

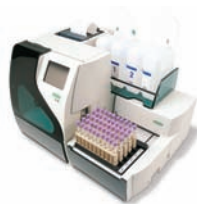


Bio-Rad provides low, mid and high-volume Hemoglobin A_{1c} testing solutions for ALL types of laboratories. Bio-Rad's platforms offer HPLC (High-Performance Liquid Chromatography), which is the Gold Standard in Hemoglobin A_{1c} testing.

High-Volume HbA_{1c} testing:

D-10™ Hemoglobin Testing System and D-10™ Rack Loader

50 sample walkaway capacity, 3 minutes per sample, fully-automated, flexible HPLC: HbA_{1c} Program, HbA₂/F/A_{1c} Dual Program, NGSP certified, excellent precision, small footprint



VARIANT™ II

100 sample walkaway capacity, 3 minutes per sample; fully-automated, flexible HPLC: HbA_{1c} Program, affinity-based total GHb program, NGSP certified

- *β-thalassemia program also available*



VARIANT™ II TURBO

100 sample walkaway capacity, 1.6 minutes per sample; fully-automated, one system runs 800 samples per month or more, NGSP certified



Mid-Volume HbA_{1c} testing:

D-10™ Hemoglobin Testing System

10 sample walkaway capacity, 3 minutes per sample; fully-automated, flexible HPLC: HbA_{1c} Program, HbA₂/F/A_{1c} Dual Program, NGSP certified, excellent precision, small footprint



Low-Volume HbA_{1c} testing:

Micromat™ II

Point-of-care instrument, 510(k) approved for prescription home use and, accordingly, is CLIA waived, NGSP certified, calibrated to the DCCT, 10 µL fingerstick sample, immediate results—5 minutes



Also Available:

Quality Control:

Lyphochek® Diabetes Control



Linearity:

Lyphochek® Hemoglobin A_{1c} Linearity Set



To order or for more information, contact your local Bio-Rad representative, call us toll free at **1-800-2BIO-RAD**, or visit us on the web at www.diabetes.bio-rad.com



World's Most Accurate Diabetes Testing

For over 27 years, Bio-Rad has been at the forefront in developing the world's most accurate and precise diabetes testing products. Today, Bio-Rad offers a full range of innovative HbA_{1c} products for clinical laboratories, hospital laboratories, physicians' offices and clinics. All Bio-Rad products are calibrated to the DCCT (Diabetes Control and Complications Trial) reference, and standardized by the NGSP (National Glycohemoglobin Standardization Program).

Ordering Information

Catalog No. Description

D-10™/D-10™ Rackloader

220-0220	D-10™ Hemoglobin Testing System	1 unit
220-0600	D-10™ Rack Loader 110/220V with Accessories	1 unit
220-0101	D-10™ Hemoglobin A _{1c} Program Reorder Pack	400 tests
220-0201	D-10™ Dual Program Reorder Pack	400 A _{1c} or 200 A ₂ /F/A _{1c} tests

VARIANT™ II

270-2000	VARIANT™ II Hemoglobin Testing System	1 unit
270-2101	VARIANT™ II HbA _{1c} Program Reorder Pack	1000 tests
270-2102	VARIANT™ II Total GHb Program Reorder Pack	500 tests

VARIANT™ II TURBO

270-2600	VARIANT™ II TURBO Hemoglobin Testing System	1 unit
270-2100	VARIANT™ II TURBO Upgrade Kit	1 unit
270-2415	VARIANT™ II TURBO Hemoglobin A _{1c} Program Reagent Kit	2000 tests
270-2416	VARIANT™ II TURBO Hemoglobin A _{1c} Wash/Diluent Set	2000 tests
270-2417	VARIANT™ II TURBO Hemoglobin A _{1c} Program Reagent Kit	1600 tests

Micromat™ II

280-0016	CLIA Waived Micromat™ II HbA _{1c} Instrument with Power Supply	1 unit
280-0007	Micromat™ II Control Cartridge	1 unit
280-0008	CLIA Waived Micromat™ II HbA _{1c} Test Cartridges	12 tests

Related Products

270-2154	VARIANT™ II β-Thalassemia Short Program Reorder Pack	500 tests
270-2103	VARIANT™ II β-Thalassemia Short Program Reorder Pack	250 tests
740	Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels)	6 x 0.5 mL
740x	Lyphochek® Diabetes Control, Bilevel MiniPak (1 each of 2 levels)	2 x 0.5 mL
120	Lyphochek® Hemoglobin A _{1c} Linearity Set (4 levels)	4 x 0.5 mL
270-0259	VARIANT™ Library of Abnormal Hemoglobins	1 unit
195-4151	CSN™ Installation Package*	1 unit

* A connection to CSN and the internet is required for access to VOL™



**Bio-Rad
Laboratories**

*For further information, please contact your Bio-Rad representative.
Call us toll free at 1-800-2BIO-RAD or visit our website
at www.diabetes.bio-rad.com.*

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics **U.S.** 1-800-2BIO-RAD **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** 36-1-455-8800 **India** 91-124-4029300 **Israel** 972-3-9636050 **Italy** +39-02-216091
Japan 81-3-5811-6290 **Korea** 82-2-3473-4460 **Mexico** +52-(555)-200-0520 **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** 65-6415-3188 **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