

AMF [4-(Aminomethyl)fluorescein]

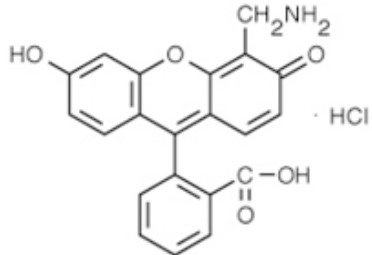
Catalog Number	Packaging Size
C301	25 mg

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

Introduction

The primary aliphatic amine of 4'-(aminomethyl)fluorescein can be reversibly coupled to aldehydes and ketones to form a Schiff base – which can be reduced to a stable amine derivative by sodium borohydride (NaBH₄) or sodium cyanoborohydride (NaCNBH₃) to form new probes. Carboxylic acids of proteins and other water-soluble biopolymers can be coupled to this molecule in aqueous solution using water-soluble carbodiimides such as EDAC.

Specifications

Label:	Fluorescein	
Ex/Em:	494/520 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₁ H ₁₆ ClNO ₅	
Molecular Weight:	397.81	
CAS Number:	91539-64-9	
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