

5-TRITC (Tetramethylrhodamine-5-isothiocyanate)

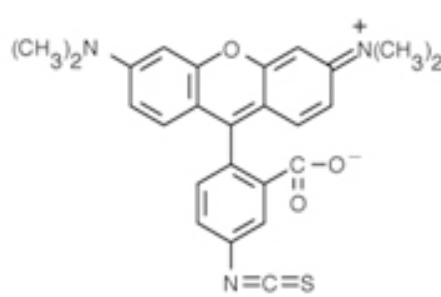
Catalog Number	Packaging Size
C131	5 mg

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

Introduction

The amine-reactive **tetramethylrhodamine-5-isothiocyanate** (5-TRITC, G isomer) can be used to create bright red-orange fluorescent bioconjugates with excitation/emission maxima ~553/575 nm.

Specifications

Label:	Tetramethylrhodamine	
Ex/Em:	553/575 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₅ H ₂₁ N ₃ O ₃ S	
Molecular Weight:	443.52	
CAS Number:	80724-19-2	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [A multistep adhesion cascade for lymphoid progenitor cell homing to the thymus.](#)
Scimone ML, Aifantis I, Apostolou I, von Boehmer H, von Andrian UH
Proc Natl Acad Sci U S A (2006) 103:7006-7011
2. [Caspase-11 regulates cell migration by promoting Aip1-Cofilin-mediated actin depolymerization.](#)
Li J, Brieher WM, Scimone ML, Kang SJ, Zhu H, Yin H, von Andrian UH, Mitchison T, Yuan J
Nat Cell Biol (2007) 9:276-286
3. [Fluorescent dyes for lymphocyte migration and proliferation studies.](#)
Parish CR
Immunol Cell Biol (1999) 77:499-508