

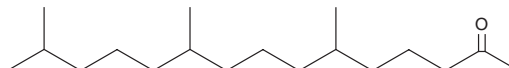
PRODUCT INFORMATION



Perhydrofarnesyl Acetone

Item No. 33825

CAS Registry No.: 502-69-2
Formal Name: 6,10,14-trimethyl-2-pentadecanone
Synonym: Hexahydrofarnesyl Acetone
MF: $C_{18}H_{36}O$
FW: 268.5
Purity: $\geq 95\%$
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Perhydrofarnesyl acetone is supplied as a liquid. A stock solution may be made by dissolving the perhydrofarnesyl acetone in the solvent of choice, which should be purged with an inert gas. Perhydrofarnesyl acetone is soluble in the organic solvent chloroform.

Description

Perhydrofarnesyl acetone is a sesquiterpenoid that has been found in orchid bees (*Euglossa*).¹

Reference

1. Zimmermann, Y., Ramírez, S.R., and Eltz, T. Chemical niche differentiation among sympatric species of orchid bees. *Ecology* **90**(11), 2994-3008 (2009).

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, 02/16/2026

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