



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, suspension	Molecules Electroporated	DNA: linearized DNA used for stable transfections.
Species Used	Mouse, SP-2, myeloma [Sp2/0-Ag14].		

Before the Pulse

Cell growth medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth phase at harvest	50 to 70% confluency
Wash solution	Wash two times in electroporation buffer	Pre-pulse incubation	4° C, 10 min.

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (5) cells/pulse, stable transfection	Voltage	0.180 kV
Volume of Cells	0.5 ml	Field Strength	0.45 kV/cm
DNA Concentration	10 µg / pulse	Capacitor	960 µF
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Resistor	(Pulse Controller) Ω none
Volume of DNA	Not given; pulse volume: 0.8 ml	Time Constant	24 msec

After the Pulse

Outgrowth Medium	DMEM, 10% Fetal Calf Serum (FCS)	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Outgrowth Temperature	37 °C	Note:	Stable transfections generally do not use carrier DNA. Also, the level of selective agent required to kill off non-transfected cells needs to be established before transfection. The level required should kill non-transfected cells in approximately 7 days.
Length of Incubation	48 to 72 hrs.		
Selection Method or Assay Used	G418 (stable transfections)		
Electroporation Efficiency	Not given		
Per Cent Survival	about 50 %		

Name of Submitter	Telephone Number
Institution Address	Fax Number
	Date Submitted 7/1/90
	Survey Number 142
	© Bio-Rad Laboratories, 1993