

PRODUCT INFORMATION



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

Item No. 28663

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

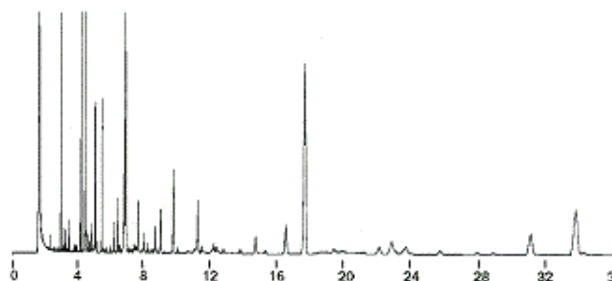
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Long-chain Fatty Acid Methyl Ester Mixture contains three saturated and 18 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 Menhaden fish oil and provided as methyl esters. This mixture is intended for use as a reference standard for the detection of these long-chain fatty acids from various sample types by GC-MS.

Included are:

Elution Order*	Item Name
1	Myristic Acid methyl ester
2	Palmitic Acid methyl ester
3	Palmitoleic Acid methyl ester
4	10,12-Hexadecadienoic Acid methyl ester
5	6,9,12-Hexadecatrienoic Acid methyl ester
6	4,8,12,16-Hexadecatetraenoic Acid methyl ester
7	Stearic Acid methyl ester
8	Oleic Acid methyl ester
9	Vaccenic Acid methyl ester
10	Linoleic Acid methyl ester
11	10,14-Octadecadienoic Acid
12	10,12,14-Octadecatrienoic Acid methyl ester
13	Linolenic Acid methyl ester
14	Stearidonic Acid methyl ester
15	11(Z)-Eicosenoic Acid methyl ester
16	Arachidonic Acid methyl ester
17	ω-3 Arachidonic Acid methyl ester
18	Eicosapentaenoic Acid methyl ester
19	Heneicosapentaenoic Acid methyl ester
20	Docosapentaenoic Acid methyl ester
21	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.

Copyright Cayman Chemical Company, 11/22/2019

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