

# Electroprotocols Species List

<b>Bacterial Cells, Gram-Negative</b>	<b>Survey Number(s)</b>	<b>Bacterial Cells, Gram-Negative</b>	<b>Survey Number(s)</b>
<i>Acetobacter xylinum</i> , ATCC 23769 .....	1	<i>Legionella pneumophila</i> , strain Nottingham N -7.....	50
<i>Agrobacterium rhizogenes</i> .....	3	<i>Legionella pneumophila (philadelphia)</i> .....	51, 203
<i>Agrobacterium tumefaciens</i> .....	3, 198	<i>Pseudomonas putida</i> ATCC 12633 .....	52
<i>Agrobacterium</i> , unspecified species .....	2	<i>Pseudomonas syringae</i> .....	53
<i>Bacteroides fragilis</i> .....	47	<i>Salmonella senftenburg</i> .....	200
<i>Bradyrhizobium japonicum</i> .....	48	<i>Salmonella typhimurium</i> .....	54, 200
Cyanobacteria, primarily filamentous, <i>Anabaena</i> species .....	4	<i>Vibrio anguillarum</i> .....	55
<i>E. coli</i> , BB4.....	5	<i>Xanthomonas campestris</i> .....	53
<i>E. coli</i> , CSR 603 maxi cells.....	6	<b>Bacterial Cells, Gram-Positive</b>	<b>Survey Number(s)</b>
<i>E. coli</i> , DH5 $\alpha$ .....	7 to 25, 35, 38, 41, 49, 198, 200, 201	<i>Bacillus sphaericus</i> 1593 .....	56
<i>E. coli</i> , DH10B.....	26, 27	<i>Brevibacterium flavum</i> , ATCC21475.....	58, 59
<i>E. coli</i> , HB101.....	10, 13, 28 to 31, 200	<i>Brevibacterium lactofermentum</i> .....	57
<i>E. coli</i> , JM83.....	12, 31	<i>Corynebacterium glutamicum</i> ATCC 13032.....	59
<i>E. coli</i> , JM105.....	28	<i>Corynebacterium</i> , unspecified strain.....	58
<i>E. coli</i> , JM109 .....	32, 35, 40	<i>Enterococcus faecalis</i> JH2-2 and UV202 .....	60
<i>E. coli</i> , K12.....	33	<i>Enterococcus hirae</i> .....	61
<i>E. coli</i> , LE 392.....	49	<i>Lactobacillus acidophilus</i> .....	62, 63
<i>E. coli</i> , MC1061 .....	34 to 39, 46	<i>Lactobacillus delbrueckii</i> .....	64
<i>E. coli</i> , MC4100 .....	201	<i>Lactobacillus fermentum</i> .....	65, 66
<i>E. coli</i> , MV1184 .....	40	<i>Lactobacillus gasseri</i> .....	67
<i>E. coli</i> , MV1190 .....	35	<i>Lactobacillus plantarum</i> .....	69
<i>E. coli</i> , N99.....	41	<i>Lactobacillus reuteri</i> .....	68, 70, 71
<i>E. coli</i> , NM522.....	35, 37, 42	<i>Lactobacillus salivarius</i> .....	72
<i>E. coli</i> , NR9162.....	46	<i>Lactobacillus sp.</i> , strain 100-33.....	73
<i>E. coli</i> , S9OC.....	202	<i>Lactobacillus sp.</i> , strain ES1 .....	74
<i>E. coli</i> , TG1 .....	5, 43, 199	<i>Lactococcus lactis</i> , subspecies <i>cremoris</i> .....	75, 76
<i>E. coli</i> , W3110 .....	8	<i>Lactococcus lactis</i> , subspecies <i>lactis</i> .....	75, 76
<i>E. coli</i> , XA90 .....	199	<i>Mycobacterium bovis</i> , BCG.....	78
<i>E. coli</i> , XL-1 Blue .....	35, 38, 199	<i>Mycobacterium smegmatis</i> .....	79, 80
<i>E. coli</i> , unspecified strain .....	44, 45, 47, 48, 50, 52	<i>Mycobacterium</i> , unspecified species .....	77
<i>Legionella longbeachae</i> .....	51	<i>Staphylococcus aureus</i> .....	81, 82, 83, 204
		<i>Streptococcus sanguis</i> .....	84

