

Cell Type Bacterial, gram negative

Molecules DNA: plasmid pISP-2, 3.3 kB, supercoiled
ElectroporatedSpecies *E. coli*, DH5 α , JM83
Used

Before the Pulse

Cell Growth Medium LB

Growth Phase at O.D. (600) = 0.65
HarvestPre-pulse 30 min at 0°C
Incubation

Wash Solution Deionized water

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation 4 °C
TemperatureElectroporation Deionized water/10% glycerol
Medium

Cuvette Gap 0.2 / 0.1 cm

Cell Density 2.9 x 10 (10) cells / ml

Voltage 2.5 / 2.5 kV

Volume of Cells 40 μ l

Field Strength 12.5 / 25 kV/cm

DNA Concentration not given

DNA Resuspension TE (made up in deionized water)
BufferCapacitor 25 / 25 μ FVolume of DNA 2 μ lResistor 200 / 100 Ω (Pulse Controller)

After the Pulse

Time Constant 4.7 / 2.1 msec

Outgrowth Medium SOC

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.**Comments:** In press, Insect Biochemistry.**SOC:** 2% Bacto tryptone, 0.5% Bacto yeast extract, 10mM NaCl, 2.5mM KCl, 10 mM MgCl₂, 10 mM MgSO₄, 20 mM glucose.**LB:** 1% Bacto tryptone, 0.5% Bacto yeast extract, 0.5% NaCl.

Outgrowth Temperature 37 °C

Length of Incubation 1 hr.

Selection Method LB + ampicillin (100 μ g / ml) + X-gal + IPTG
or Assay UsedElectroporation 1 x 10 (8) to 1 x 10 (9)
Efficiency transformants / μ g DNA

Per Cent Survival 4 - 40 %

Name of Submitter Hilary Mende

Institution CSIRO
Address Division of Entomology
Box 1700
Canberra, 2601
Australia

Survey Number

012