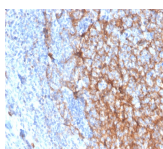


CD11b Monoclonal Mouse Antibody (ITGAM/3337)



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 α chain and β chain. Integrin α M combines with the Integrin β 2 to form a leukocyte-specific integrin referred to as macrophage receptor 1 (Mac-1), or inactivated-C3b (iC3b) receptor 3 (CR3). Integrin α M/ β 2 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 order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3337, Anti-CD11b|MAC-1 (ITGAM/3337)

Product attributes

Antibody number	#3337
Antibody reactivity (target)	CD11b, MAC-1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ITGAM/3337
Isotype	IgG2b, kappa
Molecular weight of antigen	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 polypeptide; 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)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	Monocytes & granulocytes. Lymph nodes, spleen and tonsil (IHC).
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
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
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Microglia, Monocytes/macrophages, Neutrophils

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	775/760	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 June 4, 2026 at 09:27:17 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd11b-monoclonal-mouse-antibody-itgam-3337/>