

Sulforhodamine 101 Sulfonyl Chloride (Texas Red Sulfonyl Chloride)

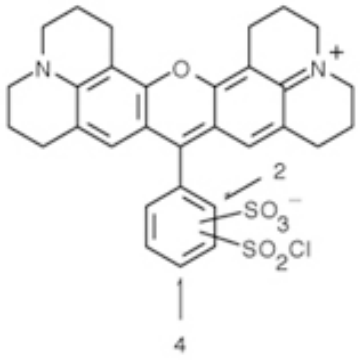
Catalog Number	Packaging Size
C152	5 mg

Storage upon receipt: -20°C, protect from light

Introduction

The amine-reactive **Sulforhodamine 101 Sulfonyl Chloride** (Known as Texas Red® sulfonyl chloride, Trade Mark of Molecular Probes) can be used to create bright red-fluorescent bioconjugates with excitation/emission maxima ~595/615 nm.

Specifications

Label:	Texas Red	
Ex/Em:	595/615 nm	
Detection Method:	Fluorescent	
Solubility:	DMF	
Molecular Formula:	C ₃₁ H ₂₉ ClN ₂ O ₆ S ₂	
Molecular Weight:	625.15	
CAS Number:	82354-19-6	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Co-expression of chemotactic ligand receptors on human peripheral blood monocytes.](#)
Ohura K, Katona IM, Wahl LM, Chenoweth DE, Wahl SM
J Immunol (1987) 138:2633-2639
2. [Texas Red, a hydrophilic, red-emitting fluorophore for use with fluorescein in dual parameter flow microfluorometric and fluorescence microscopic studies.](#)
Titus JA, Haugland R, Sharrow SO, Segal DM
J Immunol Methods (1982) 50:193-204
3. Flow Cytometry and Surface-Marker Phenotyping Using Monoclonal Antibodies: A Combined Approach to Precisely Define the State of the Immune System.
Hadam MR
New Aspects in Physiol Antitumor Substances, Gillissen G, Theurer K 1985; (na):na pp. 120-146