



Product Data Sheet

Product Name: β -Amyloid (1-42), sodium salt

Catalog Number: AS-60883-01 (0.1 mg) Lot Number: See label on vial
AS-60883 (1.0 mg)

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, Na Salt (3-letter code)
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA, Na Salt (1-letter code)

Molecular Weight: 4514.1 • 23

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use 1.0% NH₄OH as the solvent, followed by buffer (i.e. 1X PBS). Add 1.0% NH₄OH directly to the lyophilized peptide powder (add 35-40 μ l to 0.5 mg peptide or 70-80 μ l to 1 mg peptide). The peptide cannot be stored long term in 1.0% NH₄OH, and it is therefore important to immediately dilute this solution with 1X PBS or other buffer to a concentration of approximately 1mg/mL or less. Gently vortex to mix.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at -20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at -20°C or lower. Do not freeze thaw.

Description: This peptide prepared by neutralizing the TFA salt form of β -Amyloid (1-42) with a dilute sodium hydroxide solution has superior solubility and fibrillogenesis properties, and the fibrils are equally neurotoxic. Ref: Fezoui, Y. et al. *Amyloid* 7, 166 (2000).

For Research Use Only