



Lyphocheck® Quantitative Urine Control



Lyphochek® Quantitative Urine Control

Contains a comprehensive analyte menu for quantitative assays that monitor the precision for general urine chemistry, biogenic amines and hormones testing. It also contains hCG (for qualitative pregnancy tests) and Microalbumin at levels suitable for monitoring early proteinuria and associated nephropathy.



- Human urine based
- 28 month shelf life at 2–8°C for most analytes
- Comprehensive analyte menu and method listing

Analytes

3-Methoxytyramine	Calcium	Hydroxyproline (Free)*	Potassium
5-Aminolevulinic Acid	Chloride	Hydroxyproline (Total)	Pregnancy (Qualitative)
5-HIAA	Copper*	Iron	Pregnanetriol*
11-β-Hydroxy-Androsterone*	Coproporphyrin	Lead	Protein (Total)
11-β-Hydroxy-Etiocholanolone*	Cortisol (Free)	Magnesium	Sodium
11-Ketoandrosterone*	Creatine*	Mercury	Specific Gravity
11-Ketoetiocholanolone*	Creatinine	Metanephrine	Urea
17-Hydroxycorticosteroids	Cystine*	Metanephrine (Total)	Urea Nitrogen
17-Ketogenic Steroids	Dehydroepiandrosterone*	Microalbumin	Uric Acid
17-Ketosteroids	Dopamine	Norepinephrine	Uroporphyrins
Aldosterone	Epinephrine	Normetanephrine	VMA
AMP (Cyclic)	Estriol (Total)*	Osmolality	Zinc*
Androsterone*	Etiocholanolone*	pH	
Arsenic	Glucose	Phosphorus	
Barbiturates*	HVA	Porphyrins (Total)	

* No claims are made regarding performance or stability.
Refer to the package insert of currently available lots for specific analyte and stability claims

Ordering Information

Cat #	Lyphochek® Quantitative Urine Control	Quantity
376	Level 1, Normal12 x 10 mL
377	Level 2, Abnormal12 x 10 mL
375X	Bilevel MiniPak (1 of each level)2 x 10 mL



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve your laboratory quality 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. 1-800-2BIO-RAD 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 91-124-4029300 Israel 972-3-9636050
Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico 52(55)5200-0520 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 65-6415-3188 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