

PRODUCT INFORMATION



Arachidic Acid methyl ester

Item No. 26724

CAS Registry No.: 1120-28-1
Formal Name: eicosanoic acid, methyl ester
Synonyms: C20:0 methyl ester,
Eicosanoic Acid methyl ester,
Methyl Arachidate,
Methyl Eicosanoate, SFE 21:0

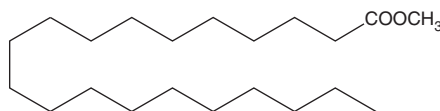
MF: C₂₁H₄₂O₂
FW: 326.6

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Arachidic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the arachidic acid methyl ester in the solvent of choice. Arachidic acid methyl ester is soluble in organic solvents such as ethanol and dimethyl formamide, which should be purged with an inert gas. The solubility of arachidic acid methyl ester in these solvents is approximately 0.1 and 2 mg/ml, respectively.

Arachidic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, arachidic acid methyl ester should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Arachidic acid methyl ester has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Arachidic acid methyl ester is an esterified form of arachidic acid (Item Nos. 9000339 | 21906). It is a minor constituent of biodiesel generated from acid-catalyzed or subcritical water-subcritical methanol (SCW-SCM) *in situ* transesterification of dry activated sludge or sludge oil, respectively.¹

Reference

1. Huynh, L.H., Tran Nguyen, P.L., Ho, Q.P., *et al.* Catalyst-free fatty acid methyl ester production from wet activated sludge under subcritical water and methanol condition. *Bioresour. Technol.* **123**, 112-116 (2012).

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