

7-Methoxycoumarin-3-carboxylic acid *CAS 20300-59-8*

Catalog Number: 560, 561

Unit Size: 1 g, 5 g

Product Details

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

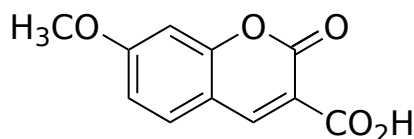
Chemical Properties

Appearance Solid off white

Molecular Weight 220.18

Soluble In DMSO

Chemical Structure

**Spectral Properties**

Excitation Wavelength 355 nm

Emission Wavelength 405 nm

Applications

7-Methoxycoumarin-3-carboxylic acid is used for HPLC derivatization. It is also used to develop FRET probes for analyzing protease activities.