

# PRODUCT INFORMATION



## Short-chain Fatty Acid Mixture 2

Item No. 28680

**Supplied as:** A solution in water  
**Storage:** -20°C  
**Stability:** ≥2 years  
**Concentration:** 1 mg/ml for each lipid

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

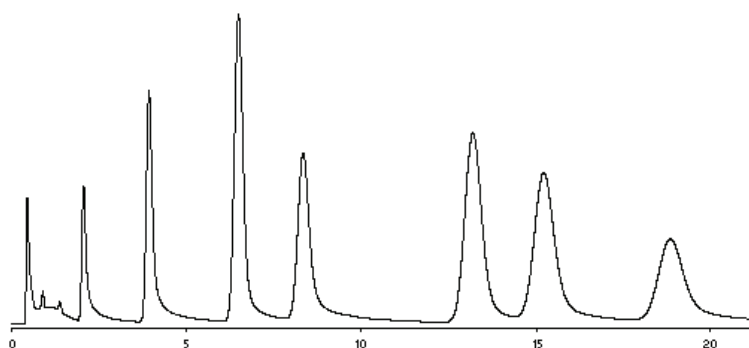
### Description and Data

Short-chain Fatty Acid Mixture 2 contains the short-chain and water-soluble fatty acids acetic, propionic, isobutyric, n-butyric, 2-methylbutyric, isovaleric, and n-valeric acids. It is intended for use as a reference standard for the detection of these fatty acids in various sample types by GC- or LC-MS.

Included are:

Elution Order*	Item Name
1	Acetic Acid
2	Propionic Acid
3	Isobutyric Acid
4	n-Butyric Acid
5	2-Methylbutyric acid
6	Isovaleric Acid
7	n-Valeric Acid

\*GC Column Used: 80/120 Carbowax B-DA/4% Carbowax 20 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, 11/13/2019

#### 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](http://WWW.CAYMANCHEM.COM)