



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 Mouse, J558-L, myeloma

Molecules Electroporated DNA: plasmids, 4.5 kB, supercoiled.

Before the Pulse

| | | | |
|---------------------------|--|--------------------------------|---------------------------|
| Cell growth medium | DMEM + 10% Fetal Calf Serum (GIBCO/BRL, Sigma) | Growth phase at harvest | Split on the previous day |
| Wash solution | None | Pre-pulse incubation | 10 min on ice |

The Pulse

| | | | |
|------------------------------------|-----------------------------|-------------------------|---|
| Electroporation Temperature | 0 °C | Instruments Used | Gene Pulser® apparatus & Capacitance Extender |
| Electroporation Medium* | DMEM + 10% Fetal Calf Serum | Cuvette Gap | 0.4 cm |
| Cell Density | 10 (7) cells in 300 µl | Voltage | 0.25 kV |
| Volume of Cells | 300 µl | Field Strength | 0.625 kV/cm |
| DNA Concentration | 4 µg DNA per 10 (7) cells | Capacitor | 960 µF |
| DNA Resuspension Buffer | Not given | Resistor | (Pulse Controller) Ω none |
| Volume of DNA | 4 µl | Time Constant | 50 msec |

After the Pulse

| | | |
|---------------------------------------|-----------------------------|--|
| Outgrowth Medium | DMEM + 10% Fetal Calf Serum | Relevant Publications and/or Comments Note: exponential values designated in parentheses. <i>PNAS</i> , 87 : 5788 (1990). Electroporation worked best for J558L cells - DEAE dextran did not; neither did lipofection. |
| Outgrowth Temperature | 37 °C | |
| Length of Incubation | 48 hours | |
| Selection Method or Assay Used | CAT assay | |
| Electroporation Efficiency | Not given | |
| Per Cent Survival | 60 % | |

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 8/23/90
Survey Number 135
© Bio-Rad Laboratories, 1993