

IVD For In Vitro Diagnostic Use

Rx Only

REF	Quantity
10015632	(3 x 18 mL Kit)
100248	(70 mL Kit)
100249	(500 mL Kit)

Intended Use

The DRI® Oxycodone Assay is intended for the qualitative and semi-quantitative determination of oxycodone in human urine at cutoffs of 100 and 300 ng/mL. The assay provides a simple and rapid analytical screening procedure to detect oxycodone in human urine.

This assay provides only a preliminary analytical test result. A more specific alternative chemical method must be used in order to obtain a confirmed analytical result. Gas chromatography/mass spectrometry (GC/MS) is the preferred confirmatory method. Clinical consideration and professional judgment should be applied to any drug of abuse test result, particularly when preliminary results are used.

Summary and Explanation of the Test

Oxycodone is a semi-synthetic opioid prescribed for pain management in patients with moderate to severe pain. It is similar to codeine and morphine in its analgesic properties but it is more potent than morphine and has higher dependence potential. The drug oxycodone is supplied as OxyContin® (Oxycodone HCl) or in combination with aspirin (Percodan®) or acetaminophen (Percocet®).¹ Drug abusers crush the pills into powder and snort them for faster effect which may result in a potentially fatal outcome. According to Drug Abuse Warning Network (DAWN), there has been a dramatic increase in oxycodone related deaths.^{2,3} Oxymorphone, noroxycodone and noroxymorphone are the only known metabolites of oxycodone.² The metabolite, oxymorphone, is a potent narcotic analgesic, while the other two metabolites are relatively inactive. From 33-61% of a single dose of oxycodone is excreted in urine within 24 hours as unconjugated oxycodone (13-19%), conjugated oxycodone (7-29%), and conjugated oxymorphone (13-14%).⁴

The DRI Oxycodone Assay is supplied as a liquid ready-to-use homogeneous enzyme immunoassay. The assay uses specific antibodies that can detect oxycodone and oxymorphone without any significant cross-reactivity to other opiate compounds. The assay is based on competition between a drug labeled with glucose-6-phosphate dehydrogenase (G6PDH), and free drug from the urine sample for a fixed amount of specific antibody binding sites. In the absence of free drug from the sample, the specific antibody binds the drug labeled with G6PDH and causes a decrease in enzyme activity. This phenomenon creates a direct relationship between the drug concentration in urine and enzyme activity. The enzyme activity is determined spectrophotometrically at 340 nm by measuring the conversion of nicotinamide adenine dinucleotide (NAD) to NADH.

Reagents

Antibody/Substrate Reagent:

Contains mouse monoclonal anti-oxycodone derivative antibody, glucose-6-phosphate (G6P), and nicotinamide adenine dinucleotide (NAD) in Tris buffer with sodium azide as a preservative.

Enzyme Conjugate Reagent:

Contains oxycodone derivative labeled with glucose-6-phosphate dehydrogenase (G6PDH) in Tris buffer with sodium azide as a preservative.

Additional Materials Required (sold separately):

REF	Kit Description
1664	DRI Negative Calibrator, 10 mL
1388	DRI Negative Calibrator, 25 mL
100250	DRI Oxycodone Calibrator 100, 10 mL
100251	DRI Oxycodone Calibrator 300, 10 mL
100252	DRI Oxycodone Calibrator 500, 10 mL
100253	DRI Oxycodone Calibrator 1000, 10 mL
DOAT-2	MAS® DOA Total – Level 2
DOAT-3	MAS® DOA Total – Level 3
DOAT-4	MAS® DOA Total – Level 4
DOAT-5	MAS® DOA Total – Level 5

⚠ Precautions and Warnings

This test is for in-vitro diagnostic use only. The reagents are harmful if swallowed.

DANGER: DRI Oxycodone Assay contains ≤0.2% bovine serum albumin (BSA) and ≤0.5% Drug-specific antibody (Mouse).

H317 - May cause allergic skin reaction.

H334 - May cause allergy or asthma symptoms or breathing difficulties if inhaled.

Avoid breathing mist or vapor. Contaminated work clothing should not be allowed out of the workplace. Wear protective gloves/eye protection/face protection. In case of inadequate ventilation wear respiratory protection. If on skin: Wash with plenty of soap and water. If INHALED: If breathing is difficult, remove victim to fresh air and keep at rest in a position comfortable for breathing. If skin irritation or rash occurs: Get medical advice/attention. If experiencing respiratory symptoms: Call a POISON CENTER or doctor/physician. Wash contaminated clothing before reuse. Dispose of contents/container to location in accordance with local/regional/national/international regulations.

