



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
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