

Profinia™ 50 ml Sample Lids

Catalog #

620-0408

Quantity

2 x 50 ml sample lids

Description

Sample lids keep dust and other debris contaminants out of the sample tubes.

Storage

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

Instructions for Use

Add sample to the 50 ml conical tubes. Place sample lids on 50 ml conical tubes. Place into the Profinia instrument by lifting the sipper tube away from the sample loading area and inserting it into the conical tube. Clean thoroughly with appropriate cleaning reagents before reusing.



Profinia 50 ml sample lids



Profinia 50 ml sample lid on conical tube