

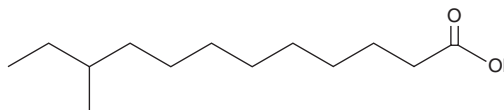
PRODUCT INFORMATION



10-methyl Lauric Acid

Item No. 36504

CAS Registry No.: 7416-57-1
Formal Name: 10-methyl-dodecanoic acid
Synonyms: 10-methyl Dodecanoic Acid,
FA 13:0
MF: C₁₃H₂₆O₂
FW: 214.3
Purity: ≥95%
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 Lauric acid 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 10-methyl lauric acid in these solvents is approximately 10 and 2 mg/ml, respectively.

Description

10-methyl Lauric acid is a methyl branched-chain fatty acid.^{1,2} It has been found in the Gram-positive bacteria *B. acetylicum*, as well as the mandibular gland secretions of carpenter ant species, including *C. herculeanus*, *C. ligniperda*, and *C. pennsylvanicus*.

References

1. Brand, J.M., Fales, H.M., Sokoloski, E.A., *et al.* Identification of mellein in the mandibular gland secretions of carpenter ants. *Life Sci.* **13(3)**, 201-211 (1973).
2. Bernard, K.A., Bellefeuille, M., and Ewan, E.P. Cellular fatty acid composition as an adjunct to the identification of asporogenous, aerobic gram-positive rods. *J. Clin. Microbiol.* **29(1)**, 83-89 (1991).

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