

Profinia™ Inline Filter Replacement Kit

Catalog #

620-0404

Quantity

12 x inline filters, 1 x 30 ml Luer-Lok syringe

Description

The inline filter is a 10 µm membrane filter that protects the system from particles in the inlet fluid stream.

Storage

Store at ambient temperature or 4°C. Keep away from direct heat sources.

Instructions for Use

The inline filter is located under the top cover of the instrument above the buffer bottles, near the left side of the instrument. The inline filter should be replaced every month, or after 20 purification runs.

Prepare filter for installation:

1. Fill the Luer-Lok syringe with 30 ml of 20% ethanol.
2. Connect 30 ml Luer-Lok syringe to the inline filter.
3. Push 30 ml of 20% ethanol through the inline filter in approximately 30 seconds.
4. Remove inline filter from syringe and follow installation instructions.



Connecting syringe to inline filter



Filter connected to syringe

Install the filter:

1. Open the top cover and locate the filter.
2. Twist both luer fittings counterclockwise to unlock them, then remove the filter.
3. Install the new filter, and tighten the fittings.



Location of the inline filter under the instrument top cover

Luer-Lok is a trademark of Becton, Dickinson and Company