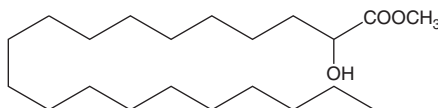


PRODUCT INFORMATION

2-hydroxy Arachidic Acid methyl ester

Item No. 24593

CAS Registry No.: 16742-49-7
Formal Name: 2-hydroxy-eicosanoic acid, methyl ester
Synonyms: α-hydroxy methyl Eicosanoate, methyl 2-hydroxy Eicosanoate
MF: C₂₁H₄₂O₃
FW: 342.6
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

2-hydroxy Arachidic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy arachidic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Arachidic acid methyl ester is soluble in organic solvents such as chloroform and ethyl ether.

Description

2-hydroxy Arachidic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in sediment samples from the Harney River and in petroleum ether extracts of *I. eriocarpa*.^{1,2} It is also a volatile flavoring compound in anhydrous milk fat (AMF) that is reduced by treatment with the lipases lipzyme-435 and TL-IM in a time-dependent manner.³

References

1. Jaffé, R., Rushdi, A.I., Medeiros, P.M., *et al.* Natural product biomarkers as indicators of sources and transport of sedimentary organic matter in a subtropical river. *Chemosphere* **64**(11), 1870-1884 (2006).
2. Das, M. and Malipeddi, H. Phytochemical screening, GC-MS analysis and biological activity of *Ipomoea eriocarpa* leaf extracts. *Int. J. Pharm. Pharm. Sci.* **6**(4), 592-594 (2014).
3. Omar, K.A., Gounga, M.E., Liu, R., *et al.* Effects of microbial lipases on hydrolyzed milk fat at different time intervals in flavour development and oxidative stability. *J. Food Sci. Technol.* **53**(2), 1035-1046 (2016).

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, 11/11/2022

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