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.

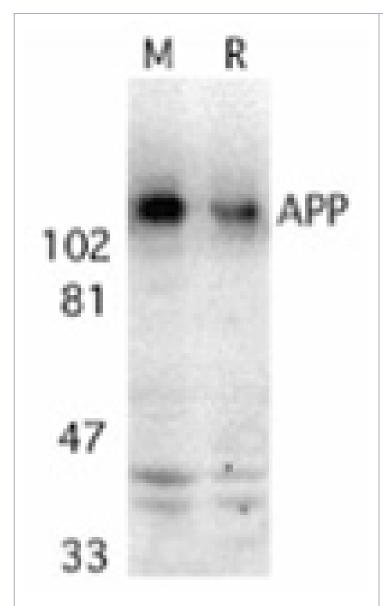
Ordering Information

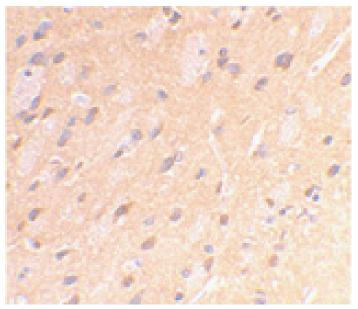
Order Online »

ADI-NBA-102-E 100μg

Manuals, SDS & CofA

View Online »





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

Western blot analysis of APP in mouse (M) and rat (R) brain tissue lysates with 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 42024 om

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com