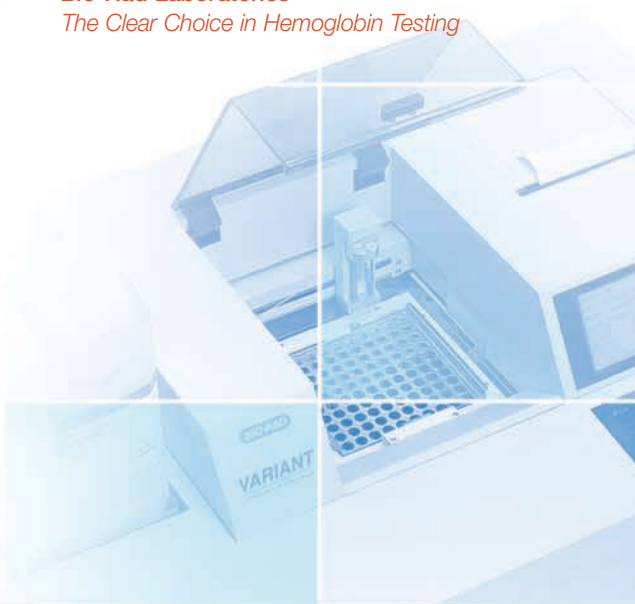




**Bio-Rad Laboratories**  
*The Clear Choice in Hemoglobin Testing*



# VARIANT™

## HbA<sub>1c</sub> Program

HEMOGLOBIN TESTING

### Hemoglobin A<sub>1c</sub> Program for Diabetes Monitoring

The VARIANT™ A<sub>1c</sub> program provides a fast, easy means to obtain gold standard HbA<sub>1c</sub> results. Laboratories worldwide have come to depend on Bio-Rad as the reference method for hemoglobin analysis. With the VARIANT™ A<sub>1c</sub> Program, laboratories can realize the added benefits of a simplified procedure and advanced automation.



### Efficient

- 3-minute analysis meets increasing throughput requirements
- 1-step sample prep improves turnaround time
- Enhanced data handling provides true walkaway operation.

### Reliable

- No interferences from Schiff base, temperature or Hemoglobins S and C

### Powerful

- Runs on a multi-analyte platform (HbA<sub>1c</sub>, β-thalassemia, Total GHb, α-thalassemia, Hemoglobinopathy and Sickle Cell Programs)
- Hemoglobins S and C detected when present in patient samples
- On-board QC program

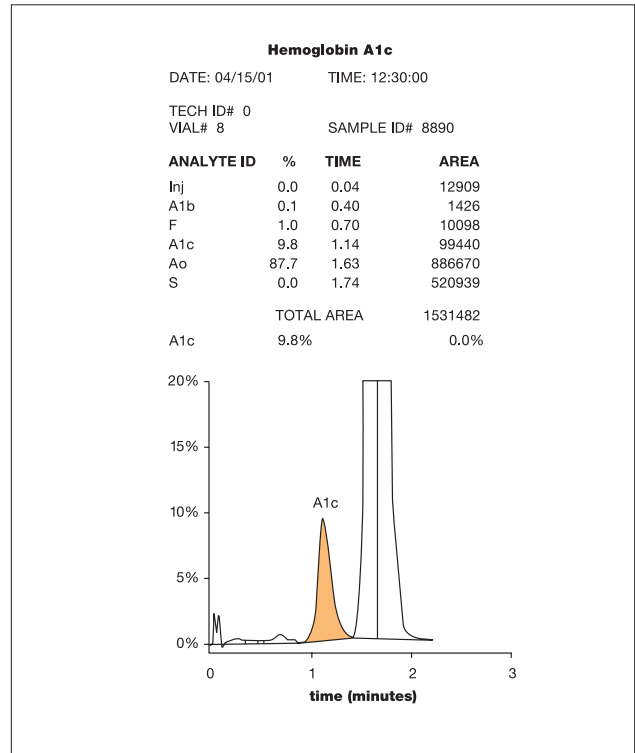




# VARIANT™ HbA<sub>1c</sub> Program



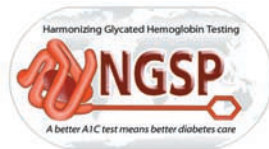
VARIANT HbA<sub>1c</sub> results: Normal patient chromatogram



VARIANT HbA<sub>1c</sub> results: Diabetic patient with sickle cell trait

## NGSP Certified

The Bio-Rad VARIANT II™ Total GHb Program has been certified by the NGSP as being traceable to the Diabetes Control and Complications Trial (DCCT) reference method using the new precision criteria of ≤ 4% CV. The mission of the NGSP Program is to harmonize glycosylated hemoglobin testing worldwide.



## Ordering Information

Catalog No.	Description
270-0003	VARIANT™ HbA <sub>1c</sub> Program Reorder Pack ..... 1000 tests
270-2167	Capillary Collection Kit* ..... 100 tests
740	Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels) ..... 6 x 0.5mL
120	Lyphochek® Hemoglobin A <sub>1c</sub> Linearity Set, (1 each of 4 levels) ..... 4 x 0.5 mL

\* Not available in the U.S.



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com](http://www.bio-rad.com).

Clinical Diagnostics Group

Website [www.bio-rad.com](http://www.bio-rad.com) U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-2507-6191 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-815-39-51 Portugal 351-21-472-7700 Russia 7-095-721-14-00 Singapore 65-6272-9877 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 44-208-328-2000