

Product Information

ExoBrite™ True EV Membrane Stains

Product list

Cat. No.	Product	Unit Size
30129	ExoBrite™ 515/540 True EV Membrane Stain	500 µL
30129-T		100 µL
30130	ExoBrite™ 555/575 True EV Membrane Stain	500 µL
30130-T		100 µL

For more information see Table 1 on page 2.

Storage and Handling

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

Spectral Properties: See Table 1 on page 2.

Product Description

Extracellular vesicles (EVs) are lipid-bound vesicles that are released from cells. EVs display specific surface proteins and can carry nucleic acids and other cargo, allowing them to transfer biological information between cells in different parts of the body. Therefore, EVs are increasingly studied for their potential use in drug delivery and medical diagnostic applications. Biotium developed the ExoBrite™ product line for fluorescent labeling and detection of EVs by flow cytometry. Other potential applications include fluorescence nanoparticle tracking analysis (fNTA), fluorescence microscopy, and other fluorescence detection platforms.

ExoBrite™ True EV Membrane Stains are the best choice for general pan-EV staining. They are a significant improvement upon classic membrane dyes like PKH or DiO/DiI/DiD-type dyes for EV staining. They have higher (near-complete) coverage of EVs as determined by fNTA, much higher than competitor membrane dyes. They also have greatly improved solubility, reducing the amount of aggregation and allowing stained EVs to be readily differentiated from background particles, which is not usually possible with other membrane dyes. ExoBrite™ True EV Membrane Stains labeled all EV types tested to date. See Table 2 on page 2 for a list of validated EV sources.

EVs are often labeled with fluorescent antibodies targeting one or more of the tetraspanin proteins CD9, CD63, and CD81. ExoBrite™ True EV Membrane staining can be combined with antibody staining, for multi-parameter analysis (see Experimental Protocols). Biotium offers a selection of fluorescent ExoBrite™ Flow Antibodies against CD9, CD63, and CD81 that are optimized for detection of EVs by flow cytometry (see Related Products).

Biotium also offers other conjugates optimized for bright and sensitive staining of EVs. This includes ExoBrite™ Annexin EV Stains, ExoBrite™ CTB EV Stains (cholera toxin B conjugates), and ExoBrite™ WGA EV Stains (wheat germ agglutinin conjugates) (see Related Products).

For more suggestions and other protocols for working with EVs see our Tech Tips: [Isolation and Staining of Extracellular Vesicles](#) and [Fluorescent Detection of EVs by Flow Cytometry](#).

Considerations for Detecting EVs by Flow Cytometry

- EVs are extremely small vesicles (~30-150 nm in diameter), a size which is near or below the detection limit of some flow cytometers. We recommend determining the size detection limit of your instrument by running sizing beads (for example, ranging from 0.02-2 µm) in SSC before attempting to detect purified EVs. We also recommend running sizing beads before each EV detection experiment and using them to set the SSC threshold.
- Consider using a 405 nm laser for the SSC instead of a 488 nm laser for improved sensitivity for small particles.
- Use a low flow rate to keep the event rate and abort rate low. This will result in reduced noise.
- For best results, buffers used for suspending and staining EVs should be filtered through a 0.2 µm filter to remove particulates.

Considerations for Staining with ExoBrite™ True EV Membrane Stains

The following are general considerations for using ExoBrite™ True EV Membrane Stains to stain EVs. See Experimental Protocols for step-by-step instructions.

- ExoBrite™ True EV Membrane Stains have been validated in flow cytometry on the CytoFLEX LX from Beckman Coulter and in fNTA on the ZetaView® QUATT NTA System from Particle Metrix. Results on other instruments may vary based on the instrument's parameters.
- ExoBrite™ True EV Membrane Stains were validated using size exclusion column (SEC) enriched EVs. Staining results may vary depending on the EV isolation method used. We do not recommend using ExoBrite™ True EV Membrane Stains to stain bead-bound EVs. For bead-bound EVs we recommend using ExoBrite™ CTB EV Stains or ExoBrite™ WGA EV Stains (see Related Products).
- Individual EVs are too small to be imaged by conventional fluorescence or confocal microscopy, but clusters of EVs taken up by cells may be visualized. ExoBrite™ True EV Membrane Stains have not been validated for labeling EVs for cellular uptake. It may be necessary to remove free stain (by ultrafiltration, for example) before attempting to apply ExoBrite™ True EV Membrane Stain-labeled EVs to cells.
- ExoBrite™ True EV Membrane Stains have not been validated for super-resolution applications. For imaging EVs by STORM, we recommend our ExoBrite™ STORM CTB EV Staining Kits (see Related Products).
- ExoBrite™ True EV Membrane Stains have been found to label EVs isolated from cell culture supernatant from every cell line tested at Biotium (see Table 2), but staining may vary for EVs from other biological fluids or sources.
- While we have found that staining with 1X ExoBrite™ True EV Membrane Stains give a bright signal and low background under our typical staining conditions, the dye concentration may need optimization for different samples and detection systems.
- ExoBrite™ True EV Membrane Stains can be used for co-staining with fluorescently labeled primary antibodies. Co-staining can be performed concurrently or sequentially (see "Antibody co-staining of purified EVs" under Experimental Protocols).

