



Product Data Sheet

Product Name: β -Amyloid (1-40), FAM-Labeled, Human

Catalog Number: AS-23514-01 (0.1 mg) Lot Number: See label on vial
AS-23513-05 (0.5 mg)

Sequence: FAM-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)
FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV
(1-letter code)

Molecular Weight: 4689.2

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized yellow powder

Peptide Reconstitution: Reconstitute by adding 35-50 μ l 1% NH_4OH to 0.5 mg β -Amyloid (1-40). 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: 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 is a fluorescent (FAM)-labeled β -Amyloid (1-40), Abs/Em=494/521 nm. FAM is preferred over FITC because of its photo- and chemical stability.

Published Citations:

Wakabayashi, M. et al. *Biochem. Biophys. Res. Comm.* **328**, 1019 (2005).
Wakabayashi, M. and K. Matsuzaki, *J. Mol. Biol.* **371**, 924 (2007).

For Research Use Only