

Cell Type Bacterial, gram negative

Molecules Electroporated DNA: plasmid, pKT240, pKT230, pUC19

Species Used *Vibrio anguillarum***Before the Pulse**

Cell Growth Medium Marine broth

Growth Phase at Harvest O.D. (600) =0.4

Pre-pulse Incubation 30°C

Wash Solution 272 mM sucrose, 15% glycerol, 7mM NaHPO₄**The Pulse**

Instruments Used Gene Pulser® apparatus

Electroporation Temperature 0 °C

Electroporation Medium 272 mM sucrose, 15% glycerol, 7mMNa HPO₂

Cuvette Gap 0.4, 0.2, 0.1 cm

Cell Density 10 (10) cells / ml

Voltage 0.65 to 2.5 kV

Volume of Cells 40 µl

Field Strength Varied

DNA Concentration 200 ng / µl

DNA Resuspension Buffer Not given

Capacitor 25 µF

Volume of DNA 1 µl

Resistor 1000 Ω (Pulse Controller)

After the Pulse

Time Constant 5 to 19 msec

Outgrowth Medium Marine broth + 1% glucose

Relevant Publications and/or Comments**Note:** exponential values designated in parentheses.Hamashima, H. *et. al.*, *Microbiol. Immunol.* **34**(f): 703-708 (1990).

Outgrowth Temperature 30 °C

Length of Incubation 1 hour

Selection Method or Assay Used Kanamycin, streptomycin, ampicillin

Electroporation Efficiency 4 x 10⁴ transformants / µg DNA

Per Cent Survival 90%

Name of Submitter Mr. Hon-a-liu Leong

Institution Address The Chinese University of Hong Kong
Biology Department
Shatin H.K.
Hong Kong

Survey Number

055