

Detect EVs, Not Dye Aggregates

Discover ExoBrite™ EV Stains & Antibodies

ExoBrite™ True EV Membrane Stains - The Best Choice for Pan-EV Labeling

Membrane dyes are frequently used to detect EVs by labeling EV membranes. However, commonly used carbocyanine dyes such as PKH, DiO, and DiI have poor solubility and thus produce poor staining or form aggregates that can be mistaken for EVs. Other membrane dyes, such as di-8-ANEPPS, simply don't generate a signal that is robust or bright enough for efficient detection.

ExoBrite™ True EV Membrane Stains are novel lipophilic dyes designed to address these challenges by offering bright signal and clear differentiation of EVs from non-specific particles by flow cytometry. The stains are offered in four colors (Fig. 1) and can be used with antibodies for multi-parameter analysis. Flow cytometry and fluorescence nanoparticle tracking analysis (fNTA) demonstrated ExoBrite™ 515/540 to be a superior choice over PKH67 with higher (near-complete) pan-EV staining (Fig. 2).

Features of ExoBrite™ True EV Membrane Stains

- Lipophilic membrane dyes designed for pan-EV labeling
- Superior alternatives to PKH, DiO, DiI, DiD, and ANEPPS
- Near-complete coverage of EVs in a sample
- Stain EVs from all sources tested (see table, next page)
- Compatible with antibody co-staining
- Available for Pacific Blue™, FITC, PE, and APC channels

ExoBrite™ True EV Membrane Stains

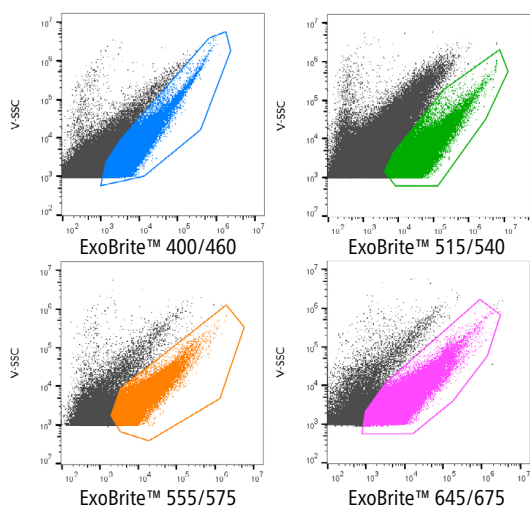
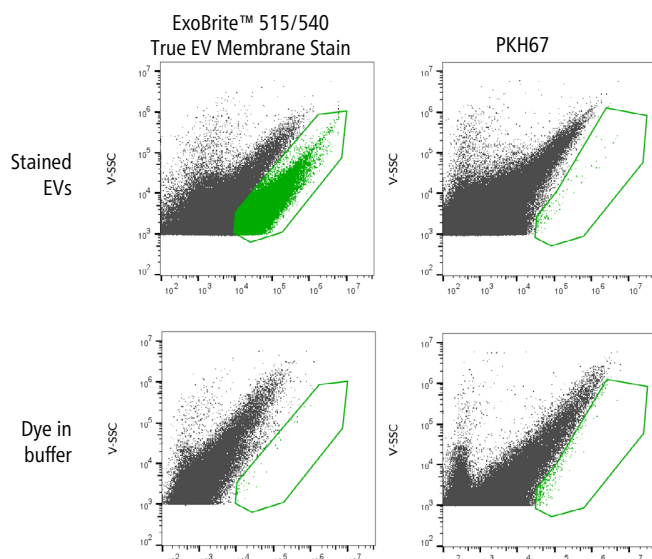


Figure 1. SEC-purified, MCF-7 cell-derived EVs stained with ExoBrite™ True EV Membrane Stains. Samples were run on a CytoFLEX® LX flow cytometer in the Pacific Blue™, FITC, PE, and APC channels.

