

Biotin PEG 4 AlkyneCatalog number: 3018
Unit size: 5 mg**Product Details**

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

Chemical Properties

Appearance	Solid
Molecular Weight	281.37
Soluble In	DMSO

Applications

Biotin and biotin derivatives bind with high affinity to both avidins and streptavidins. Biotin alkyne can be readily conjugated to many biomolecules through the well-known click chemistry.