CLONING & SYNTHETIC BIOLOGY

Application Note

Improved method for assembly of linear yeast expression cassettes using NEBuilder® HiFi DNA Assembly Master Mix

Saulius Vainauskas, Ph.D. and Christopher H. Taron, Ph.D., New England Biolabs, Inc.

Introduction

Heterologous protein production in yeast expression systems (i.e., *Kluyveromyces lactis* and *Pichia pastoris*) normally involves insertion of a linear expression cassette into a target locus in the host genome (1-3). Typically, an expression cassette is assembled in *E. coli* by first cloning a gene of interest into a circular expression vector (Figure 1A). The expression cassette comprises DNA encoding a strong yeast promoter upstream of a heterologous gene of interest, a downstream transcription terminator sequence and a selectable marker gene. The entire cassette is flanked by locus-specific targeting sequences on either end (Figure 1B). The assembled vector is then amplified by propagation of the host *E. coli* cells, isolated by standard DNA preparation techniques, and subjected to restriction digestion to create a linear expression cassette. The linear fragment is then introduced into yeast whereby it integrates into a target locus on the host chromosome (Figure 1C).

While this approach for expression strain construction has been accepted methodology for over two decades, it does have some limitations. For example, expression vector assembly using traditional cloning techniques is a multi-day process. Additionally, one is limited to the use of restriction sites that are present in the expression vector and the use of the expression machinery (i.e., promoter, terminator, locus targeting sequence, etc.) that is built into the vector. Finally, for cloning of a gene whose product is highly toxic to *E. coli*, serendipitous protein expression during vector assembly in bacteria can yield clones having deleterious mutations (1,4). The use of *in vitro* DNA assembly methods to construct linear expression cassettes suitable for direct introduction into yeast circumvents each of these limitations while shaving days off of expression strain construction time. In the presented method, we highlight the use of NEBuilder HiFi DNA Assembly Master Mix (NEB #E2621) to rapidly assemble expression cassettes for use with NEB's *K. lactis* Protein Expression Kit (NEB #E1000).

DNA CLONING

DNA AMPLIFICATION & PCR
EPIGENETICS
RNA ANALYSIS
LIBRARY PREP FOR NEXT GEN SEQUENCING
PROTEIN EXPRESSION & ANALYSIS
CELLULAR ANALYSIS

Materials

- · NEBuilder HiFi DNA Assembly Master Mix
- Q5® Hot Start High-Fidelity 2X Master Mix
- · Overlapping Primers
- NEBuilder Assembly Tool
- K. lactis Protein Expression Kit
- pKLAC2
- K. lactis GG799 Competent Cells
- Yeast Carbon Base Medium
- Biolux® Gaussia Luciferase Assay Kit

References:

- Colussi, P.A. and Taron, C.H (2005) Kluyveromyces lactis LAC4 promoter variants that lack function in bacteria but retain full function in K. lactis. Appl. Environ. Microbiol. 71:7092–7098.
- van Ooyen, A.J.J., Dekker, P., Huang, M., Olsthoorn, M.M.A., Jacobs, D.I., Colussi P.A. and Taron, C.H. (2006) Heterologous protein production in the yeast Kluyveromyces lactis. FEMS Yeast Res. 6:381–392.
- Read, J.D., Colussi, P.A., Ganatra, M.B. and Taron, C.H. (2007) Acetamide selection of Kluyveromyces lactis cells transformed with an integrative vector leads to high frequency formation of multicopy strains. Appl. Environ. Microbiol. 73:5088–5096.
- Gibbs, M., Reeves, R.A., Sunna, A., and Bergquist, P.L. (2004) A yeast intron as a translational terminator in a plasmid shuttle vector. FEMS Yeast Res. 4:573–577.



A. HindIII 1 MCS SnaBl 8812 Bael 414 ame II LACA XcmI 508 CspCI 969 BsrGI 1010 BstXI 7532 SacII 7475 **BgIII** 1701 BssHII 1748 pKLAC2 PshAl 2225 9,107 bp Eco0109I - PpuMI 2536 Kasi - Nari - PluTi - Sfoi 2614 BstEII 3146 **BspEl** 3236 Ahdl 5929 Smal - TspMI - Xmal 3360 Bsu36I 3403 SacII 4648 PfIFI - Tth1111 4366 BmgBI 4257 BstXI 4319 В. SacII SacII Recombination Arm (5´-P_{LAC4}) Selectable Target 3'-PLAC4 homologous recombination K. lactis chromosome

Figure 1. Traditional yeast expression strain construction.

