



Liquichek™ Urine Chemistry Control

BIO-RAD

Liquichek™ Urine Chemistry Control

A liquid control ideal for monitoring urine chemistry testing procedures. It provides assayed values for urine tests commonly performed on chemistry analyzers and some manual methods.

10 mL Fill Size

- Liquid, human urine based
- 2 year shelf life at 2–8°C
- 30 day open-vial stability at 2–8°C
- Assayed for quantitative microalbumin and visual pregnancy tests

2.5 mL Fill Size

- Dimension Vista® vials offer ready-to-use convenience for on-board refrigerated storage
- Offers 13 analytes to match the Dimension Vista® test menu
- Shared lot numbers and Unity™ peer groups with existing Urine Chemistry configurations
- 2D barcode technology for quick and easy product insert and vial label scanning
- 2 year shelf life at 2–8°C
- 30 day open-vial (punctured) stability for all analytes when stored 2–8°C



Analytes

Amylase	Glucose	Phosphorus	Specific Gravity*
Calcium	Magnesium	Potassium	Urea*
Chloride	Microalbumin	Pregnancy (Qualitative)*	Urea Nitrogen
Cortisol*	Osmolality*	Protein (Total)	Uric Acid
Creatinine	pH*	Sodium	

* Not included for use on the Dimension Vista® test menu. Refer to the package insert of currently available lots for specific analyte and stability claims

Ordering Information

Cat #	Liquichek™ Urine Chemistry Control	Quantity
397	Level 1	12 x 10 mL
398	Level 2	12 x 10 mL
395X	Bilevel MiniPak (1 of each level)	2 x 10 mL
195	Level 1	12 x 2.5 mL
196	Level 2	12 x 2.5 mL
198X	Bilevel MiniPak (1 of each level)	2 x 2.5 mL



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

Discover the power of Unity™ Solutions at www.QCNet.com.



**Bio-Rad
Laboratories**

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

**Clinical
Diagnostics Group**

Website www.bio-rad.com/qualitycontrol **U.S.** 1-800-224-6723 **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400 **Canada** 1-514-334-4372 **China** 86-21-64260808 **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** 1-800-180-1224 **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** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000