



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	Proteins
Species Used	Monkey, CV-1, kidney; Mouse, 3T3, embryo; Mouse, p3x63AG8, myeloma; Human: HeLa, epithelial carcinoma	Electroporated	

Before the Pulse

Cell growth medium	DMEM (GIBCO/BRL, Sigma)	Growth phase at harvest	50 to 60% confluent cells
Wash solution	(See notes)	Pre-pulse incubation	5 minutes

The Pulse

Electroporation Temperature	0 °C (See notes)	Instruments Used	Gene Pulser® apparatus
Electroporation Medium*		Cuvette Gap	0.4 cm
Cell Density	5 x 10 ⁽⁶⁾ cells / ml	Voltage	0.5 to 0.75 kV
Volume of Cells	0.2 µl	Field Strength	1.25 to 1.875 kV/cm
DNA Concentration	Not given	Capacitor	25 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	Not given	Time Constant	2 to 4 msec

After the Pulse

Outgrowth Medium	Pore-Resealing Buffer (see notes) for 10 min. ,37° C; follow with DMEM plus 10% Fetal Calf Serum.	Relevant Publications and/or Comments
Outgrowth Temperature	37° C	Note: exponential values designated in parentheses.
Length of Incubation	Not given	Ref: M.R. Mitchell et. al. (1988) <i>Experientia</i> 44 : 199-203.
Selection Method or Assay Used	Not given	M.R. Michel et. al. (1990) <i>J.Virol.</i> 64 : 5123-5131.
Electroporation Efficiency	Not given	Elgizoli, M. et. al. (1989) <i>J. Virol.</i> 63 : 2921-2928.
Per Cent Survival	Not given	Electroporation Buffer: 20 mM PIPES, pH 7; 128 mM K-glutamate, 5mM ATP, 5 mM GTP, 10 µm Ca-Acetate, 2 mM Mg-Acetate & amino acid concentration corresponding to that of DMEM media.
		Pore-Resealing Buffer: 20 mM PIPES, pH 7.0, 128 mM K-glutamate,10 µm Ca-Acetate, 2 mM Mg-Acetate

Name of Submitter	Telephone Number
Institution Address	Fax Number
	Date Submitted 3/8/91
	Survey Number 167
	© Bio-Rad Laboratories, 1993