

Biotin-beta-Amyloid (1-28) [beta-Amyloid (1-28), Biotin-Labeled]

Catalog Number: 81005

Unit Size: 100 ug

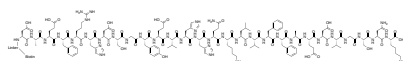
Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	4169.60
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Biotin-beta-Amyloid (1-28) 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-OH (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVG). AAT Bioquest's Biotin-beta-Amyloid (1-28) material is thoroughly analyzed to ensure its quality by HPLC, electrophoresis and cell functional analysis. While the full-length A β peptides like A β (1-40) and A β (1-42) are more commonly associated with plaque formation, A β (1-28) is often used for specific mechanistic and biochemical studies. It is used as a monomeric starting material in seeding and fibril formation experiments.