

# APP monoclonal antibody (LN27)

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

## Ordering Information

[Order Online »](#)

ENZ-ABS126-1000	1ml
ENZ-ABS126-0200	200µl
ENZ-ABS126-0500	500µl

## 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      Amyloid precursor protein, Amyloid A4 protein

Application      WB

Clone      LN27

Formulation      Liquid. In PBS. Does not contain any preservatives or carrier proteins.

Host      Mouse

Isotype      IgG1

Purity Detail      Protein G purified.

Species Reactivity      Human

Specificity      Recognizes all three known forms of APP ( $\beta$ -APP695,  $\beta$ -APP751,  $\beta$ -APP770). Epitope lies within amino acids 45-53 of amyloid precursor protein.

UniProt ID      P05067

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)