

Electroprotocols Species List

Mammalian Cells	Survey Number(s)	Mammalian Cells	Survey Number(s)
B-cell line, unspecified.....	121	Human, U373, glioblastoma	115
Hamster, CHO, ovary.....	85–91, 156, 168	Human, U937, hystiocytic lymphoma	106, 116, 117
Human, 293, kidney	108, 156	Human, UC729-6, lymphoblastoid, B-cells	118
Human, B lymphomas: BJAB, P3HR-1, B95-8	157	Human, V79, skin cells, fibroblasts.....	119
Human, C-4I, cervical carcinoma	155	Hybrid, rat/mouse, MEL cells.....	168
Human, CEMx174, T lymphoblastoid.....	92	Hybrid, mouse/human , A9 fibroblast.....	165
Human epithelial cells	158	Hybridomas	164
Human, fibroblast	160, 163	Monkey, COS, kidney	126–128, 154, 161, 195
Human, GCT, fibrous histocytoma.....	93	Monkey, CV-1, kidney	167
Human, HEL cells, eythroleukemia	158, 166, 192	Monkey, Vero, kidney.....	144, 155, 158, 160, 162
Human, HeLa, epithelial carcinoma	94–98, 158, 167, 192, 196	Mouse, 3T3, embryo	160, 167
Human, Hep3b2, hepatocytes	98, 163	Mouse, 32d, myeloma	134
Human, HepG2, hepatoma	99, 161	Mouse, A-9, derivative of mouse L cell	120
Human, HL 60, eythroleukemia	166	Mouse, BALB/c 3T3, clone A31, fibroblast, embryo	122
Human, HuT78, cutaneous T-cell lymphoma	100	Mouse, BbSutA, hematopoietic.....	166
Human, JEG-3, choriocarcinoma.....	101	Mouse, C127, fibroblast, mammary tumor.....	125, 193
Human, Jurkat.....	107	Mouse, C2, muscle myoblast	123
Human, JY, B-cell, Epstein-Barr virus transformed	102	Mouse, C2C12, muscle	124
Human, K562, chronic myeloid leukemia	103–106, 166, 168, 192	Mouse, D10.G4.1, T-cell, helper	143
Human, lymphoblast cell lines, EBV immortalized	108	Mouse, embryonic stem cells.....	129, 130
Human, lymphocytes, primary.....	194	Mouse, erythroleukemia cells	131, 132
Human, MCF-7, breast.....	110	Mouse, FDC-PI, IL-3-dependent cell line.....	133
Human, MRC-5, lung.....	111, 112	Mouse, J558-L, myeloma.....	134, 135
Human, pancreatic	113	Mouse, L-cells.....	163
Human, Raji, Burkitt lymphoma.....	107	Mouse, L929, connective tissue	109
Human, red blood cells.....	114	Mouse, LM(TK-), connective tissue.....	136, 137
Human, squamous cell carcinoma, oral & cervical lines	162	Mouse, mammary epithelial cells.....	154
Human, T-cells.....	107, 159	Mouse, NIH/3T3; embryo	139, 159
Human, TCCSUP (epithelial-like) bladder carcinoma.....	153	Mouse, NSO, myeloma cells.....	140
		Mouse, p3x63AG8; myeloma	167
		Mouse, SP2/0, [Sp-2], myeloma.....	141, 142

Mammalian Cells	Survey Number(s)
Mouse cells, unspecified	146
Mouse, WEHI-3B, myelomonocytic leukemia.....	138
Mouse, X-63, myeloma [p3 X63 - Ag8.653].....	145
Ovine (sheep), CSL503, fetal lung.....	152
Ovine (sheep), R.E., rumen	151
Rat-1	157
Rat brain	195
Rat, CA77, medullary thyroid carcinoma cell line	197
Rat, D202CC, hepatoma	153
Rat, fibroblasts.....	148
Rat, H4-11-E-C3, hepatoma	150
Rat, L-6, myoblast.....	149
Rat, N62 T cells	154
Rat, PC12, adrenal pheochromocytoma	95
Rat, submandibular acini (secretory cells).....	147

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, adherent	Molecules Electroporated	DNA: pAG-6 (pSV2-gpt derivative containing hamster APRT), 8.4 kB, linear.
Species Used	Hamster, CHO -ATS49, ovary		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Bovine Serum + Non Essential Amino Acids (NEAA)+ Pen/Strep (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	10 min., 4°C
Wash Solution	TD (analogous to Phosphate Buffered Saline)		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	4 °C		
Electroporation Medium*	Berg buffer (HEPES Buffered Saline- see notes)	Cuvette Gap	0.4 cm
Cell Density	Not given	Voltage	0.8 kV
Volume of Cells	10 (7) cells / 0.8 mls	Field Strength	2.0 kV/cm
DNA Concentration	1µg / µl		
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	25 µF
Volume of DNA	20 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	0 .9 msec

After the Pulse	
Outgrowth Medium	DMEM + 10% Fetal Bovine Serum + NEAA + Pen/Strep

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	14 days
Selection Method or Assay Used	Alanosine, Azaserine, Adenine (apt+) or HAT (gpt+)
Electroporation Efficiency	2 x10 (7) cells / 20 µg
Per Cent Survival	usually high %

Pennington, S.L., and Wilson, J.H. (1991) Gene targeting in Chinese Hamster Ovary cells is conservative. *PNAS* **88**: 9498-9502.

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

Name of Submitter Sandra Pennington

Institution Address Baylor College of Medicine
Biochemistry
1 Baylor Plaza
Houston, TX 77030

Survey Number
085

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, adherent	Molecules Electroporated	DNA: 11 to 12 kB expression vector, Rep 4.
Species Used	Hamster, CHO -K1, ovary, (requires proline)		

Before the Pulse

Cell Growth Medium	F12 + 10% serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	10 min., on ice.
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature (pulse), then ice		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	2 x 10 ⁵ (5) cells / pulse	Voltage	0.20 to 0.40 kV
Volume of Cells	0.8 ml	Field Strength	0.5 to 1.0 kV/cm
DNA Concentration	0.5 µg / pulse		
DNA Resuspension Buffer	Not given	Capacitor	500 & 960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
		Time Constant	18 to 22 msec

After the Pulse

Outgrowth Medium F12 + serum + hygromycin or neomycin

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	14 to 21 days
Selection Method or Assay Used	Hygromycin or neomycin selection
Electroporation Efficiency	Not done as yet - just beginning
Per Cent Survival	>10%

Name of Submitter Not listed

Institution Address

Survey Number

086

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections; linearized DNA used for stable transfections.
Species Used	Hamster, CHO, ovary		

Before the Pulse

Cell Growth Medium	F12, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer	Pre-pulse Incubation	4°C, 10 min. (option: add 50µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	Phosphate Buffered Saline	Voltage	1.30 kV
Cell Density	5 x 10 (6) cells/pulse for transient assay;	Field Strength	3.25 kV/cm
Volume of Cells	0.5 ml	Capacitor	25 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none. NOT
DNA Resuspension Buffer	Not given; final volume: 0.8 ml	Time Constant	0.4 msec
Volume of DNA	Not given; final volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	F12, 10% Fetal Calf Serum (FCS)		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. **It is NOT RECOMMENDED to use high voltage with out the Pulse Controller. Note: Stable transfections generally do not use carrier DNA. Also, the level of selective agent required to kill off non-transfected cells needs to be established before transfection. The level required should kill non-transfected cells in approximately 7 days. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Length of Incubation	48 to 72 hrs.	
Selection Method or Assay Used	G418 (stable transfections) and transient assays	
Electroporation Efficiency	Not given	
Per Cent Survival	about 50%	

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number
087

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, adherent, suspension	Molecules Electroporated	DNA: linearized plasmids, about 10 kB.
Species Used	Hamster, CHO, ovary		

Before the Pulse

Cell Growth Medium	Hams F12 (Gibco)	Growth Phase at Harvest	70 % confluence
		Pre-pulse Incubation	10 min. on ice
Wash Solution	Phosphate Buffered Saline, pH 7.4		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	0 to 4 °C		
Electroporation Medium*	Phosphate Buffered Saline (PBS), pH 7.4	Cuvette Gap	0.4 cm
Cell Density	5x10 ⁶ (6) cells / 700 µl	Voltage	0.75 kV
Volume of Cells	700 µl	Field Strength	1.88 kV/cm
DNA Concentration	20 to 100 µg / 700 µl		
DNA Resuspension Buffer	PBS or TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	25 µF
Volume of DNA	20 µl to 50 µl; (conc.=1µg / µl)	Resistor	(Pulse Controller) Ω none
		Time Constant	0.5 msec

After the Pulse

Outgrowth Medium Hams F12 (Gibco)

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂	
Length of Incubation	2 days before selection		
Selection Method or Assay Used	F12 - 5158 plus or minus MTX		
Electroporation Efficiency	250 to 1000 transformants / µg DNA		
Per Cent Survival	25%		

Name of Submitter Lars Adamson M.Sc

Institution Address
Kabigen AB
Cell Biology
Strandbergsgatan 49
S-11287 Stockholm, Sweden

Survey Number
088

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, adherent	Molecules Electroporated	DNA: plasmids of 8.4 kb, linearized
Species Used	Hamster, CHO, ovary		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Bovine Serum, NEAA, Pennicillin /Streptomycin	Growth Phase at Harvest	Log
		Pre-pulse Incubation	10 min on ice

Wash Solution Berg buffer (see Ref in notes)

The Pulse

Instruments Used Gene Pulser®

Electroporation Temperature	4 °C		
Electroporation Medium*	Berg buffer (see Ref. in notes)	Cuvette Gap	0.4 cm
Cell Density	10(8) cells /0.8ml	Voltage	0 .80 kV
Volume of Cells	0.8 ml	Field Strength	2.0 kV/cm
DNA Concentration	20 µg	Capacitor	25 µF
DNA Resuspension Buffer	water	Resistor	(Pulse Controller) Ω none
Volume of DNA	20 µl		

After the Pulse

Outgrowth Medium DMEM, 10% FBS, NEAA, Pen/Strep

Time Constant 0 .6 to 0.8 msec

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. Ref: Chu,G., Hayakawa, H. and Berg, P. <i>NAR</i> 15 , 1131. Berg buffer: 20 mM HEPES, pH 7.05, 137 mM NaCl, 5 mM KCl, 0.7 mM Na2HPO4, 6 mM dextrose.
Length of Incubation	2 weeks	
Selection Method or Assay Used	G418, 8-Aza adenine	
Electroporation Efficiency	10 transfectants / µg DNA	
Per Cent Survival	20 to 80% (varies)	

Name of Submitter Not given

Institution Address Not given

Survey Number

089

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, adherent	Molecules Electroporated	Proteins: restriction enzymes (<i>Eco</i> RI, <i>Sca</i> I, <i>Dra</i> I) and catalase
Species Used	Hamster, CHO, ovary		

Before the Pulse

Cell Growth Medium	McCoy's 5a (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Exponential
		Pre-pulse Incubation	Hepes Buffered Saline (HBS)
Wash Solution	Phosphate Buffered Saline (PBS) or serum-free medium		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4°C		
Electroporation Medium*	HEPES Buffered Saline (HBS)	Cuvette Gap	0.4 cm
Cell Density	2 x 10 ⁶ (6) cells / ml	Voltage	0.3 kV
Volume of Cells	0.8 ml	Field Strength	0.750 kV/cm
DNA Concentration	Not given		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	16 to 22 msec
Outgrowth Medium	McCoy's 5A		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. Cortez, F., and Ortiz, T. <i>Mutation Research</i> 246 (1):221-6 PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂	
Length of Incubation	20 hr.		
Selection Method or Assay Used	Not given		
Electroporation Efficiency	Not given		
Per Cent Survival	40 to 60%		

Name of Submitter Dr. Felipe Cortes

Institution Address Faculty of Biology
Cell Biology Avenida Reina Mercedes s.n.
E-41012 Sevilla
Spain

Survey Number

090

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Hamster, CHO, ovary		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C during and after electroporation		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.25 kV
Volume of Cells	300 µl	Field Strength	0.625 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto- chemistry)	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

091

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Human, CEMx174, T lymphoblastoid		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.200 kV
Volume of Cells	300 µl	Field Strength	0.50 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% Fetal Calf Serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto- chemistry)	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

092

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Human, GCT, fibrous histiocytoma, metastasis to lung		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.20 kV
Volume of Cells	300 µl	Field Strength	0.50 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approximately 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, β-gal, immunohisto- chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number
093

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Human, HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, used 0.3 ml	Voltage	0.300 kV
Volume of Cells	300 µl	Field Strength	0.750 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	500 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kB). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.	
Length of Incubation	48 hours		
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto-chemistry).		
Electroporation Efficiency	50 to 100%		
Per Cent Survival	20 to 75%		

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

094

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, adherent, suspension	Molecules Electroporated	DNA: plasmid
Species Used	Human, HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	Not given	Growth Phase at Harvest	50 to 80%
		Pre-pulse Incubation	Ice, 10 minutes
Wash Solution	Not given		

The Pulse

Instruments Used Not given

Electroporation Temperature	Room temperature		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁷ (7) cells / ml	Voltage	0.25 kV
Volume of Cells	800 µl	Field Strength	0.625 kV/cm
DNA Concentration	1 µg / ml		
DNA Resuspension Buffer	TE Buffer (10 mM Tris, 1 mM EDTA)	Capacitor	960 µF
Volume of DNA	5 µl	Resistor	(Pulse Controller) none Ω
		Time Constant	18 to 22 msec

