

AMCA Ethylenediamine

Catalog Number: 504

Unit Size: 5 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid light yellow
Molecular Weight	503.35
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	346 nm
Emission Wavelength	434 nm

Applications

AMCA is one of the brightest blue fluorescent labeling dyes. It can be well excited at 365 nm. It is quite photostable, and its fluorescence is pH-independent from pH 4 to 10. This AMC dye can be used for labeling carbonyl compounds (such as carboxy compounds).