

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-IIe-IIe-Gly- Leu-Met-Val-Gly-Gly-Val-Val-IIe-Ala-OH • HCI (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA • HCI (1-letter code)	
Molecular Weight:	4514.1 • 36.5	
Peptide Purity:	>95%	

Appearance: Lyophilized white powder

Peptide Reconstitution: Using basic buffer (1%NH₄OH), 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) • HCI was purchased from AnaSpec Inc. (San Jose, CA, USA). A β • HCI (1-42) at a concentration of 500 μ M was dissolved in 2 mM HCI and stored in small aliquots at -20 °C. To generate fibrils, A β • HCI (1-42)) was diluted to 100 μ M in water, NaCI and HEPES (final concentrations: 150 mM NaCI, 0.4 mM HCI 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 β • HCl (1-42) was dissolved at a concentration of 500 μ M 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 – <u>Nakagami, R. et al. *Brit. J. Pharma.* **137**, 676 (2002).</u>

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

A β • HCI (1-42) was dissolved in sterile distilled water, and allowed to aggregate for 7d at 37 °C. A β • HCI (1-42) injected daily for 7 d after cerebral ischemia – <u>Iwasaki, K. et al. *Bio. Pharma. Bull.*</u> **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) • HClÊ͡₽ັ{ æ}	ŒÙË2321F	0.5 { *
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV• HCl	ŒÙË20698	1 mg

For Research Use Only