

Autocamtide-2

CaM kinase inhibitor

Sequence, KKALRRQETVDAL, derived from Ca^{2+} /calmodulin-dependent protein kinase II autophosphorylation site RQETVD plus basic residues to facilitate phosphocellulose binding. Contains consensus phosphorylation sequence R-X-X-S/T. Highly selective CaMKII substrate ($K_m=2\text{ }\mu\text{M}$ for CaMKII; $50\text{ }\mu\text{M}$ for PKC; no activity towards PKA).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P101-0001	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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

Formula $C_{65}H_{118}N_{22}O_{20}$

Formulation Lyophilized.

MW 1527.8

Purity ≥95% (HPLC)

Sequence H-Lys-Lys-Ala-Leu-Arg-Arg-Gln-Glu-Thr-Val-Asp-Ala-Leu-OH

Solubility Soluble in water.

Source Synthetic.

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)