



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	DNA: pSU2 neo, PAG 60
Species Used	Human, pancreatic cell lines	Electroporated	

Before the Pulse

Cell growth medium	Daigo's T (10% Fetal Calf Serum)	Growth phase at harvest	Not given
		Pre-pulse incubation	Not given
Wash solution	Hanks' Balanced Salt Solution		

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus
Electroporation Medium*	Hanks' Balanced Salt Solution		
Cell Density	1.5 x 10 ⁽⁶⁾ / ml	Cuvette Gap	Not given
Volume of Cells	1 ml	Voltage	0.50 to 0.7 kV
DNA Concentration	1 to 10 µg / ml	Field Strength	Not given
DNA Resuspension Buffer	Hanks' Balanced Salt Solution	Capacitor	25 µF
Volume of DNA	1 ml	Resistor	(Pulse Controller) Ω none
		Time Constant	0.9 msec

After the Pulse

Outgrowth Medium	Daigo's T (5% FCS)	Relevant Publications and/or Comments	
		Note:	exponential values designated in parentheses.
Outgrowth Temperature	37 °C		
Length of Incubation	not given		
Selection Method or Assay Used	pSV2 Neo (G418)		
Electroporation Efficiency	1 to 20 clones / 1µl DNA; depends on the cell line		
Per Cent Survival	50 %		

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 4/10/91
Survey Number 113
 © Bio-Rad Laboratories, 1993