

Place Your JCRB Cell Line Order

| JCRB No. | Cell Name | ANIMAL | TISSUE | GENETICS | CHARACT | COMMENT | celltype | Ordering Comment |
|-----------|------------------|--------|--------|---|---------|---------|----------------------------------|------------------|
| AyuK10A10 | Cdc42se1-K1 | mouse | embryo | Cdc42se1, CDC42 pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK10C05 | Map2k1-K1 | mouse | embryo | Map2k1, mitogen a pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK10C08 | Gata2a-K1 | mouse | embryo | Gata2a, GATA zir pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK10D02 | Ubr5-K1 | mouse | embryo | Ubr5, ubiquitin proI pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK10D10 | Ppp2r2d-K1 | mouse | embryo | Ppp2r2d, protein pi pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK10H01 | Mapk1ip1-K1 | mouse | embryo | Mapk1ip1, mitiger pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK10H04 | Cbx3-K1 | mouse | embryo | Cbx3, chromobox I pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK11B06 | Atf1-K1 | mouse | embryo | Atf1, activating tr pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK11C02 | Ahsa1-K1 | mouse | embryo | Ahsa1, AHA1, acti pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK11D09 | Zfp568-K1 | mouse | embryo | Zfp568, zinc finger pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK11E04 | Csnk2a1-K1 | mouse | embryo | Csnk2a1, casein ki pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK11E11 | Wdr4-K1 | mouse | embryo | Wdr4, WD repeat pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12A07 | Saps3-K1 | mouse | embryo | Saps3, SAPS dom pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12A12 | Mipol1-K1 | mouse | embryo | Mipol1, mirror-imaq pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12C03 | Fryl-K1 | mouse | embryo | Fryl, furry homolog pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12C11 | Ctdspl2-K1 | mouse | embryo | Ctdspl2, CTD (cart pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12C12 | Igsf9b-K1 | mouse | embryo | Igsf9b, immunoglobl pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12D06 | Ccdc58-K1 | mouse | embryo | Ccdc58, coiled-coil pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12E07 | Trip10-K1 | mouse | embryo | Trip10, thyroid horr pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12E09 | Rad51ap1-K1 | mouse | embryo | Rad51ap1, RAD51 pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12E12 | Thrap3-K1 | mouse | embryo | Thrap3, thyroid hor pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12F03 | Anks3-K1 | mouse | embryo | Anks3, ankyrin rep pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK12H05 | Atad1-K1 | mouse | embryo | Atad1, ATPase fan pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13A11 | Rap1b-K1 | mouse | embryo | Rap1b, RAS relate pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13C12 | Jrk-K1 | mouse | embryo | Jrk, jerky pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13E11 | Cbx5-K1 | mouse | embryo | Cbx5, chromobox I pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13F02 | Tarbp2-K1 | mouse | embryo | Tarbp2, TAR (HIV) pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13F05 | Phf20-K1 | mouse | embryo | Phf20, PHD finger pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13G03 | Zfp206-K1 | mouse | embryo | Zfp206, zinc finger pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK13G06 | BC024479-K1 | mouse | embryo | BC024479, cDNA : pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK14A05 | Trerf1-K1 | mouse | embryo | Trerf1, transcription pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK14C04 | Fubp3-K1 | mouse | embryo | Fubp3, far upstream pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK14E01 | Kntc1-K1 | mouse | embryo | Kntc1, kinetochore pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK14E08 | Pspc1-K1 | mouse | embryo | Pspc1, paraspickei pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK14E11 | Milt6-K1 | mouse | embryo | Milt6, myeloid/lymph pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK14G01 | Cdca7-K1 | mouse | embryo | Cdca7, cell divisor pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK14G12 | Nr5a2-K1 | mouse | embryo | Nr5a2, nuclear rec pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK15A08 | Anp32e-K1 | mouse | embryo | Anp32e, acidic (leu pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK15E07 | Leo1-K1 | mouse | embryo | Leo1, Leo1, Paf1/F pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK15F02 | Deaf1-K1 | mouse | embryo | Deaf1, deformed e pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK15G02 | Kpna4-K1 | mouse | embryo | Kpna4, karyopherin pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK15G06 | Capza2-K1 | mouse | embryo | Capza2, capping p pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK15H03 | Psmc3ip-K1 | mouse | embryo | Psmc3ip, proteaseo pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16A10 | Fmn1l2-K1 | mouse | embryo | Fmn1l2, formin-like pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16C06 | Dgr8r-K1 | mouse | embryo | Dgr8r, DiGeorge's pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16C10 | Gnb2-K1 | mouse | embryo | Gnb2, guanine nuc pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16C12 | Trim71-K1 | mouse | embryo | Trim71, tripartite m pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16E07 | Lrch4-K1 | mouse | embryo | Lrch4, leucine-rich pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16E08 | Rbpms-K1 | mouse | embryo | Rbpms, RNA bindi pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK16H07 | Nmnat2-K1 | mouse | embryo | Nmnat2, nicotinan pCMT-SAHygPA-NP22 | | | mouse homozygous mutant ES cells | |
| AyuK17A06 | Trim6-K1 | mouse | embryo | Trim6, triparalite mo pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17A12 | Lnx2-K1 | mouse | embryo | Lnx2, ligand of nun pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17B02 | Tada2l-K1 | mouse | embryo | Tada2l, transcriptic pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17C03 | Hspbp1-K1 | mouse | embryo | Hspbp1, Hspb as pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17E05 | Ptnp11-K1 | mouse | embryo | Ptnp11, protein tyr pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17E10 | Myh9-K1 | mouse | embryo | Myh9, myosin, hea pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17G03 | Gtf2h1-K1 | mouse | embryo | Gtf2h1, general tra pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK17H08 | Pcbp2-K1 | mouse | embryo | Pcbp2, poly(rC) bir pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18B07 | Zfp518-K1 | mouse | embryo | Zfp518, zinc finger pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18B08 | Setd5-K1 | mouse | embryo | Setd5, SET domai pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18B12 | Epn2-K1 | mouse | embryo | Epn2, epsin 2 pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18C02 | 1110007L15Rik-K1 | mouse | embryo | 1110007L15Rik, R pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18D01 | Birc6-K1 | mouse | embryo | Birc6, baculoviral I pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18E01 | Iblk-K1 | mouse | embryo | Iblk, inhibitor of Br pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK18E04 | Thumpd3-K1 | mouse | embryo | Thumpd3, THUMP pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK19B07 | 593041619Rik-K1 | mouse | embryo | 593041619Rik, RII pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK19D05 | Fam49b-K1 | mouse | embryo | Fam49b, family wit pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK19F03 | Nmt2-K1 | mouse | embryo | Nmt2, N-myristoylti pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK19F08 | Tex19-K1 | mouse | embryo | Tex19, testis expre pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK19G11 | BC057627-K1 | mouse | embryo | BC057627, cDNA : pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK20B04 | Tjp2-K1 | mouse | embryo | Tjp2, tight junction pT2F2-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK20B09 | Srk1-K1 | mouse | embryo | Srk1, serine/argin pT2F2-SAHygPA-NP21-Rev | | | mouse homozygous mutant ES cells | |
| AyuK5F06 | Eapp-K1 | mouse | embryo | Eapp, E2F-associa pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK6A05 | Ranbp1-K1 | mouse | embryo | Ranbp1, RAN bind pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK6B03 | G3bp2-K1 | mouse | embryo | G3bp2, GTPase at pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |
| AyuK6B09 | Tbrg4-K1 | mouse | embryo | Tbrg4, transformin pCMT-SAHygPA-NP21 | | | mouse homozygous mutant ES cells | |

| | | | | | |
|----------|------------------|----------------------|-----------------------------|---|---------------------------------|
| AyuK6C01 | Jagn1-K1 | mouse | embryo | Jagn1, jagnal hon pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6C07 | 5830457O10Rik-K1 | mouse | embryo | 5830457O10Rik, R pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6D12 | Smg5-K1 | mouse | embryo | Smg5, Smg-5 hom pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6E04 | Armet-K1 | mouse | embryo | Armet, arginine-rich pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6F01 | Gpatch4-K1 | mouse | embryo | Gpatch4, G patch :pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6F07 | Bat2d-K1 | mouse | embryo | Bat2d, BAT2 doma pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6F11 | Rnf44-K1 | mouse | embryo | Rnf44, ring finger p CMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6G02 | Eif4b-K1 | mouse | embryo | Eif4b, eukaryotic tr pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK6H07 | Cbx1-K1 | mouse | embryo | Cbx1, chromobox 1 pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7A06 | Epb4.1I5-K1 | mouse | embryo | Epb4.1I5, erythroc pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7A07 | Cxxc5-K1 | mouse | embryo | Cxxc5, CXXC finger pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7B02 | Avpi1-K1 | mouse | embryo | Avpi1, arginase vas pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7C09 | Axin1-K1 | mouse | embryo | Axin1, axin 1 pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7D01 | Cfdp1-K1 | mouse | embryo | Cfdp1, craniofacial pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7G08 | Rbm5-K1 | mouse | embryo | Rbm5, RNA bindin pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7G10 | Zfp280b-K1 | mouse | embryo | Zfp280b, zinc finger pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7G11 | Rsrc2-K1 | mouse | embryo | Rsrc2, arginine/ser pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK7H06 | Cnbp-K1 | mouse | embryo | Cnbp, cellular nucl pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8A06 | Nmt1-K1 | mouse | embryo | Nmt1, N-myristyli pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8B01 | Ilf2-K1 | mouse | embryo | Ilf2, interleukin enh pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8C06 | Cnot10-K1 | mouse | embryo | Cnot10, CCR4-NO pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8C12 | Gm561-K1 | mouse | embryo | Gm561, gene mod pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8D01 | Eno1-K1 | mouse | embryo | Eno1, enolase 1, a pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8D05 | Ccdc77-K1 | mouse | embryo | Ccdc77, coiled-coil pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8E11 | Ube2c-K1 | mouse | embryo | Ube2c, ubiquitin-cc pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8F07 | Slc39a14-K1 | mouse | embryo | Slc39a14, solute c: pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8G02 | Nosip-K1 | mouse | embryo | Nosip, nitric oxide :pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK8G11 | Nedd4-K1 | mouse | embryo | Nedd4, neural prec pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK9B05 | Zfp532-K1 | mouse | embryo | Zfp532, zinc finger pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK9E04 | AU019823-K1 | mouse | embryo | AU019823, expres pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK9F02 | Pml-K1 | mouse | embryo | Pml, promyelocytic pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| AyuK9F06 | Ddx11-K1 | mouse | embryo | Ddx11, DEAD/H (A pCMT-SAhygpA-NP21 | mousehomozygous mutant ES cells |
| IFO50004 | WISH | human | HeLa-contaminant | Originally establish *Contamination of I general cells | |
| IFO50005 | J-111 | human | HeLa-cont | HeLa markers obs originally establish *Contamination of I general cells | |
| IFO50006 | M1 | mouse | hemo-lymphocytic | spontaneous myeloid leukemia, differe general cells | |
| IFO50007 | Ca Ski | human | uterine cervix | cervical epidermoic This cell line contai general cells | |
| IFO50008 | Mm1 | mouse | hemo-lym subline of M-1 mou | spontaneous myeloid leukemia, M-1-di general cells | |
| IFO50009 | G-361 | human | melanoma | malignant melanoma | general cells |
| IFO50010 | C6/36 | mosquito | hatched larvae | useful for replication of flavivirus | general cells |
| IFO50011 | HeLa S3 | human | uterine cervix | cervical epithelioid Contains DNA seq | general cells |
| IFO50012 | Pt K2 | mouse | marsupial | kidney, epithelial-like | general cells |
| IFO50013 | RPMI 8226 | human | hemo-lymphocytic | myeloma, multiple | general cells |
| IFO50015 | F2408 | rat | embryo | fibroblast-like | general cells |
| IFO50016 | Chang Liver | human | HeLa-contaminant | originally establish *Contamination of I general cells | |
| IFO50017 | lXc | rat | sarcoma | useful for assay of This cell line contai general cells | |
| IFO50018 | NIL | hamster | \$ embryo | transformed from hamster fibroblast | general cells |
| IFO50019 | NIH 3T3 p-7 cl-3 | mouse | embryo | fibroblast, contact-inhibited | general cells |
| IFO50020 | SIRC | rabbit | cornea | susceptible to rubella virus | general cells |
| IFO50021 | A31-714 C4 | mouse | embryo | fibroblast, subclone of Balb/3T3 A31. 1 | general cells |
| IFO50022 | HL-60 | human | hemo-lymphocytic | leukemia, acute prc This line is equiv | general cells |
| IFO50025 | IM-9 | human | hemo-lymphocytic | B-lymphoblastoid cells, EB virus-transf | general cells |
| IFO50026 | CCRF-SB | human | hemo-lymphocytic | leukemia, acute lymphoblastic | general cells |
| IFO50034 | RBL-1 | rat | hemo-lymphocytic | basophilic leukemia | general cells |
| IFO50036 | P388D1 | mouse | hemo-lymphocytic | lymphoid neoplasm, derived from P38 | general cells |
| IFO50037 | RPMI 1788 | human | hemo-lymphocytic | established from normal blood, IgM se | general cells |
| IFO50038 | U-937 | human | hemo-lymphocytic | lymphoma, histiocytic | general cells |
| IFO50039 | NC-37 | human | hemo-lymphocytic | lymphoblast, Conta Contamination of R | general cells |
| IFO50040 | Namalwa | human | hemo-lymphocytic | lymphoma, Burkitt's | general cells |
| IFO50041 | MBT2 | mouse | bladder | bladder carcinoma | general cells |
| IFO50042 | C3H/MCA clone 15 | mouse | embryo | 3-methylcholanthrene-transformed C31 | general cells |
| IFO50043 | WiDr | human | colon | adenocarcinoma WiDr cell line is su | general cells |
| IFO50046 | Raji | human | hemo-lymphocytic | lymphoma, Burkitt's | general cells |
| IFO50067 | LoVo | human | colon | adenocarcinoma, carcinoembryonic ar | general cells |
| IFO50069 | Alexander cells | human | liver, gallbladder | hepatoma, HBs antigen-positive | general cells |
| IFO50070 | Balb/3T3-A31-1-1 | mouse | embryo | fibroblast, subclone of Balb/3T3-A31. 1 | general cells |
| IFO50071 | MDCK (NBL-2) | dog | kidney | established from normal kidney, susce | general cells |
| IFO50072 | MRC-9 | human | lung | fetus lung normal diploid fibroblast | general cells |
| IFO50073 | MRC-5 | human | lung | fetus lung normal diploid fibroblast | general cells |
| IFO50074 | HFL1 | human | lung | fetus lung normal diploid fibroblast | general cells |
| IFO50075 | WI-38 | human | lung | fetus lung normal diploid fibroblast | general cells |
| IFO50076 | IT-45R1 | rat | thymus | T cell differentiation factor-producing | general cells |
| IFO50077 | IT-45R91 | rat | thymus | thymic fibroblast | general cells |
| IFO50078 | IT-26R21 | rat | thymus | T cell differentiation factor producing | general cells |
| IFO50079 | Flow7000 | human | skin | skin normal diploid fibroblast | general cells |
| IFO50080 | NCTC clone 1469 | mouse | liver | normal liver-derived | general cells |
| IFO50081 | Neuro-2a | mouse | neural | neuroblastoma, Td = 20 hrs. | general cells |
| IFO50082 | V79 379A | hamster | C lung | lung, fibroblast, widely used in somatic | general cells |
| IFO50089 | Flow2000 | human | lung | fetus lung normal diploid fibroblast | general cells |
| IFO50090 | 4G12 hybridoma | mouse x h antibody-p | human lymphoid | 4G12 antibody, preferentially reacting | general cells |
| IFO50098 | F2408-No.3 | rat | embryo | HGPRT-deficient F2408 cell line | general cells |

