

Autocamtide-2-related inhibitory peptide

CaM kinase II inhibitor

Highly specific and potent inhibitor of Ca^{2+} /calmodulin-dependent protein kinase II without affecting cAMP-dependent protein kinase (PKA), protein kinase C (PKC) and Ca^{2+} /calmodulin-dependent protein kinase IV.

Citations: 5

[View Online »](#)

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	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AIP
Appearance	Lyophilized.
CAS	167114-91-2
Formula	$C_{64}H_{116}N_{22}O_{19}$
MW	1497.8
Purity	≥96%
Sequence	H-Lys-Lys-Ala-Leu-Arg-Arg-Gln-Glu-Ala-Val-Asp-Ala-Leu-OH
Solubility	Soluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com