

Electroprotocols Species List

Fungal/Yeast Cells	Survey Number(s)	Plant Cells	Survey Number(s)
<i>Aspergillus</i> spp	169	<i>Hedyotis corymbosa</i>	182
<i>Aspergillus nidulans</i>	170	<i>Lactuca sativa</i> (aka chirimen chisha)	183
<i>Candida maltosa</i>	171	Maize, Black mexican sweet.....	181
<i>Colletotrichum gloeosporioides</i> (a fungal phytopathogen)	172	Maize, protoplast, DeKalb XL82	206
<i>Cryptococcus neoformans</i> , ma5 mutants	173	<i>Nicotiana plumbaginifolia</i> ; protoplasts from leaf	184
<i>Dictyostelium discoideum</i>	174, 175	<i>Oryza sativa</i> , cv. Yamahouci or cv. Nihonbare	185
<i>Pichia pastoris</i>	205		
<i>Saccharomyces cerevisiae</i> , DC5U	178	Other Cell Types	Survey Number(s)
<i>Saccharomyces cerevisiae</i> , strain S288C; a,α and a/α.....	177, 179	Chicken, HD11, macrophage.....	186
<i>Saccharomyces cerevisiae</i> , SEY6210	180	Chicken, primary hepatocytes.....	188
<i>Schizosaccharomyces pombe</i>	176	Chicken,TS34 a6 L1, [LSCC HD2], erythroblast.....	187
		Hydra cells, Cnidaria.....	189
		<i>Leishmania</i> , all species within the genus	190
		<i>Trypanosoma brucei brucei</i> , AnTat 1.3A, (blood- stream forms); EA TRO 1125 (procyclic forms).....	191

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: *Aspergillus* genomic DNA, 3 to 13 kB.

Species Used *Aspergillus* spp.

Before the Pulse

Cell Growth Medium Czapek + requirement (ATCC#312) medium, Polypeptone -Dextrin medium.

Growth Phase at Harvest Protoplast

Pre-pulse Incubation None

Wash Solution 0.8 M Sorbitol

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature Room temperature, 25 °C

Electroporation Medium 1.1 M sorbitol

Cuvette Gap 0.4 and 0.2 cm

Voltage Not given

Cell Density 2 x 10⁷ / ml

Volume of Cells 2 x 10⁷ / ml

Field Strength 4 kV/cm

DNA Concentration 5 µg

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

Capacitor 25 µF

Volume of DNA Not given

Resistor (Pulse Controller) Not given.

Time Constant 3 to 7 msec

After the Pulse

Outgrowth Medium Czapek + 0.8 M Sorbitol

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 30 °C

Length of Incubation 7 days

Selection Method or Assay Used Can grow on minimal medium

Electroporation Efficiency 10 / µg DNA

Per Cent Survival 10%

Name of Submitter Shigeomi Ushijima, Ph.D.

Institution Address Kikkoman Corp.
Research and Development Division
399 Noda
Noda City, Chiba 278 JAPAN

Survey Number

169

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: integrative plasmid

Species Used *Aspergillus nidulans*

Before the Pulse

Cell Growth Medium Not given

Growth Phase at Harvest Not given

Pre-pulse Incubation Not given

Wash Solution Not given

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Not given

Electroporation Medium 1.2 M Sorbitol, 7mM NaPO4 (pH7.2), 1 mM MgSO4

Cuvette Gap 0.2 cm

Cell Density 4 x 10 (6) protoplasts / ml

Voltage 0.400 kV, 0.700 kV

Volume of Cells Not given

Field Strength 2.0 kV/cm, 3.5 kV/cm

DNA Concentration Not given

DNA Resuspension Buffer Not given

Capacitor 25 µF

Volume of DNA Not given

Resistor (Pulse Controller) Ω none

After the Pulse

Time Constant 5.2 msec /3.3 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 Not given

Electroporation Efficiency 3.5, 4.0 transformants / µg DNA

Per Cent Survival Not given

Name of Submitter Dr. D. Sanglard

Institution Address Swiss Federal Institute of Technology
Inst. for Biotechnologie
ETH - Hoenggerberg
8093 Zurich
Switzerland

Survey Number

170

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast
Species Used *Candida maltosa*

Molecules Electroporated DNA: pTRA11, episomal plasmid

Before the Pulse

Cell Growth Medium Not given

Growth Phase at Harvest Not given

Pre-pulse Incubation Not given

Wash Solution Not given

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Not given

Electroporation Medium 0.5M sucrose, 8mM Phosphate Buffer (pH7.2), 1mM MgCl₂

Cuvette Gap 0.4 cm

Cell Density 1 x 10⁶ (6) protoplasts / ml

Voltage 0.500 kV

Volume of Cells 800 µl

Field Strength 1.25 kV/cm

DNA Concentration 100 ng DNA

DNA Resuspension Buffer Not given

Capacitor 1 µF

Volume of DNA Not given

Resistor (Pulse Controller) Ω none

After the Pulse

Outgrowth Medium Not given

Time Constant 0.2 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

Electroporation Efficiency 160 transformants/µg DNA

Per Cent Survival Not given

Name of Submitter Dr. D. Sanglard

Institution Address Swiss Federal Institute of Technology
 Inst. for Biotechnologie
 ETH - Hoenggerberg
 8093 Zurich
 Switzerland

Survey Number

171

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: pHIS (6.7 kb) circular & linear, (Hygromycin B resistance).

