

PRODUCT INFORMATION



Long-chain Fatty Acid Methyl Ester Mixture (porcine ovary)

Item No. 28662

Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years

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

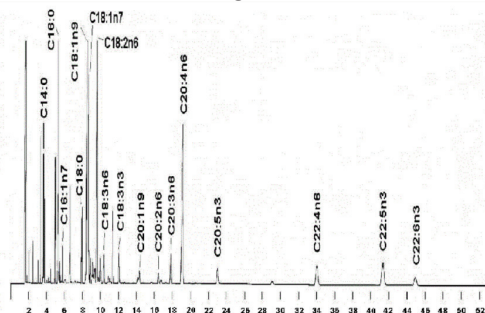
Description and Data

Long-chain fatty acid methyl ester mixture contains 14 omega fatty acids and three saturated fatty acids isolated from porcine ovaries and supplied as methyl esters. It is intended for use as a reference standard for the detection of these long-chain fatty acids from various sample types.

Included are:

Elution Order*	Item Name
1	Myristic Acid methyl ester
2	Palmitic Acid methyl ester
3	Palmitoleic Acid methyl ester
4	Stearic Acid methyl ester
5	Oleic Acid methyl ester
6	Vaccenic Acid methyl ester
7	Linoleic Acid methyl ester
8	Jacaric Acid methyl ester
9	Linolenic Acid methyl ester
10	11(Z)-Eicosenoic Acid methyl ester
11	11(Z),14(Z)-Eicosadienoic Acid methyl ester
12	Dihomo-γ-Linolenic Acid methyl ester
13	Arachidonic Acid methyl ester
14	Eicosapentaenoic Acid methyl ester
15	7(Z),10(Z),13(Z),16(Z)-Docosatetraenoic Acid methyl ester
16	Docosapentaenoic Acid methyl ester
17	Docosahexaenoic Acid methyl ester

*GC Column Used: Omegawax 250, 30 m x 0.25 mm x 0.25 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.

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