## CD32 Monoclonal Mouse Antibody (8.7)

## **Product Description**

This MAb reacts with a CD32 (FcgRII) epitope distinct from that defined by MAb 8.26 and the epitope overlaps with that of MAb 7.30 (cluster 4). It displays a stronger reaction with Daudi than with U937 cells. The epitope is located in domain 2 of FcgRIIa. Its Fab'2 fragments block immune complex binding. CD32 (FcgRII) is a type 1 transmembrane glycoprotein that mediates several functions including phagocytosis, cytotoxicity, and immunomodulation as well as platelet aggregation. Three genes (A, B, and C) encode CD32 and at least 6 isoforms are generated via alternative mRNA splicing, i.e., Ila1, Ila2, Ilb1, Ilb2, Ilb3 and Ilc. Monocytes/macrophages, placental trophoblasts and endothelial cells express all isoforms. In addition, the Ilb isoform is expressed by B cells, and the Ila isoform by platelets, granulocytes and, weakly, by B cells. NK cells and neutrophils express Isoform Ilc. CD32 binds weakly to the Fc region of monomeric IgG but more strongly to IgG aggregates and immune complexes.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1079, Anti-CD32 (8.7)

## Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	1079		
Antibody reactivity (target)	CD32		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	8.7		
Isotype	IgG1, kappa		
Molecular weight	40 kDa		
Synonyms	CD32A; Fc fragment of IgG low affinity Ila receptor; Fc gamma RIIa; Fc-gamma RII-a; Fc-gamma-RIIa; FCGR2A; FcgRII; Fcr-2; FcRII-a; IGFR2; IgG Fc receptor II-a; Low affinity immunoglobulin gamma Fc region receptor II-a; Ly-17; Ly-m20; LyM-1; Lymphocyte antigen 17		
Human gene symbol	FCGR2A		
Entrez gene ID	2212		
SwissProt	P12318		
Unigene	352642		
Immunogen	K562 and FcgRII L cells		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human		
Species reactivity  Expected antibody applications	Human Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone)		
Expected antibody	Flow, surface (published for clone), Functional studies		
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone)  Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mlL, Flow Cytometry 0.5-1 ug/mllion cells/0.1 mL, Optimal dilution for a		
Expected antibody applications  Antibody application notes	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone)  Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mlL, Flow Cytometry 0.5-1 ug/mllion cells/0.1 mL, Optimal dilution for a specific application should be determined by user		
Expected antibody applications  Antibody application notes  Positive control	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone)  Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immonfluorescence 0.5-1 ug/mlt. Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  K-562 cells. Lymph node, tonsil and placenta.		
Expected antibody applications Antibody application notes  Positive control Shipping condition	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone)  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/mll. Flow Cytometry 0.5-1 ug/mll. Thu, Optimal dilution for a specific application should be determined by user  K-562 cells. Lymph node, tonsil and placenta.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Expected antibody applications Antibody application notes  Positive control Shipping condition Storage Conditions	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immonfluorescence: 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 K-562 cells. Lymph node, tonsil and placenta. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Expected antibody applications Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immonfluorescence: 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 K-562 cells. Lymph node, tonsil and placenta. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Expected antibody applications Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immonfluorescence: 0.5-1 ug/mlt. Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user K-562 cells. Lymph node, tonsil and placenta.  Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 9.8SA-free: 1 mg/mL in PBS/0.05% BSA, Purified: 9.8SA-free: 1 mg/mL in PBS/0.05% BSA/0.05% acide, Purified, BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% acide, Purified, BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% acide, Purified, BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% acide, Purified in PBS/0.05% BSA/0.05% acide, Purified, BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% acide, Purified in PBS/0.05% BSA/0.05% acide, Purifie		
Expected antibody applications Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immonfluorescence: 0.5-1 ug/mlt. Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user K-562 cells. Lymph node, tonsil and placenta. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% 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/.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide		

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.

## References

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

- 1. J Immunol (1993) 150(5): 1794-1803. (Flow; IP; epitope mapping)
- 2. J Exp Med (2001) 194(6): 747-756. (functional studies)
- 3. J Immunol (2014) 192:792-803. (Flow)