

Profinia™ 2 ml and 10 ml Desalting Sample Loop

Catalog #

620-0402, 620-0403

Quantity

1 precut Tefzel sample loop, 1 female luer to 1/4–28 female fitting, 1 male luer to 1/4–28 female fitting, 3-way stopcock, and 10 ml syringe.

Description

The Profinia 2 ml and 10 ml desalting sample loops come preassembled with luer fittings for a quick and easy connection to the Profinia instrument. Desalting loops are required for use with desalting-only and buffer exchange-only applications on the Profinia instrument. A desalting sample loop is not required for combined affinity and desalting methods.

Storage

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

Instructions for Use

Select the desalting-only method on the Select Method Type and Options screen. Follow the touch screen instructions. The desalting sample loop connects to the instrument in the cartridge 1 position and the desalting cartridge in the cartridge 2 position.

To connect and fill the desalting loop:

1. Insert the male luer fitting from cartridge holder 1 into the top female luer fitting of the 3-way stopcock. Ensure that the second female luer fitting of the 3-way stopcock is facing out so the syringe may be easily inserted.
2. Connect the male luer of the end of the desalting loop to the female luer fitting of cartridge holder 1.
3. Install the desalting cartridge in the cartridge 2 position. Touch the Next button to proceed to the Start Run screen.
4. To apply sample, fill the syringe and insert it into the open female luer of the 3-way stopcock. Turn the 3-way stopcock lever up (toward the top inlet fitting), and inject the sample into the loop.
5. Return the 3-way stopcock lever to the middle position, facing toward the syringe. The desalting loop is now filled with sample and ready for the desalting method to begin.



Profinia desalting sample loop