

## 7-Hydroxycoumarin-4-acetic acid, succinimidyl ester

Catalog Number: 553

Unit Size: 25 mg

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid light yellow
Molecular Weight	317.25
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

7-Hydroxycoumarin-4-acetic acid is used for labeling peptides and proteins to give blue fluorescence. It can also be used for labeling other biological molecules.