

# CD48 Monoclonal Mouse Antibody (5-4.8)

## Product Description

Reacts with human CD48, a 45 kDa glycosyl phosphatidyl-inositol (GPI)-anchored cell surface protein. CD48 is strongly expressed on lymphocytes and monocytes and weakly on granulocytes but is absent on platelets, fibroblasts, epithelium and endothelium. CD48 is one of the markers for detecting the defects of GPI anchoring structure on the patients with paroxysmal nocturnal hemoglobinuria (PNH) and serves as a low affinity ligand for CD2.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 1057, Anti-CD48 (5-4.8)**

## Product attributes

<b>Antibody number</b>	#1057
<b>Antibody reactivity (target)</b>	CD48
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	5-4.8
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight of antigen</b>	45 kDa
<b>Synonyms</b>	B cell membrane protein; B-cell activation marker; BCM1 surface antigen; BLAST1; Signaling lymphocytic activation molecule 2; SLAM family member 2 (SLAMF2); TCT.1
<b>Human gene symbol</b>	CD48
<b>Entrez gene ID</b>	962
<b>SwissProt</b>	P09326
<b>Unigene</b>	243564
<b>Immunogen</b>	Human peripheral blood lymphocytes
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Expected antibody applications</b>	Flow, surface (published for clone), Functional studies (published for clone), WB (published for clone)
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	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, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
<b>Positive control</b>	Daudi, JY, Raji, Jurkat, and human lymphocytes. Human lymph node and tonsil.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Immunology
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	Lymphocytes, Monocytes/macrophages

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 Exp Med (1996) 184: 1769-1779. (functional studies)->/li>
2. J Immunol (2001) 166:5567-5577. (Flow; functional studies; WB)