| | | | | | |
|----------|---------------------|-------------------|-----------------------------|---|--|
| IFO50099 | F2408-B812 | rat | embryo | HGPRT-deficient. | HGPRT-deficient cell line from F2408. general cells |
| IFO50100 | F2408-B993 | rat | embryo | HGPRT(-), temper: | HGPRT-deficient cell line from F2408. general cells |
| IFO50101 | F2408-No.20 | rat | embryo | thymidine kinase-d | thymidine kinase-deficient F2408 cell li general cells |
| IFO50102 | F2408-No.7 | rat | embryo | HGPRT-deficient | HGPRT-deficient F2408 cell line general cells |
| IFO50103 | TuD-1 | rat | thymus | thymic stroma, keratin-positive | general cells |
| IFO50105 | GH3 | rat | pituitary | pituitary tumor, somatotrophin-secretin | general cells |
| IFO50106 | HOS | human | sarcoma | osteosarcoma | general cells |
| IFO50107 | G-292 clone A141B1 | human | sarcoma | osteosarcoma | general cells |
| IFO50108 | MG-63 | human | sarcoma | osteosarcoma | general cells |
| IFO50110 | C6 | rat | neural | glial tumor, GFAP-positive | general cells |
| IFO50111 | GRSL | mouse | hemo-lymphocytic | thymic leukemia This line was curec | general cells |
| IFO50112 | TaD-1 | rat | thymus | thymoma, keratin-positive | general cells |
| IFO50120 | TaD-1-3 | rat | thymus | thymoma, keratin-positive | general cells |
| IFO50151 | MEG-01 | human | hemo-lymphocytic | leukemia, chronic megakaryoblastic | general cells |
| IFO50152 | TuD-1-3 | rat | thymus | thymic stroma, keratin-positive | general cells |
| IFO50153 | A549 | human | lung (cancer) | adenocarcinoma | general cells |
| IFO50161 | KT-5 | mouse | neural | astrocyte cell line | general cells |
| IFO50210 | L2C | guinea pig | hemo-lymphocytic | leukemic cell, maintainable only by trar | general cells |
| IFO50221 | N18-RE-105 | mouse x r: neural | | N18TG2 x Fischer rat 18 day-embryon | general cells |
| IFO50248 | IT-79MTNC3 | mouse | thymus | thymic nurse cell | general cells |
| IFO50267 | PB-3c | mouse | hemo-lymphocytic | mast cell line | general cells |
| IFO50268 | EHS | mouse | sarcoma | chondrosarcoma, secreting basement | general cells |
| IFO50269 | GH1 | rat | pituitary | pituitary tumor, somatotrophin-secretin | general cells |
| IFO50270 | 6-23 clone 6 | rat | thyroid | medullary thyroid carcinoma, neuroten | general cells |
| IFO50271 | HUV-EC-C | human | vascular system | normal endothelial cell, umbilical cord | general cells |
| IFO50276 | A2058 | human | melanoma | malignant melanoma, metastasis to ly | general cells |
| IFO50277 | 5RP7 | rat | embryo | fibroblast transformed by c-Ha-ras onc | general cells |
| IFO50278 | PC-12 | rat | neural | pheochromocytoma, neuronal different | general cells |
| IFO50279 | NKY113 | mouse | antibody-producing hybridom | MSW113 antibody, recognizing a canc | general cells |
| IFO50280 | 77N1 | rat | kidney | transformant of NRK cells by avian sar | general cells |
| IFO50282 | RAT-2 | rat | embryo | thymidine kinase deficient RAT-1 (F24 | general cells |
| IFO50283 | IMR-32 | human | neural | N-myc amplification transplantable to nude mice | general cells |
| IFO50284 | A2B5 clone 105 | mouse | antibody-producing hybridom | A2B5 antibody, recognizing a glycolip | general cells |
| IFO50285 | U-251 MG (KO) | human | neural | astrocytoma, GFAT The DNA fingerpr | general cells |
| IFO50286 | SF126 | human | neural | astrocytoma, GFAP-negative | general cells |
| IFO50288 | U-251 MG | human | neural | GFAP-posotive, gli The DNA fingerpr | general cells |
| IFO50289 | Becker | human | neural | astrocytoma, GFAP-negative | general cells |
| IFO50293 | MA-89 | mouse | neural | normal mixed glial culture | general cells |
| IFO50294 | MEG-01SSF | human | hemo-lymj serum-free adapte | leukemia, chronic megakaryoblastic, IV | general cells |
| IFO50295 | NB-1 | human | neural | neuroblastoma, neuronal differentiatio | general cells |
| IFO50296 | WEHI-3b | mouse | hemo-lymphocytic | myelomonocyte | general cells |
| IFO50297 | HE-1 | human | embryo | Normal diploid fibrc whole fetus, normal diploid fibroblast | general cells |
| IFO50298 | Balb/c 3T3 A31-I-1 | mouse | embryo | fibroblast, subclone of Balb/3T3 A31. 1 | general cells |
| IFO50299 | Balb/c 3T3 A31-1-13 | mouse | embryo | fibroblast, subclone of Balb/3T3 A31. 1 | general cells |
| IFO50303 | T98G | human | neural | glioblastoma | general cells |
| IFO50308 | SKG-I | human | uterine cervix | cervical epidermoi This cell line contai | general cells |
| IFO50309 | SKG-II | human | uterine cervix | cervical squamous This cell line contai | general cells |
| IFO50310 | SKG-IIIa | human | uterine cervix | cervical epidermoi This cell line and S | general cells |
| IFO50311 | SKG-IIIb | human | uterine cervix | cervical epidermoi This cell line and S | general cells |
| IFO50312 | SNG-II | human | uterus | endometrial adeno This cell line does i | general cells |
| IFO50313 | SNG-M | human | uterus | endometrial adeno This cell line does i | general cells |
| IFO50314 | SKN | human | uterus | leiomyosarcoma This cell line does i | general cells |
| IFO50315 | RMG-I | human | ovary | clear cell carcinom This cell line does i | general cells |
| IFO50317 | RKN | human | ovary | leiomyosarcoma | general cells |
| IFO50320 | RMUG-S | human | ovary | mucinous cystader This cell line and R | general cells |
| IFO50321 | GAK | human | melanoma | malignant melanoma, metastasis to lyr | general cells |
| IFO50322 | NJG | human | uterus | gestational uterine This cell line does i | general cells |
| IFO50323 | BOKU | human | uterine cervix | cervical squamous This cell line contai | general cells |
| IFO50350 | Hs68 | human | skin | The enzyme aspar skin normal diploid fibroblast | general cells |
| IFO50354 | HT-1080 | human | sarcoma | acetabulum, fibrosarcoma | general cells |
| IFO50355 | ONS-76 | human | neural | medulloblastoma | general cells |
| IFO50356 | KNS-42 | human | neural | glioma, GFAP-positive, S-100 and NSI | general cells |
| IFO50357 | KNS-60 | human | neural | glioma, GFAP, S-100 and NSE-negati | general cells |
| IFO50358 | KNS-62 | human | lung (cancer) | bronchial squamous carcinoma, metas | general cells |
| IFO50359 | KNS-81 | human | neural | glioma, GFAP and S-100-positive, NSI | general cells |
| IFO50361 | B2-17 | human | neural | thymidine kinase-d astrocytoma, thymi | The DNA fingerpr |
| IFO50362 | MOLT-4 | human | hemo-lymphocytic | leukemia, acute lymphoblastic | general cells |
| IFO50363 | KU812 | human | hemo-lymphocytic | leukemia, chronic myeloid | general cells |
| IFO50364 | L6 | rat | muscular system | skeletal muscle myoblast | general cells |
| IFO50367 | SHOK | hamster, ♀ | embryo | fibroblast, contact-inhibited | general cells |
| IFO50368 | no.10 | human | neural | glioma, anaplastic, GFAP-positive | general cells |
| IFO50369 | no.11 | human | neural | glioma, anaplastic, GFAP-positive | general cells |
| IFO50373 | TKC2 | mouse | kidney | TG mouse harborir tubule cell, tsSV40 large T-immortalize | general cells |
| IFO50374 | TKD2 | mouse | vascular s | TG mouse harborir endothelial cell, tsSV40 large T-immort | general cells |
| IFO50375 | MSS31 | mouse | spleen | spleen stromal cell line | general cells |
| IFO50376 | MSS62 | mouse | spleen | spleen stromal cell line | general cells |
| IFO50377 | FLS3 | mouse | liver | liver stromal cell line | general cells |
| IFO50378 | FLS5 | mouse | liver | liver stromal cell line | general cells |
| IFO50379 | TLR2 | mouse | liver | TG mouse harborir hepatocyte immortalized with the tsSV | general cells |
| IFO50380 | TLR3 | mouse | liver | TG mouse harborir hepatocyte immortalized by the tsSV4C | general cells |

| | | | | | |
|----------|--------------------|-------------------|---|--|---|
| IFO50409 | NCTC clone 929 | mouse | skin | fibroblast, subclone of strain L | general cells |
| IFO50411 | A-431 | human | epidermoid carcinoma | epidermoid carcinoma, high expressio general cells | |
| IFO50412 | CCRF-CEM | human | hemo-lymphocytic | leukemia, acute lymphoblastic | general cells |
| IFO50414 | CHO-K1 | hamster, C ovary | | widely used for production of recombin | general cells |
| IFO50415 | C3H/10T1/2-clone8 | mouse | embryo | fibroblast, contact-inhibited | general cells |
| IFO50416 | 3T3-L1 | mouse | embryo | a substrain of 3T3- fibroblast, 3T3-Swiss derivative, adip | general cells |
| IFO50417 | 3T3-Swiss albino | mouse | embryo | fibroblast, contact-inhibited | general cells |
| IFO50418 | ASF-4-1 | human | skin | skin normal diploid fibroblast | general cells |
| IFO50419 | ASF-4-2 | human | skin | skin normal diploid fibroblast | general cells |
| IFO50421 | PKS | pig | kidney | susceptible to Vaccinia virus | general cells |
| IFO50422 | PKF | pig | kidney | susceptible to Vaccinia virus | general cells |
| IFO50423 | PKR | pig | kidney | susceptible to Vaccinia virus | general cells |
| IFO50424 | AP-16 | mouse | neural | astrocyte-progenitor cell line | general cells |
| IFO50425 | C6-SF2 | rat | neural | C6 subline adapted to serum-free med | general cells |
| IFO50426 | NIL-SF2 | hamster, S embryo | | NIL subline adapted to serum-free me | general cells |
| IFO50428 | CMK-86 | human | hemo-lym from patient with D leukemia, acute me | This cell line is use | general cells |
| IFO50430 | CMK-11-5 | human | hemo-lym from patient with D leukemia, acute me | This cell line is use | general cells |
| IFO50431 | KP-N-YN | human | neural | neuroblastoma, from lymph node, smo | general cells |
| IFO50432 | KP-N-RT-BM-1 | human | neural | neuroblastoma, from bone marrow, ne | general cells |
| IFO50433 | KP-N-SI9s | human | neural | neuroblastoma, from lymph node, smo | general cells |
| IFO50434 | KALS-1 | human | neural | glioma, GFAP, vimentin and CD13-pos | general cells |
| IFO50435 | KINGS-1 | human | neural | astrocytoma, anaplastic, GFAP, S-100 | general cells |
| IFO50436 | KS-1 | human | neural | glioblastoma | general cells |
| IFO50444 | KNS-81-FD | human | neural | serum-free adapti glioma, derivative c | The DNA fingerprint |
| IFO50466 | RC1 | mouse | antibody-p hybridoma (mouse | RC1 antibody, recognizing undifferenti | general cells |
| IFO50467 | NMC-G1 | human | neural | glioma, FGF-9-producing | general cells |
| IFO50472 | MEB5 | mouse | neural | immortalized by E7 CNS stem cell line, HPV type 16 E7-irr | general cells |
| IFO50473 | MEG-01s | human | hemo-lym suspension culture | leukemia, chronic megakaryoblastic, IV | general cells |
| IFO50474 | NOMO-1 | human | hemo-lymphocytic | leukemia, acute monocytic, promonoc | general cells |
| IFO50475 | NOMO-1s | human | hemo-lym a subline of NOMC | leukemia, acute monocytic, NOMO-1-c | general cells |
| IFO50476 | NKM-1 | human | hemo-lymphocytic | leukemia, acute myeloid, myeloid leuke | general cells |
| IFO50477 | FS-1 | human | bone marrow stroma | bone marrow strom Although this cell li | general cells |
| IFO50478 | MEG-A2 | human | hemo-lym Phl. chromosome | leukemia, chronic myelogenous, mega | general cells |
| IFO50479 | NAGL-1 | human | hemo-lymphocytic | leukemia, B cells | general cells |
| IFO50480 | NRK-52E | rat | kidney | epithelioid clone of NRK cells | general cells |
| IFO50481 | NRK-49F | rat | kidney | fibroblastic clone of NRK cells, useful f | general cells |
| IFO50483 | CCD-33Co | human | colon | established from n The ampoules pres | general cells |
| IFO50484 | DBC1.2 | mouse | olfactory epithelium | dark (horizontal) basal cell-like, keratin | general cells |
| IFO50488 | OUMS-27 | human | sarcoma | mutation of the p53 chondrosarcoma type 1, 2 and 3 colla | general cells |
| IFO50489 | GB-1 | human | neural | glioblastoma, MDR1 and P-glycoprotei | general cells |
| IFO50490 | MDM2-RNB | rat | neural | human MDM2 gen human MDM2 gene-introduced neonel | general cells |
| IFO50491 | RNB | rat | neural | neonatal rat astrocytes, GFAP-positive | general cells |
| IFO50492 | AM-38 | human | neural | glioblastoma, GFAP and S-100-positiv | general cells |
| IFO50493 | YH-13 | human | neural | glioblastoma, GFAP and S-100-positiv | general cells |
| IFO50495 | Neuro-2aTG | mouse | neural | neuroblastoma, HPRT-deficient Neuro | general cells |
| IFO50496 | PC-12TG | rat | neural | HPRT-deficient pheochromocytoma, HPRT-deficient P | general cells |
| IFO50513 | MTA | human | hemo-lymphocytic | leukemia, natural killer-like T cell line | general cells |
| IFO50516 | TMD5 | human | hemo-lymphocytic | leukemia, acute B I Mycoplasma contz | general cells |
| IFO50518 | KAI3 | human | hemo-lymphocytic | NK-like cell line from a patient with sev | general cells |
| IFO50519 | A1 | mouse | neural | Established from p p53-deficient cell li | This cell line is use |
| IFO50520 | MG5 | mouse | neural | Established from p microglial cell line from p53-deficient n | general cells |
| IFO50525 | NCO2 | human | hemo-lymphocytic | | general cells |
| JCRB0001 | F9-41 | mouse | teratoma | cells are differentia Maximum cell dens | general cells |
| JCRB0003 | C3H/10T1/2-clone 8 | mouse | embryo | contact sensitive cell line | general cells |
| JCRB0004 | A-431 | human | epidermoid epidermoid carcino | epidermoid carcinoma | general cells |
| JCRB0005 | FRSK | rat | skin | epidermal cells | Doubling time = 15 |
| JCRB0006 | HL60RG | human | hemo-lym DNA prifiling of this HL60 RG grow fast RG:rapid growth. A | general cells | |
| JCRB0007 | Pt K2 | marsupial | kidney | kidney PtK2 was used in s The cell line may b | general cells |
| JCRB0009 | P3/NS1/1-Ag4-1 | mouse | hemo-lym | resistant to 10^-4 in non-secreting mou | Used for the prepa |
| JCRB0011 | S.CET.M8.1.1 | mouse | antibody-p IgG producing hybi | Release monoclon The cell line releas | general cells |
| JCRB0013 | 16-8 | mouse | antibody-p Produce IgM. Aggl | Producing monoclon producing monoclo | general cells |
| JCRB0014 | 19-7 | mouse | antibody-p Producing IgM. Ag | Producing monoclon Production of mon | general cells |
| JCRB0016 | P388-D1 | mouse | hemo-lym ascitic form of P38 | Used by the NCI c | Release interleukin |
| JCRB0017 | P388 | mouse | hemo-lym | Low level IL1 prod This cell line is use | Target cell by cytot |
| JCRB0018 | J774.1 | mouse | hemo-lym | Produce large amo | Carry receptors for immunoglobulin am |
| JCRB0019 | K562 | human | hemo-lymphocytic | Differentiate into er | Differentiation with |
| JCRB0021 | MOLT-4F | human | hemo-lymphocytic | Rosettes formation. | High TdT. No Ig n |
| JCRB0023 | RBL-2H3 | rat | hemo-lym mast cell line | IgE mediated histai | Release IgE-medi |
| JCRB0024 | IM-9 | human | hemo-lym B lymphoblasts | syr Immunoglobulin se | The cell line has re |
| JCRB0025 | BRL-3A | rat | liver | liver | Producing MSA pr Production of some |
| JCRB0030 | CHL/IU(IVGT) | Chinese h | lung | Standard cell line f Same cell line in E | Only for researchers in Asian countries |
| JCRB0031 | CCRF-HSB2 | human | hemo-lymphocytic | Cultured in vitro fro | After CCRF-SB wa |
| JCRB0032 | CCRF-SB | human | hemo-lym | Normal human dipl | This cell line was e B-cell |
| JCRB0033 | CCRF-CEM | human | hemo-lymphocytic | Near diploid, no Ig-secreting, free of vi | general cells |
| JCRB0034 | RPMM 8226 | human | hemo-lymphocytic | hematopoietic cell line, Lambda L chal | general cells |
| JCRB0035 | RPMM 1788 | human | hemo-lym normal | IgM secreting(light chain), positive for | general cells |
| JCRB0041 | HLCL-1 | human | hemo-lymphocytic | peripheral lymphocyte | general cells |
| JCRB0042 | HEC-1 | human | uterus | STR analysis in the cell bank indicated | Mycoplasma conta |
| JCRB0044 | LU99A | human | lung (cancer) | produce colony-stimulating factor | general cells |
| JCRB0054 | LU65A | human | lung (cancer) | | general cells |

