

# PRODUCT INFORMATION



## Tricyclene

Item No. 34290

CAS Registry No.: 508-32-7  
Formal Name: 1,7,7-trimethyl-tricyclo[2.2.1.0<sup>2,6</sup>]heptane  
Synonym: NSC 86978  
MF: C<sub>10</sub>H<sub>16</sub>  
FW: 136.2  
Purity: ≥95%  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years  
Special Conditions: Low melting; temperature sensitive  
Item Origin: Synthetic



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

### Laboratory Procedures

Tricyclene is supplied as a solid. A stock solution may be made by dissolving the tricyclene in the solvent of choice, which should be purged with an inert gas. Tricyclene is slightly soluble in chloroform and methanol.

### Description

Tricyclene is a monoterpene that has been found in a variety of conifers, including *Abies*, *Pinus*, *Picea*, and *Tsuga*.<sup>1</sup>

### Reference

1. Yatagai, M. and Sato, T. Terpenes of leaf oils from conifers. *Biochem. Syst. Ecol.* **14**(5), 467-478 (1986).

#### 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, 11/02/2022

#### 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