

# Mca- RPPGFSAFK(Dnp)

## Protease substrate

Fluorogenic substrate for ECE-1 and ECE-2, ACE1 and ACE2, neprilysin, MMP-2 and MMP-9, thimet oligopeptidase, as well as cathepsin A, cathepsin X/Z, BACE, and insulinase, but not MMP-1. Mca fluorescence is quenched by the Dnp group until cleavage separates them. Ex.: 328nm, Em.: 393nm, although the following Ex/Em have also been used: 320-340/400-420. This substrate is useful for inhibitor screening and kinetic analysis.

Citations: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-P227-0001	1mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White to off-white solid.
Formulation	Lyophilized.
MW	1388.5
Purity	≥98% (HPLC)
Sequence	Mca-Arg-Pro-Pro-Gly-Phe-Ser-Ala-Phe-Lys(Dnp)-OH[Mca=(7-methoxycoumarin-4-yl)acetyl;Dnp=2,4-dinitrophenyl]
Solubility	Soluble in DMSO to at least 10mM.
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-  
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)