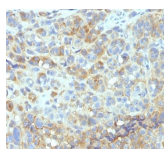


CD54 / ICAM-1 Monoclonal Mouse Antibody (Wehi-CAM-1 or W-CAM-1 or 1H4)



Product Description

Recognizes an 85-115 kDa protein (variation with cell type), identified as intercellular adhesion molecule (ICAM-1) (Workshop IV). 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. W-CAM-1 MAb blocks aggregation of cell lines mediated by the ICAM-1 and blocks homotypic binding of purified populations of activated T- and B-lymphocytes and also aggregation of mixed T- and B-cell blasts. It inhibits T-cell adhesion to normal human endothelial cells. Activation induced by cell-cell contact (mixed lymphocyte reaction, T-cell mediated B-cell activation) is significantly inhibited. This MAb blocks elements of both effector arms of immune system (cytotoxic cell function and Ig production).

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 0988, Anti-CD54 (1H4 or Wehi-CAM-1 or W-CAM-1)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#0988
Antibody reactivity (target)	CD54
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	Wehi-CAM-1 or W-CAM-1 or 1H4
Isotype	IgG2b, kappa
Molecular weight of antigen	85-115 kDa
Synonyms	Intestinal mucin-3A; intestinal mucin-like; MUC3; MUC3A; MUC3B; mucin 3, intestinal; mucin 3A, cell surface associated; mucin 3A, intestinal; mucin 3B, cell surface associated
Human gene symbol	ICAM1
Entrez gene ID	3383
SwissProt	P05362
Unigene	643447
Immunogen	Raji Burkitt lymphoma cells
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	Immunohistochemistry formalin-fixed 2-4 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Functional Studies order Ab without BSA & azide), Optimal dilution for a specific application should be determined by user
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
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 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
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	Endothelial cells, Lymphocytes, Monocytes/macrophages

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 June 15, 2026 at 09:10:46 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd54-1h4-or-wehi-cam-1-or-w-cam-1/>