

Cell Type Bacterial, gram negative

Molecules Electroporated DNA, various plasmids

Species Used *E. coli*, DH5 α

Before the Pulse

Cell Growth Medium LB

Growth Phase at Harvest O.D. (600) = 0.8 - 1.0

Pre-pulse Incubation Not done

Wash Solution Water

The Pulse

Instruments Used Gene Pulser® apparatus and Pulse Controller

Electroporation Temperature 0 °C

Electroporation Medium Water + 10% glycerol

Cuvette Gap 0.2 cm / 0.1 cm

Cell Density 3 x 10 (10) cells / ml

Voltage 2.5 kV / 1.6kV

Volume of Cells 40 μ l

Field Strength 12.5 kV/cm // 16 kV/cm

DNA Concentration 10 to 100 μ g DNA

DNA Resuspension Buffer Water or ligation buffer

Capacitor 25 μ FVolume of DNA 2 μ lResistor 200 Ω (Pulse Controller)

After the Pulse

Time Constant 4.8 msec

Outgrowth Medium LB

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
LB: 1% Bacto tryptone, 0.5% Bacto yeast extract, 0.5% NaCl.

Outgrowth Temperature 37 °C

Length of Incubation 1 hour

Selection Method or Assay Used Ampicillin resistance

Electroporation Efficiency 1 x 10 (8) to 1 X 10 (9) transformants / μ g DNA

Per Cent Survival Not given

Name of Submitter Dr. Imre Kovetsdi

Institution Address Lederle Labs
 Bldg. 205/283
 Pearl River, NY 10965

Survey Number

015