

Prescription & Illicit Drug Certified Reference Materials



Many analgesics clinically prescribed to manage chronic pain carry warranted concerns over safety and abuse liability. Further complications can arise from drug-drug interactions when stimulants or other illicit compounds are also abused. Cayman Chemical offers Certified Reference Materials (CRMs) to help identify, confirm, and quantify many of the major opioids, benzodiazepines, psychoactive drugs, and other controlled substances. Our CRMs meet ISO/IEC 17025 and ISO 17034 standards and are sold as quantitative solutions in sealed ampules.

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

Opioid CRMs

Item No.	Product Name
ISO60128	Acetyl fentanyl (hydrochloride) (CRM)
23060	Acrylfentanyl (hydrochloride) (CRM)
25936	Benzyl fentanyl (hydrochloride) (CRM)
ISO60178	Buprenorphine (hydrochloride) (CRM)
19734	Butyryl fentanyl (hydrochloride) (CRM)
19884	Carfentanil (CRM)
26637	Crotonyl fentanyl (CRM)
23603	Cyclopropyl fentanyl (hydrochloride) (CRM)
ISO60197	Fentanyl (hydrochloride) (CRM)
25615	β -methyl Fentanyl (hydrochloride) (CRM)
19795	FIBF (hydrochloride) (CRM)
19826	<i>para</i> -Fluorobutyryl fentanyl (hydrochloride) (CRM)
19633	Furanyl fentanyl (hydrochloride) (CRM)
ISO60187	Heroin (CRM)
23540	Methoxyacetyl fentanyl (hydrochloride) (CRM)
18567	Mitragynine (CRM)
19635	Ocfentanil (hydrochloride) (CRM)
25887	Phenyl fentanyl (hydrochloride) (CRM)
22429	Remifentanil (hydrochloride) (CRM)
18660	Tapentadol (hydrochloride) (CRM)
19397	U-47700 (CRM)
25522	U-48800 (hydrochloride) (CRM)
23999	U-49900 (CRM)
24264	U-51754 (hydrochloride) (CRM)
19798	Valeryl fentanyl (hydrochloride) (CRM)

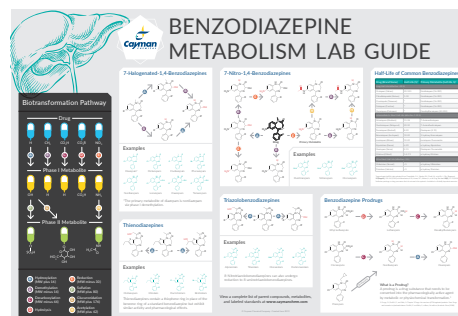
Common opioids listed, more opioids and their metabolites and internal standards available online

Benzodiazepine CRMs

Item No.	Product Name
ISO60185	Alprazolam (CRM)
ISO60181	Clonazepam (CRM)
25614	Clonazolam (CRM)
ISO60177	Diazepam (CRM)
28623	Diclazepam (CRM)
20991	Etizolam (CRM)
25577	Flualprazolam (CRM)
26352	Flubromazolam (CRM)
18160	Flurazepam (CRM)
19391	Midazolam (CRM)
19206	Nimetazepam (CRM)
ISO60172	Phenazepam (CRM)
ISO60182	Temazepam (CRM)

Common benzodiazepines listed, more benzodiazepines and their metabolites and internal standards available online

Request a physical copy of the
**Benzodiazepine Metabolism
Wall Poster for your lab!**



Visit www.caymanchem.com/literature

Stimulant and Hallucinogen CRMs

Item No.	Product Name
ISO60188	(±)-Amphetamine (hydrochloride) (CRM)
25523	BZP (hydrochloride) (CRM)
ISO60176	Cocaine (CRM)
25688	Deschloroketamine (hydrochloride) (CRM)
27239	bk-DMBDB (hydrochloride) (CRM)
29672	Eutylone (hydrochloride) (CRM)
19389	Ketamine (hydrochloride) (CRM)
ISO60189	(±)-MDA (hydrochloride) (CRM)
27240	bk-MDEA (hydrochloride) (CRM)
ISO60190	3,4-MDMA (hydrochloride) (CRM)

Common psychoactives listed, more psychoactive drugs and their metabolites and internal standards available online

Synthetic Cannabinoid CRMs

Item No.	Product Name
27237	AB-CHMINACA (CRM)
ISO60211	AB-FUBINACA (CRM)
19805	4-fluoro ADB (CRM)
25012	(S)-5-fluoro ADB (CRM)
27043	AM2201 (CRM)
28624	APP-BUTINACA (CRM)

Common synthetic cannabinoids listed, more synthetic cannabinoids and their metabolites and internal standards available online

Sedative and Muscle Relaxant CRMs

Item No.	Product Name
ISO60206	Carisoprodol (CRM)
ISO60141	Dextromethorphan (CRM)
20283	Methaqualone (CRM)
20411	Zaleplon (CRM)
20648	Zolpidem (CRM)
20397	Zopiclone (CRM)

View a full list of internal standards online

Antidepressant CRMs

Item No.	Product Name
19031	Amitriptyline (hydrochloride) (CRM)
25576	Bupropion (hydrochloride) (CRM)
23981	Quetiapine (hemifumarate) (CRM)
19049	Trazodone (hydrochloride) (CRM)
25524	Venlafaxine (hydrochloride) (CRM)

View a full list of internal standards online

Item No.	Product Name
26615	(+)-Methamphetamine (hydrochloride) (CRM)
20891	Methylone (hydrochloride) (CRM)
26279	Mexedrone (hydrochloride) (CRM)
ISO60194	PCP (hydrochloride) (CRM)
28958	3-methoxy PCP (hydrochloride) (CRM)
25670	4-methyl Pentedrone (hydrochloride) (CRM)
27504	Pentylone (hydrochloride) (CRM)
24337	N-ethyl Pentylone (hydrochloride) (CRM)
22808	α-Pyrrolidinopentiophenone (hydrochloride) (CRM)
28879	4'-chloro-α-Pyrrolidinovalerophenone (hydrochloride) (CRM)

Item No.	Product Name
17726	5-fluoro CUMYL-PINACA (CRM)
27973	JWH 018 (CRM)
22047	MAB-CHMINACA (CRM)
28186	4-fluoro MDMB-BUTINACA (CRM)
18203	PB-22 (CRM)
ISO60059	XLR11 (CRM)

Barbiturate CRMs

Item No.	Product Name
22912	Amobarbital (CRM)
22854	Barbital (CRM)
20091	Butalbital (CRM)
20966	Pentobarbital (CRM)
20987	Phenobarbital (CRM)
21707	Secobarbital (CRM)

View a full list of internal standards online

Performance-Enhancing Drug CRMs

Item No.	Product Name
ISO60161	Androstenedione (CRM)
19247	Lorcaserin (hydrochloride) (CRM)
21677	Methyldienolone (CRM)
ISO60154	Testosterone (CRM)
9002564	11-keto Testosterone (CRM)

To view a complete list of our Certified Reference Materials,
visit us online at www.caymanchem.com/forensics