

5-ROX [5-Carboxy-X-rhodamine] *CAS 216699-35-3*

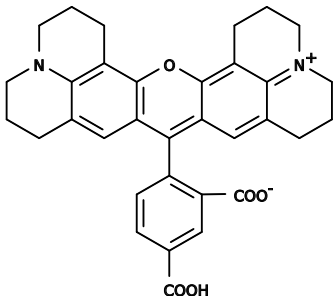
Catalog Number: 381, 385

Unit Size: 5 g, 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	534.60
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	578 nm
Emission Wavelength	604 nm

Applications

5-ROX is the purified single isomer of 5(6)-ROX. It is preferred for some complicated biological applications where reproducibility is more critical than material cost since the minor positional difference between 5-ROX and 6-ROX might significantly affect some biological properties of the underlying conjugates. 5-ROX is predominantly used to label peptides and proteins. Currently AAT Bioquest is using a proprietary process to synthesize isomer-free 5-ROX without the conventional chromatographic purification. Cautions must be exercised to prevent the deterioration during storage.