

PRODUCT INFORMATION

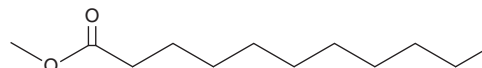


Undecanoic Acid methyl ester

Item No. 26720

CAS Registry No.: 1731-86-8
Synonyms: C11:0 Methyl ester,
Hendecanoic Acid methyl ester,
Methyl Undecanoate, SFE 12:0,
Undecylic Acid methyl ester

MF: C₁₂H₂₄O₂
FW: 200.3
Purity: ≥98%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Undecanoic acid methyl ester is supplied as a neat oil. A stock solution may be made by dissolving the undecanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Undecanoic acid methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of undecanoic acid methyl ester in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml in DMSO.

Undecanoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, undecanoic acid methyl ester should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Undecanoic acid methyl ester has a solubility of approximately 0.25 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Undecanoic acid methyl ester is an ester form of undecanoic acid (Item No. 19721). Undecanoic acid methyl ester is a minor component of biodiesel formed by the transesterification of freeze-dried municipal secondary sewage sludge.¹

Reference

1. Kumar, M., Ghosh, P., Khosla, K., *et al.* Biodiesel production from municipal secondary sludge. *Bioresour. Technol.* **216**, 165-171 (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, 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