



Examples of Electro-transfected Eukaryotic and Prokaryotic Cells

Mammalian

[†]Hamster fibroblast (CHO)
[†]Human B
[†]Human, B lymphomas: BJAB,P3HR-1, B95-8, UC729-6
[†]Human, B-cell, JY, Epstein-Barr virus transformed
[†]Human, C-41, cervical carcinoma
[†]Human epithelial cells
[†]Human erythroleukemia (K-562)
[†]Human, epithelial, cervix (HeLa)
[†]Human fibroblast
[†]Human, GCT, fibrous histiocytoma
 Human glioblastoma
[†]Human, HEL cells, eythrolekemia
[†]Human, Hep3b2, hepatocytes
[†]Human, HepG2, hepatoma
[†]Human, HL 60, eythrolekemia
[†]Human hybridoma
[†]Human, JEG-3, choriocarcinoma
[†]Human, K562, chronic myeloid leukemia
[†]Human, kidney, 293
[†]Human, lymphocytes, primary
[†]Human lymphoma EB4, Ly67, HUT-78
[†]Human, MCF-7, breast
[†]Human, MRC-5, lung
 Human medullary thyroid (MTC)
 Human melanoma (SKMEL 25, SKMEL 28)
[†]Human, pancreatic
 Human primary monocytes
[†]Human, Raji, Burkitt lymphoma
[†]Human, red blood cells
[†]Human, skin carcinoma
[†]Human, squamous cell carcinoma, oral and cervical lines
[†]Human, T lymphoma (Jurkat)
[†]Human, T lymphoblastoid, CEMx174
[†]Human, TCCSUP (epithelial-like) bladder carcinoma
[†]Human, U373, glioblastoma
[†]Human, U937, hystiocytic lymphoma
[†]Human, V79, skin cells, fibroblasts

[†]Hybrid, mouse/human, A9 fibroblast
[†]Hybrid, rat/mouse, MEL cells

[†]Monkey, kidney (COS-1, COS-7, CV-1, Vero)

[†]Mouse, 3T3, embryo
[†]Mouse, 32d, myeloma
 Mouse, B
[†]Mouse, BALB/c 3T3, clone A31, fibroblast, embryo
[†]Mouse, BbSutA, hematopoietic
[†]Mouse, C127, fibroblast, mammary tumor
[†]Mouse, C2, muscle myoblast
[†]Mouse, C2C12, muscle
[†]Mouse, D10.G4.1, T-cell, helper

Mammalian (continued)

[†]Mouse, embryonic stem cells (ES-D3, E14)
[†]Mouse, erythrolekemia cells
[†]Mouse, FDC-PI, II-3-dependent cell line
 Mouse fibroblast
 Mouse, hepatoma (Hepa 1-6)
[†]Mouse, J558-L, myeloma.
[†]Mouse, L-cells
[†]Mouse, L cell derivative, A-9
[†]Mouse, L929, connective tissue
[†]Mouse, LM(TK-), connective tissue
[†]Mouse, mammary epithelial cells
[†]Mouse, myeloma, 32d, p3x63AG8 (X-63)
 Mouse macrophage (RAW 264)
 Mouse mammary tumor (C127)
[†]Mouse, NIH/3T3; embryo
[†]Mouse, NSO, myeloma cells
[†]Mouse, p3x63AG8; myeloma
[†]Mouse, SP2/O, [Sp-2], myeloma
 Mouse T
[†]Mouse, WEHI-3B, myelomonocytic leukemia
[†]Mouse, X-63, myeloma [p3 X63 - Ag8.653]
[†]Ovine (sheep), CSL503, fetal lung
[†]Ovine (sheep), R.E., rumen
[†]Rat-1
[†]Rat brain
[†]Rat, CA77, medullary thyroid carcinoma cell line
[†]Rat, D202CC, hepatoma
[†]Rat, fibroblasts
[†]Rat, H4-11-E-C3, hepatoma
[†]Rat, L-6, myoblast
 Rat lymphoma (BW5147)
 Rat macrophage
 Rat myeloma (4B 2/O)
[†]Rat, N62 T cells
[†]Rat, PC12, adrenal pheochromocytoma
[†]Rat, submandibular acini (secretory cells)

Plant Protoplants

Rice	Sugar beet	Lettuce
Maize	Pine	Potato
Alder tree	Sugar cane	Soybean
Tobacco	Carrot	Tomato

Intact Plant Cells

Carrot
Chenopodium rubrum
[†]*Hedyotis corymbosa*
[†]*Lactuca sativa* (aka chirimen chisha)
[†]Maize
[†]*Nicotiana plumbaginifolia*; protoplasts from leaf
[†]*Oryza sativa*, cv. Yamahouci or cv. Nihonbare
Thalictrum rugosum

Fungal Cells

- † *Aspergillus* (4 species)
- † *Candida* (2 species)
- † *Colletotrichum gloeosporioides* (a fungal phytopathogen)
- † *Cryptococcus neoformans*, ma5 mutants
- † *Dictyostelium discoideum*
- Fusarium solani*
- Leptosphaeria maculans*
- Neurospora crassa*
- † *Pichia pastoris*
- Penicillium urticae*
- † *Saccharomyces cerevisiae* (6 strains)
- † *Schizosaccharomyces pombe*
- Trichoderma harzianum*

