



Urinalysis Controls





Urinalysis Controls from Bio-Rad

Versatile packaging and comprehensive analytes for today's laboratory

Urinalysis is widely documented as the first clinical laboratory test developed and dates back over 6,000 years. Even today, modern urinalysis is still one of the most widely performed tests in routine health screening.

Bio-Rad Urinalysis controls are available in a variety of convenient packaging options, including dropper bottles, flip-top squeeze bottles and centrifuge tubes.

Covering an extensive menu of analytes to monitor physical, chemical and microscopic examination of urine specimens, Bio-Rad Urinalysis controls are well suited for manual or automated methods for dipstick or microscopic analysis of urine specimens.

Participating in the Unity™ Interlaboratory Program provides valuable peer group comparisons with distinctive, easy-to-read color coded reports.

Liquichek™ Urinalysis Control

Ready to use Urine control formulated to monitor the precision of your dipstick and microscopic analyzers. Available in a 12 mL glass vial.

- Liquid, human urine based
- 30 day open-vial stability at 2–25°C
- 2.5 year shelf life at 2–8°C
- Color-coded peer group reports available

[POCT Methods Listed](#)

Analytes

Bilirubin	Nitrite
Blood	Osmolality
Clarity	pH
Color	Pregnancy (hCG)
Creatinine	Protein (Total)
Glucose	Protein-to-Creatinine Ratio
Ketones	Specific Gravity
Leukocytes	Urobilinogen
Microalbumin	
Microscopics (RBC, WBC, Crystals, Casts)	

qUAntify® Control

Try our easy to use and store qUAntify® Control that comes in a squeeze dropper 12 mL plastic bottle.

- Liquid
- 2 year shelf life at 2-8°C
- 31 day open-vial stability at 2-25°C
- Color-coded peer group reports available

[POCT Methods Listed](#)

Analytes

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

qUAntify® Plus Control

Choose from our popular, laboratory-testing friendly conical 12 mL qUAntify® Plus Control or for the cost-conscious laboratory, the 120mL flip top easy to use 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
- Color-coded peer group reports available

[POCT Methods Listed](#)

Analytes

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

qUAntify Plus Level I	
Leukocytes	
neg.	
trace	
+/small	
++/mod.	
+++/large	
Day of Month	1 2 3 4 5 6 7 8 9 1

Example Peer Group Report



*Tested and found to be negative.

Refer to myinserts.com or the package insert of currently available lots for specific analyte and stability claims.

Ordering Information

Cat #	Product Name	Configuration	Cat #	Product Name	Configuration
Liquichek™ Urinalysis Control			qUAntify® Control		
435	Bilevel (6 of each level)	12 x 12 mL	975	Bilevel (3 of each level)	.6 x 12 mL
436	Level 1	12 x 12 mL	975X	Bilevel MiniPak (1 of each level)	.2 x 12 mL
437	Level 2	12 x 12 mL			
435X	Bilevel MiniPak (1 of each level)	.2 x 12 mL	qUAntify® Plus Control		
			995	Bilevel (5 of each level)/Tube	10 x 12 mL
			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	.4 x 120 mL
			964	Level 2/Bottle	.4 x 120 mL
			962X	Bilevel MiniPak (1 of each level)/Bottle	.2 x 120 mL

Partner with the Experts in Quality Control



EQAS®

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



Independent QC

Ongoing, proactive, unbiased daily QC that identifies 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.



**Bio-Rad
Laboratories**

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

**Clinical
Diagnostics Group**

Website www.bio-rad.com/qualitycontrol **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** +32 (3)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