CD50 Monoclonal Mouse Antibody (ICAM3/1019)



Product Description

Recognizes an N-glycosylated glycoprotein of 120 kDa with intra-chain disulfide bonds, identified as CD50 or ICAM-3. CD50 is the major ligand for LFA-1 (CD11a/CD18) and may have signaling role to increase adhesion. It is expressed on thymocytes and T lymphocytes and is resistant to treatment with phosphatidylinositol (PI) phospholipase C. This MAb is excellent for staining of formalin/paraffin 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 1019, Anti-CD50 (ICAM3/1019)

Product attributes

Call us: 800-304-5357

Molecular weight 110-16	ilonal V/1019 kappa		
Antibody type Primar Host species Mouse Clonality Monoc Clone ICAM: Isotype IgG2b Molecular weight 110-16	ilonal V/1019 kappa		
Host species Mouse Clonality Monoc Clone ICAM Isotype IgG2b Molecular weight 110-16	ilonal V/1019 kappa		
Clonality Monoc Clone ICAM Isotype IgG2b Molecular weight 110-16	lonal V/1019 kappa		
Clone ICAM3 Isotype IgG2b Molecular weight 110-16	v/1019 kappa		
Isotype IgG2b Molecular weight 110-16	kappa		
Molecular weight 110-16	**		
	O LDa		
Synonyms ICAME			
	R; Intercellular adhesion molecule 3 (ICAM3)		
Human gene symbol ICAM3	ene symbol ICAM3		
Entrez gene ID 3385	ez gene ID 3385		
SwissProt P3294	rot P32942		
Unigene 35456	3		
Immunogen Recon	abinant human ICAM3 protein		
Verified antibody IHC (F applications			
Antibody target cellular Plasm localization	a membrane		
Species reactivity Human			
primar secon Immur formal citrate 20 mir Weste	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user		
Positive control HL-60 or THP-1 cells. Lymph node and tonsil			
Shipping condition Room	temperature		
	e Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
	life Guaranteed for at least 24 months from date of receipt when stored as recommended		
stored			
	search use only (RUO)		
Regulatory status For res Antibody/conjugate formulation conjug in PBS	search use only (RUO) jates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP ates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL //0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in ithout azide		
Regulatory status For res Antibody/conjugate formulation conjug in PBS	pates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP ates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL /0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in ithout azide		
Regulatory status	pates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP ates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL /0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in ithout azide		
Verified antibody applications Antibody target cellular localization Species reactivity Human Antibody application notes Higher	FPE) (verified), WB (verified) a membrane concentration may be required for direct detection using		

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 4, 2025 at 02:00:57 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd50-icam-3-icam31019/