

Product Datasheet

DC-SIGN/CD209 Antibody (109H12.03) - Azide and BSA Free DDX0207

Unit Size: 0.1 mg

Store at -20C in the dark. Avoid freeze-thaw cycles.

www.novusbio.com



technical@novusbio.com

Publications: 8

Protocols, Publications, Related Products, Reviews, Research Tools and Images at:
www.novusbio.com/DDX0207

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/DDX0207



DDX0207**DC-SIGN/CD209 Antibody (109H12.03) - Azide and BSA Free**

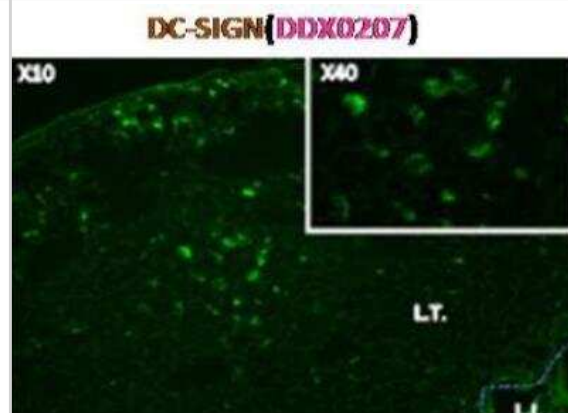
Product Information	
Unit Size	0.1 mg
Concentration	Please see the vial label for concentration. If unlisted please contact technical services.
Storage	Store at -20C in the dark. Avoid freeze-thaw cycles.
Clonality	Monoclonal
Clone	109H12.03
Preservative	No Preservative
Isotype	IgG2a
Purity	Ion exchange chromatography
Buffer	Tris-NaCl (pH 8.0)

Product Description	
Description	Novus Biologicals Mouse DC-SIGN/CD209 Antibody (109H12.03) - Azide and BSA Free (DDX0207) is a monoclonal antibody validated for use in IHC, Flow and ICC/IF. Anti-DC-SIGN/CD209 Antibody: Cited in 8 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	human DC-SIGN-transfected HeLa cells
Notes	This product is manufactured by Eurobio Scientific (formerly Dendritics) and distributed by Novus Biologicals.

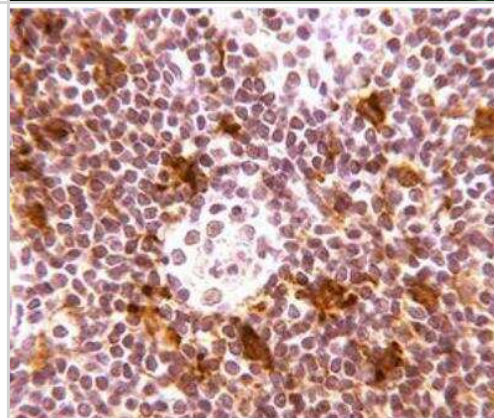
Product Application Details	
Applications	Immunohistochemistry-Paraffin, Flow Cytometry, Immunocytochemistry/ Immunofluorescence, Immunohistochemistry
Recommended Dilutions	Flow Cytometry 1:10-1:1000, Immunohistochemistry 1:10-1:500, Immunocytochemistry/ Immunofluorescence, Immunohistochemistry-Paraffin
Application Notes	Immunohistochemistry, Flow Cytometry

Images

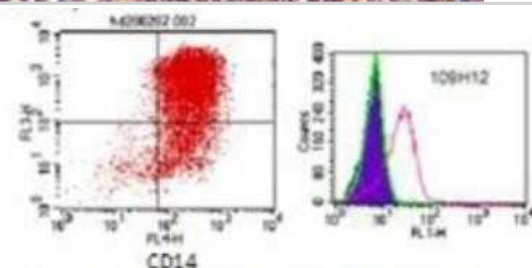
Immunocytochemistry/Immunofluorescence: DC-SIGN/CD209 Antibody (109H12.03) [DDX0207] - Murine peyer patch: DC-SIGN



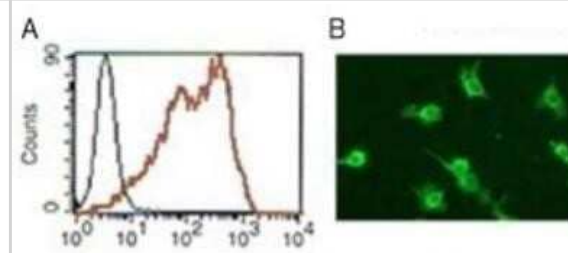
Immunohistochemistry-Paraffin: DC-SIGN/CD209 Antibody (109H12.03) [DDX0207] - Tonsil



Flow Cytometry: DC-SIGN/CD209 Antibody (109H12.03) [DDX0207] - Analysis of Alexa Fluor (R) 647 conjugate of DDX0207. FACS staining of CD14+ CD209+ cells. (CD14+ cells cultured 5 days with GM CSF and IL-10)



Flow Cytometry: DC-SIGN/CD209 Antibody (109H12.03) [DDX0207] - Analysis using the Alexa Fluor (R) 488 conjugate of DDX0207. A) Staining of DC-SIGN transfected HeLa cells. B) IF staining of DC-SIGN transfected cells with 109H12.03.



Publications

Gabriel-Segard T, Rontard J, Miny L et al. Proof-of-Concept Human Organ-on-Chip Study: First Step of Platform to Assess Neuro-Immunological Communication Involved in Inflammatory Bowel Diseases International Journal of Molecular Sciences 2023-06-24 [PMID: 37445748]

Debeer S et al. Comparative histology and immunohistochemistry of porcine versus human skin. Eur J Dermatol 2013-07-01 [PMID: 24047577]

Rochereau N et al. Phenotypic localization of distinct DC subsets in mouse Peyer Patch. Vaccine 2011-05-09 [PMID: 21439318]

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]

Kwan WH et al. Dermal-type macrophages expressing CD209/DC-SIGN show inherent resistance to dengue virus growth. PLoS Negl Trop Dis 2008-10-01 [PMID: 18827881]

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 DDX0207

HAF007	Goat anti-Mouse IgG Secondary Antibody [HRP]
NB7539	Goat anti-Mouse IgG (H+L) Secondary Antibody [HRP]
NBP1-96778	Mouse IgG2a Isotype Control (M2A)
DDX0207B	DC-SIGN/CD209 Antibody (109H12.03) [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/DDX0207

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

