

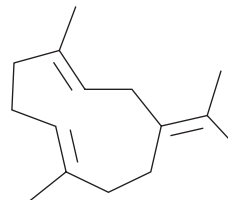
# PRODUCT INFORMATION



## Germacrene B

Item No. 34907

CAS Registry No.: 15423-57-1  
Formal Name: 1,5-dimethyl-8-(1-methylethylidene)-1E,5E-cyclodecadiene  
MF: C<sub>15</sub>H<sub>24</sub>  
FW: 204.4  
Purity: ≥90%  
Supplied as: A solution in methanol  
Storage: -20°C  
Stability: ≥2 years  
Item Origin: Synthetic



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

### Laboratory Procedures

Germacrene B is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Germacrene B is slightly soluble in chloroform.

### Description

Germacrene B is a sesquiterpene that has been found in *P. cablin* (patchouli).<sup>1</sup>

### Reference

1. Hasegawa, Y., Tajima, K., Toi, N., *et al.* An additional constituent occurring in the oil from a patchouli cultivar. *Flavour Fragr. J.* **7(6)**, 333-335 (1992).

#### 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, 11/08/2022

#### 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