

Sulforhodamine 101-X, succinimidyl ester (Texas Red-X, succinimidyl ester)

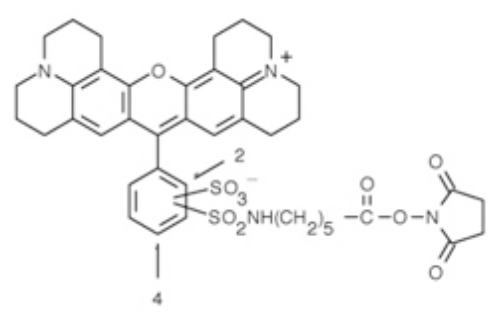
Catalog Number	Packaging Size
C154	5 mg

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

Introduction

The amine-reactive **Sulforhodamine 101-X, succinimidyl ester** (Known as Texas Red-X, succinimidyl ester, 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 ₄₁ H ₄₄ N ₄ O ₁₀ S ₂	
Molecular Weight:	816.94	
CAS Number:	216972-99-5	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

- [Influence of fluorophore and linker composition on the pharmacology of fluorescent adenosine A1 receptor ligands.](#)
Baker JG, Middleton R, Adams L, May LT, Briddon SJ, Kellam B, Hill SJ, Br J Pharmacol (2010) 159:772-786
- [A lipid receptor sorts polyomavirus from the endolysosome to the endoplasmic reticulum to cause infection.](#)
Qian M, Cai D, Verhey KJ, Tsai B, PLoS Pathog (2009) 5:e1000465-e1000465
- [H-type dimer formation of fluorophores: a mechanism for activatable, in vivo optical molecular imaging.](#)
Ogawa M, Kosaka N, Choyke PL, Kobayashi H, ACS Chem Biol (2009) 4:535-546