

Profinia™ Sipper Tube Kit

Catalog #

620-0405

Quantity

10 sipper tubes and 10 silicone couplers

Description

Profinia sipper tubes are made of polypropylene tubing. The tubing is chemically resistant to most commonly used aqueous buffers.

Storage

Store at ambient temperature. Keep away from direct heat sources.

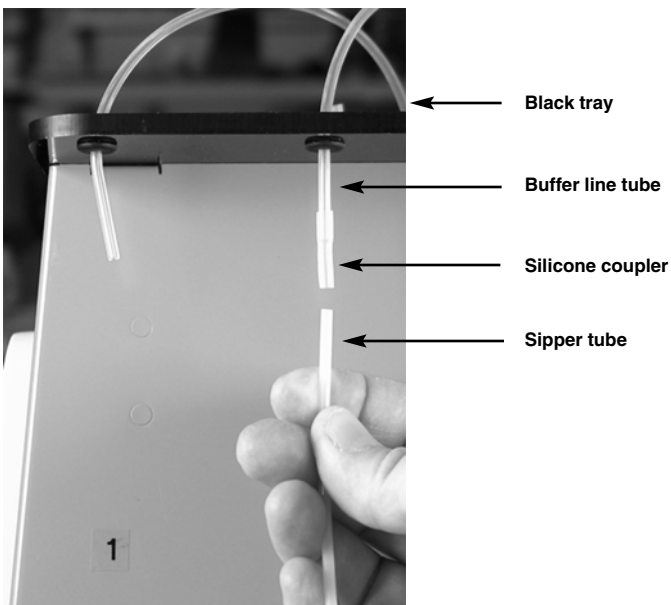
Instructions for Use

To replace silicone coupler:

1. Use one hand to hold the buffer line tube, and the other to disconnect the silicone coupler by pulling it off.
2. Install the new silicone coupler by pushing one end into the buffer line tube.
3. Make sure the buffer line tube length from the black tray is approximately 1 inch in length.

To replace sipper tube:

1. Use one hand to push up on the silicone coupler, and the other to disconnect the sipper tube by pulling it off.
2. Install the new sipper tube by pushing one end into the silicone coupler.



Sipper tube and coupler replacement