

Sulforhodamine 101 sulfonyl chloride [Texas Red®] *CAS 82354-19-6*

Catalog Number: 480

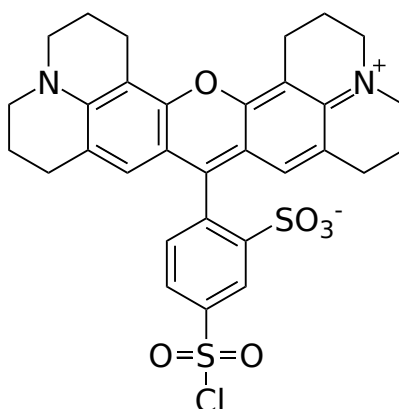
Unit Size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid purple
Molecular Weight	625.16
Soluble In	DMF
Chemical Structure	



Spectral Properties

Excitation Wavelength	586 nm
Emission Wavelength	603 nm

Applications

Sulforhodamine 101 acid chloride (also called Texas Red®) is the most popular labeling reagent of sulfonyl chloride. It is quite unstable in water, especially at the higher pH required for reaction with aliphatic amines. Protein modification by this reagent is frequently done at low temperature (preferably at 4 °C). This reagent reacts with amine compounds such as amino acids, peptides and proteins to give bright red fluorescent conjugates that are extremely stable, and resistant to protease-catalyzed hydrolysis. Unlike succinimidyl esters, sulfonyl chlorides react with both aliphatic amines and aromatic amines indiscriminately.