Table 1. ExoBrite™ True EV Membrane Stains

Cat. No.	Number of labeling reactions	Product Name	Ex/Em	Laser Line(s) (nm)	Detection Channel	Compatible Applications
30129	500	ExoBrite™ 515/540 True EV Membrane Stain	515/542 nm	488, 532	FITC	Flow cytometry, fNTA
30129-T	100					
30130	500	ExoBrite™ 555/575 True EV Membrane Stain	556/576 nm	532,561	PE	Flow cytometry, fNTA
30130-T	100					

Table 2. Validated EV Sources for ExoBrite™ True EV Membrane Stains

Staining validated with EVs from the following cell lines
MCF-7, J774, U2OS, Jurkat, HeLa, CHO, U937, A549, THP-1, RAW 264.7

Experimental Protocols

Note: Before beginning, please read "Considerations for Staining with ExoBrite™ True EV Membrane Stains".

Staining of purified EVs

This protocol was developed for staining purified EVs with ExoBrite™ True EV Membrane Stains for detection by flow cytometry.

1. Isolate or purify EVs using the procedure of your choice.
2. Aliquot 50 uL of EVs into FACS tubes or microcentrifuge tubes.
3. In addition to the EVs stained with ExoBrite™ True EV Membrane Stain, it is helpful to include the following controls (the buffer should be an appropriate negative control for the EVs, such as a mock purification or the buffer used to suspend the EVs):
 - a. Buffer alone (no EVs, no stain)
 - b. Buffer plus ExoBrite™ True EV Membrane Stains
 - c. EVs alone (no stain)
4. Prepare 1X ExoBrite™ True EV Membrane staining solution by diluting the 500X dye stock 1:500 in filtered PBS or other buffer of choice.

Notes:

- a. The 1X ExoBrite™ True EV Membrane staining solution should be used the day of preparation.
 - b. The working concentration of ExoBrite™ True EV Membrane Stain can be optimized by the user.
5. Add 450 uL of 1X ExoBrite™ staining solution to each tube containing 50 uL sample. Remember to also add the staining solution to the "buffer plus ExoBrite™" control.
 6. Incubate at room temperature for 30 minutes, protected from light.
 7. Run the samples on a flow cytometer. For tips for flow cytometry detection of purified EVs read "Considerations for Detecting EVs by Flow Cytometry" on page 1.

Staining of purified EVs for fNTA

For analysis on an fNTA device, such as ZetaView®, follow steps 1-6 in the "Staining of purified EVs" protocol above. Make sure to include the preparation of the mock sample (dye in buffer).

After step 6, prepare dilutions of the stained EVs and control dye in 5-10 mL filtered water. We recommend starting with a 1:1000 dilution, but depending on your sample, the optimum dilution could be anywhere from 1:500-1:5000. When running on the fNTA instrument, choose the particular dilution factor that falls within the recommended particle concentration range. It is desirable that for a particular concentration, the mock dye sample has fewer than 10% of the number of particles of the stained EV samples.

Antibody co-staining of purified EVs

This protocol was developed for staining purified EVs with both ExoBrite™ True EV Membrane Stains and fluorescent antibodies, and detecting them by flow cytometry.

Note: Use labeled primary antibodies at the manufacturer's recommended concentration, or try staining in the range of 0.1-5 ug/mL. Either co-incubation or sequential incubations can be performed as described below.

