CD38 Monoclonal Mouse Antibody (AT1)

Product Description

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).

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0430, Anti-CD38 (AT1)

Droduct attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#0430			
Antibody reactivity (target)	CD38			
Antibody type	rimary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	AT1			
Isotype	IgG1, kappa			
Molecular weight	~45 kDa (Glycoprotein); 35 kDa (protein core)			
Synonyms	Cell differentiation antigen 37; GP52 40; Tetraspanin-26; Tspan-26; TSPAN26			
Human gene symbol	CD38			
Entrez gene ID	951			
SwissProt	P28907			
Unigene	479214			
Immunogen	human T cell line CCRF-CEM			
Antibody target cellular localization	Plasma membrane			
Expected antibody applications	Flow, surface (published for clone), IHC (FFPE) (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)			
Species reactivity	Human			
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			
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended			
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			
Antibody research areas	Immunology			
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.			
Cell/tissue expression	B-cells, T-cells			

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- 1. Diabetologia (2002) 45:1667-1677. (IHC, FFPE; IF)
- 2. BMC Immunol (2004) 5: 21. (Flow, IF, WB)
- 3. J Cellular Biochem (2008) 103:1294-1308. (IP, WB, Flow, IF)
- 4.

This datasheet was generated on November 30, 2025 at 12:41:47 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd38-at1/