

Bio-Rad Laboratories
Results that improve lives.

VARIANT™

Sickle Cell Short Program

NEWBORN SCREENING

- Unequivocal chromatographic results
- Rapid, 3-minute analysis
- Walkaway operation
- "Smart" reports highlight detected hemoglobins, area percentages and other key sample data
- Proven, reliable HPLC technology
- Electronic data – interfaces with any lab information system

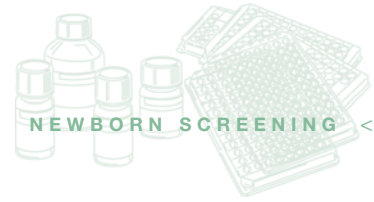
VARIANT Sickle Cell Short Program



Bio-Rad Newborn Screening Systems Approach

- Screening Assays
- DNA Confirmation Assays
- EIA Instrumentation
- HPLC Instrumentation
- Data Management Software

BIO-RAD



Sickle Cell Short Program

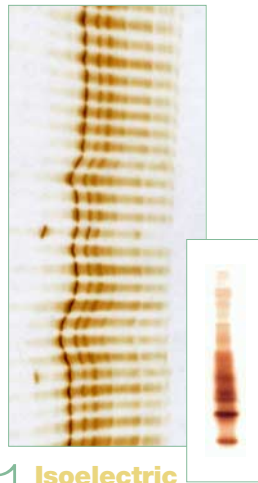
Clinical Background

Sickle Cell Anemia is a disease caused by a genetic mutation in the β -globin gene resulting in the production of hemoglobin S instead of hemoglobin A. In this condition, oxygen-carrying red blood cells quickly lose their ability to carry oxygen throughout the body. The cells tend to distort and elongate, which causes them to become rigid and assume a sickle shape. Over time, the sickle-shaped cells cause obstructions to blood microcirculation, resulting in tissue damage and other life threatening complications.

Product Highlights

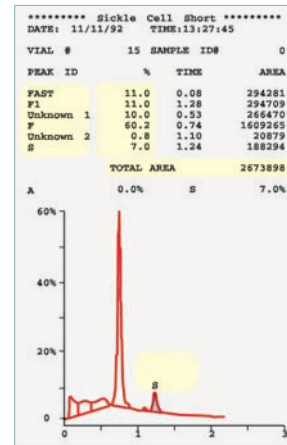
- Unmatched objectivity
- Fully automated
- No manual data entry
- Part of a full hemoglobin testing product line

Which would you rather read?



1 Isoelectric Focusing

- Labor intensive
- Subjective interpretation
- Multiple-step sample preparation
- Can require multiple reviewers for 1 result
- Manual data entry
- Hazardous chemicals



2 VARIANT Sickle Cell Short

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

Ordering Information

Catalog No.	Description	
270-0001	VARIANT™ Hemoglobin Testing System	110V
270-0021	VARIANT™ Hemoglobin Testing System	220V
270-0500	Sickle Cell Short Program Reorder Pack	1000 tests
270-0700	Alpha Thal Short Program Reagent Pack	100 tests
270-0713	Alpha Thal Short Program Cartridge/Prefilter	300 tests
270-0711	Alpha Thal Short Program Hardware Kit	1 unit
270-0450	VDMS™ 2000 System and Software complete with PC, Monitor, Printer, Rocket communication board, and necessary cables for 1 instrument connection	1 package
270-0451	VDMS™ 2000 Software	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 5521-3461-5202 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 48-22-331-99-99 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