



# 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: plasmid, closed circular, CsCl purified x2; Co-transfection with 2 plasmids; pHIV LTR-CAT and pSV40-Tat; also transfection with pCH110 alone (7.2 kB).
<b>Species Used</b>	Human, U937, hystiocytic lymphoma		

## Before the Pulse

<b>Cell growth medium</b>	RPMI 1640 + 10% fetal calf serum + penicillin/streptomycin, L-glutamine, sodium pyruvate (GIBCO/BRL, Sigma)	<b>Growth phase at harvest</b>	Log growth
<b>Wash solution</b>	Cell growth media	<b>Pre-pulse incubation</b>	10 min at 4° C.

## The Pulse

<b>Electroporation Temperature</b>	4° C	<b>Instruments Used</b>	Gene Pulser® apparatus & Capacitance Extender
<b>Electroporation Medium*</b>	Cell growth media		
<b>Cell Density</b>	1.3 x 10 <sup>(7)</sup> / ml	<b>Cuvette Gap</b>	0.4 cm
<b>Volume of Cells</b>	250 µl	<b>Voltage</b>	0.30 kV
<b>DNA Concentration</b>	Not given	<b>Field Strength</b>	0.75 kV/cm
<b>DNA Resuspension Buffer</b>	Not given	<b>Capacitor</b>	960 µF
<b>Volume of DNA</b>	Not given	<b>Resistor</b>	(Pulse Controller) Ω none
		<b>Time Constant</b>	Not given

## After the Pulse

<b>Outgrowth Medium</b>	Cell growth media	<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 to 48 hr.		
<b>Selection Method or Assay Used</b>	CAT assay, stain for β-galactosidase		
<b>Electroporation Efficiency</b>	About 500 transfectants / µg DNA		
<b>Per Cent Survival</b>	15 %		

<b>Name of Submitter</b>	<b>Telephone Number</b>
<b>Institution Address</b>	<b>Fax Number</b>
	<b>Date Submitted</b> 10/1/90
	<b>Survey Number</b> 117
	© Bio-Rad Laboratories, 1993