



# dsDNA Quantification with the Echo 525 Liquid Handler for Miniaturized Reaction Volumes, Reduced Sample Input, and Cost Savings

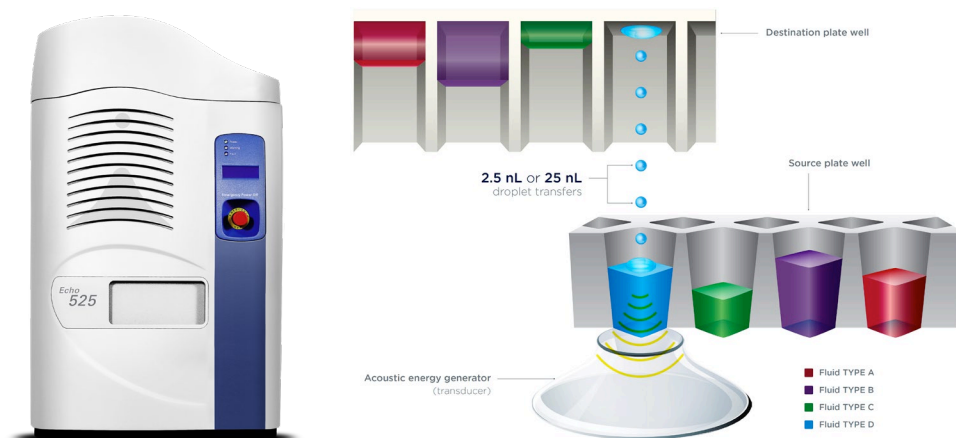
Rabia Khan, Jared Bailey, Jefferson Lai, and John Lesnick

## Abstract

Precise dsDNA quantification is a requirement for many genomics workflows including diagnostics, synthetic biology, and next-generation sequencing. A variety of quantification kits are commercially available, with fluorescence-based kits like the AccuBlue NextGen dsDNA Quantitation Kit and the Quantifluor ONE dsDNA System dominating the market. The AccuBlue NextGen dsDNA Quantitation Kit covers the lower end of dsDNA quantification (3 ng - 5 pg) and the QuantiFluor ONE dsDNA System covers the higher range of dsDNA quantification (50 ng - 20 pg). The kit manufacturer's recommended volumes are 200  $\mu$ L. The main constraints on DNA quantification of reagent cost, hands on time, and required input DNA can be substantially mitigated by miniaturizing the reaction volumes.

The Echo 525 Liquid Handler was used to miniaturize and optimize these fluorescent-based quantification assays. The DNA quantification reaction volumes can be miniaturized up to 20-fold, reducing sample input and allowing more test samples than by manual arraying reagents, with up to 200 more reactions. With optimization and miniaturization of recommended protocols, standard curves generated are equivalent to that of the regular protocols. The precision and accuracy of the miniaturized reaction volumes for the two kits in an Echo 525 Liquid Handler-enabled workflow was verified by quantifying a library of dsDNA samples in a 384-well polypropylene microplate and cross checking the quantification data against the Qubit dsDNA (HS) Assay Kit.

The Echo 525 Liquid Handler utilizes Dynamic Fluid Analysis (DFA), which enables the system to determine fluid composition, height, and impedance in the source well. The required power needed to eject a droplet from the source well is then calculated within milliseconds. The droplet ejected from the source well into an inverted microtiter destination plate is 25 nanoliters (nL) in volume. By transferring a series of droplets large volumes can be pooled in the destination plate at a rate of up to 5  $\mu$ L/sec.



**Figure 1.** (A) Echo 525 Acoustic Liquid Handler. (B) The Echo system transducer rapidly moves between wells on the source plate while the destination plate also moves, allowing rapid transfer from any well to any well for multiple fluid types.

