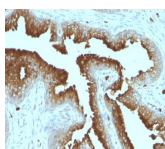


CD53 Monoclonal Mouse Antibody (161-2)



Product Description

Recognizes a protein of 33-55 kDa, identified as CD53 (Workshop VI; Code N-L033). It is expressed on monocytes and macrophages, dendritic cells, osteoblasts and osteoclasts, and on T and B cells from every stage of differentiation but is absent from platelets, red blood cells. CD53 appears to be the marker with widest reactivity as well as the marker with the strictest specificity to hematopoietic cells. CD53 is a type III membrane with both termini in the cytoplasm and two loops in the extracellular environment. This molecule, in common with other members of tetraspan family, is involved in cellular activation as part of a signal transduction complex involving other membrane glycoproteins. CD53 crosslinking induces calcium flux on human monocyte and B cells. Cross-linking of CD53 promotes activation of resting human B-lymphocytes. This MAb recognizes CD53 transfected cells and partially inhibits T-cell proliferation induced by CD3 antibody (clone: UCHT1).

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 0324, Anti-CD53 (161-2)

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

Product attributes

Antibody number	#0324
Antibody reactivity (target)	CD53
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	161-2
Isotype	IgG2a, kappa
Molecular weight	33-55 kDa
Synonyms	CD53 glycoprotein; CD53 tetraspan antigen; Cell surface glycoprotein CD53; Leukocyte surface antigen CD53; Tetraspanin-25 (Tspan-25); TSPAN25
Human gene symbol	CD53
Entrez gene ID	963
SwissProt	P19397
Unigene	443057
Immunogen	Stimulated human leukocytes
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Functional studies (published for clone)
Antibody application notes	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. Does not react with baboon or horse, others not tested, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user
Positive control	Daudi, Raji, IM9, U266, YT, HUT-78, HUT-102, Jurkat, HL-60, THP-1, KG1a, human leukocytes. Human tonsil and lymph node.
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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Hematopoietic cells, Osteoblasts/osteoclasts

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.

- Immunology (1999) 96(1): 35-47. (functional studies)
- Eur J Biochem (2002) 269: 1012-1021. (functional studies)