

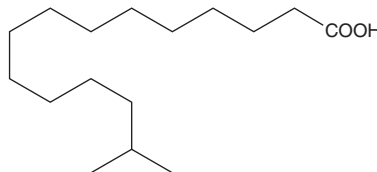
PRODUCT INFORMATION



14-methyl Pentadecanoic Acid

Item No. 36459

CAS Registry No.: 4669-02-7
Formal Name: 14-methyl-pentadecanoic acid
Synonyms: FA 16:0, *iso*-C16:0, Isohexadecanoic Acid, Isopalmitic Acid
MF: $C_{16}H_{32}O_2$
FW: 256.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

14-methyl Pentadecanoic acid is supplied as a solid. A stock solution may be made by dissolving the 14-methyl pentadecanoic acid in the solvent of choice, which should be purged with an inert gas. 14-methyl Pentadecanoic acid is soluble in ethanol, DMSO, and dimethyl formamide.

Description

14-methyl Pentadecanoic acid is a branched-chain fatty acid that has been found in soil bacteria.^{1,2}

References

1. Kim, I., Choi, J., Chhetri, G., *et al.* *Lysobacter helvus* sp. nov. and *Lysobacter xanthus* sp. nov., isolated from Soil in South Korea. *Antonie Van Leeuwenhoek* **112**(8), 1253-1262 (2019).
2. Erwin, G.S., Heikkinen, J., Halimaa, P., *et al.* *Streptomyces lasalocidi* sp. nov. (formerly '*Streptomyces lasaliensis*'), an actinomycete isolated from soil which produces the polyether antibiotic lasalocid. *Int. J. Syst. Evol. Microbiol.* **70**(5), 3076-3083 (2020).

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, 02/20/2024

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