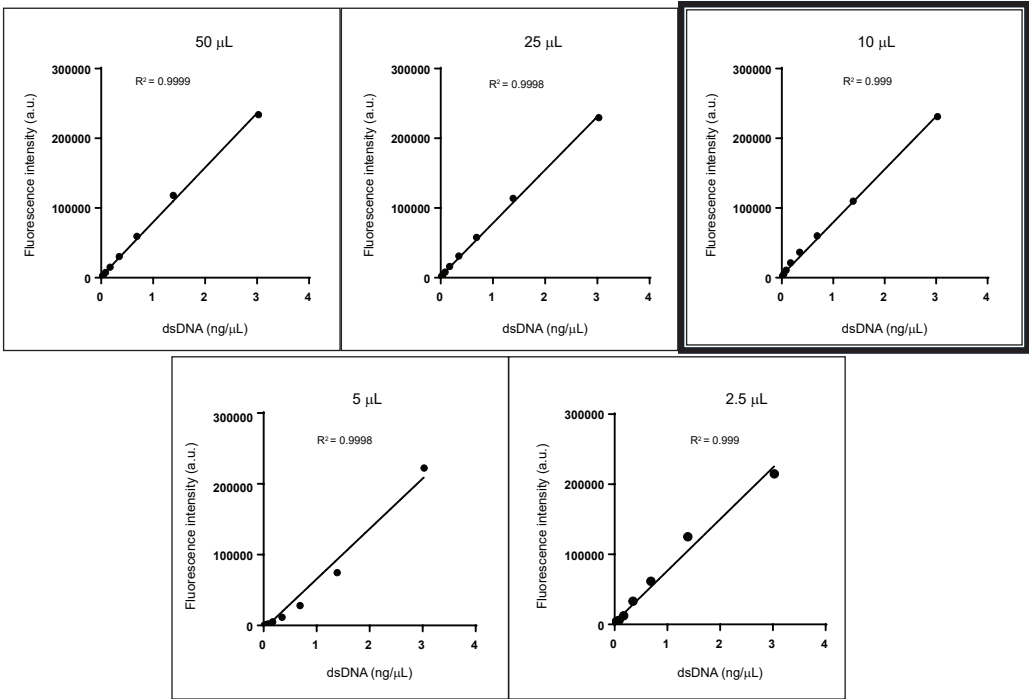
# Assay Optimization

The AccuBlue NextGen dsDNA Quantitation Kit and the QuantiFluor ONE dsDNA System are robust quantification kits used for a wide range of dsDNA with a combined linear range of 5pg - 50ng. Both the kits are fluorescence-based assays and follow a similar protocol. According to the kit-recommended protocols, the reagents are added to a 96-well microtiter plate up to the reaction volume of 200  $\mu\text{L}$ , along with dsDNA serial dilution and unknown sample. This plate is then incubated for 5 minutes in the dark and read on a microplate fluorometer. Fluorescence reads are then computed to generate a standard curve to determine the amount of dsDNA present in the unknown dsDNA sample.

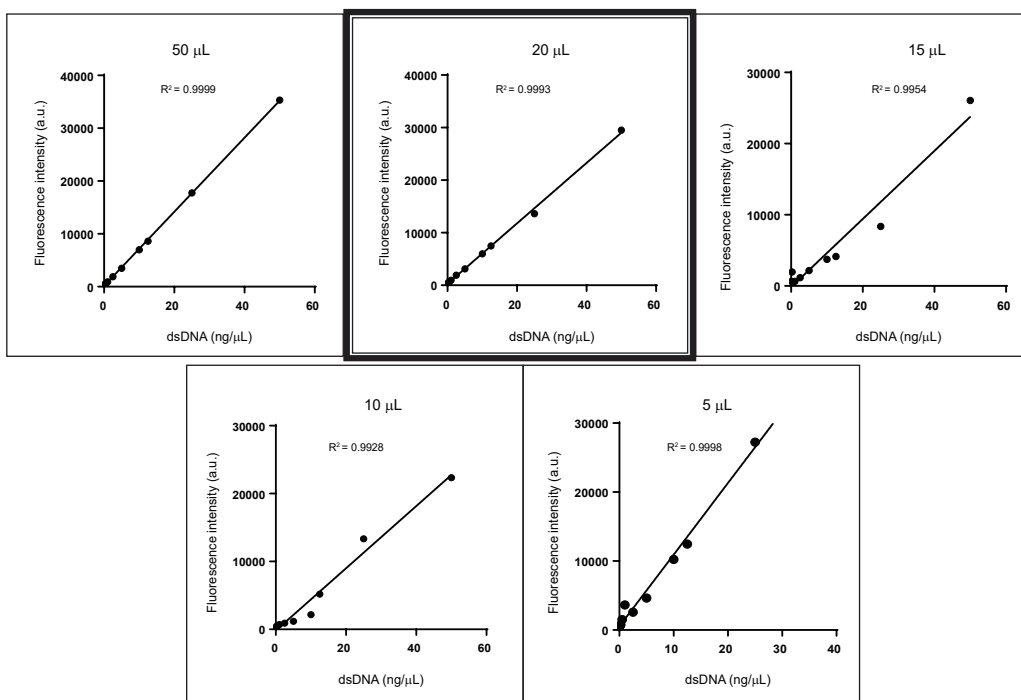
To identify the level of assay miniaturization, both the AccuBlue NextGen dsDNA Quantitation Kit and the QuantiFluor ONE dsDNA System were used to generate standard curves at variable volumes. The volumes chosen for acoustic miniaturization were 50  $\mu\text{L}$ , 25  $\mu\text{L}$ , 10  $\mu\text{L}$ , 5  $\mu\text{L}$ , and 2.5  $\mu\text{L}$  for the AccuBlue NextGen dsDNA Quantitation Kit and 50  $\mu\text{L}$ , 25  $\mu\text{L}$ , 15  $\mu\text{L}$ s, 10  $\mu\text{L}$ , and 5  $\mu\text{L}$  for the QuantiFluor ONE dsDNA System.

