

AMCA C2 Maleimide

Catalog Number: 503

Unit Size: 5 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

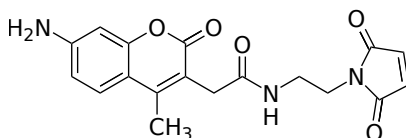
Chemical Properties

Appearance Solid yellow

Molecular Weight 355.34

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 AMCA C2 maleimide can be used for labeling thiol compounds.