

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
---------------	-------

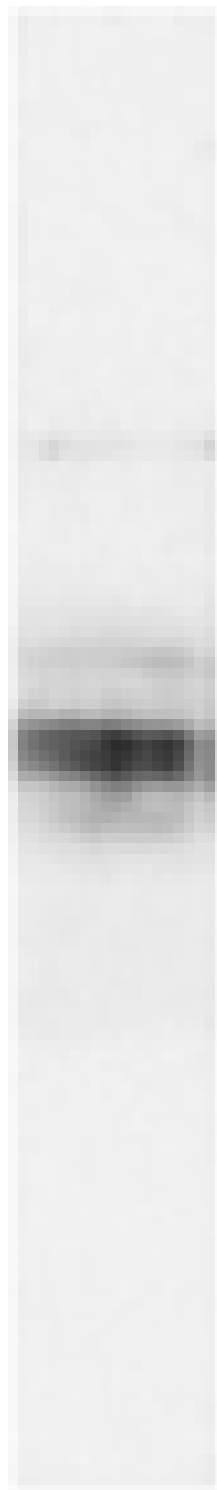
Manuals, SDS & CofA [View Online »](#)

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-
fr@enzolifesciences.com](mailto:info-fr@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)

Last modified: May 29, 2024