

PRODUCT INFORMATION



Docosanoic Acid methyl ester

Item No. 26866

CAS Registry No.: 929-77-1

Synonyms: Behenic Acid methyl ester,
C22:0 methyl ester,
Methyl Behenate,
Methyl Docosanoate,
NSC 158426, SFE 23:0

MF: C₂₃H₄₆O₂

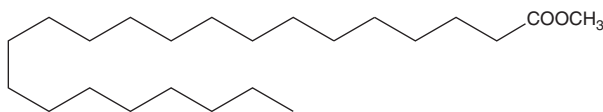
FW: 354.6

Purity: ≥95%

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

Docosanoic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the docosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Docosanoic acid methyl ester is soluble in the organic solvent ethanol at a concentration of approximately 0.5 mg/ml.

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

Description

Docosanoic acid methyl ester is an ester form of docosanoic acid (Item No. 9000338). It has been found in soybean, corn, sunflower, and canola oils, as well as beef fat, used in the production of biodiesel.¹

Reference

1. Mogollón, N.G.S., Ribeiro, F.A.L., Poppi, R.J., *et al.* Exploratory analysis of biodiesel by combining comprehensive two-dimensional gas chromatography and multiway principal component analysis. *J. Braz. Chem. Soc.* **28**(5), 740-746 (2017).

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, 02/14/2024

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