ExoBrite™ 515/540 True vs. PKH67 in flow



ExoBrite™ 515/540 True vs. PKH67 in fNTA

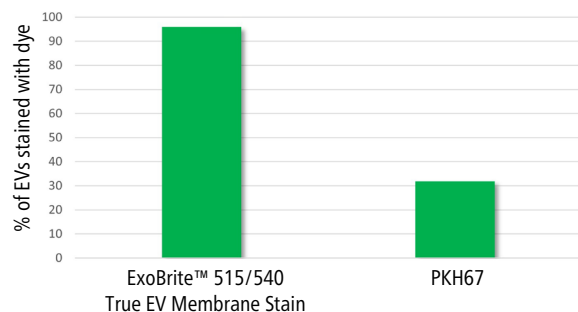


Figure 2. Flow cytometry and fNTA analysis of SEC-purified, MCF-7-derived EVs stained with ExoBrite™ 515/540 True EV Membrane Stain and PKH67. Top: Analysis on a CytoFLEX™ LX flow cytometer shows ExoBrite™ 515/540 True EV Membrane Stain was able to differentiate EVs from background particles, unlike PKH67. Bottom: In ZetaView® fNTA, ExoBrite™ 515/540 True EV Membrane Stain showed nearly complete coverage of EVs, much higher than PKH67.

ExoBrite™ EV Surface Stains

Optimized Annexin V, WGA, & CTB conjugates designed for bright staining of EVs

ExoBrite™ EV surface stains are fluorescent protein conjugates that bind to specific molecules commonly found on EV surfaces. The stains are available as Annexin V, wheat germ agglutinin (WGA), or cholera toxin B (CTB) conjugates, and were developed specifically for bright and sensitive staining of isolated EVs by flow cytometry. The conjugates display much lower background aggregation than most other EV stains and are suitable for antibody co-staining (Fig. 3 and Fig. 4). Biotium also developed ExoBrite™ STORM CTB EV Stains for examining EV morphology in fine detail with stochastic optical reconstruction microscopy (STORM) imaging (see table on page 4). ExoBrite™ EV surface stains have been found to label EVs derived from several tested cell lines, but may not stain EVs from every source. See the ExoBrite™ Stain Compatibility with EVs from Different Sources table below.

Features of ExoBrite™ EV Surface Stains

- Optimized conjugates of Annexin V, WGA, and CTB for EV labeling
- Bright staining of purified EVs with minimal dye aggregation
- Compatible with antibody co-staining
- Options for staining bead-bound EVs
- Available for Pacific Blue™, FITC, PE, and APC channels

ExoBrite™ Stain Compatibility with EVs from Different Sources

| 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 ² | ND | No ¹ | ND | ND |
| HeLa cells | Yes | No | Yes | Yes |
| HUVEC ³ | 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 | Yes | ND | ND | ND |
| Serum | ND | No | ND | Yes |
| Skeletal myoblasts | ND | Yes ¹ | ND | ND |
| THP-1 | Yes | ND | ND | ND |
| U2OS cells | Yes | No | Yes | Yes |
| U937 cells | Yes | No | Yes | Yes |

¹Customer-reported data

²Human adipose stem cells

³Human umbilical vein endothelial cells

Entry of "Yes" or "No" indicates coverage of EVs based on Biotium's internal data or customer-reported data. Value of "ND" indicates no data.

Scan QR code to view the latest information on our ExoBrite™ EV Stains



ExoBrite™ EV Surface Stain Comparison

| ExoBrite™ EV Surface Stain | Pros | Cons |
|------------------------------------|---|--|
| ExoBrite™ Annexin EV Staining Kits | <ul style="list-style-type: none"> • Broad compatibility with EVs from different sources • Validated for flow and fNTA • Low background aggregates | <ul style="list-style-type: none"> • May not stain every EV in a sample • Not recommended for bead-bound EVs |
| ExoBrite™ WGA EV Staining Kits | <ul style="list-style-type: none"> • Broad compatibility with EVs from different sources • Can be used with bead-bound EVs | <ul style="list-style-type: none"> • May not stain every EV in a sample • Not recommended for fNTA |
| ExoBrite™ CTB EV Staining Kits | <ul style="list-style-type: none"> • Validated for flow and fNTA • Extremely low background | <ul style="list-style-type: none"> • May not stain every EV in a sample • Does not stain EVs from every source |

