



Product Data Sheet

Product Name: β -Amyloid (9-42)
Catalog Number: 60084-01 (0.1 mg) Lot Number: See label on vial
60084-1 (1 mg)
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: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 100 μ l 1% NH_4OH to 1 mg β -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°C .

Storage: β -Amyloid (9-42) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°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 α -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).

Related Products:

| Name | Cat # | Size |
|--|----------|--------|
| β -Amyloid (9-42)-Lys(Biotin)-NH ₂ NEW | 62462 | 0.5 mg |
| GYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA-K(Biotin)-NH ₂ | | |
| β -Amyloid (8-42) | 60085-01 | 0.1 mg |
| SGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA | 60085-1 | 1 mg |

For Research Use Only