

AIP

CaM kinase inhibitor

AIP (autocamtide-2-related inhibitory peptide) is a highly specific and potent inhibitor of CaM kinase II. This peptide is derived from the CaM kinase II substrate autocamtide-2 (BML-P101) but alanine is substituted for threonine at the 9-position. AIP at 1 μ M completely inhibited CaM kinase II but did not affect PKC, PKA, CaMKIV and other brain extract kinases. A cell-permeable, myristoylated version is available (BML-P212).

Ordering Information

[Order Online »](#)

BML-P211-0001	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

Alternative Name	Autocamtide-2-related inhibitory peptide
MW	1498.0
Purity	≥97%
Sequence	Lys-Lys-Ala-Leu-Arg-Arg-Gln-Glu-Ala-Val-Asp-Ala-Leu



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)