

CEDIA® Buprenorphine Calibrators and Controls

Thermo
SCIENTIFIC

IVD For In Vitro Diagnostic Use

Rx Only

REF 100241 CEDIA Buprenorphine S1 Calibrator (0 ng/mL)
100242 CEDIA Buprenorphine S2 Calibrator (5 ng/mL)
100243 CEDIA Buprenorphine S3 Calibrator (20 ng/mL)
100244 CEDIA Buprenorphine S4 Calibrator (50 ng/mL)
100245 CEDIA Buprenorphine S5 Calibrator (75 ng/mL)
100246 CEDIA Buprenorphine Low (3 ng/mL) and High (7 ng/mL) Controls

Intended Use

The CEDIA® Buprenorphine calibrators are used to calibrate the CEDIA Buprenorphine Assay in human urine.

The CEDIA Buprenorphine controls are used to qualify the CEDIA Buprenorphine Assay in human urine.

Each calibrator and control set is sold separately and may be used with any reagent lot. The ready-to-use calibrators or controls contain processed human urine, buffer salts, and 0.09% sodium azide.

Precautions and Warnings

Exercise the normal precautions required for handling all laboratory reagents.



WARNING: The calibrators and controls contain 0.09% sodium azide. Avoid contact with skin and mucous membranes. Flush affected areas with copious amount of water. Get immediate medical attention for eyes, or if ingested. Azide may react with lead or copper plumbing to form potentially explosive metal azides. When disposing of such reagents, always flush with large amount of water to prevent azide build-up. Clean exposed metal surfaces with 10% sodium hydroxide.

Storage and Stability

Store calibrators or controls at 2-8°C. **Do not freeze.** For stability of the unopened calibrators or controls, refer to the box or bottle labels for the expiration date. For stability of the opened calibrators or controls, 60 days or until the printed expiration date, whichever comes first.

Instructions for Use

The calibrators or controls are liquid ready-to-use. Mix the contents of the vial before each use by gently inverting the vial 2-3 times.

Record on the vial label the date on which the calibrator or controls was opened.

Remove the cap and dispense the required quantity of the calibrators or controls into a sample cup as specified in the appropriate CEDIA assay application sheet.

Avoid cross-contamination of calibrator or control levels by matching colored caps to the appropriately colored dropper tips.

Quality Control

All quality control requirements should be performed in conformance with local, state and/or federal regulations or accreditation requirements.

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.thermoscientific.com/diagnostics

Other countries:

Please contact your local representative.

CEDIA is a registered trademark of Roche Diagnostics.

10007591-7-EN
2018 01