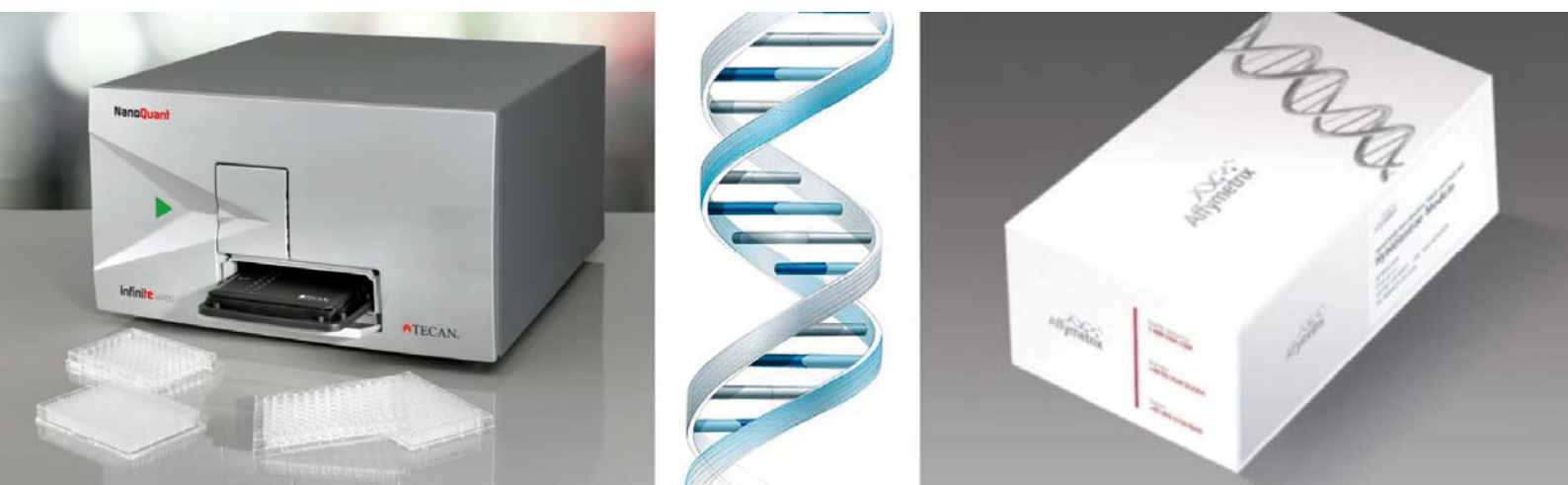




Low Volume DNA Quantification for Affymetrix® GeneChip

NanoQuant Plate™ for use on Infinite® 200 NanoQuant



Introduction

Scientists from Affymetrix® Inc., a world-leading supplier of microarray equipment and assays, evaluated the Tecan Infinite® 200 NanoQuant multimode microplate reader and two other spectrophotometers for DNA concentration analysis.

Detection, quantification and the assessment of purity of biomolecules in very small volumes are the basis of many modern assay systems.

The most used traditional method determining nucleic acid concentrations is based on measuring the absorbance at 260 nm (A_{260}). The purity of DNA or RNA is checked by comparing absorbance values from 260 and 280 nm measurements (260 / 280 ratio) and it is generally accepted that a sample of pure, double-stranded DNA (dsDNA) of 50 µg/ml has an absorbance at 260 nm of 1.0 and the ratio 260/280 nm will be greater than 1.8 [1].

The NanoQuant Plate™ from Tecan allows researchers to measure the concentration and purity of nucleic acids from up to 16 samples simultaneously.

Additionally, the efficiency of nucleic acid labeling with fluorescent dyes can be easily determined with this new tool. For these applications only 2 µl from the sample volume is needed, which makes the NanoQuant Plate an ideal tool for the latest applications in molecular biology.

In this note, data are presented from a comparison done by Affymetrix using Tecan's Infinite 200 NanoQuant, Molecular Devices SpectraMax® Plus and the Nanodrop® ND1000 instrument.

Material and Methods

NanoQuant Plate

- NanoQuant Plate with 16 (2 x 8) separate sample positions with specially designed quartz optics (Tecan Austria, Austria)



Figure 1: Tecan's NanoQuant Plate. The figure shows the NanoQuant Plate with closed lid as it will be read by the Infinite 200 readers, see also [2].

