A fluorescence-based method for accurate molar quantification

of NGS libraries in minutes.

Ashesh A. Saraiya¹, Chaitali Parikh¹, Bin Li¹, and Joe Don Heath¹ ¹ Tecan Genomics, Inc., Redwood City, CA

INTRODUCTION

Accurate quantification of NGS libraries is critical for a successful sequencing run. Currently used methods of quantification are time-consuming, costly, and can be highly variable. We have developed NuQuant[®], a novel method to accurately quantify NGS libraries, that can be performed with a simple fluorescent measurement. NuQuant is compatible with the common red fluorescence excitation/emission filter set (650/670 nm), making it compatible with a wide range of bench top fluorometers and fluorescent plate readers. We have developed a custom library quantification application for the Qubit fluorometer that directly provides the molar concentration of a library. Utilizing this application, we have demonstrated that NuQuant has excellent reproducibility across users from multiple sites. We now demonstrate the compatibility of NuQuant with standard fluorescence plate readers, enabling quantification of libraries in a high-throughput fashion. We have tested NuQuant on a variety of commonly used plate readers such as the Tecan Infinite[®] 200 Pro and the Promega GloMax. Libraries in a 96-well format can be measured in a matter of minutes, without the need for sample dilution. Molar concentration of libraries was easily determined by utilizing a standard curve. We tested libraries with various input from 10 ng to 500 ng and insert size from 200 bp to 500 bp, and found good agreement between NuQuant values and a qPCR based quantification method. Most importantly, we observe good correlation between NuQuant library concentration and total number of sequenced reads. In conclusion, scientists with access to commonly used fluorescent plate readers can now use NuQuant to achieve rapid and cost-effective quantification of NGS libraries, generating highly uniform sequence reads in multiplex runs.

High-throughput library quantification with NuQuant



Standard Standard





Figure 1: Library quantification workflow with NuQuant. After the final bead purification step of the Celero[™] DNA-Seq or Universal Plus mRNA-Seq with NuQuant library preparation kits, the libraries and standards are diluted in the NuQuant buffer. The standards and libraries are assayed on fluorometers such as Qubit or other compatible fluorescence plate readers. The molar concentration of each library is directly measured for use in pooling. The pooled libraries are ready for sequencing on Illumina NGS instruments.



re			
ence	Day	Day	Day
	1	4	7

Figure 4: Experimental design. Libraries generated with the Celero DNA-Seq kit were aliquoted into a 96well black full-skirted PCR plate. The NuQuant standard was used to generate a standard curve and test for edge effects. The plate was sealed and NuQuant fluorescence was measured using the Tecan Infinite 200 PRO F Nano+ on days 1, 4, and 7.



Figure 5: NuQuant is consistent over time and between plates. The NuQuant standard was used to generate a standard curve and aliquoted into two plates. NuQuant measurements are presented in log, of relative fluorescent units (RFU). A) Measurement of the NuQuant standard on days 1, 4, and 7 provided consistent results demonstrating that the quantification is stable even after multiple freeze/thaw cycles. **B)** The NuQuant standard curve was added to a second PCR plate and concentration was measured over time. The standard curve between the two plates is similar indicating the consistency of measurement. This also indicates that values from a standard curve can be applied to other plates. C) Replicates of the NuQuant standard were added to selected wells. Analysis of the fluorescence demonstrated no difference between samples on the edge of the plate (red text) versus samples in center of the plate. Additionally, blank wells (gray) had a consistently low fluorescence indicating minimal cross-talk across the plate.

NuQuant provides accurate quantification of library molarity







Figure 3: NuQuant accurately measures library molarity. Two users prepared eight libraries each from 10 ng of genomic DNA sheared to 200 or 300 bp using a Covaris system and amplified for 10 or 13 cycles to produce both lower and higher library concentrations. Purified libraries were quantified in duplicate by NuQuant, then equal volumes of all libraries were pooled and sequenced on one lane of an Illumina sequencer. The number of reads for each sample highly correlated with the NuQuant concentration demonstrating the





Figure 6: Consistent library quantification. The total recovered volume from twenty-four libraries were aliquoted into a 96-well, eliminating loss of sample due to standard QC methods. Library molarity was measured with NuQuant on days 1, 4, and 7. Four μ I of each library was taken from each well on day 4 for sequencing. The day 7 fluorescent readings were normalized to account for lower sample volumes. The % CV between the three measurements is 5.6%. The library molarity is consistent even after multiple freeze/thaw cycles.

www.tecan.com

For Research Use Only. Not for use in diagnostic procedures.

For disclaimer and trademarks please visit www.tecan.com · © 2019 Tecan Trading AG, Switzerland, all rights reserved.

CONCLUSIONS

- NuQuant provides a simple method to calculate library molarity eliminating the time and cost associated with current methods.
- Accurate molar concentration can be determined in minutes with a Qubit or a standard plate reader for high-throughput quantification of NGS libraries (tested with Tecan Infinite F Nano, Infinite M Plex, Promega GloMax)
- Use of 96-well PCR plates and a plate reader allows direct library quantification without the sample loss observed with other QC methods

TECAN