



## Gene Pulser® Electroprotocols

<b>Cell Type</b>	Bacterial, gram negative	<b>Molecules Electroporated</b>	DNA: various plasmids
<b>Species Used</b>	<i>Bradyrhizobium japonicum</i> ; <i>E. coli</i> , species unspecified		

### Before the Pulse

<b>Cell growth medium</b>	For <i>B. japonicum</i> , see reprint; for <i>E. coli</i> , standard protocol, see Pulse Controller or <i>E. coli</i> Pulser™ Manual.	<b>Growth phase at harvest</b>	O.D. (600) =varies - usually exponentially growing
<b>Wash solution</b>	Sterile distilled water	<b>Pre-pulse incubation</b>	5 min on ice

### The Pulse

<b>Electroporation Temperature</b>	See notes	<b>Instruments Used</b>	Gene Pulser® apparatus Pulse Controller
<b>Electroporation Medium</b>	10 % glycerol	<b>Cuvette Gap</b>	0.2 cm
<b>Cell Density</b>	10 <sup>(9)</sup> to 10 <sup>(10)</sup> cfu / ml	<b>Voltage</b>	2.5 kV
<b>Volume of Cells</b>	40 µl	<b>Field Strength</b>	12.5 kV/cm
<b>DNA Concentration</b>	Varies: 12 ng / ml to 4 µg / ml	<b>Capacitor</b>	Not given
<b>DNA Resuspension Buffer</b>	Distilled water	<b>Resistor</b>	200 Ω (Pulse Controller)
<b>Volume of DNA</b>	2 µl	<b>Time Constant</b>	5 msec

### After the Pulse

<b>Outgrowth Medium</b>	For <i>B. japonicum</i> , see reprint: YEGG	<b>Relevant Publications and/or Comments</b>	<b>Note:</b> exponential values designated in parentheses.
<b>Outgrowth Temperature</b>	30 °C		Electroporation temperature: cells and cuvette are on ice - then pulsed
<b>Length of Incubation</b>	20 hr		Publications: Gerinot, M.L., Morisseau, B.A. & T. Klapatch. 1990. Electroporation of <i>Bradyrhizobium japonicum</i> <i>Mol. Gen. Genet.</i> <b>221</b> :287
<b>Selection Method or Assay Used</b>	Various drug resistances		The info on this sheet is for <i>B. japonicum</i> . See reprint. For <i>E. coli</i> we just use standard conditions (see Pulse Controller or <i>E. coli</i> Pulser™ Manual).
<b>Electroporation Efficiency</b>	1.8 x 10 <sup>(5)</sup> ( <i>B. japonicum</i> ) transformants / µg DNA		
<b>Per Cent Survival</b>	20 to 95 %		

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

**Telephone Number**

**Fax Number**

**Date Submitted** 3/21/91

**Survey Number** 048

© Bio-Rad Laboratories, 1993