



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram positive
Species Used *Lactobacillus acidophilus* ADH;
 gastrointestinal isolate from human faeces

Molecules Electroporated DNA: pGT633, covalently closed circular form, a native 9.8kb erythromycin resistant *Lactobacillus* plasmid.

Before the Pulse

Cell growth medium Lactobacilli MRS broth (Difco)

Growth phase at harvest O.D. (600) = log phase cells, 0.8

Pre-pulse incubation 0°C for 1 min.

Wash solution 3.5X SMEB (Luchansky *et al.* 1988 BioRad Bulletin 1350:1-3)

The Pulse

Electroporation Temperature 0 °C
Electroporation Medium 3.5X SMEB (1x = 272 mM sucrose, 1 mM MgCl₂)
Cell Density 10 (9) cells / ml
Volume of Cells 800 µl
DNA Concentration 10 µg
DNA Resuspension Buffer TE (10 mM Tris, 1 mM EDTA)
Volume of DNA 5 µl

Instruments Used Gene Pulser® apparatus

Cuvette Gap 0.4 cm

Voltage 2.5 kV

Field Strength 6.25 kV/cm

Capacitor 25 µF

Resistor Pulse Controller not used.
 **See comments

Time Constant 10 to 15 msec

After the Pulse

Outgrowth Medium Lactobacilli MRS broth (Difco) 10ml

Outgrowth Temperature 37 °C

Length of Incubation 3 hrs.

Selection Method or Assay Used Erythromycin, 25 µg/ml; the Em(R) gene of pGTG33 requires a min. expression time of 3hr. to recover

Electroporation Efficiency Average 8.6 x 10 (1) transformants / µg DNA

Per Cent Survival 17 %

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
 It is NOT RECOMMENDED to use high voltage with out the Pulse Controller. **Ref: (1) H. J. Connell, "Investigation of Methods for the Transformation of Gastrointestinal Strains of Lactobacilli with Plasmid pGT633." Ph.D. Thesis, University of Otago, Dunedin, NEW ZEALAND (1990) (2) This work was carried out under the supervision of Dr. G. Tannock, Dept. of Microbiol., Univ. of Otago, P. O. Box 56, Dunedin, NEW ZEALAND; PH: +64-3-4797713; FAX: 64-3-4741607. Questions regarding the availability of strains and the plasmid pGT633 should be directed to him.

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 12/20/90

Survey Number 062

© Bio-Rad Laboratories, 1993