

# $\beta$ -Amyloid (1-40) monoclonal antibody (11A50-B10)

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

## Ordering Information

[Order Online »](#)

ENZ-ABS116-0200	200 $\mu$ l
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application	ELISA, IHC, WB
Clone	11A50-B10
Formulation	Liquid. In PBS. Does not contain any preservatives or carrier proteins.
Host	Mouse
Isotype	IgG1
Purity Detail	Protein G purified.
Species Reactivity	Human, Mouse
Specificity	Recognizes C-terminal $\beta$ -amyloid. Specific for the isoform ending at aa40.
UniProt ID	P05067
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)