

PRODUCT INFORMATION



L-threo-PPMP (hydrochloride)

Item No. 24525

Formal Name: N-[(1S,2S)-2-hydroxy-1-(4-morpholinylmethyl)-2-phenylethyl]-hexadecanamide, monohydrochloride

Synonym: L-threo-1-phenyl-2-Palmitoylamino-3-morpholino-1-propanol

MF: C₂₉H₅₀N₂O₃ • HCl

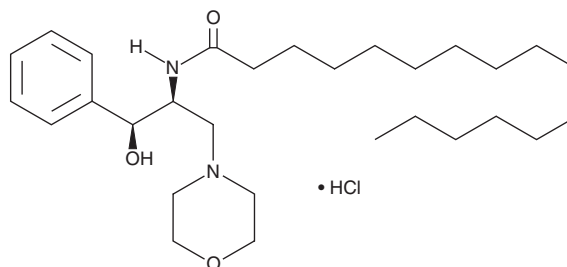
FW: 511.2

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

L-threo-PPMP (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the L-threo-PPMP (hydrochloride) in the solvent of choice, which should be purged with an inert gas. L-threo-PPMP (hydrochloride) is soluble in organic solvents such as ethanol and methanol.

Description

L-threo-PPMP is an inhibitor of glucosylceramide synthetase and isomer of D-threo-PPMP (Item No. 22677).¹ It inhibits the growth of MCF-7, MDA-MB-468, and SK-BR-3 breast cancer cells (ED₅₀s = 4, 7, and 6 μM, respectively) through damage of cell organelle membranes, increasing ceramide concentration, and induction of apoptosis.

Reference

1. Ma, R., Decker, N.M., Anilus, V., et al. Post-translational and transcriptional regulation of glycolipid glycosyltransferase genes in apoptotic breast carcinoma cells: VII. Studied by DNA-microarray after treatment with L-PPMP. *Glycoconj. J.* **26(6)**, 647-661 (2009).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/17/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM