



# Modular DNA Assembly of PIK3CA Using Acoustic Liquid Transfer in Nanoliter Volumes

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

## Abstract

Previous work has shown the ability of the Echo Liquid Handler to generate DNA constructs using various assembly chemistries with a miniaturized protocol<sup>1,2</sup>. Interchanging smaller, modular pieces of DNA is the preferred method for numerous workflows due to decreased synthesis cost and faster progression along the design, build, test, and learn cycle. In this study, NEBuilder HiFi was combined with the 25 nL increment dispensing Echo 525 Liquid Handler to assemble a modular oncogene construct encoding phosphatidylinositol-4,5-bisphosphate 3-kinase catalytic subunit alpha (PIK3CA) over five modular pieces (Figure 2A). The modules themselves were designed to contain four activating mutations: R38H, E542K, E545K, and H1047R along with the wild-type PIK3CA in a pcDNA 3.1/Zeo(+) Vector for a plasmid size of 8,940 bp. The wild-type reaction was miniaturized down with a forty-fold reduction from the NEBuilder HiFi recommended volume in a previous experiment 2. At the forty-fold reduced volume, 1.25 fmoles of DNA was all that was required for each piece in a total reaction volume of 500 nL. The goal of this experiment was to utilize the Echo 525 Liquid Handler in assembling five, five-piece assemblies by substituting modular DNA pieces at the nanoliter scale. The experiment was validated by junction colony qPCR for rapid analysis of successful assembly and Illumina MiSeq sequencing verification for nucleotide resolution. The Echo Liquid Handler enables lower-cost methods and workflows to produce high-quality synthetic DNA constructs which expands design-based testing with higher throughput and affords the scientist a broader biological landscape to interrogate.

## Introduction

Synthetic biology is an interdisciplinary science with the potential to impact academic and industrial applications including the creation of novel therapeutics and vaccines, plant science and biofuels, as well as bio-based chemical manufacturing capabilities that involve the application of engineering principles to biology. The focus is often on generating parts of natural biological systems, characterizing and isolating them, and then using them as components of an engineered biological system. A trademark of synthetic biology is the application of rational design principles to the design and assembly of biological components. The outcome of introducing a rationally designed synthetic DNA construct into a cell is difficult to predict. This creates the need to test multiple permutations to obtain the desired outcome. A greater emphasis on the modular design of DNA parts enables the assembly of a greater variety of potential constructs through combinatorial assembly of synthetic components. In addition to simplifying the overall DNA construction workflow, modular design and assembly of DNA components makes automation of the assembly process possible reducing the time, labor, and cost of generating multiple constructs to allow for an increase of throughput with an overall shortened development cycle.

Synthetic DNA plasmids are designed using computer-aided software with experimentally dependent functionality. This can range from interrogation of different domains of a single protein to the production of an entire pathway with heterologous genes. The designed DNA is then divided into synthesizable pieces. The pieces are designed with overlapping sequences to their neighbors before being chemically synthesized. The DNA fragments are then assembled together to build the designed construct using gene assembly techniques. If needed, multiple pieces can be assembled together into larger DNA assemblies. The assembled constructs are then typically cloned into an expression vector and sequence-verified. Once verified, the synthetic constructs are transferred into a production cell and the function of the designer construct is assayed. Depending on the results the constructs can then be modified or refined, and the test cycle is repeated until a DNA construct is obtained that produces the desired function.

New England Biolabs has developed the NEBuilder HiFi DNA Assembly Cloning Kit (Figure 1). The NEBuilder HiFi kit takes input DNA with 15 to 30 base pairs of terminal sequence identity and generates overhangs by using a proprietary 5' exonuclease. The complementary 3' overhangs are subsequently filled in with a DNA polymerase upon annealing. DNA ligase is used on the final product to seal nicks and create a continuous assembly. The entire reaction is incubated at a temperature of 50°C for as little as 15 minutes in a recommended volume of 20 µL.

#### DNA Preparation

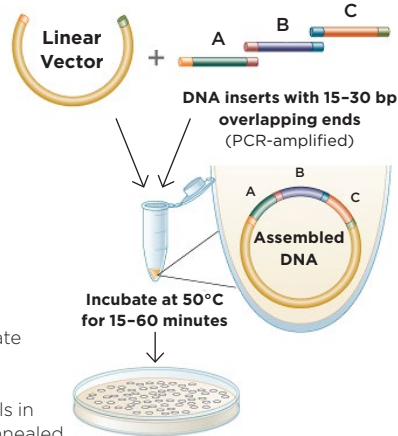
From:

- PCR
- Restriction enzyme digestion
- Synthetic DNA (e.g., gBlocks)

#### NEBuilder HiFi DNA Assembly Master Mix

Single-tube Reaction

- Exonuclease chews back 5' ends to create single-stranded 3' overhangs
- DNA polymerase fills in gaps within each annealed fragment
- DNA ligase seals nicks in the assembled DNA



**Figure 1.** The typical workflow using the NEBuilder HiFi kit showing various input DNA types, required overlaps, and general reaction conditions. Reprinted from [www.neb.com](http://www.neb.com) (2018) with permission from New England Biolabs, Inc.

