

# Thermo Scientific™ CEDIA® / DRI™ Drugs of Abuse Immunoassays

800-232-3342 | [thermoscientific.com/diagnostics](http://thermoscientific.com/diagnostics)

	Generic Names	Trade Names	Appearance	Street Names	Route of Ingestion	% Excreted Unchanged in Urine	Detection Time in Urine	Detection Time in Oral Fluid	Major Compounds / Metabolites	Progressive Symptomatology	
<b>Cannabinoids</b>	<b>Cannabis, Marijuana, Hashish</b>	Marinol, Dronabinol	Leaves, capsules, buds, liquid	Bud, Pot, Grass, Hash, Reefer, Mary Jane, Weed, Dope	Smoked, oral	Hepatic < 1% excreted unchanged in urine			THC-COOH, 11-OH Δ <sup>9</sup> THC, Δ <sup>9</sup> THC Glucuronide	Altered perception, memory loss, disorientation, psychosis	
	<b>Synthetic Cannabinoids</b>	JWH-018, UR-144, XLR-11 (5-F-UR144), AB-PINACA, 5-F-AB-PINACA, AB-CHMINACA, APP-CHMINACA, AB-FUBINACA, AB-CHMICA, ADBICA, and 5-F-ABICA to name a few	Powder, plant extracts	K2, Spice Gold, Spice, Sdiamond, Yucatan Fire, Genie, Fire N Ice, Black Mamba, Bombay Blue, Zombie World, The Moon, Bliss	Snorted, smoked, sold as incense, potpourri	< 1%	Up to 72 Hours	24-48 Hours	Acids, Hydroxylated, Glucuronides	Effects are just like marijuana	
<b>Depressants</b>	<b>BARBITURATES:</b> <b>Amobarbital (A)</b> <b>Butabarbital (B1)</b> <b>Butalbital (B2)</b> <b>Pentobarbital (P)</b> <b>Phenobarbital (Ph)</b> <b>Secobarbital (S)</b>	Amytal, Nembutal, Butisol, Luminal, Seconal, Tuinal, Fiorinal, Neodom, Immenostal, Stadolorm	Capsule, powder, tablets, solution	Blues, Blue Dolls, Rainbows, Yellows, Nebbies, Barbs, Downers, Reds, Red Devils, Seccles	Oral, injected	Hepatic (A) 1-3% (B1) 5-10% (B2) 5% (P) 1% (Ph) 20-35% (S) 5%	1-21 Days	2-3 Days	Amobarbital, Apropobarbital, Butethal, Butabarbital, Cyclopentobarbital, Butalbital, Pentobarbital, Phenobarbital, Secobarbital	Sedation, anti-anxiety, euphoria, anesthesia, decreased mental acuity and slowed speech, respiratory depression, lethargy, stupor, coma	
	<b>BENZODIAZEPINE:</b> <b>Alprazolam (A)</b> <b>Chlordiazepoxide (C1)</b> <b>Clonazepam (C2)</b> <b>Clorazepate (C3)</b> <b>Diazepam (D)</b> <b>Flurazepam (F)</b> <b>Lorazepam (L)</b> <b>Oxazepam (O)</b> <b>Temazepam (T1)</b> <b>Triazolam (T2)</b>	Xanax, Librium, Valium, Clonopin, Dalmane, Ativan, Restoril, Halcion, Tranxene, Tavor, Remestan, Rohypnol, Restoril, Lexotanil, Durayinil, Temestra, Normiston, Flunitrazepam (Rohypnol)	Capsule, powder, tablets, solution	Benzos, Bzd, Tranx, Libs, Vals, Flying V's, Ruffies, Downers, Nerve Pills, Gooftalls, Valley Girl, Heavenly Blues, Stupefy	Oral, injected	Hepatic (A) 20% (C1) < 1% (C2) < 1% (C3) 0% (D) < 1% (F) < 1% (L) < 1% (O) < 1% (T1) 0% (T2) < 1%	Therapeutic: up to 7 Days Chronic: 4-6 Weeks	1-2 Days	α-OH-alprazolam, Bromazepam, Chlorazepate, Chlordiazepoxide, Clonazepam, Delorazepam, Demoxepam, Diazepam, Flunitrazepam, Flurazepam, Halazepam, Lorazepam, Medazepam, Nitrazepam, Nordiazepam, Oxazepam, Prazepam, Temazepam, Triazolam	Hypnotic, muscle relaxation, sedation, dizziness, anti-anxiety, drowsiness, anticonvulsant, lethargy, fatigue, ataxia, coma	
