



## Quality and Confidence with Every Result

Bio-Rad offers superior confidence in reported patient results. The unparalleled lockout features of in2it™ provide administrator security and control ensuring testing is performed within established laboratory guidelines.

### in2it™ offers:

#### Operator lockout

Trained operators are critical in providing accurate results. Now you can release results with confidence knowing only trained, authorized operators have performed the test.

- Add up to 16 basic level and 16 administrator level operators.
- Operator identification provides traceability.

#### QC options

Patient care starts with controlling the performance of your testing system. Bio-Rad's in2it™ supports user specified levels of quality control testing. Now you can configure the analyzer for the desired level of QC control.

- QC prompt for simple reminders to run QC. Allows the operator flexibility in running various levels of QC or overriding the prompt.
- QC lockout requires QC testing. Patient testing cannot resume until the specified levels of QC material have been run.

Bio-Rad also provides superior results you can trust time after time. The cornerstones of Bio-Rad applications are consistent reproducibility and accuracy, providing you with reliable results for improved patient care.

### in2it™ results:

#### Reproducible

Laboratory precision produces consistent results with minimal variation.

	in2it™ 1st Result	in2it™ 2nd Result	in2it™ 3rd Result
N	5.1	5.0	5.1
D	7.1	7.1	7.0
HD	9.2	9.1	9.3

N: Normal D: Diabetic HD: High Diabetic

#### Accurate

Result comparison with NGSP reference values assigned using HPLC.

	NGSP % HbA1c	in2it™ Result % A1c
N	5.6	5.8
D	8.0	7.8
HD	10.9	10.8

N: Normal D: Diabetic HD: High Diabetic

For more information visit us on the web at [diabetes.bio-rad.com](http://diabetes.bio-rad.com)

