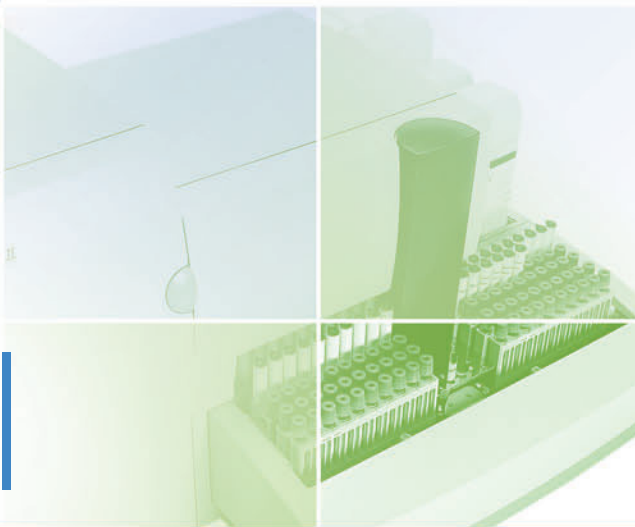




Bio-Rad Laboratories
The Clear Choice in Hemoglobin Testing



VARIANT™ II

HbA_{1c} Program

Hemoglobin A_{1c} Program for Diabetes Monitoring

The Bio-Rad VARIANT™ II Hemoglobin A_{1c} Program provides a fast, easy method for glycohemoglobin measurement that correlates to the Diabetes Control and Complications Trial (DCCT) reference study and is certified by the National Glycohemoglobin Standardization Program (NGSP).

Laboratories worldwide have come to depend on Bio-Rad for the reference method for glycohemoglobin analysis. Now, laboratories can realize the added benefits of a simplified procedure and total automation using the Bio-Rad VARIANT™ II Hemoglobin Testing System.



*VARIANT II
HbA_{1c} Program*



Efficient

- VARIANT™ II HbA_{1c} Program fits your workflow environment
- Sampling from primary whole blood tubes—no sample preparation
- Accepts pediatric tubes
- Clear, accurate results in minutes

Reliable

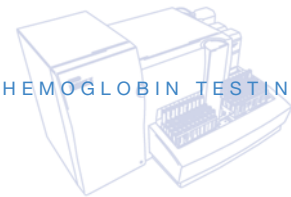
- Proven HPLC methodology ensures accuracy, minimizing repeats

Powerful

- Abnormal hemoglobins detected
- Labile A_{1c} chromatographically removed
- Positive sample identification
- LIS bi-directional interface
- Clinical Data Management (CDM™) Software
- Clinical Systems Network (CSN™) Ready – Online tools and manuals as well as quality control, instrument support and remote education

H E M O G L O B I N T E S T I N G





VARIANT™ II HbA_{1c} Program

Bio-Rad CDM System 201
Bio-Rad VARIANT II Instrument #10457-1

PATIENT REPORT
V2_A1c

Patient Data
Sample ID: Unknown-1-2230
Patient ID: 2230
Name: 84
Physician: Rack ID: 9
Sex: Tube Number: 05/14/2001 17:21:40
DOB: Report Generated: Operator ID: SW

Analysis Data
Analysis Performed: 05/14/2001 17:57:25
Injection Number: 2230
Run Number: 84
Rack ID: 9
Tube Number: 05/14/2001 17:21:40
Report Generated: Operator ID: SW

Comments:

Peak Name	Calibrated Area in	Area in	Retention Time (min)	Peak Area
A1b	—	1.2	0.16	31726
LA1c/CHb	—	2.2	0.26	59272
A1c	5.5*	—	0.89	125728
P3	—	6.1	1.68	162588
Ao	—	83.6	1.81	2225309
Total Area:				2667946

A1c Concentration = 5.5%*

Analysis comments:

VARIANT II HbA_{1c} results: Normal Patient Chromatogram

Bio-Rad CDM System 201
Bio-Rad VARIANT II Instrument #10457-1

PATIENT REPORT
V2_A1c

Patient Data
Sample ID: Unknown-1-2208
Patient ID: 2208
Name: 84
Physician: Rack ID: 0331
Sex: Tube Number: 7
DOB: Report Generated: Operator ID: SW

Analysis Data
Analysis Performed: 05/14/2001 16:45:47
Injection Number: 2208
Run Number: 84
Rack ID: 0331
Tube Number: 7
Report Generated: Operator ID: SW

Comments:

Peak Name	Calibrated Area in	Area in	Retention Time (min)	Peak Area
A1b	—	3.9	0.26	110773
LA1c/CHb	—	2.0	0.63	56673
A1c	10.7*	—	0.85	230383
Ao	—	86.0	1.81	2452540
Total Area:				28050374

A1c Concentration = 10.7%*

* Values outside of expected ranges

Analysis comments:

VARIANT II HbA_{1c} results: Diabetic Patient Chromatogram

Meets Industry Standards

The Bio-Rad VARIANT™ II HbA_{1c} Program is CE marked to meet the European IVD Directive and has been anchored to the IFCC Reference Method. The program is also certified by the NGSP as being traceable to the Diabetes Control and Complications Trial (DCCT).



IFCC Anchored



NGSP Certified



CE Marked

Ordering Information

Catalog No.	Description
270-2101	VARIANT™ II HbA _{1c} Reorder Pack 1000 tests
740	Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels) 6 x 0.5 mL
120	Lyphochek® Hemoglobin A _{1c} Linearity Set, (1 each of 4 levels) 4 x 0.5 mL
BC80	EQAS™ Hemoglobin Program (3 analytes) 12 x 0.5 mL



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com.

Clinical
Diagnostics Group

Website www.bio-rad.com U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3461-5202 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-815-39-51 Portugal 351-21-472-7700 Russia 7-095-721-14-00 Singapore 65-6272-9877 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 44-208-328-2000