

Profinia™ Bottle Starter Pack

Catalog #

620-0231

Quantity

4 x 125 ml bottles, 4 x 250 ml bottles, and 8 buffer lids

Description

The Profinia bottles are made of polypropylene and are chemically compatible for use with commonly used aqueous buffers. The starter pack includes 8 bottles for installing custom buffers and for cleaning protocols. Buffer 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 tray area of the Profinia instrument. Clean thoroughly with appropriate cleaning reagents before reusing.



Profinia bottle starter pack