

# 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 »](#)

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]

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)