## **ExoBrite™ EV Stain Enhancer (100X)**

A unique additive that can be added to extracellular vesicle (EV) stain reactions to improve the staining specificity for applications like flow cytometry.



#### Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

### **Product Description**

The ExoBrite™ EV Stain Enhancer is a unique additive that can be added to extracellular vesicle (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.

#### **Features**

- Improves signal-to-noise by reducing or eliminating aggregates of certain EV stains
- Validated with several different lectins and Annexin V
- Does not interfere with antibody staining of EVs
- Easy to use, just add directly to the staining reaction

#### Get precise, reliable results with a range of popular EV probes

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<sup>TM</sup> 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<sup>TM</sup> antibody conjugates because they are already formulated to reduce aggregation, however, it may provide benefits for certain antibodies that do show aggregates in flow.

# Learn about Biotium's optimized ExoBrite™ stains and antibodies for EV research

Biotium's ExoBrite<sup>TM</sup> True EV Membrane Stains are genuine lipophilic membrane dyes are designed for superior pan-EV labeling over other membrane dyes including PKH, DiO, DiI, and DiD. Biotium also offers

ExoBrite<sup>TM</sup> Antibody Conjugates are optimized for detection of CD9, CD63, and CD81 EV markers by flow cytometry and western blotting. For super-resolution imaging by STORM, learn about our <a href="ExoBrite\*IM"><u>ExoBrite\*IM</u></a>
STORM CTB EV Staining Kits available in four CF® Dyes validated for STORM.

## ExoBrite™ EV Stains Comparison Guide

ExoBrite™ EV Surface Stain	Pros	Cons
ExoBrite™ True EV Membrane Stains	Near-complete staining of EVs in a sample     Broad compatibility with different EV sources     Validated for flow and fNTA	Can't be used to stain bead-bound EVs     May have more aggregation than CTB & Annexin
ExoBrite™ Annexin EV Staining Kits	Broad compatibility with different EV sources     Validated for flow and fNTA     Low background aggregates	May not stain every EV in a sample     Doesn't work well on bead-bound EVs
ExoBrite™ WGA EV Staining Kits	Broad compatibility with different EV sources     Can be used with bead-bound EVs	May not stain every EV in a sample     Doesn't work well for fNTA
ExoBrite™ CTB EV Staining Kits	Validated for flow and fNTA     Extremely low background	May not stain every EV in a sample     Does not stain EVs from every source
ExoBrite™ Antibodies	Highly specific for human tetraspanins CD9, CD63, or CD81     Validated for EV flow     Broad compatibility for different EV sources     Can be used with bead-bound EVs     Can be used for WB	Depends on the expression level of the target protein on the EVs