

MGC Select DAU Control Set

IVD For In Vitro Diagnostic Use

REF 100202 Low 3 x 5 mL
High 3 x 5 mL

Intended Use

The MGC Select DAU Control Set consists of unassayed High and Low controls intended for use in the validation of drug of abuse assays in human urine.

Description of Controls

The MGC Select DAU Control Set is sold as ready-to-use liquid controls formulated in a human urine matrix containing stabilizers, preservatives and drugs in the following targeted concentrations:

Drug	Low (ng/mL)	High (ng/mL)
6-Acetylmorphine	7.5	12.5
Benzoyllecgonine	112.5	187.5
LSD	0.3	0.7
MDMA	375	625
Nitrazepam	150	250

Precautions and Warnings

Exercise the normal precautions required for handling all laboratory reagents.

DANGER: MGC Select DAU Control Set contains $\leq 0.3\%$ 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 MGC Select DAU Control Set contains $< 0.1\%$ 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 a large volume of water to prevent azide build-up.

Materials of human origin used in the manufacture of the MGC Select DAU Control Set were tested for HIV 1 and 2, Hepatitis B, and Hepatitis C using test approved by the FDA. The findings were negative. However, as no test method can rule out the potential risk of infection with absolute certainty, the material must be handled just as carefully as the patient sample. In the event of exposure, the directives of the responsible health authorities should be followed.

Storage and Stability

The MGC Select DAU Control Set should be stored refrigerated at 2-8°C and is stable until the expiration date indicated on the label.

Instructions for Use

The MGC Select DAU Control Set is liquid ready-to-use. Mix the contents of the vial before each use by gently inverting the vial 2-3 times.

Remove the cap and dispense the required quantity of the control into a sample cup as specified in the appropriate assay package insert or application sheet.

Quality Control

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

Materials provided:

MGC Select DAU Control Set

Additional materials required:

Drugs of Abuse Reagent Kits

CEDIA® is a registered trademark of ROCHE Diagnostics



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
Tel: +49 (0) 800 404 077 12
Fax: +49 (0) 800 404 077 13



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

Other countries:

Please contact your Thermo Fisher Scientific representative.