

PRODUCT INFORMATION



Cannabis Terpenes Mixture 1 (CRM)

Item No. 40526

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

Fill volume: >1 ml

Storage: -20°C

Stability: ≥1 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 1 (CRM) (Item No. 40526) is a certified reference material mixture containing terpenes found in *Cannabis*. The mixture contains α-pinene, (+)-fenchone, camphor, geraniol, and (-)-guaiaol, 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 20 terpenes and terpenoids that have been found in *Cannabis*. Included are:

Item No. 40526	Cannabis Terpenes Mixture 1 (CRM)	
Item Name	CAS Number	Concentration (µg/ml)
α-Pinene	80-56-8	2,500
Sabinene	3387-41-5	2,500
β-Myrcene	123-35-3	2,500
(1S)-(+)-3-Carene	498-15-7	2,500
(R)-(+)-Limonene	5989-27-5	2,500
p-Cymene	99-87-6	2,500
Terpinolene	586-62-9	2,500
(+)-Fenchone	4695-62-9	2,500
(1R)-endo-(+)-Fenchyl alcohol	2217-02-9	2,500
Terpineol (mixture of isomers)	8000-41-7	2,500
Camphor	76-22-2	2,500
Menthol	89-78-1	2,500
Nerol	106-25-2	2,500
Geraniol	106-24-1	2,500
(-)-trans-Caryophyllene	87-44-5	2,500
α-Humulene	6753-98-6	2,500
cis-Nerolidol	3790-78-1	2,500
(+)-Cedrol	77-53-2	2,500
(-)-Guaiaol	489-86-1	2,500
Phytol	7541-49-3	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, 08/01/2024

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