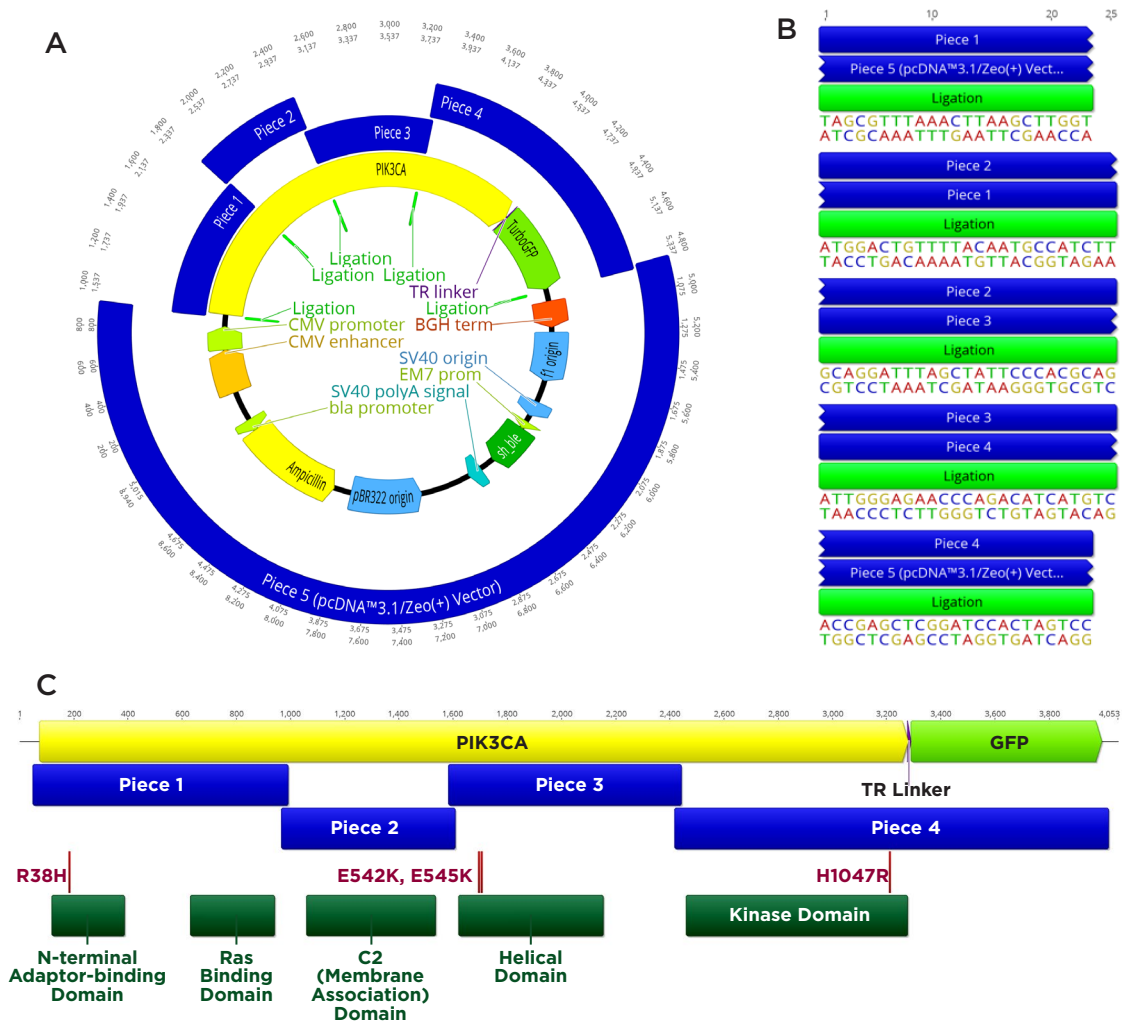
#### Transformation

Previous work has shown the ability of the Echo 550 Liquid Handler to generate two-piece assemblies using the Golden Gate and Gibson Assembly chemistries utilizing a miniaturized protocol <sup>1</sup>. The Echo 550 Liquid Handler uses a transducer to acoustically dispense in increments of 2.5 nL. In another prior experiment, NEBuilder HiFi chemistry was combined with the 25 nL increment dispensing Echo 525 Liquid Handler to generate a two and five-piece assembly at the 500 nL volume <sup>2</sup>.

Phosphatidylinositol-4,5-bisphosphate 3-kinase catalytic subunit alpha (PIK3CA) is a commonly studied gene in human oncology. The mutation of specific domains from this gene have been correlated with multiple tumor types <sup>3-7</sup>. Due to PIK3CA's importance in oncology, it was chosen to demonstrate a real-world application of an Echo 525 system-driven miniaturization of the NEBuilder HiFi kit. PIK3CA was divided into four modular pieces based on domains of interest <sup>7</sup> and was tagged with C-terminal GFP (Figure 2A). Four known oncogenic mutations, R38H, E542K, E545K, and H1047R were designed in modular pieces along with a wild-type control (Figure 2B).

The construct was also engineered to assemble into a KpnI-HF digested pcDNA 3.1/Zeo(+) Vector to allow for future transfection. The vector is interchangeably referred to as piece five after being linearized. The pieces were designed with 23 to 25 base pair DNA overlaps (Figure 2C), assembling five, five-piece assemblies by substituting modified modular DNA pieces at the nanoliter scale.

The Echo Liquid Handler enables lower-cost methods and workflows to produce high-quality synthetic DNA which expands design-based testing with higher throughput and affords the scientist a broader biological landscape to interrogate. Ultimately, the utilization of the Echo Liquid Handler for synthetic biology can help close the DNA read-write cost gap.



**Figure 2.**

**A.** The complete plasmid construct to be assembled with GFP-tagged PIK3CA being generated in a KpnI-HF Digested pcDNA 3.1/Zeo(+) backbone. gBlock Pieces ordered as well as linear vector are shown in blue, with their overlapping regions shown by a light green annotation labeled ligation. The PIK3CA gene is shown in yellow with the linker (purple) and turboGFP tag (green) also annotated. Other important parts of the plasmid machinery from pcDNA 3.1/Zeo(+) are also shown.

**B.** Nucleotide resolution for each of the designed overlaps in the assembly. This is enlarged from the Figure 2A ligation regions. Five overlaps were needed for the modular assembly.

**C.** The locations of the PIK3CA functional domains (dark green) and the planned mutations (red) within the constructs. The protein domains relationship with the synthetic piece design (blue) is also shown. The full length PIK3CA (yellow) gene is shown with its C-terminal GFP tag (light green). Each construct is designed to be either wild type or have one mutation. The planned mutations are R38H, E542K, E545K, and H1047R.

