



Product Data Sheet

Product Name: β -Amyloid (1-42)-Lys(Biotin)-NH₂
Catalog Number: AS-61484-01 (0.1 mg) **Lot Number:** See label on vial

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-Lys(Biotin)-NH₂ (3-letter code)
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA
-K(Biotin)-NH₂ (1-letter code)

Molecular Weight: 4867.6

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use 1.0% NH₄OH as the solvent, followed by buffer (i.e. 1XPBS). 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: The difference between this recently released and the previous product (Cat# 23528-01) is the change at the C terminus from a carboxylic acid to an amide. The biological activity should remain the same. Elbaum, D. et al. *Biochem. Biophys. Res. Commun.* **267**, 733 (2000).

For Research Use Only