

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... 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... P003-RPE	ExoBrite™ CD9 Flow Antibody
P018-410... P018-650	ExoBrite™ CD9 (Mouse) Flow Antibody
P004-410... P004-RPE	ExoBrite™ CD63 Flow Antibody
P022-410... P022-650	ExoBrite™ CD63 (Mouse) Flow Antibody
P005-410... P005-RPE	ExoBrite™ CD81 Flow Antibody
P019-410... P019-650	ExoBrite™ CD81 (Mouse/Rat) Flow Antibody
P008-410... P008-RPE	ExoBrite™ IgG1 Isotype Control Flow Antibody
P003-680... P003-770	ExoBrite™ CD9 Western Antibody
P004-680... P004-770	ExoBrite™ CD63 Western Antibody
P006-680... P006-770	ExoBrite™ CD81 Western Antibody
P007-770	ExoBrite™ Calnexin Western Antibody

Please visit our website at 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.