

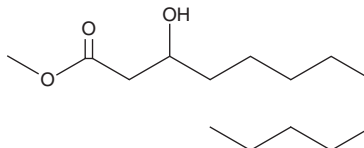
# PRODUCT INFORMATION



## 3-hydroxy Tridecanoic Acid methyl ester

Item No. 24644

**CAS Registry No.:** 150024-70-7  
**Formal Name:** 3-hydroxy-tridecanoic acid, methyl ester  
**Synonyms:** Methyl 3-hydroxytridecanoate, SFE 14:0;O  
**MF:** C<sub>14</sub>H<sub>28</sub>O<sub>3</sub>  
**FW:** 244.4  
**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

3-hydroxy Tridecanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy tridecanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Tridecanoic acid methyl ester is soluble in organic solvents such as chloroform and ethyl ether.

### Description

3-hydroxy Tridecanoic acid methyl ester is a hydroxylated fatty acid methyl ester and major fatty acid component of *A. lipolytica* and *A. glycerini*.<sup>1</sup> It has been used in the synthesis of components of the female sex pheromone of the painted apple moth (*T. anartoides*).<sup>2</sup>

### References

1. Strömpl, C., Tindall, B.J., Jarvis, G.N., et al. A re-evaluation of the taxonomy of the genus *Anaerovibrio*, with the reclassification of *Anaerovibrio glycerini* as *Anaerosinus glycerini* gen. nov., comb. nov., and *Anaerovibrio burkinabensis* as *Anaeroarcus burkinensis* [corrig.] gen. nov., comb. nov. *Int. J. Syst. Bacteriol.* **49(Pt 4)**, 1861-1872 (1999).
2. Muto, S.-e. and Mori, K. Synthesis of the four components of the female sex pheromone of the painted apple moth, *Teia anartoides*. *Biosci. Biotechnol. Biochem.* **67(7)**, 1559-1567 (2003).

#### 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