

## Product Information Sheet

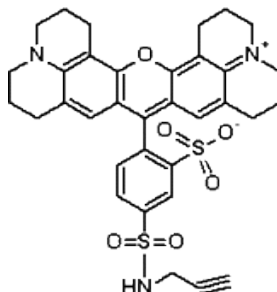
### Ordering Information

Product Number:	485
Product Name:	Texas Red® alkyne *Single Isomer*
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: 643.77

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	588 nm
Emission Wavelength:	601 nm

### Application Notes

Texas Red® 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 Texas Red® conjugates, thus all the instrument settings of the Texas Red® conjugates would be applicable to the applications of Texas Red® alkyne conjugates.