

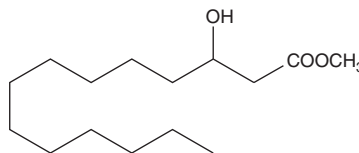
# PRODUCT INFORMATION



## 3-hydroxy Myristic Acid methyl ester

Item No. 24645

**CAS Registry No.:** 55682-83-2  
**Formal Name:** 3-hydroxy-tetradecanoic acid, methyl ester  
**Synonyms:** methyl 3-hydroxy Myristate,  
(±)-3-hydroxy Myristic Acid methyl ester,  
SFE 15:0;O,  
methyl 3-hydroxy Tetradecanoate  
**MF:** C<sub>15</sub>H<sub>30</sub>O<sub>3</sub>  
**FW:** 258.4  
**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 Myristic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy myristic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Myristic acid methyl ester is soluble in organic solvents such as chloroform, methanol, and ethyl ether.

### Description

3-hydroxy Myristic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in *E. camaldulensis* and *E. torelliana* extracts.<sup>1</sup> It is active against *M. tuberculosis* (MIC = 49.5 µg/ml) and is non-cytotoxic to Vero cells (IC<sub>50</sub> = >100 µM). 3-hydroxy Myristic acid methyl ester is also a volatile compound that contributes to the aroma of red wild strawberries (*F. pentaphylla*) but is not present in cultivated strawberries (*Fragaria x ananassa*).<sup>2</sup>

### References

1. Lawal, T.O., Adeniyi, B.A., Adegoke, A.O., *et al.* *In vitro* susceptibility of *Mycobacterium tuberculosis* to extracts of *Eucalyptus camaldulensis* and *Eucalyptus torelliana* and isolated compounds. *Pharm. Biol.* **50(1)**, 92-98 (2012).
2. Duan, W., Sun, P., Chen, L., *et al.* Comparative analysis of fruit volatiles and related gene expression between the wild strawberry *Fragaria pentaphylla* and cultivated *Fragaria x ananassa*. *Eur. Food Res. Technol.* **244(1)**, 57-72 (2018).

#### 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, 03/29/2024

#### 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