

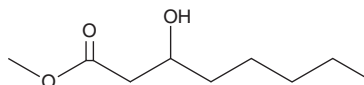
# PRODUCT INFORMATION



## 3-hydroxy Octanoic Acid methyl ester

Item No. 24610

**CAS Registry No.:** 7367-87-5  
**Formal Name:** 3-hydroxy-octanoic acid, methyl ester  
**Synonyms:** 3-hydroxy Caprylic Acid methyl ester,  
 $\beta$ -hydroxy Octanoic Acid methyl ester,  
SFE 9:0;O  
**MF:**  $C_9H_{18}O_3$   
**FW:** 174.2  
**Purity:**  $\geq 98\%$   
**Supplied as:** A liquid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 2$  years



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

### Laboratory Procedures

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

### Description

3-hydroxy Octanoic acid methyl ester is a hydroxylated fatty acid that is a volatile compound and aromatic agent that has been found in pineapple, yuzu, sudachi, and kabosu sour citrus fruits, and human axillary sweat.<sup>1-3</sup>

### References

1. Teai, T., Claude-Lafontaine, A., Schippa, C., *et al.* Volatile compounds in fresh pulp of pineapple (*Ananas comosus* [L.] Merr.) from French Polynesia. *J. Essent. Oil Res.* **13(5)**, 314-318 (2001).
2. Tomiyama, K., Aoki, H., Oikawa, T., *et al.* Characteristic volatile components of Japanese sour citrus fruits: Yuzu, Sudachi and Kabosu. *Flavour Fragr. J.* **27(5)**, 341-355 (2012).
3. Martin, A., Saathoff, M., Kuhn, F., *et al.* A functional ABCC11 allele is essential in the biochemical formation of human axillary odor. *J. Invest. Dermatol.* **130(2)**, 529-540 (2010).

#### 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, 04/01/2024

#### 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