<b>Mammalian Cells</b>	<b>Survey Number(s)</b>
B-cell line, unspecified .....	121
Hamster, CHO, ovary.....	85 to 91, 156, 168
Human, 293, kidney .....	108, 156
Human, B lymphomas: BJAB, P3HR-1, B95-8.....	157
Human, C-4I, cervical carcinoma.....	155
Human, CEMx174, T lymphoblastoid .....	92
Human epithelial cells.....	158
Human, fibroblast .....	160, 163
Human, GCT, fibrous histiocytoma.....	93
Human, HEL cells, eythroleukemia.....	158, 166, 192
Human, HeLa, epithelial carcinoma.....	94 to 98, 158, 167, 192, 196
Human, Hep3b2, hepatocytes.....	98, 163
Human, HepG2, hepatoma .....	99, 161
Human, HL 60, eythroleukemia.....	166
Human, HuT78, cutaneous T-cell lymphoma .....	100
Human, JEG-3, choriocarcinoma .....	101
Human, Jurkat .....	107
Human, JY, B-cell, Epstein-Barr virus transformed .....	102
Human, K562, chronic myeloid leukemia.....	103 to 106, 166, 168, 192
Human, lymphoblast cell lines, EBV immortalized .....	108
Human, lymphocytes, primary.....	194
Human, MCF-7, breast.....	110
Human, MRC-5, lung.....	111, 112
Human, pancreatic .....	113
Human, Raji, Burkitt lymphoma .....	107
Human, red blood cells .....	114
Human, squamous cell carcinoma, oral & cervical lines.....	162
Human, T-cells .....	107, 159
Human, TCCSUP (epithelial-like) bladder carcinoma.....	153
Human, U373, glioblastoma .....	115
Human, U937, hystiocytic lymphoma .....	106, 116, 117
Human, UC729-6, lymphoblastoid, B-cells.....	118
Human, V79, skin cells, fibroblasts .....	119
Hybrid, rat/mouse, MEL cells.....	168
Hybrid, mouse/human , A9 fibroblast .....	165
Hybridomas .....	164

<b>Mammalian Cells</b>	<b>Survey Number(s)</b>
Monkey, COS, kidney .....	126 to 128, 154, 161, 195
Monkey, CV-1, kidney .....	167
Monkey, Vero, kidney.....	144, 155, 158, 160, 162
Mouse, 3T3, embryo.....	160, 167
Mouse, 32d, myeloma.....	134
Mouse, A-9, derivative of mouse L cell.....	120
Mouse, BALB/c 3T3, clone A31, fibroblast, embryo .....	122
Mouse, BbSutA, hematopoietic.....	166
Mouse, C127, fibroblast, mammary tumor .....	125, 193
Mouse, C2, muscle myoblast.....	123
Mouse, C2C12, muscle .....	124
Mouse, D10.G4.1, T-cell, helper.....	143
Mouse, embryonic stem cells .....	129, 130
Mouse, erythroleukemia cells.....	131, 132
Mouse, FDC-PI, Il-3-dependent cell line .....	133
Mouse, J558-L, myeloma .....	134, 135
Mouse, L-cells .....	163
Mouse, L929, connective tissue.....	109
Mouse, LM(TK-), connective tissue.....	136, 137
Mouse, mammary epithelial cells .....	154
Mouse, NIH/3T3; embryo.....	139, 159
Mouse, NSO, myeloma cells.....	140
Mouse, p3x63AG8; myeloma.....	167
Mouse, SP2/0, [Sp-2], myeloma .....	141, 142
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 .....	195
Rat, submandibular acini (secretory cells) .....	147

