

## ExoBrite™ CD9/CD63/CD81 3-Color Antibody Cocktails

Antibody mixtures developed with optimal brightness and signal-to-noise for phenotyping EV samples. Available in blue/green/far-red or green/orange/far-red.



### Product Description

The ExoBrite™ CD9/CD63/CD81 3-Color Antibody Cocktails were developed for phenotyping EV samples by flow cytometry. The mixtures have been meticulously optimized for optimal brightness and signal-to-noise. The cocktails are available in blue/green/far-red or green/orange/far-red.

#### Skip the optimization, get results faster

The most commonly used markers for extracellular vesicle (EV) detection are the three tetraspanin proteins CD9, CD63, and CD81. Depending on the source of the EVs, each of these proteins will be present at varying levels. The ExoBrite CD9/CD63/CD81 3-Color Antibody Cocktails were designed for fast, reproducible phenotyping of an EV sample by flow cytometry. Each individual antibody in the mixture has been validated to have the same coverage as its respective ExoBrite™ Flow Antibody, ensuring accurate results.

In addition, like all ExoBrite™ Flow antibodies, the ExoBrite™ cocktails are provided in a proprietary buffer formulation for reduced antibody aggregation and brighter EV staining for optimal accuracy and signal-to-noise. The cocktails can be used on their own, for tetraspanin profiling, or combined with other antibodies of interest for more multi-target analyses.

#### EV antibodies you can trust

Other commercially available antibodies for human tetraspanin proteins CD9, CD63, and CD81 are generally not validated for isolated EVs and may require tedious optimization for your EV prep and staining protocol. The antibodies and dye labels of ExoBrite™ Flow Antibody Conjugates were carefully selected and validated for robust detection of isolated EVs. See the table below for a full list of available [ExoBrite™ Flow Antibody Conjugates](#). [ExoBrite™ Isotype Control Flow Antibodies](#) are also available as a negative controls.

#### Accelerate your EV research with ExoBrite™ EV stains

For general EV staining, Biotium's [ExoBrite™ True EV Membrane Stains](#) offer unparalleled coverage of EVs in a sample and address issues of dye aggregation often seen with PKH and other common membrane dyes. Biotium also offers optimized ExoBrite™ EV Surface stains conjugated to [cholera toxin B \(CTB\)](#), [wheat germ agglutinin \(WGA\)](#), and [Annexin V](#). These stains are specially formulated for bright and specific detection of isolated EVs by flow cytometry. These [ExoBrite™ EV Surface Stains](#) may also be combined with antibody staining, for multi-parameter analysis.

Biotium also offers conjugated ExoBrite™ Western Antibodies against human [CD9](#), [CD63](#), and [CD81](#) designed for optimal detection in EV extracts by fluorescent western blot.

**Note:** 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.

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

### Product attributes

Antibody number	#P028, #P029
SwissProt	P21926, P08962, P60033
Antibody type	ExoBrite™, Primary
Clonality	Monoclonal
Host species	Mouse
Isotype	IgG1, kappa
Antibody reactivity (target)	CD63, CD81, CD9
Species reactivity	Human
Human gene symbol	CD9, CD63, CD81
Entrez gene ID	928, 967, 975
Unigene	114286, 445570, 54457
Antibody target cellular localization	Exosomes/EVs, Plasma membrane
Verified antibody applications	Flow (surface) (verified)
Positive control	MCF-7 cells, MCF-7 derived EVs
Antibody application notes	Recommended concentration for flow cytometry: 5 uL per 0.1 mL exosomes
Antibody research areas	Exosomes/EVs
Antibody/conjugate formulation	Proprietary buffer containing 0.05% sodium azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light
Regulatory status	For research use only (RUO)
Shipping condition	Room temperature

## ExoBrite™ Flow Anti-Human CD9/CD63/CD81 Antibody Cocktails

Cocktail	Ex/Em	Detection Channel	Size	Catalog No.
<a href="#">ExoBrite™ CD9/CD63/CD81 Single-Color Antibody Cocktail (Green)</a>	490/516 nm	FITC	25 tests	P030-125
100 tests	P030-500			
<a href="#">ExoBrite™ CD9/CD63/CD81 Single-Color Antibody Cocktail (Orange)</a>	562/584 nm	PE	25 tests	P031-125
100 tests	P031-500			
<a href="#">ExoBrite™ CD9/CD63/CD81 Single-Color Antibody Cocktail (Far-Red)</a>	642/663 nm	APC	25 tests	P032-125
100 tests	P032-500			
<a href="#">ExoBrite™ CD9/CD63/CD81 3-Color Antibody Cocktail (Blue/Green/Far-Red)</a>	411/452 nm 490/516 nm 642/663 nm	Pacific Blue® FITC APC	25 tests	P028-125
100 tests	P028-500			
<a href="#">ExoBrite™ CD9/CD63/CD81 3-Color Antibody Cocktail (Green/Orange/Far-Red)</a>	490/516 nm 562/584 nm 642/663 nm	FITC PE APC	25 tests	P029-125
100 tests	P029-500			

