

7-Hydroxycoumarin-4-acetic acid *CAS 6950-82-9*

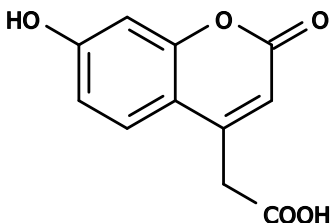
Catalog Number: 552

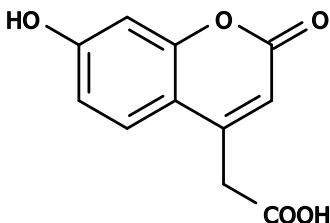
Unit Size: 100 mg

Product Details

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

Chemical Properties

| | |
|--------------------|--|
| Appearance | Solid |
| Molecular Weight | 220.18 |
| Soluble In | DMSO |
| Chemical Structure |  |



Spectral Properties

| | |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength | N/A |

Applications

7-Hydroxycoumarin-4-acetic acid is used for labeling peptides and proteins to give blue fluorescence. It can also be used for labeling other biological molecules.