

Abz-SDK(Dnp)-P

ACE substrate

Fluorogenic substrate for angiotensin converting enzyme-1 (ACE1) N-domain. The somatic form of ACE1 is a zinc-dependent peptidyl dipeptidase comprising two catalytic domains referred to as the N- and C-domains. The testicular form, due to a downstream start site, contains only the C-domain. Abz fluorescence is quenched by the Dnp group until cleavage (at Asp-Lys) separates them. Ex: 320 nm, Em: 420 nm. This quenched substrate is useful for assay of cell culture ACE1 N-domain activity, inhibitor screening and kinetic analysis.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P162-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ACE1 (N-domain) substrate (fluorogenic)

MW 730.9

Purity ≥97%

Sequence Abz-Ser-Asp-Lys(Dnp)-Pro-OH[Abz=o-aminobenzoicacid;Dnp=2,4-dinitrophenyl]



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)