

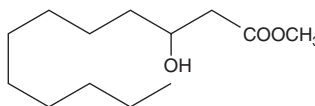
PRODUCT INFORMATION



3-hydroxy Lauric Acid methyl ester

Item No. 24643

CAS Registry No.: 72864-23-4
Formal Name: 3-hydroxy-dodecanoic acid, methyl ester
Synonyms: Methyl 3-Hydroxydodecanoate,
Methyl β -Hydroxydodecanoate
MF: C₁₃H₂₆O₃
FW: 230.3
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

3-hydroxy Lauric acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. 3-hydroxy Lauric acid methyl ester is soluble in chloroform and ether.

Description

3-hydroxy Lauric acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in methyl-branched poly(3-hydroxyalkanoate) (PHA) polymers produced by *P. putida* and *D. tsuruhatensis*.¹

Reference

1. Razaif-Mazinah, M.R.M., Abis, S.N.S., Harun, H.I., et al. Unusual poly(3-hydroxyalkanoate) (PHA) biosynthesis behavior of *Pseudomonas putida* Bet001 and *Delftia tsuruhatensis* Bet002 isolated from palm oil mill effluent. *Biotechnol. Appl. Biochem.* **64**(2), 259-269 (2017).

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, 11/04/2022

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