

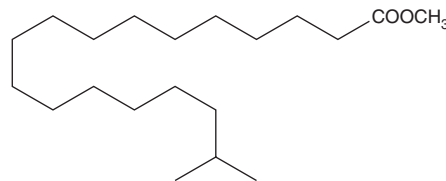
# PRODUCT INFORMATION



## 17-methyl Stearic Acid methyl ester

Item No. 24816

**CAS Registry No.:** 55124-97-5  
**Formal Name:** 17-methyl-octadecanoic acid, methyl ester  
**Synonyms** Methyl 17-methyl Octadecanoate Acid,  
SFE 20:0  
**MF:** C<sub>20</sub>H<sub>40</sub>O<sub>2</sub>  
**FW:** 312.5  
**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

17-methyl Stearic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 17-methyl stearic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 17-methyl Stearic acid methyl ester is soluble in organic solvents such as chloroform, ethanol, and ethyl ether, which should be purged with an inert gas.

### Description

17-methyl Stearic acid methyl ester is a methylated fatty acid methyl ester that has been found in mutton tallow.<sup>1</sup> It is a volatile component of roasted tigernut (*C. esculentus*) oil.<sup>2</sup>

### References

1. Chin, S.-T., Che Man, Y.B., Tan, C.P., *et al.* Rapid profiling of animal-derived fatty acids using fast GC × GC coupled to time-of-flight mass spectrometry. *J. Am. Oil Chem. Soc.* **86**(10), 949-958 (2009).
2. Lasekan, O. Volatile constituents of roasted tigernut oil (*Cyperus esculentus* L.). *J. Sci. Food Agric.* **93**(5), 1055-1061 (2013).

#### 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, 03/04/2024

#### 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