

Calcineurin inhibitory peptide

Sequence is calcineurin (CaN) autoinhibitory domain (rat brain CaN aa 457-482). It inhibits calcineurin with an IC_{50} of 10 μ M using [32 P]myosin light chain as substrate. Inhibits Mn^{2+} -stimulated calcineurin activity but has no effect on Ni^{2+} -stimulated activity. $K_i=4.4$ μ M with radiolabeled [Ala 97]-RII (81-99) peptide as the phosphatase substrate.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PR104-0500	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 CN412

Appearance Lyophilized solid.

CAS 148067-21-4

Formula $C_{124}H_{205}N_{39}O_{39}S$

MW 2930.7

Sequence Ile-Thr-Ser-Phe-Glu-Glu-Ala-Lys Gly-Leu-Asp-Arg-Ile-Asn-Glu-Arg- Met-Pro-Pro-Arg-Arg-Asp-Ala-Met-Pro

Solubility Soluble in water.

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)