



Product Data Sheet

Product Name:	Scrambled- β -Amyloid (1-40)	
Catalog Number:	AS-24625 (0.5 mg) AS-24626 (1 mg)	Lot Number: See label on vial
Sequence:	H-Ala-Glu-Gly-Asp-Ser-His-Val-Leu-Lys-Glu-Gly-Ala-Tyr-Met-Glu-Ile-Phe-Asp-Val-Gln-Gly-His-Val-Phe-Gly-Gly-Lys-Ile-Phe-Arg-Val-Val-Asp-Leu-Gly-Ser-His-Asn-Val-Ala-OH (3-letter code) AEGDSHVLKEGAYMEIFDVQGHVFGGKIFRVVDLGSNVA (1-letter code)	
Molecular Weight:	4329.9	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	
Peptide Reconstitution:	Scrambled-beta-Amyloid is freely soluble in basic buffer.	
Storage:	Scrambled-beta-Amyloid 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.	

Published Citations:

Ji, Y. et al. *J. Alzheimer's Dis.* **3**, 23 (2001).

Related Products:

Name	Cat #	Size
Scrambled- β -Amyloid (1-42)	AS-25382	0.5 mg
AIAEGDSHVLKEGAYMEIFDVQGHVFGGKIFRVVDLGSNVA	AS-25383	1 mg
β -Amyloid (1-42), Scrambled, 5-FAM labeled	AS-60892	0.1 mg
5-FAM-AIAEGDSHVLKEGAYMEIFDVQGHVFGGKIFRVVDLGSNVA		

For Research Use Only