



## Gene Pulser® Electroprotocols

<b>Cell Type</b>	Plant, protoplast	<b>Molecules Electroporated</b>	DNA: plasmids with VP1 maize gene, supercoiled; several promoter / GUS fusions that are activated by VP1.
<b>Species Used</b>	Maize cell protoplast, DeKalb XL82 (scutellum), Mpp		

### Before the Pulse

<b>Cell growth medium</b>	N6 medium	<b>Growth phase at harvest</b>	Mid-log, three days after transfer
<b>Wash solution</b>	Protoplasts made by digesting with enzymes, then washed.	<b>Pre-pulse incubation</b>	DNA plus cells held on ice for 10 minutes prior to electroporation

### The Pulse

<b>Electroporation Temperature</b>	25 °C, but sample pre-chilled	<b>Instruments Used</b>	Gene Pulser® apparatus & Capacitance Extender
<b>Electroporation Medium</b>	Not given	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	4 x 10 <sup>(6)</sup> cells / ml	<b>Voltage</b>	0.2 kV
<b>Volume of Cells</b>	1 ml <b>**</b> (SEE NOTES)	<b>Field Strength</b>	0.5 kV/cm
<b>DNA Concentration</b>	20 to 50 µg / pulse	<b>Capacitor</b>	960 µF
<b>DNA Resuspension Buffer</b>	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	<b>Resistor</b>	(Pulse Controller) none
<b>Volume of DNA</b>	20 to 50 µl / pulse	<b>Time Constant</b>	14 to 16 msec, average

### After the Pulse

<b>Outgrowth Medium</b>	KMØ medium	<b>Relevant Publications and/or Comments</b>
<b>Outgrowth Temperature</b>	25 °C	<b>Note:</b> exponential values designated in parentheses.
<b>Length of Incubation</b>	40 hours	<b>**</b> Maximum volume for 0.4 cm cuvettes is 0.8 ml; greater volumes will deliver a non-uniform pulse to sample.
<b>Selection Method or Assay Used</b>	Fluorescence assay for GUS, luminescence for luciferase	Ref: McCarty, <i>et.al.</i> , <i>Cell</i> , <b>66</b> :895-905(1991).
<b>Electroporation Efficiency</b>	Not done	
<b>Per Cent Survival</b>	70 to 80 %	

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

**Telephone Number**

**Fax Number**

**Date Submitted** 1/9/92

**Survey Number** 206

© Bio-Rad Laboratories, 1993