

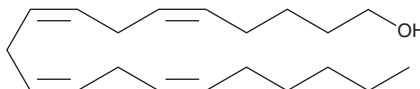
PRODUCT INFORMATION



Arachidonyl Alcohol

Item No. 32539

CAS Registry No.: 13487-46-2
Formal Name: 5Z,8Z,11Z,14Z-5,8,11,14-eicosatetraen-1-ol
Synonyms: (all-Z)-5,8,11,14-Eicosatetraen-1-ol, FOH 20:4
MF: $C_{20}H_{34}O$
FW: 290.5
Purity: $\geq 98\%$
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Arachidonyl alcohol is supplied as a neat oil. A stock solution may be made by dissolving the arachidonyl alcohol in the solvent of choice, which should be purged with an inert gas. Arachidonyl alcohol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of arachidonyl alcohol in these solvents is approximately 30 mg/ml.

Description

Arachidonyl alcohol is a polyunsaturated fatty alcohol produced by the reduction of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607).^{1,2} It has been used as a substrate in the synthesis of various ether lipids.

References

1. Natarajan, V. and Schmid, H.H. Substrate specificities in ether lipid biosynthesis: Metabolism of polyunsaturated fatty acids and alcohols by rat brain microsomes. *Biochem. Biophys. Res. Commun.* **79**(2), 411-416 (1977).
2. Nagao, T., Watanabe, Y., Tanaka, S., *et al.* Microbial conversion of arachidonic acid to arachidonyl alcohol by a new *Acinetobacter* species. *J. Am. Oil Chem. Soc.* **89**, 1663-1671 (2012).

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, 04/29/2025

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