



## Gene Pulser® Electroprotocols

<b>Cell Type</b>	Bacterial, gram positive	<b>Molecules Electroporated</b>	DNA: plasmid pCW1 (a shuttle vector between <i>E. coli</i> and <i>Coryneform</i> bacterium)
<b>Species Used</b>	<i>Corynebacterium glutamicum</i> ATCC 13032, <i>Brevibacterium flavum</i> ATCC 21475		

---

### Before the Pulse

<b>Cell growth medium</b>	ABG + 0.5% Tween 80 + 2.5% glycine	<b>Growth phase at harvest</b>	O.D. (600) = 0.25
<b>Wash solution</b>	15% glycerol	<b>Pre-pulse incubation</b>	ice

### The Pulse

<b>Electroporation Temperature</b>	4°C	<b>Instruments Used</b>	Gene Pulser® apparatus Pulse Controller
<b>Electroporation Medium</b>	15% cold glycerol	<b>Cuvette Gap</b>	0.2 cm
<b>Cell Density</b>	5 x 10 <sup>(10)</sup> / ml	<b>Voltage</b>	2.5 kV
<b>Volume of Cells</b>	50 to 60 µl	<b>Field Strength</b>	12.5 kV/cm
<b>DNA Concentration</b>	0.4 µg / µl	<b>Capacitor</b>	25 µF
<b>DNA Resuspension Buffer</b>	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	<b>Resistor</b>	200 Ω (Pulse Controller)
<b>Volume of DNA</b>	1 µl	<b>Time Constant</b>	4.6 to 4.9 msec

### After the Pulse

**Outgrowth Medium** SMMC buffer

#### Relevant Publications and/or Comments

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

<b>Outgrowth Temperature</b>	30°C
<b>Length of Incubation</b>	12 hr
<b>Selection Method or Assay Used</b>	Neomycin, 10 µg / ml
<b>Electroporation Efficiency</b>	10 (5) transformants / µg DNA
<b>Per Cent Survival</b>	15 %

---

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

**Telephone Number**

**Fax Number**

**Date Submitted** 12/4/90

**Survey Number** 059

© Bio-Rad Laboratories, 1993