Species Used *Colletotrichum gloeosporioides*
(a fungal phytopathogen)

Before the Pulse

Cell Growth Medium Clarified V8 juice

Growth Phase at Harvest Mycelium, 24 hours old.

Pre-pulse Incubation

Wash Solution Produce protoplasts per Tilbur *et. al.*, 26:205-221 (1983).

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature 0 ° C (ice)

Electroporation Medium 10% sucrose, 10 mM Tris, 2 mM MgCl₂

Cuvette Gap 0.4 cm

Cell Density 5 x 10⁷ protoplasts

Voltage 0.500-1.250 kV

Volume of Cells 0.8 ml

Field Strength 2 to 5 kV/cm

DNA Concentration 50 µg / 5 x 10⁷ protoplasts

DNA Resuspension Buffer 600 mM Sucrose, 10 mM CaCl₂, pH 7.5

Capacitor 1 µF

Volume of DNA 6.7 µl / pulse

Resistor (Pulse Controller) 200 Ω

After the Pulse

Time Constant 3 to 5 msec

Outgrowth Medium Czapeck's minerals (ATCC#312) + 10 mM sodium citrate, 1% casamino acids, 2% bacto-agar, 20% sucrose

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 26 °C

Length of Incubation 5 to 7 days

Selection Method or Assay Used Hygromycin B, 25 µg / ml

Mosel, A., Erwin, J.A.G., Manners, J.M. (1989) Transformation of the plant pathogen *Colletotrichum gloeosporioides*. Abstracts, 7th Australian Plant Pathology Society Conference, Brisbane, pg 58.

Electroporation Efficiency 0.1 to 0.3 transfectants per µg DNA

Per Cent Survival 10 to 30%

Name of Submitter Dr. J. Manners

Institution Address Dept. of Botany
Univ of Queensland
Queensland, 4072
AUSTRALIA

Survey Number

172

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: supercoiled on linear plasmids containing URA5 gene

Species Used *Cryptococcus neoformans*, ma5 mutants

Before the Pulse

Cell Growth Medium YEPD

Growth Phase at Harvest Logarithmic, O.D.(650) = ~ 1.

Pre-pulse Incubation None

Wash Solution 270 mM Sucrose, 1 mM MgCl₂, 10 mM Tris-HCl, pH 7.5, 4 mM DTT

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Room temperature

Electroporation Medium 270 mM Sucrose, 10 mM Tris HCl, pH 7.5 (no DTT)

Cuvette Gap 0.2 cm

Cell Density Cells concentrated, 100 fold

Voltage 0.470 kV

Volume of Cells 450 µl

Field Strength 2.35 kV/cm

DNA Concentration 0.1 to 1.0 µg, in 1 to 10 µl TE

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

Capacitor 25 µF

Volume of DNA 1 to 10 µl

Resistor (Pulse Controller) none Ω

After the Pulse

Time Constant 18 to 22 msec

Outgrowth Medium None

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 SD media (lacking uracil): 6.7 g yeast nitrogen base per liter without amino acids and 20 g glucose / liter

Note: the protocol described in the paper by Edman and Kwon-Chung, *Mol. & Cell. Biol.*, **10**:4538-4544 (1990) is not the one described above. The one above gives 10-100x greater efficiency.

Electroporation Efficiency 2000 transfectants / µl

Per Cent Survival Not tested

Name of Submitter Jeffrey C. Edman, Asst. Prof.

Institution Address Univ. of California
Hormone Research Institute
Box 0534
San Francisco, CA 94143-0534

Survey Number

173

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: circular, ranging from 6 kB to 20 kB.

Species Used *Dictyostelium discoideum*, strain AX4 and HUD205

Before the Pulse

Cell Growth Medium HL5 (see reference in notes)

Growth Phase at Harvest Late log phase

Pre-pulse Incubation 10 min, room temperature.

Wash Solution 10 mM NaPO₄, pH 6.1, 50 mM sucrose

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Room temperature

Electroporation Medium 10 mM NaPO₄, pH 6.1, 50 mM sucrose

Cuvette Gap 0.4 cm

Cell Density 3 x 10⁷ / ml

Voltage 0.60 kV

Volume of Cells 0.4 ml

Field Strength 1.5 kV / cm

DNA Concentration 1 µg / µl

DNA Resuspension Buffer TE (10 mM Tris, 1 mM EDTA, pH 7.4)

Capacitor 3 µF

Volume of DNA 1 to 10 µg DNA / 0.4 ml cells

Resistor (Pulse Controller) none Ω

After the Pulse

Time Constant Not given

Outgrowth Medium HL5

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 21 °C

Length of Incubation 18 to 24 hours

Selection Method or Assay Used G418

We use a modification of the technique described for this organism by Howard, Ahern, Firtel (1988) *Nucleic Acids Res.* **16**: 2613-2623.

