

PRODUCT INFORMATION



Phytocannabinoid Neutrals Mixture 9 (CRM)

Item No. 39009

Supplied as: A solution in methanol (1 mg/ml of each compound)

Fill volume: >1 ml

Storage: -20°C

Stability: ≥1 year

Special conditions: Vortex or sonicate before use

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

Description and Data

Phytocannabinoid Neutrals Mixture 9 (CRM) (Item No. 39009) is a certified reference material mixture containing phytocannabinoid neutral species. The mixture contains Δ^9 -THC, Δ^8 -THC, cannabidiol, and tetrahydrocannabivarin, 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 HPLC-UV and LC-MS instrumentation. Phytocannabinoid Neutrals Mixture 9 contains compounds that are regulated as Schedule I compounds in the United States. Phytocannabinoid Neutrals Mixture 9 (CRM) (Item No. 39009) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

This mixture contains 9 phytocannabinoid neutral species for use in method development and as a reference standard. Included are:

Item No. 39009	Phytocannabinoid Neutrals Mixture 9 (CRM)	
Item Name	CAS Number	Concentration (mg/ml)
Δ^9 -THC	1972-08-3	1.0
Δ^8 -THC	5957-75-5	1.0
Cannabinol	521-35-7	1.0
Cannabidiol	13956-29-1	1.0
(±)-Cannabichromene	20675-51-8	1.0
Cannabidivarin	24274-48-4	1.0
Cannabigerol	25654-31-3	1.0
(±)-Cannabicyclol	N/A	1.0
Tetrahydrocannabivarin	31262-37-0	1.0

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, 03/01/2024

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