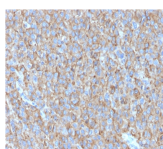


CD63 Monoclonal Mouse Antibody (LAMP3/2790)



Product Description

This MAbs recognizes protein of 26 kDa-60 kDa, which is identified as CD63. Its epitope is different from that of MAbs LAMP3/529. The tetraspanins are integral membrane proteins expressed on cell surface and granular membranes of hematopoietic cells and are components of multi-molecular complexes with specific integrins. The tetraspanin CD63 is a lysosomal membrane glycoprotein that translocates to the plasma membrane after platelet activation. CD63 is expressed on activated platelets, monocytes and macrophages, and is weakly expressed on granulocytes, T cell and B cells. It is located on the basophilic granule membranes and on the plasma membranes of lymphocytes and granulocytes. CD63 is a member of the TM4 superfamily of leukocyte glycoproteins that includes CD9, CD37 and CD53, which contain four transmembrane regions. CD63 may play a role in phagocytic and intracellular lysosome-phagosome fusion events. CD63 deficiency is associated with Hermansky-Pudlak syndrome and is strongly expressed during the early stages of melanoma progression.

Also view our [CD63 Recombinant Monoclonal Mouse Biotium Choice Antibody \(2304.63\)](#). This antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance. The antibody has been validated in flow cytometry and immunofluorescence microscopy.

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 2790, Anti-CD63 (LAMP3/2790)

Call us : [800-304-5357](tel:800-304-5357) Email: techsupport@biotium.com

Product attributes

Antibody number	#2790
Antibody reactivity (target)	CD63
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	LAMP3/2790
Isotype	IgG2a, kappa
Molecular weight	26 kDa (core protein); 30-60 kDa (glycosylated)
Synonyms	gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen ME491; MLA1; NGA; Ocular melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30
Human gene symbol	CD63
Entrez gene ID	967
SwissProt	P08962
Unigene	445570
Immunogen	Recombinant fragment (around aa 100-197) of human CD63 protein
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies
Species reactivity	Human
Positive control	U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue.
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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages
Tumor expression	Melanoma

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 September 6, 2025 at 05:01:40 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cd63-monoclonal-mouse-antibody-lamp3-2790/>