

5-ROX, SE [5-Carboxy-X-rhodamine, succinimidyl ester]

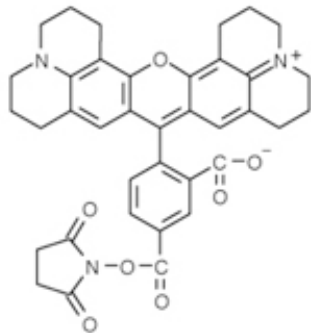
Catalog Number	Packaging Size
C150	5 mg

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

Introduction

The isomeric **5-ROX, SE** is an amine-reactive form of carboxy-X-rhodamine and is widely used for oligonucleotide labeling and automated DNA sequencing applications.

Specifications

Label:	ROX	
Ex/Em:	575/602 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₃₇ H ₃₃ N ₃ O ₇	
Molecular Weight:	631.68	
CAS Number:	344402-35-3	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Zymogen/enzyme discrimination using peptide chloromethyl ketones.](#)
Williams EB, Krishnaswamy S, Mann KG
J Biol Chem (1989) 264:7536-7545
2. [FRET and competing processes between conjugated polymer and dye substituted DNA strands: a comparative study of probe selection in DNA detection.](#)
Al Attar HA, Monkman AP,
Biomacromolecules (2009) 10:1077-1083
3. [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