

Cell Type Bacterial, gram negative

Molecules Electroporated DNA: double-stranded circular plasmids

Species Used *Salmonella typhimurium*, LT2

Before the Pulse

Cell Growth Medium L Broth

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

Pre-pulse Incubation None

Wash Solution 15% glycerol

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature 0 °C

Electroporation Medium 15% glycerol

Cuvette Gap 0.2 cm

Cell Density 1.6 x 10¹¹ (11) cells / ml

Voltage 2.4 kV

Volume of Cells 40 µl

Field Strength 12.0 kV/cm

DNA Concentration 0.3 µg / µl

Capacitor 25 µF

DNA Resuspension Buffer 10 mM Tris- Cl, EDTA, pH 8.0

Resistor 400 Ω (Pulse Controller)

Volume of DNA 1 to 2 µl

Time Constant 9 to 13 msec

After the Pulse

Outgrowth Medium SOC

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
 Outgrowth Medium: L Broth, 0.01 M MgCl₂, 0.01 M MgSO₄, 20 mM Glucose, 10mM NaCl, 2.5 mM KCl = SOC medium.
 Electroporation Efficiency: We are moving unmodified DNA from *E. coli* into restriction competent *Salmonella* species. **Ref:** Casjens, *et al.*(1991) *Genetics* **127** (4):637-647. We store washed cells (*Salmonella typhimurium*) in 15% glycerol at -80°C, with good "electro-competence" for several months.
LB: 1% Bacto tryptone, 0.5% Bacto yeast extract, 0.5% NaCl.

Outgrowth Temperature 37 °C

Length of Incubation 60 minutes

Selection Method or Assay Used Amp(R), Tet(R), Kan(R), Cam(R)

Electroporation Efficiency 10 (4) transformants / µg *E. coli* DNA

Per Cent Survival Not given

Name of Submitter Sherwood Casjens

Institution Address University of Utah Medical Center
 Cellular, Viral & Molecular Biology
 Salt Lake City, UT 84132

Survey Number

054