

New Drugs of Abuse Standards

Cayman Chemical is at the forefront of providing crime and toxicology labs with authentic reference standards for new and emerging drugs of abuse. We have more than 2,500 standards to identify, confirm, and quantify illicit opioids, cannabinoids, cathinones, amphetamines, phenethylamines, benzodiazepines, and other controlled substances. Through direct communication with forensic toxicologists, crime labs, and health labs who are tasked with the identification of unknown samples, our product line is constantly evolving to meet emergent needs. Our online drug identification tools and downloadable GC-MS library are freely available to assist in the identification of unknown compounds, and our highly trained staff of chemists is equipped to quickly synthesize currently unavailable forensic standards.



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

Opioids

Item No.	Product Name
27816	Brorphine (hydrochloride)
28070	2',5'-dimethoxy Fentanyl (hydrochloride)
26398	Metonitazene
28689	3,4-difluoro U-47700

Lysergamides

Item No.	Product Name
27818	ALD-52 (solution)
28289	LSD (D-tartrate) (solution)
27817	1B-LSD (solution)
27727	1P-LSD (solution)

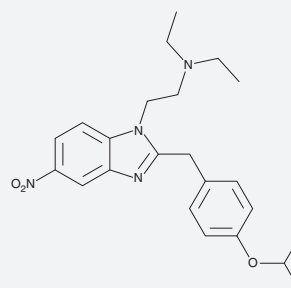
Benzodiazepines

Item No.	Product Name
27748	Bentazepam
24481	Flualprazolam
29206	α -hydroxy Flualprazolam
23524	Fluclotizolam

Stimulants

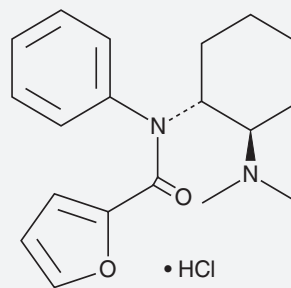
Item No.	Product Name
26596	α -ethyl 2C-D (hydrochloride)
29859	N-(1-methylpentyl)-3,4-DMA (hydrochloride)
28365	4-methyl-N-Ethylhexanophenone (hydrochloride)
27076	N-isopropyl Hexylone (hydrochloride)

Compounds showing up in forensic casework that should be on your radar



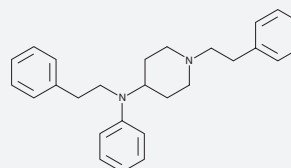
Opioid seen in USA and Canada

Isotonitazene
Item No. 27255



Utopioid seen in OH, FL, VA, and Canada

Furanyl UF-17 (hydrochloride)
Item No. 27924



Fentanyl seen in OH, VA, MD, NJ, CA, and DC.

Phenethyl 4-ANPP
Item No. 28386

Be the first to know about new emerging drug standards by subscribing to our forensic emails at www.caymanchem.com/register

Cannabinoids

Item No.	Product Name
29350	ADB-BUTINACA
28739	MDMB-BUTINACA
28508	MDMB-3en-BUTINACA
28541	MDMB-4en-PICA
29301	NE-CHMIMO

Item No.	Product Name
29210	Δ^9 -THC Acetate
25708	(6aR,9R)- Δ^{10} -THC
25707	(6aR,9S)- Δ^{10} -THC
26528	9(R)- $\Delta^{6a,10a}$ -THC
26529	9(S)- $\Delta^{6a,10a}$ -THC

TOOLS AND RESOURCES

to assist in the identification of emerging drugs of abuse

Cayman Spectral Library

Download the searchable GC-MS spectral database containing 70eV EI MS data of hundreds of Cayman's emerging forensic drug standards.

- Updated quarterly
- Available in Agilent MSD ChemStation and NIST formats
- Freely available for download at www.caymanchem.com/CSL

GC-MS Search Tool

Search unknowns by formula weight, base peak, or 2nd base peak ion at www.caymanchem.com/forensics/search



Search by FW, Base Peak, or 2nd Base Peak Ion



Select Product



View GC-MS

Can't find your unknown? Submit your GC-MS data and any relevant information (suspected chemical class, matrix, solvent, GC parameters, etc.) to techserv@caymanchem.com for our scientists to help investigate.

To view a complete list of our drugs of abuse, visit us online at www.caymanchem.com/forensics