(A) The pKLAC2 integrative expression vector. A target gene is cloned into the multiple cloning site (optionally, in-frame with the mating factor alpha secretion leader for extracellular expression) using traditional techniques. Transcription is initiated and terminated by the LAC4 promoter ($P_{\text{LAC4-PB1}}$) and LAC4 transcription terminator (TT_{LAC4}) sequences, respectively. The Saccharomyces cerevisiae ADH1 promoter (P_{ADH1}) drives expression of a fungal acetamidase gene (amdS) for the selection of transformants by growth on acetamide-containing medium. E. coli vector sequence has been inserted into a unique SacII site in P_{LAC4} to allow for propagation in bacteria. (B) Targeted integration of a linear expression cassette. The vector is linearized by digestion with SacII or BstXI to permit subsequent insertion into the LAC4 promoter locus in the K. lactis genome. In the illustrated example, the fragment becomes inserted into the LAC4 promoter locus by recombination, mediated by homologous targeting sequences at either end of the fragment. The resulting chromosomal arrangement is shown in (C). Abbreviations: $3'-P_{\text{LAC4}}$ left targeting sequence; $5'-P_{\text{LAC4}}$ right targeting sequence; SL, secretion leader; TT_{LAC4} transcription terminator.

Selectable

Marker

LAC4

promoter

LAC4

LAC4 Promoter

Target

C.

promoter



Protocol

Experimental Design - Liner expression cassette assembly

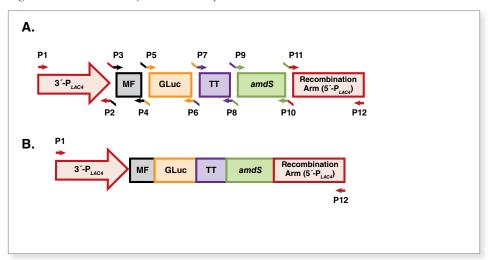
In this experiment, multiple pairs of primers were designed using the NEBuilder Assembly Tool (nebuilder.neb.com) to permit amplification of individual functional domains of an expression cassette (Table 1). Each resulting amplicon contained 22 bp overlaps with adjacent domains (Figure 2A). Purified PCR products were combined and treated with NEBuilder HiFi DNA Assembly Master Mix (NEB #E2621) and the fully assembled expression cassette was amplified using primers P1 and P12 (Figure 2B) to obtain sufficient quantities for introduction into *K. lactis* cells.

Table 1. Overlapping primers used for NEBuilder assembly of a linear *K. lactis* expression cassette.

PRIMER	SEQUENCE 5'→ 3'
P1	GATCGACTCATAAAATAGTAACC
P2	GAGAATTTCATTTTTTCAAGCTTCTCGATG
P3	GCTTGAAAAAAATGAAATTCTCTACTATATTAGCCG
P4	TCGGTGGGCTTTCTTTTCTCGAGATCATCC
P5	TCGAGAAAAGAAGCCCACCGAGAACAACG
P6	AAGGGCCTGTTTAGTCACCACCGGCCCC
P7	GTGGTGACTAAACAGGCCCCTTTTCCTTTG
P8	CACCCGGAAACAGCTTGCAAATTAAAGCCTTC
P9	ATTTGCAAGCTGTTTCCGGGTGTACAATATG
P10	ATCTGTTCCTTCTATGGAGTCACCACATTTC
P11	TGACTCCATAGAAGGAACAGATAGATAAAATTCCG
P12	CCGCGGAAA0TTTAGGAATTTTAAAC

^{*}Colors indicate DNA regions corresponding to the different functional domains shown in Figure 2A.

Figure 2. In vitro assembly of a linear expression cassette.



(A) Primer pairs P1 through P12 were used in PCR to amplify individual functional domains of the expression cassette.
(B) The domains were then assembled using the NEBuilder HiFi DNA Assembly Master Mix and the assembled linear expression cassette was amplified by PCR using primers P1 and P12. In this experiment, the CYC1 transcription terminator (TT) was used.



Fragment Preparation:

1. The following reaction conditions were used to amplify individual functional domains of the expression cassette (schematic in Figure 2A, primers in Table 1).

COMPONENT	VOLUME (μl)
Forward primer (40 µM)	1.25
Reverse primer (40 µM)	1.25
DNA template (5 ng/µl)	1
ddH_2O	46.5
Q5 Hot Start High-Fidelity 2X Master Mix	50
Total Volume	100

STEP	ТЕМР	TIME
Initial denaturation	98°C	30 seconds
30 Cycles	98°C	10 seconds
	58°C	30 seconds
	72°C	30 seconds or 1 minute
Final extension	72°C	7 minutes
Hold	4°C	∞

2. After thermocycling, each amplicon was purified by gel extraction.

Expression Cassette Assembly:

- The concentration of individual functional domain amplicons was determined by a Nanodrop™ instrument or estimated by agarose gel electrophoresis.
- 2. NEBuilder HiFi DNA Assembly Master Mix (NEB #E2621) was thawed at room temperature.
- 3. The DNA assembly reaction was set up us follows:
- 4. The reaction was incubated at 50°C for 1 hour.

