



Hemoglobin Capillary Collection System for HbA_{1c} Testing

Sample collection, preparation and transport in one convenient device!

Extended sample stability

- Sample collection and transfer from remote locations
- Laboratory testing up to 2 weeks after sample collection

Increased patient compliance, comfort and safety

- No need for venipuncture
- Plastic capillaries, reducing potential injuries

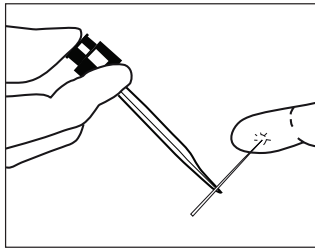
Improved workflow

- No sample preparation step
- Pierceable vials, ready for testing
- Immediate testing after sample collection
- Batch testing of capillary samples
- Multiple packaging sizes for every test volume
- For use with Bio-Rad's HPLC methods

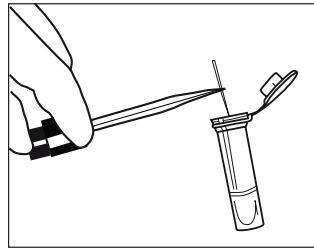


Hemoglobin Capillary Collection System for HbA_{1c} Testing

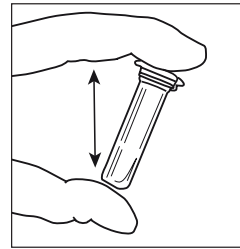
Sample Collection Procedure



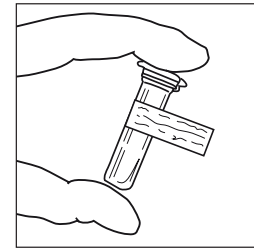
1. Collect blood sample using plastic capillary with the capillary holder.



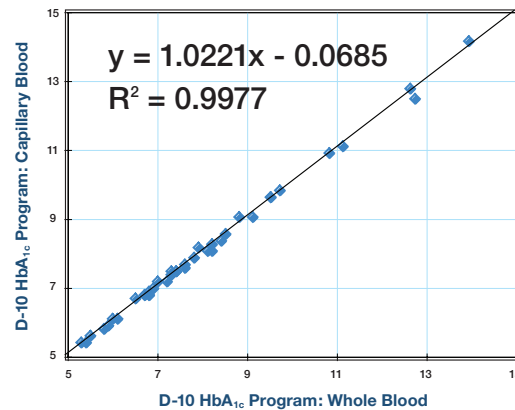
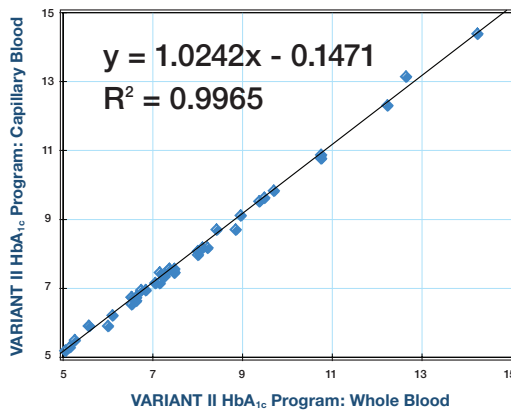
2. Place the filled capillary into the sample preparation vial.



3. Cap and shake the vial.



4. Label the vial and transport!



Reference: Study by Randie R. Little, Ph.D., University of Missouri School of Medicine

Ordering Information

Catalog No.	Description	
196-2050	Hemoglobin Capillary Collection System (HCCS).....	1 test
196-2051	Hemoglobin Capillary Collection System (HCCS).....	5 sets
196-2052	Hemoglobin Capillary Collection System (HCCS).....	100 tests
196-2053	Hemoglobin Capillary Collection System (HCCS).....	5000 tests
196-2100	Plastic Capillaries, 100 per dispenser.....	20 dispensers
196-1054	Capillary Holder.....	20 holders

Bio-Rad A1c • Be the difference



**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 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 91-124-4029300 Israel 972-3-9636050 Italy +39-02-216091
Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico 52(55)5200-0520 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