

# APP monoclonal antibody (3E9)

Deposition of APP (Amyloid A4 protein) precursor intraneuronally as neurofibrillary tangles, extracellularly as plaques, and in blood vessels, is characteristic of both Alzheimers disease and aged Downs Syndrome. There are five forms of APP: APP395, APP563, APP695, APP751, and APP770. They are produced by alternative splicing of the APP gene. APP695 is the most predominant form in neuronal tissue while APP751 is the predominant form found elsewhere.

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

Citations: 2 [View Online »](#)

Ordering Information [Order Online »](#)

ADI-NBA-100-E	100µg
---------------	-------

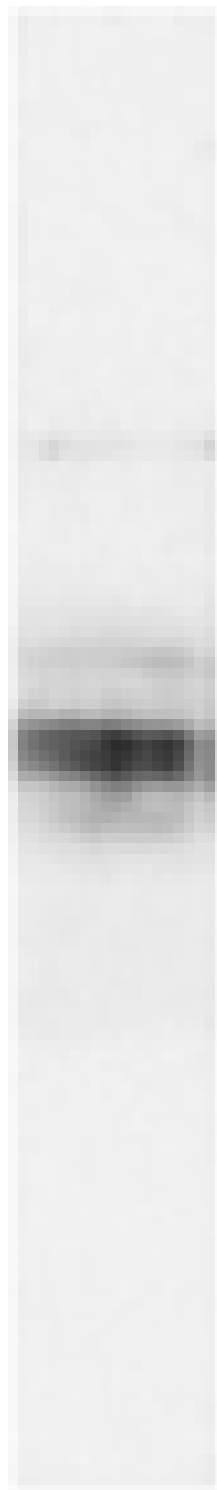
Manuals, SDS & CofA [View Online »](#)

**kDa**

**170 -**

**116 -**

**76 -**



Western blot analysis of Amyloid Precursor Protein expression in mouse brain using APP mAb (3E9).

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Amyloid A4 protein, Amyloid protein

**Application** WB

**Application Notes** Detects a band of ~102kDa by Western blot.

**Clone** 3E9

**Formulation** Liquid. In PBS pH7.2 containing 50% glycerol.

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to a portion of human APP695.

**Isotype** IgG1

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (5-10µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human, Mouse

**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)