

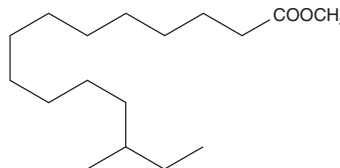
# PRODUCT INFORMATION



## 13-methyl Pentadecanoic Acid methyl ester

Item No. 24819

CAS Registry No.: 5487-50-3  
Formal Name: 13-methyl-pentadecanoic acid, methyl ester  
Synonym: Methyl 13-methyl Pentadecanoate Acid, SFE 17:0  
MF:  $C_{17}H_{34}O_2$   
FW: 270.5  
Purity:  $\geq 98\%$   
Supplied as: A solution in ethanol  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 2$  years



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

### Laboratory Procedures

13-methyl Pentadecanoic acid methyl ester 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 chloroform purged with an inert gas can be used.

### Description

13-methyl Pentadecanoic acid methyl ester is a methylated fatty acid methyl ester that has been found in the bioactive fraction of *A. malabarica* extract, *C. variabilis* microalgae, and baboon liver.<sup>1-3</sup>

### References

1. Kotha, P., Badri, K.R., Nagalapuram, R., *et al.* Anti-diabetic potential of the leaves of *Anisomeles malabarica* in streptozotocin induced diabetic rats. *Cell Physiol. Biochem.* **43**(4), 1689-1702 (2017).
2. Abdo, S.M., Ali, G.H., and El-Baz, F.K. Potential production of omega fatty acids from microalgae. *Int. J. Pharm. Sci.* **34**(2), 210-215 (2015).
3. Smith, A., Calder, A.G., Morrision, E.R., *et al.* Identification of branched chain fatty acids in baboon liver lipids. *Biomed. Mass Spectrom.* **6**(8), 345-346 (1979).

#### 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, 04/05/2024

#### 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