

Beta Amyloid 1-42 (12F4) Monoclonal Antibody, Purified (SIGNET)
Catalog Number: SIG-39142
Available Size: 0.2 mL, 0.5 mL, 1.0 mL

Description: Monoclonal Antibody against Beta Amyloid (A β) 1-42 (12F4)
(Formerly Signet Catalog Nos.: 9142-02, 9142-05, 9142-10)

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

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

Clone: 12F4

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

Host: Mouse

IsoType: IgG1

Species Reactivity: Human, Mouse

[Ab]: 0.5 mg/mL

Specificity: 12F4 is reactive to the c-terminus of beta amyloid and is specific for the isoform ending at the 42nd amino acid. This antibody was purified via Protein-G chromatography.

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:** 1:100-1:2,000

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

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.

Product Revision Date: 9/11/2009