

Technical Data Sheet

Product Name	ClearPoint™ beta - Amyloid (1 - 40), ¹³C, ¹⁵N - labeled at Arg & Lys, Human DAEF - R* - HDSGYEVHHQ - K* - LVFFAEDVGSN - K* - GAIIGLMVGGVV [R*=R(U - 13C6, U - 15N4) & K*=K(U - 13C6, U - 15N2)]
Size	50 µg
Catalog #	AS-63737
Purity	% Peak Area By HPLC ≥ 95%
Description	ClearPoint™ beta-Amyloid (1-40) is a heavy-isotope labeled peptide. All Arg and Lys residues are heavy-isotope labeled with 13C and 15N
Storage	-20°C
Molecular Weight	4355.9
Sequence (One-Letter Code)	DAEF-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVV [R*=R(U-13C6, U-15N4) & K*=K(U-13C6, U-15N2)]
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 - Asp - Val - Gly - Ser - Asn - Lys* - Gly - Ala - Ile - Ile - Gly - Leu - Met - Val - Gly - Gly - Val - Val - OH [Arg*=Arg(U - 13C6, U - 15N4) and Lys*=Lys(U - 13C6, U - 15N2)]

For Research Use Only