



VARIANT™ II TURBO
Hemoglobin Testing System

**TO USE AS AN AID IN
THE DIAGNOSIS OF DIABETES**



BIO-RAD



VARIANT™ II TURBO

Hemoglobin Testing System

Rapid
for demanding workloads
Efficient
to save you time and labor
Accurate
for better patient care

With so many patients counting on your results each day, delivering the accuracy they need as quickly as possible is a top priority. We designed the VARIANT™ II TURBO Hemoglobin Testing System with your needs—and high standards—in mind.

The VARIANT™ II TURBO Hemoglobin Testing System is a fully automated ion-exchange HPLC system for mid-to high-volume laboratories. Confidently report HbA_{1c} results to aid physicians with diagnosing those who have diabetes, identifying those who may be at risk of developing the disease, or monitoring the progress of those living with diabetes.

Let Bio-Rad help you see the full picture and make a difference in the lives of your patients.

Bio-Rad CDM System
Bio-Rad VARIANT II TURBO Instrument

PATIENT REPORT
V2 TURBO_A1c_2.0

Patient Data

Sample ID: —
Patient ID: —
Name: —
Physician: —
Sex: —
DOB: —

Analysis Data

Analysis Performed: 04/27/2017 10:06:12
Injection Number: 36
Run Number: 3
Rack ID: 18
Tube Number: 8
Report Generated: 04/27/2017 10:06:12
Operator ID: —

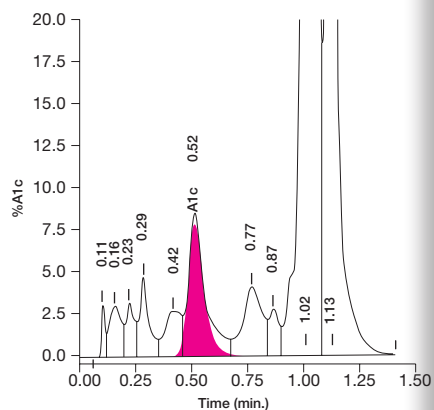
Comments:

Peak Name	IFCC mmol/mol	NGSP %	Area %	Retention Time (min)
Unknown	—	—	0.4	0.106
A1a	—	—	1.1	0.163
A1b	—	—	0.9	0.225
F	—	—	1.3	0.287
LA1c	—	—	1.0	0.424
A1c	62*	7.8*	—	0.519
P3	—	—	2.5	0.773
P4	—	—	0.9	0.869
Ao	—	—	49.3	1.018
Variant Window	—	—	38.7	1.126

*Values outside of expected range

Total Area

HbA1c (IFCC) = 62* mmol/mol HbA1c (NGSP) = 7.8*%



HbA1c reportable in the presence of hemoglobin S, C, D, and E traits*

Bio-Rad CDM System
Bio-Rad VARIANT II TURBO Instrument

PATIENT REPORT
V2 TURBO_A1c_2.0

Patient Data

Sample ID: Unknown-1-4513
Patient ID: —
Name: —
Physician: —
Sex: —
DOB: —

Analysis Data

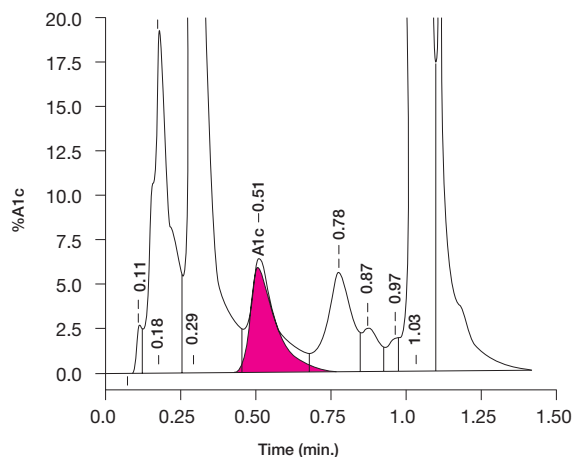
Analysis Performed: 02/14/2017 10:30:39
Injection Number: 4513
Run Number: 101
Rack ID: 0011
Tube Number: 1
Report Generated: 02/14/2017 10:44:15
Operator ID: —

Comments:

