

## ZE5™ Additive

Catalog #	Description
12004271	<b>ZE5 Additive</b> , 4 x 300 ml bottles

For research purposes only.

### Intended Use

ZE5 Additive concentrate is drip-fed into the internal sheath (deionized [DI] water) reservoir on a continuous basis while the ZE5 Cell Analyzer is running. It contains balanced salts, an antimicrobial/antifungal agent, and a surfactant. The sheath additive manages bubbles that occur during normal operation of the system and helps prevent microbial buildup. Microbes in the fluidic lines have the potential to increase background noise in the data, especially in the scatter channels, and can clog the lines or flow cell if the buildup becomes significant.

### Storage Conditions

Store at 20–30°C.

### Directions for Use

**Important:** Do not dilute.

1. Remove the flat white cap and replace it with the white cap with attached tubing, filter, and quick connector (supplied with the ZE5 Cell Analyzer). Make sure the cap is tightened.
2. Open the door to the fluidics compartment and place the additive bottle in the slot on the left below the bulk sheath (DI water) and waste containers.
3. Attach the quick connector to the additive 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 additive on a weekly basis to ensure that the remaining amount is sufficient for system usage.

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

**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.

Visit [bio-rad.com/web/ZE5Additive](http://bio-rad.com/web/ZE5Additive) for more information.

