

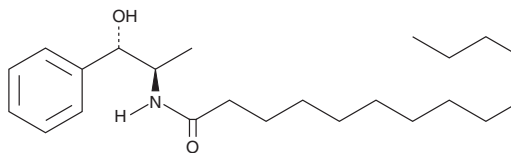
PRODUCT INFORMATION



D-erythro-MAPP

Item No. 10165

CAS Registry No.: 143492-38-0
Formal Name: N-[(1R,2S)-2-hydroxy-1-methyl-2-phenylethyl]-tetradecanamide
Synonym: 1S,2R-D-erythro-2-(N-myristoylamino)-1-phenyl-1-propanol
MF: C₂₃H₃₉NO₂
FW: 361.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

For long term storage, we suggest that D-erythro-MAPP be stored as supplied at -20°C. It should be stable for at least one year.

D-erythro-MAPP is supplied as a crystalline solid. A stock solution may be made by dissolving the D-erythro-MAPP in an organic solvent purged with an inert gas. D-erythro-MAPP is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of D-erythro-MAPP in these solvents is approximately 0.1 mg/ml in DMSO and 20 mg/ml in ethanol and DMF.

D-erythro-MAPP is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, D-erythro-MAPP should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. D-erythro-MAPP has a solubility of approximately 0.5 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

D-erythro-MAPP is an analog of ceramide that suppresses growth of HL-60 cancer cells in culture. D-erythro-MAPP inhibits alkaline ceramidase with an IC₅₀ of approximately 5 μM.¹ Treatment of cultured cells with D-erythro-MAPP elevates the levels of endogenous ceramide, resulting in growth suppression and G₀/G₁ cell cycle arrest.

Reference

1. Bielawska, A., Greenberg, M.S., Perry, D., *et al.* (1S,2R)-D-erythro-2-(N-myristoylamino)-1-phenyl-1-propanol as an inhibitor of ceramidase. *J. Biol. Chem.* **271**, 12646-12654 (1996).

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, 11/10/2017

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