Peak Name	NGSP%	Area %	Retention Time (min)	Peak Area
Unknown	—	0.4	0.112	9414
A1a	—	8.4	0.176	218589
F	—	23.9	0.292	619799
A1c	5.9	—	0.510	98995
P3	—	3.6	0.777	93134
P4	—	1.0	0.875	26310
Unknown	—	0.5	0.967	13930
Ao	—	58.3	1.035	1507942

Total Area: 2,588,114

HbA1c (NGSP) = 5.9%



HbA1c reportable in the presence of HbF up to 25%*

* See VARIANT™ II TURBO HbA1c Kit - 2.0 Instructions for Use (Catalog # 12000447)

Report both IFCC (mmol/mol) and NGSP (%) HbA1c values

Every Patient Deserves **Quality Results**



Bring speed, efficiency, and accuracy together for quality patient care

You never forget the people who count on your results – even on your busiest days. Whether physicians are ordering tests to aid in the diagnosis of diabetes, identifying patients who may be at risk for developing the disease, or monitoring their patients who are already living with diabetes, the VARIANT™ II TURBO Hemoglobin Testing System with the VARIANT™ II TURBO HbA_{1c} Kit – 2.0 is the method of choice for laboratories demanding the best for their patients. With the ability to process thousands of tests per month, this robust system lets you be the difference for the people who count on you.

Rapid

- 97 seconds per sample

Efficient

- Fully automated operation saves time and labor
- Eliminate time-consuming sample prep with primary cap-piercing tube sampling
- Flexibility in workflow through either continuous or batch sampling
- Minimize hands-on time with cartridge lifetimes of 2,500 tests
- Maintain consistent, precise HbA_{1c} results with a single calibration per cartridge
- Hemoglobin Capillary Collection System allows remote finger-stick sample collection to support outreach programs to reach more patients
- Easy-to-use tools for QC, instrument support, and education

Accurate

- No interference in the presence of:
 - Hemoglobin S, C, D, or E traits*
 - Hemoglobin F (HbF) concentrations up to 25%*
 - Carbamylated (CHb) and Labile A_{1c} (LA1c)*
 - Acetylated hemoglobins*
- Anchored to the IFCC reference method
- NGSP certified and traceable to Diabetes Control and Complications Trial (DCCT)

* See VARIANT™ II TURBO HbA_{1c} Kit - 2.0 Instructions for Use (Catalog # 12000447)

*Every Patient Deserves Quality Results. You Deserve **Efficiency and Accuracy.***



Powerful Data Management

- Results and chromatograms captured and archived by Bio-Rad's Clinical Data Management (CDM™) software
- CDM interfaces with your laboratory information system for optimal data management
- Advanced diagnostics, user alerts, and prompts reduce costs and training time

VARIANT™ II TURBO Hemoglobin Testing System



Unparalleled Support

Bio-Rad's world-class service and support team is backed by our decades-long leadership in the field of HbA_{1c} testing. The VARIANT™ II TURBO Hemoglobin Testing System represents our continued commitment to helping you be the difference for the patients who count on you.

Ordering Information

Catalog #	Description	
270-2600	VARIANT™ II TURBO Hemoglobin Testing System	1 unit
270-2601	VARIANT™ II TURBO Hemoglobin Testing System without computer	1 unit
12000447	VARIANT™ II TURBO HbA _{1c} Kit – 2.0	2500 tests
270-2730	VARIANT™ II TURBO Hemoglobin A _{1c} Wash/Diluent Set	4 x 2500 mL
740	Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels)	6 x 0.5 mL
12000070	Lyphochek® Hemoglobin A _{1c} Linearity Set (1 each of 6 levels)	6 x 0.5 mL
BC80	EQAS Hemoglobin Program (3 analytes)	12 x 0.5 mL
171	Liquichek™ Diabetes Control, Level 1	6 x 1 mL
172	Liquichek™ Diabetes Control, Level 2	6 x 1 mL
173	Liquichek™ Diabetes Control, Level 3	6 x 1 mL
196-2051	Hemoglobin Capillary Collection System (HCCS)	5 tests
196-2052	Hemoglobin Capillary Collection System (HCCS)	100 tests
196-2053	Hemoglobin Capillary Collection System (HCCS)	5000 tests
196-2100	Plastic Capillaries, 100 per dispenser	20 dispensers
196-1054	Capillary Holder	20 holders

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/diagnostics

*Clinical
Diagnostics Group*

Website www.bio-rad.com/diagnostics **U.S.** 1-800-224-6723 **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** 86-21-61698500 **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