Glowing products for science

# CD9 Recombinant Monoclonal Mouse Antibody (2310.9) - Biotium Choice

alottum

CD9 (clone 2310.9) is a validated recombinant mouse monoclonal antibody that recognizes the human tetraspanin protein CD9. The antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance.

### **Product Description**

CD9 (clone 2310.9) is a recombinant mouse monoclonal antibody that recognizes the human tetraspanin protein CD9. This antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance. This antibody has been validated in flow cytometry, immunofluorescence microscopy, and in western blot. The antibody is available unlabeled or conjugated to CF® Dyes and AstralTM tandem dyes.

CD9 is a type IV transmembrane glycoprotein with four transmembrane domains. CD9 on pre-B cells may play a role in cell-cell adhesion. In addition, CD9 may play a role in signal transduction mediated by interaction with low molecular weight GTP binding proteins. CD9 is expressed on early B cells, eosinophils, basophils and activated T cells and is a major component of the platelet cell surface. It is also expressed on most non-T acute lymphoblastic leukemia cells and on some acute myeloid and chronic lymphoid leukemia.

Tetraspanin proteins CD9, CD63, and CD81, are the most common proteins used as exosome markers. 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 exosomes. These tetraspanins are broadly expressed on many cell types and can therefore be detected on many types of exosomes, but their expression levels vary depending on the cell type of origin.

As an isotype control, we recommend <u>Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)</u> for this antibody.

#### **Exosome Stains and Antibodies**

Biotium also offers ExoBrite<sup>TM</sup> CD9 Flow Antibody and ExoBrite<sup>TM</sup> CD9 Western Antibody that are validated for optimal detection of isolated exosomes by flow cytometry or western blotting. Learn more about our products for exosome research.

## Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies

#### Powered by CF® Dyes and Astral Leap™ Tandem Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes and Astral Leap<sup>TM</sup> tandem dyes, they are our top-recommended antibodies for flow cytometry and other applications.

#### **Biotium Choice Antibody Features**

- · Robust and validated clones against common targets
- Developed and optimized for flow cytometry and other applications
- Conjugated to bright, photostable CF® Dyes for superior signal and clarity
- ◆ Also available with <u>Astral Leap™ tandem dyes</u> for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

View our full catalog of Biotium Choice antibodies

Look for this symbol next to the product name on product pages and search results

#### Call us: 800-304-5357 Email: btinfo@biotium.com

#### Product attributes

Antibody number	#P015			
SwissProt	P21926			
<b>Biotium Choice Antibody</b>	Primary			
Antibody type	Biotium Choice Flow Antibodies, Primary			
Clonality	Recombinant Monoclonal			
Host species	Mouse			
Clone	2310.9			
Isotype	IgG1, kappa			
Antibody reactivity (target)	CD9			
Synonyms	Tetraspanin-29 (TSPAN29); BA-2/p24 antigen; BA2; BTCC1; CD9; Cell growth-inhibiting gene 2 protein; DRAP27; GIG2; Leukocyte antigen MIC3; MIC3; Motility-related protein (MRP1); p24			
Species reactivity	Baboon, Bovine, Cynomolgus monkey, Dog, Horse, Human, Non-human primates, Rabbit, Sheep			
Human gene symbol	CD9			
Entrez gene ID	928			
Unigene	114286			
Molecular weight	24 kDa			
Antibody target cellular localization	Exosomes/EVs, Plasma membrane			
Cell/tissue expression	Exosomes, Platelets, Basophils, Eosinophils, Epithelial cells, Lymphocytes			
	Lymphocytes			
Verified antibody applications	Lymphocytes Flow (surface) (verified), WB (verified), IF (verified)			
	* * *			
applications	Flow (surface) (verified), WB (verified), IF (verified)			
applications Positive control	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells),			
applications  Positive control  Antibody application notes	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.05 ug/mL			
applications  Positive control  Antibody application notes  Antibody research areas  Antibody/conjugate	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.05 ug/mL  Exosomes/EVs  Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5			
applications  Positive control  Antibody application notes  Antibody research areas  Antibody/conjugate formulation	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.05 ug/mL  Exosomes/EVs  Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide  Guaranteed for at least 24 months from date of receipt when			
applications  Positive control  Antibody application notes  Antibody research areas  Antibody/conjugate formulation  Shelf life	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.05 ug/mL  Exosomes/EVs  Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide  Guaranteed for at least 24 months from date of receipt when stored as recommended  Note: store unconjugated formats at -10 °C to -35 °C, Store			
applications  Positive control  Antibody application notes  Antibody/conjugate formulation  Shelf life  Storage Conditions	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.05 ug/mL  Exosomes/EVs  Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide  Guaranteed for at least 24 months from date of receipt when stored as recommended  Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C			
applications  Positive control  Antibody application notes  Antibody/conjugate formulation  Shelf life  Storage Conditions  Volume per assay	Flow (surface) (verified), WB (verified), IF (verified)  MCF-7 cells, MCF-7 derived exosomes  Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.05 ug/mL  Exosomes/EVS  Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide  Guaranteed for at least 24 months from date of receipt when stored as recommended  Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C  5 uL/test			

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<u>CF®405M</u> 416/452 nm	416/452 nm	100 ug/mL	25 tests (125 uL)	P015-405M-125	CF®405M Features
			100 tests (500 uL)	P015-405M-500	
<u>CF®488A</u> 490/516 nm	100 ug/mL	25 tests (125 uL)	P015-488A-125	CF®488A Features	
			100 tests (500 uL)	P015-488A-500	
CF® 568	562/584 nm	100 ug/mL	25 tests (125 uL)	P015-568-125	CF®568 Features
			100 tests (500 uL)	P015-568-500	
<u>CF® 583R</u> 585/609 nm	585/609 nm	100 ug/mL	25 tests (125 uL)	P015-583R-125	CF®583R Features
			100 tests (500 uL)	P015-583R-500	
RPE-Astral <sup>TM</sup> 616	PE-Astral <sup>TM</sup> 616 496, 546, 566/617 nm	100 ug/mL	25 tests (125 uL)	P015-R616-125	RPE-Astral <sup>TM</sup> 616
			100 tests (500 uL)	P015-R616-500	<u>Features</u>
CF®647Plus	<u>8647Plus</u> 652/668 nm	100 ug/mL	25 tests (125 uL)	P015-647P-125	
			100 tests (500 uL)	P015-647P-500	
CF®700	696/721 nm	100 ug/mL	25 tests (125 uL)	P015-700-125	CF®700 Features
			100 tests (500 uL)	P015-700-500	
APC-Astral <sup>TM</sup> 813	<u>Astral™813</u> 633, 638/813 nm	100 ug/mL	25 tests (125 uL)	P015-A813-125	APC-Astral <sup>TM</sup> 813 Features
			100 tests (500 uL)	P015-A813-500	
<u>R-PE</u> 496, 546, 565/576 nm	496, 546, 565/576 nm	6 nm 50 ug/mL	25 tests (125 uL)	P015-RPE-125	
			100 tests (500 uL)	P015-RPE-500	
<u>APC</u> 651/660 nm	651/660 nm	20 ug/mL	25 tests (125 uL)	P015-APC-125	
			100 tests (500 uL)	P015-APC-500	
Purified (BSA-free)	N/A	2 mg/mL	1 mg	P015-1MG	
			5 mg	P015-5MG	
			-		

This datasheet was generated on November 4, 2025 at 03:57:36 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd9-recombinant-monoclonal-mouse-antibody-23109-biotium-choice/">https://biotium.com/product/cd9-recombinant-monoclonal-mouse-antibody-23109-biotium-choice/</a>