



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	DNA: Plasmid DNA, 5 kB -12 kB
Species Used	Human, 293, kidney cells; Hamster, CHO, ovary cells	Electroporated	

Before the Pulse

Cell growth medium	Hamms F-12 / DMEM (50:50) +10% Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth phase at harvest	Log phase
Wash solution	Phosphate Buffered Saline	Pre-pulse incubation	Change to Hamms F-12 / DMEM (50:50) media 24 hours prior to pulse.

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender, Pulse Controller
Electroporation Medium*	High glucose/DMEM (Best we tested)	Cuvette Gap	0.4 cm
Cell Density	3 x 10 ⁽⁶⁾ cells / ml	Voltage	0.245 kV
Volume of Cells	800 µl	Field Strength	0.098 kV / cm
DNA Concentration	2 µg / ml	Capacitor	960 µF
DNA Resuspension Buffer	TE	Resistor	(Pulse Controller) Ω none
Volume of DNA	2 µl	Time Constant	Not given

After the Pulse

Outgrowth Medium	Hams F-12 / DMEM (50:50) media +10% Fetal Bovine Serum	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Outgrowth Temperature	37 °C		
Length of Incubation	20 min		We are currently testing amplification procedures with linearized plasmid [See: Barsou, M., <i>DNA and Cell Biology</i> v.9 (4): 293-300].
Selection Method or Assay Used	G418		PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄
Electroporation Efficiency	5%		HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Per Cent Survival	70 to 80%		

Name of Submitter	Telephone Number
Institution Address	Fax Number
	Date Submitted 8/16/90
	Survey Number 156
	© Bio-Rad Laboratories, 1993