Strong Overlap with Tetraspanin-Positive EVs

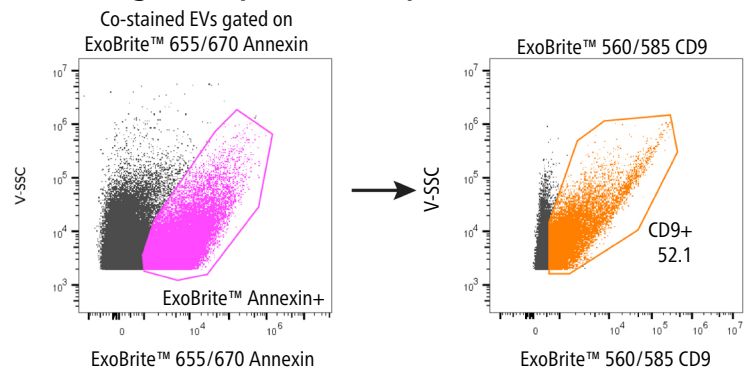


Figure 3. SEC-purified, HeLa cell-derived EVs were stained first with ExoBrite™ 560/585 CD9 Flow Antibody, followed by ExoBrite™ 655/670 Annexin EV Stain. EVs were detected on a CytoFLEX LX flow cytometer. Gating on Annexin-positive particles, ~50% were also positive for CD9.

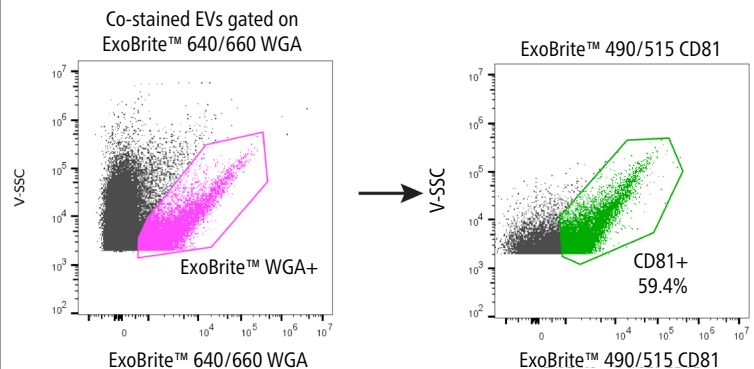


Figure 4. SEC-purified, MCF-7 cell-derived EVs were stained first with ExoBrite™ 490/515 CD81 Flow Antibody, followed by ExoBrite™ 640/660 WGA EV Stain. EVs were detected on a CytoFLEX LX flow cytometer. Gating on WGA-positive particles, ~60% were also positive for CD81.

ExoBrite™ Flow Antibody Conjugates

Validated for detection of EV markers on purified or bead-bound EVs

The tetraspanin family members CD9, CD63, and CD81 are widely used markers for EVs. While other companies may sell tetraspanin antibodies, few are validated for detection of EVs by flow cytometry. ExoBrite™ Flow Antibody Conjugates were curated and validated for flow cytometry to offer bright signal and low background for EV markers on purified and bead-bound EVs.

Features of ExoBrite™ Flow Antibody Conjugates

- Antibody conjugates against human and mouse tetraspanins
- Validated for staining purified and bead-bound EVs
- ExoBrite™ Isotype Control Flow Antibody available
- Options available for Pacific Blue™, FITC, PE, and APC channels

ExoBrite™ Western Antibody Conjugates

Validated for detection of EV markers in extracts

ExoBrite™ Western Antibodies were developed to offer bright signal and low background for EV markers CD9, CD63, and CD81 by near-IR fluorescent western blot. The antibodies are available with the near-IR dyes ExoBrite™ 680/700 and ExoBrite™ 770/800 for optimal signal-to-noise in near-IR fluorescent western detection. ExoBrite™ Calnexin Antibody Conjugates are offered as a cytoplasmic target control for EV purity.

Features of ExoBrite™ Western Antibody Conjugates

- Validated for near-IR western blot of EV extracts
- Negative control ExoBrite™ Calnexin Western Antibody available
- Available in 2 near-IR colors for the LI-COR® Odyssey and other near-IR imagers

ExoBrite™ EV Total RNA Isolation Kit

Efficient total RNA extraction from purified EVs

The ExoBrite™ EV Total RNA Isolation Kit is an optimized and easy-to-use kit for isolating total RNA, including mRNA and miRNA, from purified EVs. The isolated EV RNA can then be used for downstream analysis such as qPCR or RNAseq.

