



## Product Data Sheet

---

Product Name:	Beta - Amyloid (26 - 37)	
Catalog Number:	62920 (1mg)	Lot Number: See label on vial
Sequence:	H-Ser-Asn-Lys-Gly-Ala-Ile-Ile-Gly-Leu-Met-Val-Gly-OH (3-letter code) SNKGAIIGLMVG (1-letter code)	
Molecular Weight:	1159.4	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

**Peptide Reconstitution:** Initially try to dissolve a small amount of Beta-Amyloid (26-37) in water. If it does not dissolve, use DMSO.

**Storage:** Beta - Amyloid (26 - 37) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** This is amino acids 26 to 37 fragment of beta-amyloid peptide.

*For Research Use Only*