

## **Product Information Sheet**

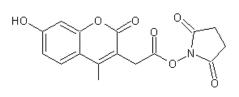
## **Ordering Information**

Product Number:	556
Product Name:	7-Hydroxy-4-methylcoumarin-3-acetic acid, succinimidyl ester
Unit Size:	25 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

331.28

## **Chemical and Spectral Properties**

Molecular Weight: Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	364 nm
Emission Wavelength:	458 nm

## **Application Notes**

7-Hydroxy-4-methylcoumarin-3-acetic acid, SE is an amine-reactive blue fluorophore that has pHdependent and environment-sensitive fluorescence. It is widely used for preparing bioconjugates of blue fluorescence.