

# Product Datasheet

## MMP-12 Antibody (706F9.01) [Biotin] DDX0281B

Unit Size: 0.1 mg

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

[www.novusbio.com](http://www.novusbio.com)



[technical@novusbio.com](mailto:technical@novusbio.com)

### Publications: 2

Protocols, Publications, Related Products, Reviews, Research Tools and Images at:  
[www.novusbio.com/DDX0281B](http://www.novusbio.com/DDX0281B)

Updated 10/23/2024 v.20.1

Earn rewards for product  
reviews and publications.

Submit a publication at [www.novusbio.com/publications](http://www.novusbio.com/publications)

Submit a review at [www.novusbio.com/reviews/destination/DDX0281B](http://www.novusbio.com/reviews/destination/DDX0281B)



**DDX0281B**

MMP-12 Antibody (706F9.01) [Biotin]

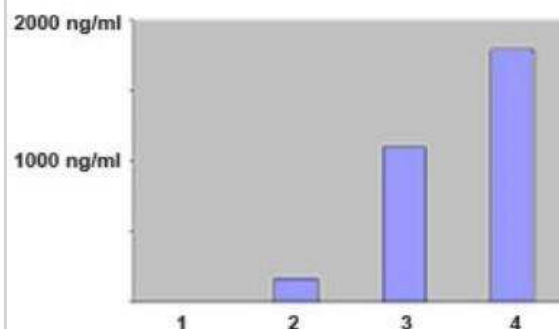
Product Information	
Unit Size	0.1 mg
Concentration	Please see the vial label for concentration. If unlisted please contact technical services.
Storage	Store at -20C in the dark. Avoid freeze-thaw cycles.
Clonality	Monoclonal
Clone	706F9.01
Preservative	No Preservative
Isotype	IgG1
Conjugate	Biotin
Purity	Ion exchange chromatography
Buffer	Antibody in Tris-NaCl (pH 8.0)

Product Description	
Host	Mouse
Gene ID	4321
Gene Symbol	MMP12
Species	Human
Reactivity Notes	Human
Immunogen	MMP-12/HME was used as immunogen.
Notes	This product is manufactured by Eurobio Scientific (formerly Dendritics) and distributed by Novus Biologicals.

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

**Images**

ELISA: MMP-12 Antibody (706F9.01) [Biotin] [DDX0281B] - 1= control; 2=HME-transfected cells; 3=CD34+ +GMCSF+TNFa+IL4; 4=patient serum of Langerhans hystiocytosis (samples were tested with DDX0281).



## Publications

Illumets H, Ryttilä P, Demedts I et al. Matrix metalloproteinases -8, -9 and -12 in smokers and patients with stage 0 COPD *Int J Chron Obstruct Pulmon Dis* 2007-01-01 [PMID: 18229576]

### Details:

This citation used the HRP version of this antibody.

Demets IK et al. Elevated MMP-12 protein levels in induced sputum from patients with COPD. *Thorax* 2006-03-01 [PMID: 16308335]





### **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 DDX0281B**

---

NBP2-29370	Streptavidin Native Protein
NBP1-96888	Mouse IgG1 Isotype Control (MG1) [Biotin]
210-TA-005	TNF-alpha [Unconjugated]
917-MPB-020	MMP-12 [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](http://www.novusbio.com/guarantee)

Earn gift cards/discounts by submitting a review: [www.novusbio.com/reviews/submit/DDX0281B](http://www.novusbio.com/reviews/submit/DDX0281B)

Earn gift cards/discounts by submitting a publication using this product:  
[www.novusbio.com/publications](http://www.novusbio.com/publications)

