



## Gene Pulser® Electroprotocols

**Cell Type** Bacterial, gram negative  
**Species Used** *E. coli*, MC1061, DH5 $\alpha$ , XL1- Blue, NM522, JM109, MV1190

**Molecules Electroporated** DNA: cDNA library in pBS-SKII+ (3 to 7 kB), Bluescript™ vectors (3 to 7 kB), (single stranded M13 phage DNA).

### Before the Pulse

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

**Growth phase at harvest** O.D. (600) = 0.9

**Pre-pulse incubation** 10% glycerol

**Wash solution** glass distilled water

### The Pulse

**Electroporation Temperature** 4 °C  
**Electroporation Medium** 10% glycerol

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

**Cell Density** 1 liter cells @ 0.9 O.D. (600) to 2 ml cell slurry

**Cuvette Gap** 0.2 cm

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

**Voltage** 2.5 kV

**Field Strength** 12.5 kV/cm

**DNA Concentration** Not given

**Capacitor** 25  $\mu$ F

**DNA Resuspension Buffer** glass distilled water

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

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

**Time Constant** 4.8 msec

### After the Pulse

**Outgrowth Medium** SOB

#### Relevant Publications and/or Comments

**Note:** exponential values designated in parentheses. I've found that I can increase my transformation efficiency by 10 - 50 fold by being ultra careful to use the purest reagents and water available, autoclaving everything that will come into contact with cells (and I mean everything - including pipet tips and microfuge tubes) in glass distilled water prior to sterilization, and by using glassware that has never been washed with soap or bleach. Also, I've found that it is important to work quickly while preparing cells and to keep them cold. **SOB:** 2% Bacto tryptone, 0.5% Bacto yeast extract, 10mM NaCl, 2.5mM KCl, 20 mM glucose.

**Outgrowth Temperature** 37 °C

**Length of Incubation** 1 hr. 225 rpm

**Selection Method or Assay Used** LB-carbampicillin 100  $\mu$ g /  $\mu$ l

**Electroporation Efficiency** Strains without F: 10 (9) to 5 x 10 (10), with F: 10 (8) to 10 (9) transformants /  $\mu$ g

**Per Cent Survival** not given

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

**Telephone Number**

**Fax Number**

**Date Submitted** 8/21/90

**Survey Number** 035

© Bio-Rad Laboratories, 1993