

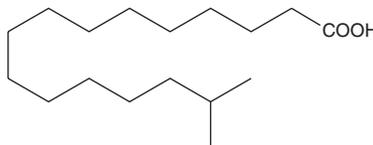
PRODUCT INFORMATION



15-methyl Palmitic Acid

Item No. 24814

CAS Registry No.: 1603-03-8
Formal Name: 15-methyl-hexadecanoic acid
Synonym: FA 17:0
MF: $C_{17}H_{34}O_2$
FW: 270.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

15-methyl Palmitic acid is supplied as a solid. A stock solution may be made by dissolving the 15-methyl palmitic acid in the solvent of choice, which should be purged with an inert gas. 15-methyl Palmitic acid is soluble in organic solvents such as ethanol, chloroform, and ethyl ether.

Description

15-methyl Palmitic acid is a methylated fatty acid that has been found in bacteria, bovine milk fat, one-humped camel (*C. dromedarius*) meat and fat, and in rabbit perirenal lipids.¹⁻⁴ It is also found in human breast milk and levels are decreased in mature breast milk compared to colostrum.⁵

References

1. Nickels, J.D., Chatterjee, S., Mostofian, B., et al. *Bacillus subtilis* lipid extract, a branched-chain fatty acid model membrane. *J. Phys. Chem. Lett.* **8(17)**, 4214-4217 (2017).
2. Henke, A., Westreicher-Kristen, E., Molkentin, J., et al. Effect of dietary quebracho tannin extract on milk fatty acid composition in cows. *J. Dairy Sci.* **100(8)**, 6229-6238 (2017).
3. Rawdah, T.N., El-Faer, M.Z., and Koreish, S.A. Fatty acid composition of the meat and fat of the one-humped camel *Camelus dromedarius*. *Meat Sci.* **37(1)**, 149-155 (1994).
4. López-Bote, C., Rey, A., Isabel, B., et al. Dietary fat reduces odd-numbered and branched-chain fatty acids in depot lipids of rabbits. *J. Sci. Food Agric.* **73(4)**, 517-524 (1999).
5. Jie, L., Qui, C., Sun, J., et al. The impact of lactation and gestational age on the composition of branched-chain fatty acids in human breast milk. *Food Funct.* **9(3)**, 1747-1754 (2018).

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, 02/29/2024

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