## ExoBrite™ Flow Antibody Conjugates

Antibody	Ex/Em	Target	Species Reactivity	Detection Channel	Catalog No.
<a href="#">ExoBrite™ 410/450 CD9 Flow Antibody</a>	416/452 nm	CD9	Human	Pacific Blue™	P003-410
<a href="#">ExoBrite™ 490/515 CD9 Flow Antibody</a>	490/516 nm	CD9	Human	FITC	P003-490
<a href="#">ExoBrite™ 560/585 CD9 Flow Antibody</a>	562/584 nm	CD9	Human	PE	P003-560
<a href="#">ExoBrite™ 640/660 CD9 Flow Antibody</a>	642/663 nm	CD9	Human	APC	P003-640
<a href="#">ExoBrite™ R-PE CD9 Flow Antibody</a>	496, 546, 565/578 nm	CD9	Human	PE	P003-RPE
<a href="#">ExoBrite™ APC CD9 Flow Antibody</a>	651/660 nm	CD9	Human	APC	P003-APC
<a href="#">ExoBrite™ 410/450 CD9 (Mouse) Flow Antibody</a>	416/452 nm	CD9	Mouse	Pacific Blue™	P018-410
<a href="#">ExoBrite™ 490/515 CD9 (Mouse) Flow Antibody</a>	490/516 nm	CD9	Mouse	FITC	P018-490
<a href="#">ExoBrite™ 560/585 CD9 (Mouse) Flow Antibody</a>	562/584 nm	CD9	Mouse	PE	P018-560
<a href="#">ExoBrite™ 640/660 CD9 (Mouse) Flow Antibody</a>	642/663 nm	CD9	Mouse	APC	P018-640
<a href="#">ExoBrite™ APC CD9 (Mouse) Flow Antibody</a>	651/660 nm	CD9	Mouse	APC	P018-APC
<a href="#">ExoBrite™ 410/450 CD47 Flow Antibody</a>	416/452 nm	CD47	Human, Mouse	Pacific Blue™	P061-410
<a href="#">ExoBrite™ 490/515 CD47 Flow Antibody</a>	490/516 nm	CD47	Human, Mouse	FITC	P061-490
<a href="#">ExoBrite™ 560/585 CD47 Flow Antibody</a>	562/584 nm	CD47	Human, Mouse	PE	P061-560
<a href="#">ExoBrite™ 640/660 CD47 Flow Antibody</a>	642/663 nm	CD47	Human, Mouse	APC	P061-640
<a href="#">ExoBrite™ 410/450 CD63 Flow Antibody</a>	416/452 nm	CD63	Human	Pacific Blue™	P004-410
<a href="#">ExoBrite™ 490/515 CD63 Flow Antibody</a>	490/516 nm	CD63	Human	FITC	P004-490
<a href="#">ExoBrite™ 560/585 CD63 Flow Antibody</a>	562/584 nm	CD63	Human	PE	P004-560
<a href="#">ExoBrite™ 640/660 CD63 Flow Antibody</a>	642/663 nm	CD63	Human	APC	P004-640
<a href="#">ExoBrite™ R-PE CD63 Flow Antibody</a>	496, 546, 565/578	CD63	Human	PE	P004-RPE