## Modular DNA Assembly

### Rational Construct Design

A modular DNA assembly was also designed to be used in the miniaturized reaction volumes to more closely simulate the challenges faced in real-world applications. The PIK3CA cancer gene was chosen to be broken up into four pieces based on domains. These pieces were ordered as either wild-type sequence or containing the known oncogenic mutations, R38H, E542K, E545K, and H1047R. The assembly was also designed to insert into a linear pcDNA 3.1/Zeo(+) mammalian vector. This vector is interchangeably referred to as piece 5 after being linearized with KpnI-HF. All the pieces were designed with 23 to 25 base pair DNA overlaps and ordered from Integrated DNA Technologies as gBlocks Gene Fragments. The vector and gBlocks were amplified to obtain the necessary amount of modular DNA required for the study.

### Modular DNA Component Production (gBlocks Pieces 1-4)

The primers listed in Table 1 were ordered from Integrated DNA Technologies (IDT) to generate additional copies of the modular DNA for assembly. The high-fidelity polymerase Q5 was used for the PCR reactions as shown in Table 2. Each reaction was set up in a 1.5 mL tube, gently mixed, and aliquoted into a 384-well Bio-Rad PCR plate. The reactions were then run on an Applied Biosystems ProFlex PCR System with the reaction conditions shown in Table 3.

**gBLOCK Amplification Primers**

Piece #	Forward Primer Sequence	Reverse Primer Sequence
Piece 1-WT	TAGCGTTTAAACTTAAGCTTGGTACATGCCTC	AAGATGGCATTGTAAAACAGTCCATTGGC
Piece 1-R38H	TAGCGTTTAAACTTAAGCTTGGTACATGCCTC	AAGATGGCATTGTAAAACAGTCCATTGGC
Piece 2-WT	AGAAAGCCTTATTCTCAACTGCCAATGGAC	CTGCGTGGAATAGCTAAATCCTGCTTC
Piece 3-WT	GCAGGATTTAGCTATTTCCACGCAG	GACATGATGTCTGGGTTCTCCCAATTCAA
Piece 3-E542K	GCAGGATTTAGCTATTTCCACGCAG	GACATGATGTCTGGGTTCTCCCAATTCAA
Piece 3-E545K	GCAGGATTTAGCTATTTCCACGCAG	GACATGATGTCTGGGTTCTCCCAATTCAA
Piece 4-WT	GGACTAGTGGATCCGAGCTCGGTAC	ATTGGGAGAACCCAGACATCATGTCCAGAG
Piece 4-H1047R	GGACTAGTGGATCCGAGCTCGGTAC	ATTGGGAGAACCCAGACATCATGTCCAGAG

**Table 1.** Primer sequences used to PCR amplify the gBlock template to increase concentration of necessary DNA for subsequent DNA assembly.

**Piece PCR Reaction Conditions**

Reagent	Volume ( $\mu$ L) per reaction
Forward Primer (10 $\mu$ M)	10
Reverse Primer (10 $\mu$ M)	10
NEB Q5 Polymerase 2x MM	100
IDT gBlock DNA (1ng/ $\mu$ l)	2
diH <sub>2</sub> O	78

**Table 2.** Volumes of reagents per PCR reaction used for template DNA expansion. The total reaction was dispensed in 20  $\mu$ L aliquots into a 384-well Bio-Rad PCR plate. Q5 polymerase was chosen due to the need for a low error rate with the amplified pieces for downstream sequence fidelity.

### Thermocycler Reaction Conditions

Piece #	Initial Denaturation	25 Cycles			Final Extension	Hold
Time	30 seconds	10 seconds	30 seconds	1 minutes	2 minutes	infinite
Temperature	98oC	98oC	71oC	72oC	72oC	4oC

**Table 3.** Proflex thermocycler reaction conditions used for the amplification of the insert piece DNA (Pieces 1-4). The reactions were held to 25 cycles to minimize the potential to induce mutations.

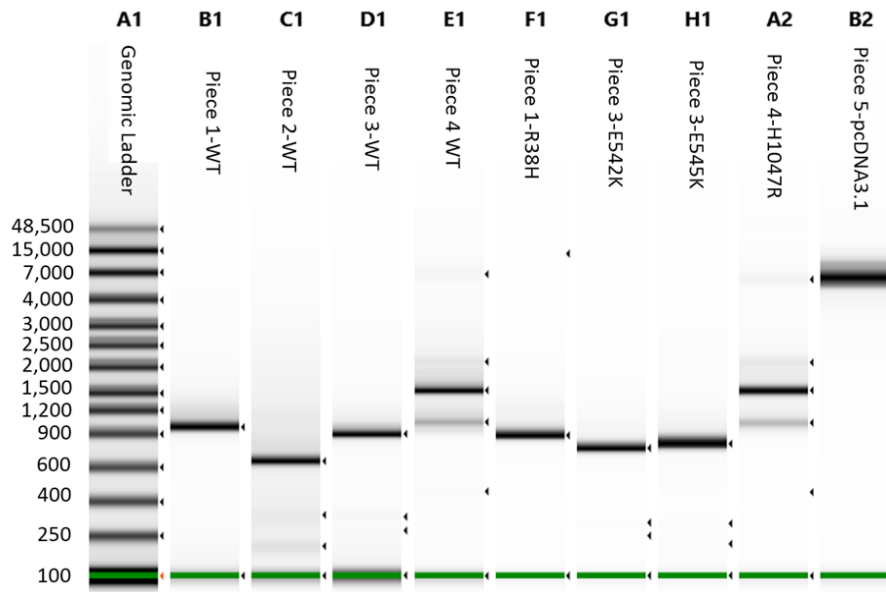
### Vector DNA Component Production (pcDNA 3.1/Zeo (+) Piece 5)

The circularized mammalian vector pcDNA 3.1/Zeo (+) was purchased from Thermo Fisher Scientific and transformed into NEB 10-beta competent E. coli. A clonal transformant was frozen in glycerol at -80°C for use in the entirety of the experiment. An overnight 25 mL culture was set up in a 250 mL Erlenmeyer Flask in LB Broth with 100 µg/mL carbenicillin and incubated at 37°C. It was split between 5 Qiaprep columns using the manufacturer recommended protocol and eluted in 250 µL. The eluted plasmid DNA was then digested overnight at 37°C with the addition of 20 µL NEB KpnI-HF and 30 µL 10x Fast Buffer.

