



Gene Pulser® Electroprotocols

Cell Type	Plant, suspension	Molecules Electroporated	DNA: pCH (pUC vector with hygromycin resistant gene)
Species Used	<i>Oryza sativa</i> , cv. Yamahouci or cv. Nihonbare		

Before the Pulse

Cell growth medium	AA medium (Amino Acid Medium)	Growth phase at harvest	Log phase
Wash solution	0.5 mM MES, 70 mM KCl, 4 mM CaCl ₂ , 0.36 M Mannitol	Pre-pulse incubation	0.5 mM MES, 70 mM KCl, 4 mM CaCl ₂ , 0.36 M Mannitol

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus, Capacitance Extender
Electroporation Medium	0.5 mM MES, 70 mM KCl, 4 mM CaCl ₂ , 0.36 M Mannitol	Cuvette Gap	0.4 cm
Cell Density	10 (6) / ml	Voltage	0.25 kV
Volume of Cells	500 µl	Field Strength	0.625 kV/cm
DNA Concentration	Not given	Capacitor	250 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) 200 Ω
Volume of DNA	50 µl	Time Constant	10 to 20 msec

After the Pulse

Outgrowth Medium	Not given	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Outgrowth Temperature	Not given		
Length of Incubation	Not given		
Selection Method or Assay Used	Not given		
Electroporation Efficiency	Not given		
Per Cent Survival	Not given		

Name of
Submittor
Institution
Address

Telephone Number

Fax Number

Date Submitted 4/22/91

Survey Number 185

© Bio-Rad Laboratories, 1993