

# PRODUCT INFORMATION



## Bacterial Fatty Acid Methyl Ester Mixture 2

Item No. 29373

**Supplied as:** A solution in methyl caproate

**Storage:** -20°C

**Stability:** ≥2 years

**Concentration:** 10 mg/ml

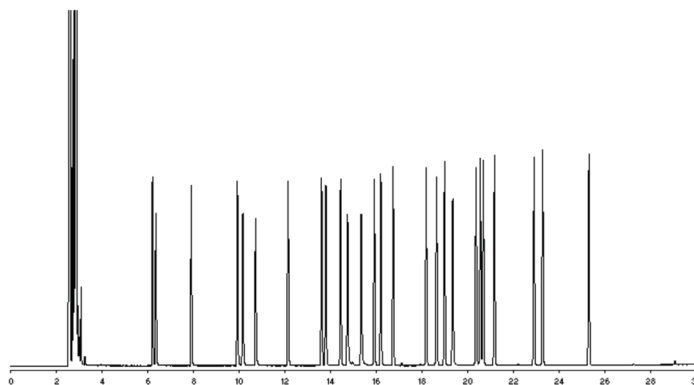
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description and Data

Bacterial Fatty Acid Methyl Ester Mixture 2 contains 26 bacterial fatty acid methyl esters. It is intended for use as a reference standard for the detection of bacterial fatty acids in various sample types.

Included are:

Component	Item Name
1	Undecanoic Acid methyl ester
2	2-hydroxy Decanoic Acid methyl ester
3	Lauric Acid methyl ester
4	Tridecanoic Acid methyl ester
5	2-hydroxy Lauric Acid methyl ester
6	3-hydroxy Lauric Acid methyl ester
7	Myristic Acid methyl ester
8	13-methyl Myristic Acid methyl ester
9	12-methyl Myristic Acid methyl ester
10	Pentadecanoic Acid methyl ester
11	2-hydroxy Myristic Acid methyl ester
12	3-hydroxy Myristic Acid methyl ester
13	14-methyl Pentadecanoic Acid methyl ester
14	Palmitoleic Acid methyl ester
15	Palmitic Acid methyl ester
16	15-methyl Palmitic Acid methyl ester
17	cis-9,10-Methylenehexadecanoic Acid methyl ester
18	Heptadecanoic Acid methyl ester
19	2-hydroxy Palmitic Acid methyl ester
20	Linoleic Acid methyl ester
21	Oleic Acid methyl ester
22	Elaidic Acid methyl ester
23	Stearic Acid methyl ester
24	cis-9,10-Methyleneoctadecanoic Acid methyl ester
25	Nonadecanoic Acid methyl ester
26	Arachidic Acid methyl ester



#### 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, 10/19/2022

#### 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