

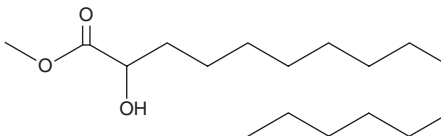
# PRODUCT INFORMATION



## 2-hydroxy Palmitic Acid methyl ester

Item No. 24589

**CAS Registry No.:** 16742-51-1  
**Formal Name:** 2-hydroxy-hexadecanoic acid, methyl ester  
**Synonyms:** Methyl  $\alpha$ -hydroxypalmitate, Methyl 2-hydroxyhexadecanoate,  $\alpha$ -hydroxy Palmitic Acid ethyl ester, SFE 17:0;O  
**MF:** C<sub>17</sub>H<sub>34</sub>O<sub>3</sub>  
**FW:** 286.5  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

2-hydroxy Palmitic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy palmitic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Palmitic acid methyl ester is soluble in methanol, chloroform, and ethyl ether.

### Description

2-hydroxy Palmitic acid methyl ester is a naturally occurring hydroxy fatty acid methyl ester. It is a minor component in essential oils of *P. sylvestris* and *B. javanica* extracts which have larvicidal and cytotoxic activities, respectively.<sup>1,2</sup>

### References

1. Fayemiwo, K.A., Adeleke, M.A., Okoro, O.P., *et al.* Larvicidal efficacies and chemical composition of essential oils of *Pinus sylvestris* and *Syzygium aromaticum* against mosquitoes. *Asian Pac. J. Trop. Biomed.* **4(1)**, 30-34 (2014).
2. Sou, Z., Huang, H., Li, J., *et al.* Chemical composition and cytotoxic activities of petroleum ether fruit extract of fruits of *Brucea javanica* (Simarubaceae). *Trop. J. of Pharma. Research* **12(5)**, 735-742 (2013).

#### 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