

PRODUCT INFORMATION



AOCS Long-chain Fatty Acid Methyl Ester Standard Mixture 3

Item No. 29366

Supplied as: A solution in dichloromethane
Storage: -20°C
Stability: ≥2 years
Concentration: 50 mg/ml
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.

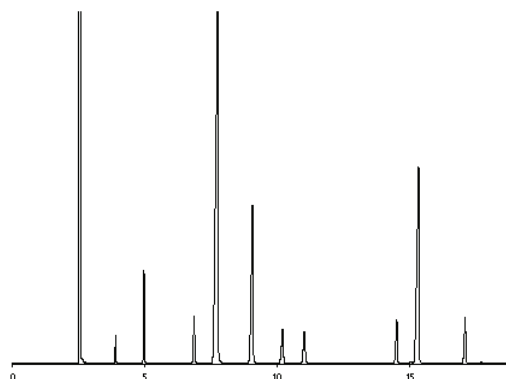
Description and Data

AOCS Long-Chain Fatty Acid Methyl Ester Standard Mixture 3 contains ten 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 peanut, rapeseed, and mustard seed 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	% Concentrations by Weight
1	Myristic Acid methyl ester	1
2	Palmitic Acid methyl ester	4
3	Stearic Acid methyl ester	3
4	Oleic Acid methyl ester	45
5	Linoleic Acid methyl ester	15
6	Arachidic Acid methyl ester	3
7	α-Linolenic Acid methyl ester	3
8	Docosanoic Acid methyl ester	3
9	13(Z)-Docosenoic Acid methyl ester	20
10	Lignoceric 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, 10/04/2022

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