

PRODUCT INFORMATION

Cannabis Terpenes Mixture 2 (CRM)

Item No. 40527

Supplied as: A solution in isopropanol:toluene (1:1) (2,500 µg/ml of each compound)

Fill volume: >1 ml

Storage: -20°C

Stability: ≥2 years

Special conditions: Vortex or sonicate before use

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

Description

Cannabis Terpenes Mixture 2 (CRM) (Item No. 40527) is a certified reference material mixture containing terpenes found in Cannabis. The mixture contains camphene, ocimene, sabinene hydrate, and geranyl acetate among others. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and as a reference standard for these substances during routine analytical checks using LC- or GC-MS instrumentation. This product is intended for research and forensic applications.

This mixture contains 21 terpenes. Included are:

Item No. 40527	Cannabis Terpenes Mixture 2 (CRM)	
Item Name	CAS Number	Concentration (µg/ml)
(±)-Camphene	79-92-5	2,500
(-)-β-Pinene	127-91-3	2,500
α-Phellandrene	99-83-2	2,500
α-Terpinene	99-86-5	2,500
β-Ocimene	13877-91-3	2,500
Eucalyptol	470-82-6	2,500
γ-Terpinene	99-85-4	2,500
Sabinene hydrate	546-79-2	2,500
(±)-Linalool	78-70-6	2,500
(-)-Isopulegol	89-79-2	2,500
Isoborneol	124-76-5	2,500
(+)-Borneol	464-43-7	2,500
(R)-(+)-Pulegone	89-82-7	2,500
Geranyl Acetate	105-87-3	2,500
α-Cedrene	469-61-4	2,500
trans-β-Farnesene	18794-84-8	2,500
(+)-Valencene	4630-07-3	2,500
trans-Nerolidol	40716-66-3	2,500
(-)-Caryophyllene oxide	1139-30-6	2,500
(-)-α-Bisabolol	23089-26-1	2,500
α-Terpineol	98-55-5	2,500

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, 05/22/2025

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