

Product Data Sheet

Product Name: β -Amyloid (3-42)

Catalog Number: AS-63715-01 (0.1 mg) Lot Number: See label on vial

AS-63715-1 (1 mg)

Sequence: H-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)

EFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA

(1-letter code)

Molecular Weight: 4328.9

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using DMSO, reconstitute by adding 100 μl to 1 mg β-Amyloid (3-42)

peptide. This peptide is also soluble in 1%NH₄OH, but not as clear as DMSO.

Storage: β -Amyloid (3-42) 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.

Related Products:

..

Name	Cat #	Size
[Pyr ³]- β-Amyloid (3-42)	AS-29907-01	0.1 mg
(Pyr-FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA)	AS-29907-1	1 mg
[Pyr³]- β-Amyloid (3-40)	AS-29906-01	0.1 mg
(Pyr-FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-29906-1	1 mg
β-Amyloid (3-40)	AS-61029	0.1 mg

For Research Use Only