

# ProFlow™ Sort Grade 8x Sheath Fluid

Catalog #	Description
145-1082	<b>ProFlow Sort Grade 8x Sheath Fluid, 20 L (5 x 4 L)</b>

**For research purposes only.**

## Intended Use

ProFlow Sort Grade 8x Sheath Fluid is a sterile, endotoxin-free, and preservative-free phosphate buffered saline solution for use with Bio-Rad's S3™ or S3e™ Cell Sorter.

## Contents

5 x 4 L containers.

## Composition

The S3 and S3e Cell Sorters automatically dilute the 8x Sheath Fluid with sterile water to create 32 L of 1x sheath fluid from each container.

Disodium EDTA	1.0 mM
Potassium phosphate, monobasic	1.9 mM
Potassium chloride	3.8 mM
Sodium phosphate, dibasic	16.6 mM
Sodium chloride	139.0 mM
Preservative	None

## Storage

Store at 15–30°C.

**Important:** Do not freeze.

## Warning and Precautions

As with all chemical solutions, avoid eye, skin, and clothing contact with the concentrated or diluted reagent. No special precautions are recommended. See Material Safety Data Sheet for more details.

The 8x concentrate is clear, colorless, and particulate free. If there is evidence of turbidity, sediment, discoloration, or bacterial/fungal contamination, discard the reagent.

## Directions for Use after S3 or S3e Cell Sorter Shutdown

**Important:** Use sterile laboratory techniques.

1. Detach the quick disconnect from the sheath fluid cap assembly (Figure 1) by pushing down the metal button and lifting up the quick disconnect until it engages the magnetic holder. This will hold the quick disconnect away from the exchange area.
2. Slightly lift and pull out the empty sheath fluid container.
3. Uncap and remove the seal from a new 4 L sheath fluid container.
4. Uncap the sheath fluid cap assembly from the empty sheath fluid container.
5. Carefully place the sheath fluid cap assembly into the new 4 L sheath fluid container and tighten the cap.
6. Place the container into position and attach the quick disconnect to the sheath fluid cap assembly. An audible click should be heard.

**Caution:** This reagent is a preservative-free sheath fluid so bacterial contamination may occur.

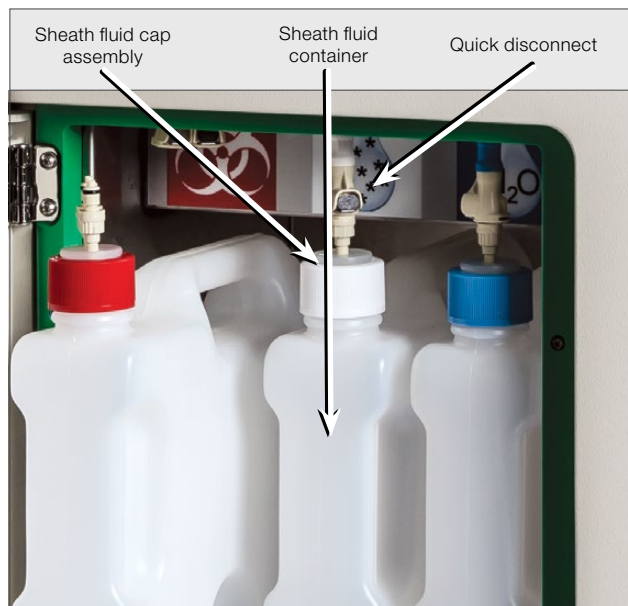


Fig. 1. S3e Cell Sorter fluidics with sheath fluid components noted.

## Directions for Performing a Hot Swap

Perform a hot swap to change the sheath fluid container during a run.

1. Select **Setup and Maintenance > Swap Fluidics** in the ProSort™ Software window. A countdown timer will automatically start (Figure 2). This is the amount of time left to swap the container before the system automatically shuts down.

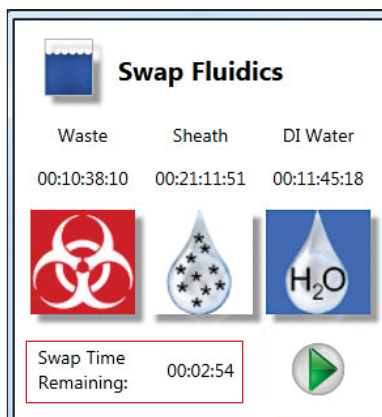


Fig. 2. Countdown timer highlighted in the Swap Fluidics window.

2. Open the fluidics door and detach the quick disconnect from the sheath fluid cap assembly by pushing down the metal button.
3. Lift the quick disconnect until it engages the magnetic holder. This will hold the quick disconnect away from the exchange area.
4. Replace the empty container with a new container.
5. Attach the quick disconnect to the sheath fluid cap assembly. An audible click should be heard.
6. Click **Play** in the Swap Fluidics window to finish the swapping procedure (Figure 3).

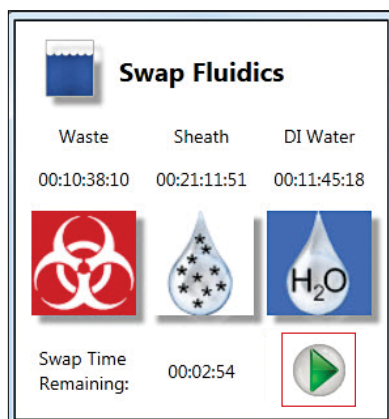


Fig. 3. Play button highlighted in the Swap Fluidics window.

Visit [bio-rad.com/web/Sheath](http://bio-rad.com/web/Sheath) for more information.