

Ac-DMQD-AMC

Caspase-3 substrate

Fluorogenic substrate for caspase-3.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-078-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-3 substrate (fluorogenic)
Appearance	White powder.
Formula	$C_{30}H_{38}N_6O_{12}S$
MW	706.7
Peptide Content	75-95%.
Purity	≥98% (HPLC)
Sequence	Ac-Asp-Met-Gln-Asp-AMC (AMC=7-Amino-4-methylcoumarin)
Solubility	Soluble in DMSO. Dilute with distilled water or buffer. Also soluble in phosphate buffer, pH 7.5
Technical Info / Product Notes	AMC has an excitation maximum of 340-360nm and an emission maximum of 440-460nm.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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