After the Pulse

Outgrowth Medium Not given

Outgrowth Temperature	Not given	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄
Length of Incubation	Not given	
Selection Method or Assay Used	Not given	
Electroporation Efficiency	Not given	
Per Cent Survival	Not given	

Name of Submitter Phang Seng Meng, Graduate Student

Institution Address Dr. Thomas Leong Laboratory
National University of Singapore
Institute Molecular Cell & Biology
Kent Ridge Crescent
Singapore 0511

Survey Number
095

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: pBLCAT2 series, 4.5 kB
Species Used	Human, HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	MEM + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	5 minutes on ice
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4 °C (from ice to chamber at 25 °C)		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	2 x 10 ⁶ (6) cells in cuvette	Voltage	0.250 kV
Volume of Cells	0.8 ml	Field Strength	0.625 kV/cm
DNA Concentration	2 to 20 µg		
DNA Resuspension Buffer	TE Buffer (10 mM Tris, 1 mM EDTA)	Capacitor	960 µF
Volume of DNA	2 to 20 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	18 to 20 msec

After the Pulse

Outgrowth Medium MEM + 10% Fetal Calf Serum

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. <i>Nucleic Acid Research</i> , 18 (3): 465-470 (1990). PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄
Length of Incubation	40 hours	
Selection Method or Assay Used	CAT assay	
Electroporation Efficiency	Not determined	
Per Cent Survival	20 to 60%	

Name of Submitter Mr. Terence Chong

Institution Address National University of Singapore
Institute Molecular Cell & Biology
Kent Ridge Crecent
Singapore 0511

Survey Number
096

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections; linearized DNA used for stable transfections.
Species Used	Human, HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum [FCS] (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
		Pre-pulse Incubation	4° C, 10 min.
Wash Solution	Wash two times in electroporation buffer		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose	Cuvette Gap	0.4 cm
Cell Density	5 x 10(6) cells/pulse for transient assay;	Voltage	0.170 kV
Volume of Cells	0.5 ml	Field Strength	0.425 kV/cm
DNA Concentration	10 µg / pulse		
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Capacitor	960 µF
Volume of DNA	Not given; pulse volume: 0.8 ml	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	30 msec
Outgrowth Medium	DMEM, 10% Fetal Calf Serum (FCS)		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Note: Stable transfections generally do not use carrier DNA. Also, the level of selective agent required to kill off non-transfected cells needs to be established before transfection. The level required should kill non-transfected cells in approximately 7 days.

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

Outgrowth Temperature	37 °C
Length of Incubation	48 to 72 hrs.
Selection Method or Assay Used	G418 (stable transfections) and transient assays
Electroporation Efficiency	Not given
Per Cent Survival	about 50%

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

097

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, adherent	Molecules Electroporated	DNA: various plasmids, about 3kB.
Species Used	Human, Hep3b2, hepatocytes; HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	DMEM, MAB87 /3 + 10% serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
Wash Solution	Not given	Pre-pulse Incubation	5 min.

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C	Cuvette Gap	0.4 cm
Electroporation Medium*	Cell growth medium	Voltage	0. 220 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.55 kV/cm
Volume of Cells	400 µl	Capacitor	960 µF
DNA Concentration	75 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Sterile deionized water	Time Constant	30 msec
Volume of DNA	5 to10 µl		
After the Pulse			
Outgrowth Medium	Same as growth medium		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	Overnight
Selection Method or Assay Used	b-gal staining (X-gal) luciferase
Electroporation Efficiency	52%
Per Cent Survival	90%

Name of Submitter Margie Wilde, Research Assistant

Institution Address Texas Children's Hospital
Pathology Department
6621 Fannin
Houston, Texas 77030

Survey Number

098

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Human, HepG2, hepatoma		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4°C, 10 min. (option: add 50µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.220 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.55 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; final volume: 0.8 ml	Time Constant	20.0 msec
Volume of DNA	Not given; final volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	F12, 10% Fetal Calf Serum (FCS)		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	HBS:	10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Length of Incubation	48 to 72 hrs.		
Selection Method or Assay Used	Transient assays		
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

099

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cyto-megalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Human, HuT78, cutaneous T-cell lymphoma		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml	Voltage	0.250 kV
Volume of Cells	300 µl	Field Strength	0.625 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto- chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	25 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number
100

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, adherent

Molecules Electroporated DNA: growth hormone reporter construct with a mammalian promoter.

Species Used Human, JEG-3, choriocarcinoma cells

Before the Pulse

Cell Growth Medium MEM + 10% Fetal Bovine Serum + 1% glutamine + 1% penicillin / streptomycin (GIBCO/BRL, Sigma)

Growth Phase at Harvest 70% confluency

Pre-pulse Incubation Not given

Wash Solution MEM + 10% Fetal Bovine Serum

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature 0 °C (ice)

Electroporation Medium* MEM + 10% FBS

Cuvette Gap 0.4cm

Cell Density 2.5 x 10⁸ (8) cells / ml

Voltage 0.140 kV

Volume of Cells 0.4 ml

Field Strength 0.35 kV/cm

DNA Concentration 4 µg / µl

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

Capacitor 960 µF

Volume of DNA 12 µl

Resistor (Pulse Controller) Ω none

After the Pulse

Time Constant 50 msec

Outgrowth Medium Not given

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature ice °C

Length of Incubation 10 min

Selection Method or Assay Used Growth hormone RIA or ELISA

Electroporation Efficiency Not given

Per Cent Survival 50 %

Name of Submitter Rich Bond, Principal Scientist

Institution Address Schering Research
Molecular Pharmacology Bldg. B-9-1
60 Orange St.
Bloomfield, NJ 07003

Survey Number

101

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, plasmid (10 kB)
Species Used	Human, JY cell (human B cell, Epstein-Barr virus transformed)		

Before the Pulse

Cell Growth Medium	DMEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	Not given
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	Room temperature		
Electroporation Medium*	Not given	Cuvette Gap	0.4 cm
Cell Density	2 x 10 ⁷ (7) cells / 0.8 ml	Voltage	Not given
Volume of Cells	Not given	Field Strength	Not given
DNA Concentration	Not given	Capacitor	Not given
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) none Ω
Volume of DNA	Not given	Time Constant	Not given
After the Pulse			
Outgrowth Medium	DMEM		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	10 min.
Selection Method or Assay Used	Not given
Electroporation Efficiency	5 x 10 ⁵ (5) transfectants / μ g DNA
Per Cent Survival	1%

Name of Submitter Mr. Junji Takeda

Institution Address Osaka University
Microphysiology Section
Immun. Deficiency Research
Japan

Survey Number

102

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: pRSVneo, linearized, 5.6 kB
Species Used	Human, K562, chronic myeloid leukemia		

Before the Pulse

Cell Growth Medium	RPMI + 20% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log
		Pre-pulse Incubation	Not given
Wash Solution	RPMI		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	RPMI	Cuvette Gap	0.4cm
Cell Density	5 x10 (6) cells / ml	Voltage	up to 2 kV
Volume of Cells	0.4 to 0.8 ml	Field Strength	up to 5 kV/cm
DNA Concentration	5 µg / 800 µl		
DNA Resuspension Buffer	Not given	Capacitor	up to 960 µF
Volume of DNA	up to 50 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	0.4 msec

After the Pulse	
Outgrowth Medium	RPMI + 20% Fetal Calf Serum

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	48 hr.
Selection Method or Assay Used	G418
Electroporation Efficiency	0.3% clonogenic cells / µg
Per Cent Survival	Not given

Name of Submitter Ms. Genevieve M. Croaker

Institution Address Royal Prince Alfred Hospital
Kanematsu Laboratories
Missenden Road
Camperdown, NSW 2050
Australia

Survey Number
103

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	Molecules Electroporated	DNA: pZ189
Species Used	Human, K562, chronic myeloid leukemia		

Before the Pulse

Cell Growth Medium	RPMI 1640 (GIBCO/BRL,Sigma)	Growth Phase at Harvest	10 (6) cells / ml
		Pre-pulse Incubation	Not given
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Not given

Electroporation Temperature	20 to 25 °C (but pre-cooled on ice)		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	1 x 10 (7) cells / ml	Voltage	0.6 kV
Volume of Cells	0.4 ml	Field Strength	1.50 V/cm
DNA Concentration	400 µg / ml		
DNA Resuspension Buffer	TE	Capacitor	25 µF
Volume of DNA	40 µg	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	1.5 to 1.9 msec
Outgrowth Medium	RPMI 1640		

Outgrowth Temperature	37 °C, 5% CO2	Relevant Publications and/or Comments
Length of Incubation	1 day	Note: exponential values designated in parentheses.
Selection Method or Assay Used	Not given	PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH2PO4, 1.15g Na2HPO4
Electroporation Efficiency	Not given	
Per Cent Survival	50%	

Name of Submitter Fumio Yatagai, Scientist

Institution Address
RIKEN (The Inst. of Phy. & Chem. Res.)
Radiation Biology
2-1 Hirosawa
Wako-shi, Saitama 351-01
JAPAN

Survey Number
104

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: linearized (5kB) n-rasCAT plasmids
Species Used	Human, K562, chronic myeloid leukemia.		

Before the Pulse

Cell Growth Medium	RPMI + 20% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	log phase
Wash Solution	RPMI	Pre-pulse Incubation	None

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	RPMI	Voltage	2.0 kV
Cell Density	10 (7) cells / cuvette	Field Strength	5 kV/cm
Volume of Cells	850 µl	Capacitor	25 µF
DNA Concentration	50 µg DNA / cuvette	Resistor	(Pulse Controller) Ω none. NOT
DNA Resuspension Buffer	Not given	Time Constant	0.4 msec
Volume of DNA	50 to 300 µl		
After the Pulse			
Outgrowth Medium	RPMI + 20% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

It is NOT RECOMMENDED to use high voltage without the Pulse Controller.

Outgrowth Temperature	37 °C
Length of Incubation	2 days
Selection Method or Assay Used	G418
Electroporation Efficiency	Not given
Per Cent Survival	Not given

Name of Submitter Ms. Jacqueline Thorn

Institution Address Royal Prince Alfred Hospital
Missenden Road,
Camperdown, NSW 2050
AUSTRALIA

Survey Number

105

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: linearized plasmid DNA; RSV LTR promoting human HLA Class II genes, along with HTK promoter (neo r gene) containing plasmid (co-transfection)
Species Used	Human, K562, chronic myeloid leukemia; U937, histiocytic lymphoma		

Before the Pulse

Cell Growth Medium	RPMI 1640 + 10% Fetal Calf Serum (GIBCO/ BRL, Sigma)	Growth Phase at Harvest	Early log phase growth
		Pre-pulse Incubation	10 min. on ice
Wash Solution	RPMI without serum		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4 °C	Cuvette Gap	0.4 cm
Electroporation Medium*	RPMI without serum	Voltage	0.2 kV
Cell Density	10 (7) cells / ml	Field Strength	0.5 kV/cm
Volume of Cells	0.4 ml	Capacitor	960 µF
DNA Concentration	1 µg / µl	Resistor	(Pulse Controller) none Ω
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Time Constant	15 to 20 msec
Volume of DNA	20 µl		
After the Pulse			
Outgrowth Medium	RPMI 1640 + 10% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. Cells resistant to G418 (1µg/ml) grew out of post-pulse cell culture after about 15-30 days, with low % of survival. G418 resistant cells were then dilution-cloned in preparation for screening a Class II expressing clone. However, before the majority of cells could be screened, the clones were lost. I don't know how efficient electroporation procedure was since we were trying to get expression of a two-chain molecule on the cell surface - many factors could go wrong post-electroporation but prior to protein expression on the cell surface.