Each NanoQuant Plate is delivered with an optimized plate definition file (pdfx) for accurate positioning within the reader.

Instruments

- Infinite M200 NanoQuant Quad4 Monochromator detection system (Tecan Austria)
- SpectraMax Plus, Molecular Devices USA
- Nanodrop ND 1000, Thermo Scientific, NanoDrop Products, USA

Reagents

- Human genomic DNA was selected to use to quantify the concentration. DNA sample preparation and DNA quantification before and after whole genome amplification were done at Affymetrix.
- A Promega Human genomic DNA was selected to use to quantify the concentration.

Measurement Parameters and Settings

For DNA quantification the measurement parameters and instrument settings for Infinite 200 NanoQuant are shown below.

Measurements on Infinite 200 instruments	
Parameter	Setting
Absorbance	260 nm
Bandwidth	5 nm
Absorbance	280 nm
Bandwidth	5 nm
Settle time	0 ms
Number of reads	25

Table 1: Measurement parameters and settings for the Infinite 200 NanoQuant using Tecan’s i-control™ software.

For easy handling with Tecan’s i-control an application tab was introduced to select for example. ‘Quantify Nucleic Acid’ in the control bar. In Figure 2 the strip is shown as it appears in the workflow pane of the i-control software after blanking followed by the sample measurement.

For measuring with MD’s SpectraMax Plus a 100-fold dilution has to be prepared and 200 µl pipetted in a 96-well UV-transparent microtiter plate before DNA quantification.

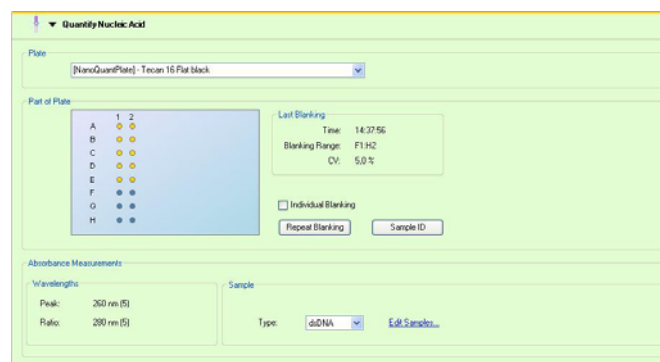


Figure 2: Quantify Nucleic Acid strip as it appears in the i-control software after blanking.

Samples were pipetted after blanking onto the NanoQuant Plate quartz positions. For sample application a multi channel pipette was used.

Two independent experiments were performed. In the first experiment genomic DNA from different origins and concentrations were used. In the second experiment a Promega Human genomic DNA was diluted to 1:1 dilution and 1:10 dilution. All samples were measured 3 to 6 independent times for each on Tecan Infinite 200 and the same DNA samples were then measured on the other two instruments in the lab ND1000 and SpectraMax Plus.

Sample	Tecan NanoQuant		MD SpectraMax Plus		NanoDrop ND1000	
	Average	CV [%]	Average	CV [%]	Average	CV [%]
Purified ProMale (n=6)	1080.3	1	1141.2	4	1009.4	0.8
Purified ProFemale (n=6)	999.8	0	1073	2	946.3	0.7
NA 19219 (n=6)	869.5	1	893.2	3	815.4	1.6
NA 19193 (n=6)	1107.5	1	893.2	4	1069.6	1.7
Average CV [%]		1		3		1.2

Table 2: Absorbance measurements of DNA samples using Tecan’s NanoQuant, MD’s SpectraMax Plus and the NanoDrop ND1000 instrument.

Sample	Tecan NanoQuant		MD SpectraMax Plus		NanoDrop ND1000	
	Average	CV [%]	Average	CV [%]	Average	CV [%]
ProMale (n=3)	40.3	5	37.0	3	39.9	2
ProFemale (n=3)	52.8	3	44.3	9	N/A	-
HapMapMale (n=3)	46.0	1	38.3	8	46.3	4
HapMapFemale (n=3)	46.3	2	36.7	3	46.0	2
Purified ProMale (n=3)	1054.3	1	1001.3	1	1110.6	5
Purified ProFemale (n=3)	974.8	1	964.0	0	995.0	1
Average [%]		2		4		3

Table 3: Absorbance measurements of PCR amplified DNA samples using Tecan’s NanoQuant, MD’s SpectraMax Plus and the NanoDrop ND1000 instrument.