1. Follow steps 1-4 in the "Staining of purified EVs" protocol. In addition to the antibody and ExoBrite™ True EV Membrane Stain co-stained EV samples, it is helpful to include the following controls (if using multiple antibodies, include "buffer plus antibody" and single-stain controls for each antibody).
 - Buffer controls
 - a. Buffer alone (no EVs, no stain)
 - b. Buffer plus ExoBrite™ True EV Membrane Stain
 - c. Buffer plus antibody
 - EV controls
 - a. Unstained EVs
 - b. Single-stain ExoBrite™ True EV Membrane Stain
 - c. Single-stain antibody
2. Choose whether to co-stain by co-incubation (proceed to step 3) or sequential incubation (proceed to step 4).
3. Co-incubation of antibodies and ExoBrite™ True EV Membrane Stain:
 - a. Add 450 uL of 1X ExoBrite™ staining solution to each tube containing 50 uL of EVs. Remember to also add the staining solution to the "buffer plus ExoBrite™" control and the ExoBrite™ single-stain control tubes.
 - b. Add fluorescent antibody conjugate to the samples at the desired concentration. For example, to the 500 uL staining reaction, add 0.5 ug antibody for 1 ug/mL. Remember to also add the antibody to the "buffer plus antibody" control and the antibody single-stain control tubes.
 - c. Continue to steps 6-7 in the "Staining of purified EVs" protocol.
4. Sequential incubation of antibodies and ExoBrite™ True EV Membrane Stain:
 - a. Add fluorescent antibody conjugate to the samples at the desired concentration. For example, to a 50 uL EV sample, add 0.05 ug antibody for 1 ug/mL. Remember to also add the antibody to the "buffer plus antibody" control and the antibody single-stain control tubes.
 - b. Incubate at room temperature for 30 minutes, protected from light.
 - c. Add 450 uL of 1X ExoBrite™ True EV Membrane staining solution to each sample tube. Remember to also add the staining solution to the "buffer plus ExoBrite™" control and the ExoBrite™ True EV Membrane Stain single-stain control tubes.
 - d. Continue to steps 6-7 in the "Staining of purified EVs" protocol.

ExoBrite™ 515/540 True EV Membrane Stain

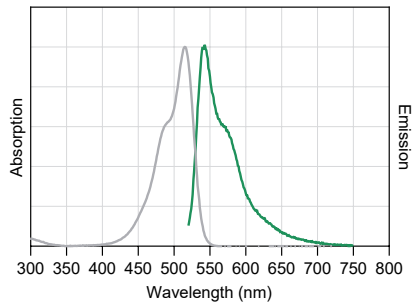


Figure 1. Absorption and emission spectra of ExoBrite™ 515/540 True EV Membrane Stain.

ExoBrite™ 555/575 True EV Membrane Stain

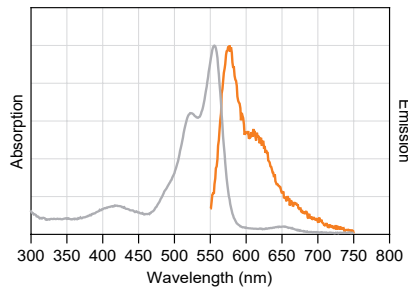


Figure 2. Absorption and emission spectra of ExoBrite™ 555/575 True EV Membrane Stain.

Related Products

Cat. No.	Product
30111-30114	ExoBrite™ CTB EV Staining Kits
30119-30122	ExoBrite™ Annexin EV Staining Kits
30123-30126	ExoBrite™ WGA EV Staining Kits
30115-30118	ExoBrite™ STORM CTB EV Staining Kits
30127	ExoBrite™ EV Surface Stain Sampler Kit, Green
28000	ExoBrite™ Streptavidin Magnetic Beads
P003-410	ExoBrite™ 410/450 CD9 Flow Antibody
P003-490	ExoBrite™ 490/515 CD9 Flow Antibody
P003-560	ExoBrite™ 560/585 CD9 Flow Antibody
P003-650	ExoBrite™ 650/665 CD9 Flow Antibody
P003-RPE	ExoBrite™ R-PE CD9 Flow Antibody
P004-410	ExoBrite™ 410/450 CD63 Flow Antibody
P004-490	ExoBrite™ 490/515 CD63 Flow Antibody
P004-560	ExoBrite™ 560/585 CD63 Flow Antibody
P004-RPE	ExoBrite™ R-PE CD63 Flow Antibody
P005-410	ExoBrite™ 410/450 CD81 Flow Antibody
P005-490	ExoBrite™ 490/515 CD81 Flow Antibody
P005-560	ExoBrite™ 560/585 CD81 Flow Antibody
P005-RPE	ExoBrite™ R-PE CD81 Flow Antibody
P008-410	ExoBrite™ 410/450 IgG1 Isotype Control Flow Antibody
P008-490	ExoBrite™ 490/515 IgG1 Isotype Control Flow Antibody
P008-560	ExoBrite™ 560/585 IgG1 Isotype Control Flow Antibody
P008-650	ExoBrite™ 650/665 IgG1 Isotype Control Flow Antibody
P008-RPE	ExoBrite™ R-PE IgG1 Isotype Control Flow Antibody
P003-680	ExoBrite™ 680/700 CD9 Western Antibody
P003-770	ExoBrite™ 770/800 CD9 Western Antibody
P004-680	ExoBrite™ 680/700 CD63 Western Antibody
P004-770	ExoBrite™ 770/800 CD63 Western Antibody
P006-680	ExoBrite™ 680/700 CD81 Western Antibody
P006-770	ExoBrite™ 770/800 CD81 Western Antibody
P007-770	ExoBrite™ 770/800 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.

Cy dyes are a registered trademark of Cytiva. CytoFLEX is a registered trademark of Beckman Coulter, Inc. ZetaView is a registered trademark of Particle Metrix.

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