

PRODUCT INFORMATION



Terpene Mixture 2 (CRM)

Item No. 25929

Supplied as: A solution in methanol

Storage: -20°C

Stability: Refer to the Certificate of Analysis

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

Description and Data

Terpene Mixture 2 (CRM) (Item No. 25929) is a certified reference material standard mixture that contains 21 terpenes, including (+)-limonene, α -humulene, α -pinene, and (-)- β -pinene, among others, that have been found in *Cannabis*. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use as a reference standard for these 21 terpenes by GC- or LC-MS. This product is intended for research and forensic applications.

This mixture contains 21 terpenes that have been found in *Cannabis*. Included are:

Item Name	CAS Number	Nominal Concentration ($\mu\text{g/ml}$)
(+)-Borneol	464-43-7	1,000
(-)-Borneol	464-45-9	1,000
(1R)-(+)-Camphor	464-49-3	1,000
(1R)-Endo-(+)-Fenchyl Alcohol	2217-02-9	1,000
(1S)-(-)-Camphor	464-48-2	1,000
(R)-(+)-Limonene	5989-27-5	1,000
α -Pinene	80-56-8	1,000
α -Cedrene	469-61-4	1,000
α -Humulene	6753-98-6	1,000
α -Terpinene	99-86-5	1,000
β -Pinene	127-91-3	1,000
γ -Terpinene	99-85-4	1,000
Geraniol	106-24-1	1,000
L(-)-Fenchone	7787-20-4	1,000
Sabinene	3387-41-5	1,000
Sabinene Hydrate	546-79-2	1,000
Terpineol (mixture of isomers)	8000-41-7	1,000
Terpinolene	586-62-9	1,000
<i>trans</i> -Nerolidol	40716-66-3	1,000
(+)-Pulegone	89-82-7	1,000
(-)-Guaial	489-86-1	1,000

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, 07/19/2022

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