### DNA Cleanup, Quantitation, and Visualization (Pieces 1-5)

Upon reaction completion, both the vector restriction digest and gBlock PCRs were cleaned and concentrated using an NEB Monarch PCR & DNA Cleanup Kit (5 µg). The reactions were processed according to the manufacturer’s protocol and eluted in 11 µL of elution buffer.

A 1 µL aliquot for each reaction was run on an Agilent 2200 Tapestation loaded with a Genomic DNA Analysis ScreenTape using their given protocol (Figure 3). The DNA was visually checked for quality, and the concentration of the expected band size was used as input for DNA normalization in DNA assembly.



**Figure 3.** Agilent 2200 Tapestation with Genomic tape run of the cleaned DNA assembly pieces. Tapestation concentration value for each correct band was used to calculate the proper reaction addition of each piece.

## DNA and NEBuilder HiFi Master Mix Arraying

Using the DNA concentrations from the Agilent 2200 Tapestation run in Figure 3, the DNA was diluted to a desirable range for assembly (Table 4). The diluted DNA was then transferred into an Echo Qualified 384-Well Low Dead Volume Plus Microplate (384 LDV Plus). NEBuilder HiFi 2x Master Mix was also separately added to the 384 LDV Plus plate to allow for a low dead volume of enzyme master mix. The NEBuilder HiFi protocol has a recommended assembly volume of 20  $\mu\text{L}$  with an input of 0.5 pmol per piece for complex assemblies. Prior work has demonstrated this reaction to be robust at the 500 nL volume with 1.25 fmol of each piece. After arraying and sealing, the reaction was conducted for 1 hour at 50°C in an Applied Biosystems ProFlex PCR System with a heated lid. The PCR plate was cooled to 4°C before being transformed into NEB 10B E. coli cells.

NEBuilder Reaction Conditions								Echo Calibration
Reagent	DNA Concentration (ng/ $\mu\text{L}$ )	ng DNA/1.25 fmol	WT Reaction (nL)	R38H Reaction (nL)	E542K Reaction (nL)	E545K Reaction (nL)	H1047R Reaction (nL)	
Piece 1 WT	23.6	16	25		25	25	25	384LDV_Plus_AQ_GP
Piece 1 R38H	25	16		25				
Piece 2 WT	14.6	11	50	50	50	50	50	
Piece 3 WT	7.69	14	75	75			75	
Piece 3 E542K	10	14			75			
Piece 3 E545K	10	14				75		
Piece 4 WT	29.6	26	50	50	50	50		
Piece 4 H1047R	30	26					50	
Piece 5	83.9	83	50	50	50	50	50	
NEBuilder Master Mix			250	250	250	250	250	

- 1 hr 50oC Incubation (Reaction in Thermocycler with Heated Lid)

**Table 4.** Volumes of reagents used for the modular DNA assembly. The total reaction volumes were transferred by the Echo 525 Liquid Handler. These reagents were all dispensed from a 384 LDV Plus plate. These reactions were all transformed into E. coli.

## Transformation/Construct Production

The modular DNA assemblies were transformed into 12.5  $\mu\text{L}$  of chemically competent NEB 10B E. coli following the conditions outlined in Table 5. In addition to the five outlined DNA constructs, pUC19 was also transformed as a plasmid control. The volume of cells listed in the table was pipetted into the 384 PCR plate containing the reactions that had been chilled to 4°C. After a 30-minute incubation for the mixture at 4°C, the cells were heat shocked at 42°C for 30 seconds then immediately transferred back to the 4°C cold block.

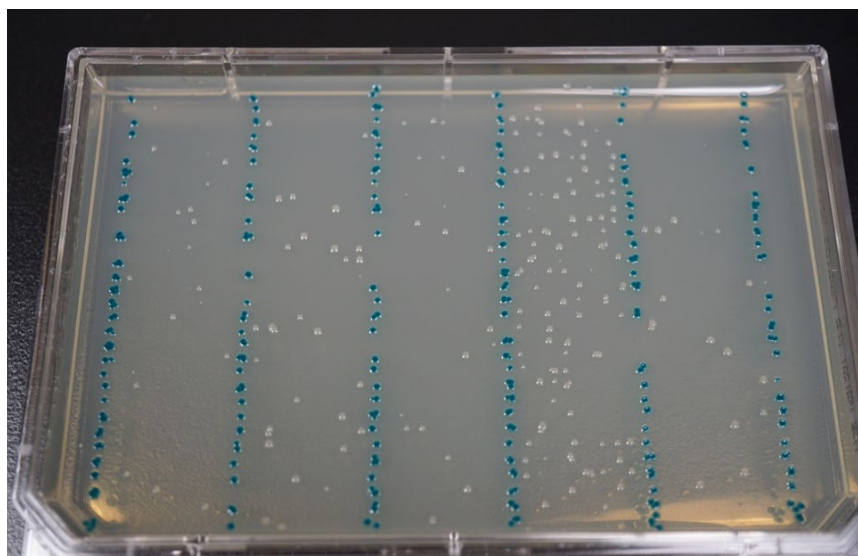
Modular DNA E. Coli Transformations					
	WT Reaction (nL)	R38H Reaction (nL)	E542K Reaction (nL)	E545K Reaction (nL)	H1047R Reaction (nL)
NEB 10B cells	12500	12500	12500	12500	12500

- 30 Minutes 4oC incubation (Cold Block)
- 30 Seconds 42oC (Heat Shock in Thermocycler)
- 2 minutes 4oC (Cold Block)

**Table 5.** Transformation conditions for NEB 10B E. coli cells into the various modular DNA assembly reactions. The cells were all added directly to the assembly reaction contained in the 384-well PCR plate.

