



# 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: 6.5 kB, linearized plamid.
<b>Species Used</b>	Mouse, FDC-PI, Il-3-dependent cell line	<b>Electroporated</b>	

## Before the Pulse

<b>Cell growth medium</b>	RPMI 1640, 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	<b>Growth phase at harvest</b>	Not given
<b>Wash solution</b>	HEPES buffered saline (see notes)	<b>Pre-pulse incubation</b>	Ice, 10 min.

## The Pulse

<b>Electroporation Temperature</b>	0 °C	<b>Instruments Used</b>	Gene Pulser® apparatus
<b>Electroporation Medium*</b>	HEPES buffered saline	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	1 x 10 <sup>(7)</sup> cells / ml	<b>Voltage</b>	1.5 kV
<b>Volume of Cells</b>	0.5 ml	<b>Field Strength</b>	3.75 kV/cm
<b>DNA Concentration</b>	1 mg / ml	<b>Capacitor</b>	25 µF
<b>DNA Resuspension Buffer</b>	TE Buffer (10 mM Tris, 1 mM EDTA, pH 8.0)	<b>Resistor</b>	(Pulse Controller) Ω none. NOT RECOMMENDED*** (see notes)
<b>Volume of DNA</b>	10 µl (10 µg) / pulse	<b>Time Constant</b>	0.8 msec

## After the Pulse

<b>Outgrowth Medium</b>	RPMI 1640, 10 % Fetal Calf Serum	<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>	Not given	<b>Ref:</b> <i>EMBO J.</i> <b>9</b> : 4367-4374 (1990) Molecular cloning and expression of the murine interlenkin -5 receptor.
<b>Selection Method or Assay Used</b>	G418, 400 µg / ml	HEPES Buffered Saline: 140 mM NaCl, 5 mM KCl, 0.75 mM Na <sub>2</sub> HPO <sub>4</sub> , 6 mM dextrose, 25 mM HEPES, pH 7.2
<b>Electroporation Efficiency</b>	10 to 20 transfectants / µg DNA	**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.
<b>Per Cent Survival</b>	60% at 10 min. after electroporation	

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

**Telephone Number**  
**Fax Number**  
**Date Submitted** 4/26/91  
**Survey Number** 133  
© Bio-Rad Laboratories, 1993