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 =0.28nM), 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 (IC₅₀=15nM).

Citations: 13

View Online »

Ordering Information

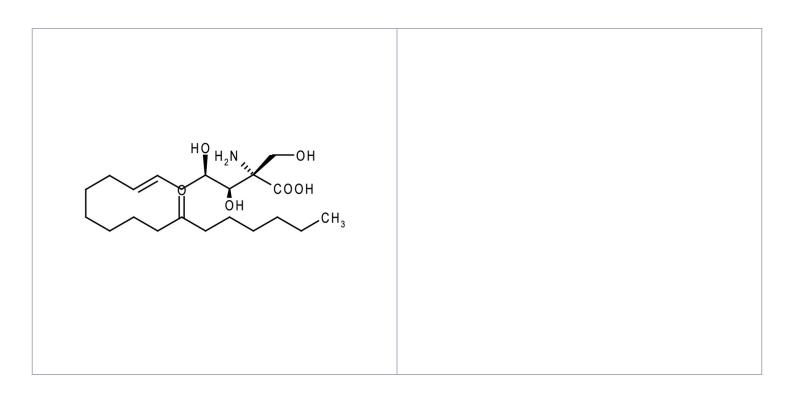
Order Online »

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, Thermozymocidin, 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).

Last modified: May 29, 2024

