

6-FAM, SE [6-Carboxyfluorescein, succinimidyl ester]

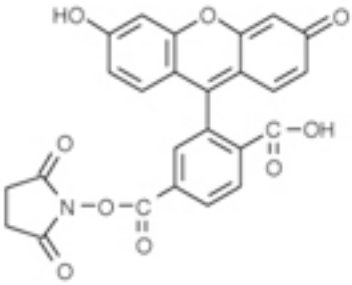
Catalog Number	Packaging Size
C112	10 mg

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

Introduction

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:	92557-81-8	
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. [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
3. [Single-cell analysis of T-cell responses in vitro and in vivo: role of costimulatory signals.](#)
Wells A, Gudmundsdottir H, Walsh M, Turka LA
Transplant Proc (1999) 31:814-815