PRODUCT INFORMATION



2-hydroxy Lignoceric Acid methyl ester

Item No. 24599

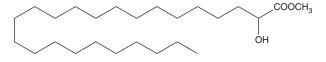
CAS Registry No.: 2433-95-6

Formal Name: 2-hydroxy-tetracosanoic acid, methyl ester

Synonyms: SFE 25:0;O,

methyl 2-hydroxy Tetracosanoate

MF: $C_{25}H_{50}O_3$ FW: 398.7 **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 Lignoceric acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy lignoceric acid methyl ester in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Lignoceric acid methyl ester is soluble in organic solvents such as chloroform and ethyl ether.

Description

2-hydroxy Lignoceric acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in ripe and unripe strawberry homogenates, Pseudosuberites and S. massa sea sponges, sediment samples from the Harney River and Lake Kivu, and the aerial parts of E. helioscopia. 1-5

References

- 1. Gorst-Allman, C.P. and Spiteller, G. Investigation of lipoxygenase-like activity in strawberry homogenates. Z. Lebensm. Unters. Forsch. 187(4), 330-333 (1988).
- 2. Barnathan, G., Kornprobst, J.-M., Doumeng, P., et al. Sponge fatty acids, 5. Characterization of complete series of 2-hydroxy long-chain fatty acids in phospholipids of two Senegalese marine sponges from the family suberitidae: Pseudosuberites sp. and Suberites massa. J. Nat. Prod. 56(12), 2104-2113 (2004).
- 3. Jaffé, R., Rushdi, A.I., Medeiros, P.M., et al. Natural product biomarkers as indicators of sources and transport of sedimentary organic matter in a subtropical river. Chemosphere 64(11), 1870-1884 (2006).
- 4. Al-Mutlaq, K., Standley, L.J., and Simoneit, B.R. Composition and sources of extractable organic matter from a sediment core in Lake Kivu, East African rift valley. Appl. Geochem. 23(5), 1023-1040 (2008).
- Cateni, F., Zilic, J., Altieri, T., et al. Lipid metabolites with free-radical scavenging activity from Euphorbia helioscopia L. Chem. Phys. Lipids 181, 90-98 (2014).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

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