

## 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.

### 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.](#)

### 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™ 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 [Astral Leap™ tandem dyes](#) for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)

### Product attributes

Antibody number	#P014
SwissProt	P60033
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice, Primary
Clonality	Monoclonal, Recombinant Monoclonal
Host species	Mouse
Clone	2304.81
Isotype	IgG1, kappa
Antibody reactivity (target)	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 of antigen	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/0.05% azide
Volume per assay	5 uL/test
Storage Conditions	Store at 2 to 8 °C
Regulatory status	For research use only (RUO)
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF@405M</a> 100 tests (500 uL)	416/452 nm <a href="#">P014-405M-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-405M-125</a>	<a href="#">CF@405M Features</a>
<a href="#">CF@488A</a> 100 tests (500 uL)	490/516 nm <a href="#">P014-488A-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-488A-125</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@ 568</a> 100 tests (500 uL)	562/584 nm <a href="#">P014-568-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-568-125</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@ 583R</a> 100 tests (500 uL)	585/609 nm <a href="#">P014-583R-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-583R-125</a>	<a href="#">CF@583R Features</a>
<a href="#">RPE-Astral™616</a> 100 tests (500 uL)	496, 546, 566/617 nm <a href="#">P014-R616-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-R616-125</a>	<a href="#">RPE-Astral™616 Features</a>
<a href="#">CF@ 647Plus</a> 100 tests (500 uL)	652/668 nm <a href="#">P014-647P-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-647P-125</a>	<a href="#">CF@647Plus Features</a>
<a href="#">CF@700</a> 100 tests (500 uL)	696/721 nm <a href="#">P014-700-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-700-125</a>	<a href="#">CF@700 Features</a>
<a href="#">RPE-Astral™775</a> 100 tests (500 uL)	496, 546, 565/774 nm <a href="#">P014-R775-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-R775-125</a>	<a href="#">RPE-Astral™775 Features</a>
<a href="#">R-PE</a> 100 tests (500 uL)	496, 546, 565/576 nm <a href="#">P014-RPE-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-RPE-125</a>	
<a href="#">APC</a> 100 tests (500 uL)	651/660 nm <a href="#">P014-APC-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P014-APC-125</a>	
<a href="#">Purified (BSA-free)</a> 5 mg (1 mL)	N/A <a href="#">P014-5MG</a>	5 mg/mL	1 mg (200 uL)	<a href="#">P014-1MG</a>	

This datasheet was generated on June 7, 2026 at 07:16:21 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cd81-recombinant-monoclonal-mouse-antibody-230481-biotium-choice/>