

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

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

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

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.

Isotype IgG1

Recommendation Dilutions/Conditions Western 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](#)



ENZO LIFE SCIENCES,
Last modified: May 29, 2024
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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