	<b>Methaqualone</b>	Rouquatone, Tualone, Normi-Nox, Mandraz, Quaalude	Capsule, powder, tablets, solution	714, Ropers, Ludes, Soape, Sopors, Love Pill, Buttons	Oral, injected	< 1%	Up to 14 Days	1-10 Days	Methaqualone, Mecloqualone, 3'-Hydroxy Methaqualone, 4'-Hydroxy Methaqualone	Euphoria, sedation, hypnosis, increased sociability and self-confidence. Sexual arousal results from depression of emotional inhibitory centers in the brain	
<b>Hallucinogens</b>	<b>Ethyl Alcohol</b>	Beer, Wine, Liqueur	Color or colorless liquid	Booze, Hooch, Brew	Oral	2-10%	4-6 Hours	6-12 Hours	Ethylglucuronide	Relaxation, impaired motor function, impaired judgement	
	<b>Dextromethorphan</b>	Benlyn, Creomulsion, Delsym, Tussein Pediatric, Vicks 44	Syrup or pills	Dex, Dxm, Robo, Skittles, Syrup	Injection, oral	2.5%	2-4 Days	2-4 Days	O-demethylation (DXO)	Trouble walking straight, the appearance of being drunk	
	<b>Phencyclidine (PCP)</b>	Sernyl, Sernylan	Powder, capsules, colorless crystals	Angel Dust, Space Cadet, Pcp, Hog Dust, White Devil	Smoked, injected, oral, nasal	30-50%			PCP, PCHP, PHP, PCM, TCP, PCE	Dissociative, anesthesia, depression, psychosis, stupor, coma, seizures	
	<b>Lysergic Acid Diethylamide (LSD)</b>	Delysid	White, odorless powder, tablets, capsules, solids, liquids, sugar cubes, gelatin, blotting paper	Acid, Purple Micro Dots, Window Panes, Sugar Cubes, Blotters, Tabs, Wedges, Frisco Special Or Speedball	Oral	1%	8-24 Hours	1-2 Days	d-LSD	Hallucinations, flashbacks, psychosis, vomiting, paralysis, respiratory depression, pupillary dilation, tachycardia, hyperthermia, elevated glucose and fatty acid levels	
	<b>Mescaline (Peyote)</b>	None	Natural-looking "buttons", tablets or capsules	Mesc, Cactus, Buttons	Smoked, oral, injected, nasal	50-60%			Mescaline, TMPEA, Mezcaline, 3,4,5-Trimethoxyphenethylamine, Mezcalin	Hallucinations, nausea, vomiting, abdominal cramps, tremors	
	<b>Ketamine</b>	Ketalar	White powder, clear or cream colored liquid	K, Special K, Super K, Kit-Kat, Kitty	Injected, mixed in beverage, sniffed	2.3%	1-3 Days		Ketamine, Nor-Ketamine	Dream-like state, euphoria, vivid hallucinations, bizarre distortions of reality, irrational behavior, major convulsions, muscle rigidity, respiratory failure	
<b>Narcotics</b>	<b>OPIATES:</b> <b>Morphine (M)</b> <b>Morphine Sulfate</b> <b>Hydromorphone</b>	Heroin (H), Codeine (C), Dilaudid, Lorpham, Hydromorphone, Percodan, OxyContin, Farcodin, Hydrocodone, Oxymorphone	White powder or crystals, clear liquid, tablets, capsules	Atom Bomb (Heroin Mixed With Marijuana) Gooftall Or Speed Ball (Heroin Mixed With Cocaine), H, Horse, Junk, Smack, Skag, Cotton Brothers (Morphine Mixed With Cocaine And Heroin)	Smoked, oral, injected, nasal	(H) < 1% (C) 5-20%	2-4 Days	1-3 Days	Codeine, Morphine-3-β-D-Glucuronide, Morphine-6-β-D-Glucuronide, 6-monoacetylmorphine (6-AM), Hydrocodone, Diacetylmorphine, Hydromorphone	(H)-euphoria, drowsiness, respiratory depression, convulsions, coma, (c)-sedation, analgesia, euphoria, nausea, coma	
	<b>Buprenorphine</b>	Suboxone, Subplex, Buprenex	Tablets, injectable	Bupe	Oral, injection	1%	1-7 Days	1-3 Days	Norbuprenorphine and its glucuronides	Nausea, vomiting, hypotension, hypoventilation, dizziness, meiosis, cramps, muscle aches, irritability	
	<b>Hydrocodone/Hydromorphone</b>	Vicodin, Vicoprofen, Repraxain, ibudone	Pills	Codone, Hydro, Viko, Norco, Vikes	Snort, injection	Detectable amounts after 16 Hours would be 1/16th of dose			Hydromorphone, Hydromorphone-3-β-D-glucuronide	Agitation and delirium, withdrawal symptoms, constipation, respiratory depression	
