

CLEC9A / DNGR1 Monoclonal Rat Antibody (1F6)



Product Description

This Mab recognizes Mouse DNGR-1, which is also known as CLEC9A, which is a highly specific marker of the CD8a and the CD103 DC subsets, as well as a receptor able to recognize a preformed signal exposed on necrotic cells. In peripheral blood, it is highly restricted on the surface of BDCA31() dendritic cells and on a small subset of CD14() and CD16(-) monocytes.

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 2209, Anti-CLEC9A (1F6)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	2209		
Antibody reactivity (target)	CLEC9A		
Antibody type	Primary		
Host species	at		
Clonality	Monoclonal		
Clone	1F6		
Isotype	G1, kappa		
Molecular weight	27 kDa		
Synonyms	C type lectin domain family 9; member A; C-type lectin domain family 8 member A; CLEC9A; DNGR1; HEEE9341; PRO34046; UNQ9341; CD370		
Entrez gene ID	32414		
SwissProt	Q8RU4		
Unigene	391518		
Immunogen	RBL-2H3 cells expressing Mouse CLEC9A fused to an HA epitope.		
Antibody target cellular localization	Plasma membrane		
	Mouse		
Species reactivity	Mouse		
Species reactivity Expected antibody applications	Mouse Flow, surface (published for clone), Functional studies (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) Flow cytometry: 0.5-1 ug/mllion cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 0.5-1 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates		
Expected antibody applications Antibody application notes	Flow, surface (published for clone), Functional studies (published for clone) Flow cytometry: 0.5-1 ug/mllion cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 0.5-1 ug/mL; 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		
Expected antibody applications Antibody application notes Positive control	Flow, surface (published for clone), Functional studies (published for clone) Flow cytometry: 0.5-1 ug/mllion cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 0.5-1 ug/mL; 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 Mouse spleen.		
Expected antibody applications Antibody application notes Positive control Shipping condition	Flow, surface (published for clone), Functional studies (published for clone) Flow cytometry: 0.5-1 ug/million cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 0.5-1 ug/mL; 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 Mouse spleen. 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) Flow cytometry: 0.5-1 ug/million cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 0.5-1 ug/mL; 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 Mouse spleen. 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) Flow cytometry: 0.5-1 ug/million cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 0.5-1 ug/mL; 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 Mouse spleen. 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		

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. Nature (2009) 458(7240): 899-903. (functional studies)
- 2. J Immunol (2015) 195:5066-5076. (Flow, surface)

This datasheet was generated on November 29, 2025 at 10:05:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/clec9a-dngr1-monoclonal-rat-antibody-1f6/