

# AIP (myristoylated)

CaM kinase inhibitor

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-P212-0500	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	Autocamtide-2-related inhibitory peptide (myristoylated)
Appearance	Powder.
Formulation	Lyophilized. Trifluoroacetate salt.
Identity	Determined by MS.
MW	1708.1
Purity	≥95% (HPLC)
Sequence	N-Myr-Lys-Lys-Ala-Leu-Arg-Arg-Gln-Glu-Ala-Val-Asp-Ala-Leu
Solubility	Soluble in 30% Acetic Acid in Water (5 mg/ml).
Source	Synthetic.

## Technical Info / Product Notes

AIP (autocamtide-2-related inhibitory peptide, Prod. No. BML-P211) is a highly-specific and potent inhibitor of CaM kinase II. The sequence is derived from autocamtide-2 (Prod. No. BML-P101) but alanine is substituted for threonine at the 9-position. AIP at 1 μM completely inhibited CaM kinase II but did not affect PKC, PKA, CaMKIV and other brain extract kinases. Myristoylated AIP (Prod. No. BML-P212) is the same sequence as AIP but is N-terminal myristoylated to increase cell-permeability.



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-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)