

Biotin-beta-Amyloid (1-42) [beta-Amyloid (1-42), Biotin-Labeled]

Catalog Number: 81110

Unit Size: 100 ug

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Solid

Molecular Weight 5421.19

Soluble In DMSO

Chemical Structure 

Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

Applications

Biotin-beta-Amyloid (1-42) is a biotin-labeled peptide used extensively in neuroscience research, particularly in studies related to Alzheimer's disease. Biotin tag is conjugated to the N-termini of beta-Amyloid (1-42) peptide with sequence: Biotin-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-Ile-Ala-OH (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA). AAT Bioquest's Biotin-beta-Amyloid (1-42) material is thoroughly analyzed to ensure its quality by HPLC, electrophoresis and cell functional analysis. Beta-Amyloid (1-42) is a peptide fragment derived from the amyloid precursor protein (APP). It is the primary component of amyloid plaques found in the brains of individuals with Alzheimer's disease. Beta-Amyloid (1-42) is used in neurotoxicity, neuroinflammation and drug screening studies. A β (1-42) aggregates into fibrils and oligomers that are neurotoxic.