



Goat (polyclonal) Anti-Human Presenilin-1 NH₂-Terminal Peptide Antiserum

PRODUCT ANALYSIS SHEET

Catalog Number:	AHB0181
Lot Number:	See product label
Quantity/Volume:	100 µL
Form of Antibody:	Processed whole goat serum
Preservation:	0.1% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
Immunogen:	A synthetic peptide corresponding to aa sequence 14-33 (₁₄ AQMSEDNHLSTVRSQNDNR ₃₃) of the N-terminal of human presenilin-1 protein.
Target:	This antibody recognizes the intact 48 kDa presenilin-1 (PS-1) protein, as well as the N-terminal fragment of the cleaved form of PS-1. PS-1 is observed in brain and cells of neuronal origin. The PS-1 gene is located on chromosome 14. Mutations in PS-1 may be responsible for an early onset form of Alzheimer's disease (AD). PS-1 has now been identified as one of the γ -secretase proteases involved in cleaving the transmembrane domain of Amyloid Precursor Protein (APP) following α and β secretase cleavage. Along with its role in β -amyloid protein formation from APP, PS-1 also is responsible for proteolytic cleavage of the C-terminus of the intracellular protein, Notch1. This antiserum can routinely be used without further purification.
Species Reactivity:	Human. Other species were not tested.
Applications:	This antiserum is suitable for use in ELISA, immunohistochemistry and Western blot analysis.
Suggested Working Dilutions:	For ELISA, a dilution of $\geq 1:5,000$ is recommended. For immunohistochemistry using formalin-fixed, paraffin-embedded sections, a dilution of $\geq 1:100$ and paraformaldehyde-fixed, frozen sections, a dilution of $\geq 1:1,000$ is recommended. For Western blots, a dilution of $\geq 1:1,000$ is recommended. The optimal concentration should be determined for each specific application.
Recommended Positive Control:	Human SH-SY5Y neuroblastoma cell lysates.
Storage:	Store at 2-8°C. For long term storage, apportion into working aliquots and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.
References:	Johnsingh, A. A., <i>et al.</i> (2000) Altered binding of mutated presenilin with cytoskeleton-interacting proteins. FEBS Letters 465:53-58.

This product is for research use only. Not for use in diagnostic procedures.

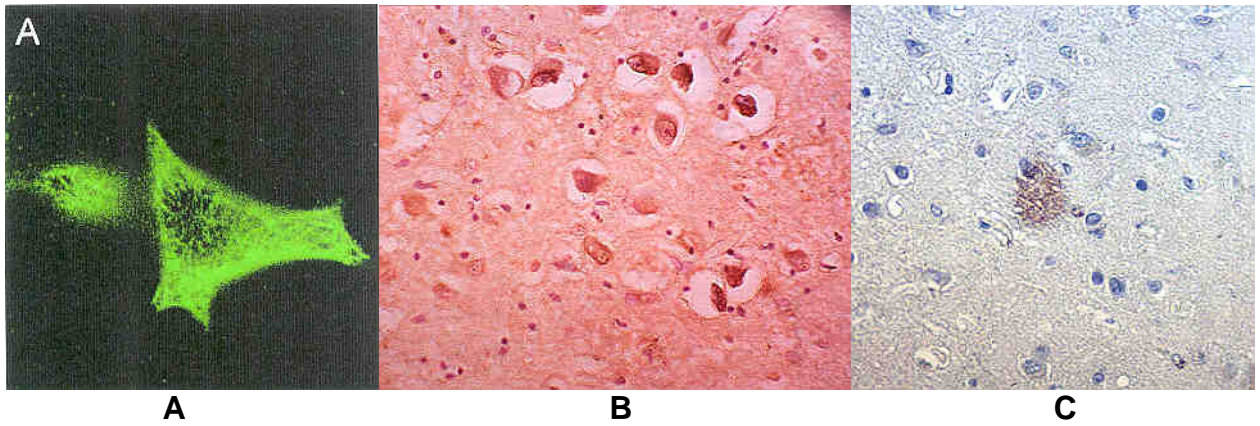
www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

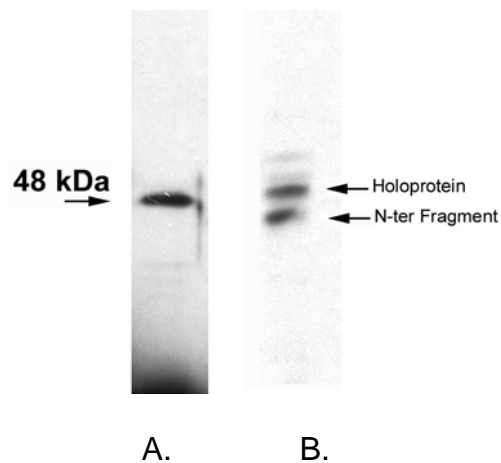
PI AHB0181

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



Immunofluorescent staining of cultured human SH-SY5Y neuroblastoma cells using the presenilin-1 antiserum (Figure A). Alzheimer's disease brain showing immunohistochemical staining of neurons (Figure B) and senile plaque (Figure C) in formalin-fixed, paraffin-embedded sections using the presenilin-1 antiserum at 1:300 dilution.



Western blot analysis using the presenilin-1 antiserum at 1:1,000 dilution. Figure A. shows SH-SY5Y human neuroblastoma cells where the presenilin-1 protein is detected. Figure B. shows human brain hippocampus where the intact presenilin-1 holoprotein and the N-terminal cleavage fragment are detected.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI AHB0181

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.