



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram negative

Species Used *E. coli*, MC1061, modified to contain F' (host for M13)

Molecules Electroported DNA: M13 recombinant RF (double stranded) and single-stranded DNA

Before the Pulse

Cell growth medium LB

Growth phase at harvest O.D. (600) = 0.4 - 0.6

Pre-pulse incubation none; done as fast as possible

Wash solution 10% glycerol, 4 washes at 4°C

The Pulse

Electroporation Temperature 0 to 4 °C

Instruments Used Gene Pulser® apparatus
Pulse Controller

Electroporation Medium 10% glycerol

Cell Density As high as possible: just vortexed drained pellet.

Cuvette Gap 0.2 cm

Volume of Cells 40 µl

Voltage 2.450 kV

Field Strength 12.2 kV/cm

DNA Concentration 7.5 µg/µl

Capacitor 25 µF

DNA Resuspension Buffer 10% glycerol

Resistor 400 Ω (Pulse Controller)

Volume of DNA 5 µl + 40µl cells

Time Constant 9.1 to 9.6 msec

After the Pulse

Outgrowth Medium SOC

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 25 °C

Transformation efficiency for M13 is about 3x as high as pUC plasmid.

Length of Incubation 1 to 10 min., plated immediately

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

Selection Method or Assay Used M13 plaques, no drugs

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

Electroporation Efficiency 1-5x10⁽¹⁰⁾ (RF); 1-50x10⁽⁸⁾ (single-stranded) transformants / µg DNA

Per Cent Survival Not measured

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 10/16/90

Survey Number 036

© Bio-Rad Laboratories, 1993