

Electroporation of YACs

Maren Bell and Robert Mortimer, Human Genome Center, Division of Cell and Molecular Biology, Lawrence Berkeley Laboratory, Berkeley, California 94720

Introduction

Transformation of yeast with yeast artificial chromosomes (YACs) has traditionally been performed by a PEG-spheroplast procedure.^{1,2} However, the procedure is complicated, often unreliable, and transformation efficiencies are relatively low when compared to more recently developed transformation methods. An additional concern is that the smaller YACs of any given ligation mixture are cloned preferentially, therefore prior size fractionation with sucrose gradients to enrich for larger molecules, or treatment with spermidine to improve transformation efficiency of larger molecules, is required to obtain reasonable efficiencies with larger YACs. Several electroporation protocols have recently been reported for the transformation of *Saccharomyces cerevisiae* with supercoiled plasmids.³⁻⁶ In an attempt to improve the currently used YAC cloning procedure, we investigated the electroporation of *S. cerevisiae* strain AB1380 (Mat a, *ade2-1*, *can1-100*, *lys2-1*, *trp1*, *ura3*, *his5*, *psi⁺*) with a 14.4 kb "mini-YAC". Initial experiments showed transformation efficiencies two-fold better than those reported in a previously published YAC electroporation protocol.⁷

Materials and Methods

The YAC used in our experiments was constructed by the insertion of a 4.6 kb lambda *Mlu*I fragment into the *Mlu*I-*SUP4* cloning site of pYAC-RC.⁸ *E. coli* was transformed by the resulting 16.1 kb plasmid, the DNA was isolated, purified over cesium chloride, and restriction digested with *Bam*HI, thereby freeing the telomeres and resulting in a 14.4 kb linear DNA molecule. The restriction-digested DNA was microdialyzed (Millipore VS, 0.025 μ m) against TE (10 mM Tris-HCl, 1 mM EDTA, pH 7.5) for 20 minutes, ethanol precipitated, and suspended in TE. The electroporation protocol used for our experiments was an adaptation of the one devised by Becker and Guarente.⁵ Exponentially growing cells were harvested by centrifugation, washed twice in water (4 °C), once in 1 M sorbitol (4 °C), and finally resuspended in a 2/3 pellet volume of 1 M sorbitol (4 °C). A 40 μ l aliquot of cells was gently mixed with 1-5 μ l of DNA (0.1 μ g) just before electroporation. A Gene Pulser[®] apparatus connected to the Pulse Controller accessory was used to pulse the sample in a chilled 0.2 cm cuvette, generating a pulse with a field strength of 7.5 kV/cm (1.5 kV), and a time constant (τ) of approximately 4.2 msec (25 μ F and 200 Ω). Ice cold 1 M sorbitol (1 ml) was

