

Synthetic Cannabinoid NPS

Stay current on emerging synthetic cannabinoid NPS with Cayman's reference standards. Cayman works closely with the forensic community to identify illicit synthetic cannabinoids and produce reference standards for the early detection of NPS. Cayman offers additional tools for the identification of unknown compounds, including GC-MS unknown analysis assistance, mass spectrometry resources like the Cayman Spectral Library and GC-MS Drug Identification Tool, as well as our lab wall poster guides.



Indole/Indazole/Pyrazole Synthetic Cannabinoids

Item No.	Product Name
29350	ADB-BUTINACA
33205	ADB-4en-PINACA
37057	CUMYL-CHSINACA
36834	CUMYL-TsINACA
36141	3,5-ADB-4en-PFUPPYCA
36146	5,3-ADB-4en-PFUPPYCA
40205	MDMB-BUTICA
31075	4-fluoro MDMB-BUTICA
28739	MDMB-BUTINACA
26645	4-fluoro MDMB-BUTINACA
39729	MDMB-PICA
20803	5-fluoro MDMB-PICA
26097	MDMB-4en-PINACA
23631	(S)-5-fluoro ADB ((S)-5-fluoro MDMB-PINACA)

OXIZID Synthetic Cannabinoids

Item No.	Product Name
34956	BZO-CHMOXIZID
36946	BZO-HEPOXIZID
34795	BZO-POXIZID
34586	5-fluoro BZO-POXIZID
35503	BZO-4en-POXIZID
10563	MDA 19

See all OXIZIDs at www.caymanchem.com

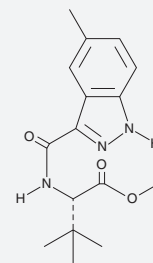
Synthetic Cannabinoid Precursors to Watch

Featured Compound

MDMB-5Me-INACA

Item No. 39658

An indazole precursor reported by the CFSRE in Nov. 2023



Additional Precursors Identified in Casework

Item No.	Product Name
39498	2-(2-Chlorophenyl)-1-(1H-indol-3-yl)ethanone
37729	ADB-INACA
36263	ADB-5Br-INACA
27249	MDMB-ICA
27248	MDMB-INACA
36757	MDMB-5Br-INACA
36883	MMB-5Br-INACA

See all synthetic cannabinoid precursors at www.caymanchem.com

New 2024 Webinar

What's New on the Designer Drug Scene: Synthetic Cannabinoids, Precursors, & More

Learn more about recent NPS developments, including the emergence of synthetic precursors for the illegitimate production of synthetic cannabinoids.

www.caymanchem.com/nps-webinar

IATA Synthetic Cannabinoids

Item No.	Product Name
38136	AB-CHMIATA
36014	ADB-BUTINAATA
35059	ADB-FUBIATA
36853	ADB-FUBINAATA
36992	AFUBIATA
37054	CH-FUBIATA
38283	CH-HEXIATA
36558	CH-PIATA
38205	CHM-FUBIATA

See all IATAs at www.caymanchem.com

Bromo Synthetic Cannabinoids

Item No.	Product Name
37741	ADB-5'Br-4en-PINACA
37963	5-fluoro ADB-5'Br-PINACA
36770	ADB-5'Br-BUTINACA
37721	ADB-5'Br-DECINACA
37758	ADB-5'Br-PINACA
37698	MDMB-5'Br-BUTINACA
39488	MDMB-5'Br-4en-PINACA

See all bromo synthetic cannabinoids at www.caymanchem.com

Sulfamoyl Synthetic Cannabinoids

Item No.	Product Name
37742	A-PONASA
38886	NMDMSB
27325	2-fluoro QMPSPB

OPYRA Synthetic Cannabinoid

Item No.	Product Name
37468	CHO-4'Me-5'Br-FUBOXPYRA



Discover 850+ Synthetic Cannabinoid Standards

www.caymanchem.com/syn-cann-standards

Synthetic & Semi-Synthetic Phytocannabinoids

Item No.	Product Name
37913	9(R)-Hexahydrocannabinol (CRM)
37942	9(S)-Hexahydrocannabinol (CRM)
36346	9(R)-Hexahydrocannabiphorol
36347	9(S)-Hexahydrocannabiphorol
37914	Δ^9 -THC Acetate (CRM)
33011	(6aR,9R)- Δ^{10} -THC (CRM)
33012	(6aR,9S)- Δ^{10} -THC (CRM)
33013	9(R)- $\Delta^{6a,10a}$ -THC (CRM)
33014	9(S)- $\Delta^{6a,10a}$ -THC (CRM)

Visit our *Cannabis & Hemp* Resource Center to Learn More

www.caymanchem.com/cannabistesting

Tools & Resources

ITEM NO: ISO60210

View Item in Cayman's Database

Search Your Findings

Go Back to AD-PROGCA

Request a Quote

Analogous Substances

This compound has at least two head/core/tail pharmacophores that are common

Synthetic Cannabinoid Flipbook Tool

This tool allows you to quickly identify drugs of abuse that share common head/core/tail pharmacophores with known synthetic cannabinoids.

www.caymanchem.com/flipbook



Cayman Spectral Library

Our free, searchable, GC-MS spectral database contains 70eV EI mass spectral data of 2,000+ of Cayman's forensic drug standards.

www.caymanchem.com/CSL