



Product Data Sheet

Product Name:	β -Amyloid (1-42) • HCl, Human	
Catalog Number:	AS-21791 (0.5 mg) AS-21793 (1.0 mg)	Lot Number: See label on vial
Sequence:	H-Asp-Ala-Glu-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 • HCl (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA • HCl (1-letter code)	
Molecular Weight:	4514.1 • 36.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	
Peptide Reconstitution:	Using basic buffer (1%NH ₄ OH), reconstitute by adding 50 μ l to 0.5 mg or 100 μ l to 1 mg β -Amyloid (1-42) • HCl peptide. Dilute this stock solution with working buffer, such as PBS or other buffers.	
Storage:	β -Amyloid (1-42) • HCl 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.	

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.*

β -Amyloid (1-42) • HCl was purchased from AnaSpec Inc. (San Jose, CA, USA). $A\beta$ • HCl (1-42) at a concentration of 500 μM was dissolved in 2 mM HCl and stored in small aliquots at -20°C . To generate fibrils, $A\beta$ • HCl (1-42) was diluted to 100 μM in water, NaCl and HEPES (final concentrations: 150 mM NaCl, 0.4 mM HCl and 20 mM HEPES) and then incubated at room temperature for 1 – 2 days. The peptide was vortexed and then diluted to desired final concentrations in artificial cerebrospinal fluid (ACSF) immediately before pretreatment in each experiment – [Nakagami, Y. and T. Oda, Japanese J. Pharma. 88, 223 \(2002\).](#)

β -Amyloid (1-42) • HCl was purchased from AnaSpec Inc. All of $A\beta$ • HCl (1-42) was dissolved at a concentration of 500 μM in phosphate-buffered saline (PBS), and stored in small aliquots at -20°C . RS-0406 was synthesized in our laboratory, and dissolved in dimethyl sulphoxide (DMSO) at a concentration of 30 mg ml^{-1} or in ethanol at a concentration of 5 mg ml^{-1} – [Nakagami, R. et al. Brit. J. Pharma. 137, 676 \(2002\).](#)

β -Amyloid (1-42) • HCl peptide was purchased from AnaSpec (San Jose, CA, USA). A 1 mg sample of peptide was dissolved in 200 μL hexafluoroisopropanol (HFIP) and aliquoted to obtain 0.1 mg stocks (handle HFIP in a chemical fume hood taking the necessary precautions) -

A β • HCl (1-42) was dissolved in sterile distilled water, and allowed to aggregate for 7d at 37 °C.
A β • HCl (1-42) injected daily for 7 d after cerebral ischemia – [Iwasaki, K. et al. *Bio. Pharma. Bull.* **30**, 698 \(2007\).](#)

Published Citations:

Kakimura, J-I. et al. *FASEB J.* **10**.1096/fj.01-0530fje (2002).
Nakagami, Y. and T. Oda, *Japanese J. Pharma.* **88**, 223 (2002).
Nakagami, Y. et al. *Eur. J. Pharma.* **457**, 11 (2002).
Nakagami, R. et al. *Brit. J. Pharma.* **137**, 676 (2002).
Kubo, T. et al. *Pharma. Toxicol.* **93**, 264 (2003).
Iwasaki, K. et al. *Neurotox. Res.* **6**, 299 (2004).
Hirata, K. et al. *J. Pharmacol. Exp. Ther.* **314**, 252 (2005).
Iwasaki, K. et al. *Brain Res.* **1097**, 216 (2006).
Iwasaki, K. et al. *Bio. Pharma. Bull.* **30**, 698 (2007).

Related Products:

Name	Cat #	Size
β -Amyloid (1-40) • HCl	AE2321F	0.5 mg
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV • HCl	AE20698	1 mg

For Research Use Only