

# MCA succinimidyl ester [7-Methoxycoumarin-4-acetic acid, succinimidyl ester]

Catalog Number: 558

Unit Size: 25 mg

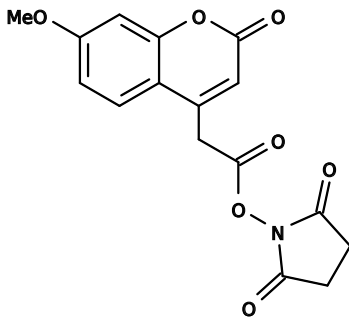
## Product Details

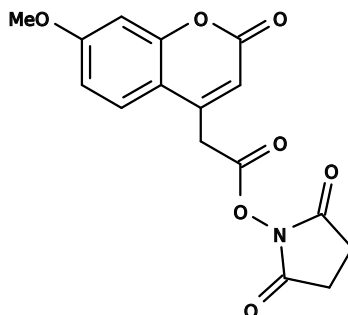
---

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

## Chemical Properties

---

Appearance	Solid off-white
Molecular Weight	331.28
Soluble In	DMSO
Chemical Structure	



## Spectral Properties

---

Excitation Wavelength	322 nm
Emission Wavelength	381 nm

## Applications

---

7-Methoxycoumarin-4-acetic acid (MCA) is widely used for developing FRET peptide substrates for analyzing protease activities. It is commonly paired with dinitrophenyl (DNP) group. It is also used for HPLC derivatization by fluorescence detection. MCA succinimidyl ester selectively reacts with aliphatic amines.