

Beta Amyloid Recombinant Peptide (1-42), Ultra Pure, TFA
Catalog Number: RPEP-590P
Available Size: 1.0 mg

Description: Beta-Amyloid (1-42) Recombinant Ultra Pure Peptide

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-Ile-Ala-OH

Molecular Weight: 4514.5 Da by Mass Spec Analysis (theoretical 4514.10)

Suggested Working Dilution: Resuspend in 1 % NH₄OH, at a concentration of 1 mg/mL. Sonicate for 30 seconds 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

prior written approval.

Related Products:

- Beta Amyloid Scrambled Recombinant Peptide (1-42)
Catalog Number RPEP-598P
- Beta Amyloid (A β) Monoclonal Antibody
Catalog Number MMS-532R

Product Revision Date: 9/11/2009