

Cell Type Bacterial, gram negative

Molecules Electroporated DNA: pRK415, pDSK519, (both about 10kB), pLARR5 (20 kB) supercoiled.

Species Used *Pseudomonas syringae*; *Xanthomonas campestris*

## Before the Pulse

Cell Growth Medium KMB

Growth Phase at Harvest O.D. (600) =0.5

Pre-pulse Incubation None

Wash Solution 0.5 M sucrose

## The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Room temperature

Electroporation Medium 0.5 M sucrose

Cuvette Gap 0.2 cm

Cell Density O.D. (600) = 1.0

Voltage 2.5 kV

Volume of Cells 100 µl

Field Strength 12.5 kV/cm

DNA Concentration 1 to 100 ng / µl

DNA Resuspension Buffer TE (10 mM Tris, 1mM EDTA, pH 8.0)

Capacitor 25 µF

Volume of DNA 1 µl

Resistor 200 Ω (Pulse Controller)

## After the Pulse

Time Constant 4 to 5 msec

Outgrowth Medium KMB

## Relevant Publications and/or Comments

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

N. T. Keen, H. Shen and D. A. Cooksey (1990) Introduction of cloned DNA into Plant pathogenic bacteria. In "Molecular Plant Pathology, a practical approach" ed. D. M. Glover (in press).

Outgrowth Temperature 28 °C

Length of Incubation 2 hours

Selection Method or Assay Used Tetracycline or Kanamycin

Electroporation Efficiency 5 x 10<sup>4</sup> transformants / µg DNA

Per Cent Survival Not known

Name of Submitter Hao Shen

Institution Address University of California  
Department of Plant Pathology  
Riverside, CA 92507

Survey Number

053