

# 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 and suspension	<b>Molecules Electroporated</b>	DNA: 2 to 4 kB, supercoiled
<b>Species Used</b>	Rat, PC12,adrenal pheochromocytoma; Rat brain; Simian (monkey) COS, kidney cells;		

## Before the Pulse

<b>Cell Growth Medium</b>	DMEM (GIBCO/BRL, Sigma)	<b>Growth Phase at Harvest</b>	Stationary growth
		<b>Pre-pulse Incubation</b>	10 min.
<b>Wash Solution</b>	Dulbecco's PBS		

## The Pulse

**Instruments Used** Gene Pulser® apparatus & Capacitance

<b>Electroporation Temperature</b>	4 °C		
<b>Electroporation Medium*</b>	Dulbecco's PBS	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	5 x 10 (6) cells/ ml	<b>Voltage</b>	0.250 kV
<b>Volume of Cells</b>	400 to 800 µl	<b>Field Strength</b>	0.625 kV/cm
<b>DNA Concentration</b>	20 to 200 µg / ml		
<b>DNA Resuspension Buffer</b>	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	<b>Capacitor</b>	25 to 500 µF
<b>Volume of DNA</b>	10 to 20 µl	<b>Resistor</b>	(Pulse Controller) Ω none
<b>After the Pulse</b>		<b>Time Constant</b>	0.6 to 16 msec
<b>Outgrowth Medium</b>	DMEM		

## Relevant Publications and/or Comments

**Note:** exponential values designated in parentheses.

<b>Outgrowth Temperature</b>	37 °C
<b>Length of Incubation</b>	days
<b>Selection Method or Assay Used</b>	Binding assays / transients; G418 selection / stable
<b>Electroporation Efficiency</b>	varies
<b>Per Cent Survival</b>	varies

**Name of Submitter** Dr.Catherine Bitler

**Institution Address** Molecular Pharmacologist  
SRI International  
Neuroscience Department  
333 Ravenswood  
Menlo Park, CA 94025

**Survey Number**

195