

CLEC9A / DNGR1 Monoclonal Mouse Antibody (8F9)



Product Description

This MAb recognizes a protein of 27 kDa, identified as CLEC9A (C-type lectin domain family 9 member A). Its epitopes maps in aa 50-110. The C-type lectin/C-type lectin-like domain (CTL/CTLD) superfamily consists of a variety of proteins that share a common protein fold and have diverse functions, including cell-cell signaling, cell adhesion, glycoprotein turnover and immune responses. CLEC-9A, also known as DNGR1 (dendritic cell natural killer lectin group receptor 1), is a 241 amino acid single-pass type II membrane protein that contains one C-type lectin domain and belongs to the CTL/CTLD superfamily. Expressed in myeloid lineage cells, brain, spleen and thymus, CLEC-9A is a group V C-type lectin-like receptor (CTLR) that acts as an activation receptor.

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 3063, Anti-CLEC9A|DNGR1 (8F9)

Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#3063			
Antibody reactivity (target)	CLEC9A, DNGR1			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	8F9			
Isotype	IgG2a, kappa			
Molecular weight	27 kDa			
Synonyms	C-type lectin domain family 9 member A (CLEC9A); DNGR1			
Human gene symbol	CLEC9A			
Entrez gene ID	283420			
SwissProt	Q6UXN8			
Unigene	531189			
Immunogen	RBL-2H3 cells expressing human DNGR-1 fused to an HA epitope			
Antibody target cellular localization	Plasma membrane			
Species reactivity	Human			
Expected antibody applications	Flow, surface (published for clone), IF (published for clone)			
Antibody application notes	ELISA: For coating, order antibody without BSA; Flow cytometry: 0.5-1 ug/million cells; Immunofluorescence: 0.5-1 ug/mil. Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody			
Positive control	MCF-7 cells (FACS/IF).			
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			
Shelf life Product origin				

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,	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 Clin Invest (2008) 118(6):2098-2110. (IF, Flow)
- 2. Clinical Translational Immunol (2016) 5:e57. (Flow)

This datasheet was generated on November 30, 2025 at 03:34:25 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/clec9a-dngr1-monoclonal-mouse-antibody-8f9/