

5(6)-FAM, SE [5-(and-6)-Carboxyfluorescein, succinimidyl ester]

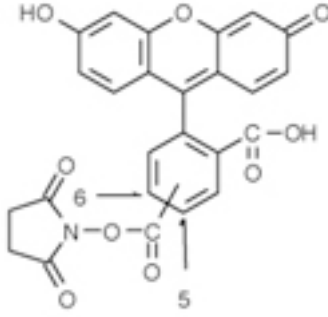
Catalog Number	Packaging Size
C109	100 mg

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

Introduction

5(6)-FAM, SE is a amine-reactive fluorescent probe for preparation of bioconjugates.

Specifications

Label:	Fluorescein	
Ex/Em:	495/520 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₅ H ₁₅ NO ₉	
Molecular Weight:	473.39	
CAS Number:	117548-22-8	
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. [Tyramide signal amplification for analysis of kinase activity by intracellular flow cytometry.](#)
Clutter MR, Heffner GC, Krutzik PO, Sacken KL, Nolan GP,
Cytometry A (2010) 77:1020-1031
3. [Stress renders T cell blasts sensitive to killing by activated syngeneic NK cells.](#)
Rabinovich BA, Shannon J, Su RC, Miller RG
J Immunol (2000) 165:2390-2397