



## Product Data Sheet

**Product Name:**  $\beta$ -Amyloid (9-42)  
**Catalog Number:** AS-60084-1 (1 mg) **Lot Number:** See label on vial

**Sequence:** H-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)  
GYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code)

**Molecular Weight:** 3556.2

**% Peak Area by HPLC:**  $\geq 95$

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Reconstitute by adding 100  $\mu$ l 1%  $\text{NH}_4\text{OH}$  to 1 mg  $\beta$ -Amyloid (9-42) peptide. Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at  $-20^\circ\text{C}$ .

**Storage:**  $\beta$ -Amyloid (9-42) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^\circ\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^\circ\text{C}$  or lower.

**Additional Information:** *Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.*

Two wash steps, as described above, followed by 2 additional washes with 50 mmol/L HEPES (pH 7.0) completed the reaction. After the arrays had dried, a 20% saturated solution of  $\alpha$ -cyano-4-hydroxycinnamic acid (CIPHERGEN Biosystems) in 5 mL/L trifluoroacetic acid–500 mL/L acetonitrile–495 mL/L water was applied to each spot. Mass analysis was performed on a ProteinChip reader (Model PBS II; CIPHERGEN). For calibration purposes, 7 fmol of  $\text{A}\beta_{9-42}$  peptide (AnaSpec) and 6 fmol of bovine insulin (CIPHERGEN) were applied and used for data calibration-  
[Vanderstichele, H. et al. \*Clin. Chem.\* \*\*51\*\*, 1650 \(2005\).](#)

**Published Citations:**

Vanderstichele, H. et al. *Clin. Chem.* **51**, 1650 (2005).

*For Research Use Only*