

## Technical Data Sheet

<b>Product Name</b>	<b><i>Beta - Amyloid (13 - 28), Human</i></b> HHQKLVFFAEDVGSNK
<b>Size</b>	1 mg
<b>Catalog #</b>	AS-61802
<b>Purity</b>	% Peak Area By HPLC $\geq$ 95%
<b>Description</b>	This $\beta$ -Amyloid peptide 13 to 27 amino acid residues was used to study the kinetics of $\beta$ -amyloid formation.
<b>Storage</b>	-20°C
<b>References</b>	Inouye, H. and DA. Kirschner <i>Subcell Biochem.</i> <b>38</b> , 203 (2005).
<b>Molecular Weight</b>	1856.1
<b>Sequence (One-Letter Code)</b>	HHQKLVFFAEDVGSNK
<b>Sequence (Three-Letter Code)</b>	H - His - His - Gln - Lys - Leu - Val - Phe - Phe - Ala - Glu - Asp - Val - Gly - Ser - Asn - Lys - OH

*For Research Use Only*