

6-TAMRA, SE [6-Carboxy-tetramethylrhodamine, succinimidyl ester]

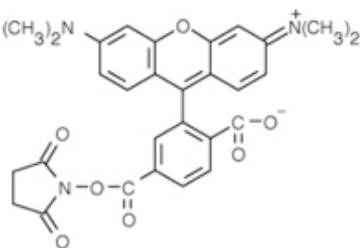
Catalog Number	Packaging Size
C142	5 mg

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

Introduction

6-TAMRA, SE is an amine-reactive form of tetramethylrhodamine. The single-isomer preparations of TAMRA are most important for high-resolution techniques such as DNA sequencing and capillary electrophoresis. 6-TAMRA is one of the traditional fluorophores used for automated DNA sequencing.

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:	150810-69-8	
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. [High sensitivity detection of cancer in vivo using a dual-controlled activation fluorescent imaging probe based on H-dimer formation and pH activation.](#)
Ogawa M, Kosaka N, Regino CA, Mitsunaga M, Choyke PL, Kobayashi H,
Mol Biosyst (2010) 6:888-893