

Traditional Drugs of Abuse & Analytical Standards Panels



The identification of traditional and emerging drugs of abuse is fundamental in forensic and clinical toxicological casework. Cayman provides a wide range of analytical standards for illicit substances and misused medications. We now offer nine comprehensive Analytical Standards Panels that provide coverage of 1,274 traditional and emerging drugs of abuse.

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

Stimulants

Item No.	Product Name
14203	(±)-Amphetamine (hydrochloride)
14204	D-Amphetamine (hydrochloride)
33764	Cocaine (hydrochloride)
11554	(±)-MDA (hydrochloride)
9001123	bk-MDEA (hydrochloride)
34480	3,4-MDMA (hydrochloride)
36566	(±)-Methamphetamine (hydrochloride)
13997	(+)-Methamphetamine (hydrochloride)
13998	(-)-Methamphetamine (hydrochloride)
11639	(±)- <i>threo</i> -Methylphenidate (hydrochloride)

See all stimulants at www.caymanchem.com

Hallucinogens

Item No.	Product Name
11734	2C-B (hydrochloride)
11157	Dextromethorphan
13959	N,N-DMT
33572	Ketamine (hydrochloride)
28289	LSD (D-tartrate) (solution)
11950	Mescaline (hydrochloride)
14276	PCP (hydrochloride)
14041	Psilocybin

See all hallucinogens at www.caymanchem.com

Opioids

Item No.	Product Name
14025	Buprenorphine (hydrochloride)
15459	Codeine
23580	Fentanyl
9001543	Heroin
15461	Hydrocodone
15463	Hydromorphone
37451	(±)-Methadone (hydrochloride)
15464	Morphine
15465	Oxycodone
15911	Oxymorphone

See all opioids at www.caymanchem.com

Anticonvulsants & Analgesics

Item No.	Product Name
22910	Gabapentin
13663	Pregabalin

See all anticonvulsants & analgesics at www.caymanchem.com

Analytical Standards Panels

View on page 2



Fit for screening method development, in-house library building, & sample testing applications.

Bulk Quantities for Traditional and Emerging Drugs of Abuse Available
Email sales@caymanchem.com to Request a Quote

Benzodiazepines

Item No.	Product Name
14255	Alprazolam
14263	Clonazepam
15887	Diazepam
15891	Lorazepam
22577	Triazolam

See all benzodiazepines at www.caymanchem.com

Phytocannabinoids

Item No.	Product Name
12068	Δ^9 -THC
30748	Tetrahydrocannabivarin
33448	THCA-A

See all phytocannabinoids at www.caymanchem.com

Sedatives & Muscle Relaxers

Item No.	Product Name
11152	Etaqualone
9002506	GHB (sodium salt)
9000636	Methaqualone
22450	Promethazine (hydrochloride)
15792	Zolpidem

See all sedatives & muscle relaxers at www.caymanchem.com

Synthetic Cannabinoids

Item No.	Product Name
10900	JWH 018
10904	JWH 073
13636	JWH 398

See all synthetic cannabinoids at www.caymanchem.com

Expand Testing with Analytical Standards Panels

New Synthetic Cannabinoid Standards Panels

Comprises a collection of 571 synthetic cannabinoid standards, including precursors and metabolites, spanning three generations of designer drugs:

Synthetic Cannabinoid Analytical Standards

Panel 3 - Item No. 38770

2017-2025 | 204 Standards

Synthetic Cannabinoid Analytical Standards

Panel 2 - Item No. 38769

2013-2016 | 207 Standards

Synthetic Cannabinoid Analytical Standards

Panel 1 - Item No. 38768

Early 2000s-2012 | 160 Standards

Convenient, cost-effective, & comprehensive

Amphetamine Analytical Standards Panel

Item No. 31615

150 analytical reference materials and standards categorized as stimulants

Benzodiazepine Analytical Standards Panel

Item No. 26260

53 analytical reference materials and standards categorized as benzodiazepines and benzodiazepine metabolites

Phenethylamine Analytical Standards Panel

Item No. 31617

71 analytical reference materials and standards categorized as phenethylamines

Atypical Stimulants Analytical Standards Panel

Item No. 31618

126 analytical reference materials and standards categorized as atypical stimulants

Cathinone Analytical Standards Panel

Item No. 31616

239 analytical reference materials and standards categorized as cathinones and cathinone metabolites

Tryptamine Analytical Standards Panel

Item No. 33685

64 analytical reference materials and standards categorized as tryptamines and tryptamine metabolites

Explore the features of Cayman's Analytical Standards Panels

www.caymanchem.com/forensicpanels