



Liquichek™ Pediatric Control



Liquichek™ Pediatric Control

Designed to monitor the precision of testing procedures for analytes commonly run for neonates. This control monitors four Bilirubin parameters at two clinically significant levels.



- Liquid
- 3 year shelf life at -20°C to -70°C
- 3 month stability when stored unopened at 2–8°C
- 14 day open-vial stability at 2–8°C

Analytes

Bilirubin (Direct)	Calcium	Potassium
Bilirubin (Indirect)	Chloride	Sodium
Bilirubin (Neonatal)	Glucose	Theophylline
Bilirubin (Total)	Magnesium	
Caffeine	Phenylalanine	

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

Ordering Information

Cat #	Liquichek™ Pediatric Control	Quantity
354	Level 1	.6 x 4 mL
355	Level 2	.6 x 4 mL
353X	Bilevel MiniPak (1 of each level)	.2 x 4 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/chemserumcontrols

*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