



# Liquichek™ Immunology Control



## Liquichek™ Immunology Control

A liquid, human serum based control available in a choice of 1mL or 3 mL fill sizes to best suit the needs of your laboratory. Each of the three levels may be purchased separately, or used in combination, to allow monitoring of assay performance at desirable levels. This comprehensive serum protein control includes an impressive menu of 32 analytes with values provided for an extensive selection of widely used test methods.



- Liquid convenience
- Human serum
- 32 serum protein related analytes
- Three distinct levels
- Values provided for wide range of methods
- 2 year shelf life at -20 to -70°C
- 30 day open-vial stability at 2-8°C
- Available in 1mL and 3 mL vials

### Analytes

α-1-Acid Glycoprotein	Apolipoprotein B	Haptoglobin	Lipoprotein (a)
α-1-Antitrypsin	C1 Inhibitor	Hemopexin*	Prealbumin
α-2-Macroglobulin	Ceruloplasmin	IgA	Properdin Factor B
β-2-Microglobulin	CH50*	IgE	Protein (Total)
ADNase B	Complement C3	IgG	Retinol Binding Protein (RBP)
Albumin	Complement C4	IgG Subclasses 1-4	Rheumatoid Factor
Antistreptolysin O	CRP	IgM	Soluble Transferrin Receptor (sTfR)*
Antithrombin III	Cystatin C	Kappa Light Chain	Transferrin
Apolipoprotein A-1	Ferritin	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 #	Liquichek™ Immunology Control	Quantity
591	Level 1 . . . . .	6 x 1 mL
592	Level 2 . . . . .	6 x 1 mL
593	Level 3 . . . . .	6 x 1 mL
590X	Trilevel MiniPak (1 of each level) . . . . .	3 x 1 mL
594	Level 1 . . . . .	6 x 3 mL
595	Level 2 . . . . .	6 x 3 mL
596	Level 3 . . . . .	6 x 3 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** +36-1-459-6100 **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