

## Technical Data Sheet

<b>Product Name</b>	<b>Beta - Amyloid (11 - 40), FAM - labeled, Human</b> 5 - FAM - EVHHQKLVFFAEDVGSNKGAIIGLMVGGVV
<b>Size</b>	0.1 mg
<b>Catalog #</b>	AS-62950
<b>Purity</b>	% Peak Area By HPLC $\geq$ 95%
<b>Detailed Information</b>	This Abeta peptide (11-40) is FAM-labeled (Abs/Em=494/521 nm). FAM is preferred over FITC because of its photo- and chemical stability. Post-mortem Alzheimer's diseased brain specimens reveal significant levels of A $\beta$ (11-40/42) within insoluble amyloid pools. The $\beta$ -secretase enzyme or $\beta$ -amyloid precursor protein-cleaving enzyme (BACE) generates the N terminus of A $\beta$ , ultimately leading to the production of full-length A $\beta$ (1-40/42) or truncated A $\beta$ (11-40/42). The abundance of A $\beta$ (11-40/42) produced by BACE suggests that their roles in AD pathogenesis may be important.
<b>Storage</b>	-20°C
<b>Molecular Weight</b>	3510.0
<b>Sequence (One-Letter Code)</b>	5-FAM-EVHHQKLVFFAEDVGSNKGAIIGLMVGGVV
<b>Sequence (Three-Letter Code)</b>	5 - FAM - 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*