

PRODUCT INFORMATION



10(Z)-Heptadecenoic Acid methyl ester

Item No. 26869

CAS Registry No.: 75190-82-8

Formal Name: 10Z-heptadecenoic acid, methyl ester

Synonyms: C17:1 methyl ester, Methyl Heptadecenoate, cis-10-Heptadecenoic Acid methyl ester, SFE 18:1

MF: $C_{18}H_{34}O_2$

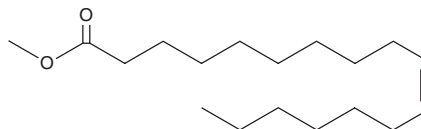
FW: 282.5

Purity: $\geq 98\%$

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

Description

10(Z)-Heptadecenoic acid methyl ester is an ester form of cis-10-heptadecenoic acid (Item No. 19748). It is a minor fatty acid methyl ester (FAME) constituent of biodiesel.¹

Reference

1. Khoury, R.R., Ebrahimi, D., Hejazi, L., *et al.* Degradation of fatty acid methyl esters in biodiesels exposed to sunlight and seawater. *Fuel* **90**(8), 2677-2683 (2011).

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, 01/16/2026

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