



## Gene Pulser® Electroprotocols

<b>Cell Type</b>	Other Cell Types	<b>Molecules Electroporated</b>	DNA: plasmid constructs (containing VSG- gene, promoter <i>T. brucei</i> )
<b>Species Used</b>	<i>Trypanosoma brucei brucei</i> , AnTat 1.3A, (blood- stream forms); EATRO 1125 (procyclic forms).		

### Before the Pulse

<b>Cell growth medium</b>	Cunningham's medium + 15% Fetal Calf Serum (See references in notes)	<b>Growth phase at harvest</b>	Mid- log phase
<b>Wash solution</b>	Not given	<b>Pre-pulse incubation</b>	Not given

### The Pulse

<b>Electroporation Temperature</b>	Ambient temperature	<b>Instruments Used</b>	Gene Pulser® apparatus
<b>Electroporation Medium</b>	Zimmerman's post-fusion medium	<b>Cuvette Gap</b>	0.4 cm
<b>Cell Density</b>	2 x 10 <sup>(7)</sup> / ml	<b>Voltage</b>	1.5 kV
<b>Volume of Cells</b>	500 µl	<b>Field Strength</b>	3.75 kV/cm
<b>DNA Concentration</b>	1 mg / ml	<b>Capacitor</b>	25 µF
<b>DNA Resuspension Buffer</b>	TES buffer	<b>Resistor</b>	(Pulse Controller) none Ω NOT RECOMMENDED*** (see notes)
<b>Volume of DNA</b>	20 to 50 µl	<b>Time Constant</b>	Not given

### After the Pulse

<b>Outgrowth Medium</b>	Cunningham's medium/Baltes Medium (procyclics) / (Bloodstream)	<b>Relevant Publications and/or Comments</b>
<b>Outgrowth Temperature</b>	27 °C procyclics/ 37° C bloodstream	<b>Note:</b> exponential values designated in parentheses.
<b>Length of Incubation</b>	12 to 18 hours	The conditions described are not those used in my own publication, Jefferies, <i>et. al.</i> 1991 <i>Mol. Cell. Biol.</i> <b>11</b> : 338-341, but were used by Clayton <i>et. al.</i> , 1990 <i>Mol. Cell. Biol.</i> <b>10</b> : 3036-3047, to transfect procyclic trypanosomes.
<b>Selection Method or Assay Used</b>	CAT assay	**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.
<b>Electroporation Efficiency</b>	Not done	
<b>Per Cent Survival</b>	30 to 60 %	

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

**Telephone Number**

**Fax Number**

**Date Submitted** 3/12/90

**Survey Number** 191

© Bio-Rad Laboratories, 1993