

## **Product Data Sheet**

Product Name: Biotin-β-Amyloid (1-40)

Catalog Number: AS-23512-01 (0.1 mg) Lot Number: See label on vial

AS-23512 (1 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-OH (3-letter code)

Biotin-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV

(1-letter code)

Molecular Weight: 4557.2 % Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 50-70 µl 1%NH<sub>4</sub>OH to

0.5 mg Biotin-β-Amyloid 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 -20C.

Storage: Biotin-β-Amyloid 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.

Reference: 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.

Biotin-A $\beta$ (1–40) were commercially available from AnaSpec, Inc. (San Jose, CA, USA). When stained with FITC-streptavidine, IMR-32 cells (1 × 10<sup>6</sup>) were incubated on ice for 30 min with biotin-A $\beta$ (1–40) (5  $\mu$ m) and washed three times with PBS containing 0.1% (v/v) BSA and 0.1% (v/v) NaN<sub>3</sub>. The cells were incubated on ice for 30 min together with FITC-streptavidin (1/100; Molecular Probes, Inc., Eugene, OR, USA). The cells were then washed three times and analyzed by an Epics Elite Flow Cytometer (Beckman Coulter, Fullerton, CA, USA)- Inaba S. et al. *J. Pept. Res.* **65**, 485 (2005).

## **Published Citations:**

Inaba S. et al. J. Pept. Res. 65, 485 (2005).

## **Related Products:**

Name	Cat #	Size
Biotin-LC-beta-Amyloid (1-40)	AS-24645-01	0.1 mg
(Biotin-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-24648 AS-24645	0.5 mg 1 mg
Beta-Amyloid (1-40)-Lys(Biotin-LC)	AS-23518-01	0.1 mg
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin-LC)	AS-23517	0.5 mg
Beta-Amyloid (1-40) Binding Peptide, Biotin-labeled	10.00407	4
(Biotin-DWGKGGRWRLWPGASGKTEA)	AS-62427	1 mg
Beta-Amyloid (1-40)-Lys(Biotin)-NH2	AS-61483-01	0.1 mg
(DAEFRHDSGYÈVHHQKLVFFAÉDVGSNKGAIIGLMVGGVV-K(Biotin)-NH2)	AS-61483-05	0.5 mg
Beta-Amyloid (1-40)-Lys(LC-biotin)-NH2, FAM-labeled, Ex/Em=490 nm/ 520 nm		

For Research Use Only

AS-61962-01

0.1 mg

(5-FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(LC-BIOTIN)-NH2)