

PRODUCT INFORMATION

Saturated/Monounsaturated Fatty Acid MaxSpec® LC-MS Mixture Item No. 17942

Supplied as: A solution in ethanol (10 µg/ml of each compound)
Fill Volume: >1 ml
Purity: ≥95%
Storage: -20°C
Stability: ≥3 years

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

Description

This mixture contains various saturated or unsaturated fatty acids. 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 saturated or unsaturated fatty acids represented in this mixture are composed of 12-24 carbons. These compounds are found naturally in plants and mammals and have important biological roles, including serving as precursors in the production of various other lipids. Some have also been used for industrial production of soaps, cosmetics, lubricants, releasing agents, and surfactants.

Contents

Item Number	Item Name
9000339	Arachidic Acid
10006626	Lauric Acid
13353	Lignoceric Acid
13351	Myristic Acid
13940	Nervonic Acid
90260	Oleic Acid
10006627	Palmitic Acid
10009871	Palmitoleic Acid
10011298	Stearic Acid

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, 07/17/2023

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