



Gene Pulser® Electroprotocols

* We recommend adapting this protocol to use the Gene Pulser electroporation buffer (catalog #165-2676, 165-2677), which increases cell viability and transfection efficiency in mammalian cell lines.

Cell Type Mammalian, suspension
Species Used Mouse, D10.G4.1, T-cell, helper

Molecules Electroporated DNA

Before the Pulse

Cell growth medium DMEM (10% fetal calf serum) + 20% con A supplement (GIBCO/BRL, Sigma)

Growth phase at harvest log phase

Pre-pulse incubation 10 to 15 min on ice

Wash solution Phosphate Buffered Saline without Ca⁺⁺, Mg⁺⁺

The Pulse

Electroporation Temperature Room Temperature
Electroporation Medium* Ca⁺⁺, Mg⁺⁺ free

Instruments Used Gene Pulser® apparatus & Capacitance Extender

Cell Density 10 (7) cells / 800 µl

Cuvette Gap 0.4 cm

Volume of Cells 800 µl

Voltage 0.3 kV

DNA Concentration 10 µg in sterile water or TE

Field Strength 0.95 kV/cm

DNA Resuspension Buffer Not given

Capacitor 960 µF

Volume of DNA 10 to 20 µl

Resistor (Pulse Controller) Ω none

Time Constant 11 msec

After the Pulse

Outgrowth Medium DMEM (10% serum) + 20% con A supplement

Relevant Publications and/or Comments

Outgrowth Temperature 37 °C

Note: exponential values designated in parentheses.

PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH₂PO₄, 1.15g Na₂HPO₄

HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl₂

Length of Incubation 2 days

Selection Method or Assay Used Hygromycin

Electroporation Efficiency not quantified (used CAT assay)

Per Cent Survival 70 to 80 %

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 8/27/90

Survey Number 143

© Bio-Rad Laboratories, 1993