| | | | | | |
|------------|-------------------|------------------|---|--|---|
| JCRB0055 | LU65B | human | lung (cancer) | | general cells |
| JCRB0056 | LU65C | human | lung (canc lung, carcinoma | | general cells |
| JCRB0057 | LU99B | human | lung (cancer) | | general cells |
| JCRB0058 | LU99C | human | lung (cancer) | | general cells |
| JCRB0060 | LLC-PK1 | pig | kidney | susceptible to viruses, rubella,vaccinia | general cells |
| JCRB0061 | AZ-521 | human | stomach | Mycoplasma conta | general cells |
| JCRB0062 | HEL | human | hemo-lymphocytic | | general cells |
| JCRB0065 | KG-1 | human | hemo-lymphocytic | | general cells |
| JCRB0066 | Mewo | human | melanoma | | general cells |
| JCRB0068 | Flow3000 | human | embryonic normal fibroblast | sensitive to CMV virus | general cells |
| JCRB0070 | MIA PaCa-2 | human | pancreas | sensitive to asparginase. Dt=40hr, colk | general cells |
| JCRB0071 | BALL-1 | human | hemo-lymphocytic | | general cells |
| JCRB0072 | MiTen | human | skin | | general cells |
| JCRB0073 | J-111 | human | HeLa-cont HeLa marker chror | Although an origin HeLa markers obse | general cells |
| JCRB0074 | VMRC-MELG | human | melanoma | | general cells |
| JCRB0075 | SF-TY | human | skin normal | normal skin fibrobz HLA:A2,A24,B5,Bv | general cells |
| JCRB0076 | A549 | human | lung (canc adenocarcinoma | Y chr.invisible. Detected by Y specific | general cells |
| JCRB0077 | PC-3 | human | lung (canc intermediately diffe | notice:different cell | Mycoplasma conta |
| JCRB0078 | P-815 | mouse | hemo-lym mastocytoma | target cell for cytotoxic T cell assay | general cells |
| JCRB0079 | LU65 | human | lung (canc c-Ki-ras2 mut. cod many floating | cells are observed | general cells |
| JCRB0080 | LU99 | human | lung (cancer) | lung, carcinoma | general cells |
| JCRB0081 | RERF-LC-MS | human | lung (canc adenocarcinoma | lung adenocarcinoma | general cells |
| JCRB0083 | 4-8H | mouse | antibody-p hybridoma | antibody to DNA px | Everyday medium |
| JCRB0084 | 4-2D | mouse | antibody-p hybridoma, produc | antibody to primas | Everyday medium |
| JCRB0085 | HL60 | human | hemo-lym APL | DMSO, butyl.,HX.,T Doubling time is 40 | general cells |
| JCRB0086 | TALL-1 | human | hemo-lym T cell | T cell lymphoblastic leukemia, mode=9 | general cells |
| JCRB0090 | MB86 | mouse | antibody-p producing antibody | recognize IgH6-4 | This cell line produ |
| JCRB0091 | P31/FUJ | human | hemo-lym monocyte, 47:XY, I Pox+ PAS+ AcPas | Lot of FCS affects | general cells |
| JCRB0094 | P30/OHK | human | hemo-lym CD5-,CD2-,la+,CD PAS+,AcPase+,ha | CD8 antigen is indi | general cells |
| JCRB0095 | P32/ISH | human | hemo-lym B cell characteristi | translocation betw Good for the resea | general cells |
| JCRB0097 | A4/Fuk | human | hemo-lym ER-,PanT-,CD4+,F | lymphoma cell line | Mycoplasma conta |
| JCRB0098 | KURAMOCHI | human | ovary hyperdiploid | Dt=24-27hr | Mycoplasma conta |
| JCRB0101 | A3/KAW | human | hemo-lym CD4+,la-,EBNA- | Dt=20 hr | Mycoplasma was e |
| JCRB0102 | RERF-LC-FM | human | lung (cancer) | create clusters on growth | general cells |
| JCRB0103 | VMRC-LCP | human | lung (cancer) | lung squamous cell carcinoma | general cells |
| JCRB0104 | KU812 | human | hemo-lym Ph1 chr., histamin(| Dt=20-30hr, colony Ph1 chromosome | (general cells |
| JCRB0104.1 | KU812E | human | hemo-lym Ph1 chr., histamin(| Dt=20-30hr, clonal | This cell line was d |
| JCRB0104.2 | KU812F | human | hemo-lym Ph1 chr., histamin(| Dt=20-30hr, clonal | This cell line was d |
| JCRB0105 | KY821 | human | hemo-lym Double minute chr. | Dt=27-37 hr | general cells |
| JCRB0105.1 | KY821A3 | human | hemo-lym leukemia | The cell was clone: Cell ID data indicat | general cells |
| JCRB0106 | SCCH-26 | human | neural neuroblastoma | neuroblastoma | Other name was di |
| JCRB0108 | SCCH-196 | human | sarcoma t(11;22) | small round cell sa | Other name was di |
| JCRB0110 | M1-T22 | mouse | hemo-lym cytologically determ | Cells are differentiated by dexamethas | general cells |
| JCRB0112 | THP-1 | human | hemo-lym Fc,C3b receptors. | No immunoglobulin HLA-DR,My 7, My | general cells |
| JCRB0112.1 | THP-1 | human | hemo-lym Fc,C3b receptor.Hi | No immuno gloguli HLA-DR, My7, My | general cells |
| JCRB0113 | PAI | mouse | hemo-lymphocytic | non screeting type for IgE. | general cells |
| JCRB0114 | LC4-1 | human | hemo-lym positive la,B4, neg | E-rosette formation. D-J joining. | EBI general cells |
| JCRB0115 | SCC-3 | human | hemo-lym LSG type. | IL2 rece monocytic cell line. Negative for ATLA, | general cells |
| JCRB0116.1 | U937 cl1-14 | human | hemo-lym express monocyte- highly susceptible | clone for HIV virus | general cells |
| JCRB0116.2 | U937 cl1-22 | human | hemo-lym express many mon low suscep | ble clo Genetic ID was the | general cells |
| JCRB0117 | B104 | human | hemo-lym IgM and IgD are ex | The cells were killed by anti-IgM but nc | general cells |
| JCRB0118 | SKM-1 | human | hemo-lym MDS->RAEB.T->A | Myeloperoxidase (MPO) release into c | general cells |
| JCRB0119 | YG10003 | hamster, C lung | salmonella O-acety | G418-r(500ug/ml), This cells are comr | general cells |
| JCRB0120 | YG10007 | hamster, C lung | human monomorp | G418-r(500ug/ml), This cells are comr | general cells |
| JCRB0121 | YG10008 | hamster, C lung | human polymorphi | G418-r(500ug/ml), This cells are comr | general cells |
| JCRB0123 | KO52 | human | hemo-lym CD12,CD33,CD2,(N-ras mutation: | G->C in codon 13. | general cells |
| JCRB0124 | Takigawa | human | stomach | Densely packed let-like colonies with sr | general cells |
| JCRB0125 | L alpha | mouse | skin | Mouse L tk- cell lin | This cell line expr: Receptors are exa |
| JCRB0126.1 | KOSC-2 cl3-43 | human | oral | Moderately differer Cells were establis | general cells |
| JCRB0129.0 | RCR-1.P3 | rat | neural | GFAP (glial fibrillar Cells having characteristics of astrocyt | general cells |
| JCRB0129.1 | RCR-1 | rat | neural | GFAP (glial fibrillar Cells having chracil Cells obtained fror | general cells |
| JCRB0129.2 | RCR-1.P3 | rat | brain | GFAP (glial fibrillar Cells having characteristics of astrocyt | general cells |
| JCRB0130 | PC-12.P3 | rat | neural | Derivative of the P Cells adapted to grow in serum free m | general cells |
| JCRB0130.1 | PC12.P3 | rat | neural | | general cells |
| JCRB0131 | RL-33 | rabbit | lung | HSV,VSV sensitive | Cell growth inhibited by cell-cell contac |
| JCRB0132.1 | AH601(JTC27) | rat | liver | hepatoma | transplantable rat ascites tumor cell lin |
| JCRB0132.2 | AH601.P3(JTC27) | rat | liver | AH-601 cells adapl | transplantable rat ascites tumor P3 cel |
| JCRB0132.3 | AH601.P3(JTC27) | rat | liver | AH-601 cells adapl | transplantable rat ascites tumor P3 cel |
| JCRB0133 | KHM-1B | human | hemo-lym HLA-DR,CD2,CD1 | Amylase and IgA productions were po: | general cells |
| JCRB0134 | MCF-7 | human | mammary | dome formation observed | general cells |
| JCRB0136.1 | CHO(pMAM-luc) | hamster, C ovary | Plasmid pMAM-luc | Luciferase gene ur See CHO-K1 cell c | general cells |
| JCRB0136.2 | CHO(pMAM-HSluc) | hamster, C ovary | pMAM-HSluc trans | HSP 70B gene pro Refer CHO-K1 cell | general cells |
| JCRB0137 | RERF-LC-KJ | human | lung (cancer) | Transplantable to SCID mice and freq | general cells |
| JCRB0138 | KHM-3S | human | lung (canc No OKT11,OKT4,C | HTLV-1 positive small cell lung cancer | general cells |
| JCRB0139 | TASK1 | human | neural | NSE,vimentin, & M | Primitive neuroect Cells are originally |
| JCRB0140 | NCE16 | human | uterine cei | E6/E7 expressed. | Normal ecto-cervic Any comments for |
| JCRB0141 | PHK16-0b | human | skin | E6/E7 expressed a | Cells immortalized Need details of the |
| JCRB0143.0 | F2408 pZipNeo | rat | embryo | No expression of t fibroblast-like | general cells |
| JCRB0143.1 | F2408 EBNA2 cl-9 | rat | embryo | Expression of the t fibroblast-like | general cells |
| JCRB0143.2 | F2408 EBNA2 cl-16 | rat | embryo | Expression of the t fibroblast-like | general cells |

| | | | | | |
|-------------|-------------------|-------------------|-----------------|--|--|
| JCRB0143.3 | F2408 EBNA2 cl-18 | rat | embryo | Expression of the fibroblast-like | general cells |
| JCRB0143.4 | F2408 EBNA2 cl-24 | rat | embryo | Expression of the fibroblast-like | Cell culture becomes general cells |
| JCRB0143.5 | F2408 EBNA2 cl-14 | rat | embryo | Expression of the fibroblast-like | general cells |
| JCRB0143.6 | F2408 EBNA2 cl-23 | rat | embryo | Expression of the fibroblast-like | general cells |
| JCRB0143.7 | F2408 EBNA2 cl-28 | rat | embryo | Expression of the fibroblast-like | general cells |
| JCRB0144.1 | LYM-1.P3 | rat | lymph node | Sticky cell from lymph node. Serum free | general cells |
| JCRB0144.1B | LYM-1.P3B | rat | lymph nod | Subclone of the LY Culture fluid of the LYM-1 strongly inhibits | general cells |
| JCRB0144.1C | LYM-1.P3D | rat | lymph nod | Subclone of the LY Culture fluid of the LYM-1 strongly inhibits | general cells |
| JCRB0144.2C | LYM-1.P3D | rat | lymph nod | Subclone of the LY Culture fluid of the LYM-1 | general cells |
| JCRB0145 | MEB 5 | mouse | neural | cells immortalized Multipotent CNS stem cell line. Differentiation | general cells |
| JCRB0146 | MLMA | human | hemo-lym | p47 X-X (X,0) IgD, IgM positive, hairy B-cell line | general cells |
| JCRB0147 | JKT-beta-del | human | hemo-lym | Deletion of rearranged Jurkat cell clone with a lack of surface | general cells |
| JCRB0148 | KHM-5M | human | thyroid | Cells are derived from Neutrophil chemotaxis Cells are decongestant | general cells |
| JCRB0149 | Bhas 42 | mouse | embryo | H-ras gene transfectants Growth on the dish | Appearing of natural general cells |
| JCRB0150 | P53LMAC01 | mouse | vascular | s p53 deficient(-/-) | smooth muscle cell line from p53 knock-out |
| JCRB0151 | T-Ag-MOSE | mouse | ovary | pEF321-T plasmid | immortalized by T ± 5% CO ₂ , 37 C cult |
| JCRB0151.1 | p53-def-MOSE | mouse | ovary | p53 gene was homozygous non-tumorigenic | Adhesion efficiency |
| JCRB0152.1 | M | rat | liver | 4-dimethylaminoaz | Morphological characteristics Reference cell line |
| JCRB0152.2 | M.P3 | rat | liver | 4-dimethylaminoaz | Morphological characteristics Cells grow in serum |
| JCRB0152.3 | M.P3 | rat | liver | 4-dimethylaminoaz | Morphological characteristics Cells grow in serum |
| JCRB0153.1 | tsJ663 | rat | embryo | A ts mutant for cell MNNG-mutagenized | 3Y1 cell derivative |
| JCRB0153.2 | tsJT16 | rat | embryo | A ts mutant for cell MNNG-mutagenized rat | 3Y1 derivative |
| JCRB0153.3 | tsJT60 | rat | embryo | A ts mutant for cell 3Y1 derived ts mutant cell line with MN | general cells |
| JCRB0154 | NB(TU)1-10 | human | neural | N-myc aberration | c Neuroblastoma cell Cells will be available |
| JCRB0156 | KHYG-1 | human | hemo-lym | p53 point mutation | high NK/LAK activity, IFN gamma prod |
| JCRB0157 | TK | human | hemo-lym | CD10, CD19, CD22 | A rare large B-cell non-Hodgkin's lymph |
| JCRB0158.0 | MY | human | hemo-lym | Chimeric BCR/ABL Bone marrow mononuclear cells. | general cells |
| JCRB0158.1 | MY-M12 | human | hemo-lym | Chimeric BCR/ABL Bone marrow mon | Please refer NIHSC |
| JCRB0158.2 | MY-M13 | human | hemo-lym | Chimeric BCR/ABL Bone marrow mononuclear cells. | general cells |
| JCRB0159 | SLVL | human | hemo-lym | Clonal rearrangement B cell line with cyto A part of the population | general cells |
| JCRB0160 | L190 | human | liver, gallbl | Karyology: 47XX, 6q-, +22 | general cells |
| JCRB0161 | Kasumi-4 | human | blood | t(9;22:11), overexp | CD34+, CD33+, CD13+ |
| JCRB0162 | NEC14 | human | testis | testicular germ cell testicular germ cell | Mycoplasma-eliminated |
| JCRB0163 | HL60(S) | human | hemo-lym | APL | DMSO, butyl, HX, STR-PCR results |
| JCRB0164 | HTC/C3 | human | thyroid | thyroid carcinoma, thyroid carcinoma, The HTC/C3 is un | general cells |
| JCRB0165 | KHM-10B | human | hemo-lym | HLA-DR, CD19 and Burkitt's lymphoma, The KHM-10B is u | general cells |
| JCRB0166 | HMC-1-8 | human | mammary | HMC-1-8 is a target cell of the T-cell k | general cells |
| JCRB0167 | RCN-9 | rat | colon | DMH induced colo | DMH induced colo Low differentiated |
| JCRB0168 | c-WRT-7-LR | rat | hemo-lym | Surface markers: F c-WRT-7 cells were WKA/Hok rat. c-WI | general cells |
| JCRB0169 | Lu-134-B | human | lung | (cancer Activities of gamma: Derived from small cell lung carcinoma | general cells |
| JCRB0170 | Lu-135 | human | lung | (cancer specific antigen Hard to observe dense core granules | general cells |
| JCRB0172 | RMG-I | human | ovary | tumor markers = C Doubling time = 60 Cell culture initiated | general cells |
| JCRB0172.1 | RMG-II | human | ovary | Chromosomes are Doubling time is ar Cells confirmed to | general cells |
| JCRB0173 | SKN | human | uterus | Mode of chromoso | Histochemical characteristics Cell culture started |
| JCRB0174.0 | NCC16-P11 | human | uterine cervix | E6/E7 positive. Tel | Columnar epithelium The NCC16-P11 w |
| JCRB0176 | RKN | human | ovary | mode of chromoso | Doubling time=31.2 Cells were contaminated |
| JCRB0178.0 | KP-3 | human | pancreas | A metastasis of adi | Unique bat the pattern |
| JCRB0178.1 | KP-3L | human | pancreas | Tissue was classifi | Cells were established Mycoplasmas were |
| JCRB0179 | SNG-M | human | uterus | Mordal No. of chro | intermediately different The cells are resp |
| JCRB0180 | GAK | human | melanoma | Mode of chromoso | The cell line GAK has an activity of 5- |
| JCRB0181 | KP-2 | human | pancreas | Mode of chromoso | The tumor cell line Cells are observed |
| JCRB0182 | KP-4 | human | pancreas | Tumor cell line sec | Cells are observed |
| JCRB0183 | QGP-1 | human | pancreas | Modal chromosom | Cell line produces Cells are observed |
| JCRB0186 | OV3121ras4 | mouse | ovary | v-Ha-ras gene tran | Ovarian granulosa cell line with v-Ha-r |
| JCRB0187 | OV3121ras7 | mouse | ovary | v-Ha-ras gene tran | Ovarian granulosa cell line with v-Ha- |
| JCRB0191 | OCUG-1 | human | liver, gallbl | Chromosome numl Only TA-4 antigen | Cells are tumoriger |
| JCRB0192 | OCUM-1 | human | stomach | karyotype=mode 5 This cell line produces | carcinogenesis |
| JCRB0193 | WB-F344 | rat | liver | Normal diploid cell | Normal rat liver epithelial cell line grew |
| JCRB0194 | KON | human | oral | Mode of chromoso | Td=32.5 hrs. Color Cells were unique |
| JCRB0195 | RP-1 | mouse | antibody-p | Rat neutrophil spe | This cell line is producing IgG1 against |
| JCRB0196 | RP-3 | mouse | antibody-p | Hybridoma X63 as | Mouse cell line producing IgM antibody |
| JCRB0197 | OSC-20 | human | oral | Cells grew in cobbl | Serum free cell cul |
| JCRB0198 | OSC-19 | human | oral | EGF receptor. | EGF (1-100ng/ml) inhibited the cell growth |
| JCRB0199 | huH-1 | human | liver, gallbl | Modal chromosom | The huH-1 cell line produce tumors in |
| JCRB0200 | KMLS-1 | human | retroperitoneum | | general cells |
| JCRB0202 | B16 melanoma | mouse | melanoma | positive for melanin | Melanin producing |
| JCRB0203 | PCC4 Typell | mouse x r: cybrid | | PCC4/L6TG Cap-r | PCC4/L6TG Cap-r |
| JCRB0206 | PCC4 AG | mouse | teratoma | testis, AG-r | testis, AG-r. Cells resistant to 8-azac |
| JCRB0207 | CaR-1 | human | rectum | | general cells |
| JCRB0208 | CCK-81 | human | colon | large intestine | carcinoembryonic antigen producing |
| JCRB0209 | L6TG | rat | muscular | : myeloblast, TG-r | myeloblast, TG-r |
| JCRB0210 | L6TG Cap | rat | muscular | : rat skeletal muscle | rat skeletal muscle myoblast, TG-r, Ca |
| JCRB0211 | A9 | mouse | skin | HPRT and skin fibroblast, AG- | Resistant to 8-azac |
| JCRB0212 | B82 | mouse | skin | skin fibroblast, Brd | fibroblast Resistant to BrdU |
| JCRB0213 | HeLa AG | human | uterine cervix | , AG-r | cervix, AG-r |
| JCRB0214 | HeLa TG | human | uterine cervix | , TG-r | cervix, TG-r |
| JCRB0215 | HeLa TG Cap | human | uterine cervix | , TG-r, Cap-r | cervix, TG-r, Cap-r |
| JCRB0216 | PCC4 AG Cap | mouse | teratoma | testis, AG-r, Cap-r | testis, AG-r, Cap-r |
| JCRB0217 | DSPT200 | hamster, S | embryo | fetus fibroblast | trar fetus fibroblast transformed cell |
| JCRB0218 | CHO/HGPRT | hamster, Covary | | HPRT deficient | HTRT deficient mur Because of the HG |

