

# 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](http://www.novusbio.com)



[technical@novusbio.com](mailto:technical@novusbio.com)

**Publications: 8**

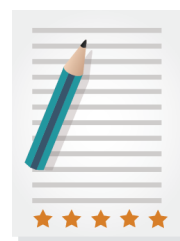
Protocols, Publications, Related Products, Reviews, Research Tools and Images at:  
[www.novusbio.com/DDX0207](http://www.novusbio.com/DDX0207)

Updated 2/21/2025 v.20.1

Earn rewards for product reviews and publications.

Submit a publication at [www.novusbio.com/publications](http://www.novusbio.com/publications)

Submit a review at [www.novusbio.com/reviews/destination/DDX0207](http://www.novusbio.com/reviews/destination/DDX0207)



**DDX0207**

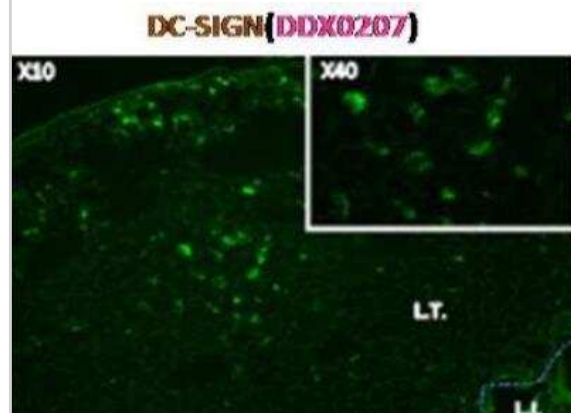
DC-SIGN/CD209 Antibody (109H12.03) - Azide and BSA Free

<b>Product Information</b>	
<b>Unit Size</b>	0.1 mg
<b>Concentration</b>	Please see the vial label for concentration. If unlisted please contact technical services.
<b>Storage</b>	Store at -20C in the dark. Avoid freeze-thaw cycles.
<b>Clonality</b>	Monoclonal
<b>Clone</b>	109H12.03
<b>Preservative</b>	No Preservative
<b>Isotype</b>	IgG2a
<b>Purity</b>	Ion exchange chromatography
<b>Buffer</b>	Tris-NaCl (pH 8.0)
<b>Product Description</b>	
<b>Host</b>	Mouse
<b>Gene ID</b>	30835
<b>Gene Symbol</b>	CD209
<b>Species</b>	Human
<b>Reactivity Notes</b>	Human
<b>Marker</b>	Dendritic Cell Marker
<b>Immunogen</b>	human DC-SIGN-transfected HeLa cells
<b>Notes</b>	This product is manufactured by Eurobio Scientific (formerly Dendritics) and distributed by Novus Biologicals.
<b>Product Application Details</b>	
<b>Applications</b>	Flow Cytometry, Immunocytochemistry/ Immunofluorescence, Immunohistochemistry, Immunohistochemistry-Paraffin
<b>Recommended Dilutions</b>	Flow Cytometry 1:10-1:1000, Immunohistochemistry 1:10-1:500, Immunocytochemistry/ Immunofluorescence, Immunohistochemistry-Paraffin
<b>Application Notes</b>	Immunohistochemistry, Flow Cytometry

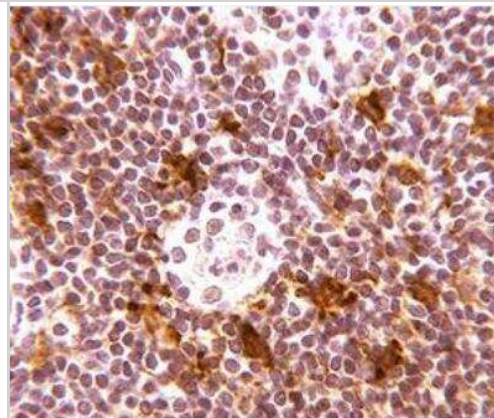


## Images

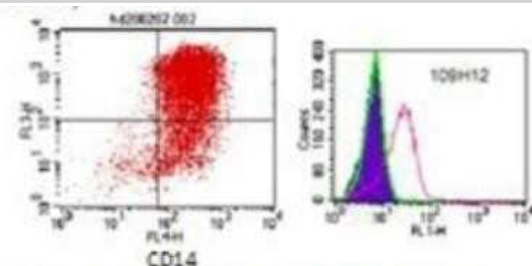
Immunocytochemistry/Immunofluorescence: DC-SIGN/CD209 Antibody (109H12.03) [DDX0207] - Murine peyer patch: DC-SIGN



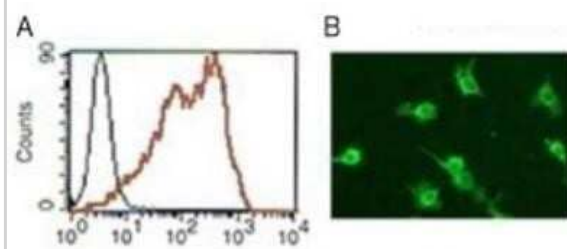
Immunohistochemistry-Paraffin: DC-SIGN/CD209 Antibody (109H12.03) [DDX0207] - Tonsil



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] (Immunocytochemistry/ Immunofluorescence)

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]
NB720-B	Rabbit anti-Mouse IgG (H+L) Secondary Antibody [Biotin]
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](http://www.novusbio.com/guarantee)

Earn gift cards/discounts by submitting a review: [www.novusbio.com/reviews/submit/DDX0207](http://www.novusbio.com/reviews/submit/DDX0207)

Earn gift cards/discounts by submitting a publication using this product:  
[www.novusbio.com/publications](http://www.novusbio.com/publications)

