

AMCA-X, SE (6-((7-amino-4-methylcoumarin-3-acetyl)amino)hexanoic acid, succinimidyl ester)

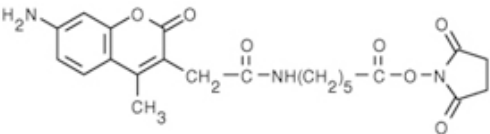
Catalog Number	Packaging Size
C161	10 mg

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

Introduction

The amine-reactive **AMCA-X, SE** (6-((7-amino-4-methylcoumarin-3-acetyl)amino)hexanoic acid, succinimidyl ester) and its conjugates yield blue-fluorescence (approximate excitation/emission maxima ~353/442) that can be used as a contrasting color in multicolor applications.

Specifications

Label:	AMCA	
Ex/Em:	350/448 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₂ H ₂₅ N ₃ O ₇	
Molecular Weight:	443.46	
CAS Number:	216309-02-3	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Identification and mapping of protein-protein interactions between gp32 and gp59 by cross-linking.](#)
Ishmael FT, Alley SC, Benkovic SJ
J Biol Chem (2001) 276:25236-25242
2. [Design of a Molecular Beacon DNA Probe with Two Fluorophores.](#)
Zhang P, Beck T, Tan W
Angew Chem Int Ed Engl (2001) 40:402-405
3. [Double-modification of lectin using two distinct chemistries for fluorescent ratiometric sensing and imaging saccharides in test tube or in cell.](#)
Nakata E, Koshi Y, Koga E, Katayama Y, Hamachi I
J Am Chem Soc (2005) 127:13253-13261