

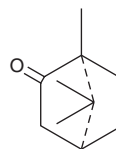
# PRODUCT INFORMATION



## (1S)-(-)-Camphor

Item No. 23176

**CAS Registry No.:** 464-48-2  
**Formal Name:** (4S)-1S,7,7-trimethyl-bicyclo[2.2.1]heptan-2-one  
**Synonyms:** (-)-Camphor, (S)-Camphor, NSC 26351  
**MF:** C<sub>10</sub>H<sub>16</sub>O  
**FW:** 152.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 291 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

(1S)-(-)-Camphor is supplied as a crystalline solid. A stock solution may be made by dissolving the (1S)-(-)-camphor in the solvent of choice. (1S)-(-)-Camphor is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of (1S)-(-)-camphor in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

(1S)-(-)-Camphor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (1S)-(-)-camphor should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (1S)-(-)-Camphor has a solubility of approximately 0.33 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

(1S)-(-)-Camphor is a monoterpene found in various plant extracts, including *P. suffruticosa* with acaricidal activity.<sup>1</sup> It exhibits fumigant toxicity against *T. putrescentiae* when used at a concentration of 10.45 mg.

### Reference

1. Tak, J.H., Kim, H.K., Lee, S.H., *et al.* Acaricidal activities of paeonol and benzoic acid from *Paeonia suffruticosa* root bark and monoterpenoids against *Tyrophagus putrescentiae* (Acari: Acaridae). *Pest Manag. Sci.* **62**(6), 551-557 (2006).

#### 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, 10/05/2022

#### 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