Outgrowth Temperature	37 °C
Length of Incubation	15 to 30 days
Selection Method or Assay Used	G418 resistance
Electroporation Efficiency	Unknown (see notes)
Per Cent Survival	<1%

Name of Submitter W. Scott Goebal

Institution Address Indiana Univ. School of Medicine
Riley Hospital Room S-09
702 Barnhill Dr.
Indianapolis, IN 46202-5200

Survey Number

106

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: linear, 5 kB
Species Used	Human, Raji, Burkitt lymphoma; Jurkat, acute T cell leukemia; T-cells		

Before the Pulse

Cell Growth Medium	RPMI 1640 + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	None
Wash Solution	Hepes Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4 °C		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁷ (7) cells / pulse	Voltage	0.40 kV
Volume of Cells	0.5 to 0.75 ml	Field Strength	1.0 kV/cm
DNA Concentration	Not given		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	11 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	Not given
Outgrowth Medium	Not given		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES,pH 7.2,150 mM NaCl, 5 mM CaCl ₂	
Length of Incubation	2 days		
Selection Method or Assay Used	Preparation of RNA or CAT assay		
Electroporation Efficiency	Not quantified		
Per Cent Survival	50%		

Name of Submitter Jeremy M. Boss, Ph.D. Asst. Prof

Institution Address Emory University
Microbiology / Immunology
Rollins Research Center
Atlanta, GA 03022

Survey Number
107

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, adherent, suspension	Molecules Electroporated	DNA, p220LTR, 11.2 kb, supercoiled.
Species Used	Human, 293s, transformed primary embryonic kidney; Lymphoblast cell lines, EBV immortalized		

Before the Pulse

Cell Growth Medium	RPMI, 10% Fetal Calf Serum, Penicillin, Streptomycin, L-Glutamine, or DMEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	80% confluent or 1 x 10 ⁶ cells / ml
		Pre-pulse Incubation	None
Wash Solution	Electroporation Buffer (see notes)		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	See notes	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁷ cells / µl	Voltage	0.260 kV
Volume of Cells	800 µl	Field Strength	0.65 kV/cm
DNA Concentration	1 to 4 µg / µl		
DNA Resuspension Buffer	TE Buffer (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	550 or 960 µF
Volume of DNA	<20 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	6.4 or 11.2 msec

After the Pulse

Outgrowth Medium RPMI, 10% Fetal Calf Serum, Penicillin, Streptomycin, L-Glutamine, or DMEM

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37°C		
Length of Incubation	Infinite		
Selection Method or Assay Used	Hygromycin resistance		
Electroporation Efficiency	4000 transformants / µg DNA		
Per Cent Survival	10%		

Name of Submitter David Van Der Berg

Institution Address Stanford University
Genetics Department
School of Medicine, S-337
Stanford, CA 94305

Survey Number

108

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Mouse, L929, connective tissue, clone of strain L		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum, +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.350 kV
Volume of Cells	300 µl	Field Strength	0.875 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	500 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto-chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

109

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Human, MCF-7, breast		

Before the Pulse

Cell Growth Medium	RPMI, 10% Fetal Calf Serum (FCS), insulin 10µg / ml (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4° C, 10 min. (optional: add 50 µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections)

The Pulse	Instruments Used	Gene Pulser® apparatus & Capacitance
------------------	-------------------------	--------------------------------------

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.22 to 0.23 kV
Cell Density	5 x 10 (6) cells/pulse	Field Strength	0.55 to 0.575 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; final pulse volume: 0.8 ml	Time Constant	22.0 msec
Volume of DNA	Not given; final pulse volume: 0.8 ml		

After the Pulse	
Outgrowth Medium	RPMI, 10% Fetal Calf Serum (FCS), insulin, 10µg/ml

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	48 to 72 hrs.
Selection Method or Assay Used	Transient assays
Electroporation Efficiency	Not given
Per Cent Survival	about 50%

Name of Submitter	Jackie Beall
--------------------------	--------------

Institution Address	University of Adelaide Department of Biochemistry P.O. Box 498 Adelaide 5001 Australia
----------------------------	--

Survey Number	110
----------------------	-----

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Human, MRC-5, lung, diploid		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.3 kV
Volume of Cells	300 µl	Field Strength	0.75 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto- chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	25 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

111

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, adherent	Molecules Electroporated	DNA: neo gene.
Species Used	Human, MRC-5 / V1, lung fibroblasts, transformed		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	37° C, 1/2 hour
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	25 °C	Cuvette Gap	0.4 cm
Electroporation Medium*	DMEM	Voltage	2.0 kV
Cell Density	5 x 10 (6) cells	Field Strength	5.0 kV/cm
Volume of Cells	1.4 ml (**see notes)	Capacitor	25 µF
DNA Concentration	20 µg	Resistor	(Pulse Controller) Ω none; NOT
DNA Resuspension Buffer	water	Time Constant	0.4 msec
Volume of DNA	20 µl		
After the Pulse			
Outgrowth Medium	DMEM, 10% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	1 month for selection
Selection Method or Assay Used	1 mg / ml geneticin
Electroporation Efficiency	Not given
Per Cent Survival	Not given

**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller. It is not recommended to use more than 0.8 ml in the 0.4 cm cuvette; greater volumes may create non-uniform field strengths during the pulse.

Name of Submitter Dr. Ann M. Simpson

Institution Address University of Sydney
Medicine Dept
N.S.W. 2006
Sydney, AUSTRALIA

Survey Number

112

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: pSU2 neo, PAG 60
Species Used	Human, pancreatic cell lines		

Before the Pulse

Cell Growth Medium	Daigo's T (10% Fetal Calf Serum)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	Not given
Wash Solution	Hanks' Balanced Salt Solution		

The Pulse

Instruments Used Gene Pulser ® apparatus

Electroporation Temperature	Room temperature		
Electroporation Medium*	Hanks' Balanced Salt Solution	Cuvette Gap	Not given
Cell Density	1.5 x 10 ⁶ / ml	Voltage	0.50 to 0.7 kV
Volume of Cells	1 ml	Field Strength	Not given
DNA Concentration	1 to 10 µg / ml		
DNA Resuspension Buffer	Hanks' Balanced Salt Solution	Capacitor	25 µF
Volume of DNA	1 ml	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	0.9 msec
Outgrowth Medium	Daigo's T (5% FCS)		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	not given
Selection Method or Assay Used	pSV2 Neo (G418)
Electroporation Efficiency	1 to 20 clones /1µl DNA; depends on the cell line
Per Cent Survival	50%

Name of Submitter Akira Kono, Ph.D.

Institution Address Research Institute, Kyushu Cancer Center
Division of Chrmotherapy
3-1-1 Notamo, Minami-ku
Fukuoka, 815 JAPAN

Survey Number

113

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	Dextrans, various proteins
Species Used	Human, red blood cells		

Before the Pulse

Cell Growth Medium	Harvested from whole blood	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	Held on ice in Phosphate Buffered Saline
Wash Solution	Isotonic Phosphate Buffered Saline (PBS)		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	0 or 25 °C		
Electroporation Medium*	20 mM Phosphate Buffered Saline	Cuvette Gap	0.2 cm
Cell Density	10 (6) to 10 (7) cells / ml	Voltage	0 to 2.5 kV
Volume of Cells	0.4 ml	Field Strength	0 to 12.5 kV/cm
DNA Concentration	10 (-5) to 10 (-4) M dextran		
DNA Resuspension Buffer	Protein in Phosphate Buffered Saline	Capacitor	25 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) not used.
		Time Constant	1 to 2 msec

After the Pulse

Outgrowth Medium Phosphate Buffered Saline

Outgrowth Temperature	0 °C	Relevant Publications and/or Comments	
Length of Incubation	0 to 5 hr.	Note: exponential values designated in parentheses.	
Selection Method or Assay Used	Flow cytometry, detection of fluorescent molecules	**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.	
Electroporation Efficiency	up to 90% of cells exhibit uptake	PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄	
Per Cent Survival	25 to 100%		

Name of Submitter Mark Prausnitz

Institution Address MIT
Chemical Engineering
Room E25-342, 77 Massachusetts Avenue
Cambridge, MA 02139

Survey Number

114

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Human, U373, glioblastoma, astrocytoma, grade III		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.3 kV
Volume of Cells	300 µl	Field Strength	0.75 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto-chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	25 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

115

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: plasmid
Species Used	Human,U937, histiocytic lymphoma		

Before the Pulse

Cell Growth Medium	RPMI 1640, 10 % Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	5 to 10 x10 (5) / ml
		Pre-pulse Incubation	Phosphate Buffered Saline (-)
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	Phosphate Buffered Saline	Voltage	0.2 kV
Cell Density	1 x 10 (8) / ml	Field Strength	0.5 kV/cm
Volume of Cells	2 x 10 (7) / 200 µl	Capacitor	960 µF
DNA Concentration	50 µg / 200 µl	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given	Time Constant	50 to 70 msec
Volume of DNA	50 µg		
After the Pulse			
Outgrowth Medium	RPMI 1640, 10% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	1 day
Selection Method or Assay Used	G418: 700 µg / ml 3 days; then 400 µg / ml 3 days.
Electroporation Efficiency	Not assayed
Per Cent Survival	Not assayed

Name of Submitter Shumsuke Mori

Institution Address Institute for Medical Genetics
Experimental Genetics
Kuhanji 4-24-1
Kumamoto, JAPAN

Survey Number
116

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: plasmid, closed circular, CsCl purified x2; Co-transfection with 2 plasmids; pHIV LTR-CAT and pSV40-Tat; also transfection with pCH110 alone (7.2 kB).
Species Used	Human, U937, hystiocytic lymphoma		

Before the Pulse

Cell Growth Medium	RPMI 1640 + 10% fetal calf serum + penicillin/streptomycin, L-glutamine, sodium pyruvate (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log growth
		Pre-pulse Incubation	10 min at 4° C.
Wash Solution	Cell growth media		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4 °C		
Electroporation Medium*	Cell growth media	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ / ml	Voltage	0.30 kV
Volume of Cells	250 µl	Field Strength	0.75 kV/cm
DNA Concentration	Not given		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	Not given
Outgrowth Medium	Cell growth media		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	24 to 48 hr.
Selection Method or Assay Used	CAT assay, stain for β-galactosidase
Electroporation Efficiency	About 500 transfectants / µg DNA
Per Cent Survival	15%

Name of Submitter Michael S. Bernstein, M.D. Asst. Prof.

Institution Address Univ of California, San Francisco
Dept. of Medicine
Box 0654
San Francisco, CA 94143

Survey Number

117

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: EBV based vectors, 12-20 kB, closed circular
Species Used	Human, UC729-6, lymphoblastoid, B-cells		

Before the Pulse

Cell Growth Medium	RPMI 1640	Growth Phase at Harvest	Mid log at 1 x 10 ⁶ (6) cells / ml
		Pre-pulse Incubation	Hepes buffered saline (HBS)
Wash Solution	Phosphate Buffered Saline (PBS)		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	Hepes Buffered Saline (HBS)	Cuvette Gap	0.4 cm
Cell Density	5 x 10 ⁷ (7) cells / ml	Voltage	0.32 kV
Volume of Cells	1.0 ml total volume (see notes)*	Field Strength	0.8 kV/cm
DNA Concentration	500 µg / ml total		
DNA Resuspension Buffer	450 µg / ml carrier + 50 µg / ml plasmid	Capacitor	960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	12 msec
Outgrowth Medium	RPMI 1640		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. (1.) Margolshee <i>et. al.</i> , <i>Mol. Cell. Biol.</i> 8 : 2837-2847 (1988). (2.) Canfield <i>et. al.</i> , <i>Mol. Cell. Biol.</i> 10 : 1367-1372 (1990). (3.) Spickofsky <i>et. al.</i> , <i>DNA and Prot. Eng. Techniques</i> , 2 : 14-18 (1990). *Maximum recommended volume for 0.4 cm cuvettes is 0.8 ml for uniform field strengths. PBS: 1x = 8g NaCl ,0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂	
Length of Incubation	24 hours		
Selection Method or Assay Used	Hygromycin		
Electroporation Efficiency	2 x 10 ⁴ (4) to 5 x 10 ⁴ (4)		
Per Cent Survival	20 to 30%		

Name of Submitter Robert Margolskee

Institution Address Roche Institute of Molecular Biology
Neuroscience
Nutley, NJ 07110

Survey Number
118

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	Molecules Electroporated	DNA: plasmid.
Species Used	Human, V79, skin cells, fibroblasts		

Before the Pulse

Cell Growth Medium	Modified MEM, 5%Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Actively growing 70% confluent
		Pre-pulse Incubation	10 min / ice
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	Room temperature		
Electroporation Medium*	10 mM HEPES	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / ml	Voltage	0.450 kV
Volume of Cells	0.4 ml	Field Strength	1.125 kV/cm
DNA Concentration	20 µg		
DNA Resuspension Buffer	Not given	Capacitor	25 µF
Volume of DNA	10 µl	Resistor	5 Ω
After the Pulse		Time Constant	10 msec
Outgrowth Medium	Not given		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	1 week
Selection Method or Assay Used	G418 or ampicillin
Electroporation Efficiency	10 (6) transformants / µg
Per Cent Survival	2%

Name of Submitter Phillip K. Liu, Ph.D., Assistant Professor

Institution Address
M.D. Anderson Cancer Center
Laboratory Medicine Box 72
1515 Halcombe Blvd.
Houston, Texas 77030

Survey Number

119

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	Molecules Electroporated	DNA, various linear constructs of human sequences & selectable markers.
Species Used	Mouse, A-9, derivative of mouse L cell (contains human chromosomes)		

Before the Pulse

Cell Growth Medium	DMEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	10 min. ice
Wash Solution	Phosphate Buffered Saline, without Ca++, Mg++		

The Pulse

Instruments Used Not given

Electroporation Temperature	Cuvette on ice just prior to pulse		
Electroporation Medium*	Phosphate Buffered Saline, without Ca++, Mg++	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / ml	Voltage	0.45 kV
Volume of Cells	0.8 ml	Field Strength	1.125 kV/cm
DNA Concentration	1 µg / 1 µl		
DNA Resuspension Buffer	water	Capacitor	500 µF and 960 µF
Volume of DNA	10 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	Not given
Outgrowth Medium	DMEM		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	48 hours
Selection Method or Assay Used	Hygromycin, G418, or MX
Electroporation Efficiency	Low
Per Cent Survival	Not given

Name of Submitter Keith Rosenbach

Institution Address UMDNJ
Microbiology Department
185 South Orange Ave
Newark, NJ 07103

Survey Number
120

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	Electrofusion
Species Used	B-cell line, unspecified, and heteromyeloma		

Before the Pulse

Cell Growth Medium	RPMI+ 10% Fetal Calf Serum+ 2 mM L-Glutamine (GIBCO/ BRL, Sigma)	Growth Phase at Harvest	Exponential growth
Wash Solution	RPMI	Pre-pulse Incubation	No

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.2 cm
Electroporation Medium*	RPMI	Voltage	0.160 kV
Cell Density	10 (8) cells / μ l	Field Strength	0.800 kV/cm
Volume of Cells	200 μ l	Capacitor	960 μ F
DNA Concentration	Not given	Resistor	(Pulse Controller) none Ω
DNA Resuspension Buffer	Not given	Time Constant	10 msec
Volume of DNA	10 (8) cells / ml		
After the Pulse			
Outgrowth Medium	HAT medium and azaserine & ouabain		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	Not given
Selection Method or Assay Used	Not given
Electroporation Efficiency	Not given
Per Cent Survival	No fusion

Name of Submitter Berta Sanchez, Ph.D.

