

# CEDIA® Multi-Drug OFT Control Set B

This test kit is intended solely for use in employment and insurance testing, and does not include test systems intended for Federal drug testing programs (e.g., programs run by the Substance Abuse and Mental Health Services Administration (SAMHSA), the Department of Transportation (DOT), and the U.S. military.)

**REF** 10022356 CEDIA® Multi-Drug OFT Control Set B (2 x 15 mL)

## Intended Use

The Thermo Scientific CEDIA Multi-Drug OFT Control Set B is unassayed quality control material intended for use in the CEDIA Barbiturates, Benzodiazepines, Methadone, and Oxycodone OFT Assays to detect and monitor systematic deviations from accuracy resulting from reagent or instrument defects.

## Description of CEDIA Multi-Drug OFT Control Set B

The CEDIA Multi-Drug OFT Control Set B is liquid ready-to-use. The two controls (low and high) are prepared by spiking secobarbital, nordiazepam, methadone, and oxycodone into buffer matrix.

Low control	-50% of cutoff
High control	+50% of cutoff

The controls are packaged as a kit and may be used with any reagent lot.

## Precautions and Warning

The CEDIA Multi-Drug OFT Control Set B is for Criminal Justice and Forensic use only. They are harmful if swallowed. Do not use the controls beyond their expiration dates printed on their respective labels.

**DANGER:** The controls contain  $\leq 0.22\%$  Bovine Serum Albumin (BSA).

H317 - May cause allergic skin reaction.

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

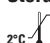
Avoid breathing mist or vapor. Contaminated work clothing should not be allowed out of the workplace. Wear protective gloves/eye protection/face protection. In case of inadequate ventilation wear respiratory protection. If on skin: Wash with plenty of soap and water. IF INHALED: If breathing is difficult, 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.

The controls contain  $\leq 0.09\%$  Sodium azide. Avoid contact with skin and mucous membranes. Refer to SDS for additional precautions, handling instructions, and accidental exposure treatment.

In the case of accidental spill, clean and dispose of material according to your laboratory's SOP, local, and state regulations.

In the case of damaged packaging on arrival, contact your technical support representative (refer to contact number listed in this Package Insert).

## Storage

 The controls should be stored refrigerated at 2° to 8°C when not in use. They are stable until the expiration date indicated on the box label as long as they remain unopened. Once opened, they are stable for 60 days from the date of opening.

## Instruction For Use

The CEDIA Multi-Drug OFT Control Set B is liquid ready-to-use. Mix the contents of the vial before each use by gentle inversion. Remove the cap and dispense. After each use, tightly close each cap and return to refrigerated storage (2-8°C).

## Assay Procedure

For instructions, refer to the instrument-specific application sheet for the CEDIA OFT Assays for Barbiturates, Benzodiazepines, Methadone, and Oxycodone.

## Interpretation of Results

The results of the controls should be within the range established by each laboratory.

## Limitations

The CEDIA Multi-Drug OFT Control Set B is unassayed and designed for use with the CEDIA OFT Assays. Values are not assigned by the manufacturer. Ensure that control results are within the established ranges determined by laboratory practices and guidelines. All quality control requirements should be performed in conformance with local, state, and/or federal regulations or accreditation requirements.

## Bibliography

1. CEDIA Barbiturates OFT Assay Package Insert
2. CEDIA Benzodiazepines OFT Assay Package Insert
3. CEDIA Methadone OFT Assay Package Insert
4. CEDIA Oxycodone OFT Assay Package Insert

## Glossary:

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



Microgenics Corporation  
46500 Kato Road  
Fremont, CA 94538 USA  
1-800-232-3342

© 2018 Thermo Fisher Scientific Inc. All rights reserved.

CEDIA® is a registered trademark of Roche Diagnostics.



For insert updates go to:  
[www.thermoscientific.com/diagnostics](http://www.thermoscientific.com/diagnostics)