

Beta Amyloid, 1-16 (6E10) Monoclonal Antibody (SIGNET) Exclusively Provided by Covance

Catalog Number: SIG-39300

Available Size: 0.2 mL, 0.5 mL, 1.0 mL

Description: Beta Amyloid (A β) Monoclonal Antibody, 6E10
(Formerly Signet Catalog Nos: 9300-02, 9300-05, 9300-10)

Covance is the Exclusive Provider of this Antibody Product

Intended Use: **Research Use Only (RUO)**

This product is sold for laboratory research use only, not for human or in-vivo use.

Clone: 6E10

Form: Ascites

Host: Mouse

IsoType: IgG1

Species Reactivity: Human

Specificity: This antibody is reactive to amino acid residues 1-16 of beta amyloid. The epitope lies within amino acids 3-8 of beta amyloid (EFRHDS). Beta amyloid forms are deposited in the CNS of patients with Alzheimer's disease and Down's syndrome. Signet's beta amyloid antibody 6E10 reacts to the abnormally processed isoforms, as well as precursor forms.

Uses: This antibody is effective in immunoblotting, immunohistochemistry (IHC),

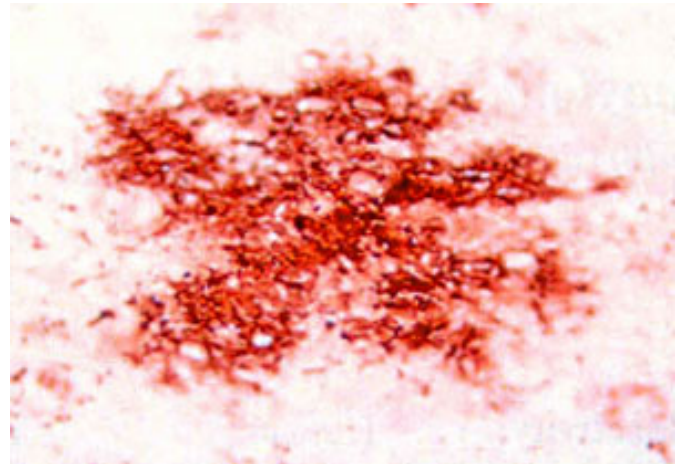


Image Description: 6E10 immunostaining on brain section of Tg-SweArc transgenic mice.

immunoaffinity purification, immunoprecipitation, and ELISA.

Suggested Working Dilution: The optimal working dilution should be determined for each specific assay condition.

- **IHC:** 1:100 - 1:1,000

Tissue Sections: Formalin-fixed human and animal paraffin-embedded brain sections

Pretreatment: Formic acid (70%) for 10-30 minutes at room temperature.

- **ELISA:** 1:1,000
- 1:10,000

Storage: Store at -20°C or below. Upon initial thawing, apportion into working aliquots and store at -20°C or below. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody. Do not store in frost-free freezers.

References: Thakker DR, et al. Intracerebroventricular amyloid-beta antibodies reduce cerebral amyloid angiopathy and associated micro-hemorrhages in aged Tg2576 mice. *Proc Natl Acad Sci USA Feb 25, 2009.*

Kimura N, et al. Age-related changes of intracellular Abeta in cynomolgus monkey brains. *Neuropath. Appl. Neurobiol.* 31(2):170-80, 2005.

Klyubin I, et al. Amyloid beta protein immunotherapy neutralizes Abeta oligomers that disrupt synaptic plasticity in vivo. *Nat. Med.* 11(5):556-61, 2005.

Iijima K, et al. Dissecting pathological effects of human Abeta40 and Abeta42 in Drosophila: a potential model for Alzheimer's disease. *PNAS* 101(17):6623-6628, 2004.

Li M, et al. Macrophage colony stimulatory factor and interferon-gamma trigger distinct mechanisms for augmentation of beta-amyloid-induced microgliamediated neurotoxicity. *J. Neurochem.* 91:623-33, 2004.

Poduloso JF, et al. Design and chemical synthesis of a magneticresonance contrast agent with enhanced in vitro binding, high blood-brain barrier permeability, and in vivo targeting to Alzheimer's disease amyloid plaques. *Biochem.* 43:6064-6075, 2004.

Venezia V, et al. Apoptotic cell death influences the signaling activityof the amyloid precursor protein through SchA and Grb2 adaptor proteins inneuroblastoma SH-SY5Y cells. *J. Neurochem.* 90:1359-1370, 2004. Kim KS, et al. Production and characterization of monoclonalantibodies reactive to synthetic cerebrovascular amyloid peptide. *Neurosci. Res. Comm.* 2130, 1988.

Kim KS, et al. *Neurosci. Res. Comm.* 7 1988.

Warranty/Conditions: Covance products may not be resold or modified for resale without prior written approval.

Related Products:

- BetaMark™ Beta-Amyloid x-40 Chemiluminescent ELISA Kit
Catalog Number SIG-38950
- BetaMark™ Beta-Amyloid x-42 Chemiluminescent ELISA Kit
Catalog Number SIG-38952
- Colorimetric BetaMark™ Beta-Amyloid x-40 ELISA Kit
Catalog Number SIG-38954
- Colorimetric BetaMark™ Beta-Amyloid x-42 ELISA Kit
Catalog Number SIG-38956
- Beta Amyloid, 1-16 (6E10) Monoclonal Antibody, Purified (SIGNET) *Exclusively Provided by Covance*
Catalog Number SIG-39320
- Beta Amyloid, 1-16 (6E10) Monoclonal Antibody, Biotin Labeled (SIGNET) *Exclusively Provided by Covance*
Catalog Number SIG-39340
- Beta Amyloid, 1-16 (6E10) Monoclonal Antibody, F(Ab')₂ Fragments (SIGNET) *Exclusively Provided by Covance*
Catalog Number SIG-39330
- Beta Amyloid, 1-16 (6E10) Monoclonal Antibody, HRP Labeled (SIGNET) *Exclusively Provided by Covance*
Catalog Number SIG-39345

Product Revision Date: 9/11/2009