

Technical Data Sheet

Product Name	<i>[Asn23] - beta - Amyloid (1 - 40), Iowa Mutation, Human</i> <i>DAEFRHDSGYEVHHQKLVFFAENVGSNKGAIIGLMVGGVV</i>
Size	0.5 mg
Catalog #	AS-62145
Purity	% Peak Area By HPLC \geq 95%
Description	This peptide is naturally occurring mutant within the beta-amyloid region of b-amyloid protein precursor (APP). This mutation is associated with severe cerebral amyloid beta-protein angiopathy (CAA) in Iowa kindred. The affected individuals share a missense mutation in APP at position 694. This site corresponds to residue 23 of beta-amyloid peptide resulting in substitution of asparagine for aspartic acid.
Storage	-20°C
References	Van Nostrand, W. et al. J. Biol. Chem. 276, 32860 (2001); Grabowski, T. et al. Ann. Neurol. 49, 697 (2001).
Molecular Weight	4328.9
Sequence (One-Letter Code)	DAEFRHDSGYEVHHQKLVFFAENVGSNKGAIIGLMVGGVV
Sequence (Three-Letter Code)	H - Asp - Ala - Glu - Phe - Arg - His - Asp - Ser - Gly - Tyr - Glu - Val - His - His - Gln - Lys - Leu - Val - Phe - Phe - Ala - Glu - Asn - Val - Gly - Ser - Asn - Lys - Gly - Ala - Ile - Ile - Gly - Leu - Met - Val - Gly - Gly - Val - Val - OH

For Research Use Only