



Product Data Sheet

Product Name: Biotin-LC- β -Amyloid (1-40)

Catalog Number: AS-24645-01(0.1 mg) Lot Number: See label on vial
AS-24648 (0.5 mg)
AS-24645 (1 mg)

Sequence: Biotin-LC-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-OH (3-letter code)
Biotin-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV (1-letter code)

Molecular Weight: 4670.3

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 50 μ l 1% NH_4OH to 0.5 mg Biotin-LC- β -Amyloid (1-40) 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: Biotin-LC- β -Amyloid (1-40) 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: This biotinylated $\text{A}\beta$ (1-40) contains a 6-carbon long chain (LC) to provide more accessibility for avidin attachment. Ref: Saito, Y. et al. *Proc. Natl. Acad. Sci. USA* **92**, 10227 (1995).

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.*

5.6 μ g biotinylated (long-chain) $\text{A}\beta$ 1-40 peptide (AnaSpec Incorporated, San Jose, CA) was reacted overnight at 4 degrees with 22.5 μ g streptavidin (Zymed, S. San Francisco, CA). This material was added to 7.5 μ g biotinylated bovine papillomavirus (BPV) L1-VLPs by incubation at 4 degrees overnight to generate $\text{A}\beta$ -conjugated VLPs. For immunization, 130 μ l complete Freund's adjuvant was added to 170 μ l VLP preparation containing 5.6 μ g $\text{A}\beta$, then emulsified and injected as 300 μ l into each mouse-[Li, Q. et al. BMC Neurosci. 5,21 \(2004\).](#)

Published Citations:

Li, Q. et al. *BMC Neurosci.* **5**, 21 (2004).

Related Products:

Name	Cat #	Size
Biotin- β -Amyloid (1-40) (Biotin-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-23512-01 AS-23512	0.1 mg 1 mg
β -Amyloid (1-40)-Lys(Biotin-LC) (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin-LC))	AS-23518-01 AS-23517	0.1 mg 0.5 mg
β -Amyloid (1-40)-Lys(Biotin)-NH ₂ (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂)	AS-61483-01 AS-61483-05	0.1 mg 0.5 mg
β -Amyloid (1-40)-Lys(Biotin)-NH ₂ , mouse, rat (DAEFGHDSGFVRRHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂)	AS-63356	0.1 mg
β -Amyloid (1-40)-Lys(LC-biotin)-NH ₂ , FAM-labeled (5-FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(LC-BIOTIN)-NH ₂)	AS-61962-01	0.1 mg
β -Amyloid (1-40) Binding Peptide, Biotin-labeled (Biotin-DWGKGGRWRLWPGASGKTEA)	AS-62427	1 mg
β -Amyloid (1-40) (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-24235 AS-24236	0.5 mg 1 mg
Biotin-LC- β -Amyloid (1-42) (Biotin-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-24641-01 AS-24640	0.1 mg 0.5 mg

For Research Use Only