Reagents used in the assay components contain ≤0.09% sodium azide. Avoid contact with skin and mucous membranes. Flush affected areas with copious amounts of water. Get immediate medical attention for eyes, or if ingested. Sodium azide may react with lead or copper plumbing to form potentially explosive metal azides. When disposing of such reagents, always flush with large volumes of water to prevent azide build-up. Clean exposed metal surfaces with 10% sodium hydroxide.

Do not use the reagents beyond their expiration dates.

Reagent Preparation and Storage

The reagents are ready-to-use. No additional reagent preparation is required. The reagents should be stored refrigerated (2-8°C). All assay components, opened or unopened, are stable until the expiration date indicated on their respective labels. Do not use the reagents beyond their expiration dates.

Specimen Collection and Handling

Collect urine specimens in plastic or glass containers. Testing of fresh urine specimens is suggested.

Specimens kept at room temperature that do not receive initial test within 7 days⁵ of arrival at the laboratory may be placed into a secure refrigeration unit at 2 to 8°C for up to two months.⁶ For longer storage prior to analysis or for sample retention after analysis, urine specimens may be stored at -20°C.^{6,7}

Laboratories following the SAMHSA mandatory guidelines should refer to SAMHSA “Short-Term Refrigerated Storage” and “Long-Term Storage” requirements.⁸

To protect the integrity of the sample, do not induce foaming and avoid repeated freezing and thawing. An effort should be made to keep pipetted samples free of gross debris. It is recommended that grossly turbid specimens be centrifuged before analysis. Frozen samples should be thawed and mixed prior to analysis. Adulteration of the urine sample may cause erroneous results. If adulteration is suspected, obtain another sample and forward both specimens to the laboratory for testing.

Handle all urine specimens as if they were potentially infectious.

Assay Procedure

Analyzers capable of maintaining a constant temperature, pipetting samples, mixing reagents, measuring enzymatic rates at 340 nm and timing the reaction accurately can be used to perform this assay.

Refer to the specific application instructions of each analyzer for chemistry parameters before performing the assay.

Quality Control and Calibration

Qualitative analysis

For qualitative analysis of samples, use either the Oxycodone 100 Calibrator, or the Oxycodone 300 calibrator, as a cutoff level.

Semi-quantitative analysis

For semiquantitative analysis, use all calibrators.

Good laboratory practice suggests the use of control specimens to ensure proper assay performance. Use controls near the cutoff calibrator to validate the calibration. Control results must fall within established ranges, as determined by laboratory procedures and guidelines. If results fall outside of established ranges, assay results are invalid. All quality control requirements should be performed in conformance with local, state and/or federal regulations or accreditation requirements.

Calibration Frequency

Recalibration is recommended

- After calibrator or reagent lot change
- After instrument maintenance is performed
- As required following quality control procedures

See below for calibration frequency recommendations for Hitachi analyzers. For other analyzers, refer to the instrument-specific application sheet.

NOTE: Reassess control targets and ranges following a change of reagent lot.

The potential cross-reactivity posed by drugs commonly coadministered with oxycodone was evaluated by adding each substance to oxycodone free urine at the concentration indicated. A drug was considered to cross-react if the observed oxycodone concentration exceeded 100 ng/mL, the lowest cutoff for the DRI Oxycodone Assay. As shown in the tables below, all of the pharmacologic compounds evaluated, including a number of the opiate compounds, exhibited no cross-reactivity at the concentrations listed. Please note that some substances including 6-acetyl morphine, codeine, dihydrocodeine, heroin, hydrocodone, hydromorphone, levorphanol, naloxone, naltrexone gave results between 75 and 99 ng/mL i.e., within 25% of the cutoff of 100 ng/mL.

Structurally related opiate compounds that tested negative at 100 ng/mL cutoff.

Compound	Concentrations (µg/mL)
6-Acetyl Morphine	75
Codeine	500
Dihydrocodeine	200
Heroin	300
Hydrocodone	200
Hydromorphone	40
Levorphanol	200
Morphine	350
Morphine-3-glucuronide	950
Naloxone	300
Norcodeine	1,000
Normorphine	1,000

Structurally unrelated compounds that tested negative at 100 ng/mL cutoff.

