

Technical Data Sheet

Product Name	<i>[Asn7] - beta - Amyloid (1 - 40), Tottori - Japanese Mutation, Human</i> <i>DAEFRHNSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV</i>
Size	0.5 mg
Catalog #	AS-63320
Purity	% Peak Area By HPLC \geq 95%
Detailed Information	This Beta-Amyloid 1-40 fragments has and Asn at position 7 instead of the normal Asp
Storage	-20°C
Molecular Weight	4328.9
Sequence (One-Letter Code)	DAEFRHNSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV
Sequence (Three-Letter Code)	H - Asp - Ala - Glu - Phe - Arg - His - Asn - Ser - Gly - Tyr - Glu - Val - His - His - Gln - Lys - Leu - Val - Phe - Phe - Ala - Glu - Asp - Val - Gly - Ser - Asn - Lys - Gly - Ala - Ile - Ile - Gly - Leu - Met - Val - Gly - Gly - Val - Val - OH

For Research Use Only