

CD11a Monoclonal Mouse Antibody (Cris-3)



Product Description

Recognizes a protein of 180 kDa, identified as CD11a (Leucocyte Workshop II; Code N202). 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.

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 0151, Anti-CD11a (Cris-3)

Call us: 800-304-5357

Product attributes

Antibody number	#0151			
Antibody reactivity (target)	CD11a			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	Cris-3			
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	Stimulated human leukocytes			
Antibody target cellular localization	Plasma membrane			
Verified antibody applications	Flow (verified)			
Species reactivity	Human			
Expected antibody applications	Functional studies (published for clone)			
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				
Positive control Shipping condition	Cytometry 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL60, U937, and human leukocytes.			
	Cytometry 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL60, U937, and human leukocytes. Human lymph nodes and tonsils.			
Shipping condition	Cytometry 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL60, U937, and human leukocytes. Human lymph nodes and tonsils. Room temperature Store at 2 to 8 ° C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	Cytometrý 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL80, U937, and human leukocytes. Human lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when			
Shipping condition Storage Conditions Shelf life	Cytometrý 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL80, U937, and human leukocytes. Human lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	Cytometrý 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL60, U937, and human leukocytes. Human lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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; 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.2 mg/mL in PBS/0.05% azide, Purified; 0			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Cytometrý 0.5-1 ug/million cells/0.1 mL JY, HUT-78, MOLT-4, HL80, U937, and human leukocytes. Human lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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			

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.

- Knapp W. et al. Leukocyte Typing IV, Oxford Unioversity Press, Oxford, 1989.
 Reinherz EL et al. eds. Leukocyte Typing II, Oxford University Press, Oxford, 1985.
 J Exp Med (1990) 172: 335-345. (Flow, surface)
 Am J Physiol Cell Physiol (1998) 274(6): C1634-C1644. (receptor blockade).

This datasheet was generated on November 30, 2025 at 03:39:16 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd11a-cris-3/