

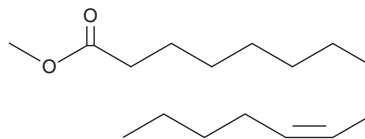
PRODUCT INFORMATION



10(Z)-Pentadecenoic Acid methyl ester

Item No. 26729

CAS Registry No.: 90176-52-6
Formal Name: 10Z-pentadecenoic acid, methyl ester
Synonyms: C15:1 (cis-10) methyl ester, cis-10-Pentadecenoic Acid methyl ester, SFE 16:1
MF: C₁₆H₃₀O₂
FW: 254.4
Purity: ≥99%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

10(Z)-Pentadecenoic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 10(Z)-pentadecenoic acid methyl ester in the solvent of choice. 10(Z)-Pentadecenoic acid methyl ester is soluble in organic solvents such as chloroform, hexane, and ethyl ether, which should be purged with an inert gas.

Description

10(Z)-Pentadecenoic acid methyl ester is a methyl ester form of the fatty acid 10(Z)-pentadecenoic acid (Item No. 22362). It has been used as a standard for the quantification of pentadecenoic acid in vernix caseosa.¹

Reference

1. Hauff, S. and Vetter, W. Exploring the fatty acids of vernix caseosa in form of their methyl esters by off-line coupling of non-aqueous reversed phase high performance liquid chromatography and gas chromatography coupled to mass spectrometry. *J. Chromatogr. A* **1217(52)**, 8270-8278 (2010).

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, 02/13/2024

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