



Meter Trax™ Control



Meter Trax™ Control

Quality Control

Meter Trax™ is an ideal quality control for point-of-care meters. It is available in easy-to-use dropper bottles and has assigned values for many meters, such as Hemoglobin on the HemoCue® systems and Hematocrit on the i-STAT®. A great compliment to Meter Trax™, the Unity™ Interlaboratory Program is designed for peer group comparisons within your laboratory or with others.

Operator Competency

Meter Trax™ can be used as part of a laboratory program to assess operator competency with glucose meters. Its whole blood matrix makes it similar to a patient sample, and is ideally suited for use as an internal blind testing sample. Meter Trax™ can help assess competency in technique, adherence to procedure and interpretation.

Analytes

Glucose
Hematocrit (HCT)
Hemoglobin (Total)

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

Ordering Information

Cat #	Meter Trax™ Control	Quantity
970	Trilevel (2 of each level)	6 x 2 mL
971	Low Level	6 x 2 mL
972	Mid Level	6 x 2 mL
973	High Level	6 x 2 mL
970X	Trilevel MiniPak (1 of each level)	



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** 1-800-180-1224 **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** 7-495-721-14-04 **Singapore** 65-6415-3170 **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



Printed on recycled paper with soy-based inks

HemoCue is a trademark of Quest Diagnostics Company. i-STAT is a trademark of Abbott Laboratories.

© 2011 Bio-Rad Laboratories, Inc. Printed in the USA 02/11 QSD11-019 Q-1531