

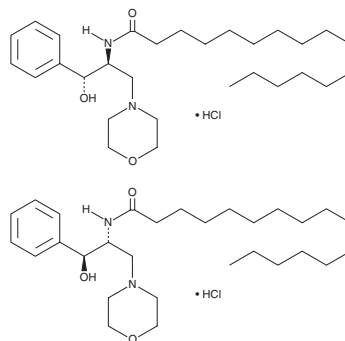
PRODUCT INFORMATION



DL-erythro-PPMP (hydrochloride)

Item No. 24524

Synonym: DL-erythro-1-phenyl-2-Palmitoylamino-3-morpholino-1-propanol
MF: $C_{29}H_{50}N_2O_3 \cdot HCl$
FW: 511.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

DL-erythro-PPMP is an inactive stereoisomer of DL-threo-PPMP (Item No. 17236).¹ It has no effect on glucosylceramide synthetase activity and is used as a negative control for DL-threo-PPMP, D-threo-PPMP (Item No. 22677), and L-threo-PPMP (Item No. 24525) inhibitory activity.

Reference

1. Cabot, M.C. A dual-action approach to multidrug-resistant breast cancer: Prophylaxis to ensure therapeutic effect. (2007).

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, 10/13/2022

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