



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 Human, lymphocytes, primary

Molecules Electroporated DNA:

Before the Pulse

Cell growth medium	RPMI-1640 + 10 % Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth phase at harvest	Not given
Wash solution	Not given	Pre-pulse incubation	Not given

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender
Electroporation Medium*	RPMI-1640 + 10 % Fetal Bovine Serum	Cuvette Gap	0.4 cm
Cell Density	5 x 10 ⁶ (6) cells/ ml	Voltage	0.250 kV
Volume of Cells	250 µl / pulse	Field Strength	0.625 kV/cm
DNA Concentration	25 µg / pulse	Capacitor	960 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω
Volume of DNA	Not given	Time Constant	60 msec

After the Pulse

Outgrowth Medium	Not given
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

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
Used conditions as published by Chen, *et al.*, Bio-Rad Technical Bulletin No.1348.

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 9/30/92
Survey Number 194
© Bio-Rad Laboratories, 1993