

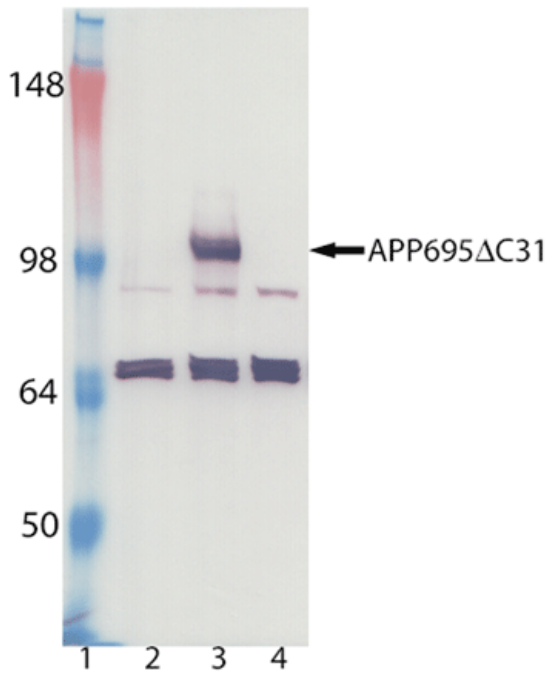
APP Δ C31 polyclonal antibody

Alzheimer's disease (AD) is a progressive neurodegenerative disease characterized by the senile plaques, neurofibrillary tangles and loss of synapses and neurons. AD has been largely viewed as a disease of toxicity being mediated by the accumulation of the amyloid beta ($A\beta$) peptide as plaques within the brain resulting in damage to brain cells from the binding of damaging metals, reactive oxygen species production and direct damage to cellular membranes. Recent research has suggested that the $A\beta$ peptide is a multifunctional peptide with non-pathological effects and that its association with AD is in conjunction with its roles in combination with other proteins such as the amyloid precursor protein (APP) resulting in the imbalance between the processes of memory formation and normal forgetting. It is through the interactions of the $A\beta$ peptide with APP that the $A\beta$ peptide itself can affect normal modulation and signaling of APP resulting in its indicated role in the pathogenesis of AD via signaling effects rather than chemical or physical effects. There are three major APP isoforms (APP695, APP751 and APP770) that are formed through alternative splicing of precursor mRNA. APP770 represents the canonical sequence. The APP695 isoform is preferentially expressed in the central nervous system, while APP770 and APP751 are more highly expressed in peripheral tissues. **It has been demonstrated that the full length APP695 can be cleaved via caspase at an intracellular site (Asp664) resulting in the release of a 31 amino acid C-terminal peptide (C31) from the remaining larger neo-APP fragment (APP Δ C31) with both of these entities being pro-apoptotic.** Immunohistochemical analysis of human brain tissue demonstrated that this cytoplasmic cleavage occurs 4-fold greater in patients with AD versus normal patients and that these cleavage products are localized to plaques and tangles in key areas of the brain affected by the disease. A single genetic mutation of aspartic acid residue 664 to alanine of APP695 led to the complete blockage of the C-terminal cleavage *in vivo* reversing many characteristics of the AD phenotype in a transgenic mouse model. Additionally, in cell culture it has been suggested that the neurotoxicity of $A\beta$ is dependent on the cleavage of APP at Asp664 and the resulting $A\beta$ -facilitated APP multimerization.

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

Citations: 2

[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-ABS445, APP Δ C31, 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 only detected 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 C31 chain, ABPP, APPI

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](#)

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-
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)