Electroporation Efficiency 6 x 10² / µg DNA

*We also electroporate *E. coli* (strains JM109 and Sure™ cells) using the procedure more or less as described in the literature which accompanied the Pulse Controller: 2.5 kV, 200 Ω, 25 µF, 0.2 cm cuvettes.

Per Cent Survival 70 to 80%

Name of Submitter Dr. Joanne E. Hughes

Institution Address Utah State University
Biology Dept
Logan, Utah 84322-5500

Survey Number

174

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: supercoiled and linear

Species Used *Dictyostelium discoideum*

Before the Pulse

Cell Growth Medium HL5 (peptone, yeast extract, glucose)
(ATCC Media #671)

Growth Phase at Harvest 1×10^6 to 1×10^7 cells / ml

Pre-pulse Incubation 5 min.

Wash Solution HEPES Buffered Saline

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature 4 °C

Electroporation Medium HEPES Buffered Saline

Cuvette Gap 0.4 cm

Cell Density 5×10^6 / ml

Voltage 1.25 kV

Volume of Cells 1 ml

Field Strength 3.125 kV/cm

DNA Concentration 20 ng to 20 µg

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

Capacitor 25 µF

Volume of DNA Any volume

Resistor (Pulse Controller) Ω none. NOT

After the Pulse

Time Constant 0.5 to 0.7 msec

Outgrowth Medium HL5

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 22 °C

Length of Incubation overnight

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

Selection Method or Assay Used G418

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

Electroporation Efficiency 10^{-3} transformants / cell;
 6×10^3 transformants / µg DNA.

Per Cent Survival >90%

Name of Submitter David Knecht

Institution Address Univ of Connecticut
Dept of Molecular & Cell Biology
U-125
Storrs, CT 06269

Survey Number

175

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: plasmid, pHILD2 & D4, linearized, 8 to 10 kB.

Species Used *Pichia pastoris* GTS115

Before the Pulse

Cell Growth Medium Yeast Extract Potato Dextrose, YEPD, (DIFCO)

Growth Phase at Harvest O.D. (600) = 1.3

Pre-pulse Incubation 5 minutes, 4 °C

Wash Solution Cold water two times; then 1 M sorbitol, one time

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature 25 °C but sample & cuvette at 4 °C

Electroporation Medium 1 M sorbitol

Cuvette Gap 0.2 cm

Cell Density 300x concentration from harvest density

Voltage 1.5 kV

Volume of Cells 50 µl

Field Strength 7.5 kV/cm

DNA Concentration 0.5 to 2 µg / pulse

DNA Resuspension Buffer 1 M sorbitol

Capacitor 25 µF

Volume of DNA 1 to 5 µl

Resistor (Pulse Controller) 400 Ω

After the Pulse

Outgrowth Medium Minimal salts plus dextrose (MD)

Time Constant approximately 8.0 msec

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. This is essentially the method described for *Saccharomyces cerevisiae* by Becker and Guarente, *Methods in Enzymol.*, **194**,182-187(1991).

Outgrowth Temperature 30 °C

Length of Incubation 3 to 5 days

Selection Method or Assay Used Complimentation of histidine auxotrophy

Electroporation Efficiency approx. 1000 transformants / µg DNA

Per Cent Survival Not tested

Name of Submitter Carl W. Despreaux

Institution Address Hoffmann La Roche
Inflammation / Autoimmune Disease
Bldg. 102, Lab 509
Nutley, NJ 07110

Survey Number

205

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: YEp24, plasmid

Species Used *Saccharomyces cerevisiae*, DC5U

Before the Pulse

Cell Growth Medium YEPD (ATCC#1202/1245)

Growth Phase at Harvest OD(600) = 1.1 to 1.3

Pre-pulse Incubation 1 M Sorbitol

Wash Solution 2 x Water, 1 M Sorbitol (Becker and Guarente protocol - see notes).

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature 0 °C (ice)

Electroporation Medium 1 M Sorbitol

Cuvette Gap 0.2 cm

Cell Density 3 x 10⁸ / ml

Voltage 1.0 to 1.5 kV

Volume of Cells 50 to 100 µl

Field Strength 5.0 to 7.5 kV/cm; optimal: 6.25kV/cm

DNA Concentration 1 µg / ml

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

Capacitor 25 µF

Volume of DNA 0.5 µl (0.5 µg)

Resistor (Pulse Controller) 200 Ω

After the Pulse

Time Constant 4.0 to 5.0 msec

Outgrowth Medium Not given

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 30 °C

Length of Incubation 3 days

Reference: Becker, D., Guarante, L. *Methods in Enzymol.* **104**:182-187 (1991).

Selection Method or Assay Used Incubate in 1 M Sorbitol for ~ 15 minutes. Place on SD + histidine + leucine + 1 M Sorbitol

Electroporation Efficiency 1.2 x 10⁵ transfectants / µg DNA

Per Cent Survival 40%

Name of Submitter Terri Dunahay

Institution Address Solar Energy Research Institute
Biotechnology
1617 Cole Blvd.
Golden, CO 80401

Survey Number

178

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroported DNA: YEp351/352, pRS 303 - 316, YEp50, usually supercoiled, 6-9 kB.

