



# Liquichek Ethanol/Ammonia Control

**BIO-RAD**

## Liquichek Ethanol/Ammonia Control

A liquid control used to monitor the precision of Ethanol and Ammonia test procedures in the clinical laboratory.

- Liquid
- Normal, elevated and toxic concentrations of Ammonia
- 2 year shelf life at 2–8°C
- 20 day open-vial stability at 2–8°C



### Analytes

Ammonia  
Ethanol

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

### Ordering Information

Cat #	Description	
544	Level 1 . . . . .	.6 x 3 mL
545	Level 2 . . . . .	.6 x 3 mL
546	Level 3 . . . . .	.6 x 3 mL
545X	Trilevel MiniPak (1 of each level) . . . . .	.3 x 3 mL

#### EQAS

An independent, external assessment of laboratory performance in comparison to your peers.

[QCNet.com/eqas](http://QCNet.com/eqas)

#### Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.

[QCNet.com/independentqc](http://QCNet.com/independentqc)

#### Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

[QCNet.com/datamanagement](http://QCNet.com/datamanagement)



**Bio-Rad  
Laboratories, Inc.**

*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/diagnostics](http://www.bio-rad.com/diagnostics) **Australia** +61(2)9914-2800 **Austria** +43-1-877-8901 **Belgium** +32(3)710-53-00 **Brazil** +55(31)3689-6600 **Canada** 1-514-334-4372 **China** +86 21 6169 8500 **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** +1800-180-1224 **Israel** +972-3-9636050 **Italy** +39-02-216091 **Japan** +81-3-6361-7070 **Korea** +82-2-3473-4460 **Mexico** +52(55)5488-7670 **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** +7-495-721-1404 **Singapore** +65-6415-3170 **South Africa** +27-11-442-85-08 **Spain** +34-91-590-5200 **Sweden** +46-8-555-127-00 **Switzerland** +41(0)26-674-55-05/06 **Taiwan** +886-2-2578-7189 **Thailand** +662-651-8311 **United Kingdom** +44(0)20-8328-2000