

## 7-Hydroxycoumarin-3-carboxylic acid \*CAS 779-27-1\*

Catalog Number: 549, 550

Unit Size: 25 mg, 250 mg

### Product Details

---

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

### Chemical Properties

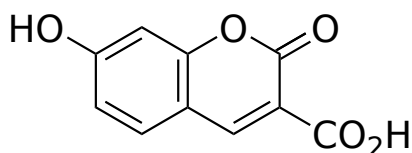
---

Appearance Solid off-white

Molecular Weight 206.15

Soluble In DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength 352 nm

Emission Wavelength 407 nm

### Applications

---

7-Hydroxycoumarin-3-carboxylic acid is a blue fluorophores for labeling proteins and nucleic acids mostly through the in situ formation of its succinimidyl ester catalyzed by EDAC.