

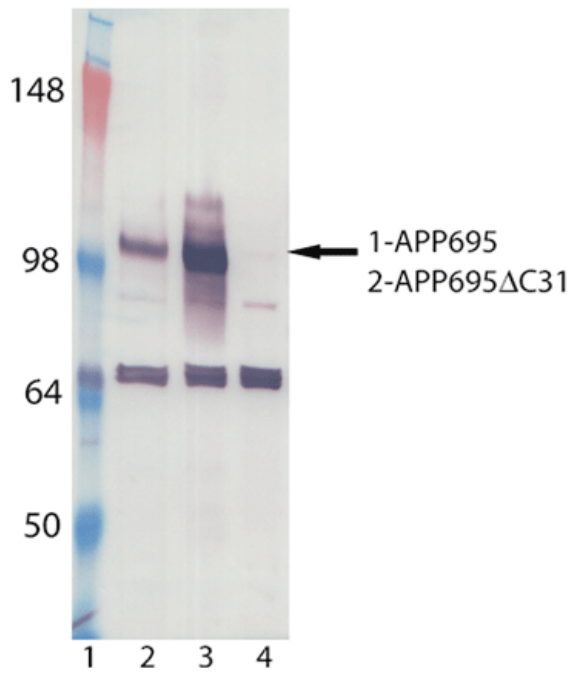
# APP polyclonal antibody

Deposition of APP (Amyloid A4 protein) precursor intraneuronally as neurofibrillary tangles, extracellularly as plaques, and in blood vessels, is characteristic of both Alzheimer's disease and aged Down's 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](#).

**Manuals, SDS & CofA**

[View Online »](#)



HEK cells were transfected with pCDNA3-APP695 (lane 2), pCDNA3-APP695  $\Delta$ C31 (lane 3) or pCDNA3 (lane 4) constructs. 10 $\mu$ g of total cell lysates were separated by 7.5% SDS-PAGE for Western Blot. The membrane was immunoblotted with primary antibody (ENZ-ABS226, APP, pAb) at 1:1000 dilution. Anti-rabbit IgG-AP (Prod. No. ADI-SAB-301) was used at 1:4000 dilution. The result showed this antibody detected both full length APP695 and APP695  $\Delta$ C31 at about 98 KDa. Pre-stained protein marker is shown in lane 1.

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Amyloid A4 protein, Amyloid protein, ABPP, APPI, Amyloid beta A4 protein
Application	ELISA, IP, WB
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human APP.
Purity Detail	Peptide affinity purified.
Source	Purified from rabbit serum.
Species Reactivity	Human, Mouse
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-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)