

**Beta Amyloid 1-40 (11A50-B10)
Monoclonal Antibody, Purified (SIGNET)
Catalog Number: SIG-39140
Available Size: 0.2 mL, 0.5 mL, 1.0 mL**

Description: Monoclonal Antibody against Beta Amyloid (A β) 1-40 (11A50-B10)
(Formerly Signet Catalog Nos.: 9140-02, 9140-05, 9140-10)

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

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

Clone: 11A50-B10

Form: Protein G Purified Antibody (in PBS; no preservatives or carrier proteins)

Host: Mouse

IsoType: IgG1

Species Reactivity: Human, Mouse

[Ab]: 0.5 mg/mL

Specificity: Clone 11A50-B10 is reactive to the c-terminus of beta amyloid and is specific for the isoform ending at the 40th amino acid.

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

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

- **Western:** 1:100-1:1,000
- **IHC:** 1:100-1:250

Tissue Sections: Formalin-fixed, paraffin-embedded tissues

Pretreatment: For optimal staining, the sections should be pretreated with formic acid (70%) for 10-30 minutes.

Incubation: For optimal staining, the primary antibody should be incubated overnight at 4°C.

- **ELISA:** 1:500-1:50,000

Storage: Do not store antibody diluted below 50 ug/mL in the absence of protein (i.e.: add 2% bovine serum albumin). It is suggested that the total volume is aliquoted upon initial thaw. The antibody is best stored undiluted at -20°C to -70°C.

References: Parvathy S, Davies P, Haroutunian V, Purohit DP, Davis KL, Mohs RC, Park H, Moran TM, Chan JY, Buxbaum JD. Correlation between Abeta₄₀-, Abeta₄₂-, and Abeta₄₃-containing amyloid plaques and cognitive decline. *Arch Neurol* 58:2025-2032, 2001.

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

Related Products:

- Beta Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Biotin Labeled (SIGNET)
Catalog Number SIG-39141

Product Revision Date: 9/11/2009