

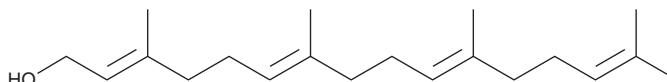
PRODUCT INFORMATION



Geranylgeranyl Alcohol

Item No. 13272

CAS Registry No.: 24034-73-9
Formal Name: 3,7,11,15-tetramethyl-2E,6E,10E,14-hexadecatetraen-1-ol
Synonyms: Geranylgeraniol, GG Alcohol
MF: C₂₀H₃₄O
FW: 290.5
Purity: ≥85%
Supplied as: A neat oil
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

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

Geranylgeranyl alcohol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, geranylgeranyl alcohol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Geranylgeranyl alcohol has a solubility of approximately 0.33 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

Geranylgeranyl alcohol is a diterpenoid produced through the isoprenoid pathway. It can be phosphorylated to give geranylgeranyl pyrophosphate (GGPP; Item No. 63330), a precursor for geranylgeranylation, a form of protein prenylation.¹ GGPP is also an intermediate in biosynthetic pathways for isoprenoids, carotenoids, plant hormones, and antibiotics.

Reference

1. McTaggart, S.J. Isoprenylated proteins. *Cell Mol. Life Sci.* **63(3)**, 255-267 (2006).

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, 06/05/2019

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