

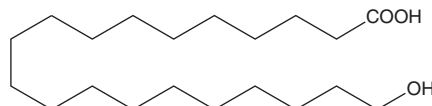
PRODUCT INFORMATION



20-hydroxy Arachidic Acid

Item No. 24652

CAS Registry No.: 62643-46-3
Formal Name: 20-hydroxy-eicosanoic acid
Synonym: ω -hydroxy Eicosanoic Acid
MF: $C_{20}H_{40}O_3$
FW: 328.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20
Stability: ≥ 4 years



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

Laboratory Procedures

20-hydroxy Arachidic acid is supplied as a solid. A stock solution may be made by dissolving the 20-hydroxy arachidic acid in the solvent of choice, which should be purged with an inert gas. 20-hydroxy Arachidic acid is soluble in chloroform.

Description

20-hydroxy Arachidic acid is a hydroxylated fatty acid that has been found in the suberin component of silver birch (*B. pendula*) outer bark as well as the leaves, roots, and wood of *Q. ilex* from Cala Violina and Colognole, Italy.^{1,2} 20-hydroxy Arachidic acid was identified as a potential inhibitor of human rhinovirus coat protein in a structure-based virtual screen.³

References

1. Karnaouri, A., Lange, H., Crestini, C., *et al.* Chemoenzymatic fractionation and characterization of pretreated birch outer bark. *ACS Sustainable Chem. Eng.* **4**(10), 5289-5302 (2016).
2. Andreetta, A., Dignac, M.-F., and Carnicelli, S. Biological and physico-chemical processes influence cutin and suberin biomarker distribution in two Mediterranean forest soil profiles. *Biogeochemistry* **112**(1-3), 41-58 (2013).
3. Rollinger, J.M., Steindl, T.M., Schuster, D., *et al.* Structure-based virtual screening for the discovery of natural inhibitors for human rhinovirus coat protein. *J. Med. Chem.* **51**(4), 842-851 (2008).

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, 12/06/2022

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