Species Used *Saccharomyces cerevisiae*, strain S288C; a, α and a/ α

Before the Pulse

Cell Growth Medium YEPD (ATCC#1202/1245) or synthetic

Growth Phase at Harvest Log phase 80 to 100 Klett units

Pre-pulse Incubation 10 mM Tris-HCl, pH 7.5, 1 M Sorbitol

Wash Solution Water

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature Room temperature

Electroporation Medium YPD, 1 M Sorbitol

Cuvette Gap 0.1 cm

Cell Density Concentrated 100 x

Voltage 0.55 kV

Volume of Cells 60 to 100 μ l

Field Strength 5.5 kV/cm

DNA Concentration 1 μ g / μ l

DNA Resuspension Buffer YEPD or SOC - yeast

Capacitor 25 μ F

Volume of DNA 7 to 10 μ l

Resistor (Pulse Controller) 600 Ω

After the Pulse

Time Constant 5 to 6 msec

Outgrowth Medium synthetic plates, - URA, -LEU, -HIS; No soft agar.

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 30 ° C, sometimes 24° C

Length of Incubation 48 hours

Selection Method or Assay Used -URA, -LEU, -HIS, or combinations of them.

Electroporation Efficiency 1000 to 3000 transfectants / μ g DNA

Per Cent Survival Not known

We are not interested in electroporation as such, but it is a very convenient method to introduce DNA into yeast cells. Electroporation is much less time consuming than the other methods available and also easier to perform. Room temperature is used because otherwise the time constant becomes too high and you get fewer transformants. This may also be used with frozen yeast cells but then the efficiency drops a lot.

Name of Submitter Stefan Astrom, BSc

Institution Address University of Umea
Dept. of Microbiology
S-90187 Umea, SWEDEN

Survey Number

177

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: pUC19 and Bluescript™ -based plasmids.

Species Used *Saccharomyces cerevisiae* - lines derived from S288C

Before the Pulse

Cell Growth Medium YEPD (ATCC#1202/1245)

Growth Phase at Harvest 0.5 to 1.0 at O.D.600; ~100 mls log phase culture concentrated 200x, in water.

Pre-pulse Incubation None required

Wash Solution Water

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature Room temperature

Electroporation Medium Water

Cuvette Gap 0.2 cm

Cell Density 6 x 10 (10) cells / ml

Voltage 0.6 kV

Volume of Cells 40 to 50 µl

Field Strength 3.0 kV/cm

DNA Concentration Total DNA = 0.2 to 1.0 µg

Capacitor 25 µF

DNA Resuspension Buffer Not given

Resistor (Pulse Controller) 200 Ω

Volume of DNA ≤5µl

Time Constant Not given

After the Pulse

Outgrowth Medium Add water to final vol. 150-200 µl. Plate on selective synthetic media (SD)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. Washed and concentrated cells can be stored by adding 80% glycerol to a final concentration of 20% and freezing at -80°C. To electroporate thawed cells: pellet cells, remove glycerol containing supernatant, wash cells in 0.5-1.0 ml water, and resuspend cells in water to original volume. Efficiency is not affected significantly by freezing but you must remove glycerol. **Comment:** removing glycerol after thawing cells may help by simply removing lysed cell contents that would alter media conductivity - this could impact efficiency (alters the time constant, t).

Outgrowth Temperature 30 °C

Length of Incubation 2 days

Selection Method or Assay Used Amino acid or nucleotide dropout

Electroporation Efficiency 2 to 10,000 transfectants / µg DNA

Per Cent Survival Not given

Name of Submitter Joachim Li

Institution Address Univ of California /San Francisco
Dept. of Biochem & Biophysics, S-960
513 Parnassus
San Francisco, CA 94143-0448

Survey Number

179

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: YCp50 plasmid DNA (yeast centrameric shuttle vector)

Species Used *Saccharomyces cerevisiae*, SEY6210

Before the Pulse

Cell Growth Medium YPDA: 10% yeast extract, 20% bacto peptone, 20% glucose, 2 µg / ml adenine

Growth Phase at Harvest 2.0 x 10⁷ cells / ml

Pre-pulse Incubation E-buffer (see notes) on ice for at least 5 minutes

Wash Solution E-buffer (see notes)

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Room temperature

Electroporation Medium E-buffer (see notes)

Cuvette Gap 0.2 cm

Cell Density 2.0 x 10⁹ cells / ml

Voltage 0.54 kV

Volume of Cells 50 µl

Field Strength 2.7 kV/cm

DNA Concentration 10 to 100 µg

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

Capacitor 25 µF

Volume of DNA 2 µl

Resistor (Pulse Controller) none Ω

After the Pulse

Time Constant 10 to 20 msec

Outgrowth Medium YPDA

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

E. Meilhoc *et. al.* Biotechnology **8**: 223-227, 1990.

Wash solution:

1. 10 mM Tris pH 8.0, 25 mM DTT in YPDA at 30° C for 10 min.
2. E-buffer: 10 mM Tris HCl pH 7.5, 270 mM Sucrose, 1mM MgCl₂

Outgrowth Temperature 30 °C

Length of Incubation 2 hours

Selection Method or Assay Used YMM - ura (4 days) [yeast introgen base without amino acids, Difco]

Electroporation Efficiency 2.5 x 10⁵ transfectants / µg DNA

Per Cent Survival 50%

Name of Submitter Martin Latterich

Institution Address University of Durham
Biological Sciences
South Road
Durham DH13LE, ENGLAND

Survey Number

180

Gene Pulser® Electroprotocol

Cell Type Fungal / Yeast

Molecules Electroporated DNA: supercoiled plasmids, linear fragments for integration.

