



Glowing products for science

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

ExoBrite™ CTB EV Staining Kits

Fluorescent cholera toxin subunit B (CTB) conjugates that are optimized for bright and clean staining of extracellular vesicles for flow cytometry.



Product attributes

Dye	ExoBrite™ 410/450, ExoBrite™ 490/515, ExoBrite™ 560/585, ExoBrite™ 640/660
Colors	Blue, Green, Orange-red, Far-red

Product Description

ExoBrite™ CTB EV Staining Kits were designed to overcome some of the challenges of EV detection, particularly in flow cytometry. ExoBrite™ CTB EV Stains bind to molecules in the EV membrane for bright, specific staining, with little to no background.

Note: The name of this product has been revised from ExoBrite™ EV Membrane Staining Kits.

ExoBrite™ CTB EV Stains are optimally formulated fluorescent conjugates of cholera toxin subunit B (CTB), which binds to GM1 gangliosides that are commonly found on the surface of mammalian lipid rafts and EVs. The stains were designed to overcome some of the challenges of EV detection, particularly in flow cytometry. Some dyes used to stain EVs can form aggregates of a similar size as exosomes or EVs, thus confounding analysis. ExoBrite™ CTB EV Stains, however, were formulated to show little to no aggregation in flow cytometry, allowing EVs to be identified with bright and specific staining. Unlike hydrophobic membrane dyes, ExoBrite™ CTB EV Stains do not bind non-specifically to polystyrene beads, meaning that they can be used to stain bead-bound EVs.

EVs are often labeled with fluorescent antibodies targeting one or more of the tetraspanin proteins CD9, CD63, and CD81. ExoBrite™ CTB staining can be combined with antibody staining, for multi-parameter analysis.

Less background and better coverage over other EV stains

ExoBrite™ CTB EV Stains were designed to offer exceptional signal:noise and more complete coverage of purified and bead-bound EVs. In the figure below, lipophilic dye DiO and plasma membrane stain CellMask™ demonstrate an unacceptable amount of dye aggregation in gated EVs. Other EV stains such as ExoFlow-ONETM and ExoGlow™ have less coverage of EVs when compared to ExoBrite™ CTB EV Stains.

Notes:

1. ExoBrite™ CTB EV Stains have been found to label EVs derived from several tested cell lines (see Validated EV Sources below), but may not stain EVs from every source.
2. In our testing, we have found that ExoBrite™ 490/515 dye may bind to streptavidin coated surfaces or beads if free biotin binding sites are not blocked. We recommend performing a biotin blocking step after binding your biotinylated capture antibody to streptavidin beads or surfaces when using ExoBrite™ 490/515 conjugates. Alternatively, consider using a different ExoBrite™ dye for staining EVs captured on streptavidin beads or surfaces.

ExoBrite™ CTB EV Staining Kits

Product	Ex/Em	Detection channels	Size	Catalog Number
ExoBrite™ 410/450 CTB EV Staining Kit	416/452 nm	Pacific Blue™	100 Labelings	30111-T
500 Labelings	30111			
ExoBrite™ 490/515 CTB EV Staining Kit	490/516 nm	FITC	100 Labelings	30112-T
500 Labelings	30112			
ExoBrite™ 560/585 CTB EV Staining Kit	562/584 nm	PE, Cy®3	100 Labelings	30113-T
500 Labelings	30113			
ExoBrite™ 640/660 CTB EV Staining Kit	642/663 nm	APC	100 Labelings	30114-T
500 Labelings	30114			

Validated EV Sources for ExoBrite™ EV Surface Stains

EV Source	ExoBrite™ True EV Membrane Stains	ExoBrite™ CTB Stains	ExoBrite™ WGA Stains	ExoBrite™ Annexin Stains
A549 cells	Yes	Yes	Yes	Yes
CHO cells	Yes	No	Yes	Yes
hASC (human adipose stem cells)	ND	No ¹	ND	ND
HEK293 cells	Yes	Yes ¹	Yes	ND
HeLa cells	Yes	No	Yes	Yes
HUVEC (human umbilical vein endothelial cells)	ND	No ¹	ND	ND
J774 cells	Yes	Yes	Yes	Yes
Jurkat cells	Yes	Yes	Yes	Yes
MCF-7 cells	Yes	Yes	Yes	Yes
Plasma	ND	No	ND	Yes
Raji cells	ND	Yes	Yes	Yes
RAW 264.7 cells	Yes	ND	ND	ND
Serum	ND	No	ND	Yes
Skeletal myoblasts	ND	Yes ¹	ND	ND
THP-1 cells	Yes	ND	ND	ND
U2OS cells	Yes	No	Yes	Yes
U937 cells	Yes	No	Yes	Yes
NIH3T3 cells	Yes	ND	ND	ND
HepG2 cells	ND	ND	Yes	ND
Yeast (S. cerevisiae)	Yes	No	Yes	Yes

¹Customer-reported data
Value of “Yes” or “No” indicates coverage of EVs based on Biotium’s internal data or customer-reported data. Value of “ND” indicates no data.

Biotium also offers other validated ExoBrite™ reagents for flow cytometry, western blotting, or super-resolution imaging.

Learn about Biotium’s new [ExoBrite™ True EV Membrane Stains](#). These genuine lipophilic membrane dyes are designed for superior pan-EV labeling over other membrane dyes including PKH, DiO, Dil, and DiD. Biotium also offers [ExoBrite™ WGA EV Stains](#) (wheat germ agglutinin) and [ExoBrite™ Annexin EV Stains](#) optimized for bright and sensitive staining of EVs. The [ExoBrite™ EV Surface Stain Sampler Kit](#) contains each of Biotium’s ExoBrite™ EV Surface Stains (CTB, WGA, and Annexin V) for assessing which stain offers the best coverage for the EV samples of interest. Biotium also offers [ExoBrite™ Antibody Conjugates](#) for optimal 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.

CellMask is a trademark of Thermo Fisher Scientific; ExoFlow-ONE and ExoGlow are trademarks of System Biosciences.