

Cytometer Cleaner

Catalog #	Description
12004272	Cytometer Cleaner , 1 L bottle

Accessories

12004395	ZE5™ Cleaner Bottle with Tubing , blue cap, includes 450 ml bottle, cap/tubing assembly with filter and quick connector
12004444	ZE5 Cleaner/Additive Bottle , includes 450 ml bottle

For research purposes only.

Intended Use

The cytometer cleaner contains a system-cleaning reagent. During the shutdown process, cleaner is delivered to portions of the fluidics system that are in contact with sample, such as the flow cell and sample line. This minimizes sample buildup over time and preserves the integrity of the system.

Storage Conditions

Store the cytometer cleaner at room temperature.

Directions for Use

Important: Do not dilute.

Important: The volume must not exceed the Max Fill line on the bottle label, otherwise the liquid will spill into the ZE5 Cell Analyzer fluidics compartment.

1. Fill the cytometer cleaner bottle (catalog #12004395) supplied with the ZE5 Cell Analyzer to the Max Fill line indicated on the bottle label.
2. Close the bottle with the supplied blue cap, which has attached tubing, filter, and quick connector. Make sure the cap is tightened.
3. Open the door to the fluidics compartment and place the cleaner bottle in the slot on the right below the bulk sheath and waste containers.
4. Attach the quick connector to the system cleaner connection. An audible click indicates that the connection is tight.

The weight of each bottle is continuously monitored so that the system can calculate the approximate remaining run time. Inspect the level of the cleaner on a weekly basis to ensure that the remaining amount is sufficient for system usage.

To refill the cytometer cleaner bottle, push the quick connector tab to release the line and remove the bottle.

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

