

**Bio-Rad Laboratories**  
Results that improve lives.

# VDMS™ 2000

## System and Software

N  
E  
W  
B  
O  
R  
N  
  
S  
C  
R  
E  
E  
N  
I  
N  
G

### **Bio-Rad VDMS™ 2000**

#### *Customized Results for Your Hemoglobin Testing*

The VDMS™ 2000 is a flexible, turnkey solution for managing large amounts of chromatography data.

- User-defined rules determine hemoglobin phenotypes based on established laboratory criteria
- Presumptively identifies abnormal hemoglobin samples
- Manages data from up to 8 chromatographic stations
- Stores records electronically, reducing transcription errors
- Unlimited customized reporting options
- Open communication protocol permits data transmission to a variety of different lab information systems

### Bio-Rad Newborn Screening Systems Approach

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

**BIO-RAD**



# VDMS™ 2000

## Flexibility for Your Laboratory Needs

**VDMS** Convenient tabs simplify software navigation

VDMS Report Tools

**BIO-RAD**

New Worklists

Results by Worklist

Results by Punch Date

Repeated Samples

ID	Punch Date	Name	Status	1st Acc #	Last Acc #	State	Comments
3	11/9/99	Plate 2	Pass	425736	425827	Received	
4	11/9/99	Plate 3	Pass	425828	425919	Received	
5	11/9/99	Plate 4	Pass	425920	426011	Received	
6	11/9/99	Plate 5	Pass	426012	426050	Received	
17	11/10/99	Plate 1	Pass	426051	426142	Received	
18	11/10/99	Plate 2	Pass	426143	426234	Received	
19	11/10/99	Plate 3	Pass	426235	426326	Received	
20	11/10/99	Plate 4	Pass	426327	426418	Received	
30	11/10/99	Plate 7	Pass	426419	426510	Received	
31	11/10/99	Plate 8	Pass	426511	426602	Received	

Archive Restore Pass Fail

VARIANT 1 Ready  
VARIANT 2 Ready  
VARIANT 3 Ready  
VARIANT 4 Ready

Results for Worklist #4 (96 Samples) Filter: All Samples

Vial	ID	SampleType	Pattern	Final Result	HighArea	LowArea	Lo F	NoF	Confirm	Tot. Area	Dther	FAST	F1	F	A	E/A2	D
1	FAES	FAES Retention	CT-FAES	CT-FAES						1.44	1	2	8.5	57.4	19.6	5.5	0
2	FADC	FADC Retention	CT-FADC	CT-FADC						1.34	0	1.2	7.9	64	14.7	0.7	5.3
3	425828	Patient	FA	FA						2.87	3.3	6.7	17.1	59.9	13	0	0
4	425829	Patient	FA	FA						1.74	0.8	10.1	18	60.8	10.3	0	0
5	425830	Patient	FS							2.99	3.9	5.2	16	61.6	0	0	0
6			FS							3.27	3.1	6.4	19	62.2	0	0	0
7			FA	FA						2.45	4.3	6	16.9	59.5	13.3	0	0
8			FAS							2.44	3.2	6.7	20.8	63.7	3.2	0	0
9			FA	FA						2.2	4.2	4.8	15.8	57.5	17.7	0	0
10			FA	FA						2.78	3.1	5.5	18.2	64	9.2	0	0
11			FA	FA						3.81	3.1	6.2	19.2	62.9	8.6	0	0
12			FS							3.27	3.1	6.4	19	62.2	0	0	0
13	425838	Patient	FA	FA						1.97	4	6.2	19.2	61.4	9.2	0	0
14	425839	Patient	FA	FA						1.63	0.8	9.7	17.7	60.2	11.6	0	0

User defined rules determine hemoglobin phenotypes based upon your established laboratory criteria

Users can quickly find individual patient data by worklist or date through different views of the database

### Overview

The VDMS™ 2000 manages the routine tasks associated with hemoglobin screening of newborns. The VDMS™ 2000 automatically creates worklists and sends sample IDs to an instrument for analysis. The software then applies a set of predetermined rules to allow users to quickly identify abnormal results, flag samples for repeat testing, and upload final results to a lab information system.

### Product Highlights

- Customize results for your laboratory with unlimited user-defined “Decision Rules”
- Presumptively identify abnormal hemoglobin samples
- Collect data from multiple chromatographic stations

## Ordering Information

Catalog No.	Description
270-0450	VDMS™ 2000 System and Software complete with PC, monitor, printer, Rocket communication board, and necessary cables for 2 instrument connections ..... 1 package
270-0451	VDMS™ 2000 Software ..... 1 unit
270-0462	Cable – 10 foot ..... 1 unit
270-0463	Cable – 25 foot ..... 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](http://www.bio-rad.com).

Clinical Diagnostics Group

Website [www.bio-rad.com](http://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 47-23-381-41-30 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