



Product Data Sheet

Product Name:	[Pyr ¹¹]- β -Amyloid (11-40)	
Catalog Number:	AS-29904-1 (1 mg)	Lot Number: See label on vial
Sequence:	Pyr-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) Pyr-VHHQKLFFAEDVGSNKGAIIGLMVGGVV (1-letter code)	
Molecular Weight:	3133.7	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Initially, dissolve a small amount of [Pyr¹¹]- β -Amyloid (11-40) peptide in water. If peptide is insoluble in water, use DMSO.

Storage: [Pyr¹¹]- β -Amyloid (11-40) 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.

Published Citations:

Piccini, A. et al., *J. Biol. Chem.* **280**, 34186 (2005).

Related Products:

Name	Cat #	Size
β -Amyloid (11-40) EVHHQKLFFAEDVGSNKGAIIGLMVGGVV	AS-60017-1	1 mg
[Pyr ¹¹]- β -Amyloid (11-42) Pyr-VHHQKLFFAEDVGSNKGAIIGLMVGGVVIA	AS-29903-01 AS-29903-1	0.1 mg 1 mg
[Pyr ³]- β -Amyloid (3-42) Pyr-FRHDSGYEVHHQKLFFAEDVGSNKGAIIGLMVGGVVIA	AS-29907-01 AS-29907-1	0.1 mg 1 mg
[Pyr ³]- β -Amyloid (3-40) Pyr-FRHDSGYEVHHQKLFFAEDVGSNKGAIIGLMVGGVV	AS-29906-01 AS-29906-1	0.1 mg 1 mg

For Research Use Only