

7-Hydroxy-4-methylcoumarin-3-acetic acid, succinimidyl ester

Catalog Number: 556

Unit Size: 25 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

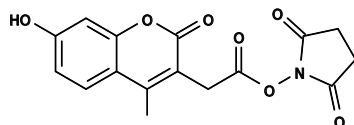
Chemical Properties

Appearance Solid

Molecular Weight 331.28

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

Applications

7-Hydroxy-4-methylcoumarin-3-acetic acid, SE is an amine-reactive blue fluorophore that has pH-dependent and environment-sensitive fluorescence. It is widely used for preparing bioconjugates of blue fluorescence.