

Sulforhodamine 101 *CAS 60311-02-6*

Catalog number: 76

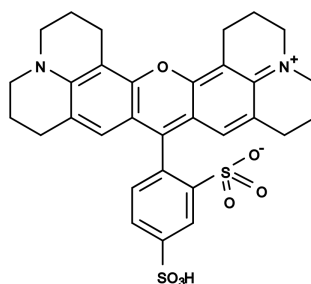
Unit size: 25 mg

Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Purple solid
Molecular Weight	606.71
Soluble In	Water
Chemical Structure	


Spectral Properties

Excitation Wavelength	578 nm
Emission Wavelength	593 nm

Applications

Sulforhodamine 101 is a fluorescence reference standard for measuring fluorescence quantum yield. It is also the key precursor used for preparing rhodamine 101 sulfonyl chloride (also called Texas Red®, a trademark of Molecular Probes).