

Product Information Sheet

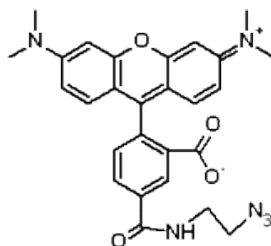
Ordering Information

Product Number:	486
Product Name:	5-TAMRA azide
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: 498.53

Chemical Structure:



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

Application Notes

5-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 5-TAMRA azide conjugates.