Species Used *Schizosaccharomyces pombe*

Before the Pulse

Cell Growth Medium YE or dropout media (recipes in paper)

Growth Phase at Harvest 1 x 10⁷ (7) cells / ml

Pre-pulse Incubation none

Wash Solution 1.2 M sorbitol (ice cold, filter sterilized)

The Pulse

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Temperature (Ice cold) 0 °C

Electroporation Medium 1.2 M sorbitol

Cuvette Gap 0.2 cm

Cell Density 1 x 10⁴ (4) cells / ml

Voltage 2.25 kV

Volume of Cells 200 µl

Field Strength 1.125 kV/cm

DNA Concentration 1 ng to 1 µg DNA per pulse

DNA Resuspension Buffer 1.2 M sorbitol

Capacitor 25 µF

Volume of DNA <10 µl

Resistor (Pulse Controller) 200 Ω

After the Pulse

Outgrowth Medium SD + necessary nutrients

Time Constant 5 msec

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Reference: Prentice, H. *Nucleic Acid Res.* **20** (3):621.

Outgrowth Temperature 30 °C

Length of Incubation 4 to 6 days

Selection Method or Assay Used auxotrophy

Electroporation Efficiency 1 x 10⁵ (5) to 1 x 10⁶ (6) for autonomous plasmids

Per Cent Survival ~50% (seeTable 1 in reference)

Name of Submitter Holly Prentice

Institution Address Mass Gen Hosp./ Harvard Med School
Molecular Biology / Genetics Dept.
Wellman 10
Boston, MA 02114

Survey Number

176

Gene Pulser® Electroprotocol

Cell Type Plant, protoplasts

Molecules Electroporated DNA: pROA93, 16.2 kB, circular.

Species Used *Hedyotis corymbosa*

Before the Pulse

Cell Growth Medium Murashige & Skoog's medium + 2% sucrose + 2 mg/l, 2,4-D, 1 g/l casein hydrolysate (GIBCO/BRL)

Growth Phase at Harvest Not given

Pre-pulse Incubation 10 min on ice

Wash Solution Not given

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature 20 °C

Electroporation Medium Phosphate HEPES buffered saline

Cuvette Gap 0.4 cm

Cell Density 1 x 10⁶ (6) protoplasts

Voltage 0.2 to 0.4 kV

Volume of Cells 1 µl

Field Strength 0.250 to 1.125 kV/cm

DNA Concentration 40 µg/ml

DNA Resuspension Buffer Not given

Capacitor 10 to 960 µF

Volume of DNA 5 to 10 µl

Resistor (Pulse Controller) none Ω

After the Pulse

Time Constant Not given

Outgrowth Medium MS medium + 6% glucose + 2 µg/l 2,4-D + 0.1% casein hydrolysate

Relevant Publications and/or Comments
Note: exponential values designated in parentheses.

Outgrowth Temperature 25 °C

Length of Incubation 2 days

Selection Method or Assay Used CAT

Electroporation Efficiency Good

Per Cent Survival Not given

Name of Submitter Dr. Pua Eng Chong

Institution Address National University of Singapore
 Inst. of Molecular and Cell Biology
 Kent Ridge
 Singapore 0511

Survey Number
 182

Gene Pulser® Electroprotocol

Cell Type Plant Molecules DNA: pCaMVneo [NPT II]
 Species *Lactuca sativa* (aka *chirimen chisha*)
 Used Electroporated

Before the Pulse

Cell Growth Medium MS (Murashige and Skoog) (GIBCO/BRL) Growth Phase at Harvest Not given
Pre-pulse Incubation Ice for 10 minutes
 Wash Solution Leaf mesophyll protoplasts

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature	0 °C		
Electroporation Medium	5 mM MES, 70 mM KCl, 0.3 M Mannitol, pH 5.8	Cuvette Gap	0.4 cm
Cell Density	2 x 10 (6) / ml	Voltage	0.250 kV
Volume of Cells	500 µl	Field Strength	0.625 kV/cm
DNA Concentration	100 µg / ml	Capacitor	250 µF
DNA Resuspension Buffer	TE (10 mM Tris, 1 mM EDTA, pH 8.0)	Resistor	(Pulse Controller) none Ω
Volume of DNA	50 µg	Time Constant	13.8 msec

After the Pulse

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

Relevant Publications and/or Comments

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

Outgrowth Temperature 24 °C
 Length of Incubation 2 months
 Selection Method or Assay Used G418 Geneticin (5 to 20 µg / ml)
 Electroporation Efficiency only one
 Per Cent Survival 10 to 30 %

