

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

Inhibits glucosylceramide synthetase. More effective than PDMP in intact cells.

Citations: 3

[View Online »](#)

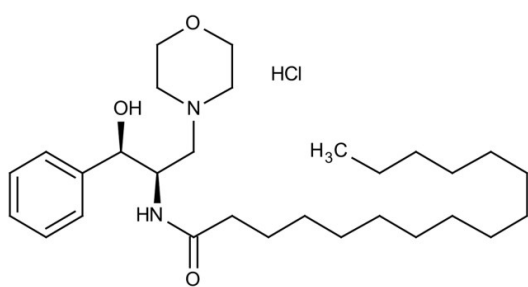
Ordering Information

[Order Online »](#)

BML-SL215-0005	5mg
BML-SL215-0025	25mg

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 DL-PPMP

Appearance White-pink solid.

CAS 149022-18-4

Couple Target Glucosylceramide synthetase

Couple Type Inhibitor

Formula $C_{29}H_{50}N_2O_3 \cdot HCl$

MW 511.2

Purity ≥98% (TLC)

Solubility Soluble in DMSO (10mg/ml).

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)