

# PRODUCT INFORMATION



## Elaidic Acid methyl ester

Item No. 26870

**CAS Registry No.:** 1937-62-8  
**Formal Name:** 9E-octadecenoic acid, methyl ester  
**Synonyms:** C18:1 (*trans*-9) methyl ester, Methyl Elaidate, Methyl 9(E)-Octadecenoate, Methyl *trans*-9-Octadecenoate, SFE 19:1, *trans*-9-Octadecenoic Acid methyl ester, *trans*-Oleic Acid methyl ester

**MF:** C<sub>19</sub>H<sub>36</sub>O<sub>2</sub>

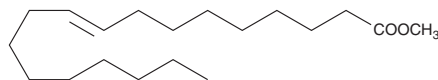
**FW:** 296.5

**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

Elaidic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the elaidic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Elaidic acid methyl ester is soluble in methanol and chloroform.

### Description

Elaidic acid methyl ester is an esterified form of elaidic acid (Item No. 90250) that has been found in biodiesel produced by the microalga *Botryococcus*.<sup>1</sup> It has also been found in Iranian olive fruit oil and certain cultivars of *O. sativa* black rice bran.<sup>2,3</sup>

### References

1. Ashokkumar, V., Agila, E., Pandian, S., *et al.* Optimization and characterization of biodiesel production from microalgae *Botryococcus* grown at semi-continuous system. *Energy Convers. Manag.* **88**, 936-946 (2014).
2. Kono, E., Abbasi, A., Parastar, H., *et al.* Analysis of olive fruit essential oil: Application of gas chromatography-mass spectrometry combined with chemometrics. *Int. J. Food Prop.* **18(2)**, 1532-2386 (2015).
3. Arjinajarn, P., Pongchaidecha, A., Chueakula, N., *et al.* Riceberry bran extract prevents renal dysfunction and impaired renal organic anion transporter 3 (Oat3) function by modulating the PKC/Nrf2 pathway in gentamicin-induced nephrotoxicity in rats. *Phytomedicine* **23(14)**, 316-331 (2016).

#### 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, 03/04/2024

#### 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