

Cyclosporin C

Immunosuppressant

Potent immunosuppressor.

Citations: 6

[View Online »](#)

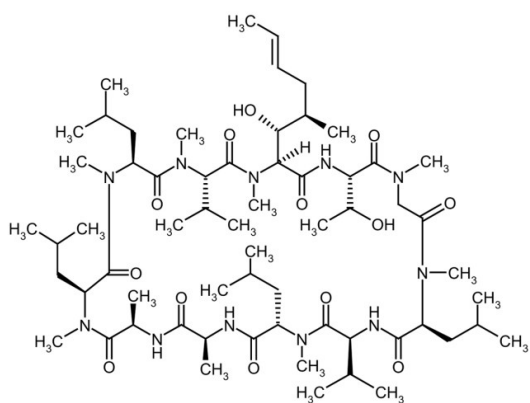
Ordering Information

[Order Online »](#)

ALX-380-282-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 -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Thr2-cyclosporin A, 7-L-Threonine-cyclosporin A

Appearance Colorless to off-white crystalline solid.

CAS 59787-61-0

Formula $C_{62}H_{111}N_{11}O_{13}$

MI **14:** 2752

MW 1218.6

Purity ≥95% (HPLC)

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

Source Isolated from *Fusarium solani*.

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)