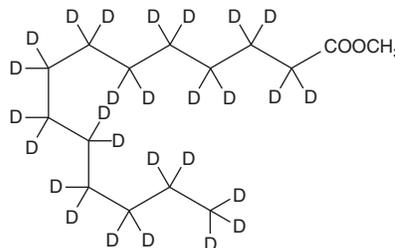


PRODUCT INFORMATION



Myristic Acid methyl ester-d₂₇ Item No. 28594

CAS Registry No.: 434349-93-6
Formal Name: tetradecanoic-d₂₇ acid, methyl ester
Synonyms: Methyl Myristate-d₂₇,
Methyl Tetradecanoate-d₂₇, SFE 15:0-d₂₇
MF: C₁₅H₃D₂₇O₂
FW: 269.6
Chemical Purity: ≥98% (Myristic Acid methyl ester)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₂₇); ≤1% d₀
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

Myristic acid methyl ester-d₂₇ is intended for use as an internal standard for the quantification of myristic acid methyl ester (Item No. 9001867) 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).

Myristic acid methyl ester-d₂₇ is supplied as a neat oil. A stock solution may be made by dissolving the myristic acid methyl ester-d₂₇ in the solvent of choice, which should be purged with an inert gas. Myristic acid methyl ester-d₂₇ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of myristic acid methyl ester-d₂₇ in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

Description

Myristic acid methyl ester is an esterified version of the free acid, which is less water soluble but more amenable for the formulation of myristate-containing diets and dietary supplements.

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