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

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	cute Lymphoblastic Leukemia Cells Antigen; ADP Ribosyl tyclase 1; cADP-ribose Hydrolase 1; CD38H; NAD(+) lucleosidase; NIM-R5 Antigen; p45; T10			
Human gene symbol	CD38			
Entrez gene ID	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				
Regulatory status Antibody/conjugate formulation	Note: store BSA-free antibodies at -10 to -35 °C			
Antibody/conjugate	Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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			
Antibody/conjugate formulation	Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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 Guaranteed for at least 24 months from date of receipt when			
Antibody/conjugate formulation Shelf life	Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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, Purified: 0.2 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when stored as recommended			

Call us: 800-304-5357 Email: btinfo@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
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 November 6, 2025 at 09:52:55 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd38-monoclonal-mouse-antibody-fs02/