



Liquichek™ Diabetes Control



Liquichek™ Diabetes Control

A liquid human whole blood based product designed to monitor the performance of hemoglobin tests associated with diabetes monitoring, including Hemoglobin A1C and Total Hemoglobin. Available in three critical levels specific to diabetes. With a 1 mL fill size and an open-vial stability of 14 days, waste is minimized. A 3 year shelf life reduces the frequency of lot crossovers, resulting in significant reagent cost savings.



- Liquid convenience
- Human whole blood based control
- Higher third level to monitor your assay range
- Convenient 1 mL fill size
- Three individual levels to fit your preference
- 3 year shelf life at -10°C to -70°C
- 6 month shelf life at 2–8°C
- 14 day open-vial stability at 2–8°C
- Value assignment listings for over 30 methods

Analytes

Hemoglobin A1C
Hemoglobin, Total

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

Ordering Information

Cat #	Liquichek™ Diabetes Control	Quantity
171	Level 1	6 x 1 mL
172	Level 2	6 x 1 mL
173	Level 3	6 x 1 mL
172X	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/diabetescontrols

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