

# Efficient Cloning and Electro-transformation of Large Eukaryotic DNA Fragments

Contributed by John M. Sedivy, Department of Molecular Biophysics and Biochemistry, Yale University, 333 Cedar Street, New Haven, CT 06510.

## Introduction

Genes of higher eukaryotes can be hundreds of kilobases in size. The need to rapidly clone and manipulate DNA molecules in that range has become acute in conjunction with the Human Genome Project. Cosmid vectors are constrained to 40–45 kb by the packaging limit of phage lambda. Recently, three systems have been developed that extend this range: yeast artificial chromosome (YAC)-based vectors (>500 kb limit),<sup>1</sup> phage P1-based vectors (100 kb limit),<sup>2</sup> and *E. coli* F factor-based vectors (limit as yet untested).<sup>3–5</sup> YAC clones have proven very successful, but are difficult to manipulate due to the extreme fragility of large linear DNA molecules. Bacterial episomes, being supercoiled, don't suffer that constraint. Unfortunately, F factor vectors have been limited by the low efficiency of transformation of large DNA molecules into *E. coli*. Recently, through the use of electroporation, efficiencies of  $10^{10}$  transformants/ $\mu\text{g}$  DNA of small plasmids have been achieved. A systematic analysis of large DNA molecules has been reported,<sup>5</sup> and is summarized here.

## Materials and Methods

*E. coli* strain JS4<sup>6</sup> was grown at 37 °C in LB broth; in mid-log phase ( $\text{OD}_{600}=0.6$ ) flasks were placed in ice water and chilled for 15 minutes. Cells were harvested, washed in ice-cold 1 mM HEPES, pH 7.0 (first wash 1x, second wash 0.5x of original culture volume), and resuspended at  $10^{11}$  cells/ml in 10% glycerol. 40  $\mu\text{l}$  aliquots were flash-frozen in dry ice/ethanol and stored at -70 °C. Cells were thawed on ice, DNA (2  $\mu\text{l}$  in TE) was added, and transferred to ice-cold 0.2 cm gap Bio-Rad electroporation cuvettes. A single pulse was delivered (Gene Pulser® II apparatus used with the Pulse Controller II) at 25  $\mu\text{F}$ , 2.50 kV and 200 ohms (time constant=4.5–4.7 msec). Cells were immediately resuspended in 1 ml of SOC (room temperature) and incubated at 37 °C for 45 minutes. Dilutions were plated on LB plates containing appropriate antibiotics. Plasmid DNA concentrations were determined by absorbance readings at  $\text{OD}_{260}$  and verified by gel electrophoresis.

## Results

We have investigated the relationship between plasmid size and electroporation efficiency in *E. coli* and found that even very large plasmids can be efficiently transfected. Six plasmids were used: pUC19 (2.6 kb),<sup>7</sup> cosmid tsA/Y/1d4 (44 kb),<sup>8</sup> and F factor clones pMBO52 (52 kb), pMBO74 (74 kb), pMBO99 (99 kb) and pMBO136 (136 kb).<sup>3</sup> The average efficiencies (number of colonies per  $\mu\text{g}$  DNA) were: pUC19,  $2 \times 10^9$ ; tsA/Y/1d4,  $6.1 \times 10^7$ ; pMBO52,  $6.6 \times 10^6$ ; pMBO74,  $3.2 \times 10^6$ ; pMBO99,  $2.7 \times 10^6$ ; pMBO136,  $1.7 \times 10^6$  (Figure 1). The data points can be fitted to a straight line when both axes are plotted on a logarithmic scale. The relation can be written as  $E=1.2 \times 10^6 M^{-0.4248}$ , where E is efficiency (colonies/fmole), and M is plasmid size (kb).

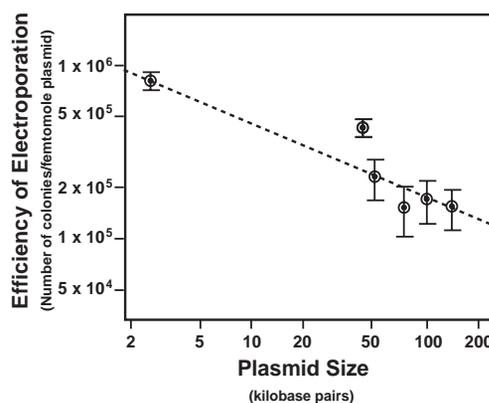
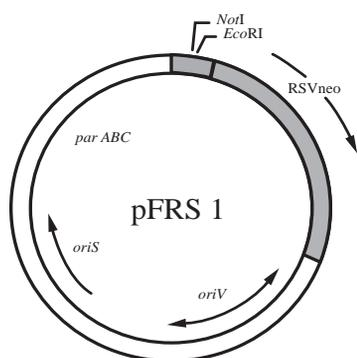


Fig. 1. Molar efficiency of electroporation as a function of plasmid size. Antibiotic selection was: pUC19 and tsA/Y/1d4, ampicillin resistance; pMBO52–pMBO136, chloramphenicol resistance. The efficiency of plating on chloramphenicol was 20–30% of that on ampicillin. Error bars are the standard deviations of six independent electroporation experiments.

