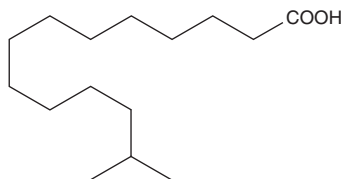


PRODUCT INFORMATION

13-methyl Myristic Acid

Item No. 24811

CAS Registry No.: 2485-71-4
Formal Name: 13-methyl-tetradecanoic acid
Synonyms: FA 15:0, Isopentadecanoic Acid, Subtilopentadecanoic Acid
MF: C₁₅H₃₀O₂
FW: 242.4
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

13-methyl Myristic acid is supplied as a solid. A stock solution may be made by dissolving the 13-methyl myristic acid in the solvent of choice, which should be purged with an inert gas. 13-methyl Myristic acid is soluble in ethanol, chloroform, and ether.

Description

13-methyl Myristic acid is a methylated fatty acid that has been found in dairy ewe rumen liquor, bovine milk, goat milk fat, the edible flesh of rainbow trout (*O. mykiss*), and *Flavobacterium*.¹⁻⁵ 13-methyl Myristic acid is also found in arable soil, and its levels decrease in response to contamination with the heavy metals cadmium, copper, nickel, and lead.⁶

References

1. Cappucci, A., Alves, S.P., Bessa, R.J.B., *et al.* Effect of increasing amounts of olive crude phenolic concentrate in the diet of dairy ewes on rumen liquor and milk fatty acid composition. *J. Dairy Sci.* **101**(6), (2018).
2. Brossillon, V., Reis, S.F., Moura, D.C., *et al.* Production, milk and plasma fatty acid profile, and nutrient utilization in Jersey cows fed flaxseed oil and corn grain with different particle size. *J. Dairy Sci.* **101**(3), 2127-2143 (2018).
3. Serment, A., Schmidely, P., Giger-Reverdin, S., *et al.* Effects of the percentage of concentrate on rumen fermentation, nutrient digestibility, plasma metabolites, and milk composition in mid-lactation goats. *J. Dairy Sci.* **94**(8), 3960-3972 (2011).
4. Calabretti, A., Cateni, F., Procida, G., *et al.* Influence of environmental temperature on composition of lipids in edible flesh of rainbow trout (*Oncorhynchus mykiss*). *J. Sci. Food Agric.* **83**(14), 1493-1498 (2003).
5. Moss, C.W. and Dees, S.B. Cellular fatty acids of *Flavobacterium meningosepticum* and *Flavobacterium* species group IIb. *J. Clin. Microbiol.* **8**(6), 772-774 (1978).
6. Frostegård, A., Tunlid, A., and Bååth, E. Phospholipid fatty acid composition, biomass, and activity of microbial communities from two soil types experimentally exposed to different heavy metals. *Appl. Environ. Microbiol.* **59**(11), 3605-3617 (1993).

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, 02/28/2024

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