



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram negative
Species Used *E. coli*, N99, DH 5 α

Molecules Electroporated DNA, various supercoiled plasmids - CsCl pure; Qigen column, after microdialysis

Before the Pulse

Cell growth medium LB

Growth phase at harvest O.D. (600) = whatever

Pre-pulse incubation Short as possible; however long it takes to set up.

Wash solution Twice in cold water; twice in 10% glycerol

The Pulse

Electroporation Temperature Not given
Electroporation Medium 10% glycerol

Instruments Used Gene Pulser® apparatus
Pulse Controller

Cell Density Not given

Cuvette Gap 0.2 cm

Volume of Cells Not given

Voltage 2.5 kV

DNA Concentration Not given

Field Strength 12.5 kV/cm

DNA Resuspension Buffer TE (10 mM Tris, 1 mM EDTA, pH 8.0)

Capacitor 25 μ F

Volume of DNA 1 to 10 μ l

Resistor 200 Ω (Pulse Controller)

Time Constant 4.0 to 5.0 msec

After the Pulse

Outgrowth Medium SOC (Add glucose after autoclaving and cooling)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. I have electroperated temperature-sensitive mu-lysogens with no problems. I was concerned, because even a slight increase in temperature will result in cell death. As long as I keep things on ice and work quickly, there seems to be no problem with temperature induction by the electroperation.

Outgrowth Temperature 30 °C (N99) 37 °C (DH5 α)

Length of Incubation 4hr (N99); 2hr (DH5 α)

Selection Method or Assay Used Ampicillin, kanamycin resistance

SOC: 2% Bacto tryptone, 0.5% Bacto yeast extract, 10mM NaCl, 2.5mM KCl, 10 mM MgCl₂, 10 mM MgSO₄, 20 mM glucose.

LB: 1% Bacto tryptone, 0.5% Bacto yeast extract, 0.5% NaCl.

Electroporation Efficiency 2 to 4 x 10⁽⁸⁾ transformants / μ g DNA

Per Cent Survival Not given

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 8/27/90

Survey Number 041

© Bio-Rad Laboratories, 1993