

## Product Information Sheet

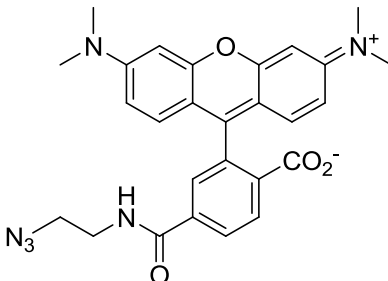
### Ordering Information

Product Number:	490
Product Name:	6-TAMRA azide
Unit Size:	5mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

### Chemical, Physical and Spectral Properties

Molecular Weight: 498.53

Chemical Structure:



Appearance:	Red Solid
Soluble in:	DMSO
Excitation Wavelength:	547 nm
Emission Wavelength:	573 nm

### Application Notes

6-TAMRA azide is an excellent alkyne-reactive fluorescent reagent for labeling alkyne-containing biological molecules through the well known click chemistry. It has the spectral properties identical to those of TRITC conjugates, thus all the instrument settings of the TRITC conjugates would be applicable to the applications of 6-TAMRA azide conjugates.