

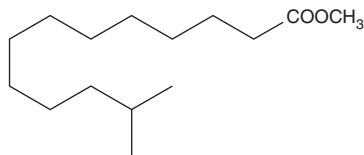
PRODUCT INFORMATION



12-methyl Tridecanoic Acid methyl ester

Item No. 24810

CAS Registry No.: 5129-58-8
Formal Name: 12-methyl-tridecanoic acid, methyl ester
Synonyms: Methyl 12-methyltridecanoate,
Methyl isomyristate, SFE 15:0
MF: C₁₅H₃₀O₂
FW: 242.4
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

12-methyl Tridecanoic 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 and ethyl ether purged with an inert gas can be used.

Description

12-methyl Tridecanoic acid methyl ester is a methylated fatty acid methyl ester that has been found in beef tallow, mutton tallow, and cod liver oil.¹ It has been used as a standard for the quantification of 12-methyl tridecanoic acid in the triglyceride fraction of skin surface lipids from human males by GC-MS.²

References

1. Chin, S.-T., Che Man, Y.B., Tan, C.P., *et al.* Rapid profiling of animal-derived fatty acids using fast GC × GC coupled to time-of-flight mass spectrometry. *J. Am. Oil Chem. Soc.* **86(10)**, 949-958 (2009).
2. Marzouki, Z.M.H., Taha, A.M., and Gomaa, K.S. Fatty acid profiles of sebaceous triglycerides by capillary gas chromatography with mass-selective detection. *J. Chromatogr.* **425(1)**, 11-24 (1988).

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, 04/05/2024

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