

PRODUCT INFORMATION



Comprehensive Fatty Acid Methyl Ester Mixture 3

Item No. 29371

Supplied as: A solution in heptane

Storage: -20°C

Stability: ≥1 year

Concentration: 15.5 mg/ml

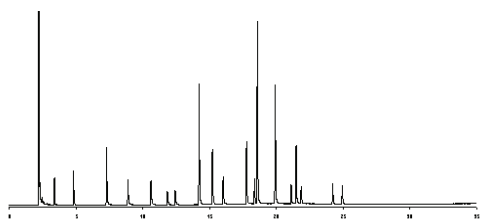
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Comprehensive Fatty Acid Methyl Ester Mixture 3 contains eleven saturated, six monounsaturated, and two polyunsaturated fatty acid methyl esters (FAMES). It is intended for use as a reference standard for the quantification of these 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	mg/ml
1	Octanoic Acid methyl ester	0.3
2	Decanoic Acid methyl ester	0.5
3	Lauric Acid methyl ester	1.0
4	Tridecanoic Acid methyl ester	0.5
5	Myristic Acid methyl ester	0.5
6	Myristoleic Acid methyl ester	0.3
7	Pentadecanoic Acid methyl ester	0.3
8	Palmitic Acid methyl ester	2.0
9	Palmitoleic Acid methyl ester	1.0
10	Heptadecanoic Acid methyl ester	0.5
11	Stearic Acid methyl ester	1.0
12	Elaidic Acid methyl ester	0.4
13	Oleic Acid methyl ester	3.0
14	Linoleic Acid methyl ester	2.0
15	Arachidic Acid methyl ester	0.3
16	α-Linolenic Acid methyl ester	1.0
17	11(Z)-Eicosenoic Acid methyl ester	0.3
18	Docosanoic Acid methyl ester	0.3
19	13(Z)-Docosenoic Acid methyl ester	0.3



*GC Column Used: 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, 04/29/2026

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