

Phytocannabinoid CRM Mixtures



Traceable reference materials are essential for measurement accuracy in *Cannabis* product quality testing and profiling. Cayman provides a collection of pre-made, multi-component CRM mixtures as well as fast turnaround of custom mixtures from our ISO 17034 labs to support quantitation workflows.

ACCREDITED
ISO/IEC 17025 #AT-1773
ISO 17034 #AR-1774

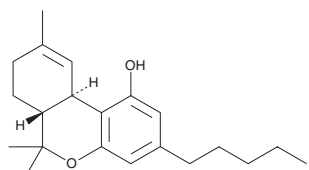
- Features:**
- Reduce time, consumable cost, and reproducibility errors associated with preparing standard curves from individual standards
 - Provides ISO 17034 traceability of measurement results
 - Intended for direct snap-and-inject analysis in HPLC-UV or LC- or GC-MS applications

Improve accuracy
& reproducibility

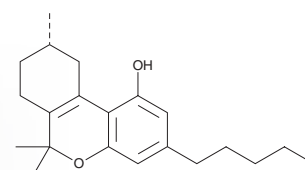
Easy to use &
cost-effective

Streamline quantitation
workflows

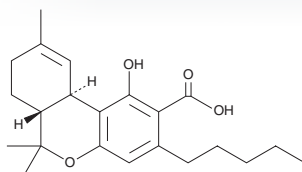
Available in alternate
batches & custom formats



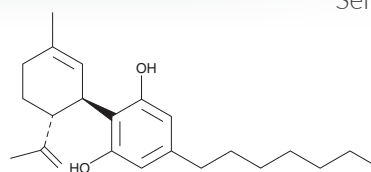
Neutral Phytocannabinoids



Semi-Synthetic Cannabinoids



Phytocannabinoid Acids



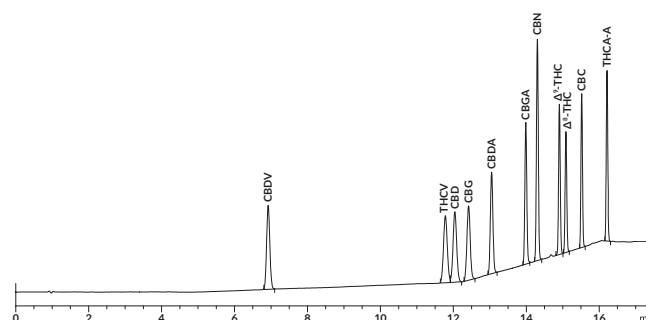
Rare Phytocannabinoids

Cannabis & Hemp Analytical Standards

Explore all products and resources for potency testing, terpene profiling, and analysis of contaminants.

- Phytocannabinoids
- Semi-Synthetic Cannabinoids
- Terpenes & Flavonoids
- Pesticides
- Mycotoxins
- Residual Solvents

Learn more at www.caymanchem.com/cannabis-standards





Phytocannabinoid CRM Mixtures

- ✓ ISO 17034 certified
- ✓ Headspace purged with argon
- ✓ Custom mixtures also available
- ✓ Designed for direct snap-and-inject analysis
- ✓ Exempt product (US)
- ✓ Available in alternate batches
- ✓ Packaged in glass ampules

Analytes Included	Quantify Up to 11 Routinely Tested Cannabinoids							Differentiate Acids & Neutrals			
	11-Mix	10-Mix	9-Mix (CA)	6-Mix	5-Mix	4-Mix	3-Mix	Neutrals 8-Mix	Neutrals 9-Mix	Acids 6-Mix	Acids 7-Mix
CBD	✓	✓	✓	✓	✓	✓	✓	✓	✓		
Δ ⁹ -THC	✓	✓	✓	✓	✓	✓	✓	✓	✓		
CBN	✓	✓	✓	✓	✓		✓	✓	✓		
THCA-A	✓	✓	✓	✓	✓	✓				✓	✓
CBDA	✓	✓	✓	✓	✓					✓	✓
CBG	✓	✓	✓	✓				✓	✓		
CBC	✓	✓	✓					✓	✓		
Δ ⁸ -THC	✓	✓	✓			✓		✓	✓		
CBDV	✓	✓						✓	✓		
THCV	✓		✓					✓	✓		
CBGA	✓	✓								✓	✓
CBDVA										✓	✓
THCVA										✓	✓
CBCA										✓	✓
CBL									✓		
CBNA											✓
Mixture Diluent	acetonitrile						methanol	acetonitrile	methanol	99:1 acetonitrile:DIPEA	
Fill Volume	0.5 ml or 1 ml	1 ml						0.5 ml	1 ml	0.5 ml	1 ml
Available Formats	View Online	View Online	View Online	View Online	View Online	View Online	View Online	View Online	View Online	View Online	View Online

Custom Phytocannabinoid CRM Mixtures

Want adjustments to formulation or components? Cayman offers fast turnaround of custom phytocannabinoid CRM mixtures from our ISO 17034 labs.

- Tailored Formulations
- Curated Component Selection
- Fast Quoting Process

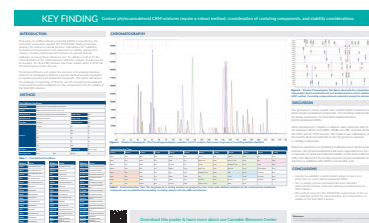
Inquire with our Sales Team at sales@caymanchem.com



Creating Custom *Cannabis* CRM Mixtures: Forty-seven Phytocannabinoids, One HPLC Method

Learn how our scientists developed a robust method to resolve coeluting components for expedited production of custom phytocannabinoid CRM mixtures.

See the method at www.caymanchem.com/cannabis-mixtures-poster

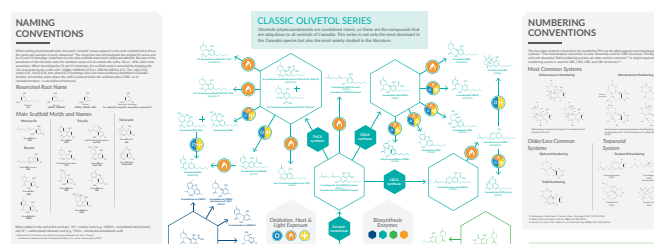


Additional *Cannabis* & Hemp Analytical Testing Resources

Lab Wall Poster

Phytocannabinoid Guide: Biosynthesis, Naming, and Numbering

Request the poster at www.caymanchem.com/phyto-poster



Application Note

GC Separation for Identification of *iso*-THC Contaminants and Accurate Quantification of Δ^8 -THC and Δ^9 -THC in *Cannabis* Samples

See the method at www.caymanchem.com/thc-contaminants



Application Note

HPLC Method to Differentiate Four THC Stereoisomers Formed from Δ^9 -THC Degradation: (6a*R*,9*R*)- Δ^{10} -THC, (6a*R*,9*S*)- Δ^{10} -THC, 9(*R*)- $\Delta^{6a,10a}$ -THC, and 9(*S*)- $\Delta^{6a,10a}$ -THC

Learn more at www.caymanchem.com/thc-stereoisomers