## E. coli Plating

The transformed E. coli were manually transferred into a 384-Well LDV Plus plate after cooling and gently mixed by pipetting. LB Agar OmniTrays with 100 µg/mL carbenicillin, 60µg/ml X-Gal, and 0.1mM IPTG were warmed and dried in the 37°C incubator for 1 hour. The Echo calibration 384LDV\_Plus\_AQ\_GP was used to array 8 columns of 25 nL per construct transformation onto OmniTrays (Figure 4). Each experimental reaction was separated by a single 25 nL column of the pUC19 control. The plates were inverted and incubated overnight at 37°C. Isolated colonies were subsequently used for verification by colony qPCR and NGS. The resulting experimental colonies plates were used in both colony qPCR and NGS sequencing.



**Figure 4.** E. coli transformations spotted by an Echo 525 Liquid Handler onto an LB Agar OmniTray with 100 µg/mL carbenicillin, 60µg/ml X-Gal, and 0.1mM IPTG. Each reaction was bracketed by a pUC19 control strain which is visible in blue. The pUC19 transformant contained an intact β-galactosidase pathway, which turns a blue color upon exposure to IPTG and X-Gal, so any cross-contamination between spots would be visible. From left to right, the reactions are: WT, R38H, E542K, E545K, and H1047R.

## Colony qPCR Quality Control

The Echo 525 Liquid Handler was used to accurately dispense 250 nL of a cell suspension into four 10  $\mu$ L qPCR reactions targeting four different junction regions of the assembled plasmid construct as a means of quality control for correct plasmid assembly (Figure 5). A qPCR reaction master mix was made for each of the four reactions containing the amplicon primer pair and LightCycler 480 SYBR Green I Master as shown in Table 6. The reagent mixture minus the colony suspension was bulk dispensed into a barcoded Bio-Rad Hard-Shell Skirted 384-Well PCR Plate. After reagent addition, the qPCR plate was centrifuged for 1 minute at 1000  $\times$  g. Isolated colonies were picked using a P10 pipette tip into 50  $\mu$ L of diH<sub>2</sub>O in an Echo Qualified 384-Well Polypropylene Microplate 2.0 (384-Well PP). The cell-containing 384-Well PP plate was sealed and briefly mixed at 2000 RPM for 2 minutes on an Eppendorf MixMate. From this, 250 nL of the cell suspension was dispensed using the Echo calibration 384PP\_AQ\_BP into each of the four qPCR reagent mixes. The qPCR plate was then sealed with optical film, vortexed for 2 minutes at 2000 RPM on the Eppendorf MixMate, and briefly centrifuged at 1000  $\times$  g for 1 minute. After centrifugation, the sealed plate was run according to the conditions specified in Table 7. Colonies with all three junction CPs closely matched to the Ampicillin control were called positive assemblies and moved to Next-Generation Sequencing (Figure 6).

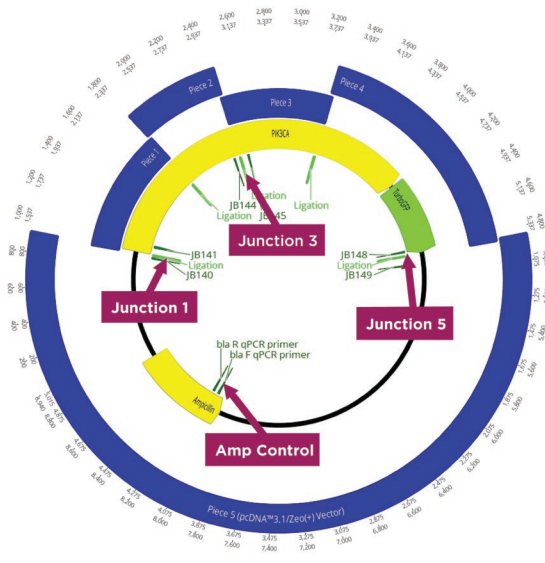
Colony qPCR Reaction Volumes		
Reagent	Volume (nL) per reaction	Echo Calibration
Forward Primer (10 $\mu$ M)	250	
Reverse Primer (10 $\mu$ M)	250	
LightCycler 480 SYBR Green I Master	5000	
diH <sub>2</sub> O	4250	
Colony Cell Suspension	250	384PP_AQ_BP

**Table 6.** Reagent additions used for the modular assembly colony qPCR reactions. Four different primer pairs were used to amplify the target regions from the colonies. Note that the Primers, SYBR Green Master Mix, and water were made as a master mix and dispensed together using a pipette. 250 nL of the cell suspension was then dispensed into each of the four different reaction conditions using the Echo 525 Liquid Handler.

