



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

# CD38 Monoclonal Mouse Antibody (FS02)

## Product Description

This antibody reacts with a type II membrane glycopeptide with a transmembrane sequence near the NH2-terminus. CD38 is a type II transmembrane glycoprotein that is present on early B- and T-cell lineages and activated B- and T-cells but is absent from most mature resting peripheral lymphocytes. CD38 is also found on thymocytes, pre-B cells, germinal center B-cells, mitogen-activated T-cells, monocytes and Ig-secreting plasma cells. CD38 is expressed on CD34 cells. The CD34 CD38-population of hematopoietic stems cells defines the most pluripotent cells (e.g. blast colony forming cells). **Catalog number key for antibody number 0171, Anti-CD38 (FS02)**

### Product attributes

Antibody number	#0171
Antibody reactivity (target)	CD38
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FS02
Isotype	IgG1, kappa
Molecular weight	~45 kDa (Glycoprotein); 35 kDa (protein core)
Synonyms	Acute Lymphoblastic Leukemia Cells Antigen; ADP Ribosyl Cyclase 1; cADP-ribose Hydrolase 1; CD38H; NAD(+) Nucleosidase; NIM-R5 Antigen; p45; T10
Human gene symbol	CD38
Entrez gene ID	952
SwissProt	P28907
Unigene	479214
Immunogen	Human CD38
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	CCRF-CEM cells, Tonsil, Spleen or Skin
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	B-cells, T-cells
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.
Antibody research areas	Immunology

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on July 12, 2025 at 02:39:40 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cd38-monoclonal-mouse-antibody-fs02/>