



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram negative
Species Used *E. coli*; *Legionella pneumophila*, strain Nottingham N -7

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

Before the Pulse

Cell growth medium Buffered charcoal yeast extract (BCYE- α) agar supplemented with L-cysteine
Wash solution Phosphate buffered saline (PBS)

Growth phase at harvest O.D. (600) =18 hr. growth on BCYE- α agar
Pre-pulse incubation on ice for 30 min.

The Pulse

Electroporation Temperature 0 °C
Electroporation Medium Electroporation buffer described by Dower: see notes
Cell Density 10 (9) cells / ml
Volume of Cells 800 μ l
DNA Concentration 5 μ g/ml
DNA Resuspension Buffer TE buffer (40 mM Tris, 2 mM EDTA (disodium), pH 7.9)
Volume of DNA 2 to 10 μ l

Instruments Used Gene Pulser® apparatus
Cuvette Gap 0.4 cm
Voltage 2.5 kV
Field Strength 6.25 kV/cm
Capacitor 25 μ F
Resistor Pulse Controller not used**. NOT RECOMMENDED (see notes).
Time Constant 4.5 to 4.8 msec

After the Pulse

Outgrowth Medium BCYE - α , supplemented with L- cysteine agar
Outgrowth Temperature 37 °C
Length of Incubation 10 days
Selection Method or Assay Used 50 μ g / ml ampicillin
Electroporation Efficiency 10 (6) transformants / μ g DNA
Per Cent Survival 60 %

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.

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 5/16/91

Survey Number 050

© Bio-Rad Laboratories, 1993