

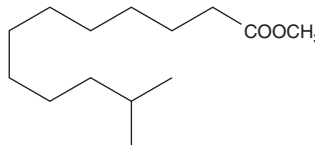
PRODUCT INFORMATION



11-methyl Lauric Acid methyl ester

Item No. 24809

CAS Registry No.: 5129-57-7
Formal Name: 11-methyl-dodecanoic acid, methyl ester
Synonyms: Methyl 11-Methyldodecanoate, SFE 14:0
MF: C₁₄H₂₈O₂
FW: 228.4
Purity: ≥98%
Supplied as: A solution in hexane
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

11-methyl Lauric acid methyl ester is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methylene chloride and ethyl ether purged with an inert gas can be used.

Description

11-methyl Lauric acid methyl ester is a methylated fatty acid methyl ester that has been used as a standard in the quantification of 11-methyl lauric acid in vernix caseosa and *M. oleifera* extract.^{1,2}

References

1. Hauff, S. and Vetter, W. Exploring the fatty acids of *vernix caseosa* in form of their methyl esters by off-line coupling of non-aqueous reversed phase high performance liquid chromatography and gas chromatography coupled to mass spectrometry. *J. Chromatogr. A* **1217(52)**, 8270-8278 (2010).
2. Surendra, T.V., Roopan, S.M., Arasu, M.V., et al. Phenolic compounds in drumstick peel for the evaluation of antibacterial, hemolytic and photocatalytic activities. *J. Photochem. Photobiol. B.* **161**, 463-471 (2016).

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