

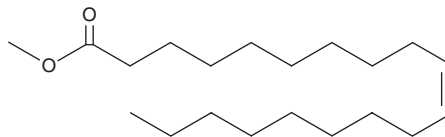
PRODUCT INFORMATION



10(Z)-Nonadecenoic Acid methyl ester

Item No. 26733

CAS Registry No.: 19788-74-0
Formal Name: 10Z-nonadecenoic acid, methyl ester
Synonyms: Methyl 10Z-Nonadecenoate,
cis-10-Nonadecenoic Acid methyl ester,
SFE 20:1
MF: C₂₀H₃₈O₂
FW: 310.5
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

10(Z)-Nonadecenoic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 10(Z)-nonadecenoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 10(Z)-Nonadecenoic acid methyl ester is soluble in chloroform.

Description

10(Z)-Nonadecenoic acid methyl ester is a fatty acid methyl ester.¹ It has been used as a standard for the quantification of *cis*-nonadecenoic acid (Item No. 19749) in chromatography applications.^{1,2}

References

1. Stránský, K., Jursík, T., Vitek, A., *et al.* An improved method of characterizing fatty acids by equivalent chain length values. *J. High Resolut. Chrom.* **15(11)**, 730-740 (1992).
2. Vine, J. Analysis of fatty acid methyl esters by high-resolution gas chromatography-chemical ionisation mass spectrometry. *J. Chromatogr. A.* **196(3)**, 415-424 (1980).

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