

## **Product Data Sheet**

Product Name:  $\beta$ -Amyloid (4-42)

Catalog Number: AS-29908-1 (1 mg) Lot Number: See label on vial

Sequence: H-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 (3-letter code)

FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA(1-letter code)

Molecular Weight: 4198.8

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Initially, dissolve a small amount of  $\beta$ -Amyloid (4-42) peptide in water. If peptide does not dissolve, then use DMSO.

Storage:  $\beta$ -Amyloid (4-42) peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at  $-20^{\circ}$ C or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}$ C or lower.

Reference: Lalowski, M. et al. J. Biol. Chem. 271, 33623 (1996).

For Research Use Only