

**Beta Amyloid Recombinant Peptide (1-40), Ultra Pure, TFA**  
**Catalog Number: RPEP-579P**  
**Available Size: 1.0 mg**

prior written approval.

**Related Products:**

- Beta Amyloid (A $\beta$ ) Monoclonal Antibody  
Catalog Number MMS-532R

**Description:** Beta Amyloid (1-40) Recombinant Ultra Pure Peptide

**Product Revision Date:** 9/11/2009

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

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

**Form:** Lyophilized Powder

**Specificity:** 99% Pure beta-Amyloid Peptide

Sequence: H-Asp-Ala-Glu-Phe-Arg-His-Asp-Ser-Gly-Tyr-Glu-Val-His-His-Gln-Lys-Leu-Val-Phe-Phe-Ala-Glu-Asp-Val-Gly-Ser-Asn-Lys-Gly-Ala-Ile-Ile-Gly-Leu-Met-Val-Gly-Gly-Val-Val-OH

**Molecular Weight:** 4330.3 Da by Mass Spec Analysis (theoretical 4329.86)

**Suggested Working Dilution:** Resuspend in 1 % NH<sub>4</sub>OH, at a concentration of 1 mg/mL. Sonicate for 30 seconds to 1 minute after it has gone into solution. To bring it into your buffer: After resuspension, add 5x or 10x buffer stock (PBS, TBS) and water, to bring to 1x buffer.

**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.

**Warranty/Conditions:** Covance products may not be resold or modified for resale without