

5-TAMRA, SE [5-Carboxy-tetramethylrhodamine, succinimidyl ester]

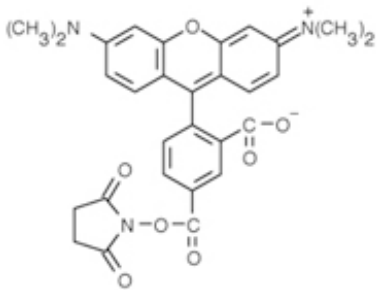
Catalog Number	Packaging Size
C140	5 mg

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

Introduction

5-TAMRA, SE is an amine-reactive fluorescent probe for preparation of bioconjugates and it is photo-stable and pH-insensitive orange-red fluorescent dye.

Specifications

Label:	Tetramethylrhodamine	
Ex/Em:	555/575 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₉ H ₂₅ N ₃ O ₇	
Molecular Weight:	527.53	
CAS Number:	150810-68-7	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Time-Resolved Fluorescence Energy Transfer DNA Helicase Assays for High Throughput Screening.](#)
Earnshaw DL, Moore KJ, Greenwood CJ, Djaballah H, Jurewicz AJ, Murray KJ, Pope AJ
J Biomol Screen (1999) 4:239-248
2. [Zymogen/enzyme discrimination using peptide chloromethyl ketones.](#)
Williams EB, Krishnaswamy S, Mann KG
J Biol Chem (1989) 264:7536-7545
3. [Multiplex detection of single-nucleotide variations using molecular beacons.](#)
Marras SA, Kramer FR, Tyagi S
Genet Anal (1999) 14:151-156