

Product Information

ExoBrite™ Flow Antibody Conjugates

Product List

Product	Cat. No.	Unit Size
ExoBrite™ 410/450 CD9 Flow Antibody	P003-410-125	25 tests
	P003-410-500	100 tests
ExoBrite™ 490/515 CD9 Flow Antibody	P003-490-125	25 tests
	P003-490-500	100 tests
ExoBrite™ 560/585 CD9 Flow Antibody	P003-560-125	25 tests
	P003-560-500	100 tests
ExoBrite™ 650/665 CD9 Flow Antibody	P003-650-125	25 tests
	P003-650-500	100 tests
ExoBrite™ R-PE CD9 Flow Antibody	P003-RPE-125	25 tests
	P003-RPE-500	100 tests
ExoBrite™ 410/450 CD9 (Mouse) Flow Antibody	P018-410-125	25 tests
	P018-410-500	100 tests
ExoBrite™ 490/515 CD9 (Mouse) Flow Antibody	P018-490-125	25 tests
	P018-490-500	100 tests
ExoBrite™ 560/585 CD9 (Mouse) Flow Antibody	P018-560-125	25 tests
	P018-560-500	100 tests
ExoBrite™ 650/665 CD9 (Mouse) Flow Antibody	P018-650-125	25 tests
	P018-650-500	100 tests
ExoBrite™ 410/450 CD63 Flow Antibody	P004-410-125	25 tests
	P004-410-500	100 tests
ExoBrite™ 490/515 CD63 Flow Antibody	P004-490-125	25 tests
	P004-490-500	100 tests
ExoBrite™ 560/585 CD63 Flow Antibody	P004-560-125	25 tests
	P004-560-500	100 tests
ExoBrite™ R-PE CD63 Flow Antibody	P004-RPE-125	25 tests
	P004-RPE-500	100 tests
ExoBrite™ 410/450 CD63 (Mouse) Flow Antibody	P022-410-125	25 tests
	P022-410-500	100 tests
ExoBrite™ 490/515 CD63 (Mouse) Flow Antibody	P022-490-125	25 tests
	P022-490-500	100 tests
ExoBrite™ 560/585 CD63 (Mouse) Flow Antibody	P022-560-125	25 tests
	P022-560-500	100 tests
ExoBrite™ 410/450 CD81 Flow Antibody	P005-410-125	25 tests
	P005-410-500	100 tests
ExoBrite™ 490/515 CD81 Flow Antibody	P005-490-125	25 tests
	P005-490-500	100 tests
ExoBrite™ 560/585 CD81 Flow Antibody	P005-560-125	25 tests
	P005-560-500	100 tests

ExoBrite™ R-PE CD81 Flow Antibody	P005-RPE-125	25 tests
	P005-RPE-500	100 tests
ExoBrite™ 410/450 CD81 (Mouse/Rat) Flow Antibody	P019-410-125	25 tests
	P019-410-500	100 tests
ExoBrite™ 490/515 CD81 (Mouse/Rat) Flow Antibody	P019-490-125	25 tests
	P019-490-500	100 tests
ExoBrite™ 560/585 CD81 (Mouse/Rat) Flow Antibody	P019-560-125	25 tests
	P019-560-500	100 tests
ExoBrite™ 410/450 IgG1 Isotype Control Flow Antibody	P008-410-125	25 tests
	P008-410-500	100 tests
ExoBrite™ 490/515 IgG1 Isotype Control Flow Antibody	P008-490-125	25 tests
	P008-490-500	100 tests
ExoBrite™ 560/585 IgG1 Isotype Control Flow Antibody	P008-560-125	25 tests
	P008-560-500	100 tests
ExoBrite™ 650/665 IgG1 Isotype Control Flow Antibody	P008-650-125	25 tests
	P008-650-500	100 tests
ExoBrite™ R-PE IgG1 Isotype Control Flow Antibody	P008-RPE-125	25 tests
	P008-RPE-500	100 tests

Storage and Handling

Store at 4°C, protected from light. Product is stable for at least 24 months from date of receipt when stored as recommended.

Note: Storage of the antibody for more than a day at final working dilution is not recommended.

Product Description

Extracellular vesicles (EVs), including exosomes, are widely studied for their potential use in drug delivery and medical diagnostic applications. The most common proteins used as EV markers are CD9, CD63, and CD81, members of the tetraspanin family. Tetraspanins are plasma membrane proteins with many proposed functions, including activation and sorting of other membrane proteins. They are also thought to play a role in the targeting of proteins to multivesicular bodies (MVBs) and EVs. These tetraspanins are broadly expressed on many cell types and can therefore be detected on many types of EVs, but their expression levels vary depending on the cell type of origin.

