

PRODUCT INFORMATION



Very Long-chain Saturated Fatty Acid Methyl Ester Mixture

Item No. 29370

Supplied as: A solution in dichloromethane

Storage: -20°C

Stability: ≥1 year

Concentration: 25 mg/ml

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

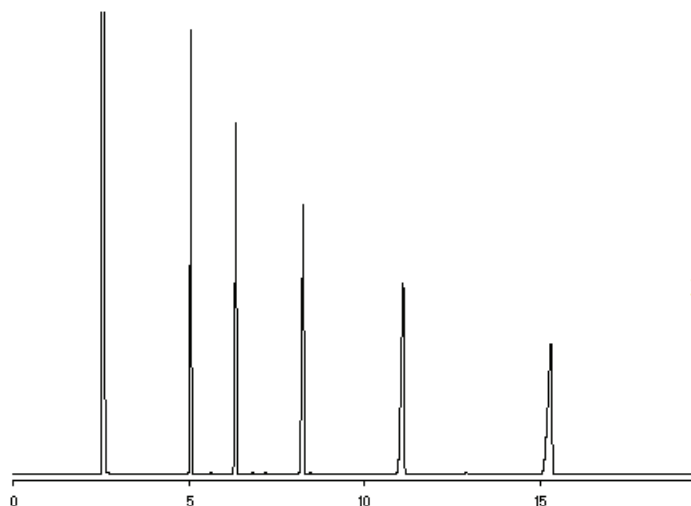
Description and Data

Very Long-chain Saturated Fatty Acid Methyl Ester Mixture contains five very long-chain saturated fatty acid methyl esters. It is intended for use as a reference standard for the quantification of very long-chain saturated fatty acids in various sample types by GC-MS.

Included are:

Elution Order*	Item Name	% Concentrations by Weight
1	Lignoceric Acid methyl ester	20
2	Hexacosanoic Acid methyl ester	20
3	Octacosanoic Acid methyl ester	20
4	Triacontanoic Acid methyl ester	20
5	Dotriacontanoic Acid methyl ester	20

*GC Column Used: RTX-1 or SPB-1, 30 m



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, 08/15/2022

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