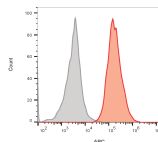


CD64 Monoclonal Mouse Antibody (10.1)



Product Description

CD64 is an Fc receptor that plays a putative role in the initiation of cell-mediated cytotoxicity. Thus far, three genes encoding four distinct CD64 transcripts have been described. CD64 has been shown to associate with signal transducing subunit of the high affinity IgE receptor. Src family kinases Hck and Lyn show increased kinase activity and will co-immunoprecipitate with CD64 subsequent to receptor cross linking. CD64 is constitutively expressed on monocytes and macrophages; exposure of granulocytes to cytokines such as IFN-gamma and G-CSF can induce CD64 expression.

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 2846, Anti-CD64 (10.1)

Product attributes

Antibody number	#2846
Antibody reactivity (target)	CD64
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	10.1
Isotype	IgG1, kappa
Molecular weight	43 kDa
Synonyms	CD64; CD64A; CD64b; CD64c; Fc fragment of IgG; high affinity Ia; receptor (CD64); FCGR1; FCGR1A; FCGR1B; FCGR1C; FcRI; FcRIB; FCRIIC; High affinity immunoglobulin gamma Fc receptor IB; IGFR1; IGFRB; IGFRIC
Human gene symbol	FCG10.1A
Entrez gene ID	2209
SwissProt	P12314
Unigene	534956; 77424
Immunogen	FCG10.1A / CD64 antibody was raised against rheumatoid synovial fluid cells and fibronectin Purified human monocytes.
Verified antibody applications	Flow (surface) (verified), IF (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Functional studies (published for clone)
Positive control	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	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.

References

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

Infect Immun (2003) 71: 5871-5880. (functional studies)

This datasheet was generated on August 1, 2025 at 07:28:13 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd64-monoclonal-mouse-antibody-10-1/>