

PRODUCT INFORMATION



EU Pharmacopoeia Fatty Acid Methyl Ester Mixture 1

Item No. 9004298

Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic

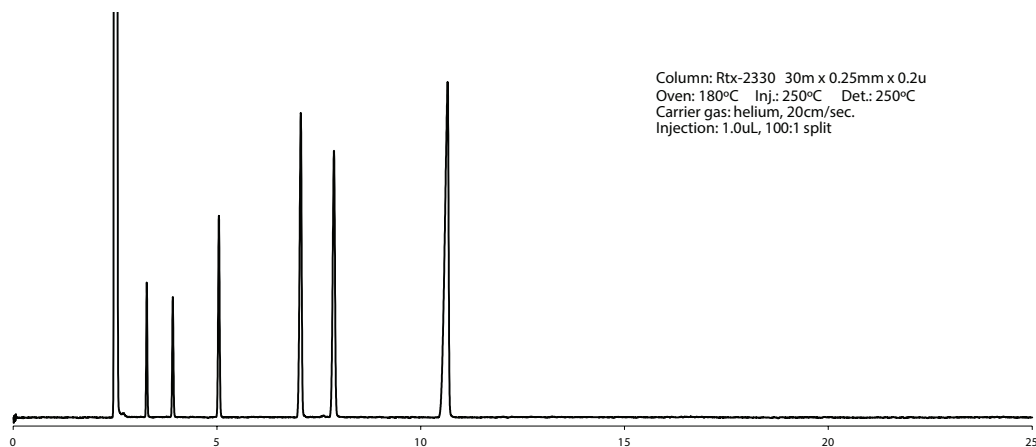
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

EU Pharmacopoeia Fatty Acid Methyl Ester Mixture 1 contains dodecanoic, tetradecanoic, hexadecanoic, octadecanoic, oleic, and eicosanoic acid methyl esters. According to the EU Pharmacopoeia 2.4.2 B, it is intended for use as a reference standard for the detection of these fatty acid methyl esters (FAMES) in various sample types by GC-MS.

This mixture contains fatty acid methyl esters. Included are:

Elution Order	Item Number	Carbon Number	Item Name	% Concentration by Weight
1	20608	C12:0	Lauric Acid methyl ester	5
2	9001867	C14:0	Myristic Acid methyl ester	5
3	10007358	C16:0	Palmitic Acid methyl ester	10
4	20609	C18:0	Stearic Acid methyl ester	20
5	20604	C18:1 [cis-9]	Oleic Acid methyl ester	20
6	26724	C20:0	Arachidic Acid methyl ester	40



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, 02/04/2026

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