



Product Data Sheet

Product Name: β -Amyloid (25-35)
Catalog Number: AS-24227 (1 mg) Lot Number: See label on vial
AS-24228 (5 mg)
Sequence: H-Gly-Ser-Asn-Lys-Gly-Ala-Ile-Ile-Gly-Leu-Met-OH (3-letter code)
GSNKGAIIGLM (1-letter code)
Molecular Weight: 1061.3
% Peak Area by HPLC: ≥ 95
Appearance: Lyophilized white powder

Peptide Reconstitution: β -Amyloid (25-35) peptide is freely soluble in H₂O.

Storage: β -Amyloid (25-35) 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 (25-35) is the main factor responsible for A β neurotoxic effects. Ref: Carvalho, K. et al. *Braz. J. Med. Biol. Res.* **3**, 1153 (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.*

The A β 25–35 peptides used in this study were purchased from AnaSpec (San Jose, CA). Purity was certified by high-performance liquid chromatography–mass spectrometry for each of the peptide. The peptides were resuspended in sterile double-deionized water, aliquoted at 5 mg/ml, and kept at -20°C -[Hashioka, S. et al. *Free Radical Bio. Med.* **42**, 945 \(2007\).](#)

Synthetic A β 25–35, obtained from AnaSpec, was dissolved in deionized distilled water at a concentration of 2.5 mM and stored at -80°C . Previous to the experiments the stock solution was diluted to the desired concentrations, maintained for 3 h at room temperature and then added to the culture medium. After treatment with the doses of A β and times indicated in the text, the cells were analyzed by optical and fluorescence microscopy to evaluate cell viability, or collected and saved for posterior RNA extraction and microarray analysis- [Martínez, T. and A. Pascual *Brain Res. Bull.* **72**, 225 \(2007\).](#)

Synthetic β -Amyloid peptides 25-35 and 35-25 (A β 25-35, A β 35-25) and FAM-labeled β -Amyloid peptide 1-40 (FAM-A β 1-40) were purchased from AnaSpec, Inc. (San Jose, CA). Aggregated A β 25-35 and A β 1-40 were prepared at 4°C for 60 h, and were then incubated at 37°C for 48 h. Oligomeric, fibrillar, and aggregated A β 1-42 and aggregated FAM-A β 1-40 were prepared as described. Briefly, A β peptides were dissolved to a final concentration of 1 mM in hexafluoroisopropanol (Sigma–Aldrich). Hexafluoroisopropanol was removed using a speed vacuum. A β aliquot was resuspended in a solution containing 10 mM HCl and 150 mM NaCl, and then incubated at 37°C for 24 h to form aggregated A β . A β aliquot was suspended in 10 mM HCl to a final concentration 100 μM and incubated at 37°C for 24 h to form fibrillar A β - [Huang, W-C. et al. *Neurosci. Res.* **63**, 280 \(2009\).](#)

Published Citations:

Yatin, SM. et al. *J. Mol. Neurosci.* **11**, 183. (1998).
Kawahara, M. and Y. Kuroda *Cell. Mol. Neurobio.* **21**, 1 (2001).
Egashira, N. et al. *Japanese J. Pharmacol.* **90**, 321 (2002).
Pu, F. et al. *J. Health Sci.* **51**, 636 (2005).
Hashioka, S. et al. *Free Radical Bio. Med.* **42**, 945 (2007).
Martínez, T. and A. Pascual *Brain Res. Bull.* **72**, 225 (2007).
Nelson, T.J. and DL. Alkon *J. Biol. Chem.* **282**, 31238 (2007).
Seyb, Kl. et al. *J. Biomol. Screen.* **13**, 870 (2008).
Huang, W-C. et al. *Neurosci. Res.* **63**, 280 (2009).

Related Products:

Name	Cat #	Size
β -Amyloid (25-35) • HCl (GSNKGAIIGLM • HCl)	AS-23212	5 mg
Biotin- β -Amyloid (25-35) (Biotin-GSNKGAIIGLM)	AS-62451	1 mg
β -Amyloid (25-35), HiLyte Fluor™ 488-labeled (HiLyte Fluor™ 488-GSNKGAIIGLM)	AS-63308	0.1 mg

For Research Use Only