



## Gene Pulser® Electroprotocols

**Cell Type** Bacterial, gram negative  
**Species Used** *E. coli*, NM522

**Molecules Electroported** DNA: plasmid DNA with inserts; pUC and pUC derivatives; relaxed circle

### Before the Pulse

<b>Cell growth medium</b>	conventional growth medium (NaCl Tryptone); LB broth; yeast extract	<b>Growth phase at harvest</b>	O.D. (600) = Log phase
<b>Wash solution</b>	Distilled water	<b>Pre-pulse incubation</b>	1 min. on ice

### The Pulse

<b>Electroporation Temperature</b>	0 °C	<b>Instruments Used</b>	Gene Pulser® apparatus Pulse Controller
<b>Electroporation Medium</b>	10% glycerol (distilled water)	<b>Cuvette Gap</b>	0.1 cm
<b>Cell Density</b>	1/500 initial volume	<b>Voltage</b>	2.0 kV
<b>Volume of Cells</b>	50 µl	<b>Field Strength</b>	20 kV/cm
<b>DNA Concentration</b>	1 to 10 ng / µl	<b>Capacitor</b>	25 µF
<b>DNA Resuspension Buffer</b>	Distilled water	<b>Resistor</b>	200 Ω (Pulse Controller)
<b>Volume of DNA</b>	10 µl	<b>Time Constant</b>	4.2 msec

### After the Pulse

<b>Outgrowth Medium</b>	LB growth medium	<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>	30 min.	<b>LB:</b>	1% Bacto tryptone, 0.5% Bacto yeast extract, 0.5% NaCl.
<b>Selection Method or Assay Used</b>	Ampicillin, 40 µg / µl		
<b>Electroporation Efficiency</b>	> 10 (8) transformants / µg DNA		
<b>Per Cent Survival</b>	Not given		

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

**Telephone Number**

**Fax Number**

**Date Submitted** 10/31/90

**Survey Number** 042

© Bio-Rad Laboratories, 1993