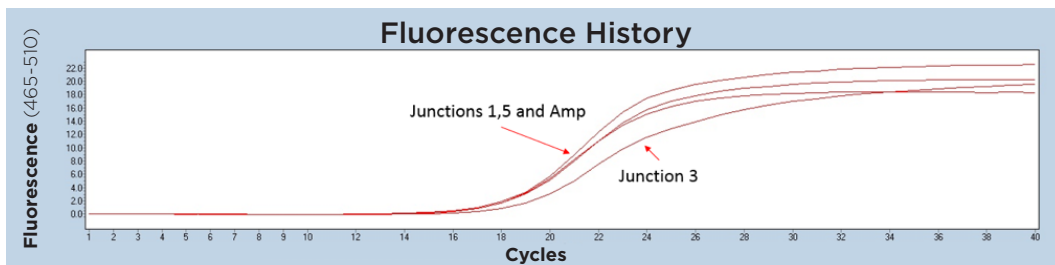
Colony qPCR Reaction Conditions					
	Temp. (oC)	Hold (sec)	Ramp Rate oC/sec	Acquisition (465 nm/520 nm)	Cycles
Pre-incubation	95	300	4.8		1
Amplification Piece 3 WT	95	10	2.5	Single	40
	50	10	2.5		
	72	10	2.5		
Melting Curve Piece 3 E545K Piece 4 WT	95	5	4.8	5/oC	1
	65	60	2.5		
	97	continuous	0.11		
Cooling	4	30	2.5		1

**Table 7.** Roche Lightcycler 480 conditions run for the DNA assembly junction qPCR looking for three generated junctions in comparison to the Ampicillin control.





**Figure 5.** Plasmid Map showing the five-piece modular assembly construct with the 4 targeted qPCR amplicons highlighted. If Junctions 1, 3, and 5 had a similar CP to the Ampicillin control, the colony was considered positive by colony qPCR. By testing these junctions, the presence of all five pieces could be seen.



**Figure 6.** Lightcycler 480 trace of a construct positive for all three experimental junctions in comparison to the Ampicillin control.

## Miniaturized Nextera XT and MiSeq Plasmid Sequencing

The same cell suspensions used in the qPCR reaction were used by the Echo 525 system to spot onto an LB+Carb100 OmniTray. These spots were then expanded into 4, 96 deep well plates with 1.0 mL of LB+Carb100 liquid media. Each colony was set up in duplicate wells (2 mL media per colony) and incubated overnight at 37°C with shaking. The cell cultures were combined the next day and run through a Macherey-Nagel NucleoSpin 96 Plasmid Kit according to the manufacturer's protocol. The DNA was eluted in diH<sub>2</sub>O and 50 µL transferred into an Echo Qualified 384-well PP plate. The Quant-iT Picogreen dsDNA Assay Kit was used to quantify the amount of plasmid present in black walled Greiner 384-well polystyrene plates with a miniaturized protocol. First, 9 µL of 1xTE was added to all sample wells through a bulk dispense. The Echo 525 Liquid Handler was then used to add 1 µL of each DNA sample. The Echo 525 Liquid Handler also dispensed a standard curve in triplicate using the provided DNA standard. The curve ranged between 0.025 and 1 ng with a 10 µL final volume in TE. 10 µL of 1:200 diluted Picogreen dye was then added to all wells. The plate was sealed and mixed on the MixMate at 2000 rpm for 2 minutes before being centrifuged at 100 xg for 1 minute. The BMG Pherastar was used for the Picogreen quantitation with the excitation and emission peaks at 485 nm and 528 nm respectively. Based on these results, the DNA was normalized to provide 0.1 ng per tagmentation reaction. The Nextera XT tagmentation reaction and indexing reactions were then dispensed as shown in Table 8.

### Nextera XT Reaction Conditions

Tagmentation	Miniaturized (μL/rxn)	Echo Calibration
Plasmid DNA Sample	0.5	384PP_AQ_BP
TD Buffer	1	384PP_AQ_GP
ATM	0.5	384PP_AQ_GPSB
<b>Total volume</b>	<b>2</b>	

- Spin @ 1500g, 1 minute -> Tagmentation reaction -> 55°C (5 min) --> 10°C (forever)

Add NT	Miniaturized (μL/rxn)	Echo Calibration
Tagmentation reaction	2	From previous (no transfer)
NT Buffer	0.5	384PP_AQ_SPHigh
<b>Total volume</b>	<b>2.5</b>	

- Spin @ 1500g, 1 minute -> Incubate 5 min RT

Indexing Amplification	Miniaturized (μL/rxn)	Echo Calibration
Tagmentation reaction	2.5	From previous (no transfer)
Indexing Primer 1 (100μM) (N7XX)	0.05	384PP_AQ_BP
Indexing Primer 2 (100μM) (55XX)	0.05	384PP_AQ_BP
diH <sub>2</sub> O	0.9	384PP_AQ_BP
NPM	1.5	384PP_AQ_GPSB
<b>Total volume</b>	<b>5</b>	

