



Product Data Sheet

| | | |
|-----------------------------|--|-------------------------------|
| Product Name: | [Pyr ¹¹]- β -Amyloid (11-42) | |
| Catalog Number: | AS-29903-01 (0.1 mg) AS-29903-1 (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-Ile-Ala-OH (3-letter code) Pyr-VHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code) | |
| Molecular Weight: | 3318.9 | |
| % Peak Area by HPLC: | ≥ 95 | |
| Appearance: | Lyophilized white powder | |

Peptide Reconstitution: Using DMSO, reconstitute by adding 30-60 μ l to 0.5 mg or 100 μ l to 1 mg [Pyr¹¹]- β -Amyloid (11-42) peptide. Dilute this stock solution with working buffer, such as PBS or other buffers.

Storage: [Pyr¹¹]- β -Amyloid (11-42) 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.

Description: A β (1-42), a major component of amyloid plaques, accumulates in neurons of Alzheimer's disease brains. Biochemical analysis of the amyloid peptides isolated from Alzheimer's disease brain indicates that A β (1-42) is the principal species associated with senile plaque amyloids, while A β (1-40) is more abundant in cerebro-vascular amyloid deposits. Ref: Nagele, R. et al. *Neurosci.* **110**, 199 (2002); Garzon-Rodriguez, W. et al. *J. Biol. Chem.* **272**, 21037 (1997).

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.

Synthetic peptides corresponding to the three major soluble A β species detected by immunoblotting (A β 1-42, py3-42, and py11-42; Anaspec) were dissolved with 1,1,1,3,3,3-hexafluoro-2-propanol (HFIP, Sigma) to produce a uniform, nonaggregated field of A β . The studies on toxicity and aggregation are all based on different lots of peptides to avoid the possibility that results could depend on a variation in the synthetic peptide batches (A β 1-42 lots 26048 and 31475, A β py3-42 lots 25479 and 28058, A β py11-42 lot 28001; Anaspec). The three synthetic A β peptides were suspended in phosphate-buffered saline at a ratio corresponding to the composition of soluble A β detected in AD (A β 1-42, 36%; A β py3-42, 48%; A β py11-42, 16%) and NA (A β 1-42, 50%; A β py3-42, 29%; A β py11-42, 21%) and kept for 24-48 h at room temperature at a final concentration of 10 μ M, pH 7.6, for subsequent analyses. For experiments in immunoprecipitation and immunoblotting, peptides were kept at 37 $^{\circ}\text{C}$ for 24 h-[Piccini, A. et al. J. Biol. Chem. 280, 34186 \(2005\).](#)

Published Citations:

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

Related Products:

| Name | Cat # | Size |
|---|---------------------------|----------------|
| β -Amyloid (11-42) (EVHHQKLFFAEDVGSNKGAIIGLMVGGVIA) | AS-63317 | 1 mg |
| β -Amyloid (11-42), HiLyte Fluor™ 488-labeled (HiLyte Fluor™ 488-EVHHQKLFFAEDVGSNKGAIIGLMVGGVIA) | AS-63327 | 0.1 mg |
| [Pyr ¹¹]- β -Amyloid (11-40) (Pyr-VHHQKLFFAEDVGSNKGAIIGLMVGGVV) | AS-29904-1 | 1 mg |
| [Pyr ³]- β -Amyloid (3-42) (Pyr-FRHDSGYEVHHQKLFFAEDVGSNKGAIIGLMVGGVIA) | 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