



## Gene Pulser® Electroprotocols

**Cell Type** Plant

**Species Used** *Lactuca sativa* (aka *chirimen chisha*)

**Molecules Electroporated** DNA: pCaMVneo [NPT II]

### Before the Pulse

**Cell growth medium** MS (Murashige and Skoog) (GIBCO/BRL)

**Growth phase at harvest** Not given

**Pre-pulse incubation** Ice for 10 minutes

**Wash solution** Leaf mesophyll protoplasts

### The Pulse

**Electroporation Temperature** 0 °C

**Instruments Used** Gene Pulser® apparatus & Capacitance Extender

**Electroporation Medium** 5 mM MES, 70 mM KCl, 0.3 M Mannitol, pH 5.8

**Cuvette Gap** 0.4 cm

**Cell Density** 2 x 10<sup>(6)</sup> / ml

**Voltage** 0.250 kV

**Volume of Cells** 500 µl

**Field Strength** 0.625 kV/cm

**DNA Concentration** 100 µg / ml

**Capacitor** 250 µF

**DNA Resuspension Buffer** TE (10 mM Tris, 1 mM EDTA, pH 8.0)

**Resistor** (Pulse Controller) none Ω

**Volume of DNA** 50 µg

**Time Constant** 13.8 msec

### After the Pulse

**Outgrowth Medium** Modified Murashige and Skoog's medium (see Comments for reference)

#### Relevant Publications and/or Comments

**Note:** exponential values designated in parentheses. Reference: Kazumi Amagasa and Toshiaki Kameya, *J. Japan Soc. Hort. Sci.* **57**(4): 620-625, 1989.

**Outgrowth Temperature** 24 °C

**Length of Incubation** 2 months

**Selection Method or Assay Used** G418 Geneticin (5 to 20 µg / ml)

**Electroporation Efficiency** only one

**Per Cent Survival** 10 to 30 %

**Name of Submitter**  
**Institution Address**

**Telephone Number**

**Fax Number**

**Date Submitted** 4/19/91

**Survey Number** 183

© Bio-Rad Laboratories, 1993