

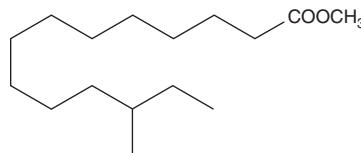
PRODUCT INFORMATION



12-methyl Myristic Acid methyl ester

Item No. 24818

CAS Registry No.: 5129-66-8
Formal Name: 12-methyl-tetradecanoic acid, methyl ester
Synonyms: Methyl 12-methyltetradecanoate, SFE 16:0
MF: $C_{16}H_{32}O_2$
FW: 256.4
Purity: $\geq 98\%$
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

12-methyl Myristic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 12-methyl myristic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 12-methyl Myristic acid methyl ester is soluble in chloroform, ethanol, and ether.

Description

12-methyl Myristic acid methyl ester is a methylated fatty acid methyl ester that has been found in vermicomposts of cattle manure, carica papaya leaves, and cuticular wax of *K. africana*.¹⁻³ It is a volatile compound in lipid-lowering granulated tea.⁴ Levels of 12-methyl myristic acid methyl ester are decreased in *T. cruzi* treated with nifurtimox (Item No. 21784) compared to non-treated controls.⁵

References

1. Balmori-Martinez, D., Spaccini, R., Aguiar, N.O., et al. Molecular characteristics of humic acids isolated from vermicomposts and their relationship to bioactivity. *J. Agric. Food Chem.* **62**(47), 11412-11419 (2014).
2. Abd El Azim, M. Fatty acid constituents and antimicrobial activities of strawberry and carica- papaya leaves. *NPAIJ* **9**(4), 143-147 (2013).
3. Olubunmi, A., Gabriel, O.A., Stephen, A.O., et al. Antioxidant and antimicrobial activity of cuticular wax from *Kigelia africana*. *FABAD J. Pharm. Sci.* **34**(4), 187-194 (2009).
4. Ding, Y., Pu, L., and Kan, J. Hypolipidemic effects of lipid-lowering granulated tea preparation from *Monascus*-fermented grains (adlay and barley bran) mixed with lotus leaves on Sprague-Dawley rats fed a high-fat diet. *J. Funct. Foods* **32**, 80-89 (2017).
5. Barreto-Bergter, E., Hogge, L., and da Cruz, F.S. Lipid alterations induced by nifurtimox in *Trypanosoma cruzi*. *Mol. Biochem. Parasitol.* **21**(3), 221-226 (1986).

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, 01/15/2026

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