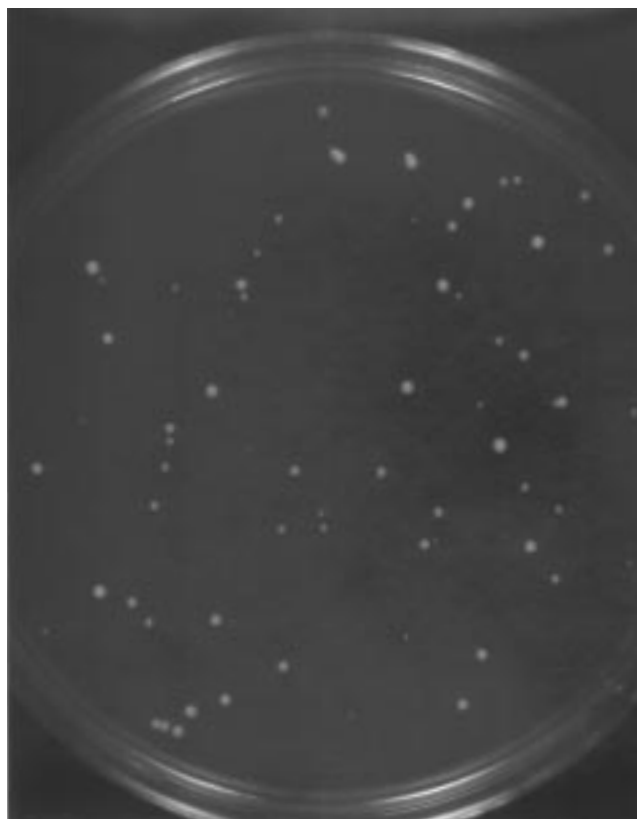


Fig. 1. Transformed yeast colonies. Shown after 3 days at 30 °C on medium lacking uracil.

added immediately after pulse delivery, and 150 μ l aliquots were plated on SD *ura⁻* plates. Transformants were visible after 3 days of incubation at 30 °C.

Results and Discussion

Transformations with the small YAC have yielded approximately 700 transformants / μ g of DNA. Transformations with a supercoiled 5 kb plasmid yielded up to 1.6×10^4 transformants / μ g of DNA. A number of parameters were investigated in attempt to increase efficiencies. We found that 0.1 μ g of DNA in 1-5 μ l of TE or water was the optimum amount for 40 μ l of cells (5×10^8 cells). More DNA had an adverse effect on transformation efficiency and less DNA did not seem to saturate the reaction (Figure 3). Spermidine-treated DNA did not yield any transformants. Exponentially growing cells gave the most consistent and highest transformation efficiencies.

A possible problem associated with the high voltage electroporation of very large DNA fragments such as YACs is that the molecules are most likely sheared before they have a chance to enter the cell. Lower voltage experiments with the Gene Pulser apparatus (which generates an exponential decay pulse) have a less disruptive effect on large DNA molecules, but they also yield lower transformation efficiencies.^{3,4} The use of alternative waveforms may improve on current results.⁶ The *S. cerevisiae* strain used here (AB1380) is traditionally employed in PEG-spheroplast transformations and we do not know whether this strain is also a good electroporation candidate. Other strains that have proven to yield high transformation rates might further improve the results presented here.

References

1. Burke, D. T., Carle, G. F. and Olson, M. V., *Science*, **236**, 806-812 (1987).
2. Burgers, P. M. J. and Percival, K. J., *Anal. Biochem.*, **163**, 391-397 (1987).
3. Delorme, E., *Appl. Environ. Microbiol.*, **55**, 2242-2246 (1989).
4. Simon, J. R. and McEntee, K., *Biochem. Biophys. Res. Comm.*, **164**, 1157-1164 (1990).
5. Becker, D. M. and Guarente, L., *Meth. Enzymol.*, **194**, 182-186 (1990).
6. Meilhoc, E., Masson, J.-M. and Teissié, J., *Bio/Technol.*, **8**, 223-227 (1990).
7. Rech, E. L., Dobson, M. J., Davey, M. R. and Mulligan, B. J., *Nucl. Acids Res.*, **18**, 1313 (1990).
8. Marchuck, D., and Collins, F. S., *Nucl. Acids Res.*, **16**, 7743 (1988).

Acknowledgements

This work was supported by Bio-Rad Laboratories and by funds administered through DOE Contract No. DE-ACO3-765F00098 to the Human Genome Center at Lawrence Berkeley Laboratory.

