

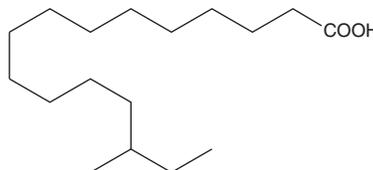
# PRODUCT INFORMATION



## 14-methyl Palmitic Acid

Item No. 24820

CAS Registry No.: 5918-29-6  
Formal Name: 14-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:  $\geq 4$  years



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

### Laboratory Procedures

14-methyl Palmitic acid is supplied as a solid. A stock solution may be made by dissolving the 14-methyl palmitic acid in the solvent of choice, which should be purged with an inert gas. 14-methyl Palmitic acid is soluble in organic solvents such as chloroform, ethanol, and ether.

### Description

14-methyl Palmitic acid is a methylated fatty acid that has been found in bacteria, bovine milk fat, Aegean jellyfish (*A. aurita*), and one-humped camel (*C. dromedarius*) meat and fat.<sup>1-4</sup> It is also found in human breast milk and levels are decreased in mature breast milk compared to colostrum.<sup>5</sup> 14-methyl Palmitic acid is produced in Florida manatee (*T. manatus latirostris*) liver in response to brevetoxin exposure.<sup>6</sup>

### 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. Kariotoglou, D.M. and Mastronicolis, S.K. Sphingophosphonolipids, phospholipids, and fatty acids from Aegean jellyfish *Aurelia aurita*. *Lipids* **36(11)**, 1255-1264 (2001).
4. 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).
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).
6. Wetzel, D.L., Reynolds, J.E., III, Sprinkle, J.M., et al. Fatty acid profiles as a potential lipidomic biomarker of exposure to brevetoxin for endangered Florida manatees (*Trichechus manatus latirostris*). *Sci. Total. Environ.* **408(24)**, 6124-6133 (2010).

#### 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