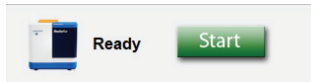
Note: Make sure you don't overfill the tank.



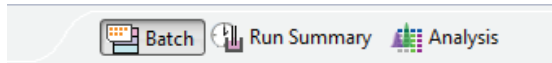
5. Seal the tank with the rubber stopper. If excess liquid comes out of the tank, make sure to wipe it with a lint-free laboratory wipe.
6. Open MauriceFlex's door.
7. Slide the cartridge into the cartridge slot. The lights around the cartridge will turn from **orange** to **blue**, and you'll feel a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



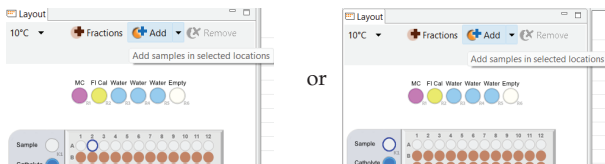
2. Click the **Batch** tab.



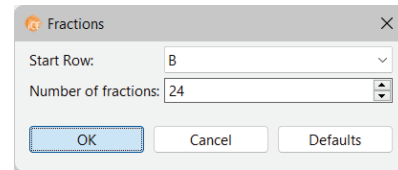
3. Set up your batch:

In the File menu, click **New Batch**. Select **MauriceFlex Fractionation**. The sample is located in well A1 and fractions are located in rows B–D in the default batch.

Changing the sample location: Highlight well A1 in the **Layout** window and click **Remove**. Then, highlight the sample location in row A or the vial location in K1 and click **Add**.



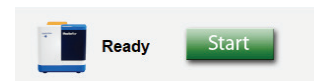
Adding fraction wells: Click **Fractions** in the **Layout** window. Choose the **Start Row** and the **Number of Fractions**, and click **OK**. Fractions must start in well 1 of a row.



Note: A minimum of 24 fractions and a maximum of 84 fractions can be collected in one batch.

Note: Fractions are collected in a serpentine pattern that cannot be adjusted.

4. Click **Start**.



At the End of the Batch

1. Open MauriceFlex's door.
2. Remove your samples.

Note: We recommend that you seal the plate to prevent evaporation if the fractions will not be used immediately.

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.