

Product Data Sheet

Product Name: β-Amyloid (1-40)-Lys(Biotin)-NH₂

Catalog Number: AS-61483-05 (0.5 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-Lys(Biotin)-NH₂ (3-letter code)

DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-

NH₂ (1-letter code)

Molecular Weight: 4685.4

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 35-40 μ l 1%NH₄OH to 0.5 mg β-Amyloid (1-40)-Lys(Biotin)-NH₂. 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 -20C.

Storage: β -Amyloid (1-40)-Lys(Biotin)-NH₂ 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.

Description: The difference between this recently released and the previous product (Cat# 23515) is the change at the C terminus from a carboxylic acid to an amide. The biological activity should remain the same.

For Research Use Only