Name of Submitter Takashi Ishibashi

Institution Address Nippon Petroleum Refining Co. Ltd.
 Petroleum Section, Bio Group
 766, Higashi-Toyoi
 Kudamatsu-shi, Yamaguchi 744
 JAPAN

Survey Number

183

Gene Pulser® Electroprotocol

Cell Type	Plant	Molecules Electroporated	DNA: ds DNA mRNA
Species Used	Maize, Black mexican sweet		

Before the Pulse

Cell Growth Medium	BMS culture medium (see reference in notes)	Growth Phase at Harvest	Log
		Pre-pulse Incubation	5 min. at 42° C
Wash Solution	Not given		

The Pulse

Instruments Used Promega X-Cell

Electroporation Temperature	0 °C		
Electroporation Medium	10 mM HEPES 1mM CaCl ₂	Cuvette Gap	0.4 cm
Cell Density	2 x 10 (6) / ml	Voltage	0. 200 kV
Volume of Cells	0.8 ml	Field Strength	0.5 kV/cm
DNA Concentration	10 µg / ml		
DNA Resuspension Buffer	Not given	Capacitor	1250 µF
Volume of DNA	Not given	Resistor	(Pulse Controller) Do not calculate
		Time Constant	Do not calculate

After the Pulse

Outgrowth Medium Plant growth media

Outgrowth Temperature	10 min. at 4 °C, then 25° C
Length of Incubation	up to 3 days
Selection Method or Assay Used	b-glucuronidase (GUS), luciferase
Electroporation Efficiency	50%
Per Cent Survival	90%

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
All of our recent work uses electroporation.
Ref: Fromm, *et. al., Meth. Enzymol.* **153:**351-366 (1987).

Name of Submitter Dr. Virginia Walbot

Institution Address Stanford University
Dept. of Biological Sciences
Stanford, CA 94305-5020

Survey Number
181

Gene Pulser® Electroprotocol

Cell Type Plant, protoplast

Molecules Electroported DNA: plasmids with VP1 maize gene, supercoiled; several promoter / GUS fusions that are activated by VP1.

Species Used Maize cell protoplast, DeKalb XL82 (scutellum), Mpp

Before the Pulse

Cell Growth Medium N6 medium

Growth Phase at Harvest Mid-log, three days after transfer

Pre-pulse Incubation DNA plus cells held on ice for 10 minutes prior to electroporation

Wash Solution Protoplasts made by digesting with enzymes, then washed.

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature 25 °C, but sample pre-chilled

Electroporation Medium Not given

Cuvette Gap 0.4 cm

Cell Density 4 x 10⁶ cells / ml

Voltage 0.2 kV

Volume of Cells 1 ml ******(SEE NOTES)

Field Strength 0.5 kV/cm

DNA Concentration 20 to 50 µg / pulse

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

Capacitor 960 µF

Volume of DNA 20 to 50 µl / pulse

Resistor (Pulse Controller) none

After the Pulse

Time Constant 14 to 16 msec, average

Outgrowth Medium KMØ medium

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
******Maximum volume for 0.4 cm cuvettes is 0.8 ml; greater volumes will deliver a non-uniform pulse to sample.
 Ref: McCarty, *et.al.*, *Cell*, **66**:895-905(1991).

Outgrowth Temperature 25 °C

Length of Incubation 40 hours

Selection Method or Assay Used Fluorescence assay for GUS, luminescence for luciferase

Electroporation Efficiency Not done

Per Cent Survival 70 to 80%

Name of Submitter Leonard Rosenkrans/ Dr. Don McCarty

Institution Address University of Florida
 Department of Vegetable Crops
 1244 Fifield Hall
 Gainesville, FL

Survey Number

206

Gene Pulser® Electroprotocol

Cell Type Plant

Molecules Electroporated DNA: plasmid pTZ 18U with a 3 kB insert (total: 6 kB), circular.

Species Used *Nicotiana plumbaginifolia*; protoplasts from leaf

Before the Pulse

Cell Growth Medium 5 mM CaCl₂; 0.5 M sucrose, 0.3% Macerozyme R-10, 0.3% cellulose "Onozuka" R-10, pH 4.7 (enzymes-Yokult Honsha Co.,Ltd.)

Growth Phase at Harvest Not given

Pre-pulse Incubation 2 hours at 4°C

Wash Solution 4 mM CaCl₂, 80mM KCl, 8.0% Mannitol, 2 mM Na₂P0₄, pH 7.2

The Pulse

Instruments Used Gene Pulser® apparatus & Capacitance

Electroporation Temperature 2 hr. at 4 °C

Electroporation Medium Same as wash solution

Cuvette Gap 0.2 cm

Cell Density 10 (6) / ml

Voltage 0.16 kV

Volume of Cells 300 µl

Field Strength 0.8 kV/cm

DNA Concentration 3 µg / µl

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

Capacitor 125 µF

Volume of DNA 10 µg DNA / pulse, 3 µl.

