



Product Data Sheet

Product Name:	β -Amyloid (42-1) • HCl	
Catalog Number:	60291-01 (0.1 mg) 60291-05 (0.5 mg)	Lot Number: See label on vial
Sequence:	H-Ala-Ile-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 • HCl (3-letter code) AIVVGGVMLGIIAGKNSGVDEAFFVLKQHHVEYGSDHRFEAD • HCl (1-letter code)	
Molecular Weight:	4515.1•36.5	
% Peak Area by HPLC:	≥ 95	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Reconstitute by adding 50-70 μ l 1%NH₄OH to 0.5 mg β -Amyloid (42-1) • HCl. Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at -20°C.

Storage: β -Amyloid (42-1) • HCl peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted A β peptide can be aliquoted and stored at -20°C or lower.

Ref: Ref: Hirata, K. et al. *J. Pharmacol. Exp. Ther.* 10.1124/jpet.105.083543 (2005).

Related Products:

Name	Cat #	Size
Beta - Amyloid (42 - 1)	27276-01	0.1 mg
(AIVVGGVMLGIIAGKNSGVDEAFFVLKQHHVEYGSDHRFEAD)	27275	0.5 mg
	27276	1 mg
Beta - Amyloid (40 - 1) • HCl	21795	0.5 mg
(VVGGVMLGIIAGKNSGVDEAFFVLKQHHVEYGSDHRFEAD • HCl)	21797	1 mg
Beta - Amyloid (40 - 1)	22817	0.5 mg
VVGGVMLGIIAGKNSGVDEAFFVLKQHHVEYGSDHRFEAD	22818	1 mg
DHL™ Fluorescent beta-Amyloid (1-42) Sampler Kit	72071	1 kit
DHL™ fluorescent beta-amyloid (1-40) sampler kit	72070	1 kit

For Research Use Only