

Product Datasheet

HIV-1 gp41 Antibody (425D11.01/ clone 3) - Azide and BSA Free DDX1306

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/DDX1306

Updated 2/21/2025 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/DDX1306



DDX1306

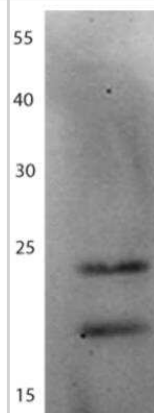
HIV-1 gp41 Antibody (425D11.01/ clone 3) - Azide and BSA Free

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	425D11.01/ clone 3
Preservative	No Preservative
Isotype	IgG1
Purity	Ion exchange chromatography
Buffer	Tris-NaCl (pH 8.0)
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	Western Blot, ELISA, Flow Cytometry, Inhibition Activity, Neutralization
Recommended Dilutions	Western Blot 1:100-1:2000, Flow Cytometry 1:10-1:1000, ELISA 1:100-1:2000, Neutralization, Inhibition Activity
Application Notes	Flow Cytometry, ELISA

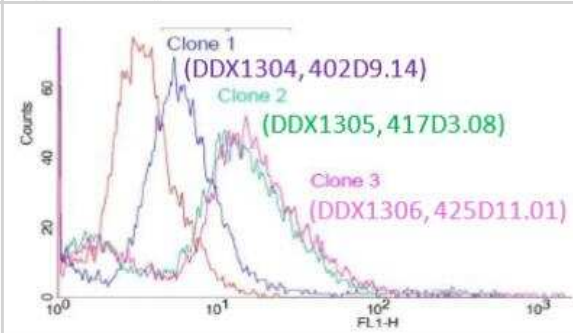


Images

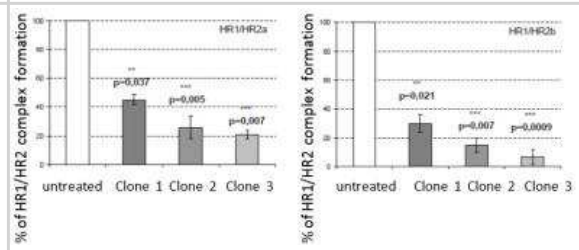
Western Blot: HIV-1 gp41 Antibody (425D11.01/ clone 3) [DDX1306] - Analysis of a purified multi-epitopic HIV antigen. The expected length of the HIV protein is 20 kDa. The second signal might represent a by-product.



Flow Cytometry: HIV-1 gp41 Antibody (425D11.01/ clone 3) [DDX1306] - FACS staining of HEK293-gp41TM cell line



Inhibition Activity: HIV-1 gp41 Antibody (425D11.01/ clone 3) [DDX1306] - Inhibition of HR1/HR2 fusion complex formation



Neutralization: HIV-1 gp41 Antibody (425D11.01/ clone 3) [DDX1306] - Neutralizing activities on laboratory strains and primary HIV-1 strains

Strain (Tropism, Clade)	Clone 1	Clone 2	Clone 3	DS	NC-1	4E10	2F5
LAU (CCR4, B)	0.78	0.16	12.5	3.1	0.53	>100	25
BAI (CCR5, B)	12.5	3.1	32.5	6.25	25	12.5	n.d.
93AG279 (CCR4, A)	0.5	0.03	1.5	0.31	0.78	0.16	50
SF162 (CCR5, B)	100	25	10	1.5	100	12.5	>100
ZMUC692 (CCR5, B)	>100	10	100	10	>100	100	n.d.
PAU640 (CCR5, B)	>100	50	>100	25	>100	50	>100
93BR825 (CCR5, B)	3.1	1.5	1.5	0.78	0.78	n.d.	n.d.
02BR825 (CCR5, C)	6.25	3.1	0.78	0.78	1.5	0.78	>100
92UK001 (CCR5, CCR4, D)	0.5	0.03	0.5	0.16	0.78	0.03	25
HIV-1 GT (CCR5, G)	12.5	3.1	32.5	3.1	6.25	1.5	50

The table contains the antibodies concentration in µg/ml of the three purified mAbs used to obtain a neutralization of HIV infection (NA50) at 50% or 90% as described in the method section. DS, NC-1, 4E10 and 2F5 mAbs were used as controls. n.d., not done. The NA50 neutralization titers have been color coded as follows: numbers with a white background indicate an K_{50} of >12.5 µg/ml, numbers with a medium grey background indicate 3.1 µg/ml, K_{50} >12.5 µg/ml and numbers with a dark grey background indicate an K_{50} of < 3.1 µg/ml.

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 DDX1306

HAF007	Goat anti-Mouse IgG Secondary Antibody [HRP]
NB720-B	Rabbit anti-Mouse IgG (H+L) Secondary Antibody [Biotin]
NBP1-97005-0.5mg	Mouse IgG1 Isotype Control (MG1)
DDX1306B	HIV-1 gp41 Antibody (425D11.01/ clone 3) [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/DDX1306

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

