

Product Name:Biotin-LC-β-Amyloid (1-42)

Catalog Number:AS-24641-01 (0.1 mg)Lot Number: See label on vialAS-24640(0.5 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-Ile-Ala-OH (3-letter code) Biotin-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code)

Molecular Weight: 4854.6

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 60-70 μ I 1%NH₄OH to 0.5 mg Biotin-LC- β -Amyloid (1-42). 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: Biotin-LC- β -Amyloid (1-42) 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.

Description: This biotinylated β -Amyloid (1-42) contains a 6-carbon long chain (LC) to provide more accessibility for avidin attachment.

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.

The detergent extracts and cytosol were divided into 4-ml aliquots and incubated in 17 x 100 mm polypropylene tubes (Falcon 2059) with 10 µg biotin-LC- -amyloid 1-42 (Anaspec) for 1h at room temperature-Nelson, TJ. & DL Aklon, *J. Biol. Chem.* 10.1074/jbc.M705558200 (2007).

Published Citations:

Nelson, TJ. & DL Aklon, J. Biol. Chem. 10.1074/jbc.M705558200 (2007).

For Research Use Only