

Mca-YVADAPK(Dnp)-OH

Substrate for caspase-1 and -4

Specific, highly fluorescent substrate for caspase-1. Also substrate for caspase-4.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-023-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 substrate (fluorogenic)
Appearance	White to yellow powder.
Formula	$C_{53}H_{64}N_{10}O_{19}$
MW	1145.1
Purity	≥97% (HPLC)
Sequence	Mca-Tyr-Val-Ala-Asp-Ala-Pro-Lys(Dnp)-OH (Mca = (7-Methoxycoumarin-4-yl)acetyl; Dnp = Dinitrophenyl)
Solubility	Soluble in DMSO or 80% acetic acid (1mg/ml, with agitation).
Technical Info / Product Notes	Mca has an excitation maximum of 325nm and an emission maximum of 392nm.

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