

Myriocin

Palmitoyltransferase inhibitor

Potent immunosuppressant ~10 to 100-fold more potent than cyclosporin A (Prod. No. BML-A195). Blocks sphingolipid biosynthesis by inhibiting serine palmitoyltransferase ($K_i=0.28\text{nM}$), therefore useful as a tool to deplete cells of sphingolipids. Induces apoptosis by depletion of cellular sphingolipids. Inhibits proliferation of an IL-2-dependent mouse cytotoxic cell line CTLL-1 ($\text{IC}_{50}=15\text{nM}$).

Citations: 13

[View Online »](#)

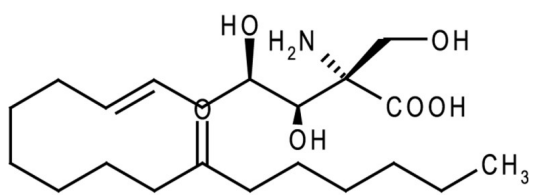
Ordering Information

[Order Online »](#)

BML-SL226-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ISP-1, Thermozytocidin, 2S,3R,4R,6E-2-Amino-3,4-dihydroxy-2-hydroxymethyl-14-oxo-6-eicosenoic acid
Appearance	White to off-white solid.
CAS	35891-70-4
Couple Target	Palmitoyltransferase
Couple Type	Inhibitor
Formula	$C_{21}H_{39}NO_6$
MW	401.5
Purity	≥98% (TLC)
RTECS	JX3890000
Solubility	Soluble in DMSO (25mg/ml; heat briefly in boiling water bath) or in dilute base (5mg/ml; 50mM NaOH).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com