



Gene Pulser® Electroprotocols

Cell Type	Bacterial, gram positive	Molecules Electroporated	DNA: pGK12, covalently closed circular form, 4.3 kB; pNZ12, covalently closed circular form, 4.3 kB
Species Used	<i>Lactobacillus plantarum</i> , strain MGD 286		

Before the Pulse

Cell growth medium	MRS (Difco) + 1% D,L - threonine	Growth phase at harvest	O.D. (600) = 0.5 to 1.0 (average = 0.7)
Wash solution	Distilled deionized water, room temperature	Pre-pulse incubation	None

The Pulse

Electroporation Temperature	23°C (room temperature)	Instruments Used	Gene Pulser® apparatus Pulse Controller
Electroporation Medium	30% PEG 1000 (Dow, Sigma, filter sterilized)	Cuvette Gap	0.2 cm
Cell Density	10 (9) cells / ml	Voltage	1.5 kV
Volume of Cells	100 µl	Field Strength	7.5 kV/cm
DNA Concentration	?	Capacitor	25 µF
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Resistor	(Pulse Controller) 400 Ω
Volume of DNA	1.5 µl	Time Constant	3.8 to 4.2 msec

After the Pulse

Outgrowth Medium	MRS + 1% D,L - threonine
Outgrowth Temperature	37 °C
Length of Incubation	1 hour
Selection Method or Assay Used	pGK12: 2 µg/ml erythromycin + lincomycin, then chloram.(10 µg/ml) pNZ12: chloramphenicol (10 µg/ml)
Electroporation Efficiency	pGK12: 2x10 ⁽³⁾ / µg DNA pNZ12: 6.5x10 ⁽²⁾ / µg DNA
Per Cent Survival	25 %

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
Strain MGD286: Hill, H.A. & Hill, J.E. 1986. *Current Microbiol.* **13**:91-94.
Method a modification of: Josson K. *et al.* 1989. *Plasmid* **21**:9-20.
Plasmid pGK12: Kok, J., *et al.* 1984. *Appl. Environ. Microbiol.* **48**: 726-731.
Plasmid pNZ12: De Vos, W. 1986. Biomolecular Engineering in the European Community (Magnien, E., ed.) pp.465-471. Martinus Nijoff, Dordrecht.

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 8/27/90

Survey Number 069

© Bio-Rad Laboratories, 1993