

5(6)-ROX [5-(and 6)-Carboxy-X-rhodamine] *CAS 198978-94-8*

Catalog Number: 380

Unit Size: 100 mg

Product Details

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

Expiration Date 6 months upon receiving

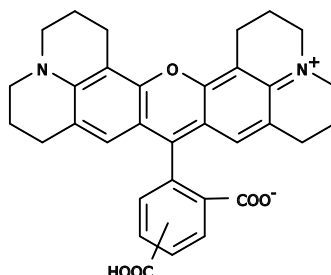
Chemical Properties

Appearance Solid purple

Molecular Weight 534.60

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength 578 nm

Emission Wavelength 604 nm

Applications

ROX dyes are strongly red fluorescent. They have longer excitation and emission wavelengths than the 'conventional' rhodamines. These dyes are used to label peptides, proteins, and other biological ligands. 5-ROX, 6-ROX and their mixture 5(6)-ROX are used to label biomolecules by EDC-mediated reactions. The ratio of 6-isomer/5-isomer is in the range 3/1 to 6/1. Compared to other rhodamines, ROX is very unstable. Cautions must be exercised to prevent the deterioration during storage.