

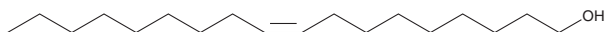
PRODUCT INFORMATION



Oleyl Alcohol

Item No. 30385

CAS Registry No.: 143-28-2
Formal Name: 9Z-octadecen-1-ol
Synonym: FOH 18:1
MF: $C_{18}H_{36}O$
FW: 268.5
Purity: $\geq 80\%$
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

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

Description

Oleyl alcohol is a monounsaturated fatty alcohol produced by the reduction of oleic acid (Item Nos. 90260 | 24659).¹ Topical application of an oleyl alcohol (5% w/w), propylene glycol, and Carbopol® 940 gel containing the non-steroidal anti-inflammatory drug (NSAID) piroxicam (Item No. 13368) enhances piroxicam absorption into isolated rat abdominal skin compared to a control gel not containing oleyl alcohol.² Oleyl alcohol also increases naloxone penetration into isolated postmortem human skin when applied at a 10% concentration.³

References

1. Klonowski, R.S., Findley, T.W., Josefson, C.M., *et al.* Oleyl alcohol from animal fats by catalytic reduction. *J. Am. Oil Chem. Soc.* **47**, 326-328 (1970).
2. Santoyo, S., Arellano, A., Ygartua, P., *et al.* Penetration enhancer effects on the in vitro percutaneous absorption of piroxicam through rat skin. *Int. J. Pharm.* **117(2)**, 219-224 (1995).
3. Aungst, B.J., Rogers, N.J., and Shefter, E. Enhancement of naloxone penetration through human skin in vitro using fatty acids, fatty alcohols, surfactants, sulfoxides and amides. *Int. J. Pharm.* **33(1-3)**, 225-234 (1986).

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