



# Multiquel<sup>®</sup> Chemistry Controls



## Multiquel® Chemistry Controls

Multiquel® Chemistry Controls are liquid, human serum based general chemistry controls that are available in both assayed and unassayed configurations and offer a wide range of claimed analytes. These trilevel controls contain more than 80 analytes, including:

- HDL, LDL Cholesterol
- Ionized Calcium
- Low Osmolality value in Level 1
- 11 of the most commonly tested TDM analytes
- 3 year shelf life at -20°C to -70°C
- Thawed and unopened stability for 30 days at 2–8°C for most analytes
- 14 day open-vial stability at 2–8°C for most analytes
- Each level sold separately



### Analytes

Acetaminophen	Ceruloplasmin	IgA	Protein (Total)
Acid Phosphatase	Chloride	IgG	Salicylate
Albumin	Cholesterol (HDL)	IgM	Sodium
Alpha-1-Antitrypsin	Cholesterol (LDL)	Iron	T3 Uptake/T-Uptake
Alpha-Hydroxybutyrate	Cholesterol (Total)	Iron (TIBC)	T3 (Total)
Dehydrogenase (αHBDH)	Cholinesterase	Iron (UIBC)	T4 (Free)
Alkaline Phosphatase	CK-MB Isoenzyme	Lactate Dehydrogenase (LDH)	T4 (Total)
ALT/SGPT	Complement C3	Lactate (Lactic Acid)	Theophylline
Amikacin	Complement C4	LAP Arylamidase	Tobramycin
Amylase	Copper	Lipase	Transferrin
Amylase (Pancreatic)	Cortisol	Lithium	Triglycerides
Apolipoprotein A-1	Creatine Kinase (CK)	Magnesium	TSH
Apolipoprotein B	Creatinine	Osmolality	Urea
AST/SGOT	Digoxin	PAP	Urea Nitrogen
Bilirubin (Direct)	Ethanol	Phenobarbital	Uric Acid
Bilirubin (Neonatal)	Ferritin	Phenytoin	Valproic Acid
Bilirubin (Total)	Gentamicin	Phospholipids	Vitamin B12
Calcium (Ionized)	GGT	Phosphorus	Zinc
Calcium (Total)	Globulin	Potassium	
Carbamazepine	Glucose	Prealbumin	
Carbon Dioxide (CO <sub>2</sub> )	Haptoglobin	Protein Electrophoresis	

Refer to the package insert of currently available lots for specific analyte and stability claims

### Ordering Information

Cat #	Liquid Assayed Multiquel®	Quantity	Cat #	Liquid Unassayed Multiquel®	Quantity
694	Level 1 .....	12 x 3 mL	697	Level 1 .....	12 x 10 mL
695	Level 2 .....	12 x 3 mL	698	Level 2 .....	12 x 10 mL
696	Level 3 .....	12 x 3 mL	699	Level 3 .....	12 x 10 mL
695X	Trilevel MiniPak (1 of each level) .....	3 x 3 mL	698X	Trilevel MiniPak (1 of each level) .....	3 x 10 mL



These controls are featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory, 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](http://www.QCNet.com).**



**Bio-Rad  
Laboratories**

*For further information, please contact your local Bio-Rad office or learn more online at [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol)*

**Clinical  
Diagnostics Group**

Website [www.bio-rad.com/qualitycontrol](http://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 +39-02-216091 Japan 81-3-6361-6902 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