

Profinia™ Cooling Accessory

Catalog #

620-0401

Quantity

2 cooling units and 4 x 15 ml conical tube holder ring inserts.

Description

The Profinia cooling accessory allows users to work with temperature-sensitive proteins for sample loading and collection on the Profinia system at room temperature environments. Each cooling unit has two slots that can hold 15 or 50 ml conical tubes.

Storage

Store at ambient temperature or 4°C. Keep away from direct heat sources.

Instructions for Use

To prepare for use with temperature-sensitive samples, fill cooling unit with water to the water line marked inside the cooling unit and freeze. Insert conical tubes into holding slots. If using 15 ml conical tubes, please use the provided black holding rings to keep tubes upright in the unit. Insert the cooling units into the sample loading area and the fraction collection compartment, located in the middle of the fraction collection area, marked as 1D and 2D.

The cooling units can maintain precooled samples and fractions at 2–8°C for 4–6 hours at 20°C ambient temperature, with the instrument powered up and operating.



Profinia cooling accessory