Resistor (Pulse Controller) none Ω

After the Pulse

Time Constant 2.8 msec

Outgrowth Medium 9% Mannitol, 3% sucrose, 100 µg/ml cefotamine (antibiotic)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature 25 °C

Length of Incubation 36 hours

I am no longer a member of this laboratory, but work is in progress there.

Selection Method or Assay Used Transient expression assays

Electroporation Efficiency Not determined

Per Cent Survival Not determined

Name of Submitter Dupriez Vincent, Student

Institution Address Universite Cottidique deLouvain
Fac. Agronomie, Unite FYSA
Place Croix du Sud No. 2 bte 20
Louvain-La-Neuve BELGIUM 1348

Survey Number

184

Gene Pulser® Electroprotocol

Cell Type Plant, suspension

Molecules Electroporated DNA: pCH (pUC vector with hygromycin resistant gene)

Species Used *Oryza sativa*, cv. Yamahouci or cv. Nihonbare

Before the Pulse

Cell Growth Medium AA medium (Amino Acid Medium)

Growth Phase at Harvest Log phase

Pre-pulse Incubation 0.5 mM MES, 70 mM KCl, 4 mM CaCl₂, 0.36 M Mannitol

Wash Solution 0.5 mM MES, 70 mM KCl, 4 mM CaCl₂, 0.36 M Mannitol

The Pulse

Instruments Used Gene Pulser® apparatus , Capacitance

Electroporation Temperature Room temperature

Electroporation Medium 0.5 mM MES, 70 mM KCl, 4 mM CaCl₂, 0.36 M Mannitol

Cuvette Gap 0.4 cm

Cell Density 10 (6) / ml

Voltage 0.25 kV

Volume of Cells 500 µl

Field Strength 0.625 kV/cm

DNA Concentration Not given

Capacitor 250 µF

DNA Resuspension Buffer Not given

Volume of DNA 50 µl

Resistor (Pulse Controller) 200 Ω

After the Pulse

Time Constant 10 to 20 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 Not given

Electroporation Efficiency Not given

Per Cent Survival Not given

Name of Submitter Dr. Seiichi Toki

Institution Address Hokkaido Univ College of Science
Biology, Plant structural studies
Sapporo-shi, kita-ku, 10 jo nishi 8 chome
JAPAN 7060

Survey Number

185

Gene Pulser® Electroprotocol

Cell Type Other Cell Types

Molecules Electroported DNA: linearized DNA used for stable transfections.

Species Used Chicken, HD11, macrophage

Before the Pulse

Cell Growth Medium DMEM, 8% Fetal Calf Serum (FCS), 2% Chicken Serum (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 0.270 kV

Volume of Cells 0.5 ml

Field Strength 0.675 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

Outgrowth Medium DMEM, 8% Fetal Calf Serum (FCS), 2% Chicken Serum

Time Constant 28 msec

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH₂PO₄, 1.15g Na₂HPO₄

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

186

Gene Pulser® Electroprotocol

Cell Type Other Cell Types

Molecules Electroported DNA: supercoiled DNA used for transient transfections.

Species Used Chicken, primary hepatocytes

Before the Pulse

Cell Growth Medium Not given

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 0.250 kV

Volume of Cells 0.5 ml

Field Strength 0.625 kV/cm

DNA Concentration 10 µg / pulse

Capacitor 960 µF

DNA Resuspension Buffer Not given; pulse volume: 0.8 ml

Resistor (Pulse Controller) Ω none

Volume of DNA Not given; pulse volume: 0.8 ml

Time Constant 25 msec

After the Pulse

Outgrowth Medium Not given

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

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 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

188

Gene Pulser® Electroprotocol

Cell Type Other Cell Types

Molecules Electroported DNA: supercoiled DNA used for transient transfections.

Species Used Chicken, TS34 a6 L1, [LSCC HD2], erythroblast

Before the Pulse

Cell Growth Medium Not given

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 (6) cells / pulse

Voltage 0.270 kV

Volume of Cells 0.5 ml

Field Strength 0.675 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 not used

After the Pulse

Outgrowth Medium Not given

Time Constant 24 msec

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
PBS: 1x = 8g NaCl, 0.2g KCl, 0.2g KH₂PO₄, 1.15g Na₂HPO₄

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

187

Gene Pulser® Electroprotocol

Cell Type	Other Cell Types	Molecules Electroporated	Analoge cytoskeletal proteins (like tubulin).
Species Used	Hydra cells, <i>Cnidaria</i>		

Before the Pulse

Cell Growth Medium	(Dissociation buffer) : 1.2 mM MgSO ₄ , 6 mM CaCl ₂ , 3.6 mM KCl, 6 mM pyruvate; 6 mM Na-Citrate, 12.5 mM TES buffer, 6 mM glucose, pH 6.9.
	Growth Phase at Harvest Not given
	Pre-pulse Incubation Not given

Wash Solution Not given

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature	Room temperature
Electroporation Medium	The same as cell growth medium
Cell Density	10 (5) cells / ml
Volume of Cells	10 (4) cells / pulse
DNA Concentration	Not given
DNA Resuspension Buffer	Not given
Volume of DNA	Not given
After the Pulse	
Outgrowth Medium	The same

Cuvette Gap 0.4 cm

Voltage 0.230 kV

Field Strength 0.575 kV/cm

Capacitor 25 µF

Resistor (Pulse Controller) none Ω

Time Constant 4.3 msec

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature	Room temperature
Length of Incubation	Not given
Selection Method or Assay Used	These cells are "primary cell cultures"
Electroporation Efficiency	Not given
Per Cent Survival	Not given

Name of Submitter Charo Gonzalez Agosti

Institution Address Zoologisches Inst.
The Univ. Zurich
Winterthurerstr. 190
8057 Zurich, Switzerland

Survey Number

189

Gene Pulser® Electroprotocol

Cell Type Other Cell Types

Molecules Electroporated DNA: supercoiled & linear; 3.3 - 33 kb pY, PR-neo.

