

**Gene Pulser® Electroprotocols**

<b>Cell Type</b>	Bacterial, gram positive	<b>Molecules</b>	DNA: plasmid, pAL 5000
<b>Species Used</b>	<i>Mycobacterium smegmatis</i> , ATCC 607	<b>Electroporated</b>	

**Before the Pulse**

<b>Cell growth medium</b>	Mycobacteria, 7H11 (Difco)	<b>Growth phase at harvest</b>	O.D. (600) = 0.5 to 1.0
<b>Wash solution</b>	272 mM sucrose, 7 mM potassium phosphate, 1 mM MgCl <sub>2</sub> , pH 7.0	<b>Pre-pulse incubation</b>	None

**The Pulse**

<b>Electroporation Temperature</b>	Room temperature	<b>Instruments Used</b>	Gene Pulser® apparatus Pulse Controller
<b>Electroporation Medium</b>	272 mM sucrose, 7 mM potassium phosphate, 1 mM MgCl <sub>2</sub> , pH 7.0	<b>Cuvette Gap</b>	0.2 cm
<b>Cell Density</b>	1 X 10 <sup>(9)</sup> cells / ml	<b>Voltage</b>	2.5 kV
<b>Volume of Cells</b>	400 µl	<b>Field Strength</b>	12.5 kV/cm
<b>DNA Concentration</b>	10 µg / 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) 200 Ω
<b>Volume of DNA</b>	10 to 20 µl	<b>Time Constant</b>	4.5 to 5.0 msec

**After the Pulse**

<b>Outgrowth Medium</b>	Mycobacteria, 7H11 (Difco)
-------------------------	----------------------------

**Relevant Publications and/or Comments**

**Note:** exponential values designated in parentheses.

<b>Outgrowth Temperature</b>	37 °C
<b>Length of Incubation</b>	15 to 21 days
<b>Selection Method or Assay Used</b>	Not given
<b>Electroporation Efficiency</b>	Not given
<b>Per Cent Survival</b>	Not given

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

**Telephone Number**

**Fax Number** 58-2-861-1258

**Date Submitted** 3/19/91

**Survey Number** 080

© Bio-Rad Laboratories, 1993