

CD9 Monoclonal Mouse Antibody (HI9a) - Biotium Choice

CD9 (clone HI9a) is a validated mouse monoclonal antibody that recognizes the human tetraspanin protein CD9. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

CD9 (clone HI9a) is a mouse monoclonal antibody that recognizes the human tetraspanin protein CD9. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. This antibody has been validated in flow cytometry, in immunofluorescence microscopy, and in western blot. The antibody is available conjugated to R-PE or Astral™ tandem dyes. They are supplied in PBS, 0.1% BSA, 0.05% azide. 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 [Isotype Control, Monoclonal Mouse IgG1 Kappa \(IGG1/1331\)](#) for this antibody. [View our full catalog of Biotium Choice antibodies](#), optimized and validated for the best results in flow cytometry. **Exosome Stains and Antibodies** Biotium also offers [ExoBrite™ CD9 Flow Antibody](#) and [ExoBrite™ 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](#).

Product attributes	
Antibody number	#P003
SwissProt	P21926
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Clonality	Monoclonal
Host species	Mouse
Clone	HI9a
Isotype	IgG1, kappa
Antibody reactivity (react)	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); p25
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	25 kDa
Antibody target cellular localization	Exosomes/EVs, Plasma membrane
Cell/tissue expression	Exosomes, Platelets, Basophils, Eosinophils, Epithelial cells, Lymphocytes
Verified antibody applications	Flow (surface) (verified), WB (verified), IF (verified)
Positive control	Human peripheral blood platelets, BT-20, MCF-7, HeLa, or ZR-75 cells. Cervix or esophagus tissue.
Antibody application notes	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
Antibody research areas	Exosomes/EVs
Antibody/conjugate formulation	Purified: 0.5 mg/mL in PBS/0.05% azide, Conjugates: PBS/0.1% BSA/0.05% 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
Volume per assay	5 uL/test
Regulatory status	For research use only (RUO)

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Availability
CF@405M	416/452 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-405M-125 P003S-405M-500	Discontinued, see replacement P015-405M
CF@488A	490/516 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-488A-125 P003S-488A-500	Discontinued, see replacement P015-488A
CF@568	562/584 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-568-125 P003S-568-500	Discontinued, see replacement P015-568
RPE-Astral™616	496, 546, 566/617 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-R616-125 P003S-R616-500	Available
CF@647Plus	652/668 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-647P-125 P003S-647P-500	Discontinued, see replacement P015-647P
CF@700	696/721 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-700-125 P003S-700-500	Discontinued, see replacement P015-700
APC-Astral™813	633, 638/813 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-A813-125 P003S-A813-500	Available
R-PE	496, 546, 565/576 nm	50 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P003S-RPE-125 P003S-RPE-500	Available
Purified (BSA-free)	N/A	0.5 mg/mL	200 uL (100 ug)	P003-200uL	Discontinued, see replacement P015