

MCA [7-Methoxycoumarin-4-acetic acid] *CAS 62935-72-2*

Catalog Number: 557

Unit Size: 1 g

Product Details

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

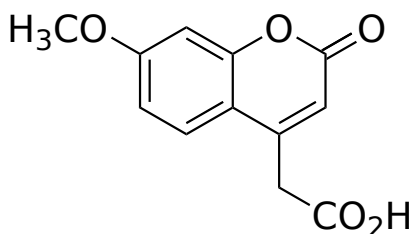
Chemical Properties

Appearance Solid off white

Molecular Weight 234.2

Soluble In DMSO

Chemical Structure

**Spectral Properties**

Excitation Wavelength 322 nm

Emission Wavelength 381 nm

Applications

7-Methoxycoumarin-4-acetic acid (MCA) is widely used for developing FRET peptide substrates for analyzing protease activities. It is commonly paired with dinitrophenyl (DNP) group. It is also used for HPLC derivatization by fluorescence detection.