



VARIANT™ II TURBO Link System

Fully-Automated HbA_{1c} Testing





VARIANT™ II TURBO Link System

**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 System— an advanced, fully-automated testing system designed to streamline operations while ensuring quality results.

The VARIANT™ II TURBO Link System integrates our proven HbA_{1c} assay performance with the Sysmex* XN-9000™ Lavender Top Management® 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.



Turn greater efficiency into greater patient care

With the VARIANT™ II TURBO Link System, 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
- Sysmex Lavender Top Management®
- Workflow optimization
- Reduced sample handling errors and delays

Save Time and Labor

- Sample ID decision logic on XN-9000 automated hematology system
- Specimen retrieval and archiving using a sorter
- Standardized automated sample data decisions using Sysmex WAM™
- Automated reflex and repeat testing using Sysmex WAM™

More than just an HbA_{1c} Result

Bio-Rad A1c provides clinically significant information by detecting

- Common hemoglobin variants
- Potentially interfering substances

Fully-integrated Lavender Top Management® solution

- CBC testing
- Integrated slide maker/stainer
- HbA_{1c}
- Automated sample sorting and archiving
- Sysmex WAM™ decision logic rules database

Lavender Top Management is a registered trademark of Sysmex

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



Patient Data

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

Analysis Data

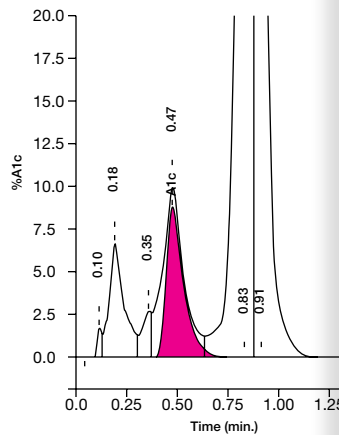
Analysis Performed:
Injection Number:
Run Number:
Rack ID:
Tube Number:
Report Generated:
Operator ID:

| 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:

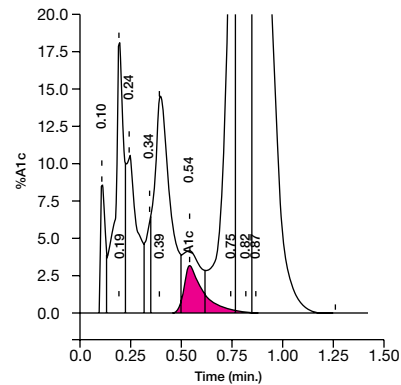
| Peak Name | Calibrated Area % | Area % | Retention Time (min) | Peak Area |
|----------------|-------------------|--------|----------------------|-----------|
| A1a | — | 0.4 | 0.104 | 9055 |
| A1b | — | 1.7 | 0.191 | 37575 |
| F | — | 1.3 | 0.242 | 28267 |
| LA1c | — | 0.4 | 0.342 | 8661 |
| CHb | — | 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

HEMOGLOBIN TESTING

VARIANT™ II TURBO Link 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 Link System 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 System 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.
577 Aptakistic Road
Lincolnshire, IL 60069
1-800-3-SYSMEX (1-800-379-7639)
www.sysmex.com/US

Lavender Top Management is a registered trademark of Sysmex

* Sysmex is a registered trademark of Sysmex Corporation.

Bio-Rad A1c • See the difference

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** 32-9-385-5511 **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