



Gene Pulser® Electroprotocols

Cell Type	Bacterial, gram negative	Molecules Electroporated	DNA: pRK415, pDSK519, (both about 10kB), pLARR5 (20 kB) supercoiled.
Species Used	<i>Pseudomonas syringae</i> ; <i>Xanthomonas campestris</i>		

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

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus Pulse Controller
Electroporation Medium	0.5 M sucrose		
Cell Density	O.D. (600) = 1.0	Cuvette Gap	0.2 cm
Volume of Cells	100 μ l	Voltage	2.5 kV
DNA Concentration	1 to 100 ng / μ l	Field Strength	12.5 kV/cm
DNA Resuspension Buffer	TE (10 mM Tris, 1mM EDTA, pH 8.0)	Capacitor	25 μ F
Volume of DNA	1 μ l	Resistor	200 Ω (Pulse Controller)
		Time Constant	4 to 5 msec

After the Pulse

Outgrowth Medium	KMB
Outgrowth Temperature	28 °C
Length of Incubation	2 hours
Selection Method or Assay Used	Tetracycline or Kanamycin
Electroporation Efficiency	5 x 10 ⁽⁴⁾ transformants / μ g DNA
Per Cent Survival	Not known

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).

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 8/20/90

Survey Number 053

© Bio-Rad Laboratories, 1993