	<b>Meperidine</b>	Demerol	White crystalline substance	Demmies	Oral, injection	17%		1-2 Days		Delirium, headache, convulsions, hyperthermic	
	<b>Fentanyl</b>	Fentora	Crushed, cake-like, crumbly	Lollipop, Perc-O-Pop, China White, Mexican Brown	Oral, injection, skin patch	0.4-6%			Fentanyl, Norfentanyl, Acetyl Fentanyl, Acryl Fentanyl, Carfentanyl, Furanyl Fentanyl, Valeryl fentanyl, Fluorofentanyl, Methylfentanyl	Euphoria, nausea, vomiting, respiratory depression, coma	
	<b>Methadone</b>	Dolophine, Amidone	Tablets, solution	Dollie, Methadose	Tablets, solution	5-50%	3-7 Days	2-3 Days	Methadone, EDDP (Methadone Metabolite)	Jitteriness, insomnia, convulsions, stupor, respiratory depression, coma	
	<b>Oxycodone/Oxymorphone</b>	Numorphan, OxyContin, Roxiphrwn, Percodan	White to off-white powder, as tablets	Blue Heaven, Blues, Mrs. O, New Blues, Octagons, Oranges, Orgasna Jr, Om, Pink, Pink Heaven, Pink Lady, Pink O, Stop Signs, The O Bomb	Chewing, snorting, injecting	< 1%			Oxymorphone-3-β-D-glucuronide, Oxymorphone	Death with overdose, erythrodermatidemia, withdrawal symptoms, respiratory depression, vomiting, constipation	
	<b>Propoxyphene</b>	Darvon, Dolene, Darvocet, Distalgic, Doloxene	Tablets, capsules, white crystalline powder	Darvon Cocktail	Oral, injected	1.1%	1-2 Days	1-2 Days	Propoxyphene, Norpropoxyphene	Analgesia, stupor, respiratory depression, seizures, coma	
	<b>Tapentadol</b>	Nucynta, Tapal, Palaxia	Oral suspension, pills	Cha Cha	Snort, oral	1%			N-desmethyl, O-desmethyl and Glucuronide	Life threatening respirator depression, hypotensive, seizures, anapylaxis, renal and hepatic impairment	
	<b>Tilidine</b>	Valoron, Valtran, Tilidin	Tablets, liquid	Lak, Tilsa, Valoron, Tildate	Oral, injection		Up to 7 Days		A 1 to 1 mixture of (1R,2S)- and (1S,2R)-Ethyl 2-(N,N-dimethylamino)-1-phenylcyclohex-3-ene-carboxylate	Respiratory depression, ventricular tachycardia	
	<b>Tramadol</b>	Ultram®, Ultracet, ConZip, Rybix ODT, Ryzoit	Capsules	Chill Pills, Ultras	Injection	30%	2-4 Days	1-4 Days	N-desmethyl, N,N-desmethyl, O-desmethyl, O,N-desmethyl and Glucuronide (60%)	Fibromyalgia, aummentation of restless leg syndrome	
	<b>Stimulants</b>	<b>Cocaine</b>	None	Powder, crystals	Coke, Snow, Crack, Flake, Blow, Toot, Cola, White Girl	Smoked, oral, injected, nasal	< 10%	2-5 Days	2-5 Days	Benzoylctgonine, Cocaethylene	Anesthesia, euphoria, confusion, depression, convulsions, cardiotoxicity
		<b>Amphetamine</b>	Adderall, Benzedrine, Dexedrine	Tablets, capsules, solution	Speed, Bennies, Uppers, Black Beauties	Smoked, oral, injected, nasal	~30%	1-4 Days	2-3 Days	4-Hydroxyamphetamine, Phenylacetone	Mood swings, malnutrition, delusion, psychosis, increased heart and breathing rates, increased blood pressure, fatal strokes and heart failure
<b>Bath Salts</b>		Cathinone / Methylendioxypropylveralone / Mephedrone	White, tan, or brown colored powdery substance	Ivory Wave, Plant Fertilizer, Vanilla Sky, Energy-1, Red Dove, White Dove, Blue Silk, Zoom, Hurricane, Aura	Ingest, smoke, snort, or inject	> 80%			Pyrolidine and its Glucuronides	Euphoria, talkative, increased motivation, severe agitation, depression, hallucinations, anxiety	
