

Gene Pulser® Electroprotocol

Cell Type Plant, protoplasts

Molecules DNA: pROA93, 16.2 kB, circular.
ElectroporatedSpecies *Hedysotis corymbosa*
Used

Before the Pulse

Cell Growth Medium Murashige & Skoog's medium + 2% sucrose + 2 mg /l, 2,4-D , 1 g/l casein hydrolysate (GIBCO/BRL)

Growth Phase at Not given
Harvest

Pre-pulse Incubation 10 min on ice

Wash Solution Not given

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature 20 °C

Electroporation Medium Phosphate HEPES buffered saline

Cuvette Gap 0.4 cm

Cell Density 1 x 10 (6) protoplasts

Voltage 0.2 to 0.4 kV

Volume of Cells 1 µl

Field Strength 0.250 to 1.125 kV/cm

DNA Concentration 40 µg/ml

Capacitor 10 to 960 µF

DNA Resuspension Buffer Not given

Resistor (Pulse Controller) none Ω

Volume of DNA 5 to 10 µl

Time Constant Not given

After the Pulse

Outgrowth Medium MS medium + 6% glucose + 2 µg/l 2,4-D + 0.1% casein hydrolysate

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 25 °C

Length of Incubation 2 days

Selection Method or Assay Used CAT

Electroporation Efficiency Good

Per Cent Survival Not given

Name of Submitter Dr. Pua Eng Chong

Institution National University of Singapore
Address Inst. of Molecular and Cell Biology
Kent Ridge
Singapore 0511

Survey Number

182