Institution Address Hospital Virgen Del Rocio
Immunologia
Avda Manuez Siurot s/n
41013 Sevilla
Spain

Survey Number

121

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Mouse, BALB/c 3T3, clone A31, fibroblast, whole embryo / fetus, normal		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase. We routinely subdivide cells 24 hours prior to electroporation.
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum, +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.4 kV
Volume of Cells	300 µl	Field Strength	1.0 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	500 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approx. 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, β-gal, immunohisto-chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology, MS1A
Davis, CA 95616

Survey Number

122

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, adherent	Molecules Electroporated	DNA: p1481D (retroviral), 11 kB, supercoiled.
Species Used	Mouse, C2 cell line, muscle myoblast		

Before the Pulse

Cell Growth Medium	DMEM (GIBCO/ BRL, Sigma)	Growth Phase at Harvest	log phase
		Pre-pulse Incubation	10 min.
Wash Solution	Hepes buffered sucrose, PBS, and phosphate buffered sucrose		

The Pulse

Instruments Used Not given

Electroporation Temperature	4 °C or Room temperature		
Electroporation Medium*	Phosphate or HEPES buffered sucrose	Cuvette Gap	0.1 cm
Cell Density	varied	Voltage	1.0 kV
Volume of Cells	200 µl to 800 µl	Field Strength	10 kV/cm
DNA Concentration	1 µg to 10 µg		
DNA Resuspension Buffer	Not given	Capacitor	1 µF
Volume of DNA	1 to 10 µl	Resistor	(Pulse Controller) infinity setting
After the Pulse		Time Constant	0.3 to 0.4 msec
Outgrowth Medium	DMEM		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PB Sucrose: 272 mM sucrose, 7 mM potassium phosphate, pH 7.4, 1 mM MgCl ₂ . HEPES Buffered Sucrose: 272 mM sucrose, 8 mM HEPES, pH 7.4.	
Length of Incubation	24 to 48 hours		
Selection Method or Assay Used	X-gal stain for LacZ		
Electroporation Efficiency	1.5 x 10 ⁽³⁾ transformants / µg DNA		
Per Cent Survival	Not known		

Name of Submitter Barb Rainish - Graduate Student

Institution Address University of Illinois
Biochemistry / Cell Biology
505 S. Goodwin - 506 Morrill Hall
Urbana, IL 61801

Survey Number
123

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Mouse, C2C12, muscle		

Before the Pulse

Cell Growth Medium	DMEM, 20% Fetal Calf Serum (FCS)	Growth Phase at Harvest	50 to 70% confluent
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4° C, 10 min. (optional: add 50 µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections)

The Pulse **Instruments Used** Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: 50 µl salmon sperm DNA).	Voltage	0.220 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.55 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Time Constant	30 msec
Volume of DNA	Not given; pulse volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	DMEM, 20% Fetal Calf Serum (FCS)		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments	
Length of Incubation	48 to 72 hrs.	Note:	exponential values designated in parentheses.
Selection Method or Assay Used	Transient assays	HBS:	10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number
124

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, adherent	Molecules Electroporated	DNA: pC59, an E2 of BPV-1 expression plasmid; p407, an E2 responsive CAT plasmid.
Species Used	Mouse, C127, fibroblast, mammary tumor		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	Room temperature
Wash Solution	1x Phosphate Buffered Saline, 3 times		

The Pulse **Instruments Used** Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	DMEM + 10% FBS + 5mM BES, pH 7.2, 50 µg salmon sperm, carrier DNA	Cuvette Gap	0.4 cm
Cell Density	1 to 5 x 10 ⁶ cells / ml	Voltage	0.21 kV
Volume of Cells	0.5 ml	Field Strength	0.525 kV/cm
DNA Concentration	5 to 10 µg DNA		
DNA Resuspension Buffer	in 5 to 10 µl 10 mM Tris Buffer, pH 8.0	Capacitor	960 µF
Volume of DNA	5 to 10 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	20 msec
Outgrowth Medium	DMEM, 10% Fetal Bovine Serum		

		Relevant Publications and/or Comments
Outgrowth Temperature	37 °C	Note: exponential values designated in parentheses.
Length of Incubation	48 hours	Similar protocol has been published by another lab. Reference is given below:
Selection Method or Assay Used	CAT assay	Ustav, M. and Stenlund, A. (1991) <i>EMBO J.</i> 10 (2): 449-457.
Electroporation Efficiency	Not given	
Per Cent Survival	40 to 60%	

Name of Submitter Shrikant Anant, Graduate Student

Institution Address	Univ of Illinois at Chicago Genetics Dept. 808 S. Wood #1404 Chicago, IL 60612	Survey Number
		125

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections; linearized DNA used for stable transfections.
Species Used	Monkey, COS-1, kidney		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer	Pre-pulse Incubation	4°C, 10 min. (option: add 50µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.3 kV
Cell Density	5 x 10 (6) cells/pulse	Field Strength	0.75 kV/cm
Volume of Cells	0.5 ml	Capacitor	250 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none.
DNA Resuspension Buffer	Not given; final volume: 0.8 ml	Time Constant	9.0 msec
Volume of DNA	Not given; final volume: 0.8 ml		

After the Pulse

Outgrowth Medium F12, 10% Fetal Calf Serum (FCS)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	HBS:	10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Length of Incubation	48 to 72 hrs.		
Selection Method or Assay Used	Transient assays		
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

126

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, adherent	Molecules Electroporated	DNA: CMV b-gal , 6 kB, supercoiled
Species Used	Monkey, COS-7, kidney, SV-40 transformed		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase, 70 to 80% confluent
		Pre-pulse Incubation	Room temperature
Wash Solution	Trypsinize		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	DMEM + 10% Fetal Bovine Serum + 5mM BES, pH 7.2	Cuvette Gap	0.4 cm
Cell Density	2 x 10 ⁷ (7) cells / ml	Voltage	0.170 kV
Volume of Cells	10 µg / 250 µl DNA	Field Strength	0.425 kV/cm
DNA Concentration	10 mM Tris, 1 mM EDTA, pH 8.0)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	45 msec
Outgrowth Medium	Not given		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	Room temperature	
Length of Incubation	10 min., then pellet& resuspend in DMEM	See reference: Ustav, M., and Stenlund, A. 1991. <i>EMBO J.</i> 10 (2):449-457.
Selection Method or Assay Used	Not given	
Electroporation Efficiency	25%	
Per Cent Survival	about 100%	

Name of Submitter Cheng-Ming Chiang

Institution Address University of Rochester Medical Center
Biochemistry, Box 607
601 Elmwood Ave
Rochester, NY 14642

Survey Number
127

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Monkey, COS-7, kidney, SV-40 transformed		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.30 kV
Volume of Cells	300 µl	Field Strength	0.75 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approximately 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, b-gal, immunohisto- chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

128

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, adherent	Molecules Electroporated	DNA: plasmid, 12 kB
Species Used	Mouse, embryonic stem cells		

Before the Pulse

Cell Growth Medium	DMEM + L Glutamine + Pen/Strep + Fetal Calf Serum + mercaptoethanol + non-essential amino acids. (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Exponential growth phase
		Pre-pulse Incubation	Not incubated pre-pulse
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	1.5 to 2.0 x 10 ⁷ cells / pulse	Voltage	0.160 kV
Volume of Cells	0.5 ml	Field Strength	0.4 kV/cm
DNA Concentration	1 µg / ml		
DNA Resuspension Buffer	Phosphate Buffered Saline	Capacitor	960 µF
Volume of DNA	100 µg / pulse	Resistor	(Pulse Controller) Ω none
		Time Constant	14.5 msec

After the Pulse

Outgrowth Medium	DMEM + L Glutamine + Pen/Strep + Fetal Calf Serum + mercaptoethanol + non-essential amino acids.
-------------------------	--

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	2 weeks and more
Selection Method or Assay Used	G418
Electroporation Efficiency	Not given
Per Cent Survival	2.38 x 10 ⁻⁵

Name of Submitter Dr. K.S.E. Cheah

Institution Address Hong Kong University
Biochemistry Department
Sassoon Road
Hong Kong

Survey Number

129

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, adherent	Molecules Electroporated	DNA: pNeo Xp 5.3TK, pHyg Xp5.3 TK, about 10 kB.
Species Used	Mouse, embryonic stem cells		

Before the Pulse

Cell Growth Medium	20% Fetal Calf Serum, DMEM + amino acids, 2- mercaptoethanol, nucleosides. (GIBCO/ BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	5 min., on ice
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4 to 10 °C		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4cm
Cell Density	1 x 10 ⁷ (7) cells / ml	Voltage	0.28 kV
Volume of Cells	0.8 ml	Field Strength	0.7 kV/cm
DNA Concentration	20 µg DNA/ pulse		
DNA Resuspension Buffer	Not given	Capacitor	500 µF
Volume of DNA	16 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	5.8 to 6.4 msec
Outgrowth Medium	Not given		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	10 days
Selection Method or Assay Used	G418, GANC
Electroporation Efficiency	20 to 30 transfectants / µg DNA
Per Cent Survival	20 to 30%

Name of Submitter Kiyoji Tanaka, Associate Professor

Institution Address Inst. Mol. Cell. Biol.
Osaka University, Cell Biology
1-3 Yamada oka
Suita, Osaka, 656
JAPAN

Survey Number
130

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: pSV2 neo, 5 kB
Species Used	Mouse, erythroleukemia cells		

Before the Pulse

Cell Growth Medium	RPMI ,10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	10 to 20 min.
Wash Solution	Not given		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	4 °C		
Electroporation Medium*	Not given	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / 50 µl	Voltage	0.4 kV
Volume of Cells	50 µl	Field Strength	1 kV/cm
DNA Concentration	Not given	Capacitor	25 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) none Ω
Volume of DNA	Not given	Time Constant	10 msec
After the Pulse			
Outgrowth Medium	RPMI,10% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

**Note: Transfection efficiemnces ofthese cells may be increased with the use of the Capacitance Extender (providing longer time constants).

Outgrowth Temperature	37 °C
Length of Incubation	10 to 30 minutes
Selection Method or Assay Used	
Electroporation Efficiency	Low (**see notes)
Per Cent Survival	20 to 70%

Name of Submitter Iliana Coccia Rese

Institution Address Istituto Superiore Sanita
Dept of Virology
Visle Regina Elena 299
Rome , ITALY 00164

Survey Number

131

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: variety of constructs, supercoiled, usually <10 kB (some pUC based)
Species Used	Mouse, erythroleukemia cells		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Calf Serum, 1 x glutamine, 1 x penicillin / streptomycin (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
Wash Solution	DMEM + 10% Fetal Calf Serum, 1 x glutamine	Pre-pulse Incubation	10 min., room temperature

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	DMEM + 1% Fetal Calf Serum	Voltage	0.250 kV
Cell Density	1.43 x 10 (7) cells / ml	Field Strength	0.625 kV/cm
Volume of Cells	700 µl	Capacitor	960 µF
DNA Concentration	20 to 30 µg /100 µl	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	10 mM Tris, pH 8.0, 1 mM EDTA	Time Constant	16 msec (some small variation)
Volume of DNA	100 µl		
After the Pulse			
Outgrowth Medium	DMEM + 10% Fetal Calf Serum, 1 x glutamine, 1 x penicillin / streptomycin		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	24 to 48 hours
Selection Method or Assay Used	CAT assay
Electroporation Efficiency	Not given
Per Cent Survival	35%

Name of Submitter Deborah L. Galson

Institution Address Center for Cancer Research, E17-536
Mass. Inst. of Technology
40 Ames St.
Cambridge, MA 02139

Survey Number
132

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: 6.5 kB, linearized plamid.
Species Used	Mouse, FDC-PI, II-3-dependent cell line		

Before the Pulse

Cell Growth Medium	RPMI 1640, 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	Ice, 10 min.

Wash Solution HEPES buffered saline (see notes)

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	0 °C		
Electroporation Medium*	HEPES buffered saline	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁷ (7) cells / ml	Voltage	1.5 kV
Volume of Cells	0.5 ml	Field Strength	3.75 kV/cm
DNA Concentration	1 mg / ml		
DNA Resuspension Buffer	TE Buffer (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	25 µF
Volume of DNA	10 µl (10 µg) / pulse	Resistor	(Pulse Controller) Ω none. NOT
After the Pulse		Time Constant	0.8 msec

Outgrowth Medium RPMI 1640, 10% Fetal Calf Serum

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	Ref: <i>EMBO J.</i> 9 : 4367-4374 (1990) Molecular cloning and expression of the murine interlenkin -5 receptor. HEPES Buffered Saline: 140 mM NaCl, 5 mM KCl, 0.75 mM Na ₂ HPO ₄ , 6 mM dextrose, 25 mM HEPES, pH 7.2 **It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.
Length of Incubation	Not given	
Selection Method or Assay Used	G418, 400 µg / ml	
Electroporation Efficiency	10 to 20 transfectants / µg DNA	
Per Cent Survival	60% at 10 min. after electroporation	

Name of Submitter Satoshi Takaki, M.D.