Supercoiled and relaxed episomes were electroporated with equal efficiency. Rearrangements incurred during electroporation were analyzed at the level of restriction digestion; none were detected. To exploit these observations we have constructed a novel mammalian-*E. coli* shuttle vector, designated pFRS1 (**F** Replicon Shuttle, Figure 2), which contains three functional units: 1) the *E. coli* F factor origin of replication, partitioning, and copy number control region, 2) a marker selectable both in *E. coli* and mammalian cells (the neo cassette from plasmid pRSVneo), and 3) the polylinker from the pBluescript SK<sup>™</sup> cloning vector. The vector stably confers resistance to kanamycin in *E. coli*, and resistance to G418 in mammalian cells. The plasmid is approximately 12 kb in size, and, like its F factor parent, is maintained at low copy number.



**Fig. 2. The pFRS1 vector.** The plasmid was assembled from three fragments: 1) a 393 bp fragment from pBluescript KS, *PvuII* (coordinate 525 nt), to *Asel* (coordinate 918 nt); 2) a 2,925 bp fragment from pRSVneoDRI (pRSVneo which had the *EcoRI* site in the LTR, coordinate 89 nt, removed by filling in), *NdeI* (coordinate 582 nt), to *BamHI* (coordinate 3,397 nt); 3) an 8,717 bp fragment from the *E. coli* F sex factor plasmid, *HindIII* (coordinate 41.1 F map units, or 847 nt in the sequence of the *EcoRI* f5 fragment, GenBank accession M12987) filled in and *BclI* linked, to *EcoRI* (coordinate 49.9 F map units, or 9,564 nt, as above) blunted with mung bean nuclease.

## Discussion

The new vector system should accept inserts well in excess of 100 kb. The efficiencies of electroporation are well above the minimum required to construct representative libraries of complex eukaryotic genomes. At this time, the only vector system available for the cloning of very large DNA fragments uses YAC vectors. *E. coli* has several advantages over yeast as an organism for the passive propagation of foreign DNA. *E. coli* grows faster, DNA can be rapidly and efficiently extracted, and yields, even of single copy plasmids, are significantly higher. In addition, circular supercoiled DNA molecules are physically quite stable and easy to handle. The most crucial issue, however, is the genetic stability of foreign DNA. Use of *E. coli* offers the distinct advantage of the very considerable body of knowledge concerning its recombination pathways. In several instances the instability of foreign DNA in *E. coli* has been overcome by the judicious choice of appropriate, multiply marked, recombination-deficient strains.<sup>9-11</sup>

## References

- Burke, D. T., Carle, G. F., and Olson, M. V., Cloning large segments of exogenous DNA into yeast by means of artificial chromosome vectors, *Science*, **236**, 806-812 (1987).
- Sternberg, N., Bacteriophage P1 cloning system for the isolation, amplification, and recovery of DNA fragments as large as 100 kilobase pairs, *Proc. Natl. Acad. Sci. USA*, **87**, 103-107 (1990).
- O'Connor, M., Peifer, M., and Bender, W., Construction of large DNA segments in *Escherichia coli*, *Science*, **244**, 1307-1312 (1989).
- Hosoda, F., Nishimura, S., Uchida, H. and Ohki, M., An F factor based cloning system for large DNA fragments, *Nucl. Acids Res.*, **18**, 3863-3869 (1990).
- Leonardo, E. D. and Sedivy, J. M., A new vector for cloning large eukaryotic DNA segments in *Escherichia coli*, *Bio/Technol.*, **8**, 841-844 (1990).
- Sedivy, J. M., Capone, J. P., RajBhandary, U. L., and Sharp, P. A., An inducible mammalian amber suppressor: propagation of a poliovirus mutant, *Cell*, **50**, 379-389 (1987).
- Yanisch-Perron, C., Vieira, J., and Messing, J., Improved M13 phage cloning vectors and host strains: nucleotide sequences of the M13mp18 and pUC19 vectors, *Gene*, **33**, 103-119 (1985).
- Schnipper, L. E., Chan, V., Sedivy, J. M., Jat, P. S., and Sharp, P. A., Gene activation by induced DNA rearrangements, *Cancer Res.*, **49**, 6640-6644 (1989).
- Wyman, A. R., Wolfe, L. B., and Botstein, D., Propagation of some human DNA sequences in bacteriophage 1 vectors requires mutant *Escherichia coli* hosts, *Proc. Natl. Acad. Sci. USA*, **82**, 2880-2884 (1985).
- Ishiura, M., Hazumi, N., Koide, T., Uchida, T., and Okada, Y., A *recB recC sbcB recJ* host prevents *recA*-independent deletions in recombinant cosmid DNA propagated in *Escherichia coli*, *J. Bacteriol.*, **171**, 1068-1074 (1989).
- Leach, D., and Lindsey, J., In vivo loss of supercoiled DNA carrying a palindromic sequence, *Molec. Gen. Genet.*, **204**, 322-327 (1986).

pBluescript is a registered trademark of Stratagene Cloning Systems.



**Bio-Rad**  
Laboratories

Life Science  
Group

Website [www.bio-rad.com](http://www.bio-rad.com) U.S. (800) 451-0894 Australia (02) 99 74-2800 Austria (7) 877 81 01 Belgium (03) 385 55 77 Canada (800) 772-2777 China (86-10) 2048822 Denmark (33) 7799 47 Finland (90) 804 2200 France (7) 43 90 45 90 Germany (089) 37804-0 Hong Kong (7533) 300 India (91-11) 467-0703 Israel (03) 957 4727 Italy (02) 2 7609 7 Japan (03) 6877-02 70 The Netherlands (0373) 76 640666 New Zealand (09) 443 3000 Singapore (65) 2 72-9877 Spain (91) 687 70 85 Sweden (46) (0) 8 62 7 50 00 Switzerland (01) 809 55 55 United Kingdom (0200) 787 734