

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](#).

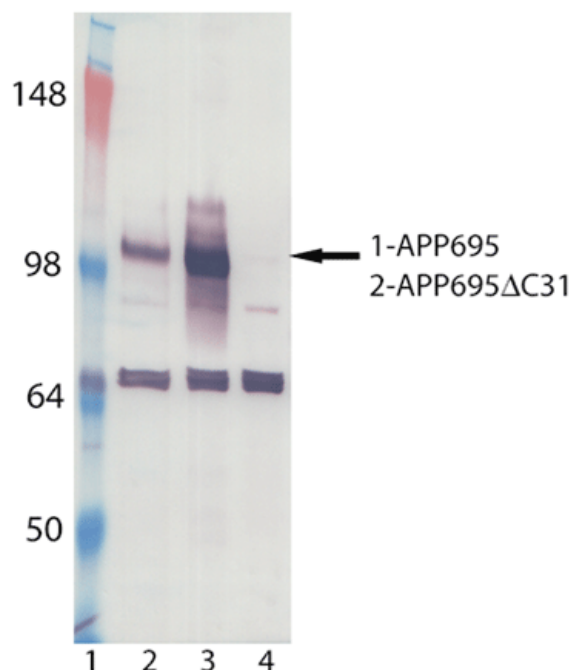
Ordering Information

[Order Online »](#)

ENZ-ABS226-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



HEK cells were transfected with pCDNA3-APP695 (lane 2), pCDNA3-APP695 Δ C31 (lane 3) or pCDNA3 (lane 4) constructs. 10 μ 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 Δ 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 [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)