

Product Information Sheet

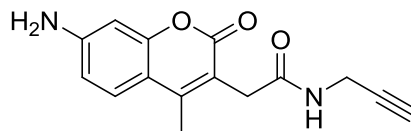
Ordering Information

Product Number: 507
Product Name: AMCA Alkyne
Unit Size: 1mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 270.28

Chemical Structure:



Appearance: Light yellow Solid
Soluble in: DMSO
Excitation Wavelength: 353 nm
Emission Wavelength: 455 nm

Application Notes

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 alkyne compounds through click chemistry.