

5-Carboxy-4'-aminomethylfluorescein

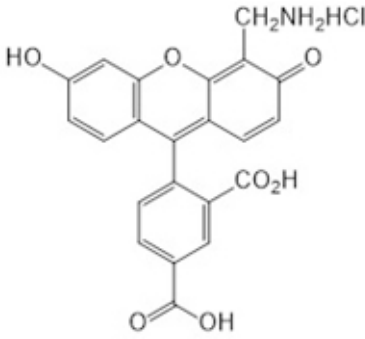
Catalog Number	Packaging Size
C302	10 mg

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

Introduction

5-Carboxy-4'-aminomethylfluorescein is a key building block for developing Big Dye DNA sequencing terminators. It can also be incorporated into a peptide sequence to give fluoresceinated peptides.

Specifications

Label:	Fluorescein	
Ex/Em:	494/520 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₂ H ₁₆ ClNO ₇	
Molecular Weight:	441.82	
CAS Number:	-	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Swelling-based method for preparing stable, functionalized polymer colloids.](#)
Kim AJ, Manoharan VN, Crocker JC
J Am Chem Soc (2005) 127:1592-1593
2. [4'-\[Aminomethyl\]fluorescein and its N-alkyl derivatives: useful reagents in immunodiagnostic techniques.](#)
Shipchandler MT, Fino JR, Klein LD, Kirkemo CL
Anal Biochem (1987) 162:89-101
3. [Redesigning an FKBP-ligand interface to generate chemical dimerizers with novel specificity.](#)
Clackson T, Yang W, Rozamus LW, Hatada M, Amara JF, Rollins CT, Stevenson LF, Magari SR, Wood SA, Courage NL, Lu X, Cerasoli F, Gilman M, Holt DA
Proc Natl Acad Sci U S A (1998) 95:10437-10442