

AMC (7-Amino-4-methylcoumarin) Standard

Fluorogenic chromophore. Ex(max): 380nm, Em(max): 460nm. Reference compound.

Citations: 3

[View Online »](#)

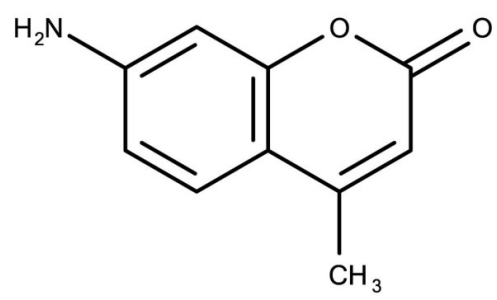
Ordering Information

[Order Online »](#)

BML-KI107-0001	1ml
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Liquid in screw-cap microfuge tube.

CAS 26093-31-2

Formula $C_{10}H_9NO_2$

Formulation 30µM in 50mM HEPES, pH 7.4, 100mM sodium chloride, 0.1% CHAPS, 10mM dithiothreitol, 1mM EDTA, 10% glycerol.

MW 175.2

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)