

Cell Type Bacterial, gram negative

Molecules DNA: pUC 19, about 2.7 kB; pLP116 (a pUC 19 derivative), about 2.8 kB

Species *E. coli*; *Legionella pneumophila*,
Used strain Nottingham N -7

Before the Pulse

Cell Growth Medium Buffered charcoal yeast extract (BCYE- α) agar supplemented with L-cysteineGrowth Phase at Harvest O.D. (600) =18 hr. growth on BCYE- α agar

Pre-pulse Incubation on ice for 30 min.

Wash Solution Phosphate buffered saline (PBS)

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature 0 °C

Electroporation Medium Electroporation buffer described by Dower: see notes

Cuvette Gap 0.4 cm

Cell Density 10 (9) cells / ml

Voltage 2.5 kV

Volume of Cells 800 μ l

Field Strength 6.25 kV/cm

DNA Concentration 5 μ g/ml

DNA Resuspension Buffer TE buffer (40 mM Tris, 2 mM EDTA (disodium). pH 7.9)

Capacitor 25 μ FVolume of DNA 2 to 10 μ l

Resistor Pulse Controller not used**. NOT

After the Pulse

Time Constant 4.5 to 4.8 msec

Outgrowth Medium BCYE - α , supplemented with L- cysteine agar

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. Electroporation medium is buffer described by Dower: 270 mM sucrose, 1 mM MgCl₂, 7 mM NaPO₄, pH 7.4, filter-sterilized.**Ref:** American Society for Microbiology Annual Meeting, Dallas, Texas, 1991, Abstract H-9.

**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.

Ref: Dower,W.J., Miller,J.F. and Ragsdale,C. 1988. *NAR* 16(13):6127-6145.

Outgrowth Temperature 37 °C

Length of Incubation 10 days

Selection Method or Assay Used 50 μ g / ml ampicillinElectroporation Efficiency 10 (6) transformants / μ g DNA

Per Cent Survival 60 %

Name of Submitter A. S. High and F. G. Rodgers - Professor

Institution Address University of New Hampshire
Microbiology Dept.
Spaulding Life Science Building
Durham, NH 03824

Survey Number

050