

PRODUCT INFORMATION



Odd-chain Fatty Acid MaxSpec® LC-MS Mixture Item No. 19227

Supplied as: A solution in ethanol (10 µg/ml of each compound)
Fill volume: >1 ml
Storage: -20°C
Stability: ≥5 years

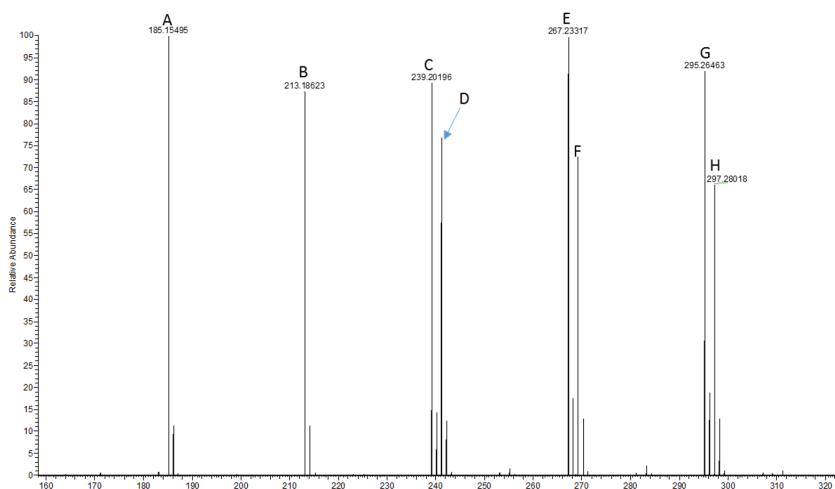
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

This mixture contains various odd-chain fatty acids (C11:0-C19:1). The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in LC-MS applications. The solution may be serially diluted for preparation of calibrators and QC standards and/or used directly as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

The odd-chain fatty acids represented in this mixture are either saturated or monounsaturated and composed of 11-19 carbons. These compounds are naturally occurring in animal fats and vegetable oils and can be used as biological markers for the intake of dietary fats in the assessment of metabolic risk factors.

Contents



Item Number: 19227 Odd-chain Fatty Acid MaxSpec® LC-MS Mixture				
Item Name	Formula	[MH] ⁻ Theoretical m/z	[MH] ⁻ Measured m/z	Mass Accuracy (ppm)
Undecanoic Acid	C ₁₁ H ₂₂ O ₂	185.15470	185.15495	1.350
Tridecanoic Acid	C ₁₃ H ₂₆ O ₂	213.18600	213.18623	1.079
14-Pentadecenoic Acid	C ₁₅ H ₃₀ O ₂	239.20165	239.20196	1.296
Pentadecanoic Acid	C ₁₅ H ₃₀ O ₂	241.21730	241.21753	0.953
cis-10-Heptadecenoic Acid	C ₁₇ H ₃₂ O ₂	267.23295	267.23317	0.823
Heptadecanoic Acid	C ₁₇ H ₃₄ O ₂	269.24860	269.24874	0.520
cis-10-Nonadecenoic Acid	C ₁₉ H ₃₆ O ₂	295.26425	295.26463	1.287
Nonadecanoic Acid	C ₁₉ H ₃₈ O ₂	297.27990	297.28018	0.942

Negative ESI-FTMS Analysis on a Q-Exactive Plus Quadrupole-Orbitrap Mass Spectrometer (Thermo Scientific)
Mass Range: 160-320 m/z Mass Resolution: 140,000 Injection Volume: 10 µl
Flow Injection Analysis in 1:1 MeOH:Water with 5 mM Ammonium Acetate at 60 µl/min

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, 07/14/2023

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