

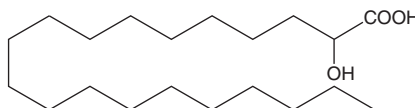
PRODUCT INFORMATION



2-hydroxy Arachidic Acid

Item No. 24592

CAS Registry No.: 16742-48-6
Formal Name: 2-hydroxy-eicosanoic acid
Synonym: α -hydroxy Arachic Acid, α -hydroxy Eicosanoic Acid, 2-Hydroxy C20:0
MF: C₂₀H₄₀O₃
FW: 328.5
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

2-hydroxy Arachidic acid is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy arachidic acid in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Arachidic acid is soluble in a 5:1 solution of chloroform:methanol.

Description

2-hydroxy Arachidic acid is a hydroxylated fatty acid that has been found in pasture grass (*S. sphacelata*) and soil from abandoned pastures in Ecuador, wool wax, and as the N-acyl chain of galactosylceramides in mouse brain.¹⁻³

References

1. Hamer, U., Rumpel, C., and Dignac, M.-F. Cutin and suberin biomarkers as tracers for the turnover of shoot and root derived organic matter along a chronosequence of Ecuadorian pasture soils. *Eur. J. Soil Sci.* **63(6)**, 808-819 (2012).
2. Jenske, R. and Vetter, W. Concentrations of medium-chain 2- and 3-hydroxy fatty acids in foodstuffs. *Food Chem.* **114(3)**, 1122-1129 (2009).
3. Alderson, N.L., Maldonado, E.N., Kern, M.J., et al. FA2H-dependent fatty acid 2-hydroxylation in postnatal mouse brain. *J. Lipid Res.* **47(12)**, 2772-2780 (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, 11/15/2022

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