



## Gene Pulser® Electroprotocols

**Cell Type** Fungal / Yeast  
**Species Used** *Cryptococcus neoformans*, ma5 mutants

**Molecules Electroporated** DNA: supercoiled on linear plasmids containing URA5 gene

### Before the Pulse

<b>Cell growth medium</b>	YEPD	<b>Growth phase at harvest</b>	Logarithmic, O.D.(650) = ~ 1.
<b>Wash solution</b>	270 mM Sucrose, 1 mM MgCl <sub>2</sub> , 10 mM Tris-HCl, pH 7.5, 4 mM DTT	<b>Pre-pulse incubation</b>	None

### The Pulse

<b>Electroporation Temperature</b>	Room temperature	<b>Instruments Used</b>	Gene Pulser® apparatus
<b>Electroporation Medium</b>	270 mM Sucrose, 10 mM Tris HCl, pH 7.5 (no DTT)	<b>Cuvette Gap</b>	0.2 cm
<b>Cell Density</b>	Cells concentrated, 100 fold	<b>Voltage</b>	0.470 kV
<b>Volume of Cells</b>	450 µl	<b>Field Strength</b>	2.35 kV/cm
<b>DNA Concentration</b>	0.1 to 1.0 µg, in 1 to 10 µl TE	<b>Capacitor</b>	25 µF
<b>DNA Resuspension Buffer</b>	TE Buffer (10 mM Tris, 1 mM EDTA, pH8.0)	<b>Resistor</b>	(Pulse Controller) none Ω
<b>Volume of DNA</b>	1 to 10 µl	<b>Time Constant</b>	18 to 22 msec

### After the Pulse

<b>Outgrowth Medium</b>	None
<b>Outgrowth Temperature</b>	Not given
<b>Length of Incubation</b>	Not given
<b>Selection Method or Assay Used</b>	SD media (lacking uracil): 6.7 g yeast nitrogen base per liter without amino acids and 20 g glucose / liter
<b>Electroporation Efficiency</b>	2000 transfectants / µl
<b>Per Cent Survival</b>	Not tested

#### Relevant Publications and/or Comments

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

**Note:** the protocol described in the paper by Edman and Kwon-Chung, *Mol. & Cell. Biol.*, **10**:4538-4544 (1990) is not the one described above. The one above gives 10-100x greater efficiency.

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

**Telephone Number**

**Fax Number**

**Date Submitted** 9/20/90

**Survey Number** 173

© Bio-Rad Laboratories, 1993