



# Gene Pulser® Electroprotocols

\* We recommend adapting this protocol to use the Gene Pulser electroporation buffer (catalog #165-2676, 165-2677), which increases cell viability and transfection efficiency in mammalian cell lines.

**Cell Type** Mammalian, adherent, suspension  
**Species Used** Mouse, NIH/3T3, embryo; Human T-cell line, (PEER)

**Molecules Electroported** DNA: pTS1-envlF, 11 kB, supercoiled

## Before the Pulse

**Cell growth medium** MDEM (GIBCO/BRL, Sigma)

**Growth phase at harvest** Not given

**Pre-pulse incubation** MDEM

**Wash solution** No

## The Pulse

**Electroporation Temperature** 25 °C  
**Electroporation Medium\*** MDEM

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

**Cell Density** 75%

**Cuvette Gap** 0.4 cm

**Volume of Cells** 0.8 ml

**Voltage** 0.270 kV

**DNA Concentration** 1 µg / µl

**Field Strength** 0.675 kV/cm

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

**Capacitor** 960 µF

**Volume of DNA** 35 µl

**Resistor** (Pulse Controller) Ω none

**Time Constant** 12.0 msec

## After the Pulse

**Outgrowth Medium** MDEM

### Relevant Publications and/or Comments

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

**Outgrowth Temperature** 25 °C

**Length of Incubation** 10 min.

**Selection Method or Assay Used** G418

**Electroporation Efficiency** Very good

**Per Cent Survival** 80 %

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

**Telephone Number**

**Fax Number**

**Date Submitted** 8/18/92

**Survey Number** 159