

Biotin PEG4 amine

Catalog Number: 3013

Unit Size: 5 mg

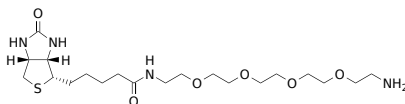
Product Details

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

Chemical Properties

Appearance	Solid white
Molecular Weight	462.61
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Biotin PEG4 amine is a biotinylation reagent with a PEG4 spacer. It is an excellent building block that can be used to make conjugates with carbonyl-containing molecules. Biotin PEG4 amine has a water-soluble and flexible spacer arm that helps to reduce aggregation and steric hindrance.