

Mca-APK(Dnp)

ACE2 substrate

Fluorogenic substrate for angiotensin converting enzyme-2 (ACE2), a zinc-dependent carboxypeptidase with 42% amino acid sequence identity to ACE1. Unlike Mca-YVADAPK(Dnp), this substrate is not cleaved by caspases. Mca fluorescence is quenched by the Dnp group until cleavage (at Pro-Lys) separates them. Ex: 328 nm, Em: 393 nm, although the following Ex/Em have also been used: 320-340/400-420. This highly quenched (99.6%) substrate is useful for inhibitor screening and kinetic analysis. Also available: fluorogenic Mca control peptide (BML-P127).

Citations: 25

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P163-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO stable at -20°C
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ACE2 substrate (fluorogenic)
Appearance	Yellow powder.
Formula	$C_{32}H_{37}N_6O_{12}$
Identity	Determined by MS.
MW	696.7
Purity	≥95% (HPLC)
Sequence	Mca-Ala-Pro-Lys(Dnp) [Mca=(7-methoxycoumarin-4-yl)acetyl;Dnp=2,4-dinitrophenyl]
Solubility	Soluble in DMSO or 10% Ammonium Hydroxide (2 mg/ml).
Source	Synthetic.



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-
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)