

Profinia™ Cleaning Tray

Catalog #

620-0400

Quantity

1 cleaning tray

Description

The Profinia cleaning tray is a chemically compatible plastic tray.

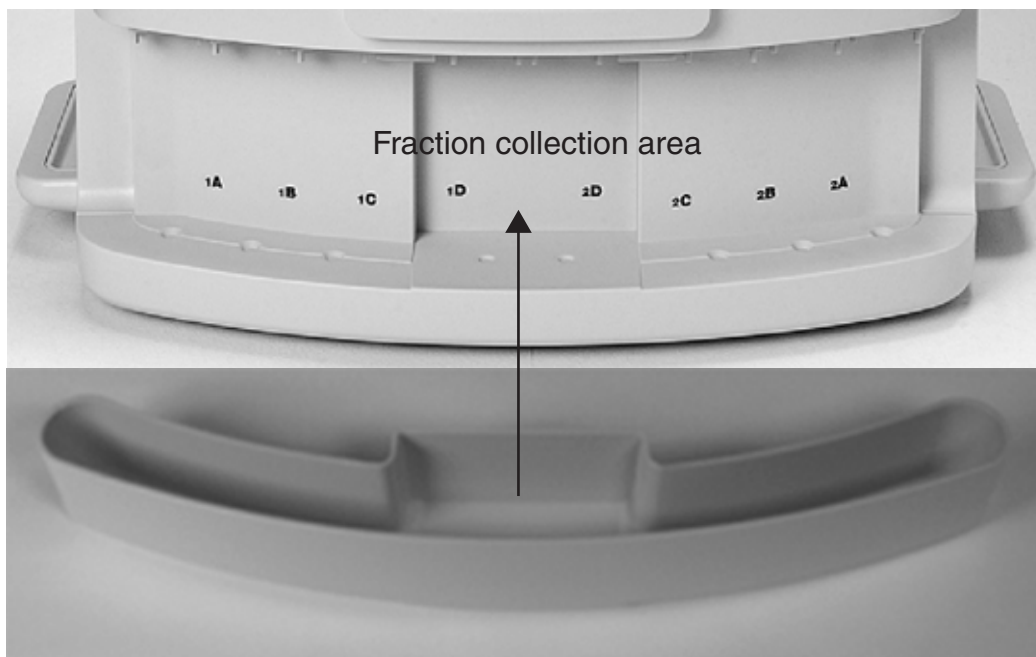
Storage

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

Instructions for Use

Insert the Profinia cleaning tray into the fraction collection area after removing the fraction tubes and before you start running the cleaning protocol.

Tray should be emptied and cleaned before storage to avoid salt buildup.



Profinia cleaning tray placement