

PKI (5-24)

PKA inhibitor

Sequence is from the heat-stable skeletal muscle inhibitor protein of PKA. Peptide binds to the catalytic subunit of PKA and displaces the regulatory subunit. Mimics protein substrate by binding to the catalytic site via the Arg-cluster basic residues. $K_i = 2.3$ nM for PKA.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P203-0500	0.5mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	cAMP-dependent protein kinase inhibitor (5-24)
Appearance	White to off-white solid
CAS	99534-03-9
Formula	$C_{94}H_{148}N_{32}O_{31}$
Formulation	Lyophilized.
MW	2222.7
Purity	≥95% (HPLC)
Sequence	Thr-Thr-Tyr-Ala-Asp-Phe-Ile-Ala-Ser-Gly-Arg-Thr-Gly-Arg-Arg-Asn-Ala-Ile-His-Asp
Solubility	Soluble in water.
Source	Synthetic.



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)