

5-ROX, SE [5-Carboxy-X-rhodamine, succinimidyl ester] *CAS 209734-74-7*

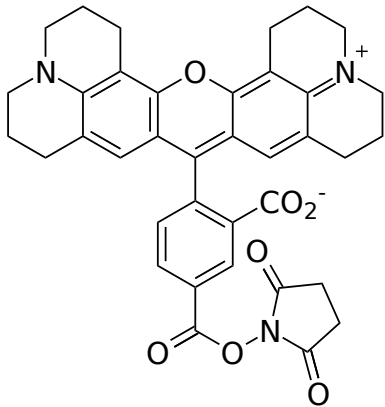
Catalog Number: 389, 391

Unit Size: 5 mg, 100 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

Single isomers of ROX are increasingly preferred for labeling peptides and nucleotides because they give better resolution in HPLC purifications that are often required in bioconjugations. Variable ratios of the 5- and 6-isomers often cause complications in the interpretation of labeling results and assay performances. 5-ROX is predominantly used to label peptides and proteins. Compared to other rhodamines, ROX is very unstable. Cautions must be exercised to prevent the deterioration during storage.