



## Gene Pulser® Electroprotocols

**Cell Type** Bacterial, gram negative  
**Species Used** *E. coli*, unspecified strain

**Molecules Electroported** DNA

### Before the Pulse

**Cell growth medium** LB

**Growth phase at harvest** O.D. (600) = 0.25

**Pre-pulse incubation** 1 min.

**Wash solution** water

### The Pulse

**Electroporation Temperature** 4 °C

**Instruments Used** Gene Pulser® apparatus  
Pulse Controller

**Electroporation Medium** water

**Cell Density** 10 (10) cells / pulse

**Cuvette Gap** 0.2 cm

**Volume of Cells** 40 µl

**Voltage** 2.5 kV

**DNA Concentration** 1 ng / µl

**Field Strength** 12.5 kV/cm

**DNA Resuspension Buffer** TE buffer

**Capacitor** 25 µF

**Volume of DNA** 1 µl

**Resistor** 300 Ω (Pulse Controller)

**Time Constant** 4.7 msec

### After the Pulse

**Outgrowth Medium** Not given

#### Relevant Publications and/or Comments

**Note:** exponential values designated in parentheses.

**Outgrowth Temperature** 37 °C

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

**Length of Incubation** 1 hour

**Selection Method or Assay Used** Ampicillin

**Electroporation Efficiency** Not given

**Per Cent Survival** Not given

**Name of Submitter**  
**Institution Address**

**Telephone Number**

**Fax Number**

**Date Submitted** 3/5/91

**Survey Number** 045

© Bio-Rad Laboratories, 1993