

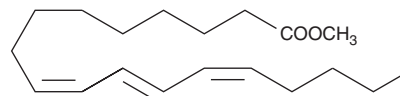
PRODUCT INFORMATION



9(Z),11(E),13(Z)-Octadecatrienoic Acid methyl ester

Item No. 24807

CAS Registry No.: 95497-55-5
Formal Name: 9Z,11E,13Z-octadecatrienoic acid, methyl ester
Synonyms: Methyl Punicate, Punicic Acid methyl ester, SFE 19:3
MF: C₁₉H₃₂O₂
FW: 292.5
Purity: ≥95%
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(Z),11(E),13(Z)-Octadecatrienoic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the 9(Z),11(E),13(Z)-octadecatrienoic acid methyl ester in the solvent of choice. 9(Z),11(E),13(Z)-Octadecatrienoic acid methyl ester is soluble in organic solvents such as hexane, ethanol, methanol, and chloroform, which should be purged with an inert gas.

Description

9(Z),11(E),13(Z)-Octadecatrienoic acid methyl ester is an isomer of 9(Z),11(E),13(E)-octadecatrienoic acid methyl ester (Item No. 10008333) and the methyl ester form of 9(Z),11(E),13(Z)-octadecatrienoic acid (Item No. 26057). It has been used as a standard for the quantification of 9(Z),11(E),13(Z)-octadecatrienoic acid in wild growing pomegranate (*P. granatum*) seed oil.¹

Reference

1. Đurđević, S., Šavikin, K., Živković, J., *et al.* Antioxidant and cytotoxic activity of fatty oil isolated by supercritical fluid extraction from microwave pretreated seeds of wild growing *Punica granatum* L. *J. Supercrit. Fluids* **133**(1), 225-232 (2018).

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