

## Andy Fluor™ 488 Alkyne

Catalog Number	Packaging Size
C317	1 µmol

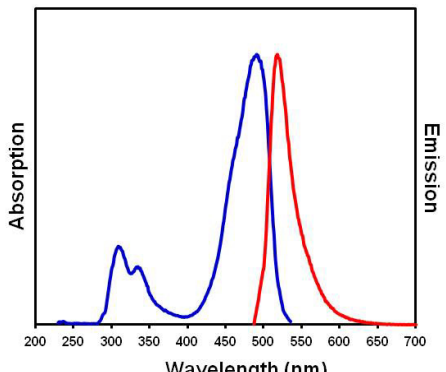
**Storage upon receipt:** -20°C, protected from light

### 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 Andy Fluor™ 488 alkyne is reactive with azide via a copper-catalyzed click reaction that allows the subsequent visualization by fluorescence spectroscopy.

### Specifications

<b>Label:</b>	Andy Fluor™ 488	
<b>Ex/Em:</b>	505/526 nm	
<b>Detection Method:</b>	Fluorescent	
<b>Solubility:</b>	DMSO, DMF	
<b>Product Size:</b>	1 µmol	
<b>Storage Conditions:</b>	-20 °C, protect from light	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Click chemistry labeling