CD11c Monoclonal Mouse Antibody (ITGAX/1284)



Product Description

This antibody recognizes a protein of 145 kDa, identified as CD11c. CD11c (ITGAX), a member of the leukointegrin family, shares the same beta subunit with other members of the leukocyte adhesion molecule family, which includes CD11a (LFA-1), CD11b (MAC-1) and CD11d (ITGAD), but has a unique alpha chain. CD11c has been shown to play a role in phagocytosis, cell migration, and cytokine production by monocytes/macrophages as well as induction of T-cell proliferation by Langerhans cells. CD11c is expressed prominently on the plasma membranes of monocytes, tissue macrophages, NK cells, and most dendritic cells (DCs). A lower level of expression is also observed on neutrophils as a result of its high level of expression on most DCs. An antibody to CD11c may aid in identification of lesions with histiocytic origin. It may also been used as a marker for hairy cell leukemia in paraffin-embedded tissues.

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 1284, Anti-CD11c (ITGAX/1284)

Product attributes

Tumor expression

Call us: 800-304-5357

Product attributes				
Antibody number	#1284			
Antibody reactivity (target)	CD11c			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	ITGAX/1284			
Isotype	lgG2b			
Molecular weight	145 kDa			
Synonyms	CD11 antigen-like family member C, Complement component 3 receptor 4 subunit, Integrin alpha-X, integrin, alpha X (antigen CD11C (p150), alpha polypeptide), Leu M5 alpha subunit, Leukocyte adhesion glycoprotein p150 95 alpha chain, Myeloid membrane antigen alpha subunit, p150/95			
Human gene symbol	ITGAX			
Entrez gene ID	3687			
SwissProt	P20702			
Unigene	248472			
Immunogen	Recombinant human ITGAX protein fragment (aa 637-827) (exact sequence is proprietary)			
Antibody target cellular localization	Plasma membrane			
Verified antibody applications	IHC (FFPE) (verified)			
Species reactivity	Human			
Antibody application notes	Higher concentration may be required for direct detection us primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology (formalin) 1-2 ug/mL, Staining of formalin-fitssues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution a specific application should be determined by user			
Positive control	Human dendritic cells. 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 PBS without azide			
Antibody research areas	Cell adhesion, Immunology			
Product origin Product may contain either bovine serum albumin (BS bovine serum (Bos taurus), or recombinant BSA produ Chinese hamster ovary cells. Inquire for the specific lo				

Leukemia/lymphoma

Email: btinfo@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
Antibody # prenx	Conjugation	EX/EIII (IIIII)	Laser inte	Detection channel	Dye i catales
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 4, 2025 at 11:03:51 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mouse-monoclonal-anti-cd11c-itgax1284/