

CaM kinase II (281-309)

CaM kinase inhibitor

Sequence is from Ca^{2+} /calmodulin-dependent kinase A subunit negative regulatory sequence containing the calmodulin binding site (290-309) plus autophosphorylation site (Thr²⁸⁶). Peptide can be phosphorylated at Thr²⁸⁶ by PKC. Useful as a CaM binding peptide and inhibits CaMKII by blocking Ca^{2+} /CaM activation and the enzyme active-site. IC_{50} =80 nM, CaM binding; IC_{50} =2 μM , active-site.

Ordering Information

[Order Online »](#)

BML-P201-0500	0.5mg
----------------------	-------

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 CaMKII (281-309), Ca²⁺/calmodulin-dependent protein kinase II (281-309)

MW 3374.5

Purity ≥97%

Sequence Met-His-Arg-Gln-Glu-Thr-Val-Asp-Cys-Leu-Lys-Lys-Phe-Asn-Ala-Arg-Arg-Lys-Leu-Lys-Gly-Ala-Ile-Leu-Thr-Thr-Met-Leu-Ala

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-
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)