

**Urinalysis Controls** 





# Comprehensive QC Solutions for Urinalysis testing

Urinalysis is one of the most widely performed tests in routine health screening. Analysis of a patient's urine can provide physicians with valuable information about their health status before symptoms appear and can be used to support the diagnosis of several diseases including Urinary Tract Infections and Diabetes. Using reliable, independent Quality Controls to monitor your Urinalysis testing can give you confidence reporting patient results.

Bio-Rad Urinalysis controls are available in a variety of convenient packaging options, including dropper bottles, flip-top squeeze bottles and centrifuge tubes. Bio-Rad Urinalysis controls provide an extensive menu of analytes to monitor physical, chemical and microscopic examination of urine specimens, all suitable for use with manual or automated methods including dipstick or microscopic analysis.

For greater confidence in the reliability of patient test results, manage your QC data and access peer group comparisons with our Unity Interlaboratory Program.

### **qUAntify Advance Control**



- Available in an easy-to-use plastic squeeze bottle
- Dipstick tests include Pregnancy and Microalbumin
- Human urine solution
- 18 month shelf life at 2-8°C
- 31 day open-vial stability at 2-25°C, including ketones

#### **Analytes**

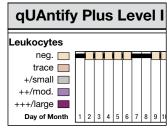
Albumin Ascorbic Acid Bilirubin Blood Creatinine Glucose Hemoglobin (HGB) Ketones Leukocytes Microalbumin Nitrite pH Pregnancy (hCG) Protein (Total) Protein-to-Creatinine Ratio Specific Gravity Urobilinogen

## **qUAntify Plus Control**



- Choose from a conical 12 mL tube or 120 mL flip top squeeze bottle
- Both options can be used for urine dipstick and microscopic testing
- Liquid
- 2 year shelf life at 2-8°C
- 31 day open-vial stability at 2-25°C for most analytes

**POCT Methods Listed** 



**Example Peer Group Report** 

#### **Analytes**

Albumin Albumin-to-Creatinine Ratio Ascorbic Acid\* Bilirubin Blood

\*Tested and found to be negative.

Creatinine Glucose Hemoglobin (HGB) Ketones Leukocytes Microalbumin Microscopics (RBC, WBC, Crystals, Casts) Nitrite pH Pregnancy (hCG) Protein (Total) Protein-to-Creatinine Ratio Specific Gravity Urobilinogen

### **Liquichek Urinalysis Control**



- Ready-to-use urine control formulated to monitor the precision of dipstick and microscopic analyzers
- Available in 12 mL glass vials
- Liquid, human urine based
- 30 day open-vial stability at 2–25°C
- 2.5 year shelf life at 2–8°C

#### Analytes

Bilirubin Blood Clarity Color Creatinine Glucose Ketones Leukocytes Microalbumin Microscopics (RBC, WBC, Crystals, Casts)

Nitrite Osmolality pH Pregnancy (hCG) Protein (Total) Protein-to-Creatinine Ratio Specific Gravity Urobilinogen



# Unity QC Data Management Solutions

Bio-Rad Quality Controls are part of the Unity Program with its industry-leading features for benchmarking laboratory performance.

- Benefit from peer data generated from more than 50,000 instruments worldwide
- Utilize real-time access to peer group comparisons to help with troubleshooting
- Consolidate your entire laboratory into one QC system
- Facilitate run-time decisions with optimized Westgard rules
- Facilitate compliance with simple tools to review rule violations and document corrective actions
- Define your own quality requirements and track key quality performance indicators

You can consolidate your entire laboratory into one QC system through our Unity Program. This sophisticated software includes best-in-class tools to support releasing results, benchmarking laboratory performance, and designing your QC system.

# **Ordering Information**

|                          | Cat #    | Description                                      | Cat #                                      | Description                         |
|--------------------------|----------|--|--|-------------------------------------|
| qUAntify Advance Control |          | Liquichek Urinalysis Control                     |  |                                     |
|                          | 12007028 | Bilevel (3 of each level)                        | 435  | Bilevel (6 of each level)12 x 12 mL |
|                          | 12007029 | Bilevel MiniPak (1 of each level)2 x 12 mL       | 436  | Level 1                             |
|                          |          |  | 437  | Level 2                             |
| qUAntify Plus Control    |          | 435X   | Bilevel MiniPak (1 of each level)2 x 12 mL |                                     |
|                          | 995      | Bilevel (5 of each level)/Tube                   |  |                                     |
|                          | 995X     | Bilevel MiniPak (1 of each level)/Tube 2 x 12 mL |  |                                     |
|                          | 962      | Bilevel (2 of each level)/Bottle 4 x 120 mL      |  |                                     |
|                          | 963      | Level 1/Bottle                                   |  |                                     |



EQA

An independent, external assessment of performance in comparison to your peers.



#### Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.



#### Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.



964

962X

Bio-Rad Laboratories, Inc.

Bilevel MiniPak (1 of each level)/Bottle .....2 x 120 mL

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/urinalysiscontrols

Clinical Diagnostics Group Web site www.bio-rad.com/qualitycontrol USA 1 800 224 6723 Australia 61 2 9914 2800 Austria 43 1 877 8901 Belgium 32 03 710 53 00 Brazil 55 31 3689 6600 Canada 1 514 334 4372 China 86 21 61698500 Czech Republic 420 241 430 532 Denmark 45 4452 1000 Finland 358 9 804 22 00 France 33 1 47 95 60 00 Germany 49 0 89 318 840 Greece 30 210 7774396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 1800 180 1224 Israel 972 3 9636050 Italy 39 02 216091 Japan 81 3 6361 7070 Korea 82 2 3473 4460 Mexico 52 55 5488 7670 The Netherlands 31 318 540666 New Zealand 64 9 415 2280 Norway 47 23 38 41 30 Poland 48 22 3319999 Portugal 351 21 472-7700 Russia 7 495 721 1404 Singapore 65 6415 3170 South Africa 27 11 442 85 08 Spain 34 91 590 5200 Sweden 46 8 555 127 00 Switzerland 41 0 26 674 55 05 06 Taiwan 886 2 2578 7189 Thailand 662 651 8311 United Kingdom 44 0 20 8328 2000