



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram negative

Species Used *E. coli*, DH5α

Molecules Electroported DNA: plasmid, pUC13 & subclones; pCDNA-1 (CDM8 variant)

Before the Pulse

Cell growth medium LB; made w/1% BACTO-tryptone, 0.5% Bacto-yeast; no NaCl added, no pH adjustment.

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

Pre-pulse incubation None

Wash solution water; best available; 4°C

The Pulse

Electroporation Temperature 25 °C

Instruments Used Gene Pulser® apparatus and Pulse Controller

Electroporation Medium water

Cell Density 2 x 10⁽¹⁰⁾ cells / ml * (see notes)

Cuvette Gap 0.2 cm

Volume of Cells 0.06 to 0.1 ml

Voltage 1.75 kV

DNA Concentration 1pg DNA

Field Strength 8.75 kV/cm

DNA Resuspension Buffer SOC

Capacitor 25 μF

Volume of DNA 1 μl

Resistor 200 Ω (Pulse Controller)

Time Constant 4.5 to 5.0 msec

After the Pulse

Outgrowth Medium SOC (1 ml added immediately)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
* prepare by resuspending pellet in 2 volumes of deionized water.

Outgrowth Temperature 37 °C

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

Length of Incubation 1 hour

Selection Method or Assay Used Ampicillin

Electroporation Efficiency 1 x 10⁽⁹⁾ transformants / μg DNA

Per Cent Survival about 25%

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 3/11/91

Survey Number 019

© Bio-Rad Laboratories, 1993