

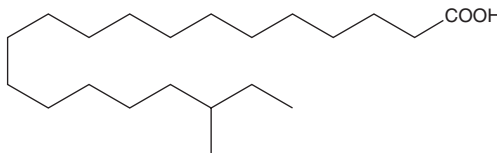
PRODUCT INFORMATION



18-methyl Eicosanoic Acid

Item No. 11009

CAS Registry No.: 36332-93-1
Formal Name: 18-methyl Eicosanoic Acid
Synonyms: Anteioheneicosanoic Acid, FA 21:0, 18-MEA
MF: $C_{21}H_{42}O_2$
FW: 326.6
Purity: $\geq 95\%$
Supplied as: A crystalline 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

18-methyl Eicosanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 18-methyl eicosanoic acid in the solvent of choice, which should be purged with an inert gas. 18-methyl Eicosanoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 18-methyl eicosanoic acid in these solvents is approximately 10, 30, and 50 mg/ml, respectively.

Description

18-methyl Eicosanoic acid is a main component of the outer layer of the epicuticle of human hair and endows hydrophobicity to the outer hair surface. It is used as an additive to hair cosmetics due to its desirable conditioning effects.¹

Reference

1. Cundy, D.J. and Gurr, P.A. An improved synthesis of (\pm) 18 methyl eicosanoic acid. *Org. Prep. Proced. Int.* **32**(5), 461-468 (2011).

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, 02/21/2024

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