



Product Data Sheet

Product Name: [Pyr3]- β -Amyloid (3-42)
Catalog Number: 29907-01 (0.1 mg) **Lot Number:** See label on vial
29907-1 (1 mg)
Sequence: Pyr-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)
Pyr-FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter
code)

Molecular Weight: 4309.9

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use 1.0% NH₄OH as the solvent, followed by buffer (i.e. 1XPBS). Add 1.0% NH₄OH directly to the lyophilized peptide powder (add 35-40 μ l to 0.5 mg peptide or 70-80 μ l to 1 mg peptide). The peptide cannot be stored long term in 1.0% NH₄OH, and it is therefore important to immediately dilute this solution with 1X PBS or other buffer to a concentration of approximately 1mg/mL or less. Gently vortex to mix.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at -20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at -20°C or lower. Do not freeze thaw.

Description: This is one of the predominant Amyloid peptide structures deposited in human brain of Alzheimer's disease and Down's syndrome patients. Ref: Russo, C. et al. *Neurobiol Nature* **405**, 531 (2000).

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-[Piccini, A. et al. J Biol Chem 280, 34186 \(2005\)](#).

Published Citations:

[Piccini, A. et al. J Biol Chem 280, 34186 \(2005\)](#).

Related Products:



Click Alzheimer's Research Brochure to view the full listing.

For Research Use Only