

Total GHb Program

GHb Affinity Method from the Leader in Diabetes Monitoring

The VARIANT II™ Hemoglobin Testing System offers a Total Glycated Hemoglobin (GHb)

answer on a fully-automated HPLC system designed specifically for the highest level of efficiency and quality of results. The Total GHb program will separate glycated hemoglobin from non-glycated hemoglobin regardless

of phenotype and without interference from carbamylated or acetylated hemoglobin.



Efficient

- Primary tube sampling
- No manual sample prep
- No primary tube stopper removal
- Fast results in minutes
- Continuous sample feeding

Reliable

- Accurate GHb results regardless of hemoglobin phenotype
- No interference
- No technique dependence
- NGSP certified

Powerful

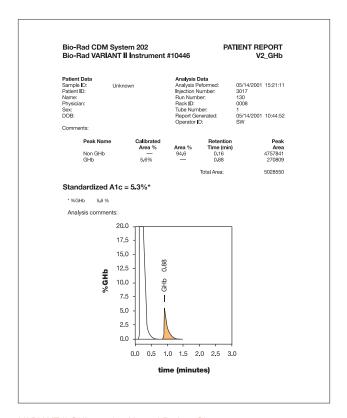
- Clinical Data Management™ (CDM) Software
- Real time, bi-directional interface with the laboratory information system
- CSN™ Ready Bio-Rad's Clinical Systems Network™
- On-board QC program
- Database archiving utility
- Positive sample identification

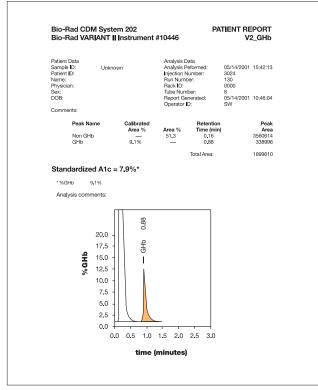


GHb Program



VARIANT II™ Total GHb Program





VARIANT II GHb results: Normal Patient Chromatogram

VARIANT II GHb results: Diabetic Patient Chromatogram

NGSP Certified

The Bio-Rad VARIANT II™ Total GHb 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 ≤ 4% CV. The mission of the NGSP Program is to harmonize glycated hemoglobin testing worldwide.

Ordering Information

| Catalog No. | Description |
|-------------|---|
| 270-2102 | VARIANT II™ Total GHb Reorder Pack500 tests |
| 740 | Lyphochek® Diabetes Control, Bilevel (3 each of 2 levels) |



Bio-Rad Laboratories For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com.