



OMNIMMP[®]

fluorogenic control

Reference peptide for OMNIMMP[®] fluorogenic substrate (BML-P126) and TACE (ADAM17) fluorogenic substrate (BML-P132). Ex.: 328nm, Em.: 393nm. The following Ex/Em have also been used: 320-340/400-420.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P127-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solution in methanol is highly stable at -20°C
Handling	Protect from light and moisture. Store under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Mca-Pro-Leu-OH
Appearance	White powder.
CAS	140430-55-3
Formula	$C_{23}H_{26}N_2O_7$
MW	444.5
Purity	≥98% (TLC)
Sequence	Mca-Pro-Leu-OH [Mca=(7-methoxycoumarin-4-yl)acetyl]
Solubility	Soluble in methanol (50 mg/ml) or DMSO (20 mg/ml).
Source	Synthetic.

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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@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)