

7-Methoxycoumarin-3-carboxylic acid, succinimidyl ester

CAS 150321-92-9

Catalog Number: 563

Unit Size: 100 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	355 nm
Emission Wavelength	405 nm

Applications

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.