

# PRODUCT INFORMATION



## 2-Hydroxy Fatty Acid Methyl Ester Mixture

Item No. 24658

**Purity:** ≥98%  
**Supplied as:** A solution in chloroform  
**Storage:** -20°C  
**Stability:** ≥2 years

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

### Description and Data

2-Hydroxy fatty acid methyl ester mixture is a mixture of fatty acid methyl esters. This mixture is supplied in an amber ampule in which the headspace has been purged with nitrogen/argon to prevent lipid oxidation. This product has been designed for direct use in GC-MS applications.

This mixture contains fatty acid methyl esters. Included are:

Elution Order	Item Number	Carbon Number	Item Name	% Concentration by Weight
1	24588	C14:0	2-hydroxy Myristic Acid methyl ester	20
2	24589	C16:0	2-hydroxy Palmitic Acid methyl ester	20
3	24591	C18:0	2-hydroxy Stearic Acid methyl ester	15
4	24593	C20:0	2-hydroxy Arachidic Acid methyl ester	15
5	24596	C22:0	2-hydroxy Docosanoic Acid methyl ester	10
6	24598	C23:0	2-hydroxy Tricosanoic Acid methyl ester	10
7	24599	C24:0	2-hydroxy Lignoceric Acid methyl ester	10

#### 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, 10/20/2022

#### 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