



# Lyphochek® Unassayed Chemistry Control (Human)

**BIO-RAD**

## Lyphochek® Unassayed Chemistry Control (Human)

Provides a comprehensive analyte menu and good stability for enzymes and CO<sub>2</sub>. It requires no special diluent and is compatible with major instruments and methodologies. Available in 5 mL and 10 mL fill sizes.



- Human serum based
- 3 year shelf life at 2–8°C
- 30 day frozen reconstituted stability for most analytes
- 7 day open-vial stability at 2–8°C for most analytes

### Analytes

|   |                      |                      |                         |                         |
|---|----------------------|----------------------|-------------------------|-------------------------|
| Acetaminophen                               | Bilirubin (Direct)   | Cortisol             | Lactate (Lactic Acid)   | T3 (Total)              |
| Acid Phosphatase (Total)                    | Bilirubin (Indirect) | Creatine Kinase (CK) | LAP-Arylamidase         | T3 Uptake/T-Uptake      |
| AFP   | Bilirubin (Total)    | Creatinine           | LDH                     | T4 (Free)               |
| Albumin                                     | Calcium              | Digoxin              | Lipase                  | T4 (Total)              |
| Alkaline Phosphatase (ALP)                  | Calcium (Ionized)    | Gentamicin           | Lithium                 | TBG                     |
| Alpha-1-Antitrypsin                         | Carbamazepine        | GGT                  | Magnesium               | Theophylline            |
| Alpha-Hydroxybutyrate Dehydrogenase (αHBDH) | CEA                  | GLDH                 | Osmolality              | Tobramycin              |
| ALT/SGPT                                    | Ceruloplasmin        | Globulin             | PAP                     | Transferrin             |
| Amylase                                     | Chloride             | Glucose              | Phenobarbital           | Triglycerides           |
| Amylase (Alpha)                             | Cholesterol (HDL)    | Haptoglobin          | Phenytoin               | TSH                     |
| Amylase (Pancreatic)                        | Cholesterol (LDL)    | hCG-Beta Subunit     | Phosphorus              | Urea                    |
| Apolipoprotein A-1                          | Cholesterol (Total)  | IgA                  | Potassium               | Urea Nitrogen           |
| Apolipoprotein B                            | Cholinesterase       | IgG                  | Protein Electrophoresis | Uric Acid               |
| AST/SGOT                                    | CO <sub>2</sub>      | IgM                  | Protein (Total)         | Valproic Acid           |
|   | Complement C3        | Iron                 | Salicylate              | Vancomycin              |
|   | Complement C4        | Iron (TIBC)          | Sodium                  | Vitamin B <sub>12</sub> |
|   | Copper               | Iron (UIBC)          | T3 (Free)               | Zinc                    |

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

### Ordering Information

| Cat #    | Lyphochek® Unassayed Chemistry Control (Human) | Quantity   |
|----------|--|------------|
| 731      | Level 1  | 25 x 5 mL  |
| 732      | Level 2  | 25 x 5 mL  |
| 730X     | Bilevel MiniPak (1 of each level)              | 2 x 5 mL   |
| C-320-10 | Level 1  | 50 x 10 mL |
| C-325-10 | Level 2  | 50 x 10 mL |
| 323X     | 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](http://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](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