

## **Product Data Sheet**

Product Name:  $\beta$ -Amyloid (1-40), mouse, rat

Catalog Number: AS-25380 (0.5 mg) Lot Number: See label on vial

AS-25230 (1 mg)

Sequence: H-Asp-Ala-Glu-Phe-Gly-His-Asp-Ser-Gly-Phe-Glu-Val-Arg-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)

DAEFGHDSGFEVRHQKLVFFAEDVGSNKGAIIGLMVGGVV

(1-letter code)

Molecular Weight: 4234.8

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 60-70 µl 1%NH<sub>4</sub>OH to

0.5 mg  $\beta$ -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 -20C.

Storage: 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.

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.

Forty-residue human or rodent  $\beta$ -Amyloid peptide and reverse human peptide were obtained from AnaSpec, Inc. Peptides were dissolved and stored in hexafluoro-isopropyl alcohol at 1 mg/ml. Samples were lyophilized by pervasion with nitrogen and then resuspended in PBS and immediately frozen as disaggregated peptide. To achieve consistent aggregation, the peptides were resuspended in cell growth medium and divided into 0.13-ml aliquots in a 96-well plate. The plate was shaken at 500 rpm for 5 h. Samples were then combined and normalized to a final A $\beta$  concentration of 50  $\mu$ M and immediately frozen-Kajkowski, EM. et al. *J. Biol. Chem.* 276, 18748 (2001).

Rat A $\beta$ 1–40 and human A $\beta$ 1–28 were purchased from AnaSpec (San Jose, CA). Synthetic A $\beta$  peptide solutions were dissolved in doubly deionized water at a concentration of 0.5–1.0 mg/mL, sonicated for 3 min, and then centrifuged for 20 min at 10000g, and the supernatant (stock A $\beta$ ) was used on the day of the experiment. The concentrations of stock A $\beta$  peptides were determined by the spectrophotometric absorbance at 214 nm or by a Micro BCA protein assay (Pierce, Rockford, IL)- Atwood, CS. et al. *Biochem.* 43, 560 (2004).

## **Published Citations:**

Sawamura, N. et al. *J. Biol. Chem.* **275**, 27901 (2000). Kajkowski, EM. et al. *J. Biol. Chem.* **276**, 18748 (2001). Atwood, CS. et al. *Biochem.* **43**, 560 (2004). Howlett, DR. et al. *Histol. Histophathol.* **23**, 67 (2008).

## Related Products:

Name	Cat #	Size
β-Amyloid (1-40)-Lys(Biotin)-NH2, mouse, rat (DAEFGHDSGFEVRHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH	AS-63356 H2)	0.1 mg
β-Amyloid Peptide (1-42), mouse, rat (DAEFGHDSGFEVRHQKLVFFAEDVGSNKGAIIGLMVGGVVIA)	AS-25381 AS-25231	0.5 mg 1 mg
β-Amyloid (1-38), mouse, rat (DAEFGHDSGFEVRHQKLVFFAEDVGSNKGAIIGLMVGG)	AS-62476	1 mg

For Research Use Only