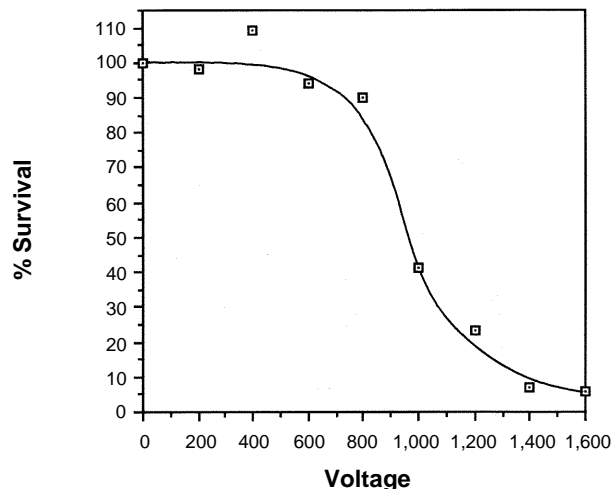


Fig. 2. Survival behavior of yeast strain AB1380. Exponentially growing cells were harvested and concentrated to 5×10^8 cells/40 μ l. Aliquots of 40 μ l were electroporated in a 0.2 cm cuvette at 25 μ F, 200 Ω , and 200 V to 1,600 V (in 200 V increments).

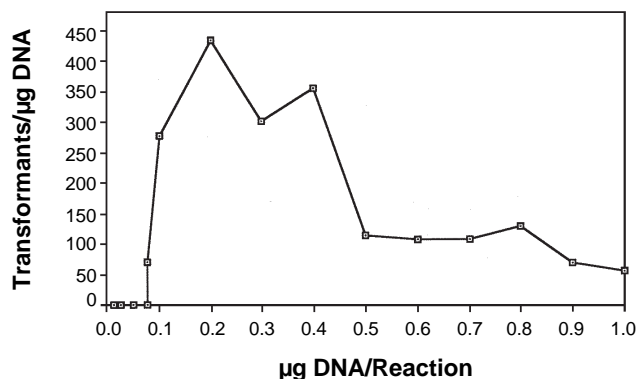


Fig. 3. Transformation efficiencies of strain AB1380 with a 14.4 kb YAC. Shown are results from several experiments performed under the same conditions: 40 μ l aliquots (5×10^8 cells in 1 M sorbitol) with 1-5 μ l of DNA were electroporated at 4 $^{\circ}$ C, 1500 V, and $\tau \sim 4.2$ msec (25 μ F and 200 Ω).

BIO-RAD

Bio-Rad
Laboratories

Life Science
Group

Bio-Rad Laboratories Main Office, 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-5800
 Also in: Reagents Park, Australia, Ph. 02-9914-2800, Fx. 02-9914-2888 Wien, Austria, Ph. (1) 877 89 01, Fx. (1) 876 56 29 Nazareth, Belgium, Ph. 09-385 55 11, Fx. 09-385 65 54
 Mississauga, Canada, Ph. (905) 712-2771, Fx. (905) 712-2990 Beijing, China, Ph. (01) 2046622, Fx. (01) 2051876 Copenhagen, Denmark, Ph. 39 17 9947, Fx. 39 27 1698
 Espoo, Finland, Ph. 90 804 2200, Fx. 90 804 1100 Ivry sur Seine Cedex, France, Ph. (1) 49 60 68 34, Fx. (1) 46 71 24 67 München, Germany, Ph. 089 31884-0, Fx. 089 31884-100
 New Delhi, India, Ph. 91-11-461-0103, Fx. 91-11-461-0765 Milano, Italy, Ph. 02-21609.1, Fx. 02-21609.399 Tokyo, Japan, Ph. 03-5811-6270 Fx. 03-5811-6272 Veendam,
 The Netherlands, Ph. 0318-540666, Fx. 0318-542216 Auckland, New Zealand, Ph. 09-443 3099, Fx. 09-443 3097 Kowloon, Hong Kong, Ph. 7893300, Fx. 7891257 Singapore,
 Ph. (65) 272-9877, Fx. (65) 273-4835 Solna, Sweden, Ph. 46 (0) 8 735 83 00, Fx. 46 (0) 735 54 60 Madrid, Spain, Ph. (91) 661 70 85, Fx. (91) 661 96 98 Glattbrugg, Switzerland,
 Ph. 01/809 55 55, Fx. 01/809 55 00 Hemel Hempstead, United Kingdom, Ph. 0800 181134, Fx. 01442 259118

SIG 020996 Printed in USA