



VARIANT™ II β -thalassemia Short Program

Fully-Automated Hemoglobin Testing

The VARIANT™ II β -thalassemia Short Program is a fully-automated method for β -thalassemia and abnormal hemoglobin testing.

Efficient

- Performs hemoglobin screening, HbA₂ quantitation and HbF quantitation in a single run
- Sampling from primary whole blood tubes—no sample preparation
- Clear, accurate results in minutes

Powerful

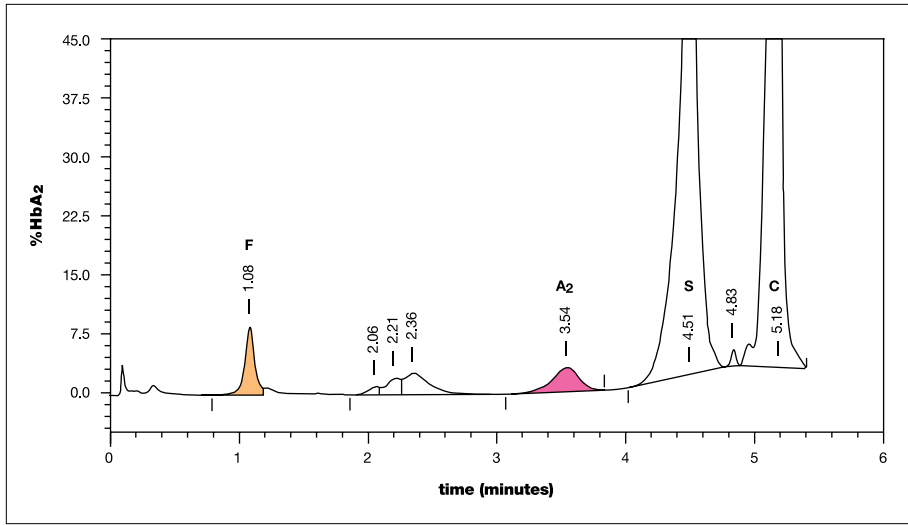
- Simultaneous quantitative results for HbA₂ and HbF
- Detection of hemoglobinopathies
- On-board QC program

Reliable

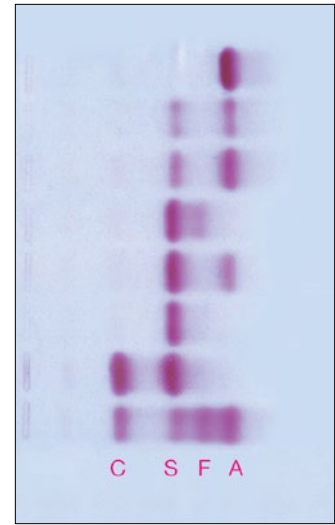
- Results are objective and operator independent
- Proven HPLC precision
- Full automation greatly reduces hands-on time
- Bar code reading of sample tubes eliminates the possibility of sample misidentification
- Results and chromatograms are stored electronically

BIO-RAD

VARIANT™ II β-thalassemia Short Program



VARIANT™ II Chromatogram



Traditional method

A typical β-thalassemia Short Program chromatogram from a patient with HbS/C phenotype. Integration of the chromatogram by the VARIANT™ II yields objective and precise results when compared to traditional electrophoresis methods.

Ordering Information

Catalog No.	Description	
270-2103	VARIANT™ II β-thalassemia Short Program.....	250 tests
270-2154	VARIANT™ II β-thalassemia Short Program.....	500 tests
553	Lyphocek® Hemoglobin A2 Control, Bilevel (2 each of 2 levels).....	4 x 0.5 mL
BC80	EQAS™ Hemoglobin Program (3 analytes).....	12 x 0.5 mL



**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