

# OMNIMMP<sup>®</sup>

## fluorogenic substrate

Highly quenched and sensitive fluorogenic  
MMP substrate

Fluorogenic substrate for nearly all matrix metalloproteinase (MMP) enzymes, and ADAM17/TACE. This highly-quenched (99.5%) substrate is useful for inhibitor screening and kinetic analysis. Ex.: 328 nm, Em.: 393 nm, although the following Ex/Em have been used: 320-340/400-420. Also available: OMNIMMP<sup>®</sup> Fluorogenic Control Peptide (BML-P127).

Citations: 35

[View Online »](#)

### Ordering Information

[Order Online »](#)

BML-P126-0001	1mg
---------------	-----

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH <sub>2</sub> . AcOH, Matrix metalloproteinase substrate
<b>MW</b>	1093.2
<b>Purity</b>	≥98%
<b>Sequence</b>	Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH <sub>2</sub> . AcOH [Mca=(7-methoxycoumarin-4-yl)acetyl; Dpa=N-3-(2,4-dinitrophenyl)-L-α,β-diaminopropionyl]
<b>Solubility</b>	Soluble in methanol (1mg/ml) or DMSO (20mg/ml).



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)