



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	Molecules Electroporated	DNA: circular plasmid recombinants (pMAM (8.8 kB + insert), pATH (9.0 kB) pGem (4.0 kB) pMSG (8.8 kB) pM6 (9.7 kB))
Species Used	Human, HeLa, cervical (C41); Human epithelial cells; Monkey, Vero, kidney.		

Before the Pulse

Cell growth medium	DMEM, MEM, M199; (10% Fetal Bovine Serum + 1% antibiotic) (GIBCO/BRL, Sigma)	Growth phase at harvest	Early log
Wash solution	Phosphate Buffered Saline (cold)	Pre-pulse incubation	Ice, 10 min.

The Pulse

Electroporation Temperature	Chilled	Instruments Used	Gene Pulser® apparatus & Capacitance Extender, Pulse Controller
Electroporation Medium*	Same as growth media	Cuvette Gap	0.4 cm
Cell Density	3 x 10 ⁽⁶⁾ cells / ml	Voltage	0.250 kV
Volume of Cells	400 µl	Field Strength	0.625 kV/cm
DNA Concentration	10 to 100 µg / sample	Capacitor	500 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	1 µl to 50 µl	Time Constant	20 to 25 msec

After the Pulse

Outgrowth Medium	Same as growth media	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Outgrowth Temperature	Brief ice, then 37°C		
Length of Incubation	16 to 48 hours prior to selection		
Selection Method or Assay Used	G418		
Electroporation Efficiency	1 x 10 ⁽⁻³⁾ to 10 ⁽⁻⁹⁾ depending on cell type and plasmids used.		
Per Cent Survival	50%		

Name of Submitter	Telephone Number
Institution Address	Fax Number
	Date Submitted 11/11/91
	Survey Number 158
	© Bio-Rad Laboratories, 1993