

Bio-Rad Laboratories
The Clear Choice in Hemoglobin Testing.

DiaSTAT™

Hemoglobin A1C Analyzer

Hemoglobin A1C Analyzer and Model 60 Dispenser

Physicians often expect HbA_{1c} determinations during the patient's routine office visit. Responding quickly and cost effectively requires a special kind of analyzer—easy to operate, adaptable to changing workloads and capable of delivering the highest quality test results.

Bio-Rad's DiaSTAT™ Hemoglobin A1C Analyzer fully satisfies the needs of low and medium volume testing environments. Compact, portable and easy-to-use, the DiaSTAT™ is also an ideal choice for the physician's office laboratory.

- Compact: ideal for small and medium-size laboratories, clinics or laboratories in physician offices
- STAT capability
- 100 and 500 test reagent kits available
- Reports HbA_{1c} results in the presence of hemoglobins S and C
- NGSP certified



Easy-to-use

- Hemoglobin A1C results in minutes
- Patient reports printed automatically
- Accommodates capillary collection application
- Simple, touch screen operation

Automated

- Intuitive, icon-driven operation
- Minimum training and maintenance
- Complete reagent kits

Efficient

- Walkaway performance
- Convenient, 15 sample capacity
- Continuous feed mechanism

BIO-RAD



DiaSTAT™ HbA_{1c} Analyzer



*DiaSTAT
HbA_{1c} Analyzer
Reagent Kit*

DiaSTAT™ Model 60 Dispenser

The Bio-Rad DiaSTAT™ Model 60 Dispenser is a dilutor/dispenser dedicated to specimen preparation for the DiaSTAT™ analyzer.

- Reduces Schiff base in five minutes
- Convenient, 1-step specimen preparation



NGSP Certified

The Bio-Rad DiaSTAT™ HbA_{1c} Program has been certified by the NGSP as being traceable to the Diabetes Control and Complications Trial (DCCT) reference method using the new precision criteria of $\leq 4\%$ CV.

The mission of the NGSP Program is to harmonize glycated hemoglobin testing worldwide.



Ordering Information

Catalog No.	Description
210-0001	DiaSTAT™ HbA _{1c} Analyzer 1 unit
210-0002	HbA _{1c} Program Reorder Pack 100 tests
210-0065	HbA _{1c} Program Reorder Pack (includes Pump Cassette) 500 tests
210-0007	Pump Cassette 500 tests
210-0064	Capillary Tubes 1000
196-1020	Printer Paper 10 rolls
210-0063	Tool Kit 1 unit
210-0097	Customer Manual (includes training guide) 1 unit
210-0075	DiaSTAT™ Model 60 Dispenser 1 unit
194-4003	Pipette Tips 10 boxes x 96 tips
210-0062	Pipette Tips (long) 200 tips
740	Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels) 6 x 0.5 mL
120	Lyphochek® Hemoglobin A1C Linearity Set (1 each of 4 levels) 4 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. 1-800-2-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 86-21-64260808 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 47-23-38-41-30 Poland 48-22-3319999 Portugal 351-21-472-7700 Russia 7-095-721-14-00 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 44-208-328-2000