



**Mouse (monoclonal)
Anti-A β [1-16]
Unconjugated**

PRODUCT ANALYSIS SHEET

Catalog Number:	44-352
Lot Number:	See product label
Quantity/Volume:	See product label
Clone:	6E10
Isotype:	IgG1
Form of Antibody:	Purified immunoglobulin in phosphate buffered saline. Preservative and carrier-free.
Immunogen:	The immunogen used in the production of this monoclonal antibody is a chemically synthesized peptide corresponding to amino acid residues 1-16 of A β (β -amyloid) peptide.
Purification:	Purified from ascites by protein G affinity chromatography.
Species Reactivity:	This antibody reacts with human A β peptide. The epitope lies within amino acid residues 3-8 of β -amyloid (A β) peptide.
Applications:	The antibody has been used in the immunohistochemical detection of A β in fixed sections of Alzheimer's Disease (AD) brain. When used in immunohistochemistry, the epitope must be re-exposed using either formic acid treatment (70% for 10-30 minutes at room temperature) or heat induced epitope retrieval. This antibody has also been used in other procedures including: ELISA, immunoprecipitation, and Western blotting.
Suggested Working Dilutions:	For immunohistochemistry, we recommend using a 1:100 to 1:1000 dilution of the antibody. The optimal antibody concentration should be determined for each specific application.
Storage:	Store at $\leq -20^{\circ}\text{C}$. Upon initial thawing, apportion into working aliquots and store at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.
Expiration Date:	Expires one year from date of receipt when stored as instructed.

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

PI44352

(Rev 11/08) 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.

References:

- Kim, K.S. et al. (1988) Production and characterization of monoclonal antibodies reactive to synthetic cerebrovascular amyloid peptide. *Neurosci. Res. Commun.* 2: 121-131.
- Kim, K.S. et al. (1990) Detection and quantitation of amyloid B-peptide with two monoclonal antibodies. *Neurosci. Res. Commun.* 7:113-123.
- Kitaguchi, T. et al. (1990) Beta-protein immunoreactivity in brains of patients with neuronal ceroid lipofuscinosis: ultrastructural and biochemical demonstration. *Neurosci. Lett.* 112(2-3):155-160.
- Melchor, J.E. and W.E. van Nostrand (2000) Fibrillar amyloid β -protein mediates the pathologic accumulation of its secreted precursor in human cerebrovascular smooth muscle cells. *J. Biol. Chem.* 275(13):9782-9791 (cites the use of this antibody in Western blotting performed under non-reducing conditions).
- Koldamova, R.P. et al. (2001) Apolipoprotein A-I directly interacts with amyloid precursor protein and inhibits A beta aggregation and toxicity. *Biochemistry* 40(12):3553-3560 (cites the use of this antibody in Western blotting performed under reducing conditions [mercaptoethanol]).
- Kusiak, J.W. et al. (2001) Expression of mutant amyloid precursor proteins decreases adhesion and delays differentiation of Hep-1 cells. *Brain Res.* 896:146-152 (cites the use of this antibody in immunoprecipitation in a system for the detection of sAPP α).
- Chishti, M.A. et al. (2001) Early-onset amyloid deposition and cognitive deficits in transgenic mice expressing a double mutant form of amyloid precursor protein 695. *J. Biol. Chem.* 276(24):21562-21570 (cites the use of this antibody in immunohistochemistry).
- Lefterov, I.M. et al. (2000) Human bleomycin hydrolase regulates the secretion of amyloid precursor protein. *FASEB J.* 14:1837-1847.

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