



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	Molecules	DNA: supercoiled plasmid; luciferase reporter gene
Species Used	Rat, CA77, medullary thyroid carcinoma cell line	Electroporated	

Before the Pulse

Cell growth medium	DMEM/F12 (1:1) + 10 % Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth phase at harvest	Active growth, about 50 to 70 % confluent
Wash solution	Dulbecco's Phosphate Buffered Saline (minus Ca ⁺⁺ , Mg ⁺⁺)	Pre-pulse incubation	None

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender
Electroporation Medium*	Dulbecco's Phosphate Buffered Saline (minus Ca ⁺⁺ , Mg ⁺⁺)	Cuvette Gap	0.4 cm
Cell Density	4 to 5 x 10 ⁽⁶⁾ cells / 800 µl	Voltage	0.22 kV
Volume of Cells	800 µl	Field Strength	0.55 kV/cm
DNA Concentration	5 to 20 µg	Capacitor	960 µF
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA)	Resistor	(Pulse Controller) Ω none
Volume of DNA	5 to 50 µl (usually less than 10 µl)	Time Constant	about 12 msec

After the Pulse

Outgrowth Medium	DMEM/F12 (1:1) + 10 % Fetal Bovine Serum
Outgrowth Temperature	37 °C
Length of Incubation	usually 24 hours
Selection Method or Assay Used	Luciferase assay, β-galactosidase assay
Electroporation Efficiency	about 1 x 10 ⁽⁴⁾ / µg [20 to 40 % cells stained blue, β-gal, transient assay]
Per Cent Survival	50 to 80 %

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
Reference: Tverberg, L.A. and Russo, A.F. 1992. *J. Biol. Chem.* **267**(5):17567-17573.

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 9/4/92
Survey Number 197
 © Bio-Rad Laboratories, 1993