



Product Data Sheet

Product Name:	β-Amyloid (40-1)	
Catalog Number:	AS-22817 (0.5 mg) AS-22818 (1 mg)	Lot Number: See label on vial
Sequence:	H-Val-Val-Gly-Gly-Val-Met-Leu-Gly-Ile-Ile-Ala-Gly-Lys-Asn-Ser-Gly-Val-Asp-Glu-Ala-Phe-Phe-Val-Leu-Lys-Gln-His-His-Val-Glu-Tyr-Gly-Ser-Asp-His-Arg-Phe-Glu-Ala-Asp-OH (3-letter code) VVGGMVLGIIAGKNSGVDEAFFVLKQHHVEYGSDHRFEAD (1-letter code)	
Molecular Weight:	4329.9	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: β-Amyloid (40-1) peptide is freely soluble in DMSO.

Storage: β-Amyloid (40-1) 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.

Aβ (1-40) (active peptide) and Aβ (40-1) (negative control with the reverse amino acid sequence) were purchased from Anaspec (San Jose, CA) and dissolved in sterile ultra-pure water obtained from ELGA Purelab Classic (Lowell, MA)-[Andras, IE. et al. Mol. Pharmacol. 10.1124/mol.107.042028 \(2008\).](#)

Published Citations:

Nakamura, S. et al. *Brain Res.* **912**, 128 (2001).
Kajkowski, EM. et al. *J. Biol. Chem.* **276**, 18748 (2001).
Ito, Y. et al. *Neurosci. Res.* **46**, 119 (2003).
Watanabe, N. et al. *The FASEB Journal.* **18**, 1013 (2004).
Andras, IE. et al. *Mol. Pharmacol.* 10.1124/mol.107.042028 (2008).

Related Products:

Name	Cat #	Size
β-Amyloid (1-40)	AS-24235	0.5 mg
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	AS-24236	1 mg
	AS-24236-5	5 mg

β -Amyloid (42-1)
AIVVGGVMLGIIAGKNSGVDEAFFVLKQHHVEYGSDHRFEAD

AS-27276-01 0.1 mg
AS-27275 0.5 mg
AS-27276 1 mg

For Research Use Only