

Cell Type Bacterial, gram positive

Molecules Electroported DNA: pGK12, covalently closed circular form, 4.3 kB; pNZ12, covalently closed circular form, 4.3 kB

Species Used *Lactobacillus plantarum*, 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)

Pre-pulse Incubation None

Wash Solution Distilled deionized water, room temperature

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature 23°C (room temperature)

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 pGK12:285 µg / ml; pNZ12:190 µg / ml

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

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 Nijhoff, Dordrecht.

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 %

Name of Submitter Dr. Russel Chan / Emily Rogers

Institution Address Pioneer Hi-Bred International
Microbial Genetics
7300 N.W. 62nd Ave.
Johnston, IA 50131

Survey Number

069