

Product Information Sheet

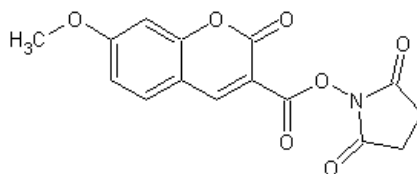
Ordering Information

Product Number:	563
Product Name:	7-Methoxycoumarin-3-carboxylic acid, succinimidyl ester
Unit Size:	100 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 317.25

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	358 nm
Emission Wavelength:	410 nm

Application Notes

7-Methoxycoumarin-3-carboxylic acid, succinimidyl ester is an excellent amino-reactive tag that has strong blue fluorescence. It is used to label peptides, nucleotides. It is also used to label cell membranes although its fluorescence wavelength is quite short.