

PRODUCT INFORMATION



AOCS Long-chain Fatty Acid Methyl Ester Standard Mixture 4 Item No. 28668

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.

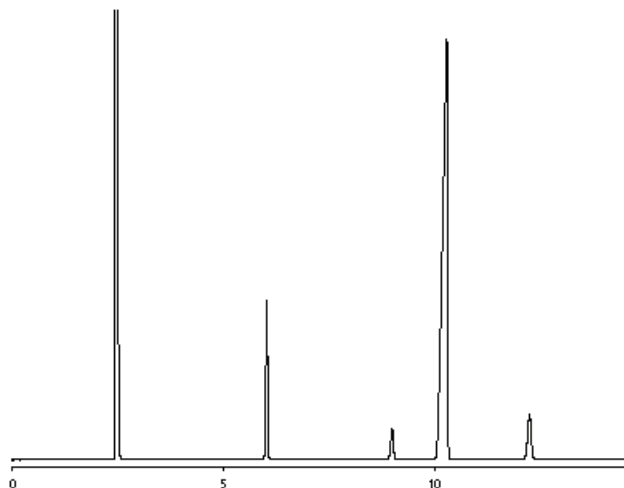
Description and Data

AOCS Long-chain Fatty Acid Methyl Ester Standard Mixture 4 contains two saturated, one monounsaturated, and one polyunsaturated fatty acid methyl ester. It is intended for use as a reference standard for the quantification of these long-chain fatty acids in olive, teaseed, and neatsfoot oils, among others, 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	Palmitic Acid methyl ester	11.0
2	Stearic Acid methyl ester	3.0
3	Oleic Acid methyl ester	80.0
4	Linoleic Acid methyl ester	6.0

*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, 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