



Gene Pulser® Electroprotocols

Cell Type Plant
Species Used *Nicotiana plumbaginifolia*; protoplasts from leaf

Molecules Electroported DNA: plasmid pTZ I8U with a 3 kB insert (total: 6 kB), circular.

Before the Pulse

Cell growth medium 5 mM CaCl₂; 0.5 M sucrose, 0.3% Macerozyme R-10, 0.3% cellulose "Onozuka" R-10, pH 4.7 (enzymes-Yokult Honsha Co.,Ltd.)
Wash solution 4 mM CaCl₂, 80mM KCl, 8.0% Mannitol, 2 mM Na₂P0₄, pH 7.2

Growth phase at harvest Not given
Pre-pulse incubation 2 hours at 4°C

The Pulse

Electroporation Temperature 2 hr. at 4°C
Electroporation Medium Same as wash solution
Cell Density 10 (6) / ml
Volume of Cells 300 µl
DNA Concentration 3 µg / µl
DNA Resuspension Buffer TE Buffer (10 mM Tris, 1 mM EDTA, pH 8.0)
Volume of DNA 10 µg DNA / pulse, 3 µl.

Instruments Used Gene Pulser® apparatus & Capacitance Extender
Cuvette Gap 0.2 cm
Voltage 0.16 kV
Field Strength 0.8 kV/cm
Capacitor 125 µF
Resistor (Pulse Controller) none Ω
Time Constant 2.8 msec

After the Pulse

Outgrowth Medium 9% Mannitol, 3% sucrose, 100 µg/ml cefotamine (antibiotic)
Outgrowth Temperature 25°C
Length of Incubation 36 hours
Selection Method or Assay Used Transient expression assays
Electroporation Efficiency Not determined
Per Cent Survival Not determined

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

I am no longer a member of this laboratory, but work is in progress there.

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 11/26/91

Survey Number 184

© Bio-Rad Laboratories, 1993