

Calyculin A

Inhibitor of PP1 and PP2A

Calyculin A is a cell-permeable and potent inhibitor with high specificity for protein phosphatase 1 (PP1) and 2A (PP2A).

May require a license for import, please [contact us](#) for more information.

Ordering Information

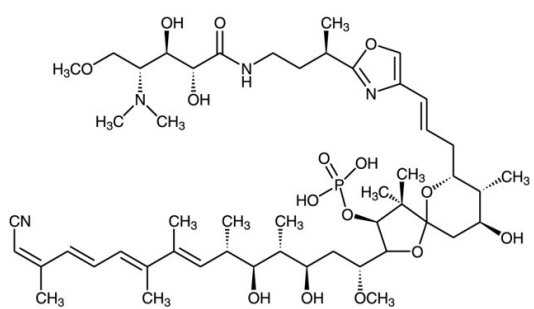
[Order Online »](#)

ENZ-CHM501-0010	10µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- Highly specific, potent, and cell-permeable inhibitor of PP1 and PP2A



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Clear film or white crystalline powder.

CAS 101932-71-2

Couple Target Serine/threonine-protein phosphatase

Couple Type Inhibitor

Formula $C_{50}H_{81}N_4O_{15}P$

MW 1009.2

Purity >98% (HPLC)

Solubility Soluble in DMSO, ethanol, acetone, or methanol.

Last modified: April 23, 2025



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)