Compound	Concentrations (µg/mL)
Acetaminophen	1,000
Acetylsalicylic acid	1,000
Amitriptyline	500
Amoxicillin	500
Amphetamine	2,000
Benzoylcegonine	2,000
Caffeine	1,000
Carbamazepine	1,000
Chlorpromazine	2,000
Clomipramine	1,000
Cimetidine	1,000
Desipramine	1,000
Dextromethorphan	200
Doxepine	200
Ephedrine	2,000
Fentanyl	200
Fluoxetine	1,000
Fluphenazine	500
Ibuprofen	1,000
Imipramine	1,000
Maprotiline	1,000
Meperidine	1,000
Methadone	1,000
Metroniazole	2,000
Nalbuphine	1,000
Nortriptyline	500
Oxazepam	500
Phencyclidine	1,000
Phenobarbital	1,000
Ranitidine	3,000
Secobarbital	1,000
Talwin	500
Thebaine	20
Thioridazine	1,000
Tramadol	500

Interference

The potential interference of pH and endogenous physiologic substances on recovery of oxycodone using the DRI Oxycodone Assay was assessed by spiking known amounts of potentially interfering substances into the low (225 ng/mL) and high (375 ng/mL) controls for the 300 ng/mL cutoff. Oxycodone concentration for each specimen (substance and final concentration noted in the table below) was determined and the percent recovery calculated as the quotient of the spiked to control value. The table below shows the substance and final concentration at which the observed oxycodone concentrations for spiked controls are within ±1% of the expected control dose. No interference was observed by the addition of the compounds upto the concentrations listed below.

Compound	Concentrations (mg/dL)
Acetone	1,000
Ascorbic Acid	1,500
Creatinine	500
Ethanol	1,000
Galactose	10
Glucose	3,000
Hemoglobin	300
Human Serum Albumin	500
Oxalic Acid	100
Riboflavin	7.5
Sodium chloride	1,000
Urea	2,000
pH	3-11

References

1. Anderson D.T., Fritz K.L., and Muto J.J. OxyContin®: The concept of a “Ghost Pill” and the Postmortem Tissue Distribution of Oxycodone in 36 Cases. *J. Anal. Toxicol.* 2002, 26: 448-459.
2. Clinical & Forensic Toxicology News, Oxycodone: Recognition and Pharmacogenomics. By Jannetto P.J. and Gock S.B. March 2003.
3. Cone E.J., et al, Oxycodone Involvement in Drug Abuse Deaths: A DAWN-Based Classification Scheme applied to an Oxycodone Postmortem Database Containing over 1000 Cases. *J. Anal. Toxicol.* 2003, 27: 57-67.
4. Oxycodone. In: Baselt R.C. and Cravey R.H. Disposition of toxic drugs and chemicals in man, 4th ed. Chemical Toxicology Institute, Foster City, California: 1995: 572-574.
5. Dixon RB, Mbeunkui F, Wiegel JV. Stability Study of Opioids and Benzodiazepines in Urine Samples by Liquid Chromatography Tandem Mass Spectrometry. *Journal of Analytical Science and Technology*, December 2015, 6:17.
6. Gonzales E, Ng G, Pesce A, West C, West R, Mikel C, Llaatyshv, S, Almazan P. Stability of pain-related medications, metabolites and illicit substances in urine. *Clinica Chimica Acta* 416: (2013) 30-35.
7. C52-A2, Toxicology and Drug Testing in the Clinical Laboratory; Approved Guideline – Second Edition, *Clinical and Laboratory Standards Institute (CLSI)* (April 2007)
8. *Notice of Mandatory Guidelines for Federal Workplace Drug Testing Program: Final Guidelines; Federal Register*, Substance Abuse and Mental Health Administration (SAMHSA), (1994) 110 (June 9):11983
9. Data on file at Microgenics, a part of Thermo Fisher Scientific.

Glossary:

<http://www.thermofisher.com/symbols-glossary>



Microgenics Corporation
46500 Kato Road
Fremont, CA 94538 USA
US Customer and
Technical Support:
1-800-232-3342



EC REP

B-R-A-H-M-S GmbH
Neuendorfstrasse 25
16761 Hennigsdorf, Germany



For insert updates go to:
www.thermofisher.com/diagnostics

Other countries:

Please contact your local Thermo Fisher Scientific representative.

10008282-11-EN
2020 11

thermo
scientific