

Cy5 Picolyl Azide

Catalog Number	Packaging Size
C338	0.5 µmol

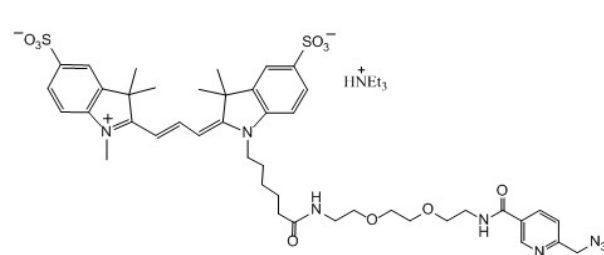
Storage upon receipt: -20°C, protected from light

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 Cy5 picolyl azide is reactive with terminal alkyne via a copper-catalyzed click reaction at a much lower copper (I) concentration without sacrificing reaction efficiency, which protects against undesired copper side reactions with proteins (e.g., GFP, RPE), nucleic acids (e.g., RNA, oligos), and even small molecules (e.g., phalloidin).

Specifications

Label:	Cy5	
Ex/Em:	650/665	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Weight:	1033.48	
Product Size:	0.5 µmol	
Storage Conditions:	-20 °C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Click chemistry labeling