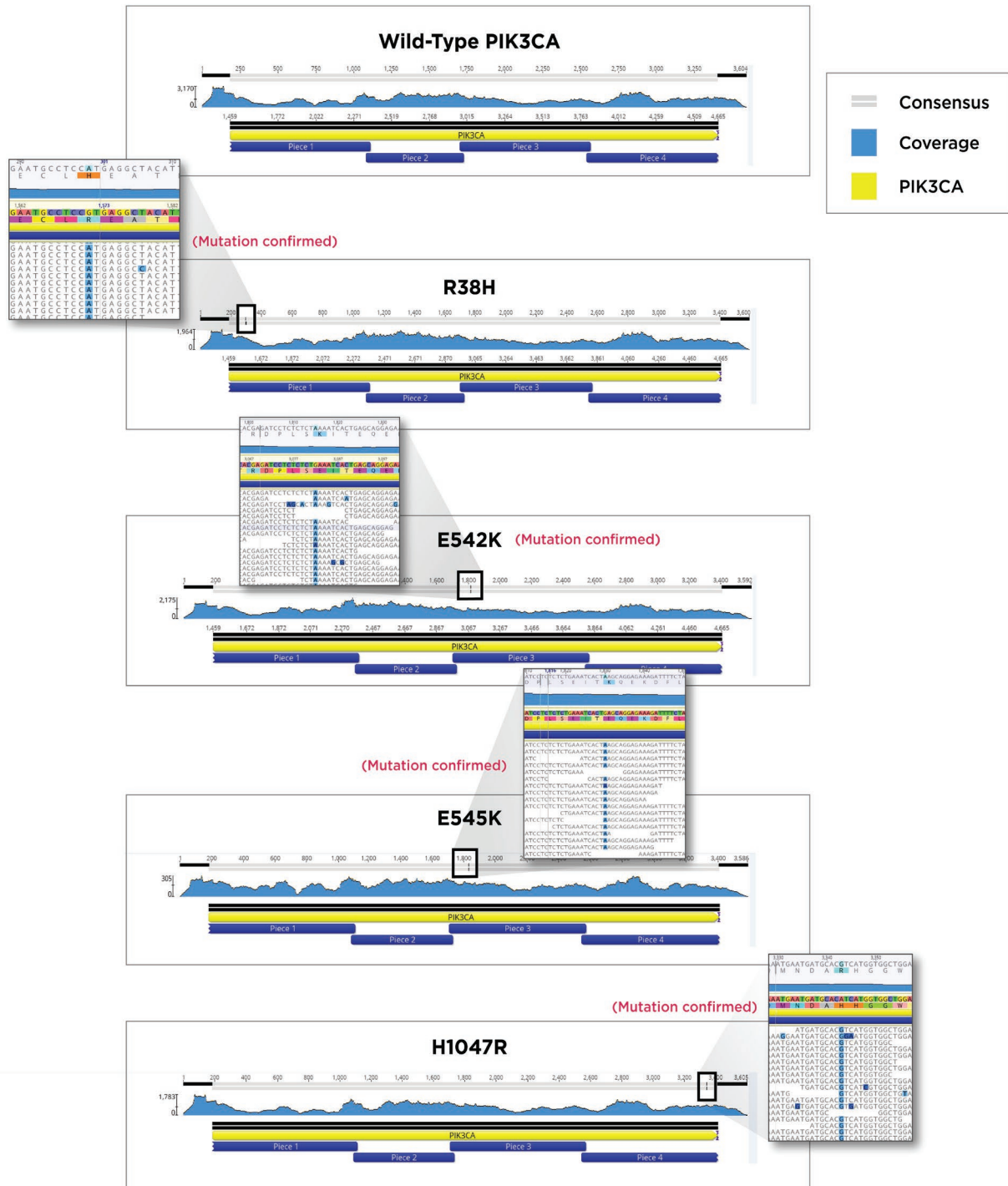
PCR Reaction	
72°C	3 min
95°C	30 sec
95°C	10 sec
55°C	30 sec
72°C	30 sec
72°C	5 min
4°C	hold

} 14 x

**Table 8.** Nextera XT Plasmid Tagmentation and Indexing.

After indexing, Agencourt AmpureXP beads were used to clean up the reactions. The SPRI bead cleanup followed the Agencourt AmpureXP PCR purification protocol for 384-well format. We utilized 9μL of beads for our 5μL indexing amplification reaction, eluted in 20μL, and removed 15 μL with care not to carry over any beads. Quantitation was performed to verify the quantity of each sample using the Picogreen quantitation assay again as described above. Fragment size analysis was performed on the Agilent TapeStation 2200 with a TapeStation D1000 HS Kit, according to the Agilent standard protocol. Using the concentration data from the Picogreen assay, as well as fragment size information from the TapeStation 2200, we built a normalization worklist in Excel to obtain equimolar pooling of samples. The Echo 525 Liquid Handler was then used to simultaneously pool and normalize the libraries. The pool was then denatured and diluted to 20pM, loaded with 1% PhiX control, and run on an Illumina MiSeq

specifying 2x75 reads. FASTQ files were generated on the MiSeq and aligned to an unmutated PIK3CA using Geneious 11.0. Correct PIK3CA constructs were obtained for each of the five desired assemblies as seen in Figure 7.



**Figure 7.** MiSeq reads from each of the five constructs aligned to an unmutated PIK3CA gene. Sequence coverage and location of the mutation in the consensus read bar are shown at the top of each alignment. A magnified region containing the mutated amino acid residues are shown next to the four mutated versions of PIK3CA.

## Discussion

Increasing the throughput of DNA construction is a goal of many research groups. The use of modular DNA pieces within a construction pipeline is a standard practice that offers the benefit of both lower synthesis cost as compared to a full-length gene and faster lead consolidation upon successful screening. The Echo 525 Liquid Handler, in combination with the NEB NEBuilder HiFi kit, can further enable this technique through the miniaturization of the reactions and workflow enhancements.

The use of a miniaturized 500 nL reaction volume while varying modular component pieces for a desired mutation was the primary aim of this experiment. The NEBuilder HiFi kit reaction volume was decreased forty-fold from the recommended volume from 20  $\mu$ L to 500 nL. At the 500 nL reduced volume, only 1.25 fmoles of DNA was used for each input piece, which would dramatically decrease the amount of DNA required from that of the recommended protocol which was 50 nmoles per piece. Obtaining sufficient high-quality input DNA is challenging, so the decrease of the required number of DNA molecules can vastly increase the number of modular assemblies possible. The use of these smaller volumes was done for PIK3CA constructs containing the mutations identified as R38H, E542K, E545K, and H1047R as well as wild-type in a mammalian expression vector. Correctly assembled clones were obtained for each of these targets as seen in the alignment from Figure 7. Successfully mutated PIK3CA constructs generated through the acoustic arraying of interchangeable modular pieces at nanoliter volumes alludes to a large scale increased-throughput DNA assembly workflow due to reduced reaction cost and DNA requirements.

Several other process improvements used in this study would help with an increase in the DNA assembly pipeline at larger scales. One such process improvement demonstrated in this experiment is the ability to acoustically plate the transformants onto solid media. OmniTray plating with the Echo 525 Liquid Handler can save operator time and incubator space as compared to traditional transformation plating. The Echo 525 Liquid Handler has technology that is uniquely positioned to make this plating possible. Tip-based systems have difficulty with poured agar plates being inherently concave and variable in height. The Echo 525 Liquid Handler is capable of transferring 25 nL increments of cell suspension very precisely onto the solid media on an X, Y grid, but with a buffered Z height due to the acoustic, tipless transfer of the droplet itself. This allows the transfer of cells onto solid media with minimal impact from the media height.

