

Product Datasheet

TLR3 Antibody (619F7.06) [Biotin] DDX0470B

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/DDX0470B

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/DDX0470B



DDX0470B

TLR3 Antibody (619F7.06) [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	619F7.06
Preservative	No Preservative
Isotype	IgM
Conjugate	Biotin
Purity	Ion exchange chromatography
Buffer	Antibody in Tris-NaCl (pH 8.0)

Product Description	
Host	Mouse
Gene ID	7098
Gene Symbol	TLR3
Species	Human, Mouse, Rat, Canine
Reactivity Notes	Human, Canine, Mouse, Rat
Immunogen	TLR3/CD283 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, Flow Cytometry
Recommended Dilutions	Western Blot 1:100-1:2000, Flow Cytometry 1:10-1:1000
Application Notes	Western Blot, Flow Cytometry





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 DDX0470B

NBP2-29370	Streptavidin Native Protein
210-TA-005	TNF-alpha [Unconjugated]
1487-TR-050	TLR3 [Unconjugated]
M6000B-1	IL-6 [HRP]

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/DDX0470B

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

