

Product Datasheet

HIV-1 gp41 Antibody (402D9.14/clone 1) - Azide and BSA Free DDX1304

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/DDX1304

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/DDX1304



DDX1304

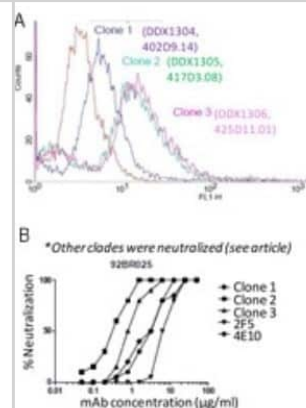
HIV-1 gp41 Antibody (402D9.14/clone 1) - Azide and BSA Free

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	402D9.14/clone 1
Preservative	No Preservative
Isotype	IgG1
Purity	Ion exchange chromatography
Buffer	Tris-NaCl (pH 8.0)
Product Description	
Description	Novus Biologicals Mouse HIV-1 gp41 Antibody (402D9.14/clone 1) - Azide and BSA Free (DDX1304) is a monoclonal antibody validated for use in ELISA and Flow. Anti-HIV-1 gp41 Antibody: Cited in 1 publication. All Novus Biologicals antibodies are covered by our 100% guarantee.
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, Block/Neutralize
Recommended Dilutions	Flow Cytometry 1:10-1:1000, ELISA 1:100-1:2000, Block/Neutralize
Application Notes	Flow Cytometry, ELISA

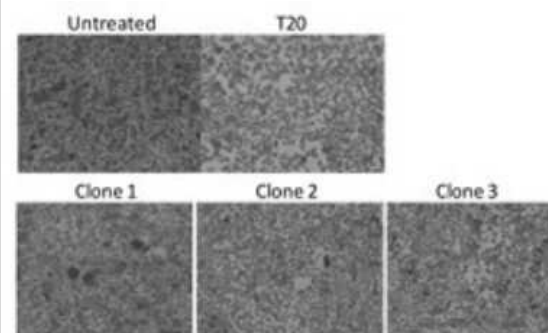


Images

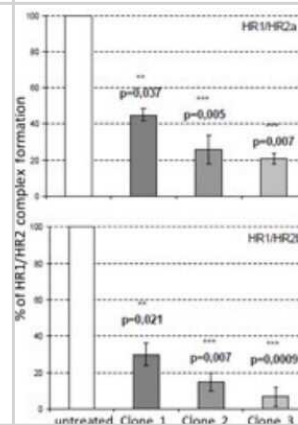
Flow Cytometry: HIV-1 gp41 Antibody (402D9.14/clone 1) [DDX1304] - Analysis using the Alexa Fluor (R) 488 conjugate of DDX1304. A) Staining of HEK293-gp41TM cell line. B) Neutralizing activity.



Blocking/Neutralizing: HIV-1 gp41 Antibody (402D9.14/clone 1) [DDX1304] - Analysis using the Alexa Fluor (R) 488 conjugate of DDX1304. Inhibition of syncytia formation.



Blocking/Neutralizing: HIV-1 gp41 Antibody (402D9.14/clone 1) [DDX1304] - Analysis using the Alexa Fluor (R) 488 conjugate of DDX1304. Inhibition of HR1/HR2 fusion complex formation.



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 DDX1304

HAF007	Goat anti-Mouse IgG Secondary Antibody [HRP]
NB7539	Goat anti-Mouse IgG (H+L) Secondary Antibody [HRP]
NBP1-97005-0.5mg	Mouse IgG1 Isotype Control (MG1)
DDX1304B	HIV-1 gp41 Antibody (402D9.14/clone 1) [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/DDX1304

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

