

PRODUCT INFORMATION

Dihydrophytol

Item No. 31692

CAS Registry No.: 645-72-7

Formal Name: 3,7,11,15-tetramethyl-1-hexadecanol

Synonyms: Octahydrogeranylgeraniol, Phytanic Alcohol

MF: $C_{20}H_{42}O$

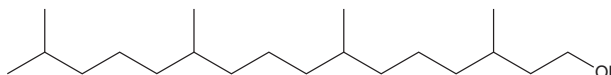
FW: 298.6

Purity: $\geq 95\%$

Supplied as: A neat oil

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Dihydrophytol is supplied as a neat oil. A stock solution may be made by dissolving the dihydrophytol in the solvent of choice, which should be purged with an inert gas. Dihydrophytol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of dihydrophytol in ethanol and DMF is approximately 10 mg/ml and approximately 5 mg/ml in DMSO.

Description

Dihydrophytol is an intermediate in the biosynthesis of phytanic acid (Item No. 90360).¹ It is produced from phytol (Item No. 17401) in the rumen of cattle.²

References

1. van den Brink, D.M. and Wanders, R.J.A. Phytanic acid: Production from phytol, its breakdown and role in human disease. *Cell Mol. Life Sci.* **63**(15), 1752-1765 (2006).
2. Patton, S. and Benson, A.A. Phytol metabolism in the bovine. *Biochim. Biophys. Acta* **125**(1), 22-32 (1966).

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, 05/07/2024

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