



# 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 Electroporated</b>	DNA: linear and circular plasmids, 4-12 kB in size
<b>Species Used</b>	Mouse, 32d, J558L, myeloma		

## Before the Pulse

<b>Cell growth medium</b>	RPMI, 10% Fetal Calf Serum GIBCO/BRL, Sigma)	<b>Growth phase at harvest</b>	log
<b>Wash solution</b>	log	<b>Pre-pulse incubation</b>	10 minutes ice

## The Pulse

<b>Electroporation Temperature</b>	(ice) 0 °C	<b>Instruments Used</b>	Gene Pulser® apparatus & Capacitance Extender, and Pulser Controller
<b>Electroporation Medium*</b>	RPMI, 10% Nu-serum	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	10 (7) cells / $\mu$ l	<b>Voltage</b>	0.20 to 0.40 kV
<b>Volume of Cells</b>	800 $\mu$ l	<b>Field Strength</b>	0.5 to 1.0 kV/cm
<b>DNA Concentration</b>	1 mg / ml	<b>Capacitor</b>	960 $\mu$ F
<b>DNA Resuspension Buffer</b>	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	<b>Resistor</b>	(Pulse Controller) none $\Omega$
<b>Volume of DNA</b>	20 to 40 $\mu$ l	<b>Time Constant</b>	17 msec

## After the Pulse

<b>Outgrowth Medium</b>	Not given	<b>Relevant Publications and/or Comments</b>	<b>Note:</b> exponential values designated in parentheses.
<b>Outgrowth Temperature</b>	Not given		
<b>Length of Incubation</b>	Not given		
<b>Selection Method or Assay Used</b>	G418; Hygromycin B		
<b>Electroporation Efficiency</b>	Not given		
<b>Per Cent Survival</b>	Not given		

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

**Telephone Number**  
**Fax Number**  
**Date Submitted** 3/5/91  
**Survey Number** 134  
© Bio-Rad Laboratories, 1993