

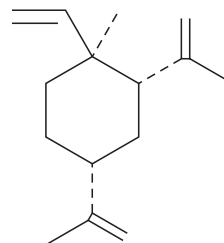
# PRODUCT INFORMATION



## β-Elemene

Item No. 19641

**CAS Registry No.:** 515-13-9  
**Formal Name:** (1S,2S,4R)-1-ethenyl-1-methyl-2,4-bis(1-methylethenyl)-cyclohexane  
**Synonyms:** (-)-β-Elemene, Levo-β-elemene  
**MF:** C<sub>15</sub>H<sub>24</sub>  
**FW:** 204.4  
**Purity:** ≥95%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

β-Elemene is supplied as a solution in ethanol. To change the solvent, simply evaporate the β-elemene under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of β-elemene in these solvents is approximately 30 mg/ml.

β-Elemene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, β-elemene, provided as a solution in ethanol, may be diluted with the aqueous buffer of choice. β-Elemene has a solubility of approximately 0.3 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

β-Elemene is a monoterpene that has been found in *C. sativa*.<sup>1</sup> It is intended for use as an analytical standard for the detection of β-elemene by GC- or LC-MS.

### Reference

1. Ross, S.A., and ElSohly, M.A. The volatile oil composition of fresh and air-dried buds of *Cannabis sativa*. *J. Nat. Prod.* **59**(1), 49-51 (1996).

#### 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, 07/02/2018

#### 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