



# 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>	Electrofusion
<b>Species</b>	B-cell line, unspecified, and	<b>Electroporated</b>	
<b>Used</b>	heteromyeloma		

## Before the Pulse

<b>Cell growth medium</b>	RPMI+ 10% Fetal Calf Serum+ 2 mM L-Glutamine (GIBCO/ BRL, Sigma)	<b>Growth phase at harvest</b>	Exponential growth
<b>Wash solution</b>	RPMI	<b>Pre-pulse incubation</b>	No

## The Pulse

<b>Electroporation Temperature</b>	Room temperature	<b>Instruments Used</b>	Gene Pulser® apparatus & Capacitance Extender
<b>Electroporation Medium*</b>	RPMI	<b>Cuvette Gap</b>	0.2 cm
<b>Cell Density</b>	10 (8) cells / $\mu$ l	<b>Voltage</b>	0.160 kV
<b>Volume of Cells</b>	200 $\mu$ l	<b>Field Strength</b>	0.800 kV/cm
<b>DNA Concentration</b>	Not given	<b>Capacitor</b>	960 $\mu$ F
<b>DNA Resuspension Buffer</b>	Not given	<b>Resistor</b>	(Pulse Controller) none $\Omega$
<b>Volume of DNA</b>	10 (8) cells / ml	<b>Time Constant</b>	10 msec

## After the Pulse

<b>Outgrowth Medium</b>	HAT medium and azaserine & ouabain	<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>	Not given		
<b>Selection Method or Assay Used</b>	Not given		
<b>Electroporation Efficiency</b>	Not given		
<b>Per Cent Survival</b>	No fusion		

<b>Name of Submitter</b>	<b>Telephone Number</b>
<b>Institution Address</b>	<b>Fax Number</b>
	<b>Date Submitted</b> 5/8/92
	<b>Survey Number</b> 121
	© Bio-Rad Laboratories, 1993