

[Ser159]-PKC ϵ (149-164)

PKC substrate

Protein kinase C ϵ (PKC ϵ) pseudosubstrate sequence (149-164) but with an Ala to Ser¹⁵⁹ substitution (ERM $\overline{R}$ PRKRQGSVRRRV) which makes it a substrate for phosphorylation. Phospholipid/phorbol ester-dependent substrate for PKC ϵ , ζ , δ and η .

Ordering Information

[Order Online »](#)

BML-P155-0500	0.5mg
----------------------	-------

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 [Ser159]-Protein kinase C ϵ (149-164), PKC ϵ substrate

MW 2067.5

Purity $\geq 97\%$

Sequence Glu-Arg-Met-Arg-Pro-Arg-Lys-Arg-Gln-Gly-Ser-Val-Arg-Arg-Arg-Val

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)