

## CD81 Monoclonal Mouse Antibody (5A6)

CD81 (clone 5A6) is a validated mouse monoclonal antibody that recognizes the human tetraspanin protein CD81.



### Product Description

CD81 (clone 5A6) is a mouse monoclonal antibody that recognizes the human tetraspanin protein CD81. This antibody has been validated in flow cytometry, in immunofluorescence microscopy, and in western blot. It is supplied conjugated to R-Phycoerythrin (R-PE or PE) at 0.1 mg/mL with 0.1% BSA, or purified at 0.5 mg/mL, BSA-free. CD81 has a very broad cellular distribution, being expressed on T- and B-lymphocytes, NK cells, thymocytes, eosinophils, fibroblasts, epithelial and endothelial cells. Neutrophils, erythrocytes and platelets are negative, while monocytes are variably positive. CD81 is a member of a family of tetraspanin transmembrane proteins, including CD9, CD37, CD53, CD63, and CD82. It associates with CD19, CD21, Leu 13, and integrins on cell membrane and is involved in signal transduction in B lymphocyte development and cell adhesion. CD81 also acts as a receptor for the envelope protein E2 of chronic hepatitis C virus. Antibodies to CD81 have anti-proliferative effects on different lymphoid cell lines, particularly those derived from large cell lymphomas. 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. **Exosome Stains and Antibodies** Biotium also offers [ExoBrite™ CD81 Flow Antibody](#) and [ExoBrite™ CD81 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	#P005
Storage Conditions	Store at 2 to 8 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody/conjugate formulation	Purified: 0.5 mg/mL in PBS/0.05% azide, Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide
Antibody research areas	Exosomes/EVs
Antibody application notes	Flow Cytometry: 0.5-1 ug/mL (or, 0.05-0.1 ug/million cells), Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL
Positive control	Ramos, Daudi, Raji, U266, HUT-78, HUT-102, Jurkat, HL-60, KG1a, A293, A431, MG63, MCF-7, and human lymphocytes. Tonsil or liver tissue.
Verified antibody applications	Flow (surface) (verified), IF (verified), WB (verified)
Cell/tissue expression	Exosomes, Hepatocytes, Lymphocytes, Monocytes/macrophages
Antibody target cellular localization	Exosomes/EVs, Plasma membrane
Molecular weight	26 kDa
SwissProt	P60033
Unigene	54457
Entrez gene ID	975
Human gene symbol	CD81
Species reactivity	Baboon, Cynomolgus monkey, Human, Non-human primates
Synonyms	CD81; Target of the antiproliferative antibody 1 (TAPA-1); Tetraspanin-28; Tspan-28
Antibody reactivity (target) isotype	CD81 IgG1, kappa
Clone	5A6
Host species	Mouse
Clonality	Monoclonal
Antibody type	Primary
Regulatory status	For research use only (RUO)

## CD81 Monoclonal Mouse Antibody (5A6)

Conjugation	Ex/Em	Size	Catalog No.
<a href="#">R-PE</a>	496, 546, 565/576 nm	25 tests 100 tests	<a href="#">P005S-RPE-125</a> <a href="#">P005S-RPE-500</a>
<a href="#">Purified (BSA-free, 0.5 mg/mL)</a>	N/A	200 uL (100 ug)	<a href="#">P005-200uL</a>