

Product Datasheet

TSLP Antibody (210H1) [Biotin]

DDX0804B

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/DDX0804B

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/DDX0804B



DDX0804B

TSLP Antibody (210H1) [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	210H1
Preservative	No Preservative
Isotype	IgG2a
Conjugate	Biotin
Purity	Ion exchange chromatography
Buffer	Antibody in PBS and 50% Glycerol

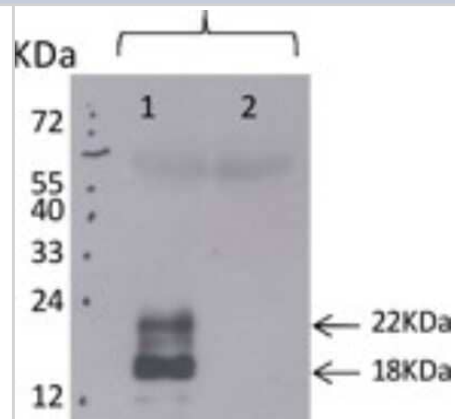
Product Description	
Host	Mouse
Gene ID	85480
Gene Symbol	TSLP
Species	Human
Reactivity Notes	Human
Immunogen	TSLP was used as immunogen.
Notes	This product is manufactured by Eurobio Scientific (formerly Dendritics) and distributed by Novus Biologicals.

Product Application Details	
Applications	Western Blot, ELISA
Recommended Dilutions	Western Blot 1:100-1:2000, ELISA 1:100-1:2000
Application Notes	Western Blot, ELISA

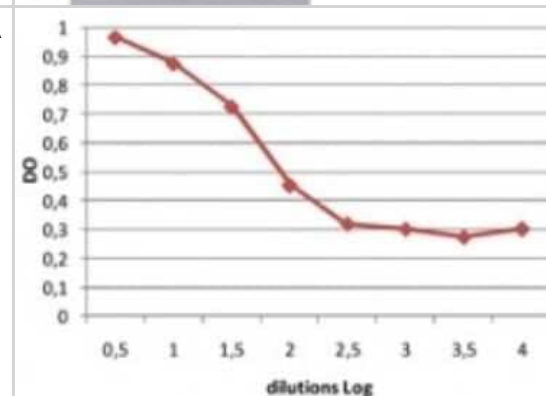


Images

Western Blot: TSLP Antibody (210H1.04) [Biotin] [DDX0804B] - Analysis of TSLP in 1) transfected 293T and 2) untransfected 293T cells using DDX0804.



ELISA: TSLP Antibody (210H1.04) [Biotin] [DDX0804B] - Capture ELISA using DDX0801 (3 ug/ml) and DDX0804B (4 ug/ml).





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 DDX0804B

NBP2-29370	Streptavidin Native Protein
IC003B	Mouse IgG2a Isotype Control (20102) [Biotin]
NBP2-35083-10ug	Recombinant Human TSLP Protein
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/DDX0804B

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

