

Cy3 Alkyne

Catalog Number	Packaging Size
C311	1 µmol

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

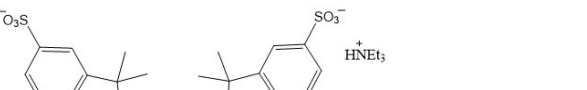
Introduction

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 alkyne is reactive with azide 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:	887.16
Product Size:	1 μ mol
Storage Conditions:	-20 $^{\circ}$ C, protect from light
Shipping Condition:	Room Temperature



The chemical structure of Cy3 dye is shown. It consists of two indole rings connected by a trimethine chain. Each indole ring has a sulfonate group (SO_3^-) at the 6-position. The nitrogen of the right indole ring is substituted with a long alkyl chain ending in a triethylammonium salt (HNEt_3^+) and a polyether chain with a terminal alkyne group.

Applications

Click chemistry labeling