

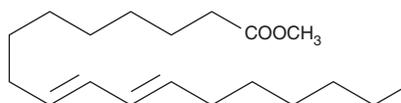
PRODUCT INFORMATION



9(E),11(E)-Conjugated Linoleic Acid methyl ester

Item No. 24580

CAS Registry No.: 13038-47-6
Formal Name: 9E,11E-octadecadienoic acid, methyl ester
Synonyms: 9E,11E-CLA methyl ester
Methyl 9(E),11(E)-Octadecadienoate,
SFE 19:2
MF: C₁₉H₃₄O₂
FW: 294.5
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 232 nm
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

9(E),11(E)-Conjugated linoleic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 9(E),11(E)-conjugated linoleic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 9(E),11(E)-Conjugated linoleic acid methyl ester is soluble in ethanol, methanol, and chloroform.

Description

9(E),11(E)-Conjugated linoleic acid methyl ester has been found in thermally stressed cooking oils and may be used as a marker of adulteration of olive oils with lower quality oils.¹

Reference

1. Saba, A., Mazzini, F., Raffaelli, A., *et al.* Identification of 9(E),11(E)-18:2 fatty acid methyl ester at trace level in thermal stressed olive oils by GC coupled to acetonitrile CI-MS and CI-MS/MS, a possible marker for adulteration by addition of deodorized olive oil. *J. Agric. Food Chem.* **53(12)**, 4867-4872 (2005).

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, 04/01/2024

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