



Product Data Sheet

Product Name: β -Amyloid (1-38)

Catalog Number: AS-24233 (0.5 mg) Lot Number: See label on vial
AS-24234 (1 mg)

Sequence: H-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-OH (3-letter code)
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGG (1-letter code)

Molecular Weight: 4132.6

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 30-50 μ l 1%NH₄OH to 0.5 mg β -Amyloid (1-38) 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 (1-38) 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.

Published Citations:

Chauhan, A. et al. *Neurochem. Res.* **25**, 1573 (2000).

Related Products:

Name	Cat #	Size
Beta-Amyloid (1-38), mouse, rat (DAEFGHDSGFVRRHQLVFFAEDVGSNKGAIIGLMVGG)	AS-62476	1 mg
Beta-Amyloid (1-37) (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGG)	AS-60251-05	0.5 mg
	AS-60251-1	1 mg

For Research Use Only