

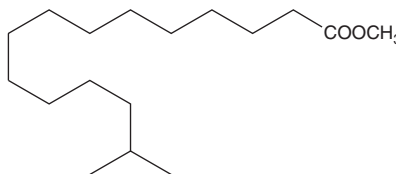
PRODUCT INFORMATION



14-methyl Pentadecanoic Acid methyl ester

Item No. 24813

CAS Registry No.: 5129-60-2
Formal Name: 14-methyl-pentadecanoic acid, methyl ester
Synonym: Methyl 14-methylpentadecanoate, SFE 17:0
MF: C₁₇H₃₄O₂
FW: 270.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

14-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

14-methyl Pentadecanoic acid methyl ester is a methylated fatty acid methyl ester that has been found in *S. zeai* sea sponges as the fatty acyl component of zeamide, *A. indica* leaf extract, and *C. vulgaris* and *H. pluvialis* microalgae.¹⁻³ It is a major component of the vancomycin-induced biofilm produced by vancomycin-resistant *S. aureus* (VRSA).⁴ 14-methyl Pentadecanoic acid methyl ester has been used as a standard for the quantification of 14-methyl pentadecanoic acid in various foods by GC-MS.⁵

References

1. Della Sala, G., Teta, R., Esposito, G., *et al.* Zeamide, a glycosylinositol phosphorylceramide with the novel core Arap(1β→6)ins motif from the marine sponge *Svenzea zeai*. *Molecules* **22(9)**, 1455 (2017).
2. Ravi, S., Shanmugam, B., Subbaiah, G.V., *et al.* Identification of food preservative, stress relief compounds by GC-MS and HR-LC/Q-TOF/MS; evaluation of antioxidant activity of *Acalypha indica* leaves methanolic extract (in vitro) and polyphenolic fraction (in vivo). *J. Food Sci. Technol.* **54(6)**, 1585-1596 (2017).
3. 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).
4. Mirani, Z.A. and Jamil, N. Role of extra-cellular fatty acids in vancomycin induced biofilm formation by vancomycin resistant *Staphylococcus aureus*. *Pak. J. Pharm. Sci.* **26(2)**, 383-389 (2013).
5. Thurnhofer, S. and Vetter, W. Application of ethyl esters and d₃-methyl esters as internal standards for the gas chromatographic quantification of transesterified fatty acid methyl esters in food. *J. Agric. Food Chem.* **54(9)**, 3209-3214 (2006).

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