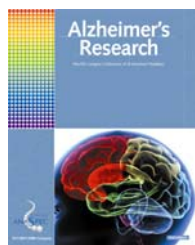




## Product Data Sheet

<b>Product Name:</b>	Beta-Amyloid (1-40) (S26C) Dimer	
<b>Catalog Number:</b>	64130-05 (0.5 mg) 64130-1 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	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-Cys-Asn-Lys-Gly-Ala-Ile-Ile-Gly-Leu-Met-Val-Gly-Gly-Val-Val-OH (dimer) (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGCNKGAIIGLMVGGVV (dimer) (1-letter code)	
<b>Molecular Weight:</b>	8689.8	
<b>% Peak Area by HPLC:</b>	≥ 90	
<b>Appearance:</b>	Lyophilized white powder	
<b>Peptide Reconstitution:</b>	Use fresh concentrated Anhydrous DMSO as the solvent. Do not use DMSO diluted with aqueous solvents, as this is not an effective solvent. Do not add aqueous solvent to the peptide prior to DMSO addition as this may prevent proper solubilization. Add concentrated DMSO directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5mg/mL to 1mg/mL. Gently vortex to mix	
<b>Storage:</b>	Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at –20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.	
<b>Related Products:</b>		



Click Alzheimer's Research Brochure to view the full listing.

***For Research Use Only***