



Lyphocheck® Assayed Chemistry Control



Lyphocek® Assayed Chemistry Control

Requires no special diluent and has good frozen reconstituted stability for enzymes and CO₂. It is assayed for more than 100 methodologies and over 30 instruments.

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

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

Ordering Information

Cat #	Lyphocek® Assayed Chemistry Control	Quantity
C-310-5	Level 1	12 x 5 mL
C-315-5	Level 2	12 x 5 mL
313X	Bilevel MiniPak (1 of each level)	2 x 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. 1-800-2BIO-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 81-3-6361-7070 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