

## CD81 Recombinant Monoclonal Mouse Antibody (2304.81) - Biotium Choice

CD81 (clone 2304.81) is a validated recombinant mouse monoclonal antibody that recognizes the human tetraspanin protein CD81. The antibody belongs to the Biotium Choice list of select antibodies that have been formulated and validated in-house for optimal performance.



### Product Description

CD81 (clone 2304.81) is a recombinant mouse monoclonal antibody that recognizes the human tetraspanin protein CD81. This antibody belongs to the Biotium Choice list of select antibodies that have been formulated and validated in-house for optimal performance. This antibody has been validated in flow cytometry and immunofluorescence microscopy. The antibody is available unlabeled or conjugated to CF® Dyes and Astral™ tandem dyes. 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. [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™ 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	#P014
SwissProt	P60033
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Clonality	Recombinant Monoclonal
Host species	Mouse
Clone	2304.81
Isotype	IgG1, kappa
Antibody reactivity (react)	CD81
Synonyms	CD81; Target of the antiproliferative antibody 1 (TAPA-1); Tetraspanin-28; Tspan-28
Species reactivity	Baboon, Cynomolgus monkey, Human, Mouse (low reactivity), Non-human primates
Human gene symbol	CD81
Entrez gene ID	975
Unigene	54457
Molecular weight	26 kDa
Antibody target cellular localization	Exosomes/EVs, Plasma membrane
Cell/tissue expression	Exosomes, Hepatocytes, Lymphocytes, Monocytes/macrophages
Verified antibody applications	Flow (surface) (verified), IF (verified)
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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.
Antibody application notes	Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells
Antibody research areas	Exosomes/EVs
Antibody/conjugate formulation	Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide
Volume per assay	5 uL/test
Storage Conditions	Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C
Regulatory status	For research use only (RUO)

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF@405M</a>	416/452 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-405M-125</a> <a href="#">P014-405M-500</a>	<a href="#">CF@405M Features</a>
<a href="#">CF@488A</a>	490/516 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-488A-125</a> <a href="#">P014-488A-500</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@ 568</a>	562/584 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-568-125</a> <a href="#">P014-568-500</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@ 583R</a>	585/609 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-583R-125</a> <a href="#">P014-583R-500</a>	<a href="#">CF@583R Features</a>
<a href="#">RPE-Astral™616</a>	496, 546, 566/617 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-R616-125</a> <a href="#">P014-R616-500</a>	
<a href="#">CF@ 647Plus</a>	652/668 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-647P-125</a> <a href="#">P014-647P-500</a>	<a href="#">CF@647Plus Features</a>
<a href="#">CF@700</a>	696/721 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-700-125</a> <a href="#">P014-700-500</a>	<a href="#">CF@700 Features</a>
<a href="#">R-PE</a>	496, 546, 565/576 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-RPE-125</a> <a href="#">P014-RPE-500</a>	
<a href="#">APC</a>	651/660 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	<a href="#">P014-APC-125</a> <a href="#">P014-APC-500</a>	
<a href="#">Purified (BSA-free)</a>	N/A	5 mg/mL	1 mg 5 mg	<a href="#">P014-1MG</a> <a href="#">P014-5MG</a>	