

## TECHNICAL DATA

**PRODUCT:** Amine Reactive Comp-Bead 2 Population Kit,  
2 Populations (Negative and High Binding), 7.0-7.9  $\mu\text{m}$ .

**DESCRIPTION:** Amine Reactive Comp-Bead 2 Population Kit is suitable for labeling with LIVE/DEAD® stains or other amine-reactive dyes to generate compensation standards for flow cytometric analysis.

**CATALOG NO.:** NBP3-00496

**LOT NO.:** AQ01

**SIZE:** 2 X 5 mL

**CONCENTRATION:**  $\sim 1 \times 10^7$  particles/mL

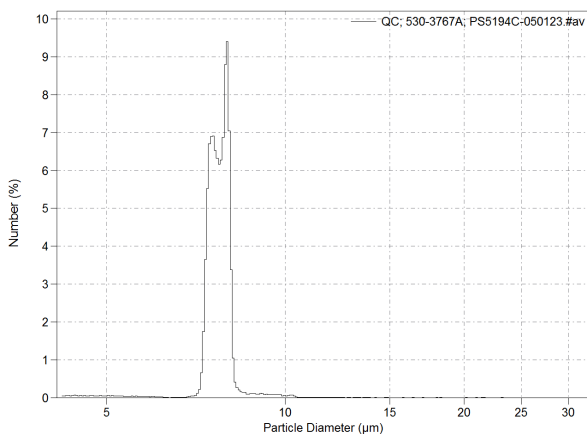
**STORAGE BUFFER:** 0.016M PBS, pH 7.4 with 0.02% Sodium Azide

**STORAGE TEMP.:** 2° to 8° C

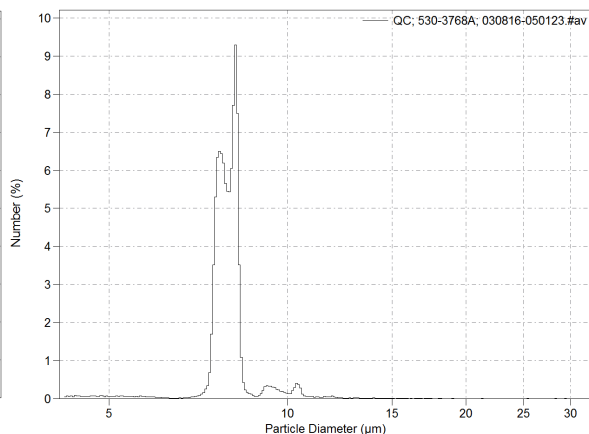
**CAUTION:** Do not freeze. Protect from light.

**NOTE:** Before use resuspend by vortexing. To achieve optimum particle suspension, sonicate the reagent after vortex mixing. Stable for 12 months from date of purchase provided the handling is in accordance with the manufacturer's recommendations.

### COULTER M3 ANALYSIS:



Negative Mean Size: 7.71  $\mu\text{m}$



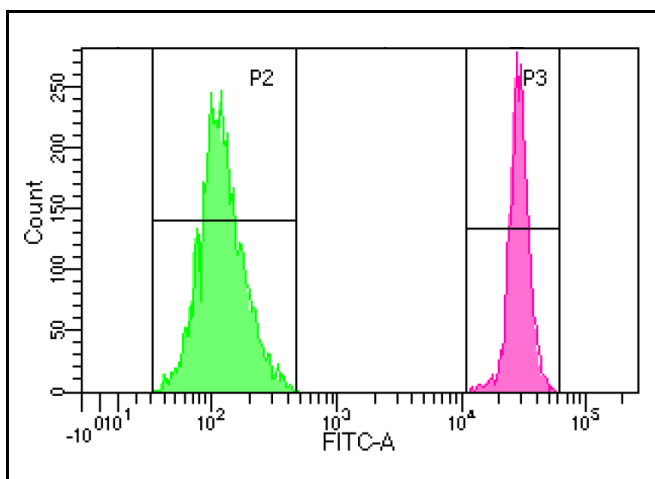
High Mean Size: 7.96  $\mu\text{m}$

**FOR RESEARCH APPLICATIONS ONLY. NOT FOR DIAGNOSTIC USE.**

### PROCEDURE FOR USE:

1. Remove Amine Reactive Comp-Beads from refrigerator and allow them to come to room temperature and briefly vortex.
2. Add 1 drop (~50 ul ) of the High Binding beads to a 1.5mL microcentrifuge tube
3. Wash the beads by adding 0.5mL of PBS that is free of surfactant and blocker to the microcentrifuge tube centrifuge at 300 x G for 5 minutes. Decant and repeat. Decant and resuspend in 50µL PBS
4. Prepare the amine reactive dye according to the manufacturer's instructions
5. Add 1-4µL of the amine reactive dye to the bead suspension and briefly vortex.  
Protect the tube from light and allow incubation for 30 minutes
6. Add 1ml of PBS to the same and briefly vortex. Centrifuge at 300 x G for 5 minutes, decant and repeat.  
Resuspend the beads in PBS containing 0.05% BSA with brief vortex
7. Add a drop (~50 ul ) of the Negative Binding beads to the labeled High Binding beads
8. Analyze on the flow cytometer using a live gate around the singlet population in the FSC/SSC dot plot
9. Create a fluorescence histogram for the appropriate detectors and perform compensation to achieve the desired results.

### HISTOGRAM:



Amine Reactive Comp-Bead 2 Population Kit stained with  
Biologend Zombie Green™

### NOTE:

1. Dye and bead concentrations may be further optimized for best results. Centrifugation force and time may need to be increased if needed.
2. Protect the beads from light after exposure to the dye. Uses immediately after staining.