

## AMCA, succinimidyl ester

Catalog Number: 502

Unit Size: 10 mg

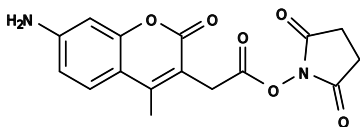
### Product Details

---

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

### Chemical Properties

---

Appearance	Solid yellow
Molecular Weight	330.29
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. AMCA, succinimidyl ester can be used for labeling aliphatic amine compounds.