



Bio-Rad Laboratories
The Clear Choice in Hemoglobin Testing



VARIANT™

Sickle Cell Short Program

H E M O G L O B I N T E S T I N G

Precise, Automated Hemoglobin Screening for Newborns

The VARIANT™ Sickle Cell Short Program provides detection and presumptive identification of hemoglobins S, F, A, C, D and E in newborn dried blood spot specimens. Rapid analysis and unattended operation maximize productivity and efficiency.

- Increased productivity: The 3-minute specimen analysis delivers high throughput for your laboratory. Operation is performed unattended.
- Objectivity: Hemoglobins S, F, A, C, D and E are detected based on quantitative thresholds. Patient data can be easily stored off-line on the laboratory's information management system or VDMS 2000™ Software Program.
- Relative ratios: Relative ratios between different hemoglobins can help distinguish sickle cell disease (FS) from sickle cell trait (FAS) and sickle/ β -thalassemia (FSA).

Efficient

- 3-minute analysis
- Walk-away operation
- 1-step sample preparation

Reliable

- Reliable HPLC technology
- Superior retention time precision
- "Smart" reports highlight hemoglobins detected
- Not technique-dependent

Powerful

- Unmatched objectivity
- Runs on a multi-analyte platform (HbA_{1c}, Total GHb, α -thalassemia, β -thalassemia and Hemoglobinopathy Programs)
- Works with VDMS 2000™ software to manage results from up to eight instruments

BIO-RAD



VARIANT™ Sickle Cell Short Program

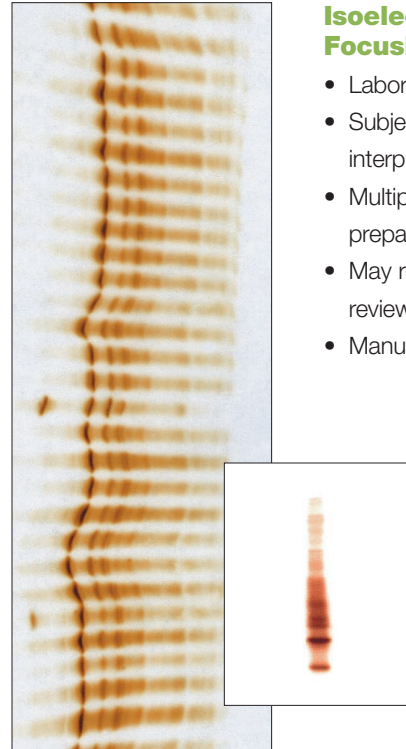
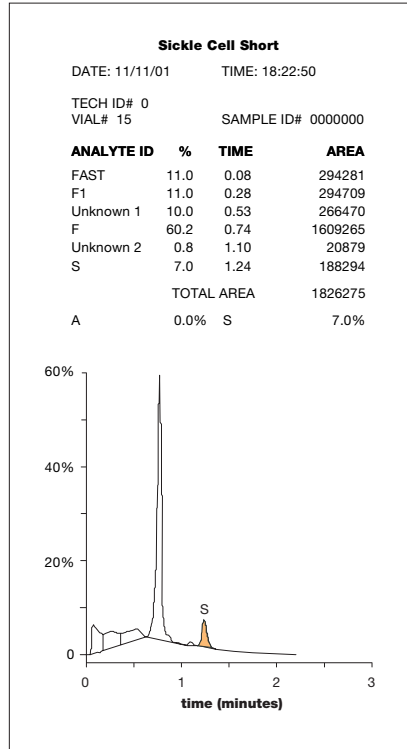
Which would you rather read?

Objective Interpretation

Subjective Interpretation

VARIANT™ Sickle Cell Short Program

- Automated processing
- Objective analysis
- 1-step sample preparation
- Printed record
- Electronic record transfer
- Better patient sample tracking



Isoelectric Focusing

- Labor intensive
- Subjective interpretation
- Multiple-step sample preparation
- May require multiple reviewers for one result
- Manual data entry

Ordering Information

Catalog No.	Description	Quantity
270-0001	VARIANT™ Hemoglobin Testing System (110V)	1 system
270-0021	VARIANT™ Hemoglobin Testing System (220V)	1 system
270-0500	VARIANT™ Sickle Cell Short Program Reorder Pack	1000 tests
270-0450	VDMS 2000™ System and Software	1 package
195-5174	Microplate Tray Insert (for microplate testing, goes with 195-5185)	1 unit
195-5185	Microplate Cover (for microplate testing, goes with 195-5174)	1 unit



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com.

Clinical
Diagnostics Group

Website 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