

Product Datasheet

DC-SIGN/CD209 Antibody (111H2.02) [Biotin] DDX0209B

Unit Size: 0.1 mg

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

www.novusbio.com



technical@novusbio.com

Protocols, Publications, Related Products, Reviews, Research Tools and Images at:
www.novusbio.com/DDX0209B

Updated 10/23/2024 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/DDX0209B



DDX0209B

DC-SIGN/CD209 Antibody (111H2.02) [Biotin]

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	111H2.02
Preservative	No Preservative
Isotype	IgG2b
Conjugate	Biotin
Purity	Ion exchange chromatography
Buffer	Antibody in PBS and 50% Glycerol

Product Description	
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	Immunohistochemistry
Recommended Dilutions	Immunohistochemistry 1:10-1:500
Application Notes	Immunohistochemistry





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 DDX0209B

NBP2-29370	Streptavidin Native Protein
NBP2-27231B	Mouse IgG2b Isotype Control (MPC-11) [Biotin]
NBP1-76787PEP	DC-SIGN/CD209 Antibody Blocking Peptide
210-TA-005	TNF-alpha [Unconjugated]

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/DDX0209B

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

