

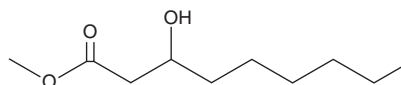
PRODUCT INFORMATION



3-hydroxy Nonanoic Acid methyl ester

Item No. 24612

CAS Registry No.: 83968-06-3
Formal Name: 3-hydroxy-nonanoic acid, methyl ester
Synonyms: Methyl 3-hydroxynonanoate, SFE 10:0;O
MF: C₁₀H₂₀O₃
FW: 188.3
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

3-hydroxy Nonanoic acid methyl ester is supplied as a liquid. 3-hydroxy Nonanoic acid methyl ester is soluble in chloroform, ethanol, and ethyl ether.

Description

3-hydroxy Nonanoic 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. oleovorans*, *P. putida*, and *D. tsuruhatensis*.^{1,2}

References

1. Hazer, B., Lenz, R.W., and Fuller, R.C. Biosynthesis of methyl-branched poly(β-hydroxyalkanoate)s by *Pseudomonas oleovorans*. *Macromol.* **27(1)**, 45-49 (1994).
2. 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, 04/02/2024

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