Institution Address Inst. for Med. Immunology
Kumamoto Univ. Medical School
Dept. of Biology , 2-2-1 Honjo
Kumamoto, JAPAN 860

Survey Number

133

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: linear and circular plasmids, 4-12 kB in size
Species Used	Mouse, 32d, J558L, myeloma		

Before the Pulse

Cell Growth Medium	RPMI, 10% Fetal Calf Serum GIBCO/BRL, Sigma)	Growth Phase at Harvest	log
		Pre-pulse Incubation	10 minutes ice
Wash Solution	log		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	(ice) 0 °C		
Electroporation Medium*	RPMI, 10% Nu-serum	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / µl	Voltage	0.20 to 0.40 kV
Volume of Cells	800 µl	Field Strength	0.5 to 1.0 kV/cm
DNA Concentration	1 mg / ml		
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	960 µF
Volume of DNA	20 to 40 µl	Resistor	(Pulse Controller) none Ω
After the Pulse		Time Constant	17 msec
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	G418; Hygromycin B
Electroporation Efficiency	Not given
Per Cent Survival	Not given

Name of Submitter Alan Matsumoto

Institution Address Johns Hopkins University
725 N. Wolfe St.
Baltimore, MD 21205

Survey Number

134

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: plasmids, 4.5 kB, supercoiled.
Species Used	Mouse, J558-L, myeloma		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Split on the previous day
		Pre-pulse Incubation	10 min on ice
Wash Solution	None		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	0 °C		
Electroporation Medium*	DMEM + 10% Fetal Calf Serum	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells in 300 µl	Voltage	0.25 kV
Volume of Cells	300 µl	Field Strength	0.625 kV/cm
DNA Concentration	4 µg DNA per 10 (7) cells		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	4 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	50 msec

After the Pulse	
Outgrowth Medium	DMEM + 10% Fetal Calf Serum

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

PNAS, **87**: 5788 (1990).

Electroporation worked best for J558L cells - DEAE dextran did not; neither did lipofection.

Outgrowth Temperature	37 °C
Length of Incubation	48 hours
Selection Method or Assay Used	CAT assay
Electroporation Efficiency	Not given
Per Cent Survival	60 %

Name of Submitter Srikumar Chellappan, Ph.D. Rsch Assoc.

Institution Address Duke University Medical Center
Microbiology and Immunology
P.O. Box 3054
Durham, NC 27710

Survey Number

135

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, adherent	Molecules Electroporated	DNA: pSV2neo, <i>Apa</i> L1 digest.
Species Used	Mouse, LM(TK-), connective tissue [L-M (TK-)].		

Before the Pulse

Cell Growth Medium	Dulbecco's MEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 75% confluent
		Pre-pulse Incubation	None
Wash Solution	Phosphate Buffered Saline, without Ca++ or Mg++		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	4 °C		
Electroporation Medium*	Phosphate Buffered Saline, without Ca++ or Mg++	Cuvette Gap	0.4 cm
Cell Density	1.25 x 10 (7) cells / ml	Voltage	0.950 kV
Volume of Cells	0.8 ml	Field Strength	2.37 kV/cm
DNA Concentration	0.5 µg / ml		
DNA Resuspension Buffer	Phosphate Buffered sucrose	Capacitor	25 µF
Volume of DNA	10 µl	Resistor	(Pulse Controller) Ω none.
After the Pulse		Time Constant	0.4msec
Outgrowth Medium	Dulbecco's MEM		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	48 hours
Selection Method or Assay Used	G-418 , 400 µg / ml after 48 hours recovery
Electroporation Efficiency	1 x 10 (-3) transformants / cell
Per Cent Survival	80 %

**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.

Name of Submitter Deborah J. Hunter, Graduate Student

Institution Address University of Utah
Dept. of Biology
Salt Lake City, Utah 84112

Survey Number

136

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Mouse, LM(TK-), connective tissue, [L-M (TK-)].		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4°C, 10 min. (option: add 50µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.300 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.75 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; final volume: 0.8 ml	Time Constant	25.0 msec
Volume of DNA	Not given; final volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	F12, 10% Fetal Calf Serum (FCS)		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	HBS:	10mM HEPES,pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Length of Incubation	48 to 72 hrs.		
Selection Method or Assay Used	Transient assays		
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

137

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: fos, jun (insert sizes are 4.2 and 1.8 kb using different vectors), linear
Species Used	Mouse, WEHI-3B, myelomonocytic leukemia		

Before the Pulse

Cell Growth Medium	McCoy's 5A (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Exponential
		Pre-pulse Incubation	10 min at room temperature
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁷ (7) cells / ml	Voltage	0.25 kV
Volume of Cells	0.8 ml	Field Strength	0.625 kV/cm
DNA Concentration	15 µg /10 µg	Capacitor	500 µF
DNA Resuspension Buffer	Phosphate Buffered Saline	Resistor	(Pulse Controller) Ω none
Volume of DNA	10 µl	Time Constant	14 msec
After the Pulse			
Outgrowth Medium	McCoy's 5A		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	2 weeks
Selection Method or Assay Used	G-418
Electroporation Efficiency	Not done
Per Cent Survival	50 to 75%

Name of Submitter Jianming Li, Postdoctoral Associate

Institution Address
Yale University School of Medicine
Pharmacology
333 Cedar Street
New Haven, CT 06510

Survey Number

138

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, adherent	Molecules Electroporated	DNA: pBR322 derived plasmids containing retroviral vectors.
Species Used	Mouse, NIH/3T3 derived retroviral vector packaging cell lines		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Bovine Serum + 1% Glutamine (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	5 min, with DNA at room temperature
Wash Solution	Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	DMEM + 10% Fetal Bovine Serum + 1% Glutamine	Cuvette Gap	0.4 cm
Cell Density	1.5 x 10 ⁶ cells / ml	Voltage	0.2 kV
Volume of Cells	0.5 ml	Field Strength	0.5 kV/cm
DNA Concentration	0.1 to 1 µg / µl		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	20 to 25 msec
Outgrowth Medium	DMEM + 10% Fetal Bovine Serum + 1% Glutamine		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	Not given
Selection Method or Assay Used	G418
Electroporation Efficiency	50 to 200 transformants / µg DNA
Per Cent Survival	Not given

Name of Submitter Bruce Sullenger - Graduate Student

Institution Address Memorial Sloan-Kettering Cancer Center
Molecular Biology Dept.
1275 York Avenue, Cost Center 6050
New York, NY 10021

Survey Number

139

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: immunoglobulin genes in pSV vector
Species Used	Mouse, NSO, myeloma cells		

Before the Pulse

Cell Growth Medium	RPMI + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	10 min. on ice
Wash Solution	Not given		

The Pulse

Instruments Used Not given

Electroporation Temperature	0 °C		
Electroporation Medium*	Not given	Cuvette Gap	0.2 cm
Cell Density	12 x 10 (6) / ml	Voltage	0.250 kV
Volume of Cells	400 µl	Field Strength	1.25 kV/cm
DNA Concentration	Not given	Capacitor	960 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	20 µl	Time Constant	0.4 to 0.8 msec
After the Pulse			
Outgrowth Medium	Not given		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	0 °C for 10 min. (ice)
Length of Incubation	Not given
Selection Method or Assay Used	mycophenolic acid
Electroporation Efficiency	None
Per Cent Survival	40 to 50 %

Name of Submitter Rosaria Orlandi

Institution Address Istituto Nazionale Tumori
Oncologia Sperimentale
Via Venezian 1
Milano, 20100
ITALY

Survey Number

140

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, 14.15 kB circular plasmid which contains a part of pSV2neo.
Species Used	Mouse, SP2/0, myeloma [Sp-2]		

Before the Pulse

Cell Growth Medium	10% Fetal Bovine Serum /RPMI 1640 (GIBCO/BRL,Sigma)	Growth Phase at Harvest	5.4 X 10 (5) cells / ml
		Pre-pulse Incubation	on ice for 10 minutes

Wash Solution Phosphate Buffered Saline, 2 times

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	21 °C		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	1.25 x 10 (7) cell / ml	Voltage	0.22 kV
Volume of Cells	800 µl	Field Strength	0.55 kV/cm
DNA Concentration	1.12 µg / µl		
DNA Resuspension Buffer	0.1 x TE buffer (1x=10 mM Tris, 1mM EDTA, pH 8.0)	Capacitor	960 µF
Volume of DNA	18 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	18.5 msec

After the Pulse
Outgrowth Medium 10% Fetal Bovine Serum / RPMI 1640

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C		
Length of Incubation	48 hours		
Selection Method or Assay Used	800 µg / ml of G418		
Electroporation Efficiency	50 µg DNA		
Per Cent Survival	50%		

- *FEBS Lett.* **244**:301 (1989). Y. Kurosawa *et. al.* Convenient plasmid vectors for construction of chimeric mouse/human antibodies.
- *Cancer Research* **50**: 3167 (1990). T. Tsuruo *et. al.* Mouse-Human Chimeric Antibody against the multidrug transporter P-Glycoprotein

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

Name of Submitter Hiroyuki Seimiya

Institution Address University of Tokyo
Institute of Applied Microbiology
1-1-1 Yayoi Bunkyo-ku, Tokyo, 113
Tokyo, Japan

Survey Number

141

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: linearized DNA used for stable transfections.
Species Used	Mouse, SP-2, myeloma [Sp2/0-Ag14].		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
		Pre-pulse Incubation	4° C, 10 min.

Wash Solution Wash two times in electroporation buffer

The Pulse **Instruments Used** Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (5) cells/pulse, stable transfection	Voltage	0.180 kV
Volume of Cells	0.5 ml	Field Strength	0.45 kV/cm
DNA Concentration	10 µg / pulse		
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Capacitor	960 µF
Volume of DNA	Not given; pulse volume: 0.8 ml	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	24 msec
Outgrowth Medium	DMEM, 10% Fetal Calf Serum (FCS)		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Note: Stable transfections generally do not use carrier DNA. Also, the level of selective agent required to kill off non-transfected cells needs to be established before transfection. The level required should kill non-transfected cells in approximately 7 days.

Outgrowth Temperature	37 °C
Length of Incubation	48 to 72 hrs.
Selection Method or Assay Used	G418 (stable transfections)
Electroporation Efficiency	Not given
Per Cent Survival	about 50%

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

142

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	Mouse, D10.G4.1, T-cell, helper		

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	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / 800 µl	Voltage	0.3 kV
Volume of Cells	800 µl	Field Strength	0.95 kV/cm
DNA Concentration	10 µg in sterile water or TE	Capacitor	960 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	10 to 20 µl	Time Constant	11 msec
After the Pulse			
Outgrowth Medium	DMEM (10% serum) + 20% con A supplement		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
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 Yang Park, Postdoc

Institution UMDNJ-RWJMS
Address Pathology
675 Hoes Lane
Piscataway, NJ 08854

Survey Number
143

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, adherent	Molecules Electroporated	DNA: multiple CAT and other vectors under different promoters including HIV, SIV, human and rhesus cytomegalovirus immediate early gene, SV40 and the simian retrovirus type 2.
Species Used	Monkey, Vero, kidney cells		

Before the Pulse

Cell Growth Medium	RPMI 1640 with 10 mM dextrose and 0.1mM dithiothreitol (Sigma, GIBCO/ BRL, Flow Labs)	Growth Phase at Harvest	Log phase (We routinely subdivide cells 24 hours prior to electroporation)
		Pre-pulse Incubation	None
Wash Solution	Trypsin harvest, washed twice in Phosphate Buffered Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	RPMI 1640 without Fetal Calf Serum , +10mM dextrose, 0.1 mM dithiothreitol	Cuvette Gap	0.4 cm
Cell Density	1.3 x 10 ⁷ (7) viable cells / ml, 0.3 ml	Voltage	0.25 kV
Volume of Cells	300 µl	Field Strength	0.625 kV/cm
DNA Concentration	500 µg / ml (5 to 10 µg per pulse)		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	33 to 38 msec
Outgrowth Medium	DMEM culture media, 10% fetal calf serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. All plasmids were prepared by alkaline lysis and banded to equilibrium twice in cesium chloride gradients. Largest DNA molecule electroporated has been the rhesus cytomegalovirus genome (approximately 225 kilobases). In general, the higher the efficiency of transfection, the higher degree of cell killing. For some cell types, > 90% of the input cells are killed, although > 90% of the survivors express very high levels of input DNA. We have found that ≥ 20 µg of DNA per electroporation resulted in greatly reduced efficiency of transfection. Similarly, 2.5 µg of DNA resulted in reduced efficiency.
Length of Incubation	48 hours	
Selection Method or Assay Used	Transient (CAT, β-gal, immunohisto- chemistry).	
Electroporation Efficiency	50 to 100%	
Per Cent Survival	20 to 75%	

Name of Submitter Peter Barry, Ph.D., Asst. Research Virologist

Institution Address Univ. of California- Davis
Department of Medical Pathology , MS1A
Davis, CA 95616

