

## **Product Data Sheet**

Product Name:	β-Amyloid (1-39)	
Catalog Number:	AS-24295 (0.5 mg) AS-24296 (1 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-OH (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGV (1-letter code)	
Molecular Weight:	4230.7	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution:  $\beta$ -Amyloid (1-39) peptide is freely soluble in basic buffer.

Storage:  $\beta$ -Amyloid (1-39) 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 number of  $A\beta$  protein variants, differing only at their carboxy terminus (1-39, 1-40, 1-42 and 1-43), are identified as the major components of the cerebral amyloid deposits in Alzheimer's disease. The length of the C-terminus is a critical determinant of the rate of amyloid formation ("kinetic solubility"), with only a minor effect on the thermodynamic solubility. Amyloid formation by the kinetically soluble peptides (e.g. 1-39) can be nucleated, or "seeded" by peptides which include the critical C-terminal residues (1-42, 26-42, 26-43, 34-42). Ref: Jarrett, JT. et al. Biochem. 32, 4693 (1993); Giacomelli, CE. and W. Norde, Macromol. Biosci. 5, 401 (2005).

## 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.

A peptide homologous to the first 39 residues of A $\beta$ ,  $\beta$ -(1-39), was purchased from AnaSpec, Inc. (San Jose, CA).  $\beta$ -(1-39), H2, or a mixture of  $\beta$ -(1-39) and H2 (1:2 molar ratio) were dissolved in 0.01 M phosphate buffer, pH 7.2. The final concentration was 0.5 mg/ml for each peptide. The solution was passed through a 0.45-µm Millipore filter to remove dust and then degassed- <u>Ghanta, J. et al. *J Biol Chem* **271**, 29525 (1996).</u>

Published Citations:

Ghanta, J. et al. *J Biol Chem* **271**, 29525 (1996). Good, TA. et al. *Biophys J* **70**, 296 (1999).. Pallitto, MM. et al. *Biochem* **38**, 3570 (1999).

For Research Use Only