

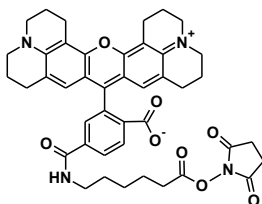
Product Information Sheet

Ordering Information

Product Number: 393
Product Name: 6-ROX-AHA, SE
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 744.83
Appearance: Purple powder
Chemical Structure:



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

Application Notes

6-ROX is predominately used for labeling oligonucleotides and sequencing nucleic acids. However, 6-ROX SE is very unstable compared to other rhodamine dyes. 6-ROX AHA has been developed to improve the stability of 6-ROX SE. The bioconjugates derived from 6-ROX AHA SE give the same spectral properties to those of the bioconjugates prepared from 6-ROX SE. In most cases 6-ROX AHA SE give higher conjugation yields than 6-ROX SE does, and in some cases the conjugates of 6-ROX AHA are brighter than those of 6-ROX probably due to the additional spacer between the ROX fluorophore and biomolecules to be labeled.