

PRODUCT INFORMATION



Phytocannabinoid Acids Mixture 7 (CRM)

Item No. 39010

Supplied as: A solution in acetonitrile:DIPEA (99:1) with 0.05% ascorbic acid (1 mg/ml each compound)

Fill volume: >1 ml

Storage: -20°C

Stability: ≥1 year

Special conditions: Vortex or sonicate before use

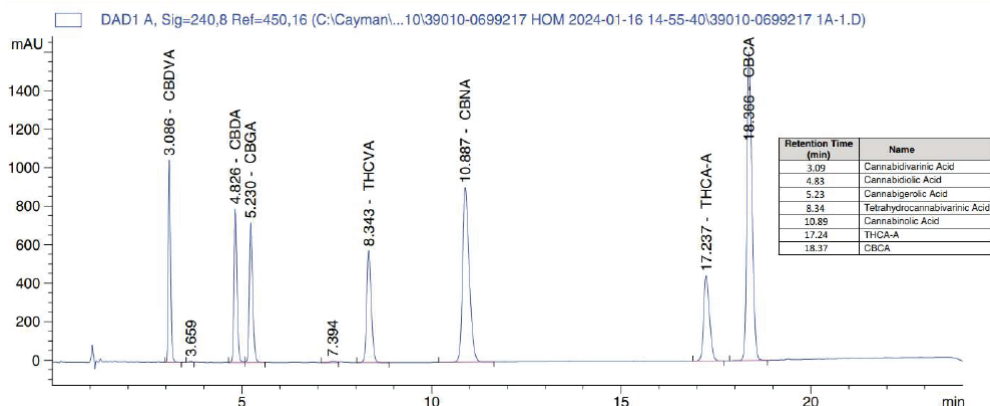
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Phytocannabinoid Acids Mixture 7 (CRM) (Item No. 39010) is a certified reference material mixture that contains compounds that can be isolated from plants in the genus *Cannabis*. The mixture contains THCA-A, cannabinolic acid, cannabidiolic acid, (±)-cannabichromenic acid, cannabidivarinic acid, and cannabigerolic acid, 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 Acids Mixture 7 (CRM) contains compounds that are regulated as Schedule I compounds in the United States. This product is intended for research and forensic applications.

This mixture contains 7 phytocannabinoid acids designed for use in method development and as a reference standard. Included are:

Item No. 39010	Phytocannabinoid Acids Mixture 7 (CRM)	
Item Name	CAS Number	Concentration (mg/ml)
THCA-A	23978-85-0	1.0
Cannabinolic Acid	2808-39-1	1.0
Cannabidiolic Acid	1244-58-2	1.0
(±)-Cannabichromenic Acid	185505-15-1	1.0
Cannabidivarinic Acid	31932-13-5	1.0
Cannabigerolic Acid	25555-57-1	1.0
Tetrahydrocannabivarinic Acid	39986-26-0	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, 03/05/2024

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