

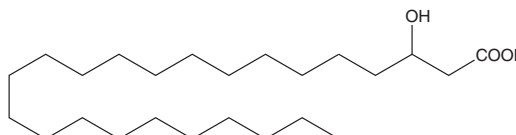
PRODUCT INFORMATION



3-hydroxy Lignoceric Acid

Item No. 22667

CAS Registry No.: 91297-89-1
Formal Name: 3-hydroxy-tetracosanoic acid
Synonym: FA 24:0;O
MF: $C_{24}H_{48}O_3$
FW: 384.6
Purity: $\geq 95\%$
Supplied as: A crystalline 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

3-hydroxy Lignoceric acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-hydroxy lignoceric acid in the solvent of choice. 3-hydroxy Lignoceric acid is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 0.3 mg/ml.

Description

3-hydroxy Lignoceric acid is a hydroxylated form of the 24-carbon saturated lignoceric acid (Item No. 13353). 3-hydroxy Lignoceric acid is found in minor amounts in *Methyloligella*, a novel Gram-negative bacteria.¹ 3-hydroxy Lignoceric acid, in the form of an acyl-CoA metabolite, is an intermediate in fatty acid chain elongation from docosanoic acid (C22:0; Item No. 9000338) to lignoceric acid (C24:0).²

References

1. Doronina, N.V., Poroshina, M.N., Kaparullina, E.N., *et al.* *Methyloligella halotolerans* gen. nov., sp. nov. and *Methyloligella solikamskensis* sp. nov., two non-pigmented halotolerant obligately methylotrophic bacteria isolated from the Ural saline environments. *Syst. Appl. Microbiol.* **36**(3), 148-154 (2013).
2. Guillou, H., Zadavec, D., Martin, P.G., *et al.* The key roles of elongases and desaturases in mammalian fatty acid metabolism: Insights from transgenic mice. *Prog. Lipid Res.* **49**(2), 186-199 (2010).

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, 03/27/2025

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