ExoBrite™ Flow Antibody Conjugates are validated by Biotium for optimal detection of EV markers CD9, CD63, and CD81 in purified or bead-bound EVs by flow cytometry. The antibodies are available in different clones for both human and mouse tetraspanins. ExoBrite™ fluorophores offer exceptional brightness and signal-to-noise over alternative fluorophores (see Table 2 on page 3 for detection settings of each ExoBrite™ conjugate).

ExoBrite™ Isotype Control Flow Antibodies are offered as a negative control for the ExoBrite™ mouse anti-human antibodies. The isotype controls have no known reactivity with any target in human cells, and have the same isotype as the mouse anti-human ExoBrite™ antibodies.

Biotium also offers other products for EV research, including ExoBrite™ conjugates of CTB, WGA, and Annexin V, as well as ExoBrite™ True EV Membrane Stains for pan-EV labeling. See Related Products and visit our [technology page](#) for more information.

Considerations for Detecting EVs by Flow Cytometry

- Obtaining a clean EV prep is crucial for obtaining robust signal and proper interpretation of results. While there are several EV isolation methods, we have found that size exclusion chromatography (SEC) is an accessible and easy-to-use method that yields a relatively pure population of EVs. For a comparison of EV isolation methods and protocols for EV isolation and staining, see the following tech tips:
[Tech Tip: Isolation and Staining of Extracellular Vesicles](#)
[Tech Tip: Fluorescent Detection of EVs by Flow Cytometry](#)
- EVs are extremely small vesicles (~30-150 nm in diameter), which is near or below the size 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 μ m) in SSC before attempting to detect purified EVs. Sizing beads should be used to set the SSC threshold before each EV detection experiment. EVs that are bound to affinity beads are large enough to detect on any instrument.
- To improve the sensitivity for detecting small particles, we recommend using the 405 nm laser for the SSC channel if it is an option on your flow cytometer.
- For best results, buffers used for suspending and staining EVs should be filtered through a 0.2 μ m filter to remove particulates.
- To reduce instrument noise, use a low flow rate to keep the event rate and abort rate low. Dilute the stained samples in filtered PBS if necessary.

Considerations for ExoBrite™ Flow Antibody Conjugates

The following are general considerations for using ExoBrite™ Flow Antibodies to stain EVs. See Experimental Protocols for step-by-step instructions for use.

- ExoBrite™ Flow Antibody Conjugates have been validated in flow cytometry on the CytoFLEX LX from Beckman Coulter. Results on other instruments may vary based on the instruments size detection limit and other parameters.
- ExoBrite™ Flow Antibody Conjugates have been validated for staining EVs isolated using several different methods, including PEG precipitation, size exclusion chromatography, and affinity bead isolation. Staining results may vary depending on the EV isolation method used.

Table 1. Antibody Attributes

Antibody	Target	Host	Species Reactivity	Target MW	Isotype	Entrez Gene ID	SwissProt	Unigene	Synonyms	Target Localization
ExoBrite™ CD9 Flow Antibody	CD9	Mouse	Human, Baboon, Bovine, Cynomolgus monkey, Dog, Horse, Rabbit, Non-human primates, Sheep	24 kDa	IgG1, kappa	928	P21926	114286	Tspan-29, MRP-1	Exosomes/EVs, Plasma membrane
ExoBrite™ CD9 (Mouse) Flow Antibody	CD9	Rat	Mouse	24 kDa	IgG2a, kappa	12527	P40240	-	DRAP-27, MRP-1, p-24	
ExoBrite™ CD63 Flow Antibody	CD63	Mouse	Human, Baboon, Cynomolgus monkey, Non-human primates	26 kDa (core protein); 30-60 kDa (glycosylated)	IgG1, kappa	967	P08962	445570	Tspan-30, LAMP-3, gp55	Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies
ExoBrite™ CD63 (Mouse) Flow Antibody	CD63	Rat	Mouse	53 kDa	IgG2a, kappa	12512	P41731	-	LIMP, LAMP-3, gp55, ME491	
ExoBrite™ CD81 Flow Antibody	CD81	Mouse	Human, Baboon, Cynomolgus monkey, Non-human primates, Mouse (low reactivity)	26 kDa	IgG1, kappa	975	P60033	54457	Tspan-28, TAPA-1	Exosomes/EVs, Plasma membrane
ExoBrite™ CD81 (Mouse/Rat) Flow Antibody	CD81	Hamster	Mouse, Rat	26 kDa	IgG1, kappa	12520 (mouse), 25621 (rat)	P35762 (mouse), Q62745 (rat)	-	TAPA-1	
ExoBrite™ IgG1 Isotype Control Flow Antibody	-	Mouse	-	-	IgG1, kappa	-	-	-	-	-

Experimental Protocols

Note: Before beginning, please read "Considerations for ExoBrite™ Flow Antibody Conjugates" section on page 2.