| | | | | | | |
|------------|-----------------|-------------------------------------|---|---|---|------------------|
| JCRB0219 | V79-6TG | hamster, C lung | lung, 6TG-r | lung, 6TG-r | general cells | |
| JCRB0220 | NTI-4 | human embryo | fetus body | fetus, half body, bo | Normal human fibr general cells | N/A via Xenotech |
| JCRB0221 | HT100 | hamster, E lung | lung, new born | lung, new born | general cells | |
| JCRB0222 | NTI-5 | human embryo | fetus, half a body | Normal human cell PDL=13 | general cells | N/A via Xenotech |
| JCRB0224 | WiDr | human colon | colon, adenocarcin colon, adenocarcinoma | colon, adenocarcin colon, adenocarcinoma | general cells | |
| JCRB0225 | COLO320 DM | human colon | colon, adenocarcin colon, adenocarcinoma | colon, adenocarcin colon, adenocarcinoma | general cells | |
| JCRB0226 | COLO201 | human colon | colon, adenocarcin colon, adenocarcinoma | colon, adenocarcin colon, adenocarcinoma | general cells | |
| JCRB0227 | C32TG | human melanoma | amelanotic melanoma | amelanotic melanoma, 6TG-r | general cells | |
| JCRB0228 | A-172 | human neural | glioblastoma | glioblastoma | general cells | |
| JCRB0229 | EGV-4T | rat lung | immortal cell line fr rat tracheal epithelial cells | immortal cell line fr rat tracheal epithelial cells | general cells | |
| JCRB0230 | SCC-131 | rat squamous | squamous cell car | squamous cell carc The cell line was ol | general cells | |
| JCRB0231 | SCC-158 | rat squamous | squamous cell car | squamous cell carc The cell line was ol | general cells | |
| JCRB0232 | SKG-IIla | human uterine cer cervix | Growth rate is slow | Tumour markers, T | general cells | |
| JCRB0232.1 | SKG-IIlb | human uterine cer tumor markers, TA | The cell line was e: The cell line was cx | general cells | general cells | |
| JCRB0233.0 | BALB-MC | mouse mammary | mammary cancer | mammary cancer | general cells | |
| JCRB0233.1 | BALB-MC.D3 | mouse mammary | mammary cancer | mammary cancer | general cells | |
| JCRB0233.2 | BALB-MC.E12 | mouse mammary | mammary cancer | mammary cancer | general cells | |
| JCRB0234.0 | TYK-nu | human ovary | ovary, undifferenti | ovary, undifferentiated carcinoma | general cells | |
| JCRB0234.1 | TYK-nu.CP-r | human ovary | ovary, undifferenti | ovary, undifferentia Sysplatin resistant. | general cells | |
| JCRB0235 | Lu-134-A-H | human lung | (canc small cell carcinom | small cell carcinom Cell number of the | general cells | |
| JCRB0236 | KG-1-C | human neural | mixed glioma | mixed glioma | general cells | |
| JCRB0238 | B8/3 | mouse hemo-lymphocytic | | Possible model sys | general cells | |
| JCRB0239 | TCO-1 | human thyroid | | non differentiated cancer | general cells | |
| JCRB0240 | MCAS | human ovary | | | general cells | |
| JCRB0241 | Shay | rat hemo-lymphocytic | | | general cells | |
| JCRB0242 | RMT-1 E4 | rat mammary | | rat mammary carci Meium Supplements | general cells | |
| JCRB0243 | B-1 | mouse testis | | estrogen dependent Leydig cell tumor | general cells | |
| JCRB0245 | RMT-1 M2 | rat mammary | rat mammary myoe | Cells were prepared from the tumor. | general cells | |
| JCRB0246 | RC-1 | rabbit cornea | cornea | cornea | general cells | |
| JCRB0247 | RL-34 | rat liver | liver | Cell growth stoppe Eliminated mycopl | general cells | |
| JCRB0248 | ARLJ301-3 | rat liver | liver epithelial cell | liver epithelial cell 10-week-old male | general cells | |
| JCRB0249 | FAA-HTC1 | rat liver | 2-AAF induced hef | 2-AAF induced hef Male, F344/DuCrj | general cells | |
| JCRB0250 | NEC8 | human testis | testicular germ cell | testicular germ cell Mycoplasma-elimir | general cells | |
| JCRB0251 | SCH | human stomach | stomach, choriocar | stomach, choriocar Mycoplasmas elimi | general cells | |
| JCRB0252 | MKN1 | human stomach | stomach, adenosq | stomach, adenosq Mycoplasma-elimir | general cells | |
| JCRB0254 | MKN45 | human stomach | stomach, 62 year-c | stomach, 62 year-c Poorly differentiate | general cells | |
| JCRB0255 | MKN74 | human stomach | | Moderately differen | general cells | |
| JCRB0256 | RCM-1 | human rectum | rectum, adenocarc | rectum, adenocarc 73 year-old, female | general cells | |
| JCRB0257 | CoCM-1 | human colon | descending colon, | Moderately differen | general cells | |
| JCRB0258 | LC-1 sq | human lung | (canc lung, squamous ca | lung, squamous ca 69 year-old, male, | general cells | |
| JCRB0259 | MSK | rat sarcoma | 32P-induced osteo | 32P-induced osteo High tumorigenecity | general cells | |
| JCRB0260 | SAS | human oral | tongue, squamous tongue, | squamous Derived from prima | general cells | |
| JCRB0261 | T.Tn | human esophagu | High alkaline phos | Squamous carcino Cells identified to b | general cells | |
| JCRB0262 | T.T | human esophagu | squamous carcinoi | 3p-, 6q- Mode=85- Derived from manc | general cells | |
| JCRB0263.1 | PC12HS.P3 | rat neural | Serum free cell cul | Cells are differentiated into sympatheti | general cells | |
| JCRB0266 | PC12HS | rat neural | NGF-highly sensitiv | Cells are differentiated into sympatheti | general cells | |
| JCRB0267 | PC12MS | rat neural | NGF-moderately si | Cells are differentiated into sympatheti | general cells | |
| JCRB0268 | PC12IS | rat neural | NGF-insensitive cl | Cells are differentiated into sympatheti | general cells | |
| JCRB0401 | HUH-6 Clone 5 | human liver, galb | hepatoblastoma, H | Produces albumin, The cell line grow ii | general cells | |
| JCRB0402 | RLN-B2 | rat liver | | normal liver | The cell line keeps | general cells |
| JCRB0403 | HuH-7 | human liver, galb | differenciated hepe | hepatoma (differen | This cell line can b | general cells |
| JCRB0404 | HLE | human liver, galb | hepatoma (non-diff | hepatoma (non-diff No protein product | general cells | |
| JCRB0405 | HLF | human liver, galb | hepatoma, non-diff hepatoma, | non-differentiated, age 68, | general cells | |
| JCRB0406 | PLC/PRF/5 | human liver, galb | Loss of D & G gro | G6PD=Type A. Tui | Grow in serum free | general cells |
| JCRB0407 | RLN-J-5-2 | rat liver | normal liver, 7 day | normal liver, 7 days | general cells | |
| JCRB0408 | Ac2F | rat liver | normal liver, diploic | normal liver, diploic | Normality is mainta | general cells |
| JCRB0409 | dRLa-74 Clone 2 | rat liver | liver adenoma | liver adenoma | general cells | |
| JCRB0410 | dRLh-84 | rat liver | hepatoma | hepatoma | Cell to cell connect | general cells |
| JCRB0411 | AH66tc | rat liver | | Yoshida ascites he | Yoshida ascites he High production of | general cells |
| JCRB0412 | AH70Btc | rat liver | | AH70B Yoshida as AH70B Yoshida as alpha-fetoprotein | a general cells | |
| JCRB0413 | HuO-3N1 | human sarcoma | osteosarcoma, age | osteosarcoma, age | Osteoblastic cell lir | general cells |
| JCRB0414 | RLN-8 | rat liver | normal liver, 9 day | normal liver, 9 days | general cells | |
| JCRB0415 | RLN-10 | rat liver | normal liver, 14 da | normal liver, 14 days | general cells | |
| JCRB0416 | Ac2F(DT) | rat liver | liver, normal | liver, normal | Cells are derived fr | general cells |
| JCRB0417 | Ac2F(ST) | rat liver | liver, normal | liver, normal | Cells were derived | general cells |
| JCRB0418 | dRLN-4 | rat liver | liver after treatmen | liver after treatment by 0.06 % | DAB | general cells |
| JCRB0419 | dRLN-9 | rat liver | liver after treatmen | liver after treatment by 0.06 % | DAB for | general cells |
| JCRB0420 | dRLN-6 | rat liver | liver after treatmen | liver after treatment by 0.06 % | DAB for | general cells |
| JCRB0421 | 3'-mRLN-30 | rat liver | liver after 3'-Me-D/ | liver after 3'-Me-DA | 3'-Me-DAB treated | general cells |
| JCRB0422 | 3'-mRLN-31 | rat liver | liver after 3'-Me-D/ | liver after 3'-Me-DAB treatment for 61 | general cells | |
| JCRB0423 | RS-A | rat spleen | spontaneo | spontaneo | Chromosome numl | general cells |
| JCRB0424 | JTC-11 | mouse ascites | JTC-11 (K strain): | Easy to get monoc | general cells | |
| JCRB0425 | HuCCT1 | human liver, galb | bile duct carcinoma | bile duct carcinoma | general cells | |
| JCRB0426 | HuH28 | human liver, galb | bile duct carcinoma | bile duct carcinoma | Doubling time is ve | general cells |
| JCRB0427 | HuO9 | human sarcoma | Liver/bone/kidney-1 | Secretes osteocalcin & | responds to vit | general cells |
| JCRB0428 | HuO9N2 | human sarcoma | Vitamin D3 respon | Tumorigenic cell line. | Growth rate is ve | general cells |
| JCRB0429 | KMS-12-BM | human | hemato-lym | late stage of B-cell Non tumorigenic, | PCA-1(+),CD38(+),C | general cells |
| JCRB0430 | KMS-12-PE | human | hemato-lym | late stage B-cell dl | non-producing, transferrin receptors p: | general cells |
| JCRB0431 | 3'-mRLh-2 | rat liver | Gamma-glutamyl tr | A dexamethasone-responsive | tyrosine | general cells |
| JCRB0432 | KMS-6 | human embryo | Normal diploid cell | fibroblast, normal d | Change all mediurr | general cells |
| | | | | | | N/A via Xenotech |

