

PRODUCT INFORMATION



10-methyl Palmitic Acid methyl ester

Item No. 24823

CAS Registry No.: 2490-51-9
Formal Name: 10-methyl-hexadecanoic acid, methyl ester
Synonyms: Methyl 10-methyl Hexadecanoic Acid, SFE 18:0
MF: C₁₈H₃₆O₂
FW: 284.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

10-methyl Palmitic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform purged with an inert gas can be used.

Description

10-methyl Palmitic acid methyl ester is a methylated fatty acid methyl ester that has been found in cyclitol-containing glycolipid analogs isolated from *P. crassa*.¹

Reference

1. Costantino, V., Fattorusso, E., and Mangoni, A. Isolation of five-membered cyclitol glycolipids, crasserides: Unique glycerides from the sponge *Pseudoceratina crassa*. *J. Org. Chem.* **58(1)**, 186-191 (1993).

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, 03/04/2024

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