



# 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, suspension	<b>Molecules</b>	DNA: variety of constructs,
<b>Species</b>	Mouse, erythroleukemia cells	<b>Electroporated</b>	supercoiled, usually <10 kB (some
<b>Used</b>			pUC based)

## Before the Pulse

<b>Cell growth medium</b>	DMEM + 10% Fetal Calf Serum, 1 x glutamine, 1 x penicillin / streptomycin (GIBCO/BRL, Sigma)	<b>Growth phase</b>	Log phase
		<b>at harvest</b>	
<b>Wash solution</b>	DMEM + 10% Fetal Calf Serum, 1 x glutamine	<b>Pre-pulse</b>	10 min., room temperature
		<b>incubation</b>	

## The Pulse

<b>Electroporation</b>	Room temperature	<b>Instruments</b>	Gene Pulser® apparatus &
<b>Temperature</b>		<b>Used</b>	Capacitance Extender
<b>Electroporation</b>	DMEM + 1% Fetal Calf Serum		
<b>Medium*</b>		<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	1.43 x 10 <sup>(7)</sup> cells / ml	<b>Voltage</b>	0.250 kV
<b>Volume of Cells</b>	700 µl	<b>Field</b>	0.625 kV/cm
		<b>Strength</b>	
<b>DNA Concentration</b>	20 to 30 µg / 100 µl	<b>Capacitor</b>	960 µF
<b>DNA Resuspension</b>	10 mM Tris, pH 8.0, 1 mM EDTA	<b>Resistor</b>	(Pulse Controller ) Ω none
<b>Buffer</b>		<b>Time</b>	16 msec (some small variation)
<b>Volume of DNA</b>	100 µl	<b>Constant</b>	

## After the Pulse

<b>Outgrowth Medium</b>	DMEM + 10% Fetal Calf Serum, 1 x glutamine, 1 x penicillin / streptomycin	<b>Relevant Publications and/or Comments</b>	
<b>Outgrowth Temperature</b>	37 °C	<b>Note:</b>	exponential values designated in parentheses.
<b>Length of Incubation</b>	24 to 48 hours		
<b>Selection Method or</b>	CAT assay		
<b>Assay Used</b>			
<b>Electroporation</b>	Not given		
<b>Efficiency</b>			
<b>Per Cent Survival</b>	35 %		

<b>Name of</b>	<b>Telephone Number</b>
<b>Submittor</b>	
<b>Institution</b>	<b>Fax Number</b>
<b>Address</b>	<b>Date Submitted</b> 8/22/90
	<b>Survey Number</b> 132
	© Bio-Rad Laboratories, 1993