

VARIANT™ II TURBO Hemoglobin Testing System for HbA_{1c}





VARIANT™ II TURBO Hemoglobin Testing System

Faster
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 System, with your needs—and high standards—in mind.

The VARIANT™II TURBO System is built for exceptional performance on even your busiest day. This state-of-the-art, fully-automated system combines Bio-Rad's HPLC precision and variant detection with the rapid throughput you need. Count on the VARIANT™II TURBO System to deliver quality results at high test volumes—with unprecedented speed, enabling you to be the difference for the many people counting on your results every day.



Bring efficiency and accuracy together for quality patient care

You never forget the people who count on your results—even on your busiest days. We kept that commitment in mind when we developed the VARIANT™II TURBO Hemoglobin Testing System. With the ability to process thousands of tests per month, this cutting-edge system lets you be the difference for the people who count on you—in a fraction of the time.

Faster

- 97 seconds per sample
- Deliver accurate results in half the time

Efficient

- Fully-automated operation saves time and labor
- Eliminate time-consuming sample prep with primary tube sampling
- Flexibility in workflow through either continuous or batch sampling
- STAT capability
- Hemoglobin Capillary Collection System (HCCS) allows remote fingerstick sample collection to support outreach programs
- Simplified data management with powerful software
- Reagents supplied in convenient kit format
- Readily viewable reagent levels in transparent buffer bottles
- Easy-to-use tools for QC, instrument support and education

*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

Patient Data

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

Analysis Data

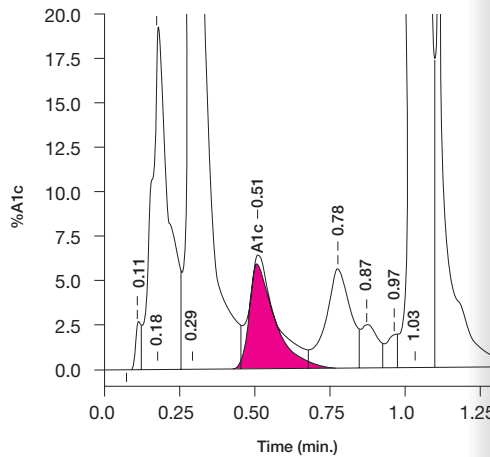
Analysis Performed: 02/14/2013 10:30:39
Injection Number: 4513
Run Number: 101
Rack ID: 001
Tube Number: 1
Report Generated: 02/14/2013 10:30:06
Operator ID:

Comments:

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

Total

HbA_{1c} (NGSP) = 5.9%



HbA_{1c} reportable in the presence of HbF up to 25%

Patient Data

Sample ID: Unknown-1-4516
Patient ID:
Name:
Physician:
Sex:
DOB:

Analysis Data

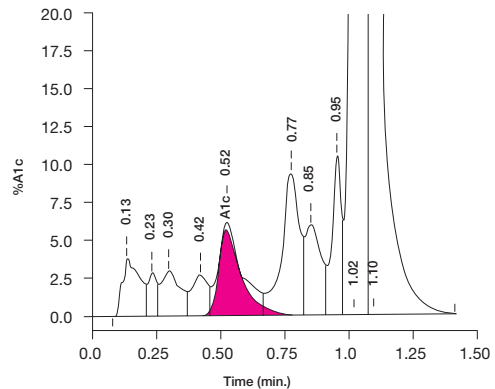
Analysis Performed: 02/14/2013 10:35:29
Injection Number: 4516
Run Number: 101
Rack ID: 0011
Tube Number: 4
Report Generated: 02/14/2013 10:39:06
Operator ID:

Comments:

Peak Name	NGSP%	Area %	Retention Time (min)	Peak Area
A1a	—	1.7	0.131	46011
A1b	—	0.6	0.229	17122
F	—	1.4	0.296	38756
LA1c	—	1.1	0.421	30277
A1c	5.6	—	0.524	90459
P3	—	4.4	0.774	117689
P4	—	2.4	0.853	64633
Unknown	—	2.5	0.953	67802
Ao	—	53.7	1.019	1441240
Variant Window	—	28.6	1.097	768180

Total Area: 2,682,168

HbA_{1c} (NGSP) = 5.6%



HbA_{1c} reportability in the presence of hemoglobin S, D and E traits*

Superior Accuracy

- Accurate and reportable HbA_{1c} results in the presence of hemoglobin S, C, D or E traits*
- Reportable HbA_{1c} results in the presence of carbamylated (CHb) and labile A1c (LA_{1c})*
- Excellent HPLC precision
- NGSP Certified and traceable to Diabetes Control and Complications Trial (DCCT)
- Anchored to the IFCC reference method
- Accurate HbA_{1c} in the presence of HbF up to 25%
- Accurate and reportable HbA_{1c} results in the presence of beta-thalassemia trait, as indicated by increased HbA₂ concentrations up to 10%.

* See VARIANT™ II TURBO HbA_{1c} Kit – 2.0 Instructions for Use.

Every patient deserves quality results. You deserve efficiency and accuracy.

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 System represents our continued commitment to helping you be the difference for the patients who count on you.



Efficiency and Accuracy

VARIANT™ II TURBO System

Rapid, fully-automated HPLC precision, and powerful software deliver quality results in half the time.

Ordering Information

Catalog No.	Description	
270-2600	VARIANT™ II TURBO Hemoglobin Testing System	1 unit
270-2601	VARIANT™ II TURBO Hemoglobin Testing System without computer	1 unit
270-2455	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
270-2100	VARIANT™ II TURBO Upgrade Kit	1 unit
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
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

Bio-Rad A_{1c} • Be the difference



**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** 32-9-385-5511 **Brazil** 5521-3237-9400
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** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200
Sweden 46-8-555-127-00 **Switzerland** +41 (0)26-674-52-01/02/03 **Taiwan** 886-2-2578-7189 **Thailand** 662-651-8311 **United Kingdom** +44 (0)20-8328-2000