<b>Methamphetamine</b>		Desoxyn, Methedrine	Tablets, capsules, solution	Doe, Crystal, Speed, Ice, Meth, Tina	Smoked, oral, injected, nasal	~10-20%	1-4 Days	Up to 24 Hours	d-Methamphetamine (S+)-Methamphetamine	Mood swings, malnutrition, delusion, psychosis, increased heart and breathing rate, increased blood pressure, fatal strokes and heart failure	
<b>Methylphenidate</b>		Ritalin	Tablets, capsules, patches, and liquid.	Kibbles And Bits, Diet Coke, Kiddy Coke, R Pop, Study Buddies, Pineapple, Poor Man's Cocaine, Skippy, Vitamin R	Snort powder, oral (dissolve in water), injected	20%			Ritalinic Acid (80%)	Agitation, psychosis, dysrhythmias, malignant hypertension, severe anxiety, tension, temporary paralysis, and hallucinations, impulsiveness, may also have learning and behavioral problems	
<b>Ecstasy</b>	None	White powder, capsules, tablets	E, Ecstasy, X-Tc, Adam, Skittles, Tabs, Pills, Molly	Oral, injected	65%			MDMA, MDEA, MDA, etc.	Muscle tension, nausea, blurred vision, fainting, alternating between chills and sweating, increased heart rate and blood pressure		
<b>Sleep Aids</b>	<b>Gamma Hydroxybutyric acid (GHB) &amp; γ-butyrolactone</b>	G; Sodium Oxybate; Xyrem, Alcover	White crude powder (hygroscopic)	G* (Most Common), Gamma-OH, Liquid E, Fantasy, Georgia Home Boy Liquid X, Harm, Grievous Bodily, Mills, Juice, Lollipops, Goop	Oral	< 5%	Up to 12 Hours	Up to 5 Hours	D-2-hydroxyglutaric acid, 4,5-dihydroxyhexanoic acid	Unrousable sleep, dizziness, nausea, and/or vomiting, fatal in an unconscious person. Sleep apnea, antiaging medicine, decreases body fat	
	<b>Zopiclone</b>	Imovane in Canada, Australia, Sweden, Finland, Norway, Zimovane in Europe, Ximovan in Germany	Fine white or slightly cream crystalline powder	Zimmers, Zimmies, Zim Zim, Zoppies	Smoked, oral	< 7%			N-desmethyl and N-oxide	Polychondritis, menopause, menopausal symptoms, hormone replacement	
	<b>Zolpidem</b>	Zolpimist, Edluar, Stilnox, and Sublinox	White to off-white crystalline powder	Z Drugs	Oral, intranasal	< 1%			Zolpidem 4-Phenyl Carboxylic Acid (ZCA), Zolpidem 6-Carboxylic Acid	Sedative-hypnotic, anxiety, aggression, agitation, confusion, unusual thoughts, hallucinations, memory problems, changes in personality	
	<b>Zaleplon</b>	Starnoc and Andante	Capsule, pill	Z Drugs	Intranasal, oral	< 1%			Glucuronides	Management/treatment of insomnia, unsteadiness, severe dizziness severe, weak muscle tone, complex sleep related behaviors	
<b>Muscle Relaxant</b>	<b>Carisoprodol</b>	Soma, Prosoma	Pills	Ds, Dance, Las Vegas Cocktail	Bumped, injected, or smoked	1%	2-7 Days	1-7 Days	Meprobamate	Visual disturbances, anxiety, fatigue	

\*Approximation of detection time. Actual detection time is dependent on dose, frequency of use and individual metabolism

Contact us at 800-232-3342 or email at [sales.diagnostics.fmt@thermofisher.com](mailto:sales.diagnostics.fmt@thermofisher.com)  
Find out more at [thermofisher.com/diagnostics](http://thermofisher.com/diagnostics)

© 2017 Thermo Fisher Scientific Inc. All rights reserved. CEDIA is a registered trademark of Thermo Scientific Diagnostics. All other trademarks are the property of Thermo Fisher Scientific or its subsidiaries.  
10023849-IMTL-DAT-POSTER-EN