

Biotin DBCO

Catalog Number	Packaging Size
C330	1 mg

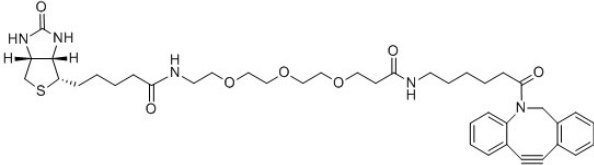
Storage upon receipt: -20°C, protected from light

Introduction

Click chemistry describes a class of chemical reactions that use bio-orthogonal or biologically unique moieties to label and detect a molecule of interest in mild, aqueous conditions. DBCO alkynes can be used to perform click reactions with azide-modified targets without the use of heavy metal catalysis. DBCO reactions are ideal for surface labeling of live cells and also minimize damage to fluorescent proteins like GFP or R-PE.

The biotin DBCO is reactive with azide via a Strain-promoted Azide-Alkyne Click Chemistry reaction (SPAAC). Biotin can be subsequently detected with streptavidin, avidin or NeutrAvidin® biotin-binding protein.

Specifications

Label:	Biotin	
Ex/Em:	-	
Detection Method:	-	
Solubility:	DMSO, DMF	
Molecular Weight:	747.95	
Product Size:	1 mg	
Storage Conditions:	-20 °C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Click chemistry labeling