

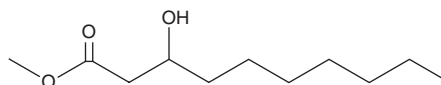
# PRODUCT INFORMATION



## 3-hydroxy Decanoic Acid methyl ester

Item No. 24614

CAS Registry No.: 62675-82-5  
Formal Name: 3-hydroxy-decanoic acid, methyl ester  
Synonyms: methyl 3-hydroxy Decanoate,  
3-Hydroxy C10:0 methyl ester, SFE 11:0;O  
MF:  $C_{11}H_{22}O_3$   
FW: 202.3  
Purity:  $\geq 98\%$   
Supplied as: A liquid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 2$  years



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

### Laboratory Procedures

3-hydroxy Decanoic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 3-hydroxy decanoic acid methyl ester in the solvent of choice. 3-hydroxy Decanoic acid methyl ester is soluble in chloroform, ethanol, and methanol.

### Description

3-hydroxy Decanoic 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*.<sup>1</sup> It has antibacterial activity against *S. faecalis* AD-4.<sup>2</sup>

### References

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).
2. Iizuka, T., Iwadare, T., and Orito, K. Antibacterial activity of myrmicacin and related compounds on pathogenic bacteria in silkworm larvae, *Streptococcus faecalis* AD-4. *J. Fac. Agric. Hokkaido Univ.* **59**(2), 262-266 (1979).

#### 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