



## Gene Pulser® Electroprotocols

**Cell Type** Other Cell Types

**Species Used** Chicken, TS34 a6 L1, [LSCC HD2], erythroblast

**Molecules Electroporated** DNA: supercoiled DNA used for transient transfections.

### Before the Pulse

**Cell growth medium** Not given

**Growth phase at harvest** 50 to 70% confluency

**Pre-pulse incubation** 4° C, 10 min.

**Wash solution** Wash two times in electroporation buffer

### The Pulse

**Electroporation Temperature** Room temperature  
**Electroporation Medium** Phosphate Buffered Saline

**Instruments Used** Gene Pulser® apparatus & Capacitance Extender

**Cell Density** 5 x 10<sup>(6)</sup> cells / pulse

**Cuvette Gap** 0.4 cm

**Volume of Cells** 0.5 ml

**Voltage** 0.270 kV

**DNA Concentration** 10 µg / pulse

**Field Strength** 0.675 kV/cm

**DNA Resuspension Buffer** Not given; pulse volume: 0.8 ml

**Capacitor** 960 µF

**Volume of DNA** Not given; pulse volume: 0.8 ml

**Resistor** (Pulse Controller) Ω not used

**Time Constant** 24 msec

### After the Pulse

**Outgrowth Medium** Not given

#### Relevant Publications and/or Comments

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

**PBS:** 1x = 8g NaCl, 0.2g KCl, 0.2g KH<sub>2</sub>PO<sub>4</sub>, 1.15g Na<sub>2</sub>HPO<sub>4</sub>

**Outgrowth Temperature** 37 °C

**Length of Incubation** 48 to 72 hrs.

**Selection Method or Assay Used** Transient assays

**Electroporation Efficiency** Not given

**Per Cent Survival** about 50 %

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

**Telephone Number**

**Fax Number**

**Date Submitted** 7/1/90

**Survey Number** 187

© Bio-Rad Laboratories, 1993