

7-Hydroxy-4-methylcoumarin-3-acetic acid *CAS 5852-10-8*

Catalog Number: 554

Unit Size: 100 mg

Product Details

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

Expiration Date 12 months upon receiving

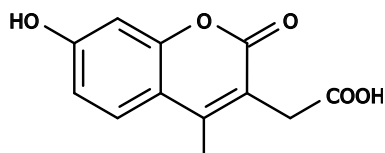
Chemical Properties

Appearance Solid off-white

Molecular Weight 234.21

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

Applications

7-Hydroxy-4-methylcoumarin-3-acetic acid is a blue fluorophore that has pH-dependent and environment-sensitive fluorescence. This coumarin is increasingly used to label peptides, nucleotides and carbohydrates.