



# Liquichek™ Spinal Fluid Control



## Liquichek™ Spinal Fluid Control

A liquid, bilevel product designed to monitor the performance of automated and manual procedures for the analysis of proteins and other critical analytes assayed in cerebrospinal fluid.



- Liquid
- Oligoclonal banding in Level 2
- 2 year shelf life at 2–8°C
- 30 day open-vial stability at 2–8°C

### Analytes

Albumin	IgG	Protein Electrophoresis*
Chloride	IgM	Protein (Total)
Glucose	Lactate (Lactic Acid)	Sodium
IgA	LDH	

\* Prealbumin, Albumin, Alpha-1-Globulin, Alpha-2-Globulin, Beta-Globulin, Gamma-Globulin.  
Refer to the package insert of currently available lots for specific analyte and stability claims

### Ordering Information

Cat #	Liquichek™ Spinal Fluid Control	Quantity
751	Level 1 . . . . .	.6 x 3 mL
752	Level 2 . . . . .	.6 x 3 mL
750X	Bilevel MiniPak (1 of each level) . . . . .	.2 x 3 mL



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve your laboratory quality 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