

# Product Information

## ExoBrite™ EV Stain Enhancer (100X)

**Catalog Number:** 28002-T, 28002

### Unit Size

28002-T: 1 x 0.5 mL

28002: 5 x 0.5 mL

### Storage and Handling

Store at 4°C. Product is stable for at least 12 months from date of receipt when stored as recommended.

### Product Description

The ExoBrite™ EV Stain Enhancer is an additive that can be added to some EV stain reactions to improve the staining specificity for applications like flow cytometry. The ExoBrite™ Stain Enhancer works by reducing the aggregation of certain EV stains, which allows the conjugate to stain the EVs more efficiently, resulting in a better signal-to-noise ratio and fewer false positives. Enhancer has been shown to be beneficial for staining EVs with WGA, other lectins, and Annexin V, but it is not recommended for lipophilic EV stains like ExoBrite™ True EV Membrane Stains or PKH dyes. Enhancer can be used to decrease aggregation of Cholera Toxin B (CTB), but generally is not required due to the intrinsically low aggregation of CTB conjugates. Enhancer is generally not required for use with ExoBrite™ antibody conjugates because they are already formulated to reduce aggregation, however, it may provide benefits for certain antibodies that do show aggregates in flow.

### Experimental Protocol

The following is a general protocol for incorporating the ExoBrite™ EV Stain Enhancer into your EV staining protocol.

1. Isolate or purify EVs or exosomes using the procedure of your choice.
2. Aliquot 50 uL of EVs into FACS tubes or microcentrifuge tubes. Also, prepare a control tube of 50 uL of PBS or the buffer the EVs are stored in.
3. Prepare your staining solution of choice. You will want 450 uL per EV sample.
4. Add ExoBrite™ EV Stain Enhancer to the stain solution prepared in step 3 by diluting the 100X Enhancer stock 1:100. For example, add 4.5 uL of 100X Enhancer stock to 450 uL of stain solution.

**Note:** You may want to empirically determine the optimal concentration of Enhancer for your particular stain or dye. We recommend trying a range between 0.5X-2X.

5. Mix the stain + Enhancer mixture from step 4 well.
6. Add 450 uL of the stain + Enhancer mixture to each tube of EVs and to the buffer-only control tube.
7. Incubate as desired and detect EVs as usual.

### Related Products

| Cat. No.                | Product                                       |
|-------------------------|---|
| 30123-30126             | ExoBrite™ WGA EV Staining Kits                |
| 30119-30122             | ExoBrite™ Annexin EV Staining Kits            |
| 30111-30114             | ExoBrite™ CTB EV Staining Kits                |
| 30127                   | ExoBrite™ EV Surface Stain Sampler Kit, Green |
| 30129...<br>30137       | ExoBrite™ True EV Membrane Stains             |
| 30115-30118             | ExoBrite™ STORM CTB EV Staining Kits          |
| 28001                   | ExoBrite™ EV Total RNA Isolation Kit          |
| 28000                   | ExoBrite™ Streptavidin Magnetic Beads         |
| P003-410...<br>P003-RPE | ExoBrite™ CD9 Flow Antibody                   |
| P018-410...<br>P018-650 | ExoBrite™ CD9 (Mouse) Flow Antibody           |
| P004-410...<br>P004-RPE | ExoBrite™ CD63 Flow Antibody                  |
| P022-410...<br>P022-650 | ExoBrite™ CD63 (Mouse) Flow Antibody          |
| P005-410...<br>P005-RPE | ExoBrite™ CD81 Flow Antibody                  |
| P019-410...<br>P019-650 | ExoBrite™ CD81 (Mouse/Rat) Flow Antibody      |
| P008-410...<br>P008-RPE | ExoBrite™ IgG1 Isotype Control Flow Antibody  |
| P003-680...<br>P003-770 | ExoBrite™ CD9 Western Antibody                |
| P004-680...<br>P004-770 | ExoBrite™ CD63 Western Antibody               |
| P006-680...<br>P006-770 | ExoBrite™ CD81 Western Antibody               |
| P007-770                | ExoBrite™ Calnexin Western Antibody           |

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.