

6-ROX C2 Maleimide

Catalog Number: 426

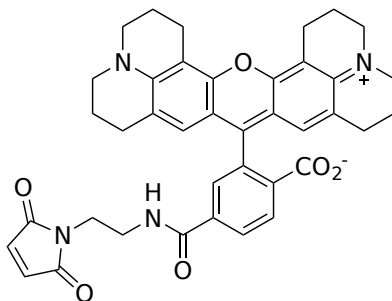
Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid purple
Molecular Weight	656.73
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	578 nm
Emission Wavelength	604 nm

Applications

6-ROX C2 maleimide is an excellent thiol-reactive ROX probe that can be used to develop new ROX bioconjugates. It selectively reacts with thiol groups of peptides, proteins and other biological molecules (such as thiol-modified oligonucleotides and small thiol-containing organic molecules). 6-ROX C2 maleimide has the spectral properties identical to those of 6-ROX conjugates, thus all the instrument settings of the 6-ROX probes would be applicable to the applications of 6-ROX C2 maleimide conjugates. Compared to the commonly used 6-ROX acid and 6-ROX SE, 6-ROX C2 maleimide is much more stable.