

## Andy Fluor™ 647 DBCO

Catalog Number	Packaging Size
C334	0.5 µmol

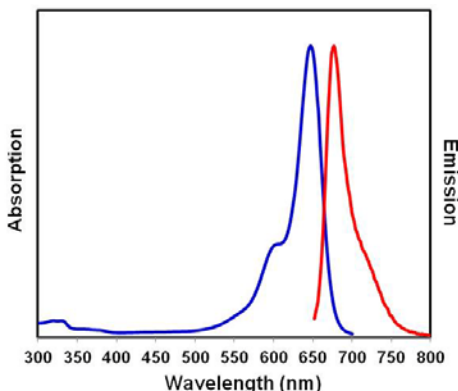
**Storage upon receipt:** -20°C, protected from light

### 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™ 647 DBCO is reactive with azide via a Strain-promoted Azide-Alkyne Click Chemistry reaction (SPAAC).

### Specifications

<b>Label:</b>	Andy Fluor™ 647	
<b>Ex/Em:</b>	650/665 nm	
<b>Detection Method:</b>	Fluorescent	
<b>Solubility:</b>	DMSO, DMF	
<b>Product Size:</b>	0.5 µmol	
<b>Storage Conditions:</b>	-20 °C, protect from light	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Click chemistry labeling