| | | | | | | | |
|------------|--------------------|------------|----------------|---|--|-----------------------|------------------|
| JCRB0433 | KMST-6 | human | embryo | fibroblast, immortal | indefinite growth and non-tumorigenic | general cells | N/A via Xenotech |
| JCRB0435 | JHH-4 | human | liver, gallb | HBV integration is | Alpha-photoprotein Cell Bank will hand | general cells | |
| JCRB0501 | TIG-1-20 | human | fetus lung | fetal lung, (F) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0502 | TIG-1-30 | human | fetus lung | fetal lung, (F) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0503 | TIG-1-40 | human | fetus lung | fetal lung, (F) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0504 | TIG-1-50 | human | fetus lung | fetal lung, (F) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0505 | TIG-1-60 | human | fetus lung | fetal lung, (F) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0506 | TIG-3-20 | human | fetus lung | fetal lung, (M) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0506.1 | SVts-8 | human | lung | TIG-3 cells(JCRB0 | Human normal fibr Culture at 34 C be | general cells | |
| JCRB0511 | TIG-7-20 | human | fetus lung | fetal lung, (M) | fetal lung, normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0518 | WI-38-40 | human | lung | lung, normal diploid | lung, normal diploid fibroblast | general cells | |
| JCRB0519 | IMR-90-40 | human | lung | | PDL=Ca 50 | general cells | |
| JCRB0521 | MRC-5-30 | human | fetus lung | lung,normal diploid fetal lung, | normal diploid fibroblast, TIC general | cells | N/A via Xenotech |
| JCRB0525 | TIG-2M-30 | human | muscular | : muscle, normal dip muscle, normal dip | Refer attached inst | general cells | |
| JCRB0526 | TIG-101 | human | skin | skin, normal diploic skin, normal diploid | fibroblast | general cells | |
| JCRB0527 | TIG-102 | human | skin | skin, normal diploic skin, normal diploid | Life span = PD41 | general cells | |
| JCRB0528 | TIG-103 | human | skin | skin,normal diploid | skin,normal diploid fibroblast | general cells | |
| JCRB0532 | TIG-107 | human | skin | skin, normal diploic skin, normal diploid | fibroblast | general cells | |
| JCRB0533 | TIG-112 | human | skin | Normal diploid, | 2N: Normal skin fibroblast cells. PD at distr | general cells | |
| JCRB0534 | TIG-114 | human | skin | Normal diploid | 2N= Normal human skin fibroblast cells. | general cells | |
| JCRB0535 | TIG-118 | human | skin | | normal | general cells | |
| JCRB0536 | TIG-121 | human | skin | | normal diploid | general cells | |
| JCRB0537 | TIG-108 | human | skin | normal diploid, 2N | : normal fibroblast | general cells | |
| JCRB0538 | TIG-109 | human | skin | normal fibroblast, 2 | normal diploid human skin | general cells | |
| JCRB0539 | TIG-113 | human | skin | 2n=46, XX | Maximum PD = 53. Cells may be distri | general cells | |
| JCRB0540 | TIG-119 | human | skin | 2N=46, XY | Maximum PD = 48. Cells may be distri | general cells | |
| JCRB0541 | TIG-111 | human | skin | normal diploid fibro | normal diploid | general cells | |
| JCRB0542 | TIG-120 | human | skin | normal diploid (2N | = normal diploid fibro NT | general cells | |
| JCRB0543 | TIG-110 | human | skin | 2n=46, XX | normal diploid fibroblast, PD35 | general cells | |
| JCRB0544 | TIG-3S | human | skin | 2n=46, XY | lifespan PD67 | general cells | |
| JCRB0601 | BALB/3T3 A31-1-1 | mouse | embryo | whole embryo | contact inhibited m The FCS requires | general cells | |
| JCRB0603 | V79 | hamster, C | lung | lung | lung | Spontaneous muta | general cells |
| JCRB0605 | L.P3(S) | mouse | skin | | Serum free cell line Culturing in the me | general cells | |
| JCRB0606 | HeLa.P3(S) | human | uterine cervix | cervix, grow in serum free medium | but general | cells | |
| JCRB0608 | JTC19 | rat | lung | lung | lung | general cells | |
| JCRB0609.1 | Mm2T.P3 | Indian Mu | thymus | Chromosome ana | Cells grew when started at 1977 but di | general cells | |
| JCRB0610 | IAR-20 | rat | liver | adult liver | Non-transformed rat liver epithelial cel | general cells | |
| JCRB0611 | KATOIII | human | stomach | stomach cancer, si | stomach cancer, signet ring cell carcin | general cells | |
| JCRB0612 | GOTO | human | neural | neuroblastoma | neuroblastoma | Mycoplasma conta | general cells |
| JCRB0612.1 | GOTO.P3 | human | neural | neroblastoma | GOTO cells cultured with serum free n | general cells | |
| JCRB0613 | ITO-II | human | testis | | F9-like antigen and Collagen-coated s | general cells | |
| JCRB0614 | NY | human | sarcoma | osteosarcoma | osteosarcoma | general cells | |
| JCRB0615 | NIH/3T3 clone 5611 | mouse | embryo | whole embryo | whole embryo, contact inhibited | general cells | |
| JCRB0618 | TGW | human | neural | | neuroblastoma | general cells | |
| JCRB0618.1 | TGW.P3 | human | neural | | | general cells | |
| JCRB0621 | NB-1 | human | neural | neuroblastoma (M) | Differentiate by (Bu | Mycoplasma conta | general cells |
| JCRB0622 | HSC-2 | human | oral | | This cell line has neither invasive nor r | general cells | |
| JCRB0623 | HSC-3 | human | oral | squamous carcin | or squamous carcinoma, tongue, age 64, | general cells | |
| JCRB0624 | HSC-4 | human | oral | squamous carcin | or squamous carcinoma | general cells | |
| JCRB0625 | Ca9-22 | human | oral | gingival carcinoma | gingival carcinoma EGF receptor is pr | general cells | |
| JCRB0626 | TG1 | hamster, C | lung | V79 deliver | lung, V79 deliver | Use for metabolic c | general cells |
| JCRB0627 | KYM-1 | human | sarcoma | neck tumor, rhabd | neck tumor, rhabd | Plasminogen activ | general cells |
| JCRB0628 | MRK-nu-1 | human | mammary | mammary carcin | mammary carcin | his is one of the lir | general cells |
| JCRB0648 | L.P3 | mouse | skin | skin | This cell line was s Using serum free n | general cells | |
| JCRB0649 | HeLa.P3 | human | uterine cervix | | This cell line is supplied from Riken G | general cells | |
| JCRB0649.1 | HeLa.P3 | human | uterine cervix | | This cell line was d Definitely no serum : | general cells | |
| JCRB0650.1 | Mm2T12.P3 | Indian Mu | thymus | | The Mm2T12.P3 h | This culture has be | general cells |
| JCRB0701 | FM3A | mouse | mammary | mammary carcin | mammary carcin | High growth rate. F | general cells |
| JCRB0702 | FM3A HPRT~ | mouse | mammary | mammary carcin | mammary carcin | Good use for gene | general cells |
| JCRB0703 | FM3A APRT~ | mouse | mammary | mammary carcin | mammary carcin | 4-carbamoylimidaz | general cells |
| JCRB0704 | FM3A TK~ | mouse | mammary | mammary carcin | mammary carcinoma | general cells | |
| JCRB0705 | ts85 | mouse | mammary | | FM3A derivative,ut | Good to analyze pr | general cells |
| JCRB0706 | tsFT101 | mouse | mammary | FM3A derivative | Cytokinesis is bloc | Nucleus separation | general cells |
| JCRB0707 | V79 HPRT~ | hamster, C | lung | lung (M) | HPRT~ cells. | Mycoplasma was e | general cells |
| JCRB0708 | P3U1 | mouse | hemo-lym | myeloma | myeloma | general cells | |
| JCRB0709 | L5178Y | mouse | hemo-lym | leukemia | culturable derivative of L5178+ | general cells | |
| JCRB0709.1 | L5178Y TK+/-3.7.2c | mouse | hemo-lym | leukemia | TK+/- heterozygot | Incomplete descrip | general cells |
| JCRB0710 | EJ-1 | human | bladder | STR-PCR data ind | Ha-ras mutation re | Decontaminated by | general cells |
| JCRB0711 | T24 | human | bladder | urinary bladder car | urinary bladder car | Ha-ras mutation | general cells |
| JCRB0712 | MCA-TCL16 | mouse | embryo | C3H10T1/2 cl-8, tr: | C3H10T1/2 cl-8, tr: | Clone 16 of C3H10 | general cells |
| JCRB0713 | HeLa S3 (sc) | human | uterine cervix | , can grow in cervix, | can grow in medium + 0.1 % | pI | general cells |
| JCRB0714 | JTC-16 P3 | rat | liver | | ascites hepatoma | JTC16 is an in vitr | general cells |
| JCRB0714.1 | AH-7974.P3 | rat | liver | | Ascites hepatoma, | This cells were def | general cells |
| JCRB0718 | RLC-10.P3 | rat | liver | | liver, no serum required | general cells | |
| JCRB0718.1 | RLC-10.P3 | rat | liver | Telomelase was negative and a length of telomere was 2. | (general cells | | |
| JCRB0719 | FSthy21 | mouse | thymidylat | complete defective | complete defective Thymidylate synth | general cells | |
| JCRB0720 | OTT6050 | mouse | teratoma | teratocarcinoma | teratocarcinoma | general cells | |
| JCRB0721 | F9 | mouse | teratoma | teratocarcinoma | teratocarcinoma | Mycoplasma was e | general cells |
| JCRB0722 | L cl 1D | mouse | skin | TK~ | skin fibroblast, TK~,(M) | general cells | |
| JCRB0723 | L5178Y ts2 | mouse | hemo-lym | Abnormal Cell divis | Abnormal Cell division observed at noi | general cells | |

| | | | |
|-------------|-----------------|---|--|
| JCRB0724 | L5178Y ts39 | mouse | hemo-lymp A ts cell line. Perm Micronucleation observed when cells < general cells |
| JCRB0726 | A9(NT18)-2 | human chi human chi The cell line contai contains human chromosome #7 | general cells |
| JCRB0727 | A9(An4)-2 | human chi human chi contains human ch contains human chromosome #1 | general cells |
| JCRB0728 | A9(7151)-3 | human chi human chi contains human chromosome 17pter>17q11::Xq11>Xqter | general cells |
| JCRB0729 | A9(3884)-1 | human chi human chi contains human ch contains human chromosome 16pter> | general cells |
| JCRB0730 | A9(2621)-4 | human chi human chi contains human ch HAT selection required | general cells |
| JCRB0731 | A9(3552)-2 | human chi human chi contains human chromosome 11pter>11q23::Xq26>Xqter | general cells |
| JCRB0732 | A9(0439)-1 | human chi human chi contains human ch contains human chromosome Xpter>X | general cells |
| JCRB0733 | PC-12 | rat neural | Cells are differenti relating to research general cells |
| JCRB0734 | 3Y1-B Clone 1-6 | rat embryo | Ref(4007)showed i whole embryo |
| JCRB0735 | SV-3Y1-C66 | rat embryo | whole embryo, SV< SV40 transformed subline of 3Y1 |
| JCRB0736 | Py-3Y1-S2 | rat embryo | whole embryo, mot whole embryo, mouse polyoma virus-tr |
| JCRB0737 | Ad12-3Y1-Z19 | rat embryo | whole embryo, Ad1 whole embryo, Ad1 Prepare many amp |
| JCRB0738 | E1A-3Y1-1 | rat embryo | whole embryo, tran whole embryo, transformed 3Y1 by E1 |
| JCRB0739 | NG-3Y1-T15L | rat embryo | whole embryo, tran whole embryo, transformed 3Y1 by MN |
| JCRB0740 | NG-3Y1-T308 | rat embryo | whole embryo,tran whole embryo,transformed 3Y1 by MN |
| JCRB0741 | NG-3Y1-T6R | rat embryo | whole embryo, tran whole embryo, transformed 3Y1 by MN |
| JCRB0742 | SR-3Y1-2 | rat embryo | whole embryo,tran whole embryo,transformed 3Y1 by Rov |
| JCRB0743 | HR-3Y1-2 | rat embryo | whole embryo, tran whole embryo, transformed 3Y1 by v-f |
| JCRB0746 | YKG-1 | human neural | GFA and S-100 pr: human glioma cell line established fror |
| JCRB0801 | Caki-1 | human kidney | renal cancer renal cancer |
| JCRB0810 | ME-180 | human uterine cei uterus, cervical car uterus, cervical cancer | general cells |
| JCRB0813 | VMRC-RCW | human kidney | renal cell carcinoma renal cell carcinoma |
| JCRB0814 | VMRC-LCD | human lung (canc adenocarcinoma | adenocarcinoma |
| JCRB0815 | ABC-1 | human lung (canc adenocarcinoma | adenocarcinoma |
| JCRB0818 | SBC-3 | human lung (canc small cell carcinom small cell carcinoma (Oat cell type) | general cells |
| JCRB0819 | SBC-5 | human lung (canc small cell carcinom small cell carcinoma (intermediate) | general cells |
| JCRB0820 | EBC-1 | human lung (canc squamous cell car | squamous cell carcinoma |
| JCRB0821 | NUGC-2 | human stomach | gastric cancer |
| JCRB0822 | NUGC-3 | human stomach | gastric cancer |
| JCRB0823 | YMB-1 | human mammary | breast cancer |
| JCRB0824 | 8305C | human thyroid | C:G->T:A mutation thyroid anaplastic carcinoma, undiffere |
| JCRB0825 | YMB-1-E | human mammary | breast cancer |
| JCRB0826 | 8505C | human thyroid | C:G->G:C transversion at codon 248 of p53 gene |
| JCRB0827 | VMRC-RCZ | human kidney | renal cancer |
| JCRB0828 | HO-1-u-1 | human oral | squamous cell carcinoma |
| JCRB0829 | LK-2 | human lung (canc low molecular weight thiol protease inhibitor | general cells |
| JCRB0830 | PEER | human hemo-lymp | Negative for B cell acute lymphocytic I Cells originally con |
| JCRB0831 | HO-1-N-1 | human oral | squamous cell carcinoma |
| JCRB0832 | NH-6 | human neural | neuroblastoma cell line from primary tu |
| JCRB0833 | NH-12 | human neural | neroblastoma cell line |
| JCRB0834 | NUGC-4 | human stomach | signet ring cell carcinoma |
| JCRB1001.1 | MS-653-A | rat sarcoma | MS-653-A is highly tumorigenic clone f |
| JCRB1001.2 | MS-653-G | rat sarcoma | Rat fibrosarcoma cell line induced by n |
| JCRB1002 | K562/ADM | human hemo-lymp | chronic myelogeno Chronic myelogenc Mycoplasma decor |
| JCRB1003 | Kasumi-1 | human hemo-lymp | t(8;21), AML-ETO I MPO+, CD34+, CC The kasumi-1 is un |
| JCRB1004 | Kasumi-3 | human hemo-lymp | t(3;7) Overexpression of Evi-1 protein, CD34 |
| JCRB1005 | NCC-IT-A3 | human teratocarc | Cells are expressir teratocarcinoma fr Cells were obtaine |
| JCRB1006.0 | OUMS-36 | human embryo | normal cell normal fibroblast Normal human emt |
| JCRB1006.1 | OUMS-36T-1 | human embryo | 10^5 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.1F | OUMS-36T-1F | human embryo | P201 fibroblast, immortal |
| JCRB1006.2 | OUMS-36T-2 | human embryo | 10^7 cells/well wer normal human emt hTRT gene transfer |
| JCRB1006.2F | OUMS-36T-2F | human embryo | OUMS-36T series P200 fibroblast, immortal |
| JCRB1006.3 | OUMS-36T-3 | human embryo | 10^6 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.3F | OUMS-36T-3F | human embryo | P206 Human normal dipl |
| JCRB1006.4 | OUMS-36T-4 | human embryo | 10^6 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.4F | OUMS-36T-4F | human embryo | P202 Immortalized norm |
| JCRB1006.5 | OUMS-36T-5 | human embryo | 10^6 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.5F | OUMS-36T-5F | human embryo | P196 fibroblast, immortal |
| JCRB1006.6 | OUMS-36T-6 | human embryo | 10^6 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.6F | OUMS-36T-6F | human embryo | P210 fibroblast, immortal |
| JCRB1006.7 | OUMS-36T-7 | human embryo | 10^6 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.7F | OUMS-36T-7F | human embryo | Cells deposited at I Human normal fibr |
| JCRB1006.8 | OUMS-36T-8 | human embryo | 10^7 cells/well wer normal human emt hTRT gene transfe |
| JCRB1006.8F | OUMS-36T-8F | human embryo | OUMS-36T-8F/5F/ P203 fibroblast, immortal |
| JCRB1008 | HFb16d | human skin | Cells transfected b Normal fibroblast h Cells line is unique |
| JCRB1009 | RERF-GC-1B | human stomach | Gastric adenocarci No reference infor |
| JCRB1010 | KMRC-1 | human kidney | VHL gene alteratio Clear cell renal car P16 at deposit. |
| JCRB1011 | KMRC-2 | human kidney | VHL gene alteratio Clear cell renal car Cells were identifie |
| JCRB1012 | KMRC-3 | human kidney | VHL gene alteratio Clear cell renal carcinoma cell line |
| JCRB1014 | KMRM-M1 | human kidney | general cells |
| JCRB1015 | HSC-1 | human skin | excess expression Skin squamous car Mouse 3T3 irradiat |
| JCRB1016 | HSC-5 | human skin | Modal number of c The cells lost their ability to differentiate |
| JCRB1017 | HKA-1 | human skin | Chromosome is b Giant keratocan Mycoplasmas dec |
| JCRB1018 | HYM-1 | human melanoma | Modal number of c Melanoma cell line established from m |
| JCRB1019 | RERF-LC-Sq1 | human lung (cancer) | Established from human lung squamot |
| JCRB1020 | RERF-LC-Ad1 | human lung (cancer) | lung cancer (adeno carcinoma) cell lin |
| JCRB1021 | RERF-LC-Ad2 | human lung (cancer) | human lung cancer, adenocarcinoma c |
| JCRB1022 | OUMS-23 | human colon | Expression of the / Cells express heat Mycoplasmas dec |
| JCRB1024 | Kasumi-6 | human hemo-lymp | 45,XY,-9,add(12)(c Mononuclear cells Cell culture started |

