



Product Data Sheet

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

Catalog Number: AS-23524-01 (0.1 mg) Lot Number: See label on vial
AS-23523-05 (0.5 mg)

Sequence: Biotin-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-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code)

Molecular Weight: 4740.7

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 30-50 μ l 1% NH_4OH to 0.5 mg Biotin- 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 -20°C .

Storage: Biotin- 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.

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.*

Under pentobarbital (50 mg/kg, i.p.) anesthesia, biotin-labeled A β (1–42) (100 μ M; AnaSpec, San Jose, CA) was infused at a rate of 0.66 μ l/min for 60 min into the lateral ventricle (coordinates relative to bregma: anterior–posterior 0.0-mm, lateral 2.0-mm, and 2.3-mm depth) of 4-month-old male L-PGDS^{-/-} mice and human L-PGDS-Tg mice-[Kanekiyo, T. et al. Proc. Natl. Acad. Sci. USA 104, 6412 \(2007\).](#)

Published Citations:

Le Vine III, H. *Anal. Biochem.* **356**, 265 (2006).
Kanekiyo, T. et al. *Proc. Natl. Acad. Sci. USA* **104**, 6412 (2007).

For Research Use Only