



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 Electroporated	DNA: pBR322 derived plasmids containing retroviral vectors.
Species Used	Mouse, NIH/3T3 derived retroviral vector packaging cell lines		

Before the Pulse

Cell growth medium	DMEM + 10% Fetal Bovine Serum + 1% Glutamine (GIBCO/BRL, Sigma)	Growth phase at harvest	Log phase
Wash solution	Phosphate Buffered Saline	Pre-pulse incubation	5 min, with DNA at room temperature

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender
Electroporation Medium*	DMEM + 10% Fetal Bovine Serum + 1% Glutamine	Cuvette Gap	0.4 cm
Cell Density	1.5 x 10 ⁽⁶⁾ cells / ml	Voltage	0.2 kV
Volume of Cells	0.5 ml	Field Strength	0.5 kV/cm
DNA Concentration	0.1 to 1 µg / µl	Capacitor	960 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	10 µl	Time Constant	20 to 25 msec

After the Pulse

Outgrowth Medium	DMEM + 10% Fetal Bovine Serum + 1% Glutamine	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Outgrowth Temperature	37 °C		
Length of Incubation	Not given		
Selection Method or Assay Used	G418		
Electroporation Efficiency	50 to 200 transformants / µg DNA		
Per Cent Survival	Not given		

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 8/21/90
Survey Number 139
 © Bio-Rad Laboratories, 1993