

# PRODUCT INFORMATION



## Long-chain Monounsaturated Fatty Acid Methyl Ester Mixture

Item No. 29353

**Supplied as:** A solution in dichloromethane

**Storage:** -20°C

**Stability:** ≥1 year

**Concentration:** 50 mg/ml

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

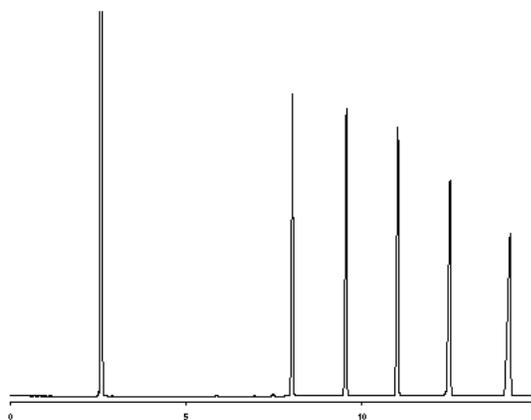
### Description and Data

Long-chain monounsaturated fatty acid methyl ester mixture contains five long-chain monounsaturated fatty acid methyl esters. It is intended for use as a reference standard for the detection of long-chain monounsaturated fatty acids in various sample types by GC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with nitrogen/argon to prevent lipid oxidation.

Included are:

Elution Order*	Item Name	% Concentration by Weight
1	Palmitoleic Acid methyl ester	20
2	Oleic Acid methyl ester	20
3	11(Z)-Eicosenoic Acid methyl ester	20
4	13(Z)-Docosenoic Acid methyl ester	20
5	Nervonic Acid methyl ester	20

\*GC Column Used: SP2330 or RTX2330, 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, 08/19/2022

#### 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