

Profinia™ Instrument Accessory Kit

Catalog No.

620-0410

Quantity

Accessory kit comes with all the necessary components to work with the Profinia instrument. This kit includes a cleaning tray, 12 inline filters, 8 buffer lids, 2 x 15 ml sample lids, 2 x 50 ml sample lids, 4 x 125 ml buffer bottles, 4 x 250 ml buffer bottles, and a waste/diluent bottle set.

Description

The Profinia bottles are made of polypropylene and are chemically compatible for use with commonly used aqueous buffers. The kit includes 8 bottles for installing custom buffers and for cleaning protocols. Buffer and sample lids are provided to prevent dust and other contaminants from entering the system.

Storage

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

Instructions for Use

Fill bottles with aqueous solutions of choice. Place green buffer lids on each bottle. Insert each buffer bottle into the corresponding buffer sipper tube located in the buffer placement tray of the Profinia instrument. Clean thoroughly with appropriate cleaning reagents before reusing.



Profinia Instrument Accessory Kit