

PRODUCT INFORMATION



C17 D-erythro/L-threo Sphingomyelin (d18:1/17:0)

Item No. 24354

Formal Name: (E)-2-heptadecanamido-3-hydroxyoctadec-4-en-1-yl
(2-(trimethylammonio)ethyl) phosphate

Synonyms: D-erythro/L-threo Heptadecanoyl Sphingomyelin,
D-erythro/L-threo N-Heptadecanoyl
Sphingosylphosphorylcholine,
D-erythro/L-threo SM (18:1/17:0);
D-erythro/L-threo (18:1/17:0)

MF: C₄₀H₈₁N₂O₆P

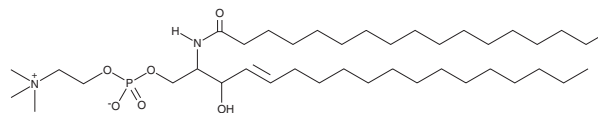
FW: 717.1

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

C17 D-erythro/L-threo Sphingomyelin (d18:1/17:0) is supplied as a solid. A stock solution may be made by dissolving the C17 D-erythro/L-threo Sphingomyelin (d18:1/17:0) in the solvent of choice, which should be purged with an inert gas. C17 D-erythro/L-threo Sphingomyelin (d18:1/17:0) is soluble in organic solvents such as chloroform, ethanol, and methanol.

Description

C17 Sphingomyelin is a synthetic derivative of sphingomyelin (Item Nos. 22674 | 24345). It has been used as an internal standard for the quantification of sphingomyelin.^{1,2} C17 D-erythro L-threo Sphingomyelin is a mixture of D-erythro and L-threo isomers. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Kirkegaard, T., Roth, A.G., Peterson, N.H.T., et al. Hsp70 stabilizes lysosomes and reverts Niemann-Pick disease-associated lysosomal pathology. *Nature* **463(7280)**, 549-553 (2010).
2. Schilling, J.D., Machkovech, H.M., He, L., et al. Palmitate and lipopolysaccharide trigger synergistic ceramide production in primary macrophages. *J. Biol. Chem.* **288(5)**, 2923-2932 (2013).

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, 11/03/2022

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