Species Used *Leishmania*, all species within the genus

Before the Pulse

Cell Growth Medium M199 (see Kapler *et. al.* 1990)

Growth Phase at Harvest Late log

Pre-pulse Incubation up to 2 hours

Wash Solution 21mM HEPES, pH 7.5, 0.7mM Na₂PO₄, 137mM NaCl, 6 mM glucose, 5mM KCl

The Pulse

Instruments Used Not given

Electroporation Temperature Cells, cuvettes, DNA on ice; 0 °C

Electroporation Medium 21 mM HEPES, pH 7.5, 0.7 mM Na₂PO₄, 137 mM NaCl, 6 mM glucose, 5 mM KCl

Cuvette Gap 0.2 cm

Cell Density 10 (8) / ml

Voltage 0.45 kV

Volume of Cells 0.4 ml

Field Strength 2.25 kV/cm

DNA Concentration 300 to 1000 µg / ml

DNA Resuspension Buffer TE

Capacitor 500 µF

Volume of DNA 1 to 100 µg

Resistor (Pulse Controller) none Ω

After the Pulse

Time Constant ~ 4 msec

Outgrowth Medium M199 medium

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.
 Kapler, *et. al.*, 1990 *Molec. Biol.* **10**:1087 (G418).
 Cruz & Beverly, 1990 *Nature* **348**:171. Gene replacement.
 LeBowitz, *et. al.*, 1990 *PNAS* **87**: 9736. Expression vector.
 Coburn, *et. al.*, 1991 *Molec. Bioch. Parasitology* **46**:169 (diverse species).
 LeBowitz, *et. al.*, 1991 *Gene* **103**:119-123. b-gal, b-gluc reporters.
 Cruz, *et. al.*, 1991 *PNAS* **88**:7170-7174. Hygromycin & gene replacement.

Outgrowth Temperature 26 °C

Length of Incubation Overnight

Selection Method or Assay Used G418, hygromycin, gancyclovir β-galactosidase, β-glucuronidase

Electroporation Efficiency 10 to 60 transformants / µg DNA, up to 10⁽⁻⁴⁾ / cell

Per Cent Survival 50%

Name of Submitter Stephen M. Beverley

Institution Address Biological Chem. & Molec. Pharmacology Dept.
 Harvard Medical School
 250 Longwood Ave
 Boston, MA 02115

Survey Number

190

Gene Pulser® Electroprotocol

Cell Type Other Cell Types

Molecules Electroported DNA: plasmid constructs (containing VSG- gene, promoter *T. brucei*)

Species Used *Trypanosoma brucei brucei*, AnTat 1.3A, (blood- stream forms); EATRO 1125 (procyclic forms).

Before the Pulse

Cell Growth Medium Cunningham's medium + 15% Fetal Calf Serum
(See references in notes)

Growth Phase at Harvest Mid- log phase

Pre-pulse Incubation Not given

Wash Solution Not given

The Pulse

Instruments Used Gene Pulser® apparatus

Electroporation Temperature Ambient temperature

Electroporation Medium Zimmerman's post-fusion medium

Cuvette Gap 0.4 cm

Cell Density 2 x 10⁷ / ml

Volume of Cells 500 µl

Voltage 1.5 kV

DNA Concentration 1 mg / ml

Field Strength 3.75 kV/cm

DNA Resuspension Buffer TES buffer

Capacitor 25 µF

Volume of DNA 20 to 50 µl

Resistor (Pulse Controller) none Ω NOT

After the Pulse

Time Constant Not given

Outgrowth Medium Cunningham's medium/Baltes Medium (procyclics) / (Bloodstream forms)

Outgrowth Temperature 27 °C procyclics/ 37° C bloodstream forms

Length of Incubation 12 to 18 hours

Selection Method or Assay Used CAT assay

Electroporation Efficiency Not done

Per Cent Survival 30 to 60%

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

The conditions described are not those used in my own publication, Jefferies, *et. al.* 1991 *Mol. Cell. Biol.* **11**: 338-341, but were used by Clayton *et. al.*, 1990 *Mol. Cell. Biol.* **10**: 3036-3047, to transfect procyclic trypanosomes.

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

Name of Submitter Dr. David Jefferies

Institution Address Univ. Brussels
Molecular Biology
67, rue dar cheveaux
B1640 Rhode-Saint-- Genese
Brussels, BELGIUM

Survey Number

191



**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
