CD54 / ICAM-1 Monoclonal Mouse Antibody (15.2)



Product Description

Recognizes an 85-115 kDa protein (variation with cell type), identified as intercellular adhesion molecule (ICAM-1). It has 7 potential N-linked glycosylation sites. ICAM-1 is a single chain glycoprotein of Ig supergene family, present on unstimulated endothelial cells (EC) and on a variety of other cell types including activated fibroblasts, EC, macrophages, and lymphocytes. ICAM-1 mediates cell adhesion by binding to integrins CD11a/CD18 (leukocyte adhesion molecule, LFA-1) and to CD11b/CD18 (Mac-1). This interaction enhances antigen-specific T-cell activation. ICAM-1 also binds to CD43 and to Plasmodium falciparum infected RBCs. ICAM-1 may also be related to progression and metastasis of tumors.

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 2853, Anti-CD54|ICAM-1 (15.2)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#2853		
Antibody reactivity (target)	CD54, ICAM-1		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	15.2		
Isotype	IgG1, kappa		
Molecular weight	85-115 kDa		
Synonyms	CD54; Cell surface glycoprotein P3.58; Human rhinovirus receptor; ICAM-1; Intercellular adhesion molecule 1; Ly 47; Major group rhinovirus receptor; MALA2; MyD10; Surface antigen of activated B cells		
Human gene symbol	ICAM1		
Entrez gene ID	3383		
SwissProt	P05362		
Unigene	643447		
Immunogen	Human monocytes.		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human, Pig		
Expected antibody applications	IHC (frozen) (published for clone), Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone)		
Positive control	Raji cells, MOLT 4, HeLa or SW480 cells. Human tonsil.		
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		
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	Endothelial cells, Lymphocytes, Monocytes/macrophages		
Antibody research areas	Immunology		

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- 1. J Leukocyte Biol (1998) 64: 519. (Flow, surface)
- 2. Inv Ophthalmol Visual Sci (2003) 44: 1888-1896. (IHC, frozen)
- 3. Blood (1999) 93(11): 3723-3735. (IF)

J Cell Biol (1992) 116(6): 1527-1535. (functional studies)

This datasheet was generated on November 30, 2025 at 04:23:09 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd54-icam-1-monoclonal-mouse-antibody-15-2/