

CERTIFICATE of ANALYSIS



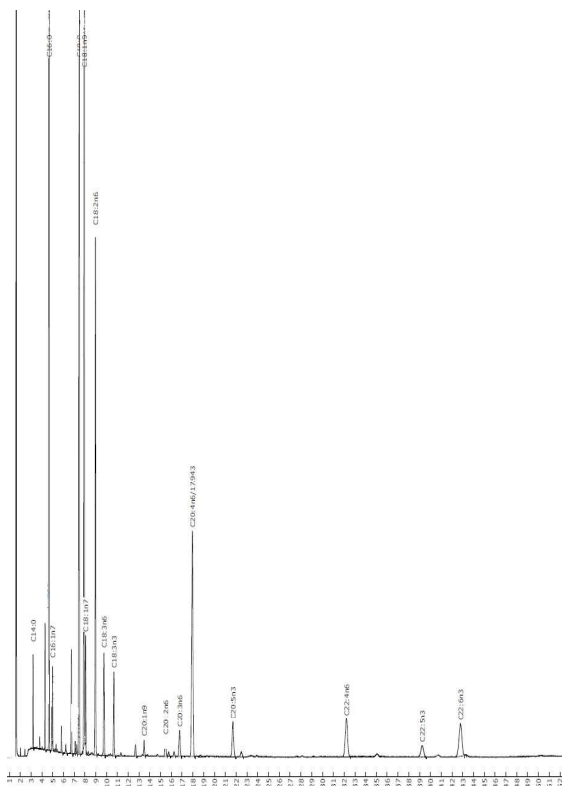
Long-chain Fatty Acid Methyl Ester Mixture (porcine ovary)

Item No.:	28662
Batch No.:	0569031
Expiry Date:	04SEP2027 (valid from date of certification)
Supplied as:	A mixture of 14 omega fatty acids and three saturated fatty acids.
Contents:	See table
Volume per Ampule:	Not less than 1 ml. Ampules are overfilled. It is advised that labs use measured volumes.
Storage & Handling:	Store upright and unopened at -20°C. Warm to room temperature prior to opening.

Description

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

Quality Information



Approval: *J. Albertson*

Certification Date: 17MAY2019

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiry date when stored unopened as recommended.

CERTIFICATE of ANALYSIS



Quality Control Table

Elution Order	Retention Time	Carbon Number	Item Name
1	3.186	C14:0	Myristic Acid methyl ester
2	4.660	C16:0	Palmitic Acid methyl ester
3	4.980	C16:1n7	Palmitoleic Acid methyl ester
4	7.453	C18:0	Stearic Acid methyl ester
5	7.926	C18:1n9	Oleic Acid methyl ester
6	8.043	C18:1n7	Vaccenic Acid methyl ester
7	8.960	C18:2n6	Linoleic Acid methyl ester
8	9.750	C18:3n6	Jacaric Acid methyl ester
9	10.670	C18:3n3	Linolenic Acid methyl ester
10	13.470	C20:1n9	11(Z)-Eicosenoic Acid methyl ester
11	15.453	C20:2n6	11(Z),14(Z)-Eicosadienoic Acid methyl ester
12	16.756	C20:3n6	Dihomo-γ-Linolenic Acid methyl ester
13	17.943	C20:4n6	Arachidonic Acid methyl ester
14	21.700	C20:5n3	Eicosapentaenoic Acid methyl ester
15	32.210	C22:4n6	7(Z),10(Z),13(Z),16(Z)-Docosatetraenoic Acid methyl ester
16	39.233	C22:5n3	Docosapentaenoic Acid methyl ester
17	42.790	C22:6n3	Docosahexaenoic Acid methyl ester
Column:	Omegawax 250, 30 m x 0.25 mm x 0.25 μm		

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	17MAY2019	Initial version
02	08SEP2025	Updated quality information, quality control table, and expiry date

Disclaimers

Material Safety Data

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some but not all of the information required for the safe and proper use of this material. Before use, review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

Warranty and Limitation of Remedy

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specification.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver of Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitations of Remedy located on our website and in our catalog.

This Certificate shall not be reproduced except in full, without written approval from the Cayman Chemical Company.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com