

CD11a Monoclonal Mouse Antibody (DF1524)

Product Description

Recognizes a protein of 180 kDa, identified as CD11a (Leucocyte Workshop IV; Code 1524). CD11a complex with the 2 subunit of the integrin family, CD18, to form the cell surface heterodimer, LFA-1 or CD11a /C18 (aLbL). LFA-1 is expressed on all leukocytes including lymphocytes, monocytes, and granulocytes. It is involved in leukocyte adhesion to its ligands including intercellular adhesion molecule-1 (ICAM-1 or CD54), ICAM-2 (CD102), ICAM-3 (CD50) and Telencephalin (TLN) and play a role in most immune/inflammatory responses. This MAb potently blocks LFA-1 dependent homotypic cell aggregation.

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 0312, Anti-CD11a (DF1524)

Product attributes

Antibody number	#0312
Antibody reactivity (target)	CD11a
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	DF1524
Isotype	IgG2b, kappa
Molecular weight	180 kDa
Synonyms	Antigen CD11A (p180), lymphocyte function associated antigen 1, alpha polypeptide; Integrin alpha-L; ITGAL; Leukocyte Adhesion Glycoprotein LFA1 Alpha Chain; LFA 1 alpha (LFA1A); Ly15; Ly21; p180
Human gene symbol	ITGAL
Entrez gene ID	3683
SwissProt	P20701
Unigene	174103
Immunogen	Human CD11a protein
Antibody target cellular localization	Plasma membrane
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL
Positive control	JY, HUT-78, MOLT-4, HL60, U937, and human leukocytes. Human lymph nodes and tonsils.
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	Hematopoietic 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. J Neuropathol Exp Neurol (2000) 59(1): 18-28. (functional studies)
2. FASEB J (2003) 17(11), 1-21. (Flow)
3. Blood (2006) 107(8): 3212-3220. (IF)

This datasheet was generated on December 12, 2025 at 03:15:35 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd11a-df1524/>