



# 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, suspension  
**Species Used** Human, Raji, Burkitt lymphoma; Jurkat, acute T cell leukemia; T-cells

**Molecules Electroporated** DNA: linear, 5 kB

## Before the Pulse

**Cell growth medium** RPMI 1640 + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)

**Growth phase at harvest** Not given

**Pre-pulse incubation** None

**Wash solution** Hepes Buffered Saline

## The Pulse

**Electroporation Temperature** 4°C  
**Electroporation Medium\*** Phosphate Buffered Saline

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

**Cell Density** 1 x 10<sup>(7)</sup> cells / pulse

**Cuvette Gap** 0.4 cm

**Volume of Cells** 0.5 to 0.75 ml

**Voltage** 0.40 kV

**DNA Concentration** Not given

**Field Strength** 1.0 kV/cm

**DNA Resuspension Buffer** Not given

**Capacitor** 960 µF

**Volume of DNA** 11 µl

**Resistor** (Pulse Controller) Ω none

**Time Constant** Not given

## 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>

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

**Outgrowth Temperature** 37°C

**Length of Incubation** 2 days

**Selection Method or Assay Used** Preparation of RNA or CAT assay

**Electroporation Efficiency** Not quantified

**Per Cent Survival** 50 %

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

**Telephone Number**

**Fax Number**

**Date Submitted** 8/17/90

**Survey Number** 107