

5(6)-TAMRA, SE [5-(and-6)-Carboxy-tetramethylrhodamine, succinimidyl ester]

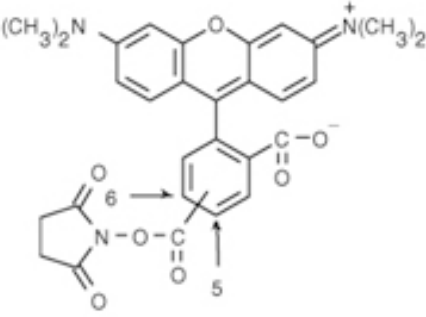
Catalog Number	Packaging Size
C138	25 mg

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

Introduction

Tetramethylrhodamine (TMR) is an important fluorophore for preparing protein conjugates. 5-(and-6)-Carboxytetramethylrhodamine, succinimidyl ester is the amine-reactive, mixed isomer form of TAMRA.

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:	246256-50-8	
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. [Optical properties and application of a reactive and bio reducible thiol-containing tetramethylrhodamine dimer.](#)
Christie RJ, Tadiello CJ, Chamberlain LM, Grainger DW,
Bioconj Chem (2009) 20:476-480