

Biotin Alkyne

Catalog Number	Packaging Size
C303	5 mg

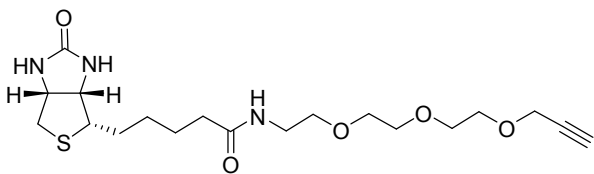
Storage upon receipt: -20°C

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. The click reaction involves a copper-catalyzed triazole formation from an azide and an alkyne. The azide and alkyne moieties can be used interchangeably; either one can be used to tag the molecule of interest, while the other is used for subsequent detection.

The biotin alkyne is reactive with azide via a copper-catalyzed click reaction. 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:	413.53	
Product Size:	5 mg	
Storage Conditions:	-20 °C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Click chemistry labeling