

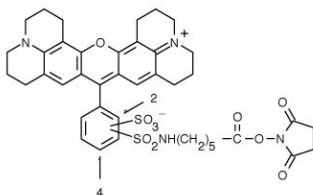
## Product Information Sheet

### Ordering Information

Product Number: 475  
Product Name: Texas Red-X, succinimidyl ester \*Mixed isomers\*  
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: 816.94



Chemical Structure:

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

### Application Notes

The amine-reactive Texas Red®-X, succinimidyl ester is a superior replacement to Texas Red. It has the same spectral properties to Texas Red, but it is much easier conjugated to a biomolecule with much higher conjugation yield. It can be used to create bright red-fluorescent bioconjugates with excitation/emission maxima ~595/615 nm. Compared to Texas Red®, this reactive dye contains an additional seven-atom aminohexanoyl spacer ("X") between the fluorophore and the succinimidyl ester group. This spacer helps to separate the fluorophore from its point of attachment, potentially reducing the interaction of the fluorophore with the biomolecule to which it is conjugated.