

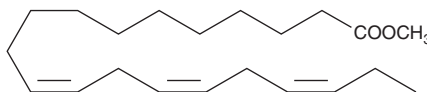
PRODUCT INFORMATION



11(Z),14(Z),17(Z)-Eicosatrienoic Acid methyl ester

Item No. 17622

CAS Registry No.: 62472-96-2
Formal Name: 11Z,14Z,17Z-eicosatrienoic acid, methyl ester
Synonyms: C21:3 (cis-11,cis-14,cis-17) methyl ester, SFE 21:3
MF: C₂₁H₃₆O₂
FW: 320.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

11(Z),14(Z),17(Z)-Eicosatrienoic 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, dimethyl formamide (DMF), and 0.1 M Na₂CO₃ purged with an inert gas can be used. The solubility of 11(Z),14(Z),17(Z)-eicosatrienoic acid methyl ester in DMSO and DMF is approximately 100 mg/ml and approximately 1 mg/ml in 0.1 M Na₂CO₃.

Description

11(Z),14(Z),17(Z)-Eicosatrienoic acid methyl ester is an esterified form of 11(Z),14(Z),17(Z)-eicosatrienoic acid (Item No. 90192).¹ It has been found in the lipid-producing microalgae *Nannochloropsis*. 11(Z),14(Z),17(Z)-Eicosatrienoic acid methyl ester also has been found in diesel-contaminated soils and residential wastewater.^{2,3}

References

1. Ahamed, T.S., Brindhadevi, K., Krishnan, R., *et al.* *In vivo* detection of triacylglycerols through Nile red staining and quantification of fatty acids in hyper lipid producer *Nannochloropsis* sp. cultured under adequate nitrogen and deficient nitrogen condition. *Fuel* **322**, 124179 (2022).
2. Ko, M.-S. and Lee, S. Discrimination methods for diesel origin by analyzing fatty acid methyl ester (FAME) composition in diesel-contaminated soil. *Sci. Rep.* **11(1)**, 16245 (2021).
3. Katam, K. and Bhattacharyya, D. Comparative study on treatment of kitchen wastewater using a mixed microalgal culture and an aerobic bacterial culture: Kinetic evaluation and FAME analysis. *Environ. Sci. Pollut. Res. Int.* **25(21)**, 20732-20742 (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