

## Biotin Azide

Catalog Number	Packaging Size
C304	5 mg

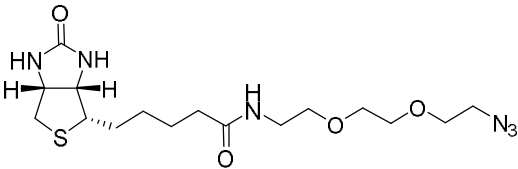
Storage upon receipt: -20°C

### 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 azide is reactive with terminal alkyne via a copper-catalyzed click reaction. Biotin can be subsequently detected with streptavidin, avidin or NeutrAvidin® biotin-binding protein.

### Specifications

<b>Label:</b>	Biotin	
<b>Ex/Em:</b>	—	
<b>Detection Method:</b>	—	
<b>Solubility:</b>	DMSO, DMF	
<b>Molecular Weight:</b>	400.50	
<b>Product Size:</b>	5 mg	
<b>Storage Conditions:</b>	-20 °C, protect from light	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Click chemistry labeling