

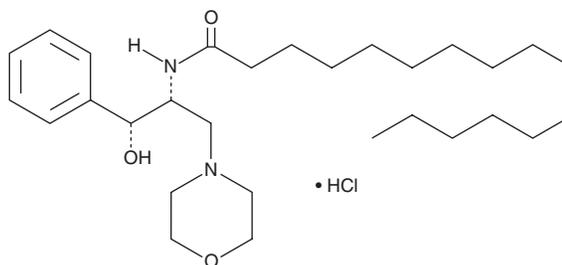
PRODUCT INFORMATION



D-threo-PPMP (hydrochloride)

Item No. 22677

CAS Registry No.: 139889-65-9
Formal Name: [R-(R*,R*)]-N-[2-hydroxy-1-(4-morpholinylmethyl)-2-phenylethyl]-hexadecanamide, monohydrochloride
Synonym: D-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

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

Description

D-threo-PPMP is a glucosylceramide (GlyCer) synthetase inhibitor.^{1,2} It is the active enantiomer and enzymatic inhibitory component of the racemic DL-threo-PPMP (Item No. 17236). In MDCK kidney epithelial cells, D-threo-PPMP induces a 70% reduction in cell growth *in vitro* at 20 μM and significantly inhibits DNA synthesis at 3 μM.³

References

1. Shen, W., Henry, A.G., Paumier, K.L., *et al.* Inhibition of glucosylceramide synthase stimulates autophagy flux in neurons. *J. Neurochem.* **129**(5), 884-894 (2014).
2. Lee, L., Abe, A., and Shayman, J.A. Improved inhibitors of glucosylceramide synthase. *J. Biol. Chem.* **274**(21), 14662-14669 (1999).
3. Abe, A., Inokuchi, J.-i., Jimbo, M., *et al.* Improved inhibitors of glucosylceramide synthase. *J. Biochem.* **111**, 191-196 (1992).

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, 11/08/2022

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