



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 Hamster, CHO, ovary

Molecules Electroporated DNA: linearized plasmids, about 10 kB.

Before the Pulse

Cell growth medium Hams F12 (Gibco)

Growth phase at harvest 70 % confluence

Pre-pulse incubation 10 min. on ice

Wash solution Phosphate Buffered Saline, pH 7.4

The Pulse

Electroporation Temperature 0 to 4 °C

Instruments Used Gene Pulser® apparatus & Capacitance Extender

Electroporation Medium* Phosphate Buffered Saline (PBS), pH 7.4

Cuvette Gap 0.4 cm

Cell Density 5x10⁽⁶⁾ cells / 700 µl

Voltage 0.75 kV

Volume of Cells 700 µl

Field Strength 1.88 kV/cm

DNA Concentration 20 to 100 µg / 700 µl

Capacitor 25 µF

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

Resistor (Pulse Controller) Ω none

Volume of DNA 20 µl to 50 µl; (conc.=1µg / µl)

Time Constant 0.5 msec

After the Pulse

Outgrowth Medium Hams F12 (Gibco)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 37 °C

PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH₂PO₄, 1.15g Na₂HPO₄

Length of Incubation 2 days before selection

HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl₂

Selection Method or Assay Used F12 - 5158 plus or minus MTX

Electroporation Efficiency 250 to 1000 transformants / µg DNA

Per Cent Survival 25 %

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 5/7/92

Survey Number 088

© Bio-Rad Laboratories, 1993