

Cy3 Azide

Catalog Number	Packaging Size
C312	1 µmol

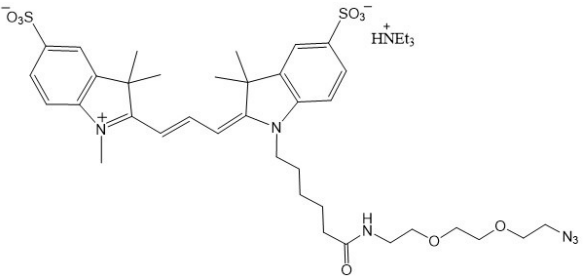
Storage upon receipt: -20°C, protected from light

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. 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 Cy3 azide is reactive with terminal alkyne via a copper-catalyzed click reaction that allows the subsequent visualization by fluorescence spectroscopy.

Specifications

Label:	Cy3	
Ex/Em:	555/565 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Weight:	874.13	
Product Size:	1 µmol	
Storage Conditions:	-20 °C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Click chemistry labeling