

PKC ζ

pseudosubstrate (myristoylated)

Selective cell permeable inhibitor of PKC ζ . Activates eNOS (NOS III) through Akt phosphorylation and can be considered as a nitric oxide production activator in endothelial cells.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-155-M001	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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Protein kinase Cζ pseudosubstrate, PKCζ (113-125) (myristoylated)

Appearance White to off-white powder.

Formula $C_{90}H_{154}N_{30}O_{17}$

Formulation Counter Ion: TFA

MW 1928.4

Peptide Content ≥60%

Purity ≥95% (HPLC)

Sequence N-Myristoyl-Ser-Ile-Tyr-Arg-Arg-Gly-Ala-Arg-Arg-Trp-Arg-Lys-Leu-OH

Solubility Soluble in water, DMSO or water-methanol (50:50) .

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)