

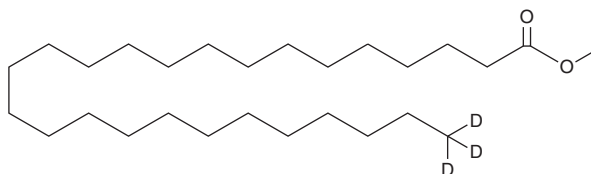
PRODUCT INFORMATION



Hexacosanoic Acid methyl ester-d₃

Item No. 28085

CAS Registry No.: 2743078-84-2
Formal Name: hexacosanoic-26,26,26-d₃ acid, methyl ester
Synonyms: C26:0 methyl ester-d₃,
Ceric Acid methyl ester-d₃,
Cerotic Acid methyl ester-d₃, SFE 27:0-d₃
MF: C₂₇H₅₁D₃O₂
FW: 413.7
Chemical Purity: ≥98% (Hexacosanoic Acid methyl ester)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A 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-d₃ is intended for use as an internal standard for the quantification of hexacosanoic acid methyl ester (Item No. 26725) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Hexacosanoic acid methyl ester-d₃ is supplied as a solid. A stock solution may be made by dissolving the hexacosanoic acid methyl ester-d₃ in the solvent of choice, which should be purged with an inert gas. Hexacosanoic acid methyl ester-d₃ is soluble in the organic solvent ethanol at a concentration of approximately 0.5 mg/ml.

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, 02/14/2024

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