The kit's reagents were thawed, equilibrated to room temperature and vortexed to ensure uniform mixing prior to being arrayed in an Echo Qualified 384-well source plate. Standard dsDNA dilutions were prepared afresh with the kit provided dsDNA. For the AccuBlue NextGen dsDNA Quantitation Kit, the working solution was made with AccuBlue NextGen Buffer, AccuBlue NextGen Dye (1:400), and AccuBlue NextGen Enhancer (1:100). The QuantiFluor ONE dsDNA System has only 2 components; fluorescein-based dye and standard dsDNA. A dilution was prepared with standard stock dsDNA for both the kits. The Echo 525 Liquid Handler was used to dispense the reagents and direct dilute standards into the assay plate to generate equivolume reaction mixtures with varying concentrations of dsDNA. The final reaction volumes were incubated for 5 minutes in the assay plates prior to being read on a BMG PHERAstar microplate fluorometer.

## A. AccuBlue NextGen dsDNA Quantitation Kit



## B. QuantiFluor ONE dsDNA System



**Figure 2.** Comparison of multiple standard curves generated via the Echo 525 Liquid Handler for (A) the AccuBlue NextGen dsDNA Quantitation Kit with reaction volume miniaturized up to 4-fold (50  $\mu$ L), 8-fold (25  $\mu$ L), 20-fold (10  $\mu$ L), 40-fold (5  $\mu$ L), and 80-fold (2.5  $\mu$ L), and (B) the QuantiFluor ONE dsDNA System with reaction volume miniaturized up to 4-fold (50  $\mu$ L), 10-fold (20  $\mu$ L), 13-fold (15  $\mu$ L), 20-fold (10  $\mu$ L), and 40-fold (5  $\mu$ L).

## Assay Validation

As shown in Figure 2, each of the reaction volumes in both the kits tested in the initial miniaturization displayed a high level of linearity. However, the reactions at 10  $\mu$ L for the AccuBlue NextGen dsDNA Quantitation Kit and 20  $\mu$ L for the QuantiFluor ONE dsDNA System and above had the most consistent R<sup>2</sup> value. For this reason, these volumes were chosen to be the reaction volume of choice for dsDNA quantification in an Echo 525 system-incorporated workflow.

