



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: linear & supercoiled: Rsv neo, Rsv gal, myc CAT, IRF CAT, fos CAT, hsp CAT, etc.
Species Used	Monkey, COS, kidney cells; Rat, N62 T cells; Mouse, mammary epithelial cells		

Before the Pulse

Cell growth medium	Fischer's + 10% Horse Serum + 10% Fetal Calf Serum (GIBCO/BRL)	Growth phase at harvest	None
Wash solution		Pre-pulse incubation	0 to 10 min, room temperature

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender
Electroporation Medium*	Hepes buffered saline or Fischer's + 10% Horse Serum + 10% Fetal Calf Serum	Cuvette Gap	0.4 and 0.2 cm
Cell Density	1 x 10 ⁽⁷⁾ cells / 0.4 ml	Voltage	0.25, 0.35, 0.45 kV
Volume of Cells	0.4 ml to 0.8 ml	Field Strength	625, 875, 1125kV/cm
DNA Concentration	80 µg/0.3 ml Hepes Buffered Saline	Capacitor	125, 250, 960 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	< 50 µl	Time Constant	2 to 30 msec

After the Pulse

Outgrowth Medium	Fischer's + 10% Horse Serum + 10% Fetal Calf Serum	Relevant Publications and/or Comments
Outgrowth Temperature	Not given	Note: exponential values designated in parentheses.
Length of Incubation	Not given	PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄
Selection Method or Assay Used	G418, 400 µg / ml	HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Electroporation Efficiency	poor	
Per Cent Survival	40 to 50 %	

Name of Submitter	Telephone Number
Institution Address	Fax Number
	Date Submitted 3/6/91
	Survey Number 154
	© Bio-Rad Laboratories, 1993