



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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

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™ 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.

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

Biotium's [ExoBrite™ 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™ EV Surface Stains which are fluorescent conjugates of probes for labeling EV membrane surface targets. ExoBrite™ EV Surface Stains are available as [cholera toxin subunit B \(CTB\)](#), [wheat germ agglutinin \(WGA\)](#), or [Annexin V](#) conjugates. A convenient [ExoBrite™ EV Surface Stain Sampler Kit](#) is also available, the kit includes each ExoBrite™ EV Surface Stains (CTB, WGA, and Annexin V) for assessing which stain offers the best coverage for the EV samples of interest.

[ExoBrite™ 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 [ExoBrite™ 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	<ul style="list-style-type: none">• Near-complete staining of EVs in a sample• Broad compatibility with different EV sources• Validated for flow and fNTA	<ul style="list-style-type: none">• Can't be used to stain bead-bound EVs• May have more aggregation than CTB & Annexin
ExoBrite™ Annexin EV Staining Kits	<ul style="list-style-type: none">• Broad compatibility with different EV sources• Validated for flow and fNTA• Low background aggregates	<ul style="list-style-type: none">• May not stain every EV in a sample• Doesn't work well on bead-bound EVs
ExoBrite™ WGA EV Staining Kits	<ul style="list-style-type: none">• Broad compatibility with different EV sources• Can be used with bead-bound EVs	<ul style="list-style-type: none">• May not stain every EV in a sample• Doesn't work well for fNTA
ExoBrite™ CTB EV Staining Kits	<ul style="list-style-type: none">• Validated for flow and fNTA• Extremely low background• Can be used with bead-bound EVs	<ul style="list-style-type: none">• May not stain every EV in a sample• Does not stain EVs from every source
ExoBrite™ Antibodies	<ul style="list-style-type: none">• Highly specific for human tetraspanins CD9, CD63, CD81, and other EV markers• Validated for EV flow• Broad compatibility for different EV sources• Can be used with bead-bound EVs• Can be used for WB	<ul style="list-style-type: none">• Depends on the expression level of the target protein on the EVs
ExoBrite™ EV Stain Enhancer	<ul style="list-style-type: none">• 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	<ul style="list-style-type: none">• Not recommended for use with lipophilic EV stains

This datasheet was generated on June 10, 2026 at 03:09:32 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/exobrite-ev-stain-enhancer-100x/>