

Profinia™ Buffer Lids

Catalog #

620-0407

Quantity

8 buffer lids

Description

Buffer lids keep dust and other debris contaminants out of the Profinia buffer bottles used in the Profinia instrument.

Storage

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

Instructions for Use

Place green buffer lids on buffer bottle. Insert each buffer bottle into buffer bottle slots in the buffer compartment of the Profinia instrument. Clean thoroughly with appropriate cleaning reagents before reusing.



Profinia buffer lid placed on buffer bottle