

Product Information Sheet

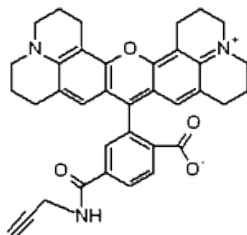
Ordering Information

Product Number:	495
Product Name:	6-ROX 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: 571.66

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	575 nm
Emission Wavelength:	602 nm

Application Notes

6-ROX 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 other 6-ROX conjugates, thus all the instrument settings of the 6-ROX conjugates would be applicable to the applications of 6-ROX alkyne conjugates.