

Cyclosporin H

TPA/PMA inhibitor

Immunologically inactive. Does not bind to immunophilin. Potent inhibitor of tumor promoting phorbol ester TPA/PMA (Prod. No. BML-PE160) in mouse skin *in vivo*. Potent inhibitor of Ca^{2+} /calmodulin dependent EF-2 phosphorylation *in vitro*. Potent and selective antagonist of formyl peptide receptor. Inhibits formyl peptide-induced superoxide formation.

Citations: 11

[View Online »](#)

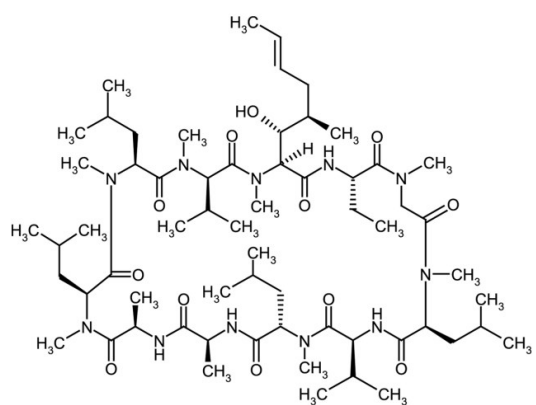
Ordering Information

[Order Online »](#)

ALX-380-286-M001	1mg
ALX-380-286-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Csh-cyclosporin, 5-(N-Methyl-D-valine)-cyclosporin A

Appearance White to off-white powder.

CAS 83602-39-5

Couple Target Formyl peptide receptor, PMA

Couple Type Inhibitor

Formula $C_{62}H_{111}N_{11}O_{12}$

MW 1202.6

Purity ≥95% (HPLC)

Solubility Soluble in 100% ethanol, acetone or ethyl acetate; insoluble in water.

Source Isolated from *Tolypocladium inflatum*.

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)