

PRODUCT INFORMATION

AOCS Medium- and Long-chain Fatty Acid Methyl Ester Standard Mixture 1

Item No. 28669

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.

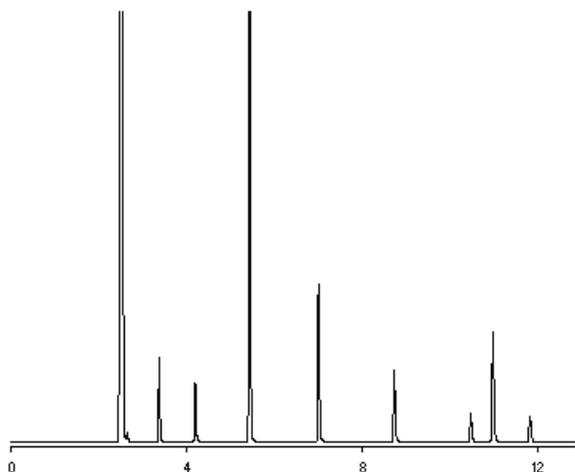
Description and Data

AOCS Medium- and Long-chain Fatty Acid Methyl Ester Standard Mixture 1 contains eight saturated and mono- and poly-unsaturated fatty acid methyl esters. It is intended for use as a reference standard for the detection of fatty acids in coconut, palm kernel, babassu, and ouri-ouri oils, among others, by GC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with nitrogen/argon to prevent lipid oxidation.

Included are:

Elution Order*	Item Name	% Concentration by Weight
1	Octanoic Acid methyl ester	7
2	Decanoic Acid methyl ester	5
3	Lauric Acid methyl ester	48
4	Myristic Acid methyl ester	15
5	Palmitic Acid methyl ester	7
6	Stearic Acid methyl ester	3
7	Oleic Acid methyl ester	12
8	Linoleic Acid methyl ester	3

*GC Column Used: SP-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, 08/16/2022

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