

PKCζ (113-125) (myristoylated)

PKC inhibitor

Pseudosubstrate sequence from human PKCζ (113-125) completely conserved in rabbit, rat, mouse and *Xenopus*. Sequence is also identical in human PKCι; (114-126) and mouse PKCλ. The N-terminus is myristoylated to allow cell permeability. Peptide inhibitor of atypical PKC's (calcium/phorbol ester/diacylglycerol insensitive). Useful in conjunction with the Myristoylated PKC (20-28), Cell-permeable PKC Inhibitor (BML-P205).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P219-0500	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	Protein kinase C ζ (113-125) (myristoylated), PKC ζ pseudosubstrate
Formula	$C_{90}H_{154}N_{30}O_{17}$
Formulation	Lyophilized.
MW	1928.2
Purity	≥97% (HPLC)
Sequence	N-Myristoyl-Ser-Ile-Tyr-Arg-Arg-Gly-Ala-Arg-Arg-Trp-Arg-Lys-Leu-OH
Solubility	Soluble in water or DMSO.

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-
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)