

HA.11 monoclonal antibody (16B12)

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS118-0500	0.5ml
ENZ-ABS118-1000	1ml
ENZ-ABS118-0200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°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. Ascites.

Host Mouse

Immunogen Twelve amino acid peptide CYPYDVPDYASL.

Isotype IgG1

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,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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