



Testing for Thalassemia and Sickle Cell

- Fully-automated HPLC methods
- Most accurate results, high reproducibility
- On-board QC program
- Worldwide references

The choice of experts for over 25 years

BIO-RAD

Thalassemia, Sickle Cell and Hemoglobinopathy Testing Systems

VARIANT™ II Hemoglobin Testing System

For full automation and breadth of menu, choose the VARIANT™ II Hemoglobin Testing System. This multi-analyte system offers a β -thalassemia Program, as well as a Dual Program with short and extended modes of operation requiring no change of column or buffers. The VARIANT™ II menu also includes HbA_{1c} for diabetes testing.



VARIANT™ nbs Newborn Screening System

The VARIANT™ nbs System is the next generation automated system for screening newborns for abnormal hemoglobins. With the VARIANT™ nbs System, dried blood spot specimens in microplates are analyzed for the detection and presumptive identification of hemoglobins S, F, A, C, D and E. The VARIANT™ nbs System delivers maximum efficiency through fully-automated analysis and advanced results reporting.



D-10™ Hemoglobin Testing System

For full automation, comprehensive menu and a small foot print, choose the D-10™ Hemoglobin Testing System. The D-10™ System offers a Dual Program for HbA_{1c} and β -thalassemia testing with short and extended modes of operation requiring no change of column or buffers. For added throughput, the D-10™ Rack Loader expands the D-10™ System sample capacity from 10 to 50 samples per run. With primary tube sampling, the D-10™ System is the easiest system available for thalassemia and diabetes testing.



Ordering Information

Catalog No.	Description	Unit
270-2000	VARIANT™ II Hemoglobin Testing System (110V/220V)	1 unit
250-3010	VARIANT™ nbs Newborn Screening System (110V/220V)	1 unit
250-3016	VARIANT™ nbs GDM Workstation Computer (110V/220V)	1 unit
220-0220	D-10™ Hemoglobin Testing System (110V/220V)	1 unit
220-0600	D-10™ Rack Loader (110V/220V)	1 unit

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. 1-800-2BIO-RAD (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 86-21-64260808 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 1-800-180-1224
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 41-61-717-95-55 Thailand 662-651-8311
United Kingdom +44-(0)20-8328-2000

Printed on recycled paper with soy-based inks

© 2011 Bio-Rad Laboratories, Inc. Printed in USA 10/11 DG11-1058 B-222