Features of ExoBrite™ EV Total RNA Isolation Kit

- Optimized for total RNA extraction from purified EVs
- Compatible with downstream applications such as qPCR or RNAseq
- Simple column-based purification
- No phenol/chloroform or ethanol precipitation steps

Robust Detection of EVs by Flow

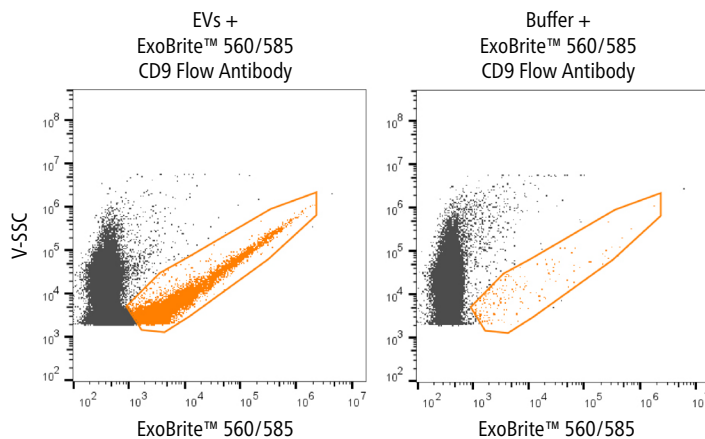


Figure 5. SEC-purified, MCF-7 cell-derived EVs were stained with ExoBrite™ CD9 560/585 Flow Antibody. Specific staining of EVs was seen (left), compared with the same antibody in buffer alone (right). EVs were detected on a CytoFLEX LX flow cytometer in the PE channel.

Conjugates for Near-IR WB of EV Extracts

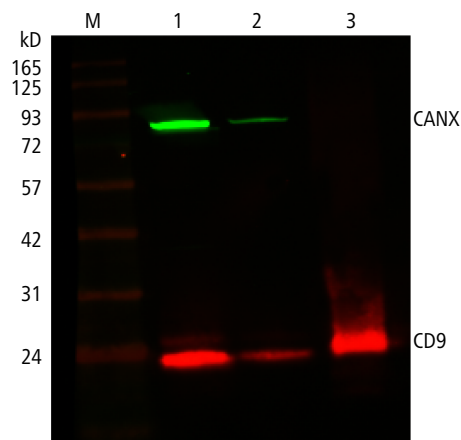


Figure 6. Western detection of human CD9 in MCF-7 cell and EV lysates using ExoBrite™ 680/700 CD9 Western Antibody (red) and ExoBrite™ 770/800 Calnexin Antibody (CANX; EV negative control) (green). Lane M: Protein marker. Lane 1: 10 ug cell lysate. Lane 2: 1 ug cell lysate. Lane 3: 1 ug EV lysate. The blots were imaged on a LI-COR® Odyssey® infrared imaging system.

ExoBrite™ Streptavidin Magnetic Beads

Convenient & optimized capture of EVs

ExoBrite™ Streptavidin Magnetic Beads were developed to offer lower background and higher sensitivity for EV capture and fluorescence detection than similar products.

Features of ExoBrite™ Streptavidin Magnetic Beads

- Streptavidin-coated and magnetic polystyrene beads (4.5 um)
- Combine with a biotinylated antibody (not included) for convenient magnetic bead-based isolation of EVs
- Isolate EVs from biofluids without an overnight precipitation step
- Less autofluorescence than competitor beads
- Compatible with downstream EV analysis methods such as flow cytometry and western blot

Available Colors for ExoBrite™ EV Surface Stains & ExoBrite™ Flow Antibodies

| Conjugates | Ex/Em (nm) | Laser Line(s) (nm) | Detection Channels | Applications |
|---------------------|-------------------|--------------------|--------------------|--------------------------|
| • ExoBrite™ 410/450 | 416/452 | 405 | Pacific Blue™ | Flow |
| • ExoBrite™ 490/515 | 490/516 | 488 | FITC | Flow, fNTA |
| • ExoBrite™ 560/585 | 562/584 | 532 or 561 | PE | Flow, fNTA |
| • ExoBrite™ 640/660 | 642/663 | 633-640 | APC | fNTA (recommended), Flow |
| • ExoBrite™ 650/665 | 652/668 | 633-640 | APC | Flow (recommended), fNTA |
| • ExoBrite™ R-PE | 496, 546, 565/578 | 488, 532, or 561 | PE | Flow |

