

Cell Type Bacterial, gram negative

Molecules Electroported DNA: plasmid (pBluescript, pSVCAT, pBR322)

Species Used *E. coli*, DH5 α **Before the Pulse**

Cell Growth Medium 2 x YT

Growth Phase at Harvest O.D. (600) =1.0

Pre-pulse Incubation on ice for 20 minutes

Wash Solution ice-cold water

The Pulse

Instruments Used Gene Pulser® apparatus, Pulse Controller,

Electroporation Temperature 0 °C

Electroporation Medium 10% glycerol

Cuvette Gap 0.1 cm

Cell Density 5 x 10 (7) cfu / μ l

Voltage 1.5 kV

Volume of Cells 20 μ l

Field Strength 15kV/cm

DNA Concentration 1 μ g / ml

DNA Resuspension Buffer TE buffer (pH 8.0)

Capacitor 25 μ FVolume of DNA 1 μ lResistor 200 Ω (Pulse Controller)**After the Pulse**

Time Constant 4.6 msec

Outgrowth Medium SOC

Relevant Publications and/or Comments**Note:** exponential values designated in parentheses.**Reference:** T. Tsuji, *et al.* (1990)PNAS. **87**:8835-8839.**SOC:** 2% Bacto tryptone, 0.5% Bacto yeast extract, 10mM NaCl, 2.5mM KCl, 10 mM MgCl₂, 10 mM MgSO₄, 20 mM glucose.**2xYT:** 1.6% Bacto tryptone, 1.0% Bacto yeast extract, 0.5% NaCl.

Outgrowth Temperature 37 °C

Length of Incubation 1 hour

Selection Method or Assay Used Ampicillin

Electroporation Efficiency 0.5 to 1.0 x 10 (9) cfu / μ g DNA

Per Cent Survival 38 %

Name of Submitter Takashi Tsuji

Institution Address Kyushu University
Science, Biology
Hakozaki 6-10-1 Fukuoka, 812
Japan

Survey Number

021