| | | | | | | |
|------------|-------------------|------------|--------------------------------|--|--|----------------------|
| JCRB1025 | MKN7 | human | stomach | stomach, well differi stomach,well differ | 39 year-old, male. | general cells |
| JCRB1026 | OV3121 | mouse | ovary | Non-metastatic epithelial tumor cells fr | general cells | |
| JCRB1027 | SAT | human | oral | MMP and TIMP-1 Cells grow in monolayer with doubling | general cells | |
| JCRB1028 | JHH-2 | human | liver, gallb | Integration of the H hepatoma (Ed.II). + Cell Bank will hand | general cells | |
| JCRB1029 | JHH-5 | human | liver, gallb | Integration of the H Hepatocellular carc Cell Bank will hand | general cells | |
| JCRB1030 | JHH-6 | human | liver, gallb | HBV-DNA was not Undifferentiated m Cell Bank will hand | general cells | |
| JCRB1031 | JHH-7 | human | liver, gallb | HBV-DNA integration observed. 6.0/2, Cell Bank will hand | general cells | |
| JCRB1032 | OZ | human | liver, gallb | No. of chromosom Cells were establis Cell Bank will hand | general cells | |
| JCRB1033 | NOZ | human | liver, gallb | No. of chromosom Dt=48 hrs. Platting + Cell Bank will hand | general cells | |
| JCRB1034 | OUS-11 | human | lung | Cells are derived fr No references ava | general cells | |
| JCRB1039 | SKN-3 | human | oral | Cells were established through xenotra | general cells | |
| JCRB1040 | FF101 | rat | liver | Cells have charact The cells are exper | Cells have been m | general cells |
| JCRB1041 | SEKI | human | melanoma | Cells are consisted from both melanin | general cells | |
| JCRB1043 | OVISE | human | ovary | ER(-), PR(-) / tumo | Cells are transplantable on mouse hyp | general cells |
| JCRB1044 | OVKATE | human | ovary | ER(6.2 fmol/mg), P | Cells are transplan Td=18.3 hours | general cells |
| JCRB1045 | OVMANA | human | ovary | ER(-), PR(-) / tumo | Cells are transplant Td=26.1 hours | general cells |
| JCRB1046 | OVSAHO | human | ovary | PR(-) / tumour mar | Cells are transplan Td=18.3 hours | general cells |
| JCRB1047 | OVSYAO | human | ovary | ER(-), PR(-) / tumo | Cells were not tran The STR pattern of | general cells |
| JCRB1048 | OVTOKO | human | ovary | ER(-), PR(-) / tumo | Cells are transplant Td=26.1 hours | general cells |
| JCRB1049 | OVMIU | human | ovary | ER(-) / tumor mark | Cells are not trans The STR pattern of | general cells |
| JCRB1050 | OVMIU-II | human | ovary | tumour markers, C | Experiments on m The STR pattern w | general cells |
| JCRB1051 | OUMS-24/P61X | human | embryo | Human normal cell Life span of the OL | Cells were transfor | general cells |
| JCRB1052 | PA-1/6TG-r | human | teratocarc | 6TG resistant (HGI Teratocarcinoma cell line with 6TG-r, t) | general cells | |
| JCRB1053 | STC 1 | human | lung (cancer) | | general cells | |
| JCRB1054 | Hep G2 | human | liver, gallbladder | Hepatoma cell line. Mycoplasmas were | general cells | |
| JCRB1065 | Ki-JK | human | hemo-lymph | Positive markers, C | Mononuclear cells | |
| JCRB1066 | KMH-2 | human | thyroid | Chromosome No. 5 | doubling time is 58 hours and multidru | general cells |
| JCRB1067 | CPT-K5 | human | hemo-lymphocytic | Camptothecin resistant cell line. | Purific | general cells |
| JCRB1069 | NOMO-1/ADM | human | hemo-lymphocytic | Adriamycin-resistant derived from NO! | general cells | |
| JCRB1070 | HSGc-C5 | human | HeLa-contaminant | Radiation sensitiv | Although the STR-I | general cells |
| JCRB1071 | KMRC-20 | human | kidney | | general cells | |
| JCRB1073 | ASH-3 | human | thyroid | Chromosome numl | Doubling time is 84 Cells were identifie | general cells |
| JCRB1074 | FU97 | human | stomach | Td = 95 hrs. AFP production and tumo | general cells | |
| JCRB1075.0 | IM95 | human | stomach | Modal chromosom | Td=25 hrs. CEA related antigen obser | general cells |
| JCRB1075.1 | IM95m | human | stomach | Clone from the IM95. | A clone from the | general cells |
| JCRB1077 | DD-762 | mouse | mammary | MuMTV is positive. | DD-762 was isolated directly from DD- | general cells |
| JCRB1078 | TMH-1 | human | thyroid | Chromosome No. Doubling time is 74 Cross | contaminati | general cells |
| JCRB1079 | IHH-4 | human | thyroid | Doubling time = 42 Original tumor appeared in thyroid. | Ce | general cells |
| JCRB1080.0 | HARA | human | lung (cancer) | This cell line expre | STR analysis indic | general cells |
| JCRB1080.1 | HARA-B | human | lung (canc originally human ce | Delivertive of the HARA cell. This cell | I general cells | |
| JCRB1081 | MS-1-L | human | lung (canc | Chromosome distri | Cells are produc | Original name of th |
| JCRB1088 | RI-T | rat | liver | SV40 antigen posit | The cells were est | Early subculture is |
| JCRB1090 | NCE 16IIA | human | uterine cer | teromerase positiv HPV16 transformed noraml | human cel | general cells |
| JCRB1091 | NCE SVIIA3 | human | cervix | SV40 ori(-), terome | SV40 transformed human keratinocyte | general cells |
| JCRB1092 | NCE SVIA6 | human | cervix | teromerase positiv | SV40 transformed | Cell line was uniqu |
| JCRB1093 | PSVK1 | human | foreskin | Cells transfected b | Human primary cell line immortalized | b general cells |
| JCRB1094 | SUIT-2 | human | pancreas | Mode of chromoso | PDL=38.2 hrs | general cells |
| JCRB1096 | hs-103-3 | hamster, C | lung | Temperature sensi | anchorage independent cell line | transf general cells |
| JCRB1097 | hs-164-2 | hamster, C | lung | Temperature sensi | Anchorage independent cell line | transf general cells |
| JCRB1098 | hs-171-1 | hamster, C | lung | Temperature sensitive mutant, which are transformed | bele | general cells |
| JCRB1099 | hs-172-3 | hamster, C | lung | Temperature sensitive mutant, which are transformed | bele | general cells |
| JCRB1100 | hs-211 | hamster, C | lung | Temperature sensi | Anchorage independent cell line | transf general cells |
| JCRB1101 | hs-222-3 | hamster, C | lung | temperature sensit | Anchorage independent cell line | transf general cells |
| JCRB1102 | cs-17-25 | hamster, C | lung | Temperature sensi | Anchorage independent cell line | transf general cells |
| JCRB1103 | cs-19-36 | hamster, C | lung | Temperature sensi | Anchorage independent cell line | transf general cells |
| JCRB1104 | cs-20-5 | hamster, C | lung | Temperature sensi | Anchorage independent cell line | transf general cells |
| JCRB1105 | AIG | hamster, C | lung | Anchorage independent cell line | transf | general cells |
| JCRB1107 | UCBTERT-21 | human | mesenchy | Human umbilical cx ** | | general cells |
| JCRB1108 | UCB408E6E7-31 | human | mesenchy | Human umbilical cx ** | | general cells |
| JCRB1109 | UCB408E7-32 | human | mesenchy | Human umbilical cx ** | | general cells |
| JCRB1110 | UCB408E6E7TERT-33 | human | mesenchy | Human umbilical cx ** | | general cells |
| JCRB1117 | HEC-1-A | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1118 | HEC-6 | human | uterus | endometrial adeno | endometrial adenocarcinoma | general cells |
| JCRB1119 | KUSA-A1 | mouse | bone marrow | | | general cells |
| JCRB1120 | HEC-59 | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1121 | HEC-88nu | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1122 | HEC-151 | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1123 | HEC-108 | human | uterus | | endometrial adenocarcinoma | general cells |
| JCRB1124 | HEC-116 | human | uterus | | | general cells |
| JCRB1129 | KUSA-H1 | mouse | bone marrow | | | general cells |
| JCRB1131 | UE6E7T-1 | human | bone marr | Human bone marrow-derived mesenchymal stem cells | wh | general cells |
| JCRB1132 | KUSA-0 | mouse | bone marrow | | | general cells |
| JCRB1133 | UE6E7T-2 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells | wh | general cells |
| JCRB1134 | KUM3 | mouse | bone marrow | | | general cells |
| JCRB1135 | KUM4 | mouse | bone marrow | | | general cells |
| JCRB1136 | UE6E7T-3 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells | wh | general cells |
| JCRB1137 | UE6E7TC-4 | human | bone marr | Human bone marrow-derived mesenchymal stem cells | wh | general cells |
| JCRB1138 | NRG | mouse | bone marrow | | | general cells |
| JCRB1139 | KUM9 | mouse | bone marrow | | | general cells |
| JCRB1140 | UBE6T-6 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells | wh | general cells |

| | | | | | | |
|-------------|-----------------|----------|-----------------------------|--|--|---------------------|
| JCRB1141 | HEC-251 | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1142 | HEC-265 | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1143 | UBE6T-7 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1145 | HEC-50B | human | uterus | endometrial carcin | endometrial carcinoma | general cells |
| JCRB1146 | RSMG-1 | rat | submandibular gland | Growth of the RSM Co-operative rese | general cells | |
| JCRB1147 | UE7T-9 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1148 | KMBC-2 | human | bladder | ***** | Doubling time for about 60 hours. A ro | general cells |
| JCRB1149 | UE6E7T-11 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1150 | KMPC-3 | human | kidney | ***** | Doubling time for about 40 hours. The | general cells |
| JCRB1151 | UE6E7T-12 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1152 | RSMG-2 | rat | submandibular gland | Growth of the RSM Co-operative rese | general cells | |
| JCRB1153 | HCA-1 | human | uterine cei | endocervical aden | endocervical adenocarcinoma | general cells |
| JCRB1154 | UE7T-13 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1159 | UBE6T-15 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1160 | UE6E7-16 | human | mesenchy | Human bone marrow-derived mesenchymal stem cells wh | general cells | |
| JCRB1166 | NCR-G1 | human | embryoni | c testis | Complex type germ cell tumor of human embryonic testis | c general cells |
| JCRB1167 | NCR-G2 | human | embryoni | c testis | Complex type germ cell tumor of human embryonic testis | c general cells |
| JCRB1168 | NCR-G3 | human | embryoni | c testis | Complex type germ cell tumor of human embryonic testis | c general cells |
| JCRB1170 | SKNO-1 | human | hemo-lym | t(8;21), monosomy MPO+, CAE+, NBE- | | general cells |
| JCRB1175 | NCR-G4 | human | embryoni | c testis | Complex type germ cell tumor of human embryonic testis | c general cells |
| JCRB1176 | KUM10 | mouse | bone marrow | | | general cells |
| JCRB1178 | KUM5 | mouse | bone marrow | | | general cells |
| JCRB1179 | KMS-11 | human | hemo-lym | t(4;14)(p16.3;q32.3 small round-cell morphology, part is mi | general cells | |
| JCRB1180 | KMM-1 | human | hemo-lym | t(8;14)(q24;q32), t(small round-cell morphology, part is mi | general cells | |
| JCRB1181 | 9-15c | mouse | bone marrow | | | general cells |
| JCRB1182 | NCC-RbC-51 | human | neural | familial bilateral retinoblastoma | | general cells |
| JCRB1183 | NCC-RbC-54 | human | neural | Bilateral retinoblastoma | | general cells |
| JCRB1185 | KMS-21BM | human | hemo-lym | t(11;14)(q13;q32) small round-cell morphology, part is mi | general cells | |
| JCRB1186 | NCC-RbC-59 | human | retina | Bilateral retinoblastoma | | general cells |
| JCRB1187 | KMS-26 | human | hemo-lym | t(4;14)(p16.3;32.3) small round-cell morphology, part is mi | general cells | |
| JCRB1188 | KMS-27 | human | hemo-lym | t(11;14)(q16.3;q32 small round-cell morphology, part is mi | general cells | |
| JCRB1189 | F56 | mouse | antibody-p | Monoclonal antibody Producing IgG1 | | general cells |
| JCRB1190 | NCC-RbC-57 | human | neural | Unilateral retinoblastoma | | general cells |
| JCRB1191 | KMS-28PE | human | hemo-lym | t(4;14)(p16.3;q32.3 small round-cell morphology, part is mi | general cells | |
| JCRB1192 | KMS-28BM | human | hemo-lym | t(4;14)(p16.3;q32.3 small round-cell morphology, part is mi | general cells | |
| JCRB1193 | HEC-1-B | human | uterus | endometrial carcin | endometrial carcinoma HEC-1-B is the sar | general cells |
| JCRB1195 | KMS-34 | human | hemo-lym | t(4;14)(p16.3;32.3) small round-cell morphology, part is mi | general cells | |
| JCRB1196 | KMS-20 | human | hemo-lymphocytic | | small round-cell morphology, part is mi | general cells |
| JCRB1198.01 | GP8 | mouse | skin | cell line having the human beta-hgalactosidase gene | | general cells |
| JCRB1199 | R201C | mouse | skin | | | general cells |
| JCRB1200 | NCC-RbC-53 | human | retina | Unilateral retinoblastoma | | general cells |
| JCRB1201 | GCH-nu-YS | human | uterus | choriocarcinoma | hCG production | general cells |
| JCRB1202 | KUM6 | mouse | bone marrow | | | general cells |
| JCRB1203 | HCS-2 | human | uterine cei | squamous cell car | squamous cell carcinoma | general cells |
| JCRB1204 | KUM7 | mouse | bone marrow | | | general cells |
| JCRB1205 | HCSC-1 | human | uterine cei | small cell carcinom | small cell carcinoma | general cells |
| JCRB1206 | TK | human | hemo-lym | 45,X,-Y,del(2)(q21 TK cells grow in su Both NIH | NIHS0245(TK | general cells |
| JCRB1207 | SV | mouse | skin | beta-hgal gene deficient | | general cells |
| JCRB1208 | HM-1 | hamster, | hemo-lym | hamster cell | macrophage | Cell growth is slow |
| JCRB1212 | NCC-RbC-56 | human | neural | Bilateral retinoblastoma | | general cells |
| JCRB1213.2 | DIG222.4D.5 | mouse | antibody-producing hybridom | Producing anti Digoxin IgG2b (k) isotyp | general cells | |
| JCRB1213.3 | DIG104.10H.1 | mouse | antibody-producing hybridom | Producing anti Digoxin IgG1 (k) isotyp | general cells | |
| JCRB1214 | NCC-RbC-67 | human | neural | Unilateral retinoblastoma | | general cells |
| JCRB1221 | NCC-RbC-92 | human | neural | Unilateral retinoblastoma | | general cells |
| JCRB1222.01 | HGH1.1.3 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.02 | HGH1.6.5 | mouse | antibody-producing hybridom | Producing IgG gamma2a, and k. | | general cells |
| JCRB1222.03 | HGH1.17.2 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.04 | HGH1.35.2 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.05 | HGH2.6.1 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.06 | HGH1.9.3 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.07 | HGH1.8.1 | mouse | antibody-producing hybridom | Producing IgG gamma2b and k | | general cells |
| JCRB1222.08 | HGH1.11.1 | mouse | antibody-producing hybridom | Producing IgG gamma2b and k | | general cells |
| JCRB1222.09 | HGH1.15.3 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.10 | HGH1.41.1 | mouse | antibody-producing hybridom | Producing IgG gamma2b and k | | general cells |
| JCRB1222.11 | HGH1.46.1 | mouse | antibody-producing hybridom | Producing IgG gamma2b and k | | general cells |
| JCRB1222.12 | HGH1.31.3 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.13 | HGH1.32.4 | mouse | antibody-producing hybridom | Producing IgG gamma2a and k | | general cells |
| JCRB1222.14 | HGH1.37.3 | mouse | x n hybridoma | Producing IgG gamma 2a and k | | general cells |
| JCRB1222.15 | HGH2.4.2 | mouse | antibody-producing hybridom | Producing IgG gamma1 and k | | general cells |
| JCRB1222.16 | HGH2.15.4 | mouse | antibody-producing hybridom | Producing IgG gamma1 and lambda | | general cells |
| JCRB1223 | NCC-RbC-T1 | human | neural | Bilateral retinoblastoma | | general cells |
| JCRB1225 | I51T | mouse | skin | | Cells were originally | general cells |
| JCRB1226 | GHBP116 | mouse | antibody-producing hybridom | Producing IgG1(k) | | general cells |
| JCRB1228 | NCC-StC-K140 | human | stomach | | | general cells |
| JCRB1301 | SC78.H81.C81.A9 | mouse | antibody-producing hybridom | Producing IgG1(k) | | general cells |
| JCRB1303 | NCC-RbC-83 | human | retina | Unilateral retinoblastoma | | general cells |
| JCRB1304 | MT-45-5-3 | mouse | antibody-producing hybridom | Producing IgG1(K) | | general cells |

