

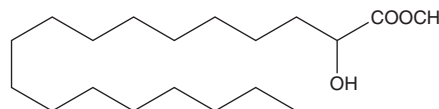
PRODUCT INFORMATION



2-hydroxy Stearic Acid methyl ester

Item No. 24591

CAS Registry No.: 2420-35-1
Formal Name: 2-hydroxy-octadecanoic acid, methyl ester
Synonyms: Methyl 2-Hydroxyoctadecanoate,
Methyl 2-Hydroxystearate,
SFE 19:0;O
MF: C₁₉H₃₈O₃
FW: 314.5
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

2-hydroxy Stearic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy stearic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Stearic acid methyl ester is soluble in organic solvents such as methanol, chloroform, and ethyl ether.

Description

2-hydroxy Stearic acid is a hydroxylated fatty acid methyl ester that broadens phase transition in dimyristoylphosphatidylcholine (DMPC) lipid membranes.¹ It has been used in the synthesis of lipid-nucleotide conjugate anti-HIV agents to increase phosphodiester bond cleavage and the amount of liberated intracellular nucleotides.²

References

1. Jenske, R., Lindstrom, F., Grobner, G., *et al.* Impact of free hydroxylated and methyl-branched fatty acids on the organization of lipid membranes. *Chem. Phys. Lipids* **154(1)**, 26-32 (2008).
2. Gangadhara, K.L., Lescrinier, E., Pannecouque, C., *et al.* Hydroxy fatty acids for the delivery of dideoxynucleosides as anti-HIV agents. *Bioorg. Med. Chem. Lett.* **24(3)**, 817-820 (2014).

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, 03/29/2024

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