

# Texas Red-X, succinimidyl ester \*Single isomer\* \*178623-11-5\*

Catalog number: 474 Unit size: 5 mg

#### **Product Details**

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

### **Chemical Properties**

Appearance Purple solid

Molecular Weight 816.94

Soluble In DMSO

**Chemical Structure** 

## **Spectral Properties**

Excitation Wavelength 586 nm

Emission Wavelength 603 nm

### **Applications**

The amine-reactive Texas Red®-X, succinimidyl ester is a superior replacement to Texas Red (Texas Red® is the trademark of Invitrogen). It has the same spectral properties to Texas Red, but it is much easier conjugated to a biomolecure 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. Compared to the mixed isomers (#474), this product is the single 5-isomer, giving more reproducible results.