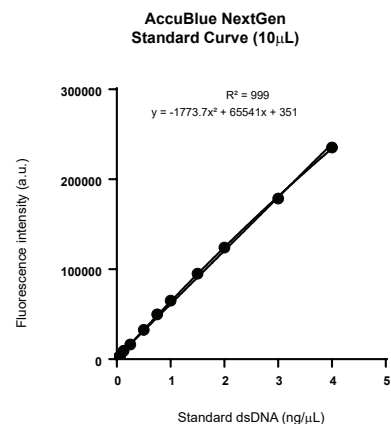
### 1. AccuBlue NextGen dsDNA Quantitation Kit

A 10  $\mu$ L reaction volume enabled us to extend the linear range beyond the AccuBlue NextGen kit-recommended linearity, without compromising the R<sup>2</sup> value. The extended standard curve also enabled us to quantify a broader range of unknown dsDNA samples. Beyond 5 ng we began to lose linearity, so the recommended extended range in an Echo 525 system-incorporated workflow is 4 ng – 0.0025 ng. The kit recommends 10 ng/ $\mu$ L stock dsDNA to generate the standard curve. This range can be achieved by using two stock dsDNA dilutions, 2 ng/ $\mu$ L and 0.1 ng/ $\mu$ L.

In 384-well PS black microtiter plates 10  $\mu$ L reaction volumes were dispensed in quadruplicate via the Echo 525 Liquid Handler, incubated for 5 minutes in the dark, and read on the PHERAstar fluorometer. The average fluorescence intensity of each reaction, standard deviation, and CV were calculated (Table 1) to generate a standard curve (Figure 3). The miniaturized reaction standard curve displays the linearity and R<sup>2</sup> values comparable to the standard curve values provided by the kit.

Source dsDNA (ng/ μL)	Source Vol. (μL)	Buffer (μL)	Enhancer (μL)	Dye (μL)	Reaction volume (μL)	dsDNA (ng)	Average FI Counts	StDev FI Counts	% CV
2	2	7.875	0.1	0.025	10	4	235320.3	3991.9	1.70
2	1.5	8.375	0.1	0.025	10	3	178717.6	3958.1	2.20
2	1	8.875	0.1	0.025	10	2	124089.8	3326.9	2.70
2	0.75	9.125	0.1	0.025	10	1.5	95020.3	1735.6	1.80
2	0.5	9.375	0.1	0.025	10	1	65221.8	1849.2	2.80
2	0.375	9.5	0.1	0.025	10	0.75	49895.6	2013.4	4.00
2	0.25	9.625	0.1	0.025	10	0.5	32683.3	1772.6	5.40
2	0.125	9.75	0.1	0.025	10	0.25	16342.8	886.5	5.40
2	0.075	9.8	0.1	0.025	10	0.125	9377.6	203.2	2.20
0.1	0.625	9.25	0.1	0.025	10	0.0625	3615.8	172.3	4.80
0.1	0.325	9.55	0.1	0.025	10	0.0313	1910.8	89.6	4.70
0.1	0.15	9.725	0.1	0.025	10	0.0150	996.1	38.4	3.90
0.1	0.075	9.8	0.1	0.025	10	0.0075	685.6	25.0	3.60
0.1	0.05	9.825	0.1	0.025	10	0.0050	544.3	23.5	4.30
0.1	0.0375	9.8375	0.1	0.025	10	0.0038	529.3	26.3	5.00
0.1	0.025	9.85	0.1	0.025	10	0.0025	453.6	18.8	4.10

**Table 1.** Protocol for preparing AccuBlue NextGen standard curve in a 10 μL reaction volume. Direct dilution is achieved by dispensing AccuBlue NextGen reagents and standard dsDNA (2 ng/μL and 0.1 ng/μL) via the Echo 525 Liquid Handler in a 384-well PS black, clear bottom microtiter plate.



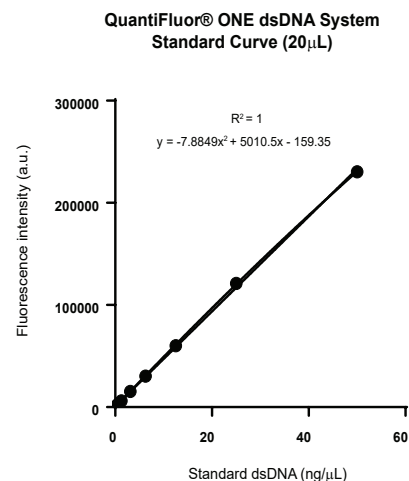
**Figure 3.** Standard curve created from the AccuBlue NextGen dsDNA Quantitation Kit in a 10 μL reaction volume (4ng – 5 pg) with R2 value of 0.999.

