

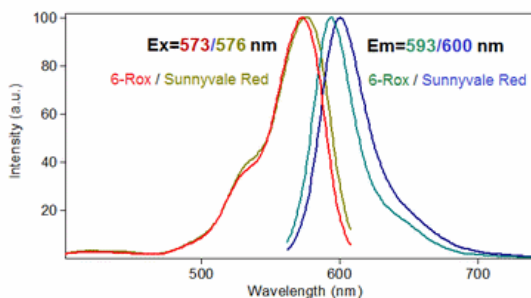
Product Information Sheet

Ordering Information

Product Number: 394
Product Name: Sunnyvale Red™ SE *Superior 6-ROX Replacement*
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: ~650
Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 576 nm
Emission Wavelength: 605 nm

Application Notes

Although ROX dyes have strong fluorescence they are notoriously unstable. Some commercial samples only have a few months shelf life. Compared to 6-ROX labeling compounds, our Sunnyvale Red™ dyes have greatly enhanced stability while they have essentially identical spectral properties to those of 6-ROX when conjugated to a biological substrate (such as oligonucleotides). Our preliminary studies indicated that Sunnyvale Red-derived oligos are more easily purified than the corresponding 6-ROX oligos in some cases.