



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram positive
Species Used *Bacillus sphaericus* 1593

Molecules Electroporated DNA: plasmid pUB110

Before the Pulse

Cell growth medium	MM2G: 3.5% antibiotic medium #3, Difco; 0.5% yeast extract; 0.5% glycerol	Growth phase at harvest	O.D. (600) =exponential growth
Wash solution	10% glycerol	Pre-pulse incubation	Not given

The Pulse

Electroporation Temperature	0 °C	Instruments Used	Gene Pulser® apparatus Pulse Controller
Electroporation Medium	10% glycerol	Cuvette Gap	0.2 cm
Cell Density	10 (8) to 10 (9) cfu / ml	Voltage	2.5 kV
Volume of Cells	50 µl	Field Strength	12.5 kV/cm
DNA Concentration	Not given	Capacitor	25 µF
DNA Resuspension Buffer	Not given	Resistor	400 Ω (Pulse Controller)
Volume of DNA	Not given	Time Constant	7.8 msec

After the Pulse

Outgrowth Medium MM2G

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 30°C
Length of Incubation 90 min.
Selection Method or Assay Used Neomycin resistance
Electroporation Efficiency 10 (6) transformants / µg DNA
Per Cent Survival 25 %

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 8/22/90

Survey Number 056

© Bio-Rad Laboratories, 1993