

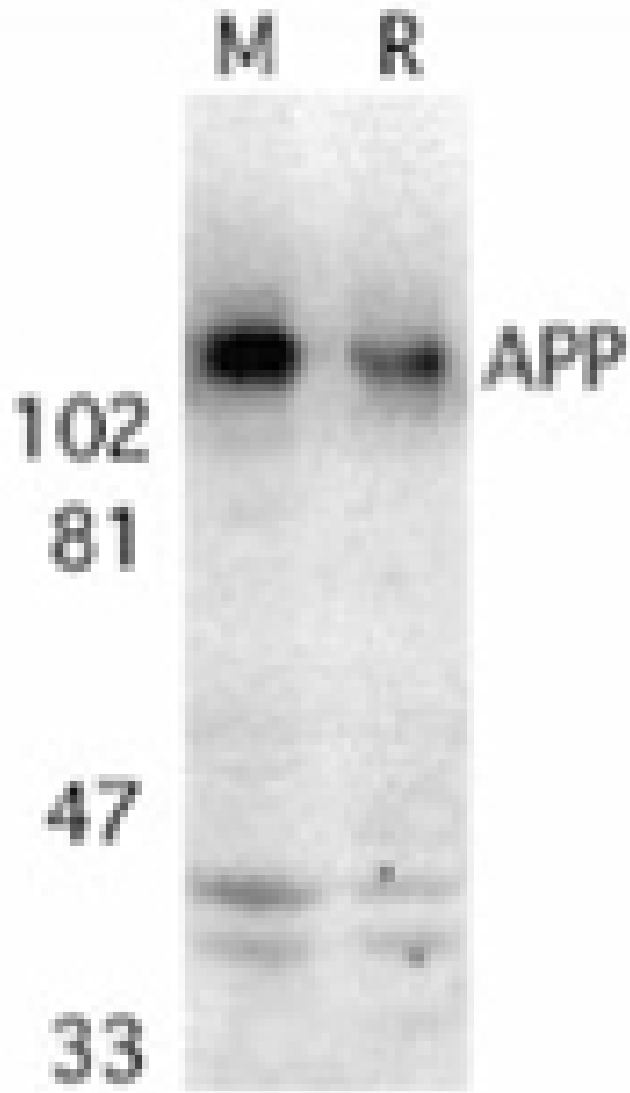
# APP polyclonal antibody

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.

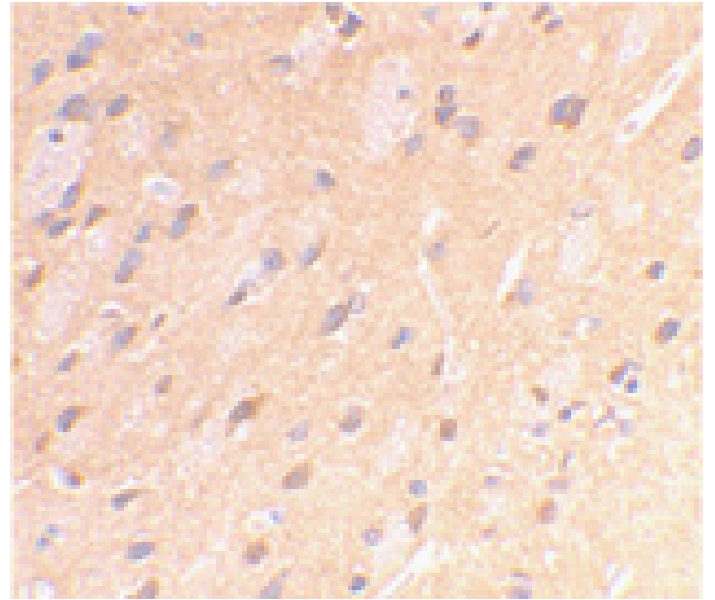
This antibody is covered by our [Worry-Free Guarantee](#).

**Manuals, SDS & CofA**

[View Online »](#)



Western blot analysis of APP in mouse (M) and rat (R) brain tissue lysates with APP antibody at 2 $\mu$ g/ml.



Immunohistochemistry analysis of rat brain using APP antibody at 2 $\mu$ g/ml.

## 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** IHC, WB

**Application Notes** Immunohistochemistry  
Western Blot (1µg/ml, ECL)  
Optimal conditions must be determined individually for each application.

**Formulation** Liquid. In PBS containing 0.02% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 737-751 of human APP or aa 85-99 of human C99 (a fragment generated by β-secretase cleavage). The sequence of this peptide is completely conserved in mouse and rat APP.

**Purity Detail** Peptide affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1µg/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, Rat

**UniProt ID** P05067

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



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430

Last modified: May 29, 2024

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

France  
Phone: +33 472 440 655

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)