

Protein kinase C (20-28) (myristoylated)

PKC inhibitor

Pseudosubstrate sequence from PKC α and β . Highly specific inhibitor, which is N-terminal myristoylated to allow membrane permeability. IC₅₀=8 μ M for TPA activation of MARCKS phosphorylation in fibroblast primary cultures, and 98% inhibition at 100 μ M. Partially reverses multidrug resistance in human breast cancer cells.

Ordering Information

[Order Online »](#)

BML-P205-0500	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 PKC (20-28) (myristoylated)

Formulation TFA salt. Lyophilized.

MW 1256.0

Purity ≥98% (HPLC)

Sequence N-Myristoyl-Phe-Ala-Arg-Lys-Gly-Ala-Leu-Arg-Gln-OH

Solubility Soluble in water or in 20mM TRIS-HCl, pH 7.5.

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)