



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