

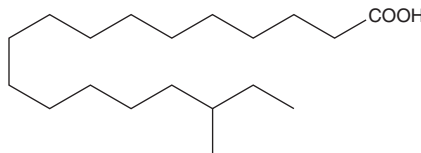
PRODUCT INFORMATION



16-methyl Stearic Acid

Item No. 36476

CAS Registry No.: 17001-28-4
Formal Name: 16-methyl-octadecanoic acid
Synonyms: Ai19:0, Anteisononadecanoic Acid, anteiso-19:0, FA 19:0
MF: $C_{19}H_{38}O_2$
FW: 298.5
Purity: $\geq 90\%$
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

16-methyl Stearic acid is supplied as a solid. A stock solution may be made by dissolving the 16-methyl stearic acid in the solvent of choice, which should be purged with an inert gas. 16-methyl Stearic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 16-methyl stearic acid in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

16-methyl Stearic acid is a methyl branched-chain fatty acid. It has been found in a variety of sea sponge species of the class Calcarea.¹

Reference

1. Schreiber, A., Wörheide, G., and Thiel, V. The fatty acids of calcareous sponges (Calcarea, Porifera). *Chem. Phys. Lipids* **143**(1-2), 29-37 (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, 04/07/2026

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