



Lyphocheck® Diabetes Control & Hemoglobin A1C Linearity Set

BIO-RAD

Lyphochek® Diabetes Control

A human whole blood based product designed to monitor the performance of hemoglobin tests used in diabetes monitoring, including Hemoglobins A1, A1C, F and Total Glycated Hemoglobin.

- Human whole blood based control
- 2 significant levels
- Convenient 0.5 mL fill size
- 3 year shelf life at 2–8°C
- 7 day open-vial stability at 2–8°C
- Extensive value assignment listings

Analytes

Hemoglobin A1
Hemoglobin A1C
Hemoglobin F
Hemoglobin, Total
Hemoglobin (Total Glycated)

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

Ordering Information

Cat #	Lyphochek® Diabetes Control	Quantity
740	Bilevel (3 of each level)	6 x 0.5 mL
740X	Bilevel MiniPak (1 of each level)	2 x 0.5 mL

Lyphochek® Hemoglobin A1C Linearity Set

A human whole blood based product designed to verify linearity throughout the patient reportable range of Hemoglobin A1C assays.

- Human whole blood based control
- 4 significant levels
- 3 year shelf life at 2–8°C
- 7 day open-vial stability at 2–8°C
- Linearity Web™ software tool is available online at www.QCNet.com

Assayed Values Available for Common Analyzers

Bio-Rad Analyzers

- D-10™
- in2it™
- Micromat II™
- DiaSTAT™
- VARIANT™
- VARIANT™ II/TURBO

Other Analyzers

- Beckman Coulter® SYNCHRON® Systems
- Abbott Architect™ Series
- Roche COBAS Integra® & Hitachi®
- Siemens ADVIA® 1650 and Dimension® Series
- TOSOH A1C 2.2 Plus, G7 and G8
- Primus® Diagnostics Series

Analyte

Hemoglobin A1C

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

Target Values for Hemoglobin A1C*

Level 1	2.5 - 4.0%
Level 2	5.0 - 7.0%
Level 3	9.0 - 12.5%
Level 4	17 - 22%

* Ranges reflect typical recovery values for this product. Exact values may vary depending on methodology. Absolute values may also be dependent on the accuracy of the instruments/ reagent system calibration, if applicable.

Ordering Information

Cat #	Hemoglobin A1C Linearity Set	Quantity
120	Set (1 of each level)	4 x 0.5 mL



These controls are 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

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