

Biotin PEG2 amine *CAS 138529-46-1*

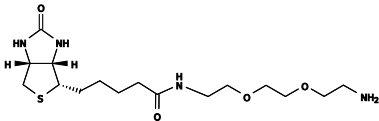
Catalog Number: 3014

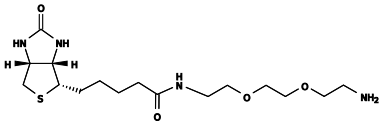
Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid white
Molecular Weight	374.50
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

This carbonyl-reactive biotin derivative contains a long arm (~20 angstrom) increase its avidin-binding affinity. It is widely used to label a variety of biological molecules and samples. Red cells are labeled with this spacers biotin, and the labeled cells can be detected in small blood samples (5 uL) with flow cytometry. Improved labeling efficiency and binding affinity allows an easy detection of positive red cells.