



Glowing products for science

CD56 (NCAM1) Recombinant Monoclonal Mouse Antibody (r5.1H11) - Biotium Choice



CD56 (clone r5.1H11) is a recombinant mouse monoclonal antibody that recognizes transmembrane glycoprotein CD56 (NCAM, neural cell adhesion molecule). The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.

Product Description

CD56 (clone r5.1H11) is a recombinant mouse monoclonal antibody that recognizes transmembrane glycoprotein CD56 (NCAM, neural cell adhesion molecule). This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to CF® Dyes, R-PE, PerCP, and APC. They are supplied in PBS, 0.1% BSA, 0.05% azide.

CD56 exists in isoforms ranging from 120 to 180 kDa and is primarily found on T cells and NK cells. It plays a crucial role in both homophilic and heterophilic adhesion by binding to itself or heparin sulfate. CD56 is also involved in key cellular processes such as migration, axonal growth, pathfinding, and synaptic plasticity. It is widely expressed in neuroectodermal-derived cell lines, tissues, and tumors, including retinoblastoma, medulloblastoma, astrocytomas, and neuroblastoma. Additionally, CD56 serves as a highly sensitive neuroendocrine marker, commonly used for detecting neuroendocrine tumors and ovarian granulosa cell tumors. Dysfunction of CD56 has been linked to diseases such as rabies and blastic plasmacytoid dendritic cell neoplasms.

As an isotype control, we recommend [Isotype Control, Monoclonal Mouse IgG1 Kappa \(IGG1/1331\)](#) for this antibody.

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](#)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	P026
Antibody reactivity (target)	CD56
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Biotium Choice Primary, Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	r5.1H11
Isotype	IgG1, kappa
Molecular weight	N/A
Synonyms	NCAM1, NKH1
Human gene symbol	CD45
Entrez gene ID	4684
SwissProt	P13591
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (surface) (verified)
Species reactivity	Human
Antibody application notes	Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL
Positive control	Human peripheral blood mononuclear cells (PBMCs)
Shipping condition	Room temperature
Storage Conditions	Note: store unconjugated formats at -10 °C to -35 °C. Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Volume per assay	5 uL/test
Antibody/conjugate formulation	Conjugates: PBS/0.1% rBSA/0.05% azide, Purified: PBS, no BSA, no azide
Antibody research areas	Cancer, Cell adhesion, Neuroscience
Product origin	Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line; recombinant BSA produced in Chinese hamster ovary cells.
Cell/tissue expression	Skeletal Muscle, Neurons, NK cells
Tumor expression	Neuroectodermal cancer

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF@405M	416/452 nm	100 ug/mL	25 tests (125 uL)	P026-405M-125	CF@405M Features
100 tests (500 uL)	P026-405M-500				
CF@488A	490/516 nm	200 ug/mL	25 tests (125 uL)	P026-488A-125	CF@488A Features
100 tests (500 uL)	P026-488A-500				
CF@ 575	575/601 nm	100 ug/mL	25 tests (125 uL)	P026-575-125	
100 tests (500 uL)	P026-575-500				
CF@647Plus	652/668 nm	25 ug/mL	25 tests (125 uL)	P026-647P-125	CF@647 Features
100 tests (500 uL)	P026-647P-500				
CF@660C	667/685 nm	100 ug/mL	25 tests (125 uL)	P026-660C-125	CF@660C Features
100 tests (500 uL)	P026-660C-500				
CF@700	696/721 nm	100 ug/mL	25 tests (125 uL)	P026-700-125	CF@700 Features
100 tests (500 uL)	P026-700-500				
CF@820	822/835 nm	20 ug/mL	25 tests (125 uL)	P026-820-125	CF@820 Features
100 tests (500 uL)	P026-820-500				
R-PE	496, 546, 565/576 nm	20 ug/mL	25 tests (20 uL)	P026-RPE-125	
100 tests (20 uL)	P026-RPE-500				
PerCP	482/675 nm	100 ug/mL	25 tests (125 uL)	P026-PCP-125	
100 tests (500 uL)	P026-PCP-500				
APC	651/660 nm	80 ug/mL	25 tests (80 uL)	P026-APC-125	
100 tests (80 uL)	P026-APC-500				
Purified (BSA-free, 5 mg/mL)	N/A	5 mg/mL	200 uL (1 mg)	P026-1MG	
1 mL (5 mg)	P026-5MG				

This datasheet was generated on January 13, 2026 at 01:20:46 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd56-recombinant-monoclonal-mouse-antibody-r5-1h11-biotium-choice/>