

Instruction Manual

Catalog Number	PK-AB718-2133
Synonyms	APP Antibody: APP AbNT , APP AbNT
Description	Accumulation of the amyloid- β peptide (A β) in the cerebral cortex is a critical event in the pathogenesis of Alzheimer's disease. The β -amyloid protein precursor (APP) is cleaved by β -secretase, producing a soluble derivative of the protein and a membrane anchored 99-amino acid carboxy-terminal fragment (C99). The C99 fragment serves as substrate for γ -secretase to generate the 4 kDa amyloid- β peptide (A β), which is deposited in the brains of all sufferers of Alzheimer's disease.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal APP antibody was raised against a peptide corresponding to amino acids of human amyloid A4 protein precursor (APP) corresponding to the amino terminus of the 4K A β peptide generated by (- and (- secretases . The peptide sequences are identical to those of rabbit, pig, bovine, guinea pig, and chicken.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ l or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. APP antibody can be used for detection of APP and A β peptide by Western blot at 1 μ g/mL. Antibody can also be used for immunohistochemistry starting at 10 μ g/mL.
Images	Available upon request.
References	Ponte P, Gonzalez-DeWhitt P, Schilling J, Miller J, Hsu D, Greenberg B, Davis K, Wallace W, Lieberburg I, Fuller F. A new A4 amyloid mRNA contains a domain homologous to serine proteinase inhibitors. Nature 1988;331:525-7. Selkoe D.J. Cell biology of the amyloid β -protein precursor and the mechanism of Alzheimer's disease. Annu Rev Cell Biol 1994;10:373-403 (RD1000)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2133P Mouse brain Tissue Lysate, Cat. No. PK-AB718-1403

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.