

PRODUCT INFORMATION

Hexacosanoic Acid methyl ester

Item No. 26725

CAS Registry No.: 5802-82-4

Synonyms: C26:0 methyl ester,
Ceric Acid methyl ester,
Cerotic Acid methyl ester,
Methyl hexacosanoate, SFE 27:0

MF: $C_{27}H_{54}O_2$

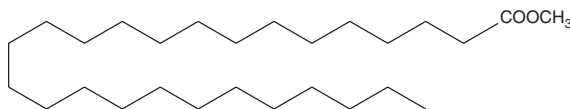
FW: 410.7

Purity: $\geq 98\%$

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

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

Hexacosanoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hexacosanoic acid methyl ester should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Hexacosanoic 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

Hexacosanoic acid methyl ester is a fatty acid methyl ester. It has been used as a standard for the detection of hexacosanoic acid (Item No. 13354) in beeswax.¹

Reference

1. Du, J., Li, C., and Qiu, F.-p. GC-MS analysis of fatty acids in three kinds of natural beeswax. *Shipin Keji* 39(8), 268-271 (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, 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