



Liquichek™ Elevated CRP Control



Liquichek™ Elevated CRP Control

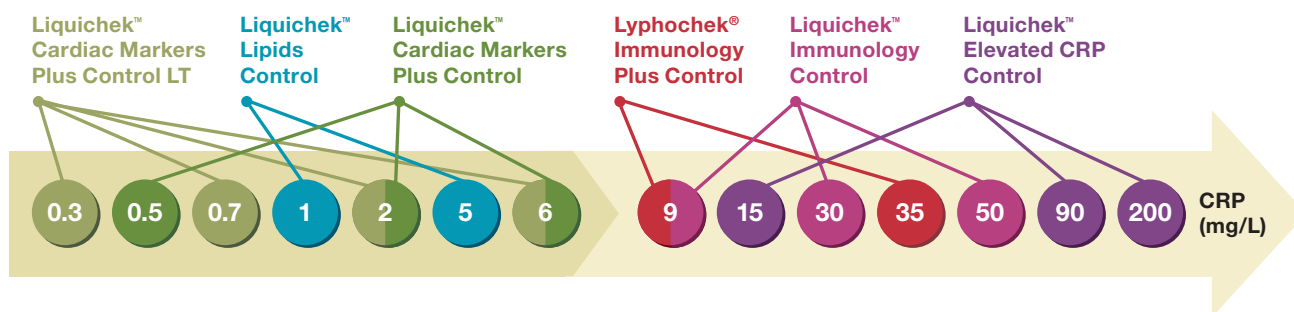
A liquid, human serum control designed to monitor the precision of C-Reactive Protein (CRP) assays as a general marker for infection and inflammation. The control is available in three individual levels to monitor the higher end of the linear range of CRP assays or CRP dilution.

- 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.



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/immunocontrols

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