

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	Molecules Electroporated	DNA:
Species Used	Human, lymphocytes, primary		

Before the Pulse

Cell Growth Medium	RPMI-1640 + 10 % Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	Not given
Wash Solution	Not given		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

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

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Used conditions as pulished by Chen, *et al.*, Bio-Rad Technical Bulletin No.1348.

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 Submitter Steve Dewhurst

Institution Address
University of Rochester Medical Center
Dept. of Microbiology
601 Elmwood Ave.
Box 672
Rochester, NY 14642

Survey Number

194