

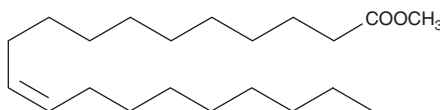
PRODUCT INFORMATION



11(Z)-Eicosenoic Acid methyl ester

Item No. 21332

CAS Registry No.: 2390-09-2
Formal Name: 11Z-eicosenoic acid, methyl ester
Synonyms: *cis*-11-Eicosenoic Acid methyl ester, Gondoic Acid methyl ester, SFE 21:1
MF: C₂₁H₄₀O₂
FW: 324.5
Purity: ≥98%
Supplied as: A 250 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

11(Z)-Eicosenoic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 11(Z)-eicosenoic acid methyl ester in these solvents is approximately 100 mg/ml.

Description

11(Z)-Eicosenoic acid methyl ester is a natural derivative of the monounsaturated 20-carbon fatty acid 11(Z)-eicosenoic acid (Item No. 20606).¹ Methyl esters of fatty acids are used in formulations for nutritional supplements.

Reference

1. Menichini, F., Losi, L., Bonesi, M., *et al.* Chemical profiling and *in vitro* biological effects of *Cardiospermum halicacabum* L. (Sapindaceae) aerial parts and seeds for applications in neurodegenerative disorders. *J. Enzyme Inhib. Med. Chem.* **29**(5), 677-685 (2014).

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, 03/25/2026

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