



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 Human, squamous cell carcinoma lines, oral & cervical; Monkey, Vero, kidney cells.

Molecules Electroporated DNA: Recombinant pMAM neo vectors (8.8 to 9.2 kB) - confirm via expression (Northern blot) immunoblotting, PCR & Southern blot

Before the Pulse

Cell growth medium M199, MEM, DMEM/F12 (GIBCO/BRL, Sigma)
Wash solution Ice-cold Phosphate Buffered Saline, pH 7.2

Growth phase at harvest Mid-log phase
Pre-pulse incubation Ice, 10 min., in media with DNA.

The Pulse

Electroporation Temperature 4 °C
Electroporation Medium* Growth media
Cell Density 1 x 10⁽⁷⁾ cells / ml
Volume of Cells 3 x 10⁽⁶⁾ cells / pulse
DNA Concentration Not given
DNA Resuspension Buffer Not given
Volume of DNA 100 µl

Instruments Used Gene Pulser® apparatus & Capacitance Extender
Cuvette Gap 0.4 cm
Voltage 0.250 kV
Field Strength 0.625 kV/cm
Capacitor 500 µF
Resistor (Pulse Controller) none Ω
Time Constant 8 to 12 msec

After the Pulse

Outgrowth Medium Growth media, plus or minus G418
Outgrowth Temperature 37°C
Length of Incubation at least 10 days
Selection Method or Assay Used G418
Electroporation Efficiency 6 x 10⁽⁻⁵⁾ to 9 x 10⁽⁻⁸⁾; depends on cell line used
Per Cent Survival 50 %

Relevant Publications and/or Comments
Note: exponential values designated in parentheses.

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 3/5/91
Survey Number 162
© Bio-Rad Laboratories, 1993