

Product Datasheet

DC-SIGN/CD209 Antibody (111H2.02) - Azide and BSA Free DDX0209

Unit Size: 0.1 mg

Store at -20C. Avoid freeze-thaw cycles.

www.novusbio.com



technical@novusbio.com

Publications: 5

Protocols, Publications, Related Products, Reviews, Research Tools and Images at:
www.novusbio.com/DDX0209

Updated 9/9/2025 v.20.1

Earn rewards for product
reviews and publications.

Submit a publication at www.novusbio.com/publications

Submit a review at www.novusbio.com/reviews/destination/DDX0209



DDX0209

DC-SIGN/CD209 Antibody (111H2.02) - Azide and BSA Free

Product Information	
Unit Size	0.1 mg
Concentration	Please see the vial label for concentration. If unlisted please contact technical services.
Storage	Store at -20C. Avoid freeze-thaw cycles.
Clonality	Monoclonal
Clone	111H2.02
Preservative	No Preservative
Isotype	IgG2b
Purity	Ion exchange chromatography
Buffer	Tris-NaCl (pH 8.0)

Product Description	
Description	Novus Biologicals Mouse DC-SIGN/CD209 Antibody (111H2.02) - Azide and BSA Free (DDX0209) is a monoclonal antibody validated for use in IHC and Flow. Anti-DC-SIGN/CD209 Antibody: Cited in 5 publications. All Novus Biologicals antibodies are covered by our 100% guarantee.
Host	Mouse
Gene ID	30835
Gene Symbol	CD209
Species	Human
Reactivity Notes	Human
Marker	Dendritic Cell Marker
Immunogen	HeLa cells stably transfected-with human DC-SIGN
Notes	This product is manufactured by Eurobio Scientific (formerly Dendritics) and distributed by Novus Biologicals.

Product Application Details	
Applications	Flow Cytometry, Immunohistochemistry
Recommended Dilutions	Flow Cytometry, Immunohistochemistry 1:10-1:500
Application Notes	Immunohistochemistry

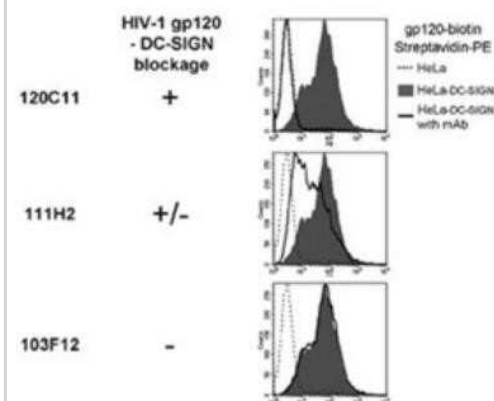


Images

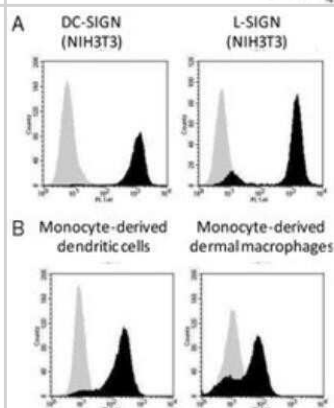
Immunohistochemistry: DC-SIGN/CD209 Antibody (111H2.02) [DDX0209] - Staining of formalin fixed and paraffin-embedded human skin sections with 111H2.02.



Flow Cytometry: DC-SIGN/CD209 Antibody (111H2.02) [DDX0209] - Analysis using the Alexa Fluor (R) 488 conjugate of DDX0209.



Flow Cytometry: DC-SIGN/CD209 Antibody (111H2.02) [DDX0209] - Analysis using the Alexa Fluor (R) 488 conjugate of DDX0209. A) Staining of NIH3T3 transfected cells with 111H2.02. B) Endogenous expression revealed by Flow staining with 111H2.02.



Publications

Schaeffer E et al. Dermal CD14(+) Dendritic Cell and Macrophage Infection by Dengue Virus Is Stimulated by Interleukin-4. *J Invest Dermatol* 2015-01-01 [PMID: 25521455]

Canard B et al. Generation of anti-DC-SIGN monoclonal antibodies capable of blocking HIV-1 gp120 binding and reactive on formalin-fixed tissue. *Immunol Lett* 2011-03-01 [PMID: 21078343]

Dakappagari N et al. Internalizing antibodies to the C-type lectins, L-SIGN and DC-SIGN, inhibit viral glycoprotein binding and deliver antigen to human dendritic cells for the induction of T cell responses. *J Immunol* 2006-01-01 [PMID: 16365436]

Soilleux EJ et al. DC-SIGN (dendritic cell-specific ICAM-grabbing non-integrin) and DC-SIGN-related (DC-SIGNR): friend or foe. *Clin Sci (Lond)* 2003-04-01 [PMID: 12653690]

Bashirova AA et al. A dendritic cell-specific intercellular adhesion molecule 3-grabbing nonintegrin (DC-SIGN)-related protein is highly expressed on human liver sinusoidal endothelial cells and promotes HIV-1 infection. *J Exp Med* 2001-03-01 [PMID: 11257134]



Novus Biologicals USA

10730 E. Briarwood Avenue
Centennial, CO 80112
USA
Phone: 303.730.1950
Toll Free: 1.888.506.6887
Fax: 303.730.1966
nb-customerservice@bio-techne.com

Bio-Techne Canada

21 Canmotor Ave
Toronto, ON M8Z 4E6
Canada
Phone: 905.827.6400
Toll Free: 855.668.8722
Fax: 905.827.6402
canada.inquires@bio-techne.com

Bio-Techne Ltd

19 Barton Lane
Abingdon Science Park
Abingdon, OX14 3NB, United Kingdom
Phone: (44) (0) 1235 529449
Free Phone: 0800 37 34 15
Fax: (44) (0) 1235 533420
info.EMEA@bio-techne.com

General Contact Information

www.novusbio.com
Technical Support: nb-technical@bio-techne.com
Orders: nb-customerservice@bio-techne.com
General: novus@novusbio.com

Products Related to DDX0209

HAF007	Goat anti-Mouse IgG Secondary Antibody [HRP]
NB7539	Goat anti-Mouse IgG (H+L) Secondary Antibody [HRP]
NBP2-27231	Mouse IgG2b Isotype Control (MPC-11)
DDX0209B	DC-SIGN/CD209 Antibody (111H2.02) [Biotin]

Limitations

This product is for research use only and is not approved for use in humans or in clinical diagnosis. Primary Antibodies are guaranteed for 1 year from date of receipt.

For more information on our 100% guarantee, please visit www.novusbio.com/guarantee

Earn gift cards/discounts by submitting a review: www.novusbio.com/reviews/submit/DDX0209

Earn gift cards/discounts by submitting a publication using this product:
www.novusbio.com/publications

