



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
Species Used Hamster, CHO, ovary

Molecules Electroporated DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.

Before the Pulse

Cell growth medium RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)

Growth phase at harvest Log phase (We routinely subdivide cells 24 hours prior to electroporation)

Pre-pulse incubation None

Wash solution Trypsin harvest, washed twice in Phosphate Buffered Saline

The Pulse

Electroporation Temperature 25 °C during and after
Electroporation Medium* RPMI 1640 without Fetal Calf Serum, +10mM dextrose, 0.1 mM dithiothreitol

Instruments Used Gene Pulser® apparatus & Capacitance Extender

Cell Density 1.3 x 10⁷ viable cells / ml, 0.3 ml

Cuvette Gap 0.4 cm

Volume of Cells 300 µl

Voltage 0.25 kV

DNA Concentration 500 µg / ml (5 to 10 µg per pulse)

Field Strength 0.625 kV/cm

DNA Resuspension Buffer Not given

Capacitor 960 µF

Volume of DNA 10 to 20 µl

Resistor (Pulse Controller) Ω none

Time Constant 33 to 38 msec

After the Pulse

Outgrowth Medium DMEM culture media, 10% fetal calf serum

Relevant Publications and/or Comments

Outgrowth Temperature 37 °C

Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.

Length of Incubation 48 hours

Selection Method or Assay Used Transient (CAT, β-gal, immunohisto-chemistry)

Electroporation Efficiency 50 to 100%

Per Cent Survival 20 to 75 %

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 3/1/91

Survey Number 091

© Bio-Rad Laboratories, 1993