

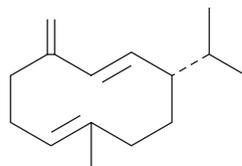
PRODUCT INFORMATION



Germacrene D

Item No. 26539

CAS Registry No.: 23986-74-5
Formal Name: (1E)-1-methyl-5-methylene-8S-(1-methylethyl)-1,6E-cyclodecadiene
Synonym: (-)-Germacrene D
MF: C₁₅H₂₄
FW: 204.4
Purity: ≥90%
Supplied as: A solution in heptane
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Cananga odorata*



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

Laboratory Procedures

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

Description

Germacrene D is a major volatile component of *Bursera* species and a precursor in sesquiterpene biosynthesis in a variety of plants, including *Solidago* species.^{1,2}

References

1. Noge, K. and Becerra, J.X. Germacrene D, a common sesquiterpene in the genus *Bursera* (Burseraceae). *Molecules* **14**(12), 5289-5297 (2009).
2. Bülow, N. and König, W.A. The role of germacrene D as a precursor in sesquiterpene biosynthesis: Investigations of acid catalyzed, photochemically and thermally induced rearrangements. *Phytochemistry* **55**(2), 141-168 (2000).

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, 01/18/2023

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