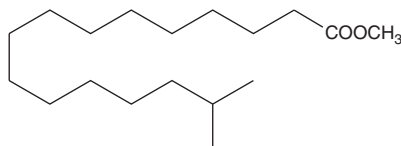


PRODUCT INFORMATION

15-methyl Palmitic Acid methyl ester

Item No. 24815

CAS Registry No.: 6929-04-0
Formal Name: 15-methyl-hexadecanoic acid, methyl ester
Synonyms: Methyl 15-methylhexadecanoate, SFE 18:0
MF: $C_{18}H_{36}O_2$
FW: 284.5
Purity: $\geq 98\%$
Supplied as: A liquid
Storage: -20°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 methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 15-methyl palmitic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 15-methyl Palmitic acid methyl ester is soluble in organic solvents such as ethanol and chloroform.

Description

15-methyl Palmitic 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, oolong tea (*C. sinensis*), and *P. sulcata* extract.¹⁻³ It has been used as a standard for the quantification of 15-methyl palmitic acid (Item No. 24814) in various foods.⁴

References

1. 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).
2. Zheng, L., Zhou, Y., Fu, X., *et al.* Does oolong tea (*Camellia sinensis*) made from a combination of leaf and stem smell more aromatic than leaf-only tea? Contribution of the stem to oolong tea aroma. *Food Chem.* **237**, 488-498 (2017).
3. Ari, F., Ulukaya, E., Oran, S., *et al.* Promising anticancer activity of a lichen, *Parmelia sulcata* Taylor, against breast cancer cell lines and genotoxic effect on human lymphocytes. *Cytotechnology* **67**(3), 531-543 (2015).
4. Thurnhofer, S. and Vetter, W. Application of ethyl esters and d_3 -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, 01/15/2026

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