## 2. QuantiFluor ONE dsDNA System

A range of 50 ng-20 pg standard curve can be generated by utilizing 12.5 ng/μL and 1 ng/μL standard stock dilutions in a 20 μL reaction volume. This encompasses a wide range for dsDNA quantification without compromising R<sup>2</sup> value. A 20 μL reaction volume was dispensed via the Echo 525 Liquid Handler into a 384-well PS black microtiter plate, incubated for 5 minutes in the dark, and read on a PHERAstar fluorometer. Average fluorescence intensity of each reaction, standard deviation and CV were calculated (Table 2) to generate a standard curve (Figure 4). The standard curve from these miniaturized reaction volumes displays the linearity and R<sup>2</sup> values comparable to the standard curve values provided by the kit.

Source dsDNA (ng/μL)	Source Vol (μL)	Buffer (μL)	Dye (μL)	Reaction volume (μL)	dsDNA (ng/μL)	Average FI Counts	StDev FI Counts	% CV
12.5	4	0	0.025	16	50	230478	1624.6	0.70
12.5	2	2	0.025	16	25	121160	2326.3	1.92
12.5	1	3	0.025	16	12.5	60209	1493.8	2.48
12.5	0.5	3.5	0.025	16	6.25	30305	1019.5	3.36
12.5	0.25	3.75	0.025	16	3.125	15422	255.8	1.66
12.5	0.1	3.9	0.025	16	1.25	6199	135.6	2.19
12.5	0.05	3.95	0.025	16	0.625	3173	82.8	2.61
12.5	0.025	3.975	0.025	16	0.3125	1500	39.4	2.63
1	0.15	3.85	0.025	16	0.15	598	63.3	5.48
1	0.075	3.925	0.025	16	0.075	266	4.8	1.81
1	0.05	3.95	0.025	16	0.05	182	3.7	2.06
1	0.025	3.975	0.025	16	0.025	76	1.8	2.37

**Table 2.** Protocol for preparing QuantiFluor ONE dsDNA System standard curve in a 20 μL reaction volume. Serial dilution is achieved by dispensing QuantiFluor ONE reagents and standard dsDNA (12.5 ng/μL and 1 ng/μL) via the Echo 525 Liquid Handler in a 384-well PS black, clear bottom microtiter plate.



**Figure 4.** Standard curve created from the QuantiFluor One dsDNA Quantitation Kit in a 20 μL reaction volume (50 ng – 20 pg) with R<sup>2</sup> value of 1.

## Assay Verification dsDNA Quantification

In an Echo system-compatible workflow, the same protocol was followed with miniaturized volumes. Reagents were dispensed in quadruplicate by the Echo 525 Liquid Handler to 384-well black microtiter plates with a range of final volumes, incubated for 5 minutes in the dark and read on the PHERAstar fluorometer. Average fluorescence intensity of each reaction, standard deviation, and CV were calculated to generate a standard curve. The miniaturized reaction standard curve displays linearity and  $R^2$  values comparable to the standard curve values provided by the kit.