Antibody	Ex/Em	Target	Species Reactivity	Detection Channel	Catalog No.
<a href="#">ExoBrite™ APC CD63 Flow Antibody</a>	651/660 nm	CD63	Human	APC	P004-APC
<a href="#">ExoBrite™ 410/450 CD63 (Mouse) Flow Antibody</a>	416/452 nm	CD63	Mouse	Pacific Blue™	P022-410
<a href="#">ExoBrite™ 490/515 CD63 (Mouse) Flow Antibody</a>	490/516 nm	CD63	Mouse	FITC	P022-490
<a href="#">ExoBrite™ 560/585 CD63 (Mouse) Flow Antibody</a>	562/584 nm	CD63	Mouse	PE	P022-560
<a href="#">ExoBrite™ 640/660 CD63 (Mouse) Flow Antibody</a>	642/663 nm	CD63	Mouse	APC	P022-640
<a href="#">ExoBrite™ APC CD63 (Mouse) Flow Antibody</a>	651/660 nm	CD63	Mouse	APC	P022-APC
<a href="#">ExoBrite™ 410/450 CD81 Flow Antibody</a>	416/452 nm	CD81	Human	Pacific Blue™	P005-410
<a href="#">ExoBrite™ 490/515 CD81 Flow Antibody</a>	490/516 nm	CD81	Human	FITC	P005-490
<a href="#">ExoBrite™ 560/585 CD81 Flow Antibody</a>	562/584 nm	CD81	Human	PE	P005-560
<a href="#">ExoBrite™ 640/660 CD81 Flow Antibody</a>	642/663 nm	CD81	Human	APC	P005-640
<a href="#">ExoBrite™ R-PE CD81 Flow Antibody</a>	496, 546, 565/578	CD81	Human	PE	P005-RPE
<a href="#">ExoBrite™ APC CD81 Flow Antibody</a>	651/660 nm	CD81	Human	APC	P005-APC
<a href="#">ExoBrite™ 410/450 CD81 (Mouse/Rat) Flow Antibody</a>	416/452 nm	CD81	Mouse, Rat	Pacific Blue™	P019-410
<a href="#">ExoBrite™ 490/515 CD81 (Mouse/Rat) Flow Antibody</a>	490/516 nm	CD81	Mouse, Rat	FITC	P019-490
<a href="#">ExoBrite™ 560/585 CD81 (Mouse/Rat) Flow Antibody</a>	562/584 nm	CD81	Mouse, Rat	PE	P019-560
<a href="#">ExoBrite™ 640/660 CD81 (Mouse/Rat) Flow Antibody</a>	642/663 nm	CD81	Mouse, Rat	APC	P019-640
<a href="#">ExoBrite™ APC CD81 (Mouse) Flow Antibody</a>	651/660 nm	CD81	Mouse, Rat	APC	P019-APC
<a href="#">ExoBrite™ 410/450 CD29 Flow Antibody</a>	416/452 nm	CD29	Human	Pacific Blue™	P040-410
<a href="#">ExoBrite™ 490/515 CD29 Flow Antibody</a>	490/516 nm	CD29	Human	FITC	P040-490
<a href="#">ExoBrite™ 560/585 CD29 Flow Antibody</a>	562/584 nm	CD29	Human	PE	P040-560
<a href="#">ExoBrite™ 640/660 CD29 Flow Antibody</a>	642/663 nm	CD29	Human	APC	P040-640
<a href="#">ExoBrite™ 410/450 IgG1 Isotype Control Flow Antibody</a>	416/452 nm	Isotype Control	None	Pacific Blue™	P008-410
<a href="#">ExoBrite™ 490/515 IgG1 Isotype Control Flow Antibody</a>	490/516 nm	Isotype Control	None	FITC	P008-490
<a href="#">ExoBrite™ 560/585 IgG1 Isotype Control Flow Antibody</a>	562/584 nm	Isotype Control	None	PE	P008-560
<a href="#">ExoBrite™ R-PE IgG1 Isotype Control Flow Antibody</a>	496, 546, 565/578	Isotype Control	None	PE	P008-RPE
<a href="#">ExoBrite™ 640/660 IgG1 Isotype Control Flow Antibody</a>	642/663 nm	Isotype Control	None	APC	P008-640
<a href="#">ExoBrite™ APC IgG1 Isotype Control Flow Antibody</a>	651/660 nm	Isotype Control	None	APC	P008-APC
<a href="#">ExoBrite™ Armenian Hamster IgG Isotype Control Flow Antibody</a>	416/452 nm	Isotype Control	None	Pacific Blue™	P056-410
<a href="#">ExoBrite™ 490/515 Armenian Hamster IgG Isotype Control Flow Antibody</a>	490/516 nm	Isotype Control	None	FITC	P056-490

Antibody	Ex/Em	Target	Species Reactivity	Detection Channel	Catalog No.
<a href="#">ExoBrite™ 560/585 Armenian Hamster IgG Isotype Control Flow Antibody</a>	562/584 nm	Isotype Control	None	PE	P056-560
<a href="#">ExoBrite™ 640/660 Armenian Hamster IgG Isotype Control Flow Antibody</a>	642/663 nm	Isotype Control	None	APC	P056-640
<a href="#">ExoBrite™ Armenian Hamster IgG Isotype Control Flow Antibody</a>	651/660 nm	Isotype Control	None	APC	P056-APC
<a href="#">ExoBrite™ 410/450 Rat IgG2a Isotype Control Flow Antibody</a>	416/452 nm	Isotype Control	None	Pacific Blue™	P057-410
<a href="#">ExoBrite™ 490/515 Rat IgG2a Isotype Control Flow Antibody</a>	490/516 nm	Isotype Control	None	FITC	P057-490
<a href="#">ExoBrite™ 560/585 Rat IgG2a Isotype Control Flow Antibody</a>	562/584 nm	Isotype Control	None	PE	P057-560
<a href="#">ExoBrite™ 640/660 Rat IgG2a Isotype Control Flow Antibody</a>	642/663 nm	Isotype Control	None	APC	P057-640
<a href="#">ExoBrite™ APC Rat IgG2a Isotype Control Flow Antibody</a>	651/660 nm	Isotype Control	None	APC	P057-APC

This datasheet was generated on June 11, 2026 at 08:13:07 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/exobrite-cd9-cd63-cd81-3-color-antibody-cocktails/>