



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	Molecules	DNA: neo gene.
Species Used	Human, MRC-5 / V1, lung fibroblasts, transformed	Electroporated	

Before the Pulse

Cell growth medium	DMEM + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth phase at harvest	Log phase
Wash solution	Phosphate Buffered Saline	Pre-pulse incubation	37° C, 1/2 hour

The Pulse

Electroporation Temperature	25°C	Instruments Used	Gene Pulser® apparatus
Electroporation Medium*	DMEM	Cuvette Gap	0.4 cm
Cell Density	5 x 10 ⁽⁶⁾ cells	Voltage	2.0 kV
Volume of Cells	1.4 ml (**see notes)	Field Strength	5.0 kV/cm
DNA Concentration	20 µg	Capacitor	25 µF
DNA Resuspension Buffer	water	Resistor	(Pulse Controller) Ω none; NOT RECOMMENDED*** (see notes)
Volume of DNA	20 µl	Time Constant	0.4 msec

After the Pulse

Outgrowth Medium	DMEM, 10% Fetal Calf Serum
Outgrowth Temperature	37 ° C
Length of Incubation	1 month for selection
Selection Method or Assay Used	1 mg / ml geneticin
Electroporation Efficiency	Not given
Per Cent Survival	Not given

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller. It is not recommended to use more than 0.8 ml in the 0.4 cm cuvette; greater volumes may create non-uniform field strengths during the pulse.

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 10/15/90
Survey Number 112
 © Bio-Rad Laboratories, 1993