

Azido-PEG4-NHS

Catalog Number: 4535

Unit Size: 25 mg

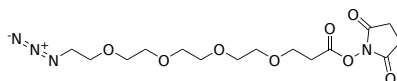
Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	388.38
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Azido-PEG4-NHS is a versatile, high-purity crosslinker designed for efficient bioconjugation and click chemistry applications. It contains an NHS (N-hydroxysuccinimide) ester group that reacts readily with primary amines on proteins, peptides, or amine-coated surfaces, forming stable amide bonds. The azido (N_3) group on the opposite end facilitates click chemistry reactions, including copper-catalyzed azide-alkyne cycloaddition (CuAAC) and the Staudinger ligation with phosphine-modified molecules. The Staudinger ligation, a bioorthogonal and catalyst-free reaction, generates a stable amide linkage, enabling specific, covalent attachment in complex biological environments.