Survey Number

144

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: linearized DNA used for stable transfections.
Species Used	Mouse, X-63, myeloma [p3 X63 - Ag8.653]		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
		Pre-pulse Incubation	4° C, 10 min.
Wash Solution	Wash two times in electroporation buffer		

The Pulse	Instruments Used	Gene Pulser® apparatus & Capacitance
------------------	-------------------------	--------------------------------------

Electroporation Temperature	Room temperature		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (5) cells/pulse, stable transfection	Voltage	0.180 kV
Volume of Cells	0.5 ml	Field Strength	0.45 kV/cm
DNA Concentration	10 µg / pulse		
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Capacitor	960 µF
Volume of DNA	Not given; pulse volume: 0.8 ml	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	24 msec
Outgrowth Medium	DMEM, 10% Fetal Calf Serum (FCS)		

		Relevant Publications and/or Comments	
Outgrowth Temperature	37 °C	Note: exponential values designated in parentheses.	
Length of Incubation	48 to 72 hrs.	Note: Stable transfections generally do not use carrier DNA. Also, the level of selective agent required to kill off non-transfected cells needs to be established before transfection. The level required should kill non-transfected cells in approximately 7 days.	
Selection Method or Assay Used	G418 (stable transfections)	PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄	
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number
145

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, adherent	Molecules Electroporated	DNA: circular
Species Used	Mouse cells, unspecified		

Before the Pulse

Cell Growth Medium	DMEM	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	Not given

Wash Solution Phosphate Buffered Saline

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Room temperature
Electroporation Medium* Not given

Cuvette Gap 0.4 cm

Cell Density 10 (8) cells /250 µl

Voltage 0.40 kV

Volume of Cells 250 µl

Field Strength 1.0 kV/cm

DNA Concentration 20 µg

DNA Resuspension Buffer Phosphate Buffered Saline

Capacitor 25 µF

Volume of DNA 250 µl

Resistor (Pulse Controller) Ω none

After the Pulse

Outgrowth Medium DMEM

Time Constant 1.2 msec

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
 Electroporation efficiencies may be enhanced by use of the Capacitance Extender (provides longer time constants).

Outgrowth Temperature 37 °C
Length of Incubation 48 hours
Selection Method or Assay Used CAT assay

Electroporation Efficiency Not known

Per Cent Survival 20 to 50%

Name of Submitter S. Matsubara

Institution Address Kagoshima University
 Dept. of Medicine
 8-35-1 Sakuragaoka
 Kagoshima City, JAPAN

Survey Number

146

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	Fluorescent dyes (SBFI and MQAE); (3)H-inositol.
Species Used	Rat, submandibular acini (secretory cells).		

Before the Pulse

Cell Growth Medium	KRH + Ca	Growth Phase at Harvest	Not applicable
		Pre-pulse Incubation	A variation of the growth medium, but high in potassium and low in sodium and calcium.
Wash Solution	Cell growth medium		

The Pulse

Instruments Used Not given

Electroporation Temperature	Room temperature		
Electroporation Medium*	Same as pre-pulse	Cuvette Gap	0.4 cm
Cell Density	10 mg protein / ml	Voltage	0.6 to 1.25 kV
Volume of Cells	10 mg protein / ml	Field Strength	1.5 to 3.125 kV/cm
DNA Concentration	1 mg protein / ml		
DNA Resuspension Buffer	Same as Pre-Pulse buffer or KRH + Ca	Capacitor	25 μ F
Volume of DNA	Not assayed	Resistor	(Pulse Controller) Ω none.
After the Pulse		Time Constant	0.4 msec
Outgrowth Medium	Not applicable		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.

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 Sheryll Barker, Research Technologist

Institution Address Lovelace Medical Foundation
Research Dept.
2425 Ridgerest Dr. S.E.
Albuquerque, NM 87108

Survey Number
147

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, adherent	Molecules Electroporated	DNA, retroviral vector, pN2, 9 kB, linearized at <i>Hind</i> III site.
Species Used	Rat, fibroblasts		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	70% confluent
		Pre-pulse Incubation	2.0 min
Wash Solution	None		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	DMEM, 10% Fetal Calf Serum	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / ml	Voltage	0.25 kV
Volume of Cells	250 µl	Field Strength	0.625 kV/cm
DNA Concentration	0.4 µg / µl		
DNA Resuspension Buffer	water	Capacitor	500 µF
Volume of DNA	5 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	18.0 msec
Outgrowth Medium	DMEM, 10% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	10 days
Selection Method or Assay Used	G418 resistance, 500 µl / ml
Electroporation Efficiency	1.21 x 10 (3) transfectants / µg DNA
Per Cent Survival	Not recorded

Name of Submitter Marc D. Corina, Graduate Student

Institution Address Baylor College of Medicine
Cell Biology
One Baylor Plaza
Houston, Texas 77030

Survey Number

148

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, adherent

Molecules Electroporated DNA: supercoiled DNA used for transient transfections.

Species Used Rat, L-6, myoblast

Before the Pulse

Cell Growth Medium DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)

Growth Phase at Harvest 50 to 70% confluency

Pre-pulse Incubation 4° C, 10 min. (optional: add 50 µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

Wash Solution Wash two times in electroporation buffer.

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature Room temperature

Electroporation Medium* HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).

Cuvette Gap 0.4 cm

Cell Density 5 x 10 (6) cells / pulse

Voltage 0.350 kV

Volume of Cells 0.5 ml

Field Strength 0.875 kV/cm

DNA Concentration 10 µg / pulse

DNA Resuspension Buffer Not given; final volume: 0.8 ml

Capacitor 960 µF

Volume of DNA Not given; final volume: 0.8 ml

Resistor (Pulse Controller) Ω none

After the Pulse

Time Constant 25.0 msec

Outgrowth Medium F12, 10% Fetal Calf Serum (FCS)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 37 °C

Length of Incubation 48 to 72 hrs.

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

Selection Method or Assay Used Transient assays

Electroporation Efficiency Not given

Per Cent Survival about 50%

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

149

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Rat, H4-11-E-C3, hepatoma		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4°C, 10 min. (option: add 50µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.240 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.6 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; final volume: 0.8 ml	Time Constant	20.0 msec
Volume of DNA	Not given; final volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	F12, 10% Fetal Calf Serum (FCS)		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	HBS:	10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Length of Incubation	48 to 72 hrs.		
Selection Method or Assay Used	Transient assays		
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number

150

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Ovine (sheep), R.E., rumen		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4° C, 10 min. (optional: add 50 µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse	Instruments Used	Gene Pulser® apparatus & Capacitance
------------------	-------------------------	--------------------------------------

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.270 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.675 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Time Constant	Not given
Volume of DNA	Not given; pulse volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	DMEM, 10% Fetal Calf Serum (FCS)		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Length of Incubation	48 to 72 hrs.	HBS:	10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂
Selection Method or Assay Used	Transient assays		
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number
151

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, adherent	Molecules Electroporated	DNA: supercoiled DNA used for transient transfections.
Species Used	Ovine (sheep), CSL503, fetal lung		

Before the Pulse

Cell Growth Medium	DMEM, 10% Fetal Calf Serum (FCS) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 70% confluency
Wash Solution	Wash two times in electroporation buffer.	Pre-pulse Incubation	4° C, 10 min. (optional: add 50 µl FCS if using HeBS as electroporation media; 50 µl salmon sperm DNA for transient transfections).

The Pulse	Instruments Used	Gene Pulser® apparatus & Capacitance
------------------	-------------------------	--------------------------------------

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	HEPES Buffered Saline, 6mM glucose, (optional: add 50 µl FCS, 50 µl salmon sperm DNA).	Voltage	0.380 kV
Cell Density	5 x 10 (6) cells / pulse	Field Strength	0.95 kV/cm
Volume of Cells	0.5 ml	Capacitor	960 µF
DNA Concentration	10 µg / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Not given; pulse volume: 0.8 ml	Time Constant	28 to 29 msec
Volume of DNA	Not given; pulse volume: 0.8 ml		
After the Pulse			
Outgrowth Medium	DMEM, 10% Fetal Calf Serum (FCS)		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments	
Length of Incubation	48 to 72 hrs.	Note:	exponential values designated in parentheses.
Selection Method or Assay Used	Transient assays	HBS:	10mM HEPES,pH 7.2,150 mM NaCl, 5 mM CaCl2
Electroporation Efficiency	Not given		
Per Cent Survival	about 50%		

Name of Submitter Jackie Beall

Institution Address University of Adelaide
Department of Biochemistry
P.O. Box 498
Adelaide 5001
Australia

Survey Number
152

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, adherent	Molecules Electroporated	DNA: linearized, (pMSG-derivative)
Species Used	Rat, D202CC, hepatoma; Human, TCCSUP (epithelial-like) bladder carcinoma		

Before the Pulse

Cell Growth Medium	MEM, 10% Fetal Calf Serum, (+aminopterin, MPA, hypoxanthin, xanthine, thymidine) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	10 min. on ice, in Phosphate Buffered Sucrose
Wash Solution	Phosphate Buffered Sucrose		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	Room temperature		
Electroporation Medium*	Phosphate Buffered Sucrose (272 mM sucrose, 7 mM potassium phosphate, pH 7.4, 1 mM MgCl ₂)	Cuvette Gap	0.4 cm
Cell Density	1 to 50 x 10 ⁵ (5) cells / ml	Voltage	0.20 to 0.35 kV
Volume of Cells	0.4 ml	Field Strength	0.5 to 0.875 kV/cm
DNA Concentration	Not given		
DNA Resuspension Buffer	Not given	Capacitor	25 µF
Volume of DNA	5 to 20 µg /ml	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	7 to 12 msec
Outgrowth Medium	As above		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	Not given
Length of Incubation	Stable transfectants
Selection Method or Assay Used	MPA
Electroporation Efficiency	1 to 10 x 10 ⁶ (6) cells
Per Cent Survival	Not given

Name of Submitter Mikael Widersten, Ph.D. Student

Institution Address BMC
Biochemistry Dept
Box 576
S-75123 Uppsala, SWEDEN

Survey Number

153

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, adherent, suspension	Molecules Electroporated	DNA: linear & supercoiled: Rsv neo, Rsv gal, myc CAT, IRF CAT, fos CAT, hsp CAT, etc.
Species Used	Monkey, COS, kidney cells; Rat, N62 T cells; Mouse, mammary epithelial cells		

Before the Pulse

Cell Growth Medium	Fischer's + 10% Horse Serum + 10% Fetal Calf Serum (GIBCO/BRL)	Growth Phase at Harvest	None
		Pre-pulse Incubation	0 to 10 min, room temperature

Wash Solution

The Pulse	Instruments Used	Gene Pulser® apparatus & Capacitance
------------------	-------------------------	--------------------------------------

Electroporation Temperature	Room temperature		
Electroporation Medium*	Hepes buffered saline or Fischer's + 10% Horse Serum + 10% Fetal Calf Serum	Cuvette Gap	0.4 and 0.2 cm
Cell Density	1 x 10 ⁷ (7) cells / 0.4 ml	Voltage	0.25, 0.35, 0.45 kV
Volume of Cells	0.4 ml to 0.8 ml	Field Strength	625, 875, 1125kV/cm
DNA Concentration	80 µg/0.3 ml Hepes Buffered Saline		
DNA Resuspension Buffer	Not given	Capacitor	125, 250, 960 µF
Volume of DNA	< 50 µl	Resistor	(Pulse Controller) Ω none

After the Pulse

Outgrowth Medium	Fischer's + 10% Horse Serum + 10% Fetal Calf Serum	Time Constant	2 to 30 msec
-------------------------	--	----------------------	--------------

Outgrowth Temperature	Not given	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES, pH 7.2, 150 mM NaCl, 5 mM CaCl ₂	
Length of Incubation	Not given		
Selection Method or Assay Used	G418, 400 µg / ml		
Electroporation Efficiency	poor		
Per Cent Survival	40 to 50%		

Name of Submitter	Li-yuan Yu-Lee, Assistant Professor
--------------------------	-------------------------------------

Institution Address	Baylor College of Medicine Medicine Department 1 Baylor Plaza Houston, Texas 77030	Survey Number	154
----------------------------	---	----------------------	-----

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, adherent, suspension,	Molecules Electroporated	DNA: circular, 9.2kB, expression vectors
Species Used	Monkey, Vero, kidney cells; Human, C-4I, cervical carcinoma cells		

Before the Pulse

Cell Growth Medium	M199 (GIBCO/BRL, Sigma)	Growth Phase at Harvest	3 x 10 ⁶ (6) cells / ml
Wash Solution	M199	Pre-pulse Incubation	None

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4cm
Electroporation Medium*	M199	Voltage	0.250 kV
Cell Density	3 x 10 ⁶ (6) cells / ml	Field Strength	0.625 kV/cm
Volume of Cells	0.3 ml / pulse	Capacitor	500 µF
DNA Concentration	100 µg DNA / pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	distilled water	Time Constant	Not given
Volume of DNA	5 µl		

After the Pulse

Outgrowth Medium M199 - DMEM /F12,5% FBS,5% serum supplement, 1% pen/ strep, 400 µg G418

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37°C
Length of Incubation	24 hours recovery before selection
Selection Method or Assay Used	G418
Electroporation Efficiency	1.3 / µg DNA
Per Cent Survival	Not given

Name of Submitter Cherrilee Steele

Institution Address UTHSC-DB
Microbiology Dept.
P.O. Box 20068
Houston, Texas 77225

Survey Number
155

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, adherent, suspension	Molecules Electroporated	DNA: Plasmid DNA, 5 kB -12 kB
Species Used	Human, 293, kidney cells; Hamster, CHO, ovary cells		

Before the Pulse

Cell Growth Medium	Hamms F-12 / DMEM (50:50) +10% Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
Wash Solution	Phosphate Buffered Saline	Pre-pulse Incubation	Change to Hamms F-12 / DMEM (50:50) media 24 hours prior to pulse.

