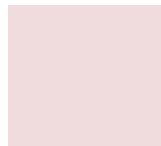


CD11b / MAC-1 Monoclonal Mouse Antibody (ITGAM/271)



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 αM combines with the Integrin β2 to form a leukocyte-specific integrin referred to as macrophage receptor 1 (Mac-1), or inactivated-C3b (iC3b) receptor 3 (CR3). Integrin αM/β2 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 0271, Anti-CD11b (ITGAM/271)**

Product attributes

Antibody number	#0271
Antibody reactivity (target)	CD11b, MAC-1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ITGAM/271
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 polypeptide; MO1A; Neutrophil adherence receptor alpha M subunit
Human gene symbol	ITGAM
Entrez gene ID	3684 (Human)
SwissProt	P11215 (Human)
Unigene	172631 (Human)
Immunogen	Recombinant human ITGAM protein
Antibody target cellular localization	Membrane/cell surface
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Human monocytes & granulocytes. 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 buffer 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, Neuroscience
Cell/tissue expression	Microglia, Monocytes/macrophages, Neutrophils

Catalog Number Key for Primary Antibody Conjugates

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	CF@405S Features
BNC05	CF@405M	408/452	405	DAPI (microscopy), Pacific Blue™	CF@405M Features
BNC06	CF@405L	395/545	405	Pacific Orange™	CF@405L Features
BNC88	CF@488A	490/515	488	GFP, FITC	CF@488A Features
BNC14	CF@514	516/548	488	GFP, FITC	CF@514 Features
BNC43	CF@543	541/560	532	RFP, TRITC	CF@543 Features
BNC55	CF@555	555/565	532, 561	RFP, TRITC	CF@555 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
BNC60	CF@660C	667/685	633-640	Cy@5	CF@660C Features
BNC61	CF@660R	663/682	633-640	Cy@5	CF@660R Features
BNC80	CF@680	681/698	633-640	Cy@5.5, IR@700	CF@680 Features
BNC81	CF@680R	680/701	633-640	Cy@5.5, IR@700	CF@680R Features
BNC00	CF@700	695/720	633-640	AF700	CF@700 Features
BNC70	CF@770	770/797	785, 808	IR@800	CF@770 Features
BNCR	R-PE (PE)	496, 546, 565/578	488-561	PE	
BNCA	APC	650/660	594-633	APC	
BNCP	PerCP	482/677	488-561	PerCP, PE-Cy@5	
BNCB	Biotin	N/A	N/A	N/A	
BNCAP	Alkaline Phosphatase	N/A	N/A	N/A	
BNCH	HRP	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 GE Healthcare; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.