

Cell Type Bacterial, gram positive

Molecules DNA: plasmid pUB110
ElectroporatedSpecies *Bacillus sphaericus* 1593
Used**Before the Pulse**Cell Growth Medium MM2G: 3.5% antibiotic medium #3, Difco;
0.5% yeast extract; 0.5% glycerolGrowth Phase at O.D. (600) =exponential growth
HarvestPre-pulse Not given
Incubation

Wash Solution 10% glycerol

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation 0 °C
TemperatureElectroporation 10% glycerol
Medium

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

DNA Resuspension Not given
Buffer

Capacitor 25 µF

Volume of DNA Not given

Resistor 400 Ω (Pulse Controller)

After the Pulse

Time Constant 7.8 msec

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 Neomycin resistance
or Assay UsedElectroporation 10 (6) transformants / µg DNA
Efficiency

Per Cent Survival 25%

Name of Submitter William F. Burke

Institution Arizona State University
Address Microbiology Dept.
Tempe, AZ 85287-2701

Survey Number

056