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature	Cuvette Gap	0.4 cm
Electroporation Medium*	High glucose/DMEM (Best we tested)	Voltage	0.245 kV
Cell Density	3 x 10 ⁶ (6) cells / ml	Field Strength	0.098 kV / cm
Volume of Cells	800 µl	Capacitor	960 µF
DNA Concentration	2 µg / ml	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	TE		
Volume of DNA	2 µl		

After the Pulse

Time Constant Not given

Outgrowth Medium	Hams F-12 / DMEM (50:50) media +10% Fetal Bovine Serum
-------------------------	--

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	<p>We are currently testing amplification procedures with linearized plasmid [See: Barsou, M., <i>DNA and Cell Biology</i> v.9 (4): 293-300]. 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</p>
Length of Incubation	20 min	
Selection Method or Assay Used	G418	
Electroporation Efficiency	5%	
Per Cent Survival	70 to 80%	

Name of Submitter George Cachiares

Institution Address Genentech Inc.
Molecular Biology
460 Pt. San Bruno
San Francisco, CA 94080

Survey Number

156

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, adherent, suspension	Molecules Electroporated	DNA: <i>ccc Bam Z</i> (5 kB), <i>EcoA</i> cosmid > 40 kB, <i>SalA</i> cosmid > 40 kB
Species Used	Human, B lymphomas: BJAB, P3HR-1, B95-8; Rat-1		

Before the Pulse

Cell Growth Medium	RPMI-10 (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log
		Pre-pulse Incubation	10 min.
Wash Solution	None		

The Pulse

Instruments Used Gene Pulser® apparatus, Capacitance

Electroporation Temperature	4°C		
Electroporation Medium*	RPMI-10 + 15% serum	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (6) or 10 (7) per 350 µl	Voltage	0.20 to 0.25 kV
Volume of Cells	350 µl	Field Strength	0.08 to 0.1 kV/cm
DNA Concentration	10 to 50 µg / cuvette		
DNA Resuspension Buffer	RPMI-10 or Phosphate Buffered Saline (PBS) or water	Capacitor	960 µF
Volume of DNA	10 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	28 to 40 msec
Outgrowth Medium	Not given		

Outgrowth Temperature	37°C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. <i>PNAS</i> 88:1546-1550 (1991). PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄	
Length of Incubation	Variable		
Selection Method or Assay Used	Protein expression by immunoblot, immunofluorescence, generation of recombinants.		
Electroporation Efficiency	> 20%		
Per Cent Survival	Variable %		

Name of Submitter Sankar Swaminathan, M.D.

Institution Address Harvard Medical School
Brigham & Women's Hospital
Dept of Infectious Diseases
20 Shattuck St.
Boston, MA 02115

Survey Number
157

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, adherent, suspension	Molecules Electroporated	DNA: circular plasmid recombinants (pMAM (8.8 kB + insert), pATH (9.0 kB) pGem (4.0 kB) pMSG (8.8 kB) pM6 (9.7 kB)
Species Used	Human, HeLa, cervical (C41); Human epithelial cells; Monkey, Vero, kidney.		

Before the Pulse

Cell Growth Medium	DMEM, MEM, M199; (10% Fetal Bovine Serum + 1% antibiotic) (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Early log
		Pre-pulse Incubation	Ice, 10 min.

Wash Solution Phosphate Buffered Saline (cold)

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Chilled		
Electroporation Medium*	Same as growth media	Cuvette Gap	0.4 cm
Cell Density	3 x 10 ⁶ (6) cells / ml	Voltage	0.250 kV
Volume of Cells	400 µl	Field Strength	0.625 kV/cm
DNA Concentration	10 to 100 µg / sample	Capacitor	500 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	1 µl to 50 µl	Time Constant	20 to 25 msec
After the Pulse			
Outgrowth Medium	Same as growth media		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	Brief ice, then 37°C
Length of Incubation	16 to 48 hours prior to selection
Selection Method or Assay Used	G418
Electroporation Efficiency	1 x 10 ⁻³ to 10 ⁻⁹ depending on cell type and plasmids used.
Per Cent Survival	50%

Name of Submitter Dr. Cherrilee Steele

Institution Address Univ. of Texas Health Sciences Center Microbiology / Dental Branch
P.O. Box 20068 / John Freeman Ave.
Houston, TX 77225

Survey Number
158

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, adherent, suspension	Molecules Electroporated	DNA: pTS1-envIF, 11 kB, supercoiled
Species Used	Mouse, NIH/3T3, embryo; Human T-cell line, (PEER)		

Before the Pulse

Cell Growth Medium	MDEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Not given
		Pre-pulse Incubation	MDEM
Wash Solution	No		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	MDEM	Cuvette Gap	0.4 cm
Cell Density	75%	Voltage	0.270 kV
Volume of Cells	0.8 ml	Field Strength	0.675 kV/cm
DNA Concentration	1 µg / µl		
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	960 µF
Volume of DNA	35 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	12.0 msec
Outgrowth Medium	MDEM		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	25 °C
Length of Incubation	10 min.
Selection Method or Assay Used	G418
Electroporation Efficiency	Very good
Per Cent Survival	80%

Name of Submitter Dr. Yuejin Yu, Postdoctoral Fellow

Institution Address University of Texas
M.D. Anderson Cancer Center
Science Park Research Division
Park Road 1C, P.O. Box 389
Smithville, Tx 78957

Survey Number

159

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, adherent	Molecules Electroporated	DNA: plasmids and cosmids, 40 kb and smaller.
Species Used	Mouse, 3T3, embryo; Human, fibroblast; primary cell lines; Monkey, Vero, kidney cells.		

Before the Pulse

Cell Growth Medium	DMEM + 10% Nu serum (GIBCO/BRL, Sigma, Flow Labs)	Growth Phase at Harvest	Passaged 24 hours prior to pulse.
		Pre-pulse Incubation	minimum
Wash Solution	Not given		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C		
Electroporation Medium*	Opti medium (BRL)	Cuvette Gap	0.4 cm
Cell Density	10 (7) cells / pulse	Voltage	0.220 kV
Volume of Cells	0.4 ml	Field Strength	0.55 kV/cm
DNA Concentration	10 to 20 µg / pulse		
DNA Resuspension Buffer	20 µl TE (10 mM Tris, 1 mM EDTA, pH 8.0) / pulse	Capacitor	960 µF
Volume of DNA	20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	30 msec
Outgrowth Medium	DMEM + 10% Nu serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	24 hours
Selection Method or Assay Used	b-gal expression: X-gal staining. Replication of Herpes origin of replication between S plasmids
Electroporation Efficiency	30 to 50% cells transfected
Per Cent Survival	60%

Name of Submitter Dr. Marie Masse

Institution Address Stanford University
Microbiology and Immunology
Stanford, CA 94305

Survey Number

160

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	cDNA's : supercoiled and linear.
Species Used	Human, HepG2, hepatoma; Monkey, COS-7, kidney.		

Before the Pulse

Cell Growth Medium	RPMI, 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Per Bio-Rad protocol
		Pre-pulse Incubation	Per Bio-Rad protocol
Wash Solution	Per Bio-Rad protocol (see Gene Pulser Instruction Manual)		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	0 °C (ice)		
Electroporation Medium*	Phosphate Buffered Saline	Cuvette Gap	0.2 cm
Cell Density	10 (7) cells / ml	Voltage	0.30 kV
Volume of Cells	0.4 ml	Field Strength	1.5 kV/cm
DNA Concentration	10 to 20 µg		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	1 µl to 10 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	Not given
Outgrowth Medium	RPMI /10% Fetal Calf Serum		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	Not given
Selection Method or Assay Used	FACS
Electroporation Efficiency	100 transfectants/ µg DNA
Per Cent Survival	30 to 50%

Name of Submitter J.S. Maltor

Institution Address Tulane Medical Center
Pathology Dept.
1430 Tulane Avenue
New Orleans, LA 70112

Survey Number

161

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, adherent	Molecules Electroporated	DNA: Recombinant pMAM neo vectors (8.8 to 9.2 kB) - confirm via expression (Northern blot) immunoblotting, PCR & Southern blot
Species Used	Human, squamous cell carcinoma lines, oral & cervical; Monkey, Vero, kidney cells.		

Before the Pulse

Cell Growth Medium	M199, MEM, DMEM/F12 (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Mid-log phase
		Pre-pulse Incubation	Ice, 10 min., in media with DNA.
Wash Solution	Ice-cold Phosphate Buffered Saline, pH 7.2		

The Pulse

Electroporation Temperature	4 °C	Instruments Used	Gene Pulser® apparatus & Capacitance
Electroporation Medium*	Growth media	Cuvette Gap	0.4 cm
Cell Density	1 x 10 ⁷ (7) cells / ml	Voltage	0.250 kV
Volume of Cells	3 x 10 ⁶ (6) cells / pulse	Field Strength	0.625 kV/cm
DNA Concentration	Not given	Capacitor	500 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) none Ω
Volume of DNA	100 µl	Time Constant	8 to 12 msec
After the Pulse			
Outgrowth Medium	Growth media, plus or minus G418		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37°C
Length of Incubation	at least 10 days
Selection Method or Assay Used	G418
Electroporation Efficiency	6 x 10 ⁻⁵ (-5) to 9 x 10 ⁻⁸ (-8); depends on cell line used
Per Cent Survival	50 %

Name of Submitter C. Steele

Institution Address UTHSC - DB
Dept of Microbiology
P.O. Box 20068
Houston, TX 77225

Survey Number

162

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, adherent	Molecules Electroporated	DNA: varies, fibronectin β -gal, genomic DNA, CMUB, etc.
Species Used	Human, fibroblast; Human Hep3b2, hepatocyte; Mouse, L-cells.		

Before the Pulse

Cell Growth Medium	MEM + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log
Wash Solution	Phosphate Buffered Saline and Trypsin	Pre-pulse Incubation	5 minutes

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25 °C	Cuvette Gap	0.4 cm
Electroporation Medium*	MEM	Voltage	0.320 kV
Cell Density	1 to 10 x 10 (6) cells / pulse	Field Strength	0.8 kV/cm
Volume of Cells	500 μ l	Capacitor	500 μ F
DNA Concentration	40 to 100 μ g DNA per pulse	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Time Constant	Not given
Volume of DNA	20 to 40 μ l		
After the Pulse			
Outgrowth Medium	MEM		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	48 hours to 1 month
Selection Method or Assay Used	GPT and Neomycin
Electroporation Efficiency	60%
Per Cent Survival	50%

Name of Submitter Ranee Taylor - Student

Institution Address Baylor College of Medicine
Molecular Genetics
1 Baylor Plaza
Houston, Texas 77030

Survey Number

163

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: Antibody genes with selectable markers.
Species Used	Hybridomas		

Before the Pulse

Cell Growth Medium	DMEM + 15% Horse Serum + L- glutamine (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Mid- log
Wash Solution	Cold Phosphate Buffered Saline	Pre-pulse Incubation	10 minutes on ice.

