

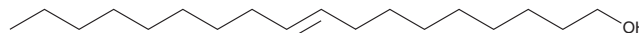
PRODUCT INFORMATION



Elaidyl Alcohol

Item No. 30459

CAS Registry No.: 506-42-3
Formal Name: 9E-octadecen-1-ol
Synonyms: NSC 72786, *trans*-9-Octadecenoyl Alcohol
MF: C₁₈H₃₆O
FW: 268.5
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

Elaidyl alcohol is supplied as a solid. A stock solution may be made by dissolving the elaidyl alcohol in the solvent of choice, which should be purged with an inert gas. Elaidyl alcohol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of elaidyl alcohol in these solvents is approximately 15, 10, and 5 mg/ml, respectively.

Elaidyl alcohol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, elaidyl alcohol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Elaidyl alcohol has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Elaidyl alcohol is a monounsaturated fatty alcohol produced by the hydrogenation of elaidic acid (Item No. 90250).¹ It is active against herpes simplex virus 2 (HSV-2) and bacteriophage $\phi 6$, but not *Pseudoalteromonas* phage PM2, virions in plaque assays (IC₅₀s = 1.9, 0.4, and >75 μ M, respectively).² Elaidyl alcohol has been used in the formation of bimolecular films to study the effect of unsaturation on model lipid membranes.³

References

- Cheah, K.Y., Tang, T.S., Mizukami, F., *et al.* Selective hydrogenation of oleic acid to 9-octadecen-1-ol: Catalyst preparation and optimum reaction conditions. *J. Am. Oil Chem. Soc.* **69(5)**, 410-416 (1992).
- Sands, J., Auperin, D., and Snipes, W. Extreme sensitivity of enveloped viruses, including herpes simplex, to long-chain unsaturated monoglycerides and alcohols. *Antimicrob. Agents Chemother.* **15(1)**, 67-73 (1979).
- Legaly, G., Weiss, A., and Stuke, E. Effect of double-bonds on bimolecular films in membrane models. *Biochim. Biophys. Acta* **470(3)**, 331-341 (1977).

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, 10/25/2022

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