

Product Datasheet

HIV-1 gp41 Antibody (417D3.08/ clone 2) [Biotin] DDX1305B

Unit Size: 0.1 mg

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

www.novusbio.com



technical@novusbio.com

Publications: 1

Protocols, Publications, Related Products, Reviews, Research Tools and Images at:
www.novusbio.com/DDX1305B

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/DDX1305B



DDX1305B

HIV-1 gp41 Antibody (417D3.08/ clone 2) [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	417D3.08/ clone 2
Preservative	No Preservative
Isotype	IgG1
Conjugate	Biotin
Purity	Ion exchange chromatography
Buffer	Antibody in 100 ug in 200 ul / 50 ug in 100 ul PBS and 50% Glycerol

Product Description	
Host	Mouse
Species	Virus
Reactivity Notes	virus
Immunogen	HR1-PID-HR2-transfected HEK 293 cells (HIV-1, 92BR025, Clade C)
Notes	This product is manufactured by Eurobio Scientific (formerly Dendritics) and distributed by Novus Biologicals.

Product Application Details	
Applications	ELISA, Flow Cytometry
Recommended Dilutions	Flow Cytometry 1:10-1:1000, ELISA 1:100-1:2000
Application Notes	Flow Cytometry, ELISA

Publications

Dawood R, Benjelloun F, Pin JJ et al. Generation of HIV-1 potent and broad neutralizing antibodies by immunization with postfusion HR1/HR2 complex. AIDS. 2013-03-13 [PMID: 23719346]





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 DDX1305B

NBP2-29370	Streptavidin Native Protein
NBP1-96888	Mouse IgG1 Isotype Control (MG1) [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/DDX1305B

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

