



# Liquichek™ Elevated CRP Control



## Liquichek™ Elevated CRP Control

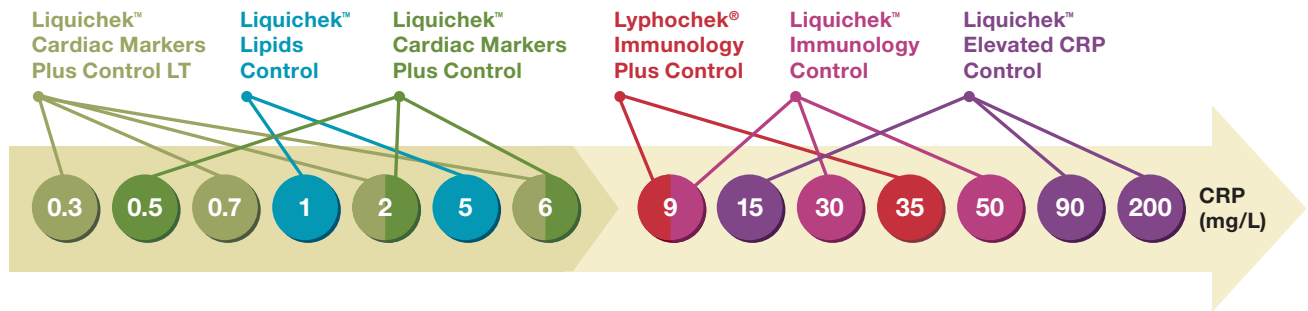
A liquid, human serum control designed to monitor the precision of C-Reactive Protein (CRP) assays as a general marker for infection and inflammation. The control is available in three individual levels to monitor the higher end of the linear range of CRP assays or CRP dilution.

- Liquid, human serum based
- 30 day open-vial stability at 2–8°C
- 3 year shelf life at -20°C to -70°C
- Target CRP levels of 15, 90, 200 mg/L



### Target CRP Level in Bio-Rad Controls

Values reflect typical target values for these products. Assay values may vary depending on methodology.



### Analyte

CRP

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

### Ordering Information

Cat #	Liquichek™ Elevated CRP Control	Quantity
254	Level 1	12 x 1 mL
255	Level 2	12 x 1 mL
256	Level 3	12 x 1 mL
255X	Trilevel MiniPak (1 of each level)	3 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](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/immunocontrols](http://www.bio-rad.com/immunocontrols)**

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