

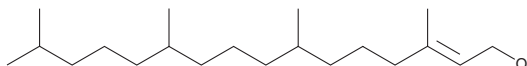
PRODUCT INFORMATION



Phytol

Item No. 17401

CAS Registry No.: 150-86-7
Formal Name: 3,7,11,15-tetramethyl-2-hexadecen-1-ol
MF: $C_{20}H_{40}O$
FW: 296.5
Purity: $\geq 85\%$ (mixture of isomers)
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

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

Description

Phytol is a diterpene alcohol obtained from the degradation of chlorophyll and has been used in the synthesis of Vitamins E and K. During the digestion process of ruminants, phytol is liberated from chlorophyll and converted to phytanic acid (Item No. 90360) to be stored in fats. While humans cannot derive phytol from chlorophyll, free phytol, obtained from the consumption of ruminant adipose tissue and dairy products, is readily absorbed in the small intestine and converted to phytanic acid.¹ Phytanic acid accumulates to toxic levels in a number of metabolic disorders, and the conversion of phytol to phytanic acid has been shown to be regulated via the activation of peroxisome proliferator-activated receptor α (PPAR α).^{1,2} Phytol and its metabolites have also been reported to activate retinoid X receptors (RXRs; K_i s range from 2.3-67.2 μM) and to promote the activity of PPAR/RXR heterodimers.^{2,3} Phytol also demonstrates sedative and anxiolytic effects through interaction with the GABA $_A$ receptor, and it has been explored as an antischistosomal agent in a mouse model of schistosomiasis.^{4,5}

References

1. Gloerich, J., van den Brink, D.M., Ruiter, J.P.N., et al. *J. Lipid Res.* **48**(1), 77-85 (2007).
2. Elmazar, M.M., El-Abhar, H.S., Schaalen, M.F., et al. *PLoS One* **8**(1), 1-10 (2013).
3. Kitareewan, S., Burka, L.T., Tomer, K.B., et al. *Mol. Biol. Cell* **7**(8), 1153-1166 (1996).
4. Costa, J.P., de Oliveira, G.A., de Almeida, A.A., et al. *Brain Res.* **1547**, 34-42 (2014).
5. de Moraes, J., de Oliveira, R.N., Costa, J.P., et al. *PLoS One* **8**(1), 1-12 (2014).

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, 11/07/2022

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