

# PRODUCT INFORMATION



## C18 Fatty Acid Methyl Ester Mixture

Item No. 29369

**Synonym:** SFE 18 Fatty Acid Methyl Ester Mixture  
**Supplied as:** A solution in dichloromethane  
**Storage:** -20°C  
**Stability:** ≥2 years  
**Concentration:** 25 mg/ml

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

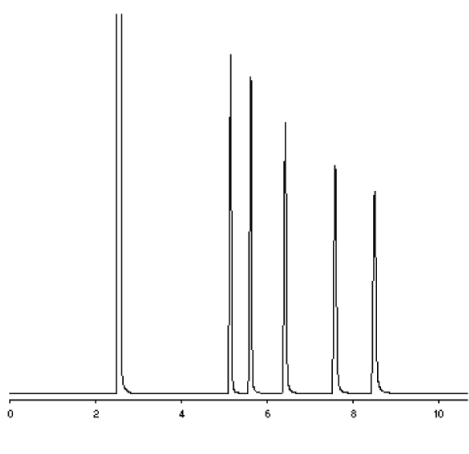
### Description and Data

C18 Fatty Acid Methyl Ester Mixture contains five C18 fatty acid methyl esters. It is intended for use as a reference standard for the detection of C18 fatty acids in various sample types by GC-MS.

Included are:

Elution Order*	Item Name	% Concentrations by Weight
1	Stearic Acid methyl ester	20
2	Oleic Acid methyl ester	20
3	Linoleic Acid methyl ester	20
4	α-Linolenic Acid methyl ester	20
5	Stearidonic Acid methyl ester	20

\*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, 03/13/2024

#### 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](http://WWW.CAYMANCHEM.COM)