



**Mouse (monoclonal)
Anti-A β [17-24]
Unconjugated**

PRODUCT ANALYSIS SHEET

Catalog Number:	44-356
Lot Number:	See product label
Quantity/Volume:	See product label
Clone:	4G8
Form of Antibody:	Purified monoclonal antibody in phosphate buffered saline. Carrier protein and preservative-free.
Immunogen:	The immunogen used in the production of this monoclonal antibody is a chemically synthesized peptide corresponding to amino acid residues 17-24 of A β (β -amyloid) peptide.
Purification:	Purified from ascites by Protein A affinity chromatography.
Isotype:	IgG2b
Species Reactivity:	Human, mouse, and rat A β .
Applications:	The antibody is suitable for use in the detection of A β by immunohistochemistry with fixed sections of Alzheimer's Disease (AD) brain. The epitope for this antibody must be re-exposed by treatment of fixed tissue sections with 70% formic acid for 10-30 minutes at room temperature or by heat-induced epitope retrieval. This antibody is also suitable for use in Western blot analysis, ELISA, immunoprecipitation, and immunoaffinity purification.
Suggested Working Dilutions:	For immunohistochemistry, the recommended working dilution is 1:500 to 1:5,000. For Western blotting, the recommended working dilution is 1:100-1:1,000. For ELISA, the recommended working dilution is 1:1,000-1:10,000. The optimal antibody concentration should be determined for each specific application.
Storage:	Store at -80°C . Upon initial thawing, apportion into working aliquots and store at -80°C . Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
References:	<p>Kim, K.S. et al. (1988) Production and characterization of monoclonal antibodies reactive to synthetic cerebrovascular amyloid peptide. <i>Neurosci. Res. Commun.</i> 2: 121-131.</p> <p>Kim, K.S. et al. (1990) Detection and quantitation of amyloid B-peptide with two monoclonal antibodies. <i>Neurosci. Res. Commun.</i> 7:113-123.</p>

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

PI44356

(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 (cont'd):

Chisti, 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.

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).

Nunomura, A. et al. (2000) Neuronal oxidative stress precedes amyloid- β deposition in Down Syndrome. *J. Neuropathol. Exp. Neurol.* 59(11):1011-1017 (cites the use of this antibody in immunohistochemistry).

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

PI44356

(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.