



Liquichek™ Unassayed Chemistry Control (Human)

BIO-RAD

Liquichek™ Unassayed Chemistry Control (Human)

A comprehensive, liquid stable control designed to monitor the precision of general and special chemistry tests.

- Liquid, human serum based
- 2 year shelf life at -20 to -70°C
- 15 day open-vial stability 2–8°C
- Suitable for use with most major chemistry instruments



Analytes

Acetaminophen	Chloride	Glucose	Osmolality	T3 Uptake/T-Uptake
Acid Phosphatase (Total)	Cholesterol (Total)	HBDH	PAP*	T4 (Free)*
Albumin	CO ₂	IgA	Phenobarbital	T4 (Total)
Aldolase	Copper	IgG	Phenytoin	Theophylline
Alkaline Phosphatase	Cortisol	IgM	Phospholipids	Tobramycin
ALT/SGPT	Creatine Kinase (CK)	Iron	Phosphorus	Transferrin
Amylase	Creatinine	Iron (TIBC)	Potassium	Triglycerides
AST/SGOT	Digoxin	Iron (UIBC)*	Protein Electrophoresis	TSH
Bilirubin (Direct)*	Ethanol	Lactate (Lactic Acid)	Protein (Total)	Urea Nitrogen
Bilirubin (Indirect)*	Folate	LAP-Arylamidase*	Pseudocholesterase	Uric Acid
Bilirubin (Total)	Gentamicin	LDH	Salicylate	Valproic Acid
Calcium	GGT	Lipase	Sodium	Vancomycin
Calcium (Ionized)*	GLDH*	Lithium	T3 (Free)*	Vitamin B ₁₂
Carbamazepine	Globulin	Magnesium	T3 (Total)	Zinc

* 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 #	Liquichek™ Unassayed Chemistry Control (Human)	Quantity
691	Level 125 x 10 mL
692	Level 225 x 10 mL
690X	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 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.** 1-800-2BIO-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** 36-1-455-8800 **India** 91-124-4029300 **Israel** 972-3-9636050 **Italy** +39-02-218091 **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** 65-6415-3188 **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