| | | | | | |
|-------------|----------------|------------|------------------------------|--|------------------------------------|
| JCRB1305 | NCC-RbC-39 | human | neural | Unilateral retinoblastoma | general cells |
| JCRB1308 | NCC-OvC-K119 | human | ovary | Ovarian carcinoma | general cells |
| JCRB1311 | ZEN | mouse | antibody-producing hybridoma | Producing IgG1 lambda. | general cells |
| JCRB1314 | HPSV178-4-3 | mouse | antibody-p | Producing IgG2b(k) producing mAb HPSV178, against the general cells | general cells |
| JCRB1316.1 | OHP57.G6.1 | mouse | antibody-producing hybridoma | Producing IgG k. | general cells |
| JCRB1316.2 | OHP101.B11.1 | mouse | antibody-producing hybridoma | Producing IgG k. | general cells |
| JCRB1316.3 | OHP138.H8.1 | mouse | antibody-producing hybridoma | Producing IgG k. | general cells |
| JCRB1316.4 | OHP1E9.3.1 | mouse | antibody-producing hybridoma | Producing IgG k. | general cells |
| JCRB1316.5 | OHP4B2.2.3 | mouse x n | hybridoma | Producing IgG k. | general cells |
| JCRB1316.6 | OHP7D7.2.3 | mouse | antibody-producing hybridoma | Producing IgG k. | general cells |
| JCRB1317 | KCL-22 | human | hem-o-lymp | Ph1-positive. | general cells |
| JCRB1318 | HeLa9903 | human | uterine cervix | HeLa clone. | genetically-modified cells |
| JCRB1319 | PL-21 | human | hem-o-lymp | 13q+, A marker chi immature cells with azurophilic granule | general cells |
| JCRB1320 | NCC-CoC-K115P | human | colon | colon carcinoma | general cells |
| JCRB1322 | mSS | mouse | skin | Near-diploid chori normal | general cells |
| JCRB1323 | F25 | mouse | antibody-producing hybridoma | Producing IgG2a | general cells |
| JCRB1324 | F119 | mouse | antibody-producing hybridoma | Producing IgG2b | general cells |
| JCRB1325 | MKPL-1 | human | hem-o-lymp | 47,X,-Y,+21,+3mar. Dt = 30 hr. Large primitive blasts with ζ | general cells |
| JCRB1326 | NCC-RbC-60 | human | retina | Bilateral retinoblastoma | general cells |
| JCRB1328 | AR-EcoScreen | hamster, C | ovary | This cell is that AR that stably express human androgen receptor. | genetically-modified cells |
| JCRB1330 | Ty-82 | human | thymus | 46XX, t(15;19)(q15;15) Thymic carcinoma. | general cells |
| JCRB1332 | NALL-1 | human | hem-o-lymp | EB virus was not d The cell line is negative for E-rosettes | general cells |
| JCRB1333 | NCC-CoC-K115L | human | colon | Metastatic colon carcinoma | general cells |
| JCRB1334 | RC-K8 | human | hem-o-lymp | 46XY, with t(11;14) Dt=48-60hr. The cell line produces pla | general cells |
| JCRB1335 | HD-70 | human | hem-o-lymp | CD30+, CD15+, C γ Dt=28 hr. | general cells |
| JCRB1337 | DL-40 | human | hem-o-lymp | Cells are positive f Cells grow without forming clumps with | general cells |
| JCRB1339 | NCC-CoC-K115B | human | colon | B-cell lymphocyte transformed by EBV. | general cells |
| JCRB1340 | RS-2 | human | brain | | general cells |
| JCRB1342.01 | MOR8.2.1 | mouse | antibody-producing hybridoma | Producing IgG1k. | general cells |
| JCRB1342.02 | MOR35.4.12 | mouse | antibody-producing hybridoma | Producing IgG1k. | general cells |
| JCRB1342.03 | MOR39.3.9 | mouse x n | hybridoma | Producing IgG1lambda. | general cells |
| JCRB1342.04 | MOR76.7F.16 | mouse x n | hybridoma | Producing IgG1 k. | general cells |
| JCRB1342.05 | MOR83.3.10 | mouse | antibody-producing hybridoma | Producing IgG1 k. | general cells |
| JCRB1342.06 | MOR115.1.13 | mouse x n | hybridoma | Producing IgG1 lambda. | general cells |
| JCRB1342.07 | MOR124.2.2 | mouse x n | hybridoma | Producing IgG1 k. | general cells |
| JCRB1342.08 | MOR131.5.13 | mouse x n | hybridoma | Producing IgG1 lambda. | general cells |
| JCRB1342.09 | MOR158.1.3 | mouse x n | hybridoma | Producing IgG1 k. | general cells |
| JCRB1342.10 | MOR33.2.9 | mouse x n | hybridoma | Producing IgG1 k. | general cells |
| JCRB1342.11 | MOR180.2.4 | mouse x n | hybridoma | Producing IgG1 lambda. | general cells |
| JCRB1342.12 | MOR44.4.1 | mouse x n | hybridoma | Producing IgG1 k. | general cells |
| JCRB1343 | OKa-C-1 | human | lung | (canc Chromosome num) The cell line was established from the | general cells |
| JCRB1344 | delta-47 | human | hem-o-lymp | The cell line expres Cells are secreting IgD. | general cells |
| JCRB1345 | PALL-2 | human | hem-o-lymp | Single Ph1 chromosome observed. | general cells |
| JCRB1346 | FLAM-76 | human | hem-o-lymp | Hypodiploid chrom kappa IgG (nonsecretory), CD38, CD4 | general cells |
| JCRB1347 | KML-1 | human | hem-o-lymp | Non-Hodgkin's lymphoma of large-cell type. Loss of chro | general cells |
| JCRB1348 | 3LL | mouse | lung | Cell line having high metastasis poten | general cells |
| JCRB1349 | Ex-3LL | mouse | lung | Lewis lung carcinoma derived cell line | general cells |
| JCRB1350 | EL-4 | mouse | lymphoma | | general cells |
| JCRB1351 | Int-3LL | mouse | lung | Lewis lung carcinoma derived cell line | general cells |
| JCRB1352 | EL-4ad | mouse | lymphoma | adhered to a tissue culture dish | general cells |
| JCRB1353 | TER-119 | mouse x r | hybridoma | Producing rat IgG 2b and kappa | general cells |
| JCRB1354 | HSC-3-M3 | human | tongue | squamous carcinori Highly metastatic cell line. | general cells |
| JCRB1355 | 1-1ras1000 | mouse | embryo | Fibroblast, subclone of Balb/3T3-A31. | general cells |
| JCRB1356 | A31-1-1 | mouse | embryo | whole embryo fibroblast, subclone The FCS requires | general cells |
| JCRB1357 | 1-src | mouse | embryo | Fibroblast, subclone of Balb/3T3-A31. | general cells |
| JCRB1358.01 | SRBC-P13-7-3 | mouse x n | hybridoma | Producing IgM(k) Recloned from P13- | general cells |
| JCRB1358.02 | SRBC-P18-3-8 | mouse x n | hybridoma | Producing IgM(k). Recloned from P18- | general cells |
| JCRB1358.03 | SRBC-P20-13-13 | mouse x n | hybridoma | Producing IgM(k). Recloned from P20- | general cells |
| JCRB1358.04 | SRBC-F12-1-19 | mouse x n | hybridoma | Producing IgM(k). Recloned from F12- | general cells |
| JCRB1358.05 | SRBC-F12-18-7 | mouse x n | hybridoma | Producing IgM(k). Recloned from F12- | general cells |
| JCRB1358.06 | SRBC-F23-20-22 | mouse x n | hybridoma | Producing IgM(k). Recloned from F23- | general cells |
| JCRB1358.07 | SRBC-F31-2-11 | mouse x n | hybridoma | Producing IgM(k). Recloned from F31- | general cells |
| JCRB1358.08 | SRBC-F31-19-1 | mouse x n | hybridoma | Producing IgM(k). Recloned from F31- | general cells |
| JCRB1358.09 | SRBC-RP8-2-4 | mouse x n | hybridoma | Producing IgG1(k). Recloned from RP1 | general cells |
| JCRB1358.10 | SRBC-RP8-18-15 | mouse x n | hybridoma | Producing IgG1(k). Recloned from RP1 | general cells |
| JCRB1358.11 | SRBC-RF29-1-3 | mouse x n | hybridoma | Producing IgG1(k). Recloned from RF1 | general cells |
| JCRB1359 | HAC-2 | human | ovary | Ovarian clear cell Clear cell carcinoma. Td = 46.8 hrs, C | general cells |
| JCRB1361 | Vn1919 | mouse | brain | GFAP (glial fibrillary acidic protein) pos | general cells |
| JCRB1362 | HCT-2 | hamster, C | hem-o-lymp | HTLV-1 positive b The cells are forming loosely aggregat | general cells |
| JCRB1365 | Vn1919 cl4 | mouse | brain | GFAP (glial fibrillary acidic protein) pos | general cells |
| JCRB1367 | Vn1919 cl49 | mouse | brain | GFAP (glial fibrillary acidic protein) pos | general cells |
| JCRB1368 | Vn1919 cl59 | mouse | brain | GFAP (glial fibrillary acidic protein) pos | general cells |
| JCRB1370 | TAS-5 | mouse x r | hybridoma | Producing rat IgG 2b and kappa | general cells |
| JCRB1371 | GS-3A4-HepG2 | human | liver | human glutamine synthetase and CYP3A4 introduced | genetically-modified cells |
| JCRB1372 | MCF-7-Luc | human | breast | pMSCV-luc transfected MCF-7 cells. Luciferase stably ex | luciferase-expressing cancer cells |
| JCRB1373 | BT-549-Luc | human | breast | pMSCV-luc transfected BT-549 cells. Luciferase stably ex | luciferase-expressing cancer cells |
| JCRB1375 | Hs578T-Luc | human | breast | pMSCV-luc transfected Hs578T cells. Luciferase stably ex | luciferase-expressing cancer cells |
| JCRB1376 | SK-MEL-28-Luc | human | melanoma | pMSCV-luc transfected SK-MEL-28 cells. Luciferase stabl | luciferase-expressing cancer cells |
| JCRB1377 | COLO 679-Luc | human | melanoma | pMSCV-luc transfected COLO 679 cells. Luciferase stably luciferase-expressing cancer cells | |
| JCRB1378 | MM-RU-Luc | human | melanoma | pMSCV-luc transfected MM-RU cells. Luciferase stably ex | luciferase-expressing cancer cells |

