

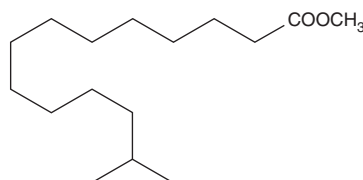
# PRODUCT INFORMATION



## 13-methyl Myristic Acid methyl ester

Item No. 24812

**CAS Registry No.:** 5129-59-9  
**Formal Name:** 13-methyl-tetradecanoic acid, methyl ester  
**Synonyms:** Methyl Isopentadecanoate,  
13-methyl Tetradecanoate,  
13-methyl Tetradecanoic Acid methyl ester,  
SFE 16:0  
**MF:** C<sub>16</sub>H<sub>32</sub>O<sub>2</sub>  
**FW:** 256.4  
**Purity:** ≥98%  
**Supplied as:** A 10 mg/ml solution in ethanol  
**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 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 13-methyl myristic acid methyl ester in these solvents is approximately 20 and 30 mg/ml, respectively.

### Description

13-methyl Myristic acid methyl ester is a methylated bacterial fatty acid methyl ester.<sup>1,2</sup> It has been used as a marker of bacteria in sewage sludges and of bacterial contaminants in pharmaceutical manufacturing by GC-MS.<sup>3,4</sup>

### References

1. Murugan, K., Prabu, R.V., Sangeetha, S., *et al.* Antiviral activity of *Cardiospermum Halicacabum* L. extract against coinfecting agents HIV and HBV. *J. Herbs Spices Med. Plants* **17(4)**, 403-418 (2011).
2. de Blas, E., Almendros, G., and Sanz, J. Molecular characterization of lipid fractions from extremely water-repellent pine and eucalyptus forest soils. *Geoderma* **206(2013)**, 75-84 (2013).
3. Jardé, E., Mansuy, L., and Faure, P. Organic markers in the lipidic fraction of sewage sludges. *Water Res.* **39(7)**, 1215-1232 (2005).
4. Olson, W.P., Groves, M.J., and Klegerman, M.E. Identifying bacterial contaminants in a pharmaceutical manufacturing facility by gas chromatographic fatty acid analysis. *Pharm. Tech.* **14(2)**, 32-36 (1990).

#### 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, 03/25/2026

#### 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