

DL-threo-1-Phenyl-2-decanoylamino-3-morpholino-1-propanol

Glucosylceramide synthetase inhibitor

Blocks glucosylation of ceramide by inhibiting UDP-glucose:ceramide glucosyltransferase (glucosylceramide synthetase). Inhibits the expression of cell surface glycolipid antigens and inhibits the growth of cultured rabbit skin fibroblasts. Possesses antitumor activity and inhibits Lewis lung carcinoma cell metastasis *in vitro*. Useful tool for studying the effects of cellular glycosphingolipid depletion. Enhances the apoptotic response to ionizing radiation by enhancing the ceramide signal.

Citations: 15

[View Online »](#)

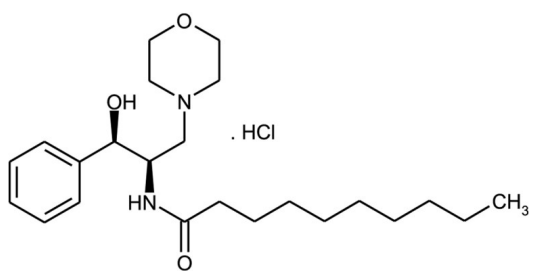
Ordering Information

[Order Online »](#)

BML-SL210-0010	10mg
BML-SL210-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name D,L-threo-PDMP . hydrochloride, DL-PDMP

Appearance White to off-white solid.

CAS 80938-69-8

Couple Target Glucosylceramide synthetase

Couple Type Inhibitor

Formula $C_{23}H_{38}N_2O_3 \cdot HCl$

MW 390.6 . 36.5

Purity ≥98% (TLC)

Solubility 25mg/ml soluble in 100% ethanol, DMSO or water (50°C, vortex or sonicate).

Technical Info / Product Notes Note: Product is not sterile.



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-
be@enzolifesciences.com](mailto:info-be@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)