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4°C	Cuvette Gap	0.4 cm
Electroporation Medium*	Phosphate Buffered Saline	Voltage	0.160 kV
Cell Density	10 (7) cells / ml	Field Strength	0.4 kV/cm
Volume of Cells	0.8 ml	Capacitor	960 µF
DNA Concentration	20 µg / ml	Resistor	(Pulse Controller) Ω none
DNA Resuspension Buffer	Phosphate Buffered Saline		
Volume of DNA	Not given	Time Constant	15 msec
After the Pulse			
Outgrowth Medium	DMEM + 15% Fetal Calf Serum + Selector		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C	
Length of Incubation	> 1 week	<i>Immunol.</i> , Jan 15, 1991.
Selection Method or Assay Used	G418 or mycophenolic acid	
Electroporation Efficiency	0.5 to 1 transfectants / µg	
Per Cent Survival	Not given	

Name of Submitter Dr. Richard Near

Institution Address Mass. General Hospital
Cardiology Dept.
Jackson 1422 Cardiology
Boston, MA 02114

Survey Number

164

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, adherent	Molecules Electroporated	DNA: linearized 8.5 kB plasmid
Species Used	Hybrid, mouse/human, A9 fibroblast, hybrid containing human chromosomes		

Before the Pulse

Cell Growth Medium	DMEM + 10% Fetal Calf Serum + mycophenolic acid + xanthine	Growth Phase at Harvest	10 (7) cells / ml
		Pre-pulse Incubation	10 min. on ice, (0°C)
Wash Solution	HEPES Buffered Saline (HBS)		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	HEPES Buffered Saline (HBS)	Cuvette Gap	0.4 cm
Cell Density	Approximately 10 (7) cells / ml	Voltage	0.250, 0.450 kV
Volume of Cells	0.8 ml	Field Strength	0.625, 1.125 kV/cm
DNA Concentration	10 µg		
DNA Resuspension Buffer	Not given	Capacitor	125, 250, 500, 960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	Not given
Outgrowth Medium	Not given		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C		
Length of Incubation	Not given		
Selection Method or Assay Used	Hygromycin B		
Electroporation Efficiency	Not given		
Per Cent Survival	Not given		

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

Name of Submitter Keith A. Rosenbach

Institution Address UMDNJ New Jersey Medical School
Microbiology & Molecular Genetics
185 South Orange Ave.
Newark, NJ 07103

Survey Number

165

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, pN2 (retroviral vector), pCDM8-CD34 (both linears and supercoil).
Species Used	Mouse, B6SutA, hematopoietic; Human, K562, HEL, HL 60 (leukemia lines)		

Before the Pulse

Cell Growth Medium	B6SutA: McCoy's 5A + supplements (15% FCS + 10% WEHI-CM) human lines: RPMI 1640 + 10% calf serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Log phase
		Pre-pulse Incubation	5 to 10 min., room temperature
Wash Solution	Dulbecco's PBS, no Ca++ or Mg++		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	Dulbecco's PBS, no Ca++ or Mg++	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (6) to 10 (7) / ml	Voltage	0.3 kV
Volume of Cells	0.8 ml	Field Strength	0.75 kV/cm
DNA Concentration	100 to 1000 ng / µl		
DNA Resuspension Buffer	TE Buffer (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	500 µF
Volume of DNA	10 to 50 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	5 to 10 msec

After the Pulse

Outgrowth Medium	B6: McCoy's 5A + supplements, Leukemia lines: RPMI 1640 + 10% calf serum
-------------------------	--

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37°C
Length of Incubation	24 to 48 hours
Selection Method or Assay Used	G418 (100 to 800 µg / ml); cell survival.
Electroporation Efficiency	Unknown
Per Cent Survival	Unknown

Name of Submitter Mayumi Yagi

Institution Address Seattle Biomed Research Inst.
4 Nickerson St.
Seattle, WA 98109

Survey Number

166

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	Proteins
Species Used	Monkey, CV-1, kidney; Mouse, 3T3, embryo; Mouse, p3x63AG8, myeloma; Human: HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	DMEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	50 to 60% confluent cells
		Pre-pulse Incubation	5 minutes
Wash Solution	(See notes)		

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	0 °C		
Electroporation Medium*	(See notes)	Cuvette Gap	0.4 cm
Cell Density	5 x 10 ⁶ (6) cells / ml	Voltage	0.5 to 0.75 kV
Volume of Cells	0.2 µl	Field Strength	1.25 to 1.875 kV/cm
DNA Concentration	Not given		
DNA Resuspension Buffer	Not given	Capacitor	25 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
		Time Constant	2 to 4 msec

After the Pulse

Outgrowth Medium Pore-Resealing Buffer (see notes) for 10 min. ,37° C; follow with DMEM plus 10% Fetal Calf Serum.

Outgrowth Temperature	37° C
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.
Ref: M.R. Mitchell et. al. (1988) *Experientia* **44**: 199-203.
M.R. Michel et. al. (1990) *J. Virol.* **64**: 5123-5131.
Elgizoli, M. et. al. (1989) *J. Virol.* **63**: 2921-2928.
Electroporation Buffer: 20 mM PIPES, pH 7; 128 mM K-glutamate, 5mM ATP, 5 mM GTP, 10 µm Ca-Acetate, 2 mM Mg-Acetate & amino acid concentration corresponding to that of DMEM media.
Pore-Resealing Buffer: 20 mM PIPES, pH 7.0, 128 mM K-glutamate, 10 µm Ca-Acetate, 2 mM Mg-Acetate

Name of Submitter Marcel R. Michel, Ph.D.

Institution Address Institute for Medical Microbiology
Molecular Biology
Friedbuehlstrasse 51
3010 Berne, Switzerland

Survey Number

167

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, adherent, suspension	Molecules Electroporated	DNA: linearized constructs
Species Used	Human, K562, chronic myeloid leukemia; Hamster, CHO, ovary; Hybrid, rat / mouse, MEL cells.		

Before the Pulse

Cell Growth Medium	RPMI 1640, 10% Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	3 x 10 ⁵ to 5 x 10 ⁵ cells / ml (suspension)
		Pre-pulse Incubation	10 min. at room temp in Hepes buffered saline with dextrose
Wash Solution	Phosphate Buffred Saline		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	HEPES buffered saline with dextrose	Cuvette Gap	0.4 cm
Cell Density	2 x 10 ⁷ cells / ml	Voltage	0.2 kV
Volume of Cells	0.4 ml	Field Strength	0.5 kV/cm
DNA Concentration	Not given		
DNA Resuspension Buffer	RPMI 1640 +10% Fetal Bovine Serum	Capacitor	960 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	25.0 msec
Outgrowth Medium	RPMI 1640+10% Fetal Bovine Serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH ₂ PO ₄ , 1.15g Na ₂ HPO ₄ HBS: 10mM HEPES,pH 7.2,150 mM NaCl, 5 mM CaCl ₂	
Length of Incubation	Not given		
Selection Method or Assay Used	G418		
Electroporation Efficiency	Not given		
Per Cent Survival	Not given		

Name of Submittor No name or address given

Institution Address

Survey Number

168

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: pGL-luciferase vector [Promega]containing b-globin promoter ; co-porated SV 40, b-gal.
Species Used	Human, K562, chronic myeloid leukemia; HeLa, epithelial carcinoma; HEL cells, eythroleukemia.		

Before the Pulse

Cell Growth Medium	RPMI + 5% Fetal Calf Serum + 5% DCS	Growth Phase at Harvest	2 to 5 x 10 (5) cells / ml
		Pre-pulse Incubation	10 min., ice
Wash Solution	Phosphate Bufferd Saline + 5% DCS		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	25°C		
Electroporation Medium*	Hepes Buffered Saline	Cuvette Gap	0.4 cm
Cell Density	4 x 10 (7) cells / ml	Voltage	0.300 kV
Volume of Cells	0.5 ml	Field Strength	0.75 kV/cm
DNA Concentration	50 µg per pulse		
DNA Resuspension Buffer	Not given	Capacitor	960 µF
Volume of DNA	50 µl	Resistor	(Pulse Controller) Ω none
		Time Constant	31 msec

After the Pulse

Outgrowth Medium	RPMI + 5% Fetal Calf Serum + 5% DCS
-------------------------	-------------------------------------

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	24 hours
Selection Method or Assay Used	luciferease, β-gal
Electroporation Efficiency	Not given
Per Cent Survival	50%

Name of Submittor Dr. C. Anthony Blau

Institution Address University of Washington- Med Center
Department of Medical Genetics
Room BB510 HSB
1959 Pacific Avenue, NE
Seattle , WA 98006

Survey Number

192

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: Bovine papilloma virus E1 in pML2d
Species Used	Mouse, C127, fibroblast, mammary tumor		

Before the Pulse

Cell Growth Medium	DMEM + 10 % Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	logarithmic
		Pre-pulse Incubation	10 minutes on ice in DMEM + 10 % Fetal Calf Serum + BES
Wash Solution	PBS (Phosphate Buffered Saline) and Electroporation Medium		

The Pulse

Electroporation Temperature	Room temperature		
Electroporation Medium*	DMEM + 10 % Fetal Calf Serum + BES	Cuvette Gap	0.4 cm
Cell Density	1.5 to 2.0 x10 ⁶ (6) cells/ pulse	Voltage	0.21 kV
Volume of Cells	0.25 ml	Field Strength	0.525 kV/cm
DNA Concentration	10 to 20 µg / pulse		
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	960 µF
Volume of DNA	25 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	70 to 80 msec
Outgrowth Medium	DMEM + 10% Fetal Calf Serum		

		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
Length of Incubation	3 days /transient; 2-3 weeks/ stable	
Selection Method or Assay Used	G418 resistance for stable transformation	
Electroporation Efficiency	20-30%	
Per Cent Survival	30-40%	

Name of Submitter Dr. Belyavski Michail

Institution Address Texas A&M University
Department of Medical Microbiology and Immunology
College Station, TX 77834

Survey Number
193

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 DNA:
Electroporated

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

Outgrowth Medium Not given

Time Constant 60 msec

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

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

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

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, adherent and suspension	Molecules Electroporated	DNA: 2 to 4 kB, supercoiled
Species Used	Rat, PC12,adrenal pheochromocytoma; Rat brain; Simian (monkey) COS, kidney cells;		

Before the Pulse

Cell Growth Medium	DMEM (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Stationary growth
		Pre-pulse Incubation	10 min.
Wash Solution	Dulbecco's PBS		

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	4 °C		
Electroporation Medium*	Dulbecco's PBS	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (6) cells/ ml	Voltage	0.250 kV
Volume of Cells	400 to 800 µl	Field Strength	0.625 kV/cm
DNA Concentration	20 to 200 µg / ml		
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Capacitor	25 to 500 µF
Volume of DNA	10 to 20 µl	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	0.6 to 16 msec
Outgrowth Medium	DMEM		

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	days
Selection Method or Assay Used	Binding assays / transients; G418 selection / stable
Electroporation Efficiency	varies
Per Cent Survival	varies

Name of Submitter Dr.Catherine Bitler

Institution Address Molecular Pharmacologist
SRI International
Neuroscience Department
333 Ravenswood
Menlo Park, CA 94025

Survey Number

195

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	Molecules Electroporated	DNA: pRC/CMV- Human 5-HT(10β) receptor construct, linearized. [See notes]
Species Used	Human: HeLa, epithelial carcinoma		

Before the Pulse

Cell Growth Medium	EMEM + 10 % Fetal Calf Serum, + Non-essential amino acids + Pen / Strep	Growth Phase at Harvest	Pre-confluent
		Pre-pulse Incubation	10 min on ice
Wash Solution	Phosphate Buffered Sucrose		

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature	about 4 °C		
Electroporation Medium*	Phosphate Buffer Sucrose (see Gene Pulser Instruction Manual)	Cuvette Gap	0.4 cm
Cell Density	5 x 10 (6) cells / 800 µl	Voltage	0.4 kV
Volume of Cells	400 µl	Field Strength	1.0 kV/cm
DNA Concentration	20 µg / pulse		
DNA Resuspension Buffer	Phosphate Buffered Sucrose	Capacitor	25 µF
Volume of DNA	400 µl	Resistor	(Pulse Controller) 200 Ω
		Time Constant	2.5 msec

After the Pulse

Outgrowth Medium Post pulse: (in cuvette) 10 min. on ice, then
EMEM + 10 % FCS,
+ non-essential a.a.+ Pen / Strep

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	37 °C
Length of Incubation	plated, split next day
Selection Method or Assay Used	G418 selection (800 µg/ml) 24 hours post pulse
Electroporation Efficiency	100 transformants / µg
Per Cent Survival	Not known

DNA: 7.2 kB linear ds DNA construct including 5.6 kB of the mammalian expression vector pRC/ CMV (Invitrogen) and 1.8 kB of human genomic DNA including the coding region of the 5-HT10b (5 HT-1B) serotonin receptor, linearized with *Bgl* II.

Phosphate Buffered Sucrose: 272 mM sucrose, 7 mM sodium phosphate, pH 7.4, 1 mM MgCl₂.

Ref: *Biophys.Biochem. Res. Comm.* **184**:752-759.

Name of Submitter Dr. Mark W. Hamblin

Institution Address GRECC 182B, Seattle VAMC &
Dept. of Psychiatry & Behavioral Science
Univ.of Washington School of Medicine
1660 S. Columbia Way
Seattle, WA 98108

Survey Number

196

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, adherent	Molecules Electroporated	DNA: supercoiled plasmid; luciferase reporter gene
Species Used	Rat, CA77, medullary thyroid carcinoma cell line		

Before the Pulse

Cell Growth Medium	DMEM/F12 (1:1) + 10 % Fetal Bovine Serum (GIBCO/BRL, Sigma)	Growth Phase at Harvest	Active growth, about 50 to 70 % confluent
Wash Solution	Dulbecco's Phosphate Buffered Saline (minus Ca++, Mg++)	Pre-pulse Incubation	None

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	Room temperature		
Electroporation Medium*	Dulbecco's Phosphate Buffered Saline (minus Ca++, Mg++)	Cuvette Gap	0.4 cm
Cell Density	4 to 5 x10 (6) cells / 800 µl	Voltage	0.22 kV
Volume of Cells	800 µl	Field Strength	0.55 kV/cm
DNA Concentration	5 to 20 µg		
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA)	Capacitor	960 µF
Volume of DNA	5 to 50 µl (usually less than 10 µl)	Resistor	(Pulse Controller) Ω none
After the Pulse		Time Constant	about 12 msec
Outgrowth Medium	DMEM/F12 (1:1) + 10 % Fetal Bovine Serum		

Outgrowth Temperature	37 °C	Relevant Publications and/or Comments Note: exponential values designated in parentheses. Reference: Tverberg, L.A. and Russo, A.F. 1992. <i>J. Biol. Chem.</i> 267 (5):17567-17573.	
Length of Incubation	usually 24 hours		
Selection Method or Assay Used	Luciferase assay, β-galactosidase assay		
Electroporation Efficiency	about 1 x 10 (4) / µg [20 to 40 % cells stained blue, b-gal, transient assay]		
Per Cent Survival	50 to 80 %		

Name of Submitter Andrew Russo

Institution Address University of Iowa
Department of Physiology
5-632 BSB
Iowa City, IA 52242

Survey Number
197



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000