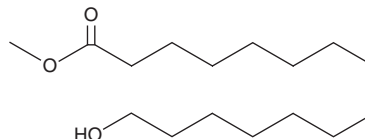


# PRODUCT INFORMATION

## 15-hydroxy Pentadecanoic Acid methyl ester

Item No. 24649

**CAS Registry No.:** 76529-42-5  
**Formal Name:** 15-hydroxy-pentadecanoic acid, methyl ester  
**Synonyms:** Methyl 15-hydroxy Pentadecanoate, SFE 16:0;O  
**MF:**  $C_{16}H_{32}O_3$   
**FW:** 272.4  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}\text{C}$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

15-hydroxy Pentadecanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 15-hydroxy pentadecanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 15-hydroxy Pentadecanoic acid methyl ester is soluble in chloroform, ethyl ether, and warm ethanol.

### Description

15-hydroxy Pentadecanoic acid methyl ester is a hydroxylated fatty acid methyl ester that is a key intermediate in the synthesis of the musk-odored macrocyclic lactones cyclopentadecanolide and exaltolide.<sup>1,2</sup>

### References

1. Lai, F.A., Liu, X., Li, W., et al. Macrolactonization of methyl 15-hydroxypentadecanoate to cyclopentadecanolide over Mo-Fe/HZSM-5 catalyst. *Reac. Kinet. Mech. Cat.* **100**(2), 407-415 (2010).
2. Ravi, S., Padmanabhan, D., and Mamdapur, V.R. Macrocyclic musk compounds: Synthetic approaches to key intermediates for exaltolide, exaltone and dilactones. *J. Indian Inst. Sci.* **81**(3), 299-312 (2001).

#### 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