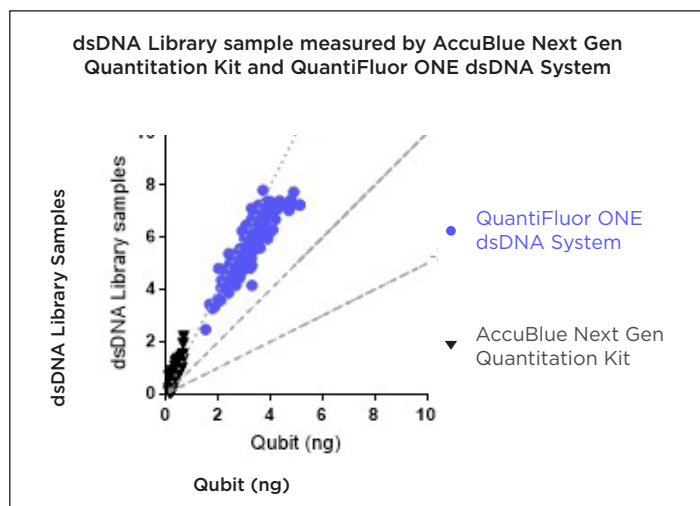
With the help of the Echo 525 Liquid Handler, a library of dsDNA assembled in a 384-well PP plate was quantified in miniaturized volumes by the AccuBlue NextGen dsDNA Quantitation Kit and the QuantiFluor ONE dsDNA System. The above mentioned miniaturized and optimized workflow was adapted for this test. The Echo Plate Reformat Application was used to dispense kit reagents and samples into 384-well PS black microtiter plates followed by a 5-minute incubation and reading the plate on the PHERAstar fluorometer. A second-order polynomial equation was derived from the standard curve, and fluorescence counts were then converted from arbitrary units to ng of dsDNA. The acquired data was plotted against the values of the same dsDNA libraries quantified by the Qubit dsDNA (HS) Assay Kit. The graph in Figure 5 displays a comparable correlation between Qubit values and the ones generated from the 2 kits mentioned above.

## Conclusion

The AccuBlue NextGen dsDNA Quantitation Kit and the QuantiFluor ONE dsDNA System can be miniaturized up to 20-fold using the Echo 525 Liquid Handler. Kit reagents can be dispensed in miniaturized volumes using the Echo Plate Reformat Application into a destination plate.

Standard curves generated for both kits in miniaturized reaction volumes display  $R^2$  values comparable to the kit-recommended protocols, which can be used to determine the amount of dsDNA present in a sample volume as low as 0.025  $\mu\text{L}$ .

These kits can be effectively used to quantify dsDNA library samples via the Echo 525 Liquid Handler. This can mitigate the constraints on DNA quantification of reagent cost, hands-on time, and required input DNA, and be highly beneficial for applications such as synthetic biology, NGS library prep, or quantifying precious low concentrations of unknown dsDNA samples.



**Figure 5:** dsDNA NGS library quantified by the AccuBlue NextGen dsDNA Quantitation Kit and the QuantiFluor ONE dsDNA System and plotted against the values of the same dsDNA libraries quantified by the Qubit dsDNA (HS) Assay Kit.

Equipment	Manufacturer	
Echo 525 Liquid Handler	Beckman Coulter Life Sciences	
Allegra X-14 Centrifuge	Beckman Coulter Life Sciences	
BMG PHERAstar	BMG Labtech	
Qubit	ThermoFisher	
<b>Materials for AccuBlue NextGen dsDNA Quantitation Kit Biotium #31060</b>	<b>Kit Recommended</b>	<b>1/10</b>
Reaction volume	200 µL	10 µL
AccuBlue NextGen Dye, 400X	0.5 µL	0.025 µL
AccuBlue NextGen Enhancer, 100X	2 µL	0.1 µL
AccuBlue NextGen Buffer, 1X	197.5 µL	>-9 µL
High Sensitivity DNA Standard	10 µL	(0.025--2) µL
Reactions per kit	1250	250,000
<b>Materials for QuantiFluor ONE dsDNA System Promega #E4871</b>	<b>Kit Recommended</b>	<b>1/20</b>
Reaction volume	200 µL	20 µL
AccuBlue NextGen Dye, 400X	200 µL	16 µL
High Sensitivity DNA Standard	1 µL	(4-0.025) µL
<b>Consumables</b>	<b>Manufacturer</b>	<b>Part Number</b>
384-well PP Microplate	Beckman Coulter Life Sciences	001-14555
384-well LDV Microplate	Beckman Coulter Life Sciences	001-12782
384-well Black Clear-Bottom Microplate	Greiner	#781096
384-well Black µClear Microplate, Small Vol, LoBase	Greiner	#788096
TapeStation Plate	Agilent	#5067-5150
Qubit dsDNA HS Assay Kit / Qubit Microtube	ThermoFisher	#Q32851 / #Q32856

This protocol is for demonstration only, and is not validated by Beckman Coulter. Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions.

Biomek Automated Workstations and Echo Liquid Handlers are not intended or validated for use in the diagnosis of disease or other conditions.



©2020 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)

AAG-6742APP03.20