

5(6)-ROX, SE [5-(and-6)-Carboxy-X-rhodamine, succinimidyl ester] *Mixed isomers*

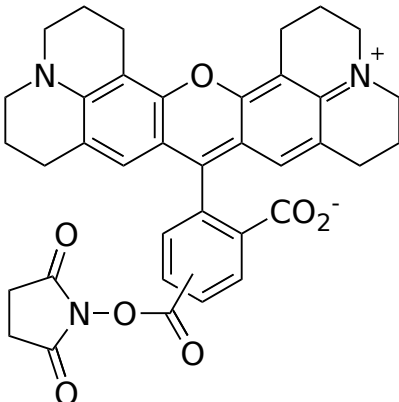
Catalog Number: 390

Unit Size: 25 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	6 months upon receiving

Chemical Properties

Appearance	Solid dark red
Molecular Weight	631.67
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	578 nm
Emission Wavelength	604 nm

Applications

5(6)-ROX, SE is the succinimidyl esters of 5(6)-ROX. It is the mixed isomers of ROX dye. The ratio of 6-isomer/5-isomer is in the range 3/1 to 6/1. This product is the primary labeling reagents for preparing various ROX-based bioconjugates. Compared to other rhodamines, ROX is very unstable. Cautions must be exercised to prevent the deterioration during storage.