



## Product Data Sheet

<b>Product Name:</b>	[Pyr <sup>11</sup> ]- $\beta$ -Amyloid (11-42)	
<b>Catalog Number:</b>	AS-29903-01 (0.1 mg) AS-29903-1 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	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)	
<b>Molecular Weight:</b>	3318.9	
<b>% Peak Area by HPLC:</b>	≥ 95	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Using DMSO, reconstitute by adding 30-60  $\mu$ l to 0.5 mg or 100  $\mu$ l to 1 mg [Pyr<sup>11</sup>]- $\beta$ -Amyloid (11-42) peptide. Dilute this stock solution with working buffer, such as PBS or other buffers.

**Storage:** [Pyr<sup>11</sup>]- $\beta$ -Amyloid (11-42) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** A $\beta$  (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 $\beta$  (1-42) is the principal species associated with senile plaque amyloids, while A $\beta$  (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 $\beta$  species detected by immunoblotting (A $\beta$ 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 $\beta$ . 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 $\beta$ 1-42 lots 26048 and 31475, A $\beta$ py3-42 lots 25479 and 28058, A $\beta$  py11-42 lot 28001; Anaspec). The three synthetic A $\beta$  peptides were suspended in phosphate-buffered saline at a ratio corresponding to the composition of soluble A $\beta$  detected in AD (A $\beta$ 1-42, 36%; A $\beta$  py3-42, 48%; A $\beta$  py11-42, 16%) and NA (A $\beta$  1-42, 50%; A $\beta$  py3-42, 29%; A $\beta$  py11-42, 21%) and kept for 24-48 h at room temperature at a final concentration of 10  $\mu$ 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:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
$\beta$ -Amyloid (11-42) (EVHHQKLVFFAEDVGSNKGAIIGLMVGGVIA)	AS-63317	1 mg
$\beta$ -Amyloid (11-42), HiLyte Fluor™ 488-labeled (HiLyte Fluor™ 488-EVHHQKLVFFAEDVGSNKGAIIGLMVGGVIA)	AS-63327	0.1 mg
[Pyr <sup>11</sup> ]- $\beta$ -Amyloid (11-40) (Pyr-VHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-29904-1	1 mg
[Pyr <sup>3</sup> ]- $\beta$ -Amyloid (3-42) (Pyr-FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVIA)	AS-29907-01 AS-29907-1	0.1 mg 1 mg
[Pyr <sup>3</sup> ]- $\beta$ -Amyloid (3-40) (Pyr-FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV)	AS-29906-01 AS-29906-1	0.1 mg 1 mg

*For Research Use Only*