## CD11b Monoclonal Mouse Antibody (ITGAM/3339)



## **Product Description**

CD11b is a cell adhesion molecule that acts as a receptor for cell surface ligands such as intracellular adhesion molecules (ICAMs) or soluble ligands. Integrins are heterodimeric proteins that contain an a chain and b chain. Integrin M2 is important in the adherence of neutrophils and monocytes to stimulated endothelium, and also in the phagocytosis of complement coated particles. The protein CD11b has been implicated in the various adhesion-related interactions of cells such as monocytes, macrophages, natural killer (NK) cells, and granulocytes. It is part of a heterodimer that consists of CD11b and CD18. It also modulates the uptake of complement-coated particles within the cell. It is commonly used as a microglial marker in tissues derived from the nervous system.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3339, Anti-CD11b|MAC-1 (ITGAM/3339)

## Product attributes

Call us: 800-304-5357

Antibody number	#3339		
Antibody reactivity (target)	CD11b, MAC-1		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	ITGAM/3339		
Isotype	IgG1, kappa		
Molecular weight	95 kDa & 170 kDa		
Synonyms	CD18; CD49d; Cell surface glycoprotein MAC-1 subunit alpha; Complement Component Receptor 3 Alpha; CR3 Alpha Chain (CR3A); Integrin alpha-M; Integrin beta 2 alpha subunit; Leukocyte adhesion receptor MO1; Ly-40; MAC1A; Macrophage antigen alpha polypeptic; MO1A; Neutrophil adherence receptor alpha M subunit		
Human gene symbol	ITGAM		
Entrez gene ID	3684		
SwissProt	P11215		
Unigene	172631		
Immunogen	Recombinant fragment (around aa941-1074) of human ITGAM protein (exact sequence is proprietary)		
	protein (exact sequence is proprietary)		
Verified antibody applications	IHC (FFPE) (verified)		
applications Antibody target cellular	IHC (FFPE) (verified)		
applications  Antibody target cellular localization	IHC (FFPE) (verified) Plasma membrane		
applications  Antibody target cellular localization  Species reactivity	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil		
applications Antibody target cellular localization Species reactivity Positive control	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).		
applications Antibody target cellular localization Species reactivity Positive control Shipping condition	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, 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: 0.2 mg/mL in PBS/0.05% azide,		
applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% aride, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 1 mg/mL in PBS without azide		
applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Validated in protein array	IHC (FFPE) (verified)  Plasma membrane  Human  Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, 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, Purffled, BSA-free: 1 mg/mL in PBS without azide  Monospecific  Guaranteed for at least 24 months from date of receipt when		

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 9, 2025 at 11:44:51 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd11b-monoclonal-mouse-antibody-itgam-3339/">https://biotium.com/product/cd11b-monoclonal-mouse-antibody-itgam-3339/</a>