

Product Information

Description:	Biotin-PEG3-azide is a biotinylation reagent that enable Click Chemistry reaction with alkyne, DBCO, BCN or prppargyl moiety to form a stable triazole linkage.
Molecular Formula:	C ₁₈ H ₃₂ N ₆ O ₅ S
Supplied As:	Solid
Solubility:	PBS (10 mg/ml) or DMSO (10 mg/ml)
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	444.6 Da
Cas No.	875770-34-6
Purity:	98%

Chemical Structure

