

Catalog Number	Packaging Size
C313	1 µmol

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

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

Label:	Cy5
Ex/Em:	650/665 nm
Detection Method:	Fluorescent
Solubility:	DMSO, DMF
Molecular Weight:	913.20
Product Size:	1 μ mol
Storage Conditions:	-20 $^{\circ}$ C, protect from light
Shipping Condition:	Room Temperature

Click chemistry labeling