

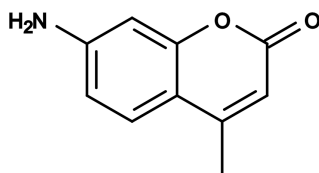
## Product Information Sheet

### Ordering Information

Product Number: 51  
Product Name: AMC [7-Amino-4-methylcoumarin] \*CAS 26093-31-2\*  
Unit Size: 1 g  
Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light  
Expiration Date: 24 months upon receiving

### Chemical and Spectral Properties

Appearance: Solid  
Molecular Weight: 175.19  
Chemical Structure:



Soluble In: DMSO  
Excitation Wavelength: 351  
Emission Wavelength: 430

### Application Notes

This compound is the precursor for a variety of AMC enzyme substrates. It is also used as a calibration standard for AMC-based enzyme substrates. Fluorogenic AMC substrates are widely used for measuring a variety of protease activities.