



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 Electroporated	DNA: plasmids of 8.4 kb, linearized
Species Used	Hamster, CHO, ovary		

Before the Pulse

Cell growth medium	DMEM, 10% Fetal Bovine Serum, NEAA, Pennicillin /Streptomycin	Growth phase at harvest	Log
Wash solution	Berg buffer (see Ref in notes)	Pre-pulse incubation	10 min on ice

The Pulse

Electroporation Temperature	4 °C	Instruments Used	Gene Pulser®
Electroporation Medium*	Berg buffer (see Ref. in notes)	Cuvette Gap	0.4 cm
Cell Density	10(8) cells /0.8ml	Voltage	0.80 kV
Volume of Cells	0.8 ml	Field Strength	2.0 kV/cm
DNA Concentration	20 µg	Capacitor	25 µF
DNA Resuspension Buffer	water	Resistor	(Pulse Controller) Ω none
Volume of DNA	20 µl	Time Constant	0.6 to 0.8 msec

After the Pulse

Outgrowth Medium	DMEM, 10% FBS, NEAA, Pen/Strep	Relevant Publications and/or Comments Note: exponential values designated in parentheses. Ref: Chu,G., Hayakawa, H. and Berg, P. <i>NAR</i> 15 , 1131. Berg buffer: 20 mM HEPES, pH 7.05, 137 mM NaCl, 5 mM KCl, 0.7 mM Na ₂ HPO ₄ , 6 mM dextrose.
Outgrowth Temperature	37 °C	
Length of Incubation	2 weeks	
Selection Method or Assay Used	G418, 8-Aza adenine	
Electroporation Efficiency	10 transfectants / µg DNA	
Per Cent Survival	20 to 80 % (varies)	

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 2/27/91
Survey Number 089
© Bio-Rad Laboratories, 1993