

PRODUCT INFORMATION



EU Pharmacopoeia Fatty Acid Methyl Ester Mixture 2

Item No. 9004297

Synonym: SFE 20 Fatty Acid Methyl Ester Mixture
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years

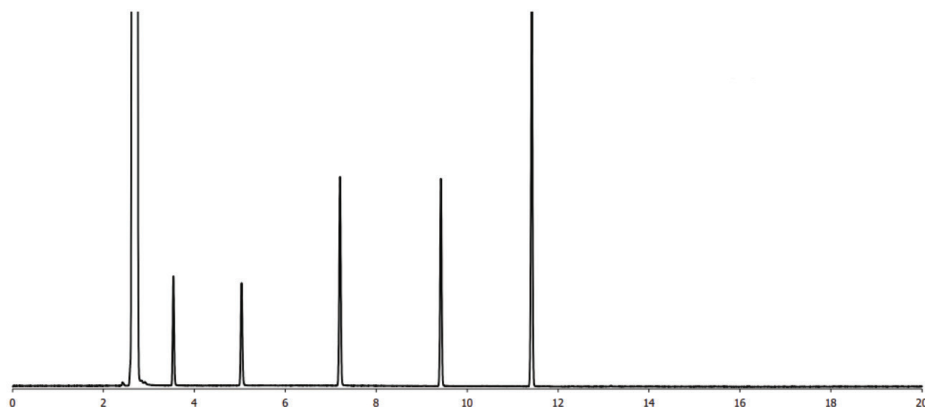
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

EU Pharmacopoeia Fatty Acid Methyl Ester Mixture 2 contains hexanoic, octanoic, decanoic, dodecanoic, and tetradecanoic acid methyl esters. According to the EU Pharmacopoeia 2.4.2 A, 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	33675	C6:0	Hexanoic Acid methyl ester	10
2	26717	C8:0	Octanoic Acid methyl ester	10
3	26864	C10:0	Decanoic Acid methyl ester	20
4	20608	C12:0	Lauric Acid methyl ester	20
5	9001867	C14:0	Myristic 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, 08/06/2026

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