Results

Measurement results for blanking and referencing, of nucleic acid quantification, purity check and results for determination of the labeling efficiency are displayed in Excel® automatically for all 16 measuring positions.

The results from experiment 1 are displayed in table 2 and 3 and the data derived from experiment 2 are shown in table 4.

The yields of the measurements are comparable between the three instruments, but the Tecan Infinite 200 NanoQuant exhibits an exceptionally high reproducibility (<3%).

DNA – Quantification

DNA samples ranging from about 40 to over 1000 ng/µl were measured using the NanoQuant Plate and for comparison with a MD’s SpectraMax Plus and the NanoDrop ND1000 instrument.

Sample (n=3)	Tecan NanoQuant			MD SpectraMax Plus			NanoDrop ND1000		
	ng/µl	ratio	CV [%]	ng/µl	ratio	CV [%]	ng/µl	ratio	CV [%]
PromegaFemaleDNA	227.6	1.8	1.0	217.9	1.93	4	215.7	1.86	2.2
PromegaFemaleDNA 1:1 dilution	112.6	1.8	0.1	118.7	1.98	2	106.9	1.88	0.7
PromegaFemaleDNA 1:10 dilution	22.1	1.8	2.3	25.3	1.96	5	20.6	1.99	3.7

Table 4: Experiment 2: Absorbance measurements of DNA samples using Tecan’s NanoQuant, MD’s SpectraMax Plus and the NanoDrop ND1000 instrument.

Conclusion

In many modern molecular biology applications sample size has become increasingly smaller and the determination of nucleic acid concentrations in these small volumes has become an even more demanding question in modern molecular biology. In forensic analysis DNA probes for genetic fingerprinting are sometimes only from a small number of cells.

With the NanoQuant Plate, Tecan provides a new tool for reproducible and sensitive quantification and purity check of nucleic acids.

The unique mathematic algorithm referencing every measurement position used for calculations makes the NanoQuant Plate an ideal tool to get reliable and highly accurate results.

For nucleic acid quantification and purity check, the Infinite 200 NanoQuant system is now listed by Affymetrix as a recommended instrument.

Literature

- [1] Maniatis T, Fritsch EF, Sambrook J (1982) Molecular Cloning. A Laboratory Manual. Cold Spring Harbor Laboratory, Cold Springs Harbor, NY
- [2] Quick Guide NanoQuant Plate™, 2008

User Statement

Vicky Huynh from the Product Development Organization at Affymetrix states:

“The Infinite M200 NanoQuant is one of my favorite quantification instruments in the lab. The software has many nice user-friendly interfaces. The data is automatically calculated and displayed in an Excel sheet after measurement. The measurement is quick and allows the measurement of 16 samples at a time in a minute or so. Also, there is no need to dilute samples, which saves time, diluting errors and plastic ware. Thus, to measure a plate, it would take less than 15 minutes. The instrument shows high reproducibility data <5%. With all these reasons, I would highly recommend the Infinite M200 NanoQuant as a quantification instrument. “

Acknowledgements

Tecan would like to thank Vicky Huynh from Product Development Organization at Affymetrix, in Santa Clara, California, USA for performing the experiments, the validation and providing the data. We would also like to thank Jim Collins, Director of Assay Development Molecular Diagnostic at Affymetrix for their constant help, the valuable comments and the excellent cooperation.

April, 2009

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.
© 2009, Tecan Trading AG, Switzerland, all rights reserved. Tecan is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland

Infinite is a registered trademark of Tecan Group Ltd. Männedorf, Switzerland.

NanoQuant Plate is trademark of Tecan Group Ltd. Männedorf, Switzerland.

Affymetrix is a registered trademark of Affymetrix, Inc. USA, SpectraMax Plus is a registered trademark of Molecular Devices USA, Nanodrop ND 1000 is a registered trademark of Thermo Scientific, NanoDrop products, USA

Austria +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 10 5869 5936 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
Italy +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 174 **Portugal** +351 21 000 82 16 **Singapore** +65 644 41 886
Spain +34 93 490 01 74 **Sweden** +46 31 75 44 000 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200
Other countries +41 44 922 8125