

PRODUCT INFORMATION



Short-chain Fatty Acid Mixture 3

Item No. 28682

Supplied as: A solution in water
Storage: 4°C
Stability: ≥2 years
Concentration: 10 mM for each lipid

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

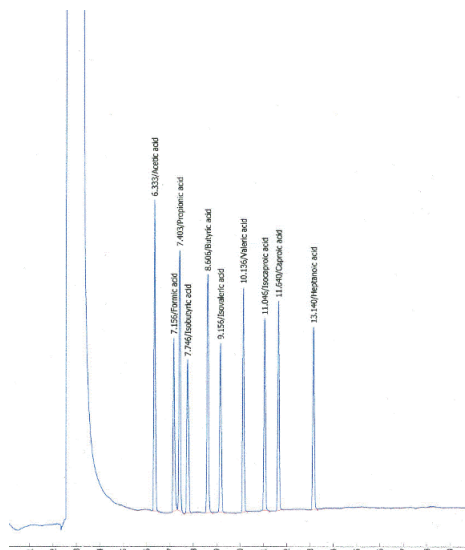
Description and Data

Short-chain Fatty Acid Mixture 3 contains ten short-chain volatile fatty acids. It is intended for use as a reference standard for the detection of these short-chain fatty acids in various sample types by GC-FID, GC-MS, or LC-MS. For some applications, derivatization may be required.

Included are:

Components*	Item Name
1	Acetic Acid
2	Formic Acid
3	Propionic Acid
4	Isobutyric Acid
5	N-Butyric Acid
6	Isovaleric Acid
7	N-Valeric Acid
8	4-Methylpentanoic Acid
9	N-Hexanoic Acid
10	Heptanoic Acid

*GC Column Used: Nukol 30 m x 0.53 mm



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.

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