

Beta-Amyloid (1-28) Azide

Catalog Number: 81001

Unit Size: 100 ug

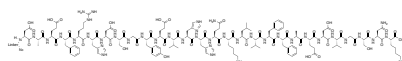
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	3888.22
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Beta-Amyloid (1-28) Azide is an excellent building block for developing Beta-Amyloid (1-28) probes through the well-known click chemistry reaction with an alkynyl group-containing biomolecule under mild reaction conditions with a robust yield. The azide functional group enables bioorthogonal conjugation through azide-alkyne cycloaddition, allowing attachment of fluorophores, affinity tags, or other functional molecules for probe development.

Beta-Amyloid (1-28) Azide is a synthetic peptide corresponding to the N-terminal 28 amino acids of the amyloid- β protein derived from amyloid precursor protein (APP). Similar to the native Beta-Amyloid (1-28) fragment, it can be used as a defined model peptide for studying amyloid peptide structure, aggregation behavior, and intermolecular interactions relevant to amyloid biology.