

PRODUCT INFORMATION



Medium-chain Saturated Fatty Acid Methyl Ester Mixture 1

Item No. 28646

Supplied as: A liquid

Storage: -20°C

Stability: ≥2 years

Special Conditions: This standard is accurately prepared by gravimetric technique (+/-0.5%)

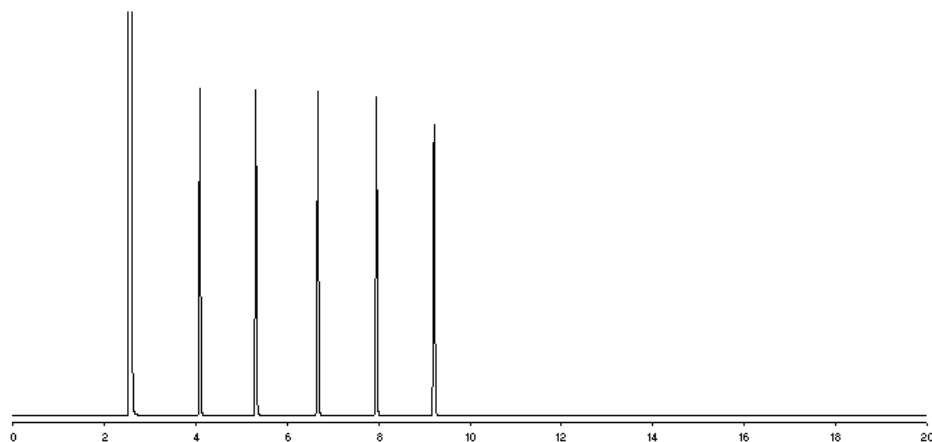
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Medium-chain Saturated Fatty Acid Methyl Ester Mixture 1 contains five saturated medium-chain fatty acid methyl esters (FAMES). It is intended for use as a reference standard for the quantification of these medium-chain fatty acids in various sample types by GC-FID or GC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation.

Included are:

Elution Order*	Item Name	% Concentration by Weight
1	Octanoic Acid methyl ester	20
2	Decanoic Acid methyl ester	20
3	Lauric Acid methyl ester	20
4	Myristic Acid methyl ester	20
5	Palmitic Acid methyl ester	20



*GC Column Used: SP-2330 or RTX-2330, 30 m x 0.25 mm x 0.2 µm

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, 04/27/2026

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