

## Product Information Sheet

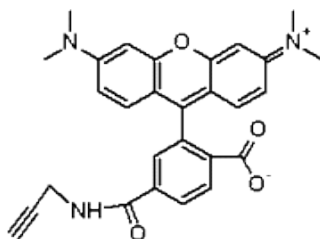
### Ordering Information

Product Number:	491
Product Name:	6-TAMRA alkyne
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 467.52

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	547 nm
Emission Wavelength:	573 nm

### Application Notes

6-TAMRA alkyne is an excellent azide-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 alkyne conjugates.