

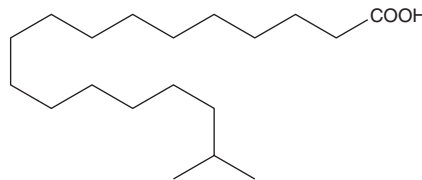
PRODUCT INFORMATION



17-methyl Stearic Acid

Item No. 26766

CAS Registry No.: 2724-59-6
Formal Name: 17-methyl-octadecanoic acid
Synonyms: iso-19:0, iso-C19, FA 19:0, Isononadecanoic Acid
MF: C₁₉H₃₈O₂
FW: 298.5
Purity: ≥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

17-methyl Stearic acid is supplied as a solid. A stock solution may be made by dissolving the 17-methyl stearic acid in the solvent of choice. 17-methyl Stearic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 17-methyl stearic acid in these solvents is approximately 20, 10, and 30 mg/ml, respectively.

Description

17-methyl Stearic acid is a methyl branched-chain fatty acid that has been found in murine meibomian glands, *C. cornucopioides* mushrooms, and *Phyllobacterium* species.¹⁻³ It has also been found in the aerial parts of *C. ladanifer* and its concentration exhibits seasonal variation.⁴

References

1. Yilmaz, N., Türkekul, I., Bulut, S., et al. Fatty acid composition in ten mushroom species collected from middle Black Sea Region of Turkey. *Asian J. Chem.* **25(3)**, 1216-1220 (2006).
2. Harvey, D.J. and Tiffany, J.M. Identification of meibomian gland lipids by gas chromatography-mass spectrometry: Application to the meibomian lipids of the mouse. *J. Chromatogr.* **301(1)**, 173-187 (1984).
3. Jiao, Y.S., Yan, H., Ji, Z.J., et al. *Phyllobacterium sophorae* sp. nov., a symbiotic bacterium isolated from root nodules of *Sophora flavescens*. *Int. J. Syst. Evol. Microbiol.* **65(Pt 2)**, 399-406 (2015).
4. Guerreiro, O., Alves, S.P., Duarte, M.F., et al. *Cistus ladanifer* L. shrub is rich in saturated and branched chain fatty acids and their concentration increases in the Mediterranean dry season. *Lipids* **50(5)**, 493-501 (2015).

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, 03/07/2024

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