

PRODUCT INFORMATION



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

Item No. 28661

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.

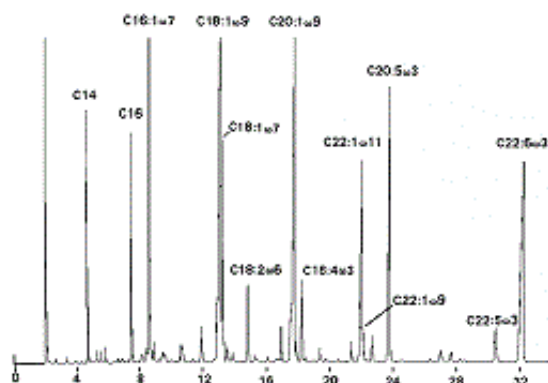
Description and Data

Long-chain Fatty Acid Methyl Ester Mixture contains 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.

Included are:

Elution Order*	Item Name
1	Myristic Acid methyl ester
2	Palmitic Acid methyl ester
3	Palmitoleic Acid methyl ester
4	Oleic Acid methyl ester
5	Vaccenic Acid methyl ester
6	Linoleic Acid methyl ester
7	11(Z)-Eicosenoic Acid methyl ester
8	Stearidonic Acid methyl ester
9	11(E)-Docosenoic Acid methyl ester
10	13(Z)-Docosenoic Acid methyl ester
11	Eicosapentaenoic Acid methyl ester
12	Docosapentaenoic Acid methyl ester
13	Docosahexaenoic Acid methyl ester

*GC Column Used: SP-2330, 30 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, 11/27/2019

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