



VARIANT™ II TURBO Link

Fully-Automated HbA_{1c} Testing





VARIANT™ II TURBO Link

**Efficient
streamlined process**

**Powerful
integrated system**

**Automated
saves time and
resources**

You work hard every day to provide accurate, reliable results in the most efficient way possible. That commitment inspired us to develop the VARIANT™ II TURBO Link—an advanced, fully-automated testing system designed to streamline operations while ensuring quality results.

The VARIANT™ II TURBO Link integrates our proven HbA_{1c} assay performance with the Sysmex* HST-N fully-automated hematology system to handle over 90% of whole blood EDTA samples while delivering maximum flexibility and accuracy. The result is a consolidated offering that lets you be the difference—for your patients and your lab.

Patient Data

Sample ID: A001899564
Patient ID:
Name:
Physician:
Sex:
DOB:
Comments:

Analysis Data

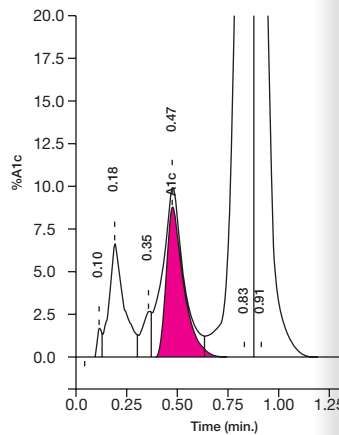
Analysis Performed:
Injection Number:
Run Number:
Rack ID:
Tube Number:
Report Generated:
Operator ID:

Peak Name	Calibrated Area %	Area %	Retention Time (min)
A1a	—	0.2	0.10
A1b	—	2.3	0.18
LA1c	—	0.6	0.35
A1c	8.8*	—	0.47
A0	—	54.7	0.82
S-Window	—	38.6	0.91

*Values outside of expected range

A1c Concentration: 8.8%*

Analysis Comments:



A1C accuracy in the presence of heterozygous hemoglobins S or C

Patient Data

Sample ID: M02
Patient ID:
Name:
Physician:
Sex:
DOB:
Comments:

Analysis Data

Analysis Performed: 06/28/2007 09:57:22
Injection Number: 2082
Run Number: 39
Rack ID: 009170
Tube Number: 7
Report Generated: 10/25/2007 10:00:12
Operator ID:

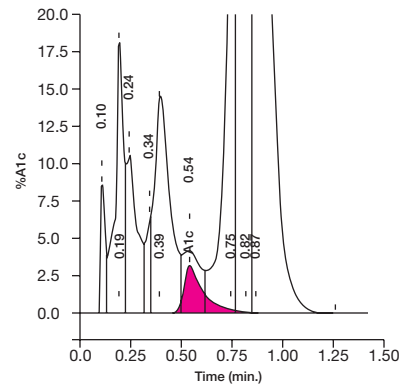
Peak Name	Calibrated Area %	Area %	Retention Time (min)	Peak Area
A1a	—	0.4	0.104	9055
A1b	—	1.7	0.191	37575
F	—	1.3	0.242	28267
LA1c	—	0.4	0.342	8661
CHb	—	2.7	0.390	60182
A1c	3.2*	—	0.537	15579
P3	—	3.5	0.745	79664
A0	—	48.7	0.818	1092831
Variant Window	—	40.7	0.867	912471

*Values outside of expected range

Total Area: 2244284

A1c Concentration: 3.2%*

Analysis Comments:



A1C accuracy in the presence of carbamylated hemoglobin (CHb) and labile A1C (LA_{1c})

Experience superior HbA_{1c} quality and accuracy

- Ion-Exchange methodology detects potential interferences
- Anchored to the IFCC reference method
- NSGP certified and traceable to the Diabetes Control and Complications Trial (DCCT)

quality results



Turn greater efficiency into greater patient care

With the VARIANT™II TURBO Link, we're combining Bio-Rad's decades of leadership in HbA_{1c} with Sysmex's proven expertise in automation to improve your laboratory's workflow. The result is a fully-automated solution that helps you be the difference with timely, accurate results.

Streamlined Operations

- On-demand testing 24/7
- Automated lavender top tube management
- Workflow optimization
- Reduced sample handling errors and delays

Save Time and Labor

- Sample ID decision logic on HST-N automation line
- Specimen retrieval and archiving using a sorter
- Standardized automated sample data decisions using MOLIS WAM™
- Automated reflex and repeat testing using MOLIS WAM™



*Every patient deserves quality results. You deserve **operational efficiency.***



More than just an HbA_{1c} Result

Bio-Rad A1c provides clinically significant information by detecting

- common hemoglobin variants
- potentially interfering substances

Fully-automated solution with Sysmex HST-N

- CBC
- Slide Smear/Prep
- HbA_{1c}
- Sorter
- Information Systems

VARIANT™ II TURBO Link

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



Powerful testing, simplified

Combining Performance and Efficiency

VARIANT™ II TURBO Link adds greater efficiency to the combination of benefits you've come to expect from our VARIANT™ II family, including precision, accuracy and real-time result delivery.

Ordering Information

Available through:

Sysmex America, Inc.
One Nelson C. White Parkway
Mundelein, IL 60060
1-866-7-SYSMEX (1-866-779-7639)

* Sysmex is a registered trademark of Sysmex Corporation.

Bio-Rad A1c • 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. 1-800-2BIO-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 36-1-455-8800 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)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 65-6415-3188 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