



Gene Pulser® Electroprotocols

Cell Type Bacterial, gram positive
Species Used *Enterococcus faecalis* JH2-2 and UV202

Molecules Electroporated DNA: pLAR33 (an 18kB ds DNA plasmid), pAM401 (an 10.4 kB ds DNA plasmid)

Before the Pulse

Cell growth medium Brain Heart Infusion (BHI) broth (Difco)

Growth phase at harvest O.D. (600) =1.0 (exponential)

Pre-pulse incubation Buffer, cuvettes, and cells placed on ice

Wash solution 1.5x EP (see notes)

The Pulse

Electroporation Temperature 25 °C
Electroporation Medium 1.5x EP

Instruments Used Gene Pulser® apparatus
Pulse Controller

Cell Density Cells concentrated 140x

Cuvette Gap 0.2 cm

Voltage 2.5 kV

Volume of Cells 50 µl

Field Strength 12.5 kV/cm

DNA Concentration 1 ng DNA

Capacitor 25 µF

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

Resistor 200 Ω (Pulse Controller)

Volume of DNA 2 µl

Time Constant 5.6 to 5.8 msec

After the Pulse

Outgrowth Medium BHI -Difco

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 37 °C

Wash solution: EP=1mM HEPES (pH7.4), 1mM MgCl₂, 0.5M sucrose

Length of Incubation 3 hours

Selection Method or Assay Used BHI agar plates with 10 µg / ml erythromycin

Electroporation Efficiency 2 x 10⁽³⁾ transformants / ng DNA

Per Cent Survival Not determined

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 8/21/90

Survey Number 060

© Bio-Rad Laboratories, 1993