Antibody staining of purified EVs

This protocol was developed for staining purified EVs with ExoBrite™ Flow Antibody Conjugates for detection by flow cytometry.

1. Isolate or purify EVs using the procedure of your choice. Thaw EVs if they have been stored frozen.
2. Aliquot 100 µL of EV sample into FACS tubes or microcentrifuge tubes. We recommend setting up control tubes of antibody in buffer alone, as well as isotype controls (if available).

Buffer controls

- a. Buffer alone (no EVs, no antibody)
- b. Buffer plus ExoBrite™ Flow Antibody

EV samples

- a. Unstained EVs
 - b. ExoBrite™ Flow Antibody
 - c. ExoBrite™ Isotype Control (if available)
3. Add 5 µL of ExoBrite™ antibody to each 100 µL sample. Remember to also add the antibody to the buffer plus ExoBrite™ Flow Antibody control.
 4. Incubate at room temperature for 30 minutes, protected from light.
 5. Add filtered PBS to the desired volume and 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 2. See Table 2 for recommended detection settings for ExoBrite™ Flow Antibody Conjugates.

Antibody staining of bead-bound EVs

This protocol was developed for EVs bound to magnetic antibody capture beads, stained with ExoBrite™ Flow Antibody Conjugates and detected by flow cytometry.

1. Prepare EVs bound to the magnetic capture beads of your choice, according to the manufacturer's recommended procedure.
2. Prepare sample tubes and the following control tubes:

Beads controls (no EVs)

- a. Beads alone
- b. Beads plus ExoBrite™ Flow Antibody

Bead-bound EV samples

- a. Unstained bead-bound EVs
 - b. ExoBrite™ Flow Antibody
 - c. ExoBrite™ Isotype Control (if available)
3. Place the tubes with bead-bound EVs on a magnet for 1 minute, remove and discard the supernatant.
Note: If the beads are not completely recovered from the buffer after 1 minute on the magnet, leave the tubes on the magnet for a longer time (up to 4 minutes). We have found that briefly centrifuging tubes to collect the contents near the bottom before placing them on the magnet can improve bead recovery.
 4. Remove the tubes from the magnet, add 300 µL of 0.2 µm-filtered PBS and gently pipet up and down to resuspend.
 5. Repeat steps 3-4 once.
 6. Remove the tubes from the magnet and suspend in 100 µL of PBS. Add 5 µL of ExoBrite™ antibody to each sample, including applicable controls.
 7. Incubate at room temperature for 30 minutes, protected from light.
 8. Place the tube on the magnet for 1 minute and discard the supernatant.
 9. Remove the tubes from the magnet, add 300 µL of PBS and gently pipet up and down to resuspend.
 10. Place the tube on the magnet for 1 minute and discard the supernatant.
 11. Remove the tubes from the magnet, add 500-900 µL of PBS and transfer to flow tubes.
 12. Run the samples on a flow cytometer. See Table 2 for recommended detection settings for ExoBrite™ Flow Antibody Conjugates.

Table 2. Detection Settings for ExoBrite™ Flow Antibodies

Conjugate	Ex/Em (nm)	Laser Line(s) (nm)	Detection Channel
ExoBrite™ 410/450	411/452	405	Pacific Blue™
ExoBrite™ 490/515	490/516	488	FITC
ExoBrite™ 560/585	562/584	532 or 561	PE
ExoBrite™ 650/665	652/668	633-640	APC
ExoBrite™ R-PE	496, 546, 565/578	488, 532, or 561	PE

Related Products

Cat. No.	Product
30129, 30130	ExoBrite™ True EV Membrane Stains
30111-30114	ExoBrite™ CTB EV Staining Kits
30123-30126	ExoBrite™ WGA EV Staining Kits
30119-30122	ExoBrite™ Annexin EV Staining Kits
30115-30118	ExoBrite™ STORM CTB EV Staining Kits
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
28000	ExoBrite™ Streptavidin Magnetic Beads
28001	ExoBrite™ EV Total RNA Isolation Kit

Please visit our website at www.biotium.com for more information on our products for EV detection and western blotting including EV stains and antibodies for flow cytometry, western blot blocking buffers, and total protein stains.

Pacific Blue is a trademark of Thermo Fisher Scientific. CYTOFLEX is a registered trademark of Beckman Coulter, Inc.

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