



# 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.

<b>Cell Type</b>	Mammalian, adherent	<b>Molecules</b>	DNA: CMV $\beta$ -gal , 6 kB, supercoiled
<b>Species</b>	Monkey, COS-7, kidney, SV-40	<b>Electroporated</b>	
<b>Used</b>	transformed		

## Before the Pulse

<b>Cell growth medium</b>	DMEM + 10% Fetal Bovine Serum (GIBCO/BRL, Sigma)	<b>Growth phase at harvest</b>	Log phase, 70 to 80% confluent
<b>Wash solution</b>	Trypsinize	<b>Pre-pulse incubation</b>	Room temperature

## The Pulse

<b>Electroporation Temperature</b>	Room temperature	<b>Instruments Used</b>	Gene Pulser® apparatus & Capacitance Extender
<b>Electroporation Medium*</b>	DMEM + 10% Fetal Bovine Serum + 5mM BES, pH 7.2	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	2 x 10 <sup>(7)</sup> cells / ml	<b>Voltage</b>	0.170 kV
<b>Volume of Cells</b>	10 $\mu$ g / 250 $\mu$ l DNA	<b>Field Strength</b>	0.425 kV/cm
<b>DNA Concentration</b>	10 mM Tris, 1 mM EDTA, pH 8.0)	<b>Capacitor</b>	960 $\mu$ F
<b>DNA Resuspension Buffer</b>	Not given	<b>Resistor</b>	(Pulse Controller) $\Omega$ none
<b>Volume of DNA</b>	Not given	<b>Time Constant</b>	45 msec

## After the Pulse

<b>Outgrowth Medium</b>	Not given	<b>Relevant Publications and/or Comments</b>
<b>Outgrowth Temperature</b>	Room temperature	Note: exponential values designated in parentheses.
<b>Length of Incubation</b>	10 min., then pellet& resuspend in DMEM	See reference: Ustav, M., and Stenlund, A. 1991. <i>EMBO J.</i> <b>10</b> (2):449-457.
<b>Selection Method or Assay Used</b>	Not given	
<b>Electroporation Efficiency</b>	25%	
<b>Per Cent Survival</b>	about 100 %	

**Name of Submitter**  
**Institution Address**

**Telephone Number**  
**Fax Number**  
**Date Submitted** 8/12/91  
**Survey Number** 127  
© Bio-Rad Laboratories, 1993