



## Gene Pulser® Electroprotocols

**Cell Type** Bacterial, gram negative  
**Species Used** *E. coli*, DH5 $\alpha$

**Molecules Electroporated** DNA: plasmid from 2kb to 30 kb from a ligation mix, sometimes supercoiled.

### Before the Pulse

**Cell growth medium** LB

**Growth phase at harvest** O.D. (600) = between 0.4 and 0.6

**Pre-pulse incubation** 30 seconds to 1 minute

**Wash solution** double distilled water (ice cold) and (0°C, ice cold) glycerol

### The Pulse

**Electroporation Temperature** 0°C

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

**Electroporation Medium** 10% glycerol/water

**Cell Density** Not given

**Cuvette Gap** 0.2 cm

**Volume of Cells** 40  $\mu$ l

**Voltage** 2.5 kV

**Field Strength** 12.5 kV/cm

**DNA Concentration** See notes

**Capacitor** 25  $\mu$ F

**DNA Resuspension Buffer** Not given

**Resistor** 200  $\Omega$  (Pulse Controller)

**Volume of DNA** from 0.5  $\mu$ l to 2  $\mu$ l

**Time Constant** Not given

### After the Pulse

**Outgrowth Medium** SOC

#### Relevant Publications and/or Comments

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

**Outgrowth Temperature** 37°C

**DNA concentration:** Depends on ligation-usually straight out of gel slice (low melt) ligation.

**Length of Incubation** 20 min. to 1 hour

**SOC:** 2% Bacto tryptone, 0.5% Bacto yeast extract, 10mM NaCl, 2.5mM KCl, 10 mM MgCl<sub>2</sub>, 10 mM MgSO<sub>4</sub>,

**Selection Method or Assay Used** Have used carpenicillin, kanamycin, tetracycline, Xgal / IPTG (depending on plasmid)

20 mM glucose.

**Electroporation Efficiency** 10 (9) transformants /  $\mu$ g DNA

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

**Per Cent Survival** Not given

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

**Telephone Number**

**Fax Number**

**Date Submitted** 3/4/91

**Survey Number** 020

© Bio-Rad Laboratories, 1993