

PRODUCT INFORMATION



10(E),12(Z)-Conjugated Linoleic Acid methyl ester

Item No. 24584

CAS Registry No.: 21870-97-3
Formal Name: 10E,12Z-octadecadienoic acid, methyl ester
Synonyms: 10E,12Z-CLA methyl ester,
Methyl 10(E),12(Z)-Octadecadienoate methyl ester,
SFE 19:2

MF: C₁₉H₃₄O₂

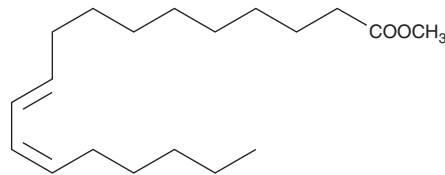
FW: 294.5

Purity: ≥98%

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(E),12(Z)-Conjugated linoleic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 10(E),12(Z)-conjugated linoleic acid methyl ester in the solvent of choice. 10(E),12(Z)-Conjugated linoleic acid methyl ester is soluble in ethanol, methanol, hexane, and chloroform.

Description

10(E),12(Z)-Conjugated linoleic acid methyl ester has been used as a standard in the quantification of 10(E),12(Z)-conjugated linoleic acid from *L. plantarum* culture samples.¹

Reference

1. Sun, C., Black, B.A., Zhao, Y.Y., *et al.* Identification of conjugated linoleic acid (CLA) isomers by silver ion-liquid chromatography/in-line ozonolysis/mass spectrometry (Ag⁺-LC/O₃-MS). *Anal. Chem.* **85(15)**, 7345-7352 (2013).

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, 10/16/2018

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