COMPONENT	AMOUNT
PCR Amplicons	0.03 pmol (each)
2X NEBuilder HiFi DNA Assembly Master Mix	10 μΙ
ddH_2O	х µІ
Total volume	20 μΙ

- 5. A 2 μ l aliquot of the reaction was used as template in a fresh PCR reaction using primers P1 and P12 to amplify the assembled cassette. (Note: occasionally, more than one 100 μ l reaction is needed to generate enough pooled assembled product for introduction into yeast).
- 6. Following thermocycling, the amplicon was purified by gel extraction.

COMPONENT	VOLUME (μl)
Forward primer (40 µM)	1.25
Reverse primer (40 µM)	1.25
DNA Template (fragment assembly mix)	2
ddH_2O	45.5
Q5 High-Fidelity 2X Master Mix	50
Total Volume	100

STEP	ТЕМР	TIME
Initial denaturation	98°C	30 seconds
	98°C	10 seconds
30 Cycles	58°C	30 seconds
	72°C	3 minutes
Final extension	72°C	7 minutes
Hold	4°C	∞

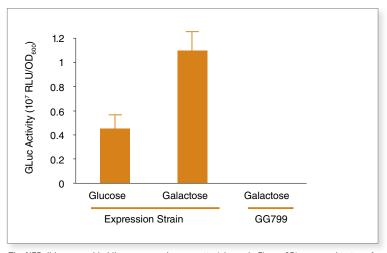
7. An aliquot containing 0.1–1 μg of the assembled linear expression cassette was used to transform *K. lactis* GG799 Competent Cells (NEB #C1001). The transformation reaction was plated on Yeast Carbon Base Medium (NEB #B9017) containing 5 mM acetamide and incubated for 3-4 days at 30°C until colonies formed. Targeted integration of the linear expression cassette into the *K. lactis* genome was confirmed by colony PCR using standard techniques described in the *K. lactis* Protein Expression Kit Manual (NEB #E1000).



Results:

A transformation efficiency of about 8 x 10³ colonies/µg DNA was obtained with the *in vitro* assembled expression construct, a similar efficiency to that observed when using a cloned expression cassette liberated from a pKLAC2 expression construct. *K. lactis* cells harboring at least one integrated copy of the *in vitro* assembled expression cassette successfully secreted *Gaussia princeps* luciferase (GLuc) into the growth medium (Figure 3).

Figure 3. Expression of Gaussia princeps luciferase (GLuc) in K. lactis GG799 cells.



The NEBuilder-assembled linear expression cassette (shown in Figure 2B) was used to transform K. lactis GG799 cells. A representative transformant was cultured in triplicate in rich medium (yeast extract/peptone) supplemented with glucose or galactose as a carbon source. Untransformed GG799 cells were grown in medium containing galactose as a negative control. GLuc enzyme activity secreted into the growth medium was measured using the BioLux Gaussia Luciferase Assay Kit (NEB #E3300). GLuc expression from the LAC4 promoter is repressed in the presence of glucose, but is de-repressed in galactose, as expected (1-3).

Conclusions:

In this application note, the use of NEBuilder HiFi DNA Assembly Master Mix (NEB #E2621) to rapidly assemble a linear expression fragment for introduction into *K. lactis* was demonstrated. The method permits fast preparation of linear expression cassettes in 1-2 days whereas the classical *E. coli* cloning approach takes at least 4-5 days. The *in vitro* assembly method is also suitable for the cloning of toxic genes because it eliminates the potential for accumulation of deleterious mutations during expression vector construction steps in bacteria. Moreover, NEBuilder HiFi DNA assembly permits more flexibility in the creation of constructs having different promoters, signal sequences, reporter genes and integration sites compared to using pKLAC2 in the *K. lactis* Protein Expression Kit (NEB #E1000). The general methodology outlined in this application note can also be applied more widely to assemble linear DNAs containing expression machinery for introduction into other yeasts (e.g. *Pichia pastoris*), fungi, bacteria, plant cells or animal cells.

NEW ENGLAND BIOLABS®, NEB®, BIOLUX®, NEBUILDER® and Q5® are registered trademarks of New England Biolabs, Inc. NANODROP® is a trademark of Thermo Fisher Scientific. Inc.







