



Lyphocheck[®] Immunology Plus Control



Lyphochek® Immunology Plus Control

With a broad menu of serum proteins and related analytes, this human serum based control provides the laboratory with an independent assessment on various test systems. Two levels are available for precision monitoring within the assay range. This lyophilized product offers an excellent 3 year shelf life to allow for long term QC monitoring and less frequent lot crossovers.



- Contains 22 serum proteins and related analytes
- Lyophilized, human serum based
- 3 year shelf life at 2–8°C
- 15 day reconstituted stability at 2–8°C

Analytes

α-1-Acid Glycoprotein	Ceruloplasmin	IgA	Prealbumin
α-1-Antitrypsin	Complement C1q*	IgE	Retinol Binding Protein (RBP)
α-2-Macroglobulin	Complement C3	IgG	Rheumatoid Factor
β-2-Microglobulin	Complement C4	IgM	Transferrin
Albumin	CRP	Kappa Light Chain	
Antistreptolysin O	Haptoglobin	Lambda Light Chain	

* 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® Immunology Plus Control	Quantity
430	Bilevel (6 of each level)	12 x 1 mL
430X	Bilevel MiniPak (1 of each level)	2 x 1 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** + 36-1-459-6100 **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