

PRODUCT INFORMATION



Short-chain Fatty Acid MaxSpec® LC-MS Mixture

Item No. 18824

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

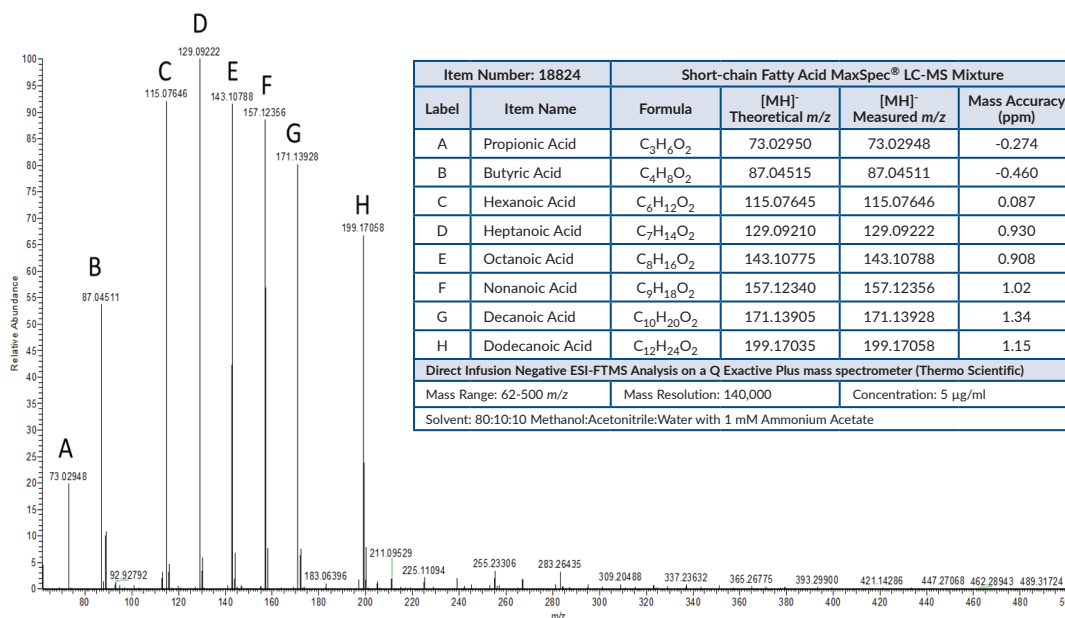
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

This mixture contains various short-chain fatty acids (≤12:0). 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 short-chain fatty acids represented in this mixture are saturated and are composed of 12 or less carbons. These compounds are naturally occurring and have been used as intermediates in chemical syntheses and in the manufacture of esters for industrial production of perfumes, lubricants, greases, rubber, dyes, plastics, food additives, herbicides, and pharmaceuticals.

Contents



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