

HDAC substrate (fluorogenic)

A fluorogenic substrate for histone deacetylase (HDAC). Ex(max): 390nm; Em(max): 460nm.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-137-M001	1mg
ALX-260-137-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Boc-Lys(Ac)-AMC, MAL, Histone deacetylase substrate (fluorogenic), N-(4-Methyl-7-aminocoumarinyl)-N α -(t-butoxycarbonyl)-N ω -acetyllysineamide
Appearance	White crystalline solid
CAS	233691-67-3
Formula	C ₂₃ H ₃₁ N ₃ O ₆
MW	445.5
Purity	≥99%
Sequence	Boc-Lys(Ac)-AMC (AMC=7-Amino-4-methylcoumarin)
Solubility	Soluble in DMSO and 100% ethanol (1mg/ml)

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-be@enzolifesciences.com)

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