

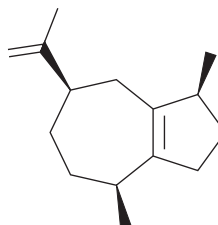
# PRODUCT INFORMATION



## $\alpha$ -Guaiene

Item No. 34686

CAS Registry No.: 3691-12-1  
Formal Name: (1S,4S,7R)-1,2,3,4,5,6,7,8-octahydro-1,4-dimethyl-7-(1-methylethenyl)-azulene  
MF:  $C_{15}H_{24}$   
FW: 204.4  
Purity:  $\geq 85\%$   
Supplied as: A solution in chloroform  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 2$  years  
Item Origin: Synthetic



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

### Laboratory Procedures

$\alpha$ -Guaiene is supplied as a solution in chloroform. To change the solvent, simply evaporate the  $\alpha$ -guaiene under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol and chloroform purged with an inert gas can be used.  $\alpha$ -Guaiene is slightly soluble in chloroform and methanol.

### Description

$\alpha$ -Guaiene is a sesquiterpene that has been found in *P. cablin*.<sup>1</sup>

### Reference

1. Kocovski, D., Du, M., Kan, J., et al. Antifungal effect of *Allium tuberosum*, *Cinnamomum cassia*, and *Pogostemon cablin* essential oils and their components against population of *Aspergillus* species. *J. Food Sci.* **78**(5), M731-M737 (2013).

#### 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, 08/03/2021

#### 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