

## Technical Data Sheet

<b>Product Name</b>	<b><i>[Lys22] - beta - Amyloid (1 - 40), Italian Mutation, Human</i></b> <i>DAEFRHDSGYEVHHQKLVFFAKDVGSNKGAIIGLMVGGVV</i>
<b>Size</b>	0.5 mg
<b>Catalog #</b>	AS-62147
<b>Purity</b>	% Peak Area By HPLC $\geq$ 95%
<b>Description</b>	This is amino acids 1 to 40 fragment of the beta-amyloid peptide with lysine substituted for glutamic acid at position 22, found in Italian families. The Italian mutation of beta-amyloid 1-40 (E22K) aggregates more rapidly and with more potent neurotoxicity than wild-type beta-amyloid 1-40. The formation of a salt bridge between Lys-22 and Asp-23 in the minor conformer might be a reason why E22K-beta-amyloid 40 is more pathogenic than wild-type beta-amyloid 40.
<b>Storage</b>	-20°C
<b>References</b>	Masuda, Y. et al. Bioorg. Med. Chem. 13, 6803 (2005).
<b>Molecular Weight</b>	4328.9
<b>Sequence (One-Letter Code)</b>	DAEFRHDSGYEVHHQKLVFFAKDVGSNKGAIIGLMVGGVV
<b>Sequence (Three-Letter Code)</b>	H - Asp - Ala - Glu - Phe - Arg - His - Asp - Ser - Gly - Tyr - Glu - Val - His - His - Gln - Lys - Leu - Val - Phe - Phe - Ala - Lys - Asp - Val - Gly - Ser - Asn - Lys - Gly - Ala - Ile - Ile - Gly - Leu - Met - Val - Gly - Gly - Val - Val - OH

*For Research Use Only*