



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, p220LTR, 11.2 kb, supercoiled.
Species Used	Human, 293s, transformed primary embryonic kidney; Lymphoblast cell lines, EBV immortalized		

Before the Pulse

Cell growth medium	RPMI, 10% Fetal Calf Serum, Penicillin, Streptomycin, L-Glutamine, or DMEM (GIBCO/BRL, Sigma)	Growth phase at harvest	80% confluent or 1 x 10 ⁽⁶⁾ cells / ml
Wash solution	Electroporation Buffer (see notes)	Pre-pulse incubation	None

The Pulse

Electroporation Temperature	25 °C	Instruments Used	Gene Pulser® apparatus & Capacitance Extender, Pulse Controller
Electroporation Medium*	See notes	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁽⁷⁾ cells / μl	Voltage	0.260 kV
Volume of Cells	800 μl	Field Strength	0.65 kV/cm
DNA Concentration	1 to 4 μg / μl	Capacitor	550 or 960 μF
DNA Resuspension Buffer	TE Buffer (10 mM Tris, 1 mM EDTA, pH 8.0)	Resistor	(Pulse Controller) Ω none
Volume of DNA	<20 μl	Time Constant	6.4 or 11.2 msec

After the Pulse

Outgrowth Medium	RPMI, 10% Fetal Calf Serum, Penicillin, Streptomycin, L-Glutamine, or DMEM	Relevant Publications and/or Comments
Outgrowth Temperature	37°C	Note: exponential values designated in parentheses.
Length of Incubation	Infinite	* Electroporation buffer from Chu, <i>et. al.</i> , <i>NAR</i> 15 (3): 1311 (1987); 1x HeBS, 20 mM HEPES, pH 7.05, 137 mM NaCl, 5 mM KCl, 0.7 mM Na ₂ HPO ₄ , 6 mM dextrose.
Selection Method or Assay Used	Hygromycin resistance	
Electroporation Efficiency	4000 transformants / μg DNA	
Per Cent Survival	10 %	

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 3/8/91
Survey Number 108
 © Bio-Rad Laboratories, 1993