

CERTIFICATE of ANALYSIS



Long-chain Fatty Acid Methyl Ester Mixture (fish oil)

Item No.: 28661
Batch No.: 0595917

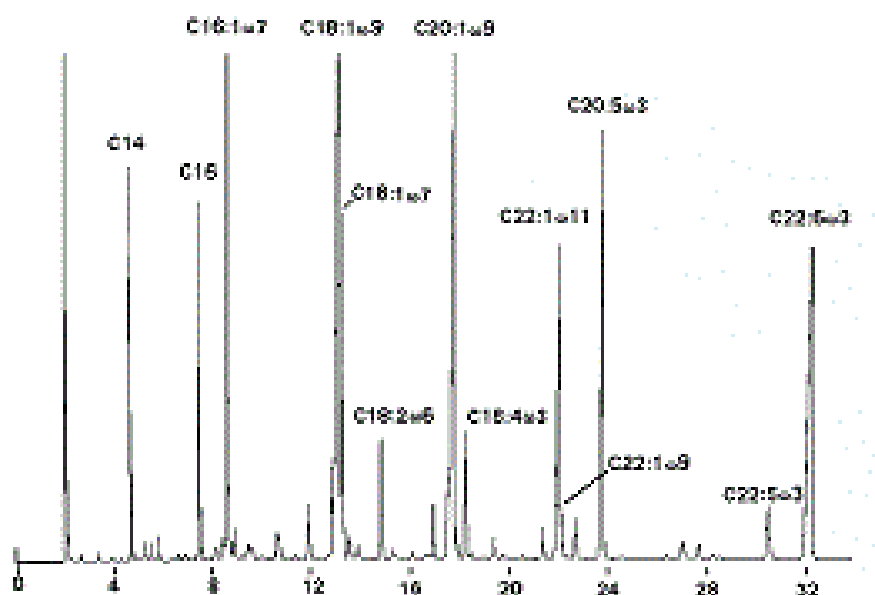
Expiry Date: 07NOV2026 (valid from date of certification)
Supplied as: A mixture of long-chain fatty acids from fish oil

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 two saturated and 11 mono- and poly-unsaturated fatty acids characterized by a terminal double bond located various distances from the omega end of the molecule. The fatty acids are isolated from fish oil and provided 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 by GC-MS.

Quality Information



Approval:

Certification Date: 23MAY2019

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	Item Name
1	6.556	Myristic Acid methyl ester
2	10.290	Palmitic Acid methyl ester
3	11.570	Palmitoleic Acid methyl ester
4	15.916	Oleic Acid methyl ester
5	16.076	Vaccenic Acid methyl ester
6	17.610	Linoleic Acid methyl ester
7	20.270	11(Z)-Eicosenoic Acid methyl ester
8	20.880	Stearidonic Acid methyl ester
9	24.600	11(E)-Docosenoic Acid methyl ester
10	24.746	13(Z)-Docosenoic Acid methyl ester
11	26.830	Eicosapentaenoic Acid methyl ester
12	34.926	Docosapentaenoic Acid methyl ester
13	36.890	Docosahexaenoic Acid methyl ester
Column:		SP-2330, 30 m x 0.25 mm x 0.2 μ
Oven:		150°C (8 min.), 3°C/min. to 190°C (29 min.)
Inj./Det.:		250°C/250°C
Carrier Gas:		helium
Injection:		1.0 μ l, 100:1 split

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	23MAY2019	Initial version
02	19APR2024	Expiry date extension
03	JAN2025	Updated expiry date and added retention times

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