

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





VARIANT™ II Hemoglobin Testing System

Superior
HbA_{1c} Accuracy

Exceptional
Ease of Use

Powerful
Data Management

Quality patient care is at the heart of what you do. And the VARIANT™ II system enables you to be the difference for the patients who count on you most—delivering high quality chemistry and providing accurate results.

The VARIANT II hemoglobin testing system is flexible, efficient and uses gold-standard, high performance liquid chromatography (HPLC) to deliver precise HbA_{1c} values in the presence of hemoglobin variants and other interfering substances.



*Your patients count on you.
You can count on the
VARIANT™ II System.*

Exceptional Ease of Use

- Fully automated operation saves time and labor
- Built-in quality control management simplifies quality assurance compliance
- Reagents supplied in convenient kit format
- Readily viewable reagent levels in transparent buffer bottles
- Primary tube sampling with cap piercing eliminates manual sample handling
- Positive sample tracking eliminates sample misidentification
- STAT capability
- Batch sampling and continuous feeding offer flexibility in workflow
- Easy-to-use online tools for QC, instrument support, remote education and communication services
- Hemoglobin Capillary Collection System (HCCS) allows remote fingerstick sample collection to support outreach programs

*Every patient deserves quality results. You deserve **ease of use.***

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



Bio-Rad CDM System
CDM 5.1 VII Instrument

Patient Report
V2_A1c_NU

Patient Data

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

Analysis Data

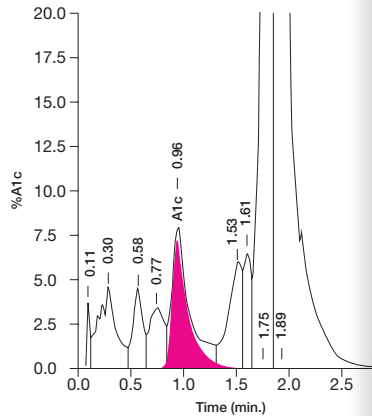
Analysis Performed: 06/21/2010 13:44:00
Injection Number:
Run Number:
Rack ID:
Tube Number:
Report Generated:
Operator ID:

Comments:

Peak Name	NGSP %	Area %	Retention Time
A1a	—	0.3	0.11
A1b	—	1.8	0.30
F	—	1.0	0.58
LA1c	—	1.2	0.77
A1c	7.2*	—	0.96
P3	—	1.8	1.53
P4	—	1.1	1.61
A0	—	47.1	1.75
E,D	—	43.0	1.89

*Values outside of expected range

HbA_{1c} (NGSP) = 7.2*%



HbA_{1c} reportable in patient sample with HbD trait

Bio-Rad CDM System
CDM 5.1 VII Instrument

Patient Report
V2_A1c_NU

Patient Data

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

Analysis Data

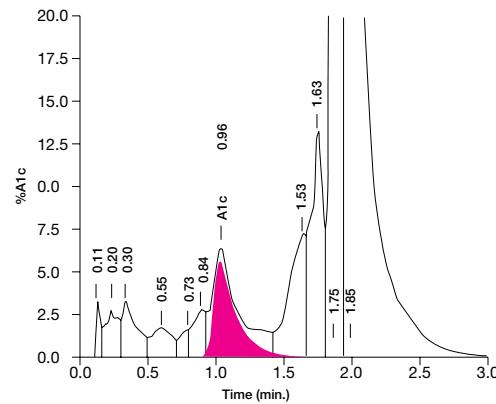
Analysis Performed: 06/21/2010 13:51:23
Injection Number: 207
Run Number: 24
Rack ID:
Tube Number: 5
Report Generated: 04/08/2011 12:22:29
Operator ID: YG-IFCC_DAY1

Comments:

Peak Name	IFCC mmol/mol	NGSP %	Area %	Retention Tim (min)	Peak Area
Unknown	—	—	0.3	0.11	6852
A1a	—	—	0.7	0.20	17722
A1b	—	—	0.9	0.30	23288
F	—	—	0.6	0.55	17159
LA1c	—	—	0.3	0.73	7405
Unknown	—	—	0.8	0.84	22534
A1c	37	5.6	—	0.96	62642
P3	—	—	2.5	1.53	66287
P4	—	—	2.7	1.63	70902
A0	—	—	53.4	1.75	1428051
E,D	—	—	35.6	1.85	950966

HbA_{1c} (IFCC) = 37 mmol/mol

Total Area: 2,673,808



HbA_{1c} reportable in patient sample with HbE trait

HbA_{1c} reportable in patient sample with CHb and LA_{1c}
(CHb is identified as the unknown peak between A1c and LA_{1c})

Superior HbA_{1c} Accuracy

- HbA_{1c} accuracy in the presence of hemoglobins D, E, S or C
- HbA_{1c} accuracy in the presence of carbamylated hemoglobin (CHb) and labile A_{1c} (LA_{1c})
- Anchored to the IFCC Reference Method
- NGSP certified and traceable to the Diabetes Control and Complications Trial (DCCT)

With both IFCC (mmol/mol)
and NGSP (%) HbA_{1c} Values

quality results

VARIANT™ II 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 system represents our continued commitment to helping you be the difference for the patients who count on you.

**Uncompromising Performance**

VARIANT™ II system combines superior HbA_{1c} results, exceptional ease of use and powerful data management in one powerful automated solution offering HbA_{1c}, β -thalassemia and dual assays*.

*Not available in the US.

Ordering Information

Catalog No.	Description	
270-2000	VARIANT™ II Hemoglobin Testing System	1 unit
270-2101NU	VARIANT™ II Hemoglobin A1c Program Reorder Pack	1000 tests
740	Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels)	6 x 0.5 mL
120	Lyphochek® Hemoglobin A1c Linearity Set, (1 each of 4 levels)	4 x 0.5 mL
BC80	EQAS™ Hemoglobin Program (3 analytes)	12 x 0.5 mL
196-2052	Hemoglobin Capillary Collection System (HCCS)	100 tests

Bio-Rad A1c • See 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 (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