

Beta-Amyloid (1-40) Alkyne

Catalog Number: 81052

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	4924.56
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Beta-Amyloid (1–40) Alkyne is an excellent building block for developing Beta-Amyloid (1–40) probes through the well-known click chemistry reaction with an azido group-containing biomolecule under mild reaction conditions with a robust yield. The alkyne functional group enables bioorthogonal conjugation through azide–alkyne cycloaddition, allowing attachment of fluorophores, affinity tags, or other functional molecules for probe development.

Beta-Amyloid (1–40) Alkyne is a synthetic peptide corresponding to the 40 amino acid sequence of the amyloid- β protein derived from amyloid precursor protein (APP). As one of the most abundant amyloid- β isoforms with moderate aggregation propensity, it is commonly used in aggregation kinetics studies, comparative fibrillization assays, and in vitro model systems under defined experimental conditions.