

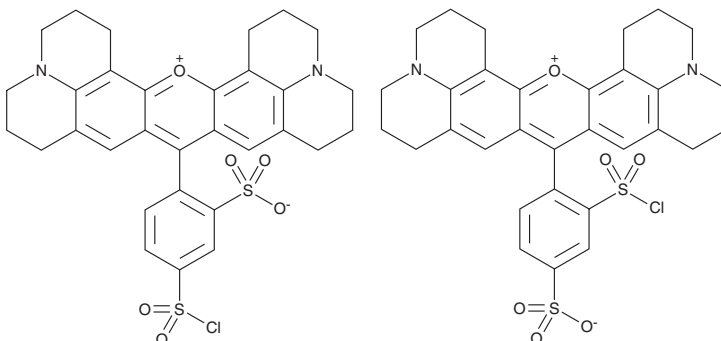
# PRODUCT INFORMATION



## Sulforhodamine 101 Sulfonyl Chloride

Item No. 28547

**CAS Registry No.:** 82354-19-6  
**Formal Name:** 9-[2(or 4)-(chlorosulfonyl)-4(or 2)-sulfophenyl]-2,3,6,7,12,13,16,17-octahydro-1H,5H,11H,15H-xantheno[2,3,4-ij:5,6,7-i'j']diquinolizin-18-ium, inner salt  
**MF:** C<sub>31</sub>H<sub>29</sub>ClN<sub>2</sub>O<sub>6</sub>S<sub>2</sub>  
**FW:** 625.2  
**Purity:** ≥85%  
**Ex./Em.:** 585/602 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Sulforhodamine 101 sulfonyl chloride is supplied as a solid. A stock solution may be made by dissolving the sulforhodamine 101 sulfonyl chloride in the solvent of choice, which should be purged with an inert gas. Sulforhodamine 101 sulfonyl chloride is soluble in the organic solvent dimethyl formamide.

### Description

Sulforhodamine 101 sulfonyl chloride is a fluorescent probe that binds to free amino groups and is a derivative of sulforhodamine 101 (Item No. 16953).<sup>1</sup> It displays excitation/emission maxima of 585/602 nm, respectively. Sulforhodamine 101 sulfonyl chloride has commonly been used as a fluorescent conjugate on antibodies or proteins for the detection of proteins via fluorescent microscopy and flow cytometry applications.<sup>1</sup>

### Reference

1. Titus, J.A., Haugland, R., Sharrow, S.O., *et al.* Texas Red, a hydrophilic, red-emitting fluorophore for use with fluorescein in dual parameter flow microfluorometric and fluorescence microscopic studies. *J. Immunol. Methods* **50**(2), 193-204 (1982).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/19/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)