ExoBrite™ EV Surface Staining Kits

| Product | Size | Cat. No. |
|---|----------------------|---------------|
| ExoBrite™ Annexin EV Staining Kits | 100 or 500 labelings | 30119 - 30122 |
| ExoBrite™ WGA EV Staining Kits | 100 or 500 labelings | 30123 - 30126 |
| ExoBrite™ CTB EV Staining Kits | 100 or 500 labelings | 30111- 30114 |
| ExoBrite™ EV Surface Stain Sampler Kit, Green | 100 labelings | 30127 |

ExoBrite™ Flow Antibodies

| Product | Host | Target | Cat. No. |
|--|---------|-------------------|----------|
| ExoBrite™ CD9 Flow Antibody | Mouse | Human CD9 | P003 |
| ExoBrite™ CD9 (Mouse) Flow Antibody | Rat | Mouse CD9 | P018 |
| ExoBrite™ CD63 Flow Antibody | Mouse | Human CD63 | P004 |
| ExoBrite™ CD63 (Mouse) Flow Antibody | Rat | Mouse CD63 | P022 |
| ExoBrite™ CD81 Flow Antibody | Mouse | Human CD81 | P005 |
| ExoBrite™ CD81 (Mouse/Rat) Flow Antibody | Hamster | Mouse or Rat CD63 | P019 |
| ExoBrite™ IgG1 Isotype Control Flow Antibody | Mouse | Negative control | P008 |

Available in 25 or 100 test sizes.

ExoBrite™ True EV Membrane Stains

| Product | Ex/Em (nm) | Detection Channels | Cat. No. |
|--|------------|--------------------|----------|
| ExoBrite™ 400/460 True EV Membrane Stain | 402/460 | Pacific Blue™ | 30136 |
| ExoBrite™ 515/540 True EV Membrane Stain | 515/542 | FITC | 30129 |
| ExoBrite™ 555/575 True EV Membrane Stain | 556/576 | PE | 30130 |
| ExoBrite™ 645/675 True EV Membrane Stain | 644/671 | APC and Cy®5 | 30137 |

Available in 100 or 500 labeling sizes.

ExoBrite™ Western Antibodies

| Product | Conjugates | Cat. No. |
|-------------------------------------|-------------------|----------|
| ExoBrite™ CD9 Western Antibody | ExoBrite™ 680/700 | P003-680 |
| | ExoBrite™ 770/800 | P003-770 |
| ExoBrite™ CD63 Western Antibody | ExoBrite™ 680/700 | P004-680 |
| | ExoBrite™ 770/800 | P004-770 |
| ExoBrite™ CD81 Western Antibody | ExoBrite™ 680/700 | P006-680 |
| | ExoBrite™ 770/800 | P006-770 |
| ExoBrite™ Calnexin Western Antibody | ExoBrite™ 770/800 | P007-770 |

Available in 25 or 100 test sizes.

ExoBrite™ STORM CTB EV Staining Kits

| Product | Ex/Em (nm) | Cat. No. |
|---|------------|----------|
| ExoBrite™ STORM CF®505 CTB EV Staining Kit | 505/519 | 30115 |
| ExoBrite™ STORM CF®583R CTB EV Staining Kit | 583/609 | 30116 |
| ExoBrite™ STORM CF®647 CTB EV Staining Kit | 652/668 | 30117 |
| ExoBrite™ STORM CF®680 CTB EV Staining Kit | 681/698 | 30118 |

Available in 100 or 500 labeling sizes.

Other Products for EV Research

| Product | Size | Cat. No. |
|---------------------------------------|----------|----------|
| ExoBrite™ Streptavidin Magnetic Beads | 5 mL | 28000 |
| ExoBrite™ EV Total RNA Isolation Kit | 50 preps | 28001 |

Pacific Blue is a trademark of Thermo Fisher Scientific; ZetaView is a registered trademark of Particle-Matrix GmbH; CytoFLEX is a registered trademark of Beckman Coulter, Inc.; LI-COR and Odyssey are registered trademarks of LI-COR Inc; Cy Dye is a registered trademark of Cytiva.