

PRODUCT INFORMATION

Comprehensive Fatty Acid Methyl Ester Mixture 2

Item No. 29335

Supplied as: A solution in dichloromethane

Storage: -20°C

Stability: ≥1 year

Concentration: 30 mg/ml

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

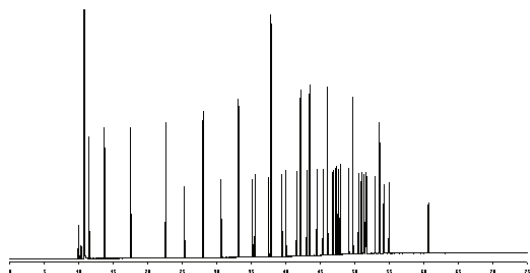
Description and Data

Comprehensive Fatty Acid Methyl Ester Mixture 2 contains 17 saturated, nine monounsaturated, and eleven 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	% Concentration by Weight	Elution Order*	Item Name	% Concentration by Weight
1	Butyric Acid methyl ester	4	20	Linoleic Acid methyl ester	2
2	Hexanoic Acid methyl ester	4	21	Arachidic Acid methyl ester	4
3	Octanoic Acid methyl ester	4	22	γ-Linolenic Acid methyl ester	2
4	Decanoic Acid methyl ester	4	23	11(Z)-Eicosenoic Acid methyl ester	2
5	Undecanoic Acid methyl ester	2	24	α-Linolenic Acid methyl ester	2
6	Lauric Acid methyl ester	4	25	Heneicosanoic Acid methyl ester	2
7	Tridecanoic Acid methyl ester	2	26	11(Z),14(Z)-Eicosadienoic Acid methyl ester	2
8	Myristic Acid methyl ester	4	27	Docosanoic Acid methyl ester	4
9	Myristoleic Acid methyl ester	2	28	Dihomo-γ-Linolenic Acid methyl ester	2
10	Pentadecanoic Acid methyl ester	2	29	13(Z)-Docosanoic Acid methyl ester	2
11	10(Z)-Pentadecenoic Acid methyl ester	2	30	11(Z),14(Z),17(Z)-Eicosatrienoic Acid methyl ester	2
12	Palmitic Acid methyl ester	6	31	Tricosanoic Acid methyl ester	2
13	Palmitoleic Acid methyl ester	2	32	Arachidonic Acid methyl ester	2
14	Heptadecanoic Acid methyl ester	2	33	13(Z),16(Z)-Docosadienoic Acid methyl ester	2
15	10(Z)-Heptadecenoic Acid methyl ester	2	34	Lignoceric Acid methyl ester	4
16	Stearic Acid methyl ester	4	35	Eicosapentaenoic Acid methyl ester	2
17	Elaidic Acid methyl ester	2	36	Nervonic Acid methyl ester	2
18	Oleic Acid methyl ester	4	37	Docosahexaenoic Acid methyl ester	2
19	Linoelaidic Acid methyl ester	2			

*GC Column Used: SP-2560, 100 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