

Cyclosporin A

Immunosuppressant

Cyclosporin A (CsA) is an immunosuppressant that has revolutionized organ transplantation through its use in the prevention of graft rejection. CsA binds to cyclophilin and suppresses its proline rotamase activity. The CsA-cyclophilin complex inhibits calcineurin, a type 2B phosphatase. It is an important diagnostic tool for the characterization of permeability transition-related phenomena. It inhibits mitochondrial permeability transition pore opening however, pore inhibition becomes less efficient at increasing Ca^{2+} loads and inhibition is transient. Inhibits staurosporin-induced apoptosis in cultured neural cells and rotenone-induced apoptosis in PC-12 cells. Inhibits cytochrome c release from mitochondria. Inhibits nitric oxide (NO) synthesis.

Citations: 36

[View Online »](#)

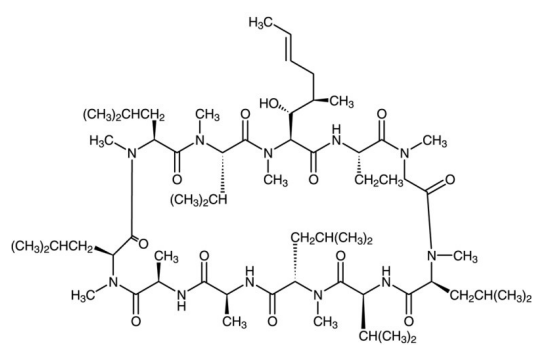
Ordering Information

[Order Online »](#)

BML-A195-0100	100mg
BML-A195-0500	500mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CsA
Appearance	White to off-white powder.
CAS	59865-13-3
Couple Target	NOS, Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{62}H_{111}N_{11}O_{12}$
MW	1202.6
Purity	≥98% (Assay)
Solubility	Soluble in DMSO (25mg/ml), methanol, 100% ethanol (25mg/ml), acetone or petroleum ether; insoluble in water.
Source	Isolated from <i>Fusarium solani</i>

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)