Another major process improvement was seen using colony qPCR for the transformed NEBuilder reactions. One of the major difficulties with any DNA construction technique is the removal of false positives, which can exceed 50% of clones upon Next-Generation Sequencing. Background vector is especially problematic whether it is re-circularized or uncut from the initial preparation. Using colony qPCR on the Echo 525 Liquid Handler, the background colonies can be removed prior to the time-consuming plasmid recovery step. The colony qPCR is advantageous for scale-up when compared to end-point colony PCR due to the larger number of screened clones possible due to the removal of the laborious, low-throughput use of agarose gels. The quantitative results provided by the colony qPCR also remove the subjective nature of agarose gel reading. The colony qPCR was able to correctly predict 100% of vector background present, as well as 74% of the insertion deletion constructs from 192 samples sequenced allowing for a vastly enriched sequencing sample set. This process will be further optimized in future work. Finally, the sequencing of the passed plasmids can be done in a cost-effective way using a miniaturized Nextera XT protocol in conjunction with the Illumina MiSeq. Prior work had demonstrated the ability to miniaturize the Nextera XT kit ten to one hundredfold. In this experiment, tenfold Nextera XT reaction volumes were confirmed to also work for plasmid validation. Acoustic liquid transfers with the Echo 525 Liquid Handler allows the prospective user to decrease the overall reagent needs for NGS plasmid validation. Beyond overall DNA assembly reaction volume reduction, use of the Echo 525 Liquid Handler dramatically improves the DNA construction pipeline.

Working at the nanoliter scale in synthetic biology would have been inconceivable in the recent past, but the Echo Acoustic Technology makes these volumes a reality. Complex, modular DNA assemblies were generated using the NEBuilder HiFi kit at a forty-fold reduction in volume, which will significantly

decrease the cost of DNA assembly allowing for substantial increases in throughput. The integration of a Echo 525 Liquid Handler into a DNA construction pipeline, such as one using the NEB NEBuilder HiFi kit, can allow for significant volume reductions as well as workflow enhancements with both complex and simple plasmid designs.

## Materials

Equipment	Manufacturer	
Echo 525 Liquid Handler	Beckman Coulter Life Sciences	
Allegra X-14 Centrifuge	Beckman Coulter Life Sciences	
MixMate	Eppendorf	
TapeStation 2200	Agilent	
Incu-Shaker Mini, Shaking Incubator	Benchmark Scientific	
ProFlex PCR System	Thermo Fisher	
LightCycler 480	Roche	
PHERASTAR FS	BMG Labtech	
MiSeq	Illumina	
Reagents	Manufacturer	Part Number
DNA Assembly Cloning Kit	New England BioLabs	#E5520
KpnI-HF	New England BioLabs	#R3142
Q5 High-Fidelity 2X Master Mix	New England BioLabs	#M0492
QIAprep Spin Miniprep Kit	QIAGEN	#27104
Genomic DNA Analysis ScreenTape Kit	Agilent	#5067-5365, #5067-5366
NEB 10-beta Competent E. coli (High Efficiency)	New England BioLabs	#C2019I
T7 Express Competent E. coli (High Efficiency)	New England BioLabs	#C2566I
LB Broth with 100 µg/mL carbenicillin	Teknova	#L8185
LB Agar OmniTrays with 100 µg/mL carbenicillin	Teknova	#L2010
LB Agar OmniTrays with 100 µg/mL carbenicillin, 60µg/ml X-Gal, and 0.1mM IPTG	Teknova	#L2906
gBlocks Gene Fragments	Integrated DNA Technologies	Custom DNA
Junction qPCR primers	Integrated DNA Technologies	Custom Oligos
Piece amplification primers	Integrated DNA Technologies	Custom Oligos
LightCycler 480 SYBR Green I Master	Roche	#04707516001
pcDNA 3.1/Zeo (+) Mammalian Expression Vector	Thermo Fisher Scientific	#V86020
Monarch PCR & DNA Cleanup Kit (5 µg)	New England BioLabs	#T1030S
Nextera XT DNA 96-Sample Prep Kit	Illumina	#FC-131-1096
Nextera XT Index Kit v2 Set A,B,C,D	Illumina	#FC-131-2001
PhiX Control v3	Illumina	#FC-110-3001
MiSeq Reagent Kit v3 (150-cycle)	Illumina	#MS-102-3003
Quant-iT Picogreen dsDNA Assay Kit	Thermo Fisher Scientific	#P11496
Agencourt AMPure Beads	Beckman Coulter Life Sciences	#A63881
200 Proof Ethanol	Sigma Aldrich	#E7023

## Materials (continued)

Consumables	Manufacturer	Part Number
Echo Qualified 384-well PP Microplate	Beckman Coulter Life Sciences	#001-14555
Echo Qualified 384-well LDV Plus Microplate	Beckman Coulter Life Sciences	#001-12782
TapeStation Plate	Agilent	#5067-5150
14 mL Round Bottom Snap Cap Polypropylene Tubes	EK Scientific	#EK-62261
Hard-Shell Skirted 384-Well PCR Plate	Bio-Rad	#HSP3805
Hard-Shell 384-Well PCR Plates, Barcoded	Bio-Rad	#HSP3901
Microseal 'C' PCR Plate Sealing Film	Bio-Rad	#MSC1001
384 well plates, polystyrene black (with micro-clear bottom)	Greiner	#781096
Nalgene Single-Use PETG Erlenmeyer Flask (250 mL)	Thermo Fisher	#4112-0250
1.5 mL DNA LoBind Tubes	Eppendorf	#022431048

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.

Beckman Coulter Life Sciences 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. ECHO is a trademark or registered trademark of Labcyte 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-6623APP02.20