

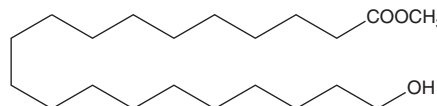
# PRODUCT INFORMATION



## 20-hydroxy Arachidic Acid methyl ester

Item No. 24653

CAS Registry No.: 37477-29-5  
Formal Name: 20-hydroxy-eicosanoic acid, methyl ester  
Synonym: Methyl 20-hydroxyeicosanoate  
MF:  $C_{21}H_{42}O_3$   
FW: 342.6  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

20-hydroxy Arachidic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 20-hydroxy arachidic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 20-hydroxy Arachidic acid methyl ester is soluble in ethanol (warm), chloroform, and ethyl ether.

### Description

20-hydroxy Arachidic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in *P. reticulata* cork extracts as well as peat samples from inland Florida.<sup>1,2</sup>

### References

1. Mota, G.S., Sartori, C.J., Ferreira, J., et al. Cellular structure and chemical composition of cork from *Plathymenia reticulata* occurring in the Brazilian Cerrado. *Ind. Crops Prod.* **90**, 65-75 (2016).
2. 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).

#### 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, 12/17/2022

#### 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