| | | | | | | |
|----------|---------------------|-------|----------------|---|------------------------------------|---|
| JCRB1379 | MKN45-Luc | human | stomach | pMSCV-luc transfected MKN45 cells. Luciferase stably ex luciferase-expressing cancer cells | | |
| JCRB1380 | HBs320 | mouse | x n hybridoma | hybridoma producing monoclonal antib general cells | | |
| JCRB1381 | Hbs128 | mouse | x n hybridoma | hybridoma producing monoclonal antib general cells | | |
| JCRB1382 | DLD-1 clone#1-Luc | human | colon | pMSCV-luc transfected DLD-1 clone#1 cells. Luciferase s luciferase-expressing cancer cells | | |
| JCRB1383 | HT-29-Luc | human | colon | pMSCV-luc transfected HT-29 cells. Luciferase stably exp luciferase-expressing cancer cells | | |
| JCRB1384 | STR-428 | human | ascites | IgH, c-Myc, BCL2 : Immunophenotypic Contact to Dr. Tair: general cells | | |
| JCRB1385 | CAS-9 | mouse | x r; hybridoma | Producing rat IgG 2b and kappa | general cells | |
| JCRB1386 | U-251 MG-Luc | human | glioblastor | pMSCV-luc transfe astrocytoma, GFAP-positive | luciferase-expressing cancer cells | |
| JCRB1387 | DU-145-Luc | human | prostate | pMSCV-luc transfected DU-145 cells. Luciferase stably ex luciferase-expressing cancer cells | | |
| JCRB1388 | A-498-Luc | human | renal | pMSCV-luc transfected A-498 cells. Luciferase stably exp luciferase-expressing cancer cells | | |
| JCRB1389 | KM12-Luc | human | colon | pMSCV-luc transfected KM12 cells. Luciferase stably expi luciferase-expressing cancer cells | | |
| JCRB1390 | G-401-Luc | human | renal | pMSCV-luc transfected G-401 cells. Luciferase stably exp luciferase-expressing cancer cells | | |
| JCRB1391 | KMS-33 | human | melanoma | small round-cell morphology, part is m general cells | | |
| JCRB1392 | AS17 | mouse | x n hybridoma | Producing mouse IgG 2b and kappa | general cells | |
| JCRB1393 | SK-MEL-2-Luc | human | melanoma | pMSCV-luc transfected SK-MEL-2 cells. Luciferase stably luciferase-expressing cancer cells | | |
| JCRB1394 | KHM-2B | human | leukemia | t(8;14)(q24;q32) ar HLA-DR, CD10, CD20, and CD19 wer general cells | | |
| JCRB1395 | Kasumi-2 | human | leukemia | t(1;19)(q23;p13), E2A-PBX fusion gene | general cells | |
| JCRB1396 | Mewo-Luc | human | melanoma | pMSCV-luc transfected Mewo cells. Luciferase stably exp luciferase-expressing cancer cells | | |
| JCRB1397 | 786-O-Luc | human | renal | pMSCV-luc transfected 786-O cells. Luciferase stably expi luciferase-expressing cancer cells | | |
| JCRB1398 | Kasumi-5 | human | leukemia | t(4;11)(q21;p15), R CD7+, CD10+, CD56+, CD38+, CD3-, general cells | | |
| JCRB1399 | SW620-Luc | human | colon | pMSCV-luc transfected SW620 cells. Luciferase stably ex luciferase-expressing cancer cells | | |
| JCRB1400 | Caki-1-Luc | human | renal | pMSCV-luc transfected Caki-1 cells. Luciferase stably exp luciferase-expressing cancer cells | | |
| JCRB1401 | Kasumi-7 | human | leukemia | del (9p), [der(9)(q, CD10+, CD19+, CD20+, HLA-DR+, CI general cells | | |
| JCRB1402 | K562-Luc | human | leukemia | pMSCV-luc transfected K562 cells. Luciferase stably expr luciferase-expressing cancer cells | | |
| JCRB1403 | Kasumi-8 | human | leukemia | t(9;22)(q32;q11), b Ph1 positive | general cells | |
| JCRB1404 | Minami-1 | human | lymphoma | t(14;18)(q32;q21), bcl gene rearrangement | general cells | |
| JCRB1405 | Minami-2 | human | lymphoma | del(13q), add(18)(c CD10+, CD19+, CD20+, CD22+, HLA-I general cells | | |
| JCRB1406 | PC-3-Luc | human | prostate | pMSCV-luc transfected PC-3 cells. Luciferase stably expr luciferase-expressing cancer cells | | |
| JCRB1407 | NCI-H460-Luc | human | lung | pMSCV-luc transfected NCI-H460 cells. Luciferase stably luciferase-expressing cancer cells | | |
| JCRB1408 | HCT-116 clone#2-Luc | human | colon | pMSCV-luc transfected HCT-116 clone#2 cells. Luciferast luciferase-expressing cancer cells | | |
| JCRB1409 | Kasumi-9 | human | leukemia | normal karyotype (no EBV) | general cells | |
| JCRB1410 | Kasumi-10 | human | leukemia | t(11;19)(q23;p13.3), MLL gene rearrangement | general cells | |
| JCRB1411 | KasumiA-541 | human | leukemia | nomal karyotype (n CD34+, CD33+, CD10+, CD19+, HLA- general cells | | |
| JCRB1413 | ST 13 | mouse | mammary | differential to adipocyte | general cells | |
| JCRB1414 | A549-Luc | human | lung | pMSCV-luc transfected A549 cells. Luciferase stably expr luciferase-expressing cancer cells | | |
| JCRB1415 | KasumiA-568 | human | leukemia | 45, XY, del(2)(q7), CD13+, CD33+, c-kit+, GPA+ | general cells | |
| JCRB1416 | KasumiA-554 | human | leukemia | t(11;19)(q23;p13.1 CD33+, CD15+, CD4+, CD56+, HLA-C general cells | | |
| JCRB1433 | MKN-1/CMV-Luc | human | stomach | Luciferase stably expressing cell line (MKN1; Human storr luciferase-expressing cancer cells | | |
| JCRB1435 | TK6(IVGT) | human | spleen | This line is derivative of the WIL-2 cell Same cell line in E1 general cells | | Only for researchers in Asian countries |
| JCRB1438 | BT-20/CMV-Luc | human | mammary | Luciferase stably expressing cell line (BT-20; Human brea luciferase-expressing cancer cells | | |
| JCRB1447 | 4T1-Luc | mouse | mammary | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1448 | BxPC-3-Luc#2 | human | pancrea | pMSCV-luc transfe Cellular Products: mucin: pancreas ca luciferase-expressing cancer cells | | |
| JCRB1450 | BT-474/CMV-Luc | human | mammary | Luciferase stably expressing cell line (BT-474; Human bre luciferase-expressing cancer cells | | |
| JCRB1451 | COLO 205/CMV-Luc | human | colon | pLL3.7-luc transfet Cellular Products: (The line was deriv: luciferase-expressing cancer cells | | |
| JCRB1453 | CLAC | dog | lung | This cell line including cancer stem cel general cells | | |
| JCRB1454 | AsPC-1/CMV-Luc | human | pancreas | Luciferase stably expressing cell line (AsPC-1; human par luciferase-expressing cancer cells | | |
| JCRB1458 | HUEhT-1 | human | umbilical v | HUEhT-1 was esta Cell culture in collagen I coated dish. | genetically-modified cells | |
| JCRB1459 | HUEhT-2 | human | umbilical v | HUEhT-2 was esta Cell culture in collagen I coated dish. | genetically-modified cells | |
| JCRB1466 | HCC-1419-Luc | human | breast | pMSCV-luc transfe The cells are poorly differentiated. The luciferase-expressing cancer cells | | |
| JCRB1471 | MKN-7/CMV-Luc | human | stomach | pLL3.7-luc transfected. Luciferase stably expressing cell li luciferase-expressing cancer cells | | |
| JCRB1473 | MKN-74/CMV-Luc | human | stomach | Luciferase stably expressing cell line (MKN74; Human mo luciferase-expressing cancer cells | | |
| JCRB1474 | B16-F0-Luc | mouse | melanoma | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1475 | Nalm-6-MSH+ | human | leukemia | MSH+, POL(WT), TK+/- | general cells | |
| JCRB1476 | HCC-1954-Luc | human | breast | pMSCV-luc transfe HCC1954 is positive for the epithelial c luciferase-expressing cancer cells | | |
| JCRB1478 | NCI-H1650-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1479 | HCT-15-Luc#1 | human | colon | pMSCV-luc transfe Cellular Products: carcinembryonic ai luciferase-expressing cancer cells | | |
| JCRB1484 | JMU-RTK-2 | human | kidney | highly aggressive malignancy | general cells | |
| JCRB1485 | IM95/CMV-Luc | human | stomach | Luciferase stably expressing cell line (IM95; Human gastri luciferase-expressing cancer cells | | |
| JCRB1486 | NCI-H1975-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1487 | KATOII/CMV-Luc | human | stomach | tetraploid, pLL3.7-1 Derived from stomach cancer, signet r luciferase-expressing cancer cells | | |
| JCRB1489 | POLZ D2781N | human | peripheral | MSH+, POLZ D27: Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1490 | POLZ L2618M | human | peripheral | MSH+, POLZ L261 Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1491 | KP-2/CMV-Luc | human | pancrea | pLL3.7-luc transfet The tumor cell line is derived from pan luciferase-expressing cancer cells | | |
| JCRB1492 | NCI-H2228-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1494 | POLZ KO MSH- | human | peripheral | MSH-, POLZ KO/K Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1496 | Colon26-Luc | mouse | colorectal | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1497 | NCI-H441/CMV-Luc | human | lung | pLL3.7-luc transfected. Luciferase stably expressing cell li luciferase-expressing cancer cells | | |
| JCRB1498 | POLZ CD MSH- | human | peripheral | MSH-, POLZ catal Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1502 | MDA-MB-361-Luc#1 | human | breast | Luciferase stably expressing cell line (MDA-MB-361; Hum luciferase-expressing cancer cells | | |
| JCRB1503 | NIH3T3/ATCC/CMV-Luc | mouse | embryo | Luciferase stably expressing cell line (NIH3T3; Mouse em luciferase-expressing cancer cells | | |
| JCRB1504 | hTf2-9F | mouse | x n hybridoma | This cell line produ This cell line produ general cells | | |
| JCRB1505 | Ishikawa 3-H-12 | human | Endometrium | | general cells | |
| JCRB1506 | POLK KO-comp | human | peripheral | MSH+, POLK KO- Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1508 | HCC-1937/CMV-Luc | human | breast | pLL3.7-luc transfec The cells are negative for expression c luciferase-expressing cancer cells | | |
| JCRB1509 | NCI-H23-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1510 | POLZ Y2779F | human | peripheral | MSH+, POLZ Y277 Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1514 | POLK KO homo MSH- | human | leukemia | MSH-, POLK KO/KO, TK+/- | genetically-modified cells | |
| JCRB1515 | POLZ L2618F | human | peripheral | MSH+, POLZ L261 Human cell line derived from B cell leu genetically-modified cells | | |
| JCRB1516 | HCC-827-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | | |
| JCRB1517 | POLK F171A MSH- | human | leukemia | MSH-, POLK KO/F171A, TK+/- | genetically-modified cells | |

| | | | | | |
|-------------|------------------------|------------|---------------|---|--|
| JCRB1518 | POLK KO hetero MSH- | human | leukemia | MSH-, POLK KO/+, TK+/+ | genetically-modified cells |
| JCRB1523 | POLK CD MSH- | human | peripheral | MSH-, POLK catal Human cell line derived from B cell leu genetically-modified cells | |
| JCRB1524 | POLK CD | human | peripheral | MSH-, POLK catal Human cell line derived from B cell leu genetically-modified cells | |
| JCRB1525 | NCI-H650-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | |
| JCRB1526 | POLK KO-CD comp | human | peripheral | MSH+, POLK KO-: Human cell line derived from B cell leu genetically-modified cells | |
| JCRB1529 | NCI-H2009-Luc | human | lung | pMSCV-luc transfected. Luciferase stably expressing cell luciferase-expressing cancer cells | |
| JCRB1531 | MEFs | Mouse | embryo | Mouse embryonic fibroblasts (control I general cells | |
| JCRB1533 | 293 TRE/Cont | human | kidney | G418 resistance | general cells |
| JCRB1534 | IKK-i-DEF | mouse | embryo | Mouse embryonic fibroblasts prepared from IkB kinase (IK general cells | |
| JCRB1535 | TBK1-DEF | mouse | embryo | Mouse embryonic fibroblasts prepared from TANK-binding genetically-modified cells | |
| JCRB1536 | EPC-1 | human | uterine | Human uterine endometrium vascular endothelial progenit genetically-modified cells | |
| JCRB1537 | LoVo-Luc#2 | human | colorectal | pMSCV-luc transfe Carcinoembryonic antigen-producing luciferase-expressing cancer cells | |
| JCRB1538 | EMC100 | human | uterine | Human uterine endometrium vascular endothelial progenit genetically-modified cells | |
| JCRB1539 | HAC-2 | human | ascites | Human ovary cancer cell line derived from mesonephros. general cells | |
| JCRB1542 | KP-3L-Luc#5 | human | pancreas | pMSCV-luc transfe Tissue was classified as a metastasis luciferase-expressing cancer cells | |
| JCRB1544 | EMC214 | human | uterine | Human uterine endometrium vascular endothelial progenit genetically-modified cells | |
| JCRB1545 | EM-E6/E7/hTERT-2 | human | uterine | Human uterine endometrium vascular endothelial progenit genetically-modified cells | |
| JCRB1546 | UCB408E7-TERT34 | human | umbilical c | Immortalized human umbilical cord blood-derived mesencl immortalized mesenchymal stem cell | |
| JCRB1548 | H-1A(H123) | mouse | bone marrow | | general cells |
| JCRB1549 | SK-OV-3-Luc | human | ovary | Luciferase stably expressing cell line (SK-OV-3; Human o' luciferase-expressing cancer cells | |
| JCRB1550 | HAdpc-25-Bmi-1-TERT | human | mesench | Immortalized human adipose tissue-derived mesenchymal immortalized mesenchymal stem cell | |
| JCRB1551 | KKU-055 | human | liver, gallb | Human poorly differentiated cholangiocarcinoma cell line e general cells | |
| JCRB1552 | HARA-B4 | human | human lun | originally human cc Deliveritive of the HARA cell. This cell I general cells | |
| JCRB1554 | MMNK-1 | human | liver, gallb | Expressing catalytic subunit of human Highly differentiated genetically-modified cells | |
| JCRB1555 | HAdpc-26-E6-Bmi-1-TERT | human | adipose tis | Human umbilical cord blood-derived mesenchymal stem c genetically-modified cells | |
| JCRB1557 | KKU-213 | human | biliary tract | Mixed papillary arc in publishing the re general cells | |
| JCRB1558 | RERF-LC-KJ-CMV-Luc | human | lung | Luciferase stably expressing cell line (RERF-LC-KJ; Hum: luciferase-expressing cancer cells | |
| JCRB1559 | MDA-MB-231-Luc | human | mammary | pLVSIN-luc transfected MDA-MB-231 cells. Luciferase sta luciferase-expressing cancer cells | |
| JCRB1563 | HAdpc-29-E7-TERT | human | adipose tis | Human adipocite-derived mesenchymal stem cells which v genetically-modified cells | |
| JCRB1564 | TMNK-1 | human | liver, gallb | Immortalized huma Expressed factor V Distribution restrict genetically-modified cells | |
| JCRB1566 | NP 2 | human | glioma | P53 mutation, P14) - | general cells |
| JCRB1567 | NP 3 | human | glioma | P53 mutation, P14) - | general cells |
| JCRB1568 | KKU-100 | human | biliary tract | poorly differentiate In publishing the re general cells | |
| JCRB1569 | NP 5 | human | glioma | P53 mutation, P14) - | general cells |
| JCRB1570 | H-1D(H162) | mouse | bone marr | Mouse bone marrow-derived osteoblast-like cell line. | general cells |
| JCRB1571 | F6B3 | mouse | brain | SR-RSV-induced b GFAP (glial fibrillary acidic protein pos general cells | |
| JCRB1572 | HAdpc-28-E6E7-TERT | human | adipocyte | Human adipocyte-derived mesenchymal stem cells which genetically-modified cells | |
| JCRB1574 | RSV-M | mouse | brain | SR-RSV-induced brain tumor cells | general cells |
| JCRB1575 | Onda 7 | human | glioma | P14/P15/P16 deletion, PTEN mutation | general cells |
| JCRB1576 | Onda 9 | human | glioma | P14/P15/P16 delet - | general cells |
| JCRB1577 | HSC-1-Luc | human | skin | Luciferase stably expressing cell line. (HSC-1: A cell line c luciferase-expressing cancer cells | |
| JCRB1578 | Onda 8 | human | glioma | P53 mutation, P14) - | general cells |
| JCRB1579 | Ishikawa 3-H-12-Luc | human | uterus | pLVSIN-luc transfected Ishikawa cells. | luciferase-expressing cancer cells |
| JCRB1580 | SR-CDF1 DBT | mouse | brain | SR-RSV-induced b sensitive to mouse coronavirus infect general cells | |
| JCRB1582 | TWNT-1 | human | liver | Expressing cataly Immortalized huma The distribution is l genetically-modified cells | Only for users in non-profit institute |
| JCRB1583 | Onda 9R | human | glioma | - | general cells |
| JCRB1585 | ACI-RCC | rat | renal | Semi-floating cells | general cells |
| JCRB1586 | HSC-5-Luc | human | skin | Luciferase stably expressing cell line. (HSC-5 : A cell line c luciferase-expressing cancer cells | |
| JCRB1588 | ACI-RCC-flLuc | rat | kidney | Luc expressing rat renal cell carcinoma cell line. | luciferase-expressing cancer cells |
| JCRB1591 | HLF-Luc | human | liver | Luciferase stably expressing cell line (HLF; Human Hepat luciferase-expressing cancer cells | |
| JCRB1592 | Hep G2-Luc | human | liver | Luciferase stably expressing cell line (Hep G2: Human He luciferase-expressing cancer cells | |
| JCRB1594 | SK-OV-3/CMV-Luc | human | ovary | pLL3.7-luc transf SK-OV-3 cells are resistant to tumor n luciferase-expressing cancer cells | |
| JCRB1596 | ACI-RCC-cbLuc | rat | kidney | Luc expressing rat renal cell carcinoma cell line. | luciferase-expressing cancer cells |
| JCRB1599 | hTf30-8H | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1600 | HuH-7-Luc | human | liver, gallb | pLVSIN-luc transfected HuH-7 cells. Luciferase stably exp luciferase-expressing cancer cells | |
| JCRB1601 | EMTOKA | human | Uterine bo | Human cell line derived from uterine carcinosarcoma. | general cells |
| JCRB1602 | hTf30-8H/sf | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1603 | hTf18-10C | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1605 | Onda 11 | human | brain | Human cell line derived from human glioma. | general cells |
| JCRB1606 | Onda 10 | human | brain | Human cell line derived from human glioma. | general cells |
| JCRB1607 | Onda 9R-B12 | human | brain | Human clonal cell line derived from human recurred gliom: general cells | |
| JCRB1608 | NP 8 | human | brain | Human cell line derived from human glioma. | general cells |
| JCRB1609 | OVCAR-3/CMV-Luc | human | ovary | Luciferase stably expressing cell line (OVCAR-3; human c luciferase-expressing cancer cells | |
| JCRB1610 | HLE-Luc | human | liver | Luciferase stably expressing cell line (HLE; Human Hepat luciferase-expressing cancer cells | |
| JCRB1611 | hTf18-10C/sf | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1612 | HuS-L12/10 | human | lung | Human immortalized cell line derived from MRC-5 cells. genetiically-modified cells | |
| JCRB1613 | hTf26-9B | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1614 | hTf26-9B/sf | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1616 | hTf/CHO-S #17 | hamster, C | ovary | pREP4-hTf transf This cell line produce human transferri general cells | |
| JCRB1617 | hTf9-11E | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1620 | hTf/CHO-S #29 | hamster, C | ovary | pREP4-hTf transf This cell line produce human transferri general cells | |
| JCRB1621 | hTf2-9F/sf | mouse x n | hybridoma | hybridoma This cell line produce monoclonal antit general cells | |
| JCRB1626 | pREP4 cont/CHO-S #5 | hamster, C | ovary | pREP4-cont transf This cell line produce human transferri general cells | |
| JCRB1