

Beta-Amyloid (1-42) Azide

Catalog Number: 81101

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	5139.82
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Beta-Amyloid (1–42) Azide is an excellent building block for developing Beta-Amyloid (1–42) 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 (1–42) Azide is a synthetic peptide corresponding to the 42 amino acid sequence of the amyloid- β protein derived from amyloid precursor protein (APP). As a form of amyloid- β associated with amyloid plaque formation and high aggregation propensity, it is commonly used in fibril formation studies, oligomerization research, cellular interaction assays, and in vitro or tissue-based models relevant to Alzheimer's disease research.