

Andy Fluor™ 488 DBCO

Catalog Number	Packaging Size
C331	0.5 µmol

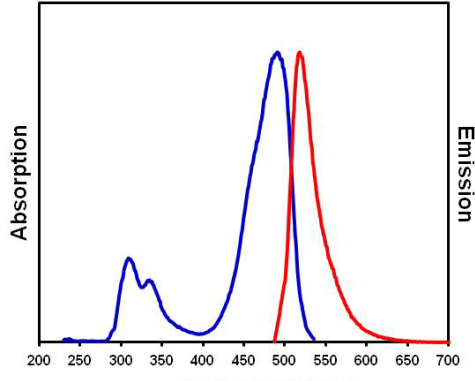
Storage upon receipt: -20°C, protected from light

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. DBCO alkynes can be used to perform click reactions with azide-modified targets without the use of heavy metal catalysis. DBCO reactions are ideal for surface labeling of live cells and also minimize damage to fluorescent proteins like GFP or R-PE.

The Andy Fluor™ 488 DBCO is reactive with azide via a Strain-promoted Azide-Alkyne Click Chemistry reaction (SPAAC).

Specifications

Label:	Andy Fluor™ 488	
Ex/Em:	495/520 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Product Size:	0.5 µmol	
Storage Conditions:	-20 °C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Click chemistry labeling