

7-Hydroxycoumarin-3-carboxylic acid, succinimidyl ester

CAS 134471-24-2

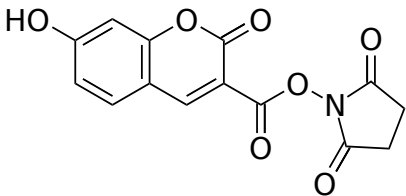
Catalog Number: 551

Unit Size: 50 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	303.23
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	352 nm
Emission Wavelength	407 nm

Applications

7-Hydroxycoumarin-3-carboxylic acid, SE is the amine-reactive succinimidyl ester of 7-Hydroxycoumarin-3-carboxylic acid. It is a blue fluorescent dyes for labeling proteins and nucleic acids. This coumarin is also increasingly used to label peptides, nucleotides and carbohydrates.