



# 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 Electroporated</b>	DNA: plasmids and cosmids, 40 kb and smaller.
<b>Species Used</b>	Mouse, 3T3, embryo; Human, fibroblast; primary cell lines; Monkey, Vero, kidney cells.		

## Before the Pulse

<b>Cell growth medium</b>	DMEM + 10% Nu serum (GIBCO/BRL, Sigma, Flow Labs)	<b>Growth phase at harvest</b>	Passaged 24 hours prior to pulse.
<b>Wash solution</b>	Not given	<b>Pre-pulse incubation</b>	minimum

## The Pulse

<b>Electroporation Temperature</b>	25 °C	<b>Instruments Used</b>	Gene Pulser® apparatus & Capacitance Extender
<b>Electroporation Medium*</b>	Opti medium (BRL)	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	10 (7) cells / pulse	<b>Voltage</b>	0.220 kV
<b>Volume of Cells</b>	0.4 ml	<b>Field Strength</b>	0.55 kV/cm
<b>DNA Concentration</b>	10 to 20 µg / pulse	<b>Capacitor</b>	960 µF
<b>DNA Resuspension Buffer</b>	20 µl TE (10 mM Tris, 1 mM EDTA, pH 8.0) / pulse	<b>Resistor</b>	(Pulse Controller) Ω none
<b>Volume of DNA</b>	20 µl	<b>Time Constant</b>	30 msec

## After the Pulse

<b>Outgrowth Medium</b>	DMEM + 10% Nu serum	<b>Relevant Publications and/or Comments</b>	<b>Note:</b> exponential values designated in parentheses.
<b>Outgrowth Temperature</b>	37 °C		
<b>Length of Incubation</b>	24 hours		
<b>Selection Method or Assay Used</b>	β-gal expression: X-gal staining. Replication of Herpes origin of replication between S plasmids		
<b>Electroporation Efficiency</b>	30 to 50% cells transfected		
<b>Per Cent Survival</b>	60%		

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

**Telephone Number**  
**Fax Number**  
**Date Submitted** 9/27/90  
**Survey Number** 160  
© Bio-Rad Laboratories, 1993