Bacteria, gram negative

- † *Acetobacter xylinum*, ATCC 23769
- Actinobacillus* (2 species)
- † *Agrobacterium* (3 species)
- Azospirillum brasilense*
- † *Bacteroides* (3 species)
- Bradyrhizobium* (2 species)
- Brucella abortus*
- Butyrivibrio fibrisolvens*
- Campylobacter* (2 species)
- Caulobacter* (3 species)
- Citrobacter freundii*
- † Cyanobacteria, primarily filamentous, *Anabaena* species
- Enterobacter aerogenes*
- † *Escherichia coli* (62 strains)
- Erwinia* (2 species)
- Francisella* (2 species)
- Haemophilus influenzae*
- Klebsiella* (3 species)
- † *Legionella* (3 species)
- Pasteurella* "aemolytica"
- Proteus* (2 species)
- Pseudomonas* (8 species)
- Rhizobium* (unspciated)
- Rhodopseudomonas viridis*
- Rhodospirillum molischanium*
- Rochalimaea quintana*
- † *Salmonella* (4 species)
- Serratia* (2 species)
- † *Vibrio* (2 species)
- Xanthomonas campestris*
- Yersinia* (3 species)

Bacteria, gram positive

- Amycolatopsis mediterranea*
- Amycolatopsis orientalis*
- † *Bacillus* (7 species)
- † *Brevibacterium* (3 species)
- Clavibacter* (4 species)
- Clostridium* (2 species)
- † *Corynebacterium* (3 species)
- Cytophaga johnsonae*
- † *Enterococcus* (4 species)
- Fremyella diplosiphon*
- † *Lactobacillus* (21 species)

Bacteria, gram positive (continued)

- Leuconostoc* (3 species)
- Listeria* (5 species)
- † *Mycobacterium* (3 species)
- Pediococcus acidilactici*
- Propionibacterium jensenii*
- Rhodococcus fascians*
- † *Staphylococcus* (7 species)
- † *Streptococcus* (11 species)

Cyanobacteria

- Anabaena*
- Fremyella diplosiphon*
- Nostoc*
- Synechococcus*

Other Bacteria

- Acholeplasma laidlawii*
- Methylophilus* (3 species)
- Mycobacterium* (4 species)
- Serpula hydrysenteriae*
- Spiroplasma citri*
- Treponema hydrysenteriae*

Other Cell Types

- Anaerobic bacteria
- † Chicken, HD11, macrophage
- † Chicken, primary hepatocytes
- † Chicken, TS34 a6 L1, [LSCC HD2], erythroblast
- Chloroplast, spinach
- Encapsulated bacteria
- Frog, retina
- Giardia lamblia*
- † Hydra cells, *Cnidaria*.
- † *Leishmania*, all species within the genus
- Marine organisms
- Photosynthetic bacteria
- Primary chicken oviduct
- Sea urchin egg
- Tetrahymena*
- Thermotolerant bacteria
- † *Trypanosoma brucei* (blood stream and procyclic forms)

Other Applications

- Cotransformation
- Cosmids
- Carbohydrates
- Direct transfer (donor to recipient)
- Fluorescent molecules
- Influence of DNA size on efficiency
- Influence of DNA conformation on efficiency
- Large DNA fragments
- Ligation mixtures
- Library construction
- M13 DNA
- Plasmid incompatibility
- RNA
- Release of cell components
- Use of more than one electroperoration unit

† Experimental conditions available as an Electroprotocol



**Bio-Rad
Laboratories**

Life Science
Group

Website: www.bio-rad.com **Bio-Rad Laboratories** **Minis Office** 2000 Alfred Nobel Drive, Hercules, CA 94547, Ph. (510) 767-7000, Fx. (510) 747-5800
Also in: **Alitalia** Ph. (02) 99 74-2000, Fx. (02) 99 74-2000 **Austria** Ph. (7) 877 69 07, Fx. (7) 876 66 29 **Belgium** Ph. (0) 308 66 77, Fx. (0) 308 66 54
Canada Ph. (605) 772-2777, Fx. (605) 772-2900 **China** Ph. (86-70) 2046622, Fx. (86-70) 2057876 **Denmark** Ph. (39) 77 9947, Fx. (39) 27 7698
Finland Ph. (9) 804 2300, Fx. (9) 804 7700 **France** Ph. (7) 43 90 46 90, Fx. (7) 46 77 24 67 **Germany** Ph. 089 37864-0, Fx. 089 37864-700
Hong Kong Ph. 7203300 Fx. 720 257 **India** Ph. 91-7-7467-0703, Fx. 91-77-467-0705 **Israel** Ph. 03 957 4 727, Fx. 03 957 4 729
Italy Ph. 02-27809-7, Fx. 02-27809-309 **Japan** Ph. 03-5877-6270, Fx. 03-5877-6272 **The Netherlands** Ph. 0373 76-540066, Fx. 0373 76-542276
New Zealand Ph. 09-443 3000, Fx. 09-443 3007 **Singapore** Ph. (65) 272-9877, Fx. (65) 273-4238 **Spain** Ph. (91) 667 70 85, Fx. (91) 667 9698
Sweden Ph. 46 (0) 8 627 80 00, Fx. 46 (0) 8 627 54 00 **Switzerland** Ph. 07-800 88 58, Fx. 07-800 99 00
United Kingdom Ph. 0800 707 734, Fx. 01442 259778