HA.11 monoclonal antibody (16B12) (purified)

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

Ordering Information

Order Online »

ENZ-ABS120-0500	0.5ml
ENZ-ABS120-1000	1ml
ENZ-ABS120-0200	200μΙ

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Hemagglutinin tag

Application IF, IHC, IP, WB

Clone 16B12

Formulation Liquid. In PBS containing 50% glycerol and 0.03%

thimerosal.

Host Mouse

Immunogen Twelve amino acid peptide CYPYDVPDYASL.

lgG1

Recommendation Dilutions/ConditionsWestern Blot (1:2,000-1:20,000).Immunoprecipitation

(1:50-1:100).Immunohistochemistry (1:1,000).Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Specificity Recognizes the influenza hemagglutinin epitope

(YPYDVPDYA) which has been used extensively as a general epitope tag in expression vectors. Recognizes HA

epitopes located in the middle of protein sequences as

well as at the N- or C-terminus.

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com