

PRODUCT INFORMATION



Fatty Acid Methyl ester GC-MS Mixture

Item No. 20503

Supplied as: A solution in dichloromethane (1 mg/ml each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥7 years

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

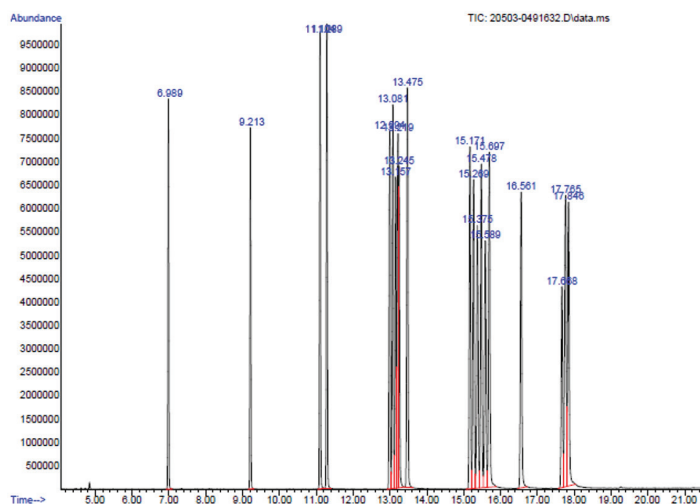
Description

This mixture contains various fatty acid methyl esters. The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in GC-MS applications. The solution may be serially diluted for preparation of calibrators and QC standards and/or used directly as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

The fatty acid methyl esters represented in this mixture include the ester versions of various important essential fatty acids, as well as endogenous fatty acids, including those that serve as precursors in the production of many other lipids. As a more lipophilic form of the free acid, these lipids are often more amenable for the formulation of fatty acid-containing diets and dietary supplements.

Contents

Each ampule contains at least 1 ml.



Item Number: 20503 Fatty Acid methyl ester GC-MS Mixture		
Item Name	Item No.	Retention Time (min)
Lauric Acid methyl ester	20608	6.99
Myristic Acid methyl ester	9001867	9.21
Palmitoleic Acid methyl ester	20605	11.10
Palmitic Acid methyl ester	10007358	11.29
γ -Linolenic Acid methyl ester	10006579	12.99
Stearidonic Acid methyl ester	10005000	13.08
Linoleic Acid methyl ester	20603	13.16
Oleic Acid methyl ester	20604	13.22
Linolenic Acid methyl ester	9000290	13.25
Stearic Acid methyl ester	20609	13.48
Arachidonic Acid methyl ester	90014	15.17
Eicosapentaenoic Acid methyl ester	9000295	15.27
Dihomo- γ -Linolenic Acid methyl ester	10006580	15.37
ω -3 Arachidonic Acid methyl ester	10006454	15.48
Eicosadienoic Acid methyl ester	20750	15.59
11(Z),14(Z),17(Z)-Eicosatrienoic Acid methyl ester	17622	15.70
Heptacosapentaenoic Acid methyl ester	11622	16.56
Docosahexaenoic Acid methyl ester	10006865	17.67
cis-4,10,13,16-Docosatetraenoic Acid methyl ester	10006866	17.77
Docosapentaenoic Acid methyl ester	9001870	17.85

Conditions:
Column: 30 m x 0.32 mm, 0.5 μ m Rtx-5MS Flow Rate: 2 ml/min Inlet Temp: 300°C
Split: 15:1 Carrier gas: He Transfer Line temp: 300°C
Oven program: Hold 120°C for 1 min, 10°C/min to 220°C, 5°C/min to 300°C, hold at 300°C for 3 min

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, 10/19/2023

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