



## Gene Pulser® Electroprotocols

**Cell Type** Fungal / Yeast

**Species Used** *Dictyostelium discoideum*

**Molecules Electroporated** DNA: supercoiled and linear

### Before the Pulse

**Cell growth medium** HL5 (peptone, yeast extract, glucose)

(ATCC Media #671)

**Growth phase at harvest**  $1 \times 10^{(6)}$  to  $1 \times 10^{(7)}$  cells / ml

**Pre-pulse incubation** 5 min.

**Wash solution** Heps Buffered Saline

### The Pulse

**Electroporation Temperature** 4 °C

**Electroporation Medium**

Heps Buffered Saline

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

**Cell Density**  $5 \times 10^{(6)}$  / ml

**Cuvette Gap** 0.4 cm

**Volume of Cells** 1 ml

**Voltage** 1.25 kV

**DNA Concentration** 20 ng to 20 µg

**Field Strength** 3.125 kV/cm

**DNA Resuspension Buffer** TE (10 mM Tris, 1 mM EDTA, pH 8.0)

**Capacitor** 25 µF

**Volume of DNA** Any volume

**Resistor** (Pulse Controller) Ω none. NOT RECOMMENDED\*\*\* (see notes)

**Time Constant** 0.5 to 0.7 msec

### After the Pulse

**Outgrowth Medium** HL5

#### Relevant Publications and/or Comments

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

**Outgrowth Temperature** 22 °C

\*\*It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.

**Length of Incubation** overnight

**Selection Method or Assay Used** G418

**HBS:** 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl<sub>2</sub>

**Electroporation Efficiency**  $10^{(-3)}$  transformants/ cell;  
 $6 \times 10^{(3)}$  transformants / µg DNA.

**Per Cent Survival** >90 %

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

**Telephone Number**

**Fax Number**

**Date Submitted** 11/13/91

**Survey Number** 175

© Bio-Rad Laboratories, 1993