

PKI (14-22) amide (myristoylated)

PKA inhibitor

Heat-stable protein kinase inhibitor (PKI) peptide sequence (14-22). N-terminus is myristoylated to increase cell membrane permeability. The non-myristoylated version of this peptide is a highly specific inhibitor of cAMP-dependent protein kinase (PKA) with $K_i=36$ nM. This product is a useful tool for studying PKA in cellular systems.

Citations: 14

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P210-0500	0.5mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	cAMP-dependent protein kinase inhibitor (14-22) amide (myristoylated)
Appearance	Solid.
CAS	201422-03-9
Formula	$C_{53}H_{100}N_{20}O_{12}$
Formulation	Lyophilized. Trifluoroacetate salt.
MW	1209.5
Purity	≥95% (HPLC)
Sequence	N-myristoyl-Gly-Arg-Thr-Gly-Arg-Arg-Asn-Ala-Ile-NH ₂
Solubility	Soluble to at least 1mg/ml in water.
Source	Synthetic.
Technical Info / Product Notes	Replacement for ADI-HPK-108

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-
de@enzolifesciences.com](mailto:info-de@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)