

ProFlow™ Sort Grade Water

Catalog #	Description
145-1083	ProFlow Sort Grade Water , 5 x 4 L

For research purposes only.

Description

Designed to work optimally with the S3™ or S3e™ Cell Sorter, ProFlow Sort Grade Water gives you confidence in your instrument fluidics.

ProFlow Sort Grade Water is prepared by reverse osmosis, passed through fine carbon, deionized through two resin beds, and serially filtered twice through 0.1 µm positively charged membranes to yield a sterile, endotoxin-free product. Each container is gamma irradiated to ensure sterility prior to filling.

Contents

5 x 4 L containers.

Storage

Store at 15–30°C.

Important: Do not freeze.

Quality Control

Each lot has been created with a careful selection of raw materials, current good manufacturing practice (cGMP) procedures, and diligent monitoring of each step to produce high-quality, sterile sort grade water for all cell sorting purposes.

Product-Specific Testing

ProFlow Sort Grade Water is tested in a cGMP environment to verify that it meets product specifications for the following parameters:

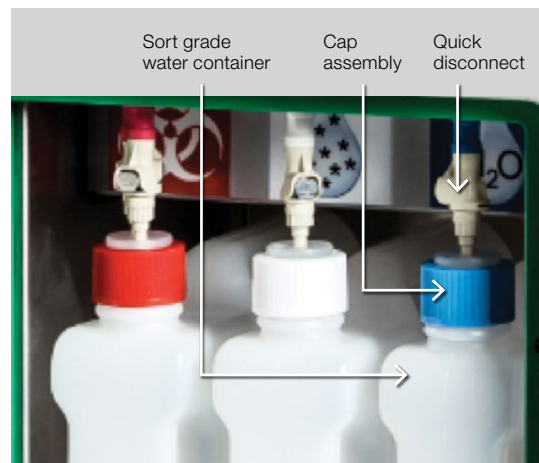
- Endotoxin level — less than 0.03 EU/ml
- Sterility — no bacterial or fungal growth
- Appearance — clear liquid

Directions for Use after S3 or S3e Cell Sorter Shutdown

Important: Use sterile laboratory techniques.

1. Detach the quick disconnect from the cap assembly (see figure).
2. Slightly lift and pull out the empty sort grade water container.
3. Uncap and remove the seal from a new 4 L ProFlow Sort Grade Water container.

4. Uncap the cap assembly from the empty sort grade water container.
5. Carefully place the cap assembly into the new 4 L sort grade water container and tighten the cap.
6. Place the bottle into position and attach the quick disconnect to the cap assembly. An audible click should be heard.



S3e Cell Sorter fluidics with sort grade water container components.

Note: Refer to the S3e Cell Sorter manual or quick guide at bio-rad.com/waterinfo to change the sort grade water container when the S3 or S3e Cell Sorter is turned on.

Visit bio-rad.com/web/WaterInfo for more information.