



# 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.

|                  |   |                       |  |
|------------------|---|-----------------------|--|
| <b>Cell Type</b> | Mammalian, suspension   | <b>Molecules</b>      | DNA: pGL-luciferase vector   |
| <b>Species</b>   | Human, K562, chronic myeloid leukemia; HeLa, epithelial carcinoma; HEL cells, eythroleukemia. | <b>Electroporated</b> | [Promega]containing $\beta$ -globin promoter ; co-porated SV 40, $\beta$ -gal. |

---

## Before the Pulse

|                           |                                     |                                |                                       |
|---------------------------|-------------------------------------|--------------------------------|---------------------------------------|
| <b>Cell growth medium</b> | RPMI + 5% Fetal Calf Serum + 5% DCS | <b>Growth phase at harvest</b> | 2 to 5 x 10 <sup>(5)</sup> cells / ml |
| <b>Wash solution</b>      | Phosphate Bufferd Saline + 5% DCS   | <b>Pre-pulse incubation</b>    | 10 min., ice                          |

## The Pulse

|                                    |                                  |                         |   |
|------------------------------------|----------------------------------|-------------------------|---|
| <b>Electroporation Temperature</b> | 25°C                             | <b>Instruments Used</b> | Gene Pulser® apparatus & Capacitance Extender |
| <b>Electroporation Medium*</b>     | Hepes Buffered Saline            | <b>Cuvette Gap</b>      | 0.4 cm  |
| <b>Cell Density</b>                | 4 x 10 <sup>(7)</sup> cells / ml | <b>Voltage</b>          | 0.300 kV                                      |
| <b>Volume of Cells</b>             | 0.5 ml                           | <b>Field Strength</b>   | 0.75 kV/cm                                    |
| <b>DNA Concentration</b>           | 50 $\mu$ g per pulse             | <b>Capacitor</b>        | 960 $\mu$ F                                   |
| <b>DNA Resuspension Buffer</b>     | Not given                        | <b>Resistor</b>         | (Pulse Controller) $\Omega$ none              |
| <b>Volume of DNA</b>               | 50 $\mu$ l                       | <b>Time Constant</b>    | 31 msec                                       |

## After the Pulse

|                                       |                                     |
|---------------------------------------|-------------------------------------|
| <b>Outgrowth Medium</b>               | RPMI + 5% Fetal Calf Serum + 5% DCS |
| <b>Outgrowth Temperature</b>          | 37 °C                               |
| <b>Length of Incubation</b>           | 24 hours                            |
| <b>Selection Method or Assay Used</b> | luciferease, $\beta$ -gal           |
| <b>Electroporation Efficiency</b>     | Not given                           |
| <b>Per Cent Survival</b>              | 50 %                                |

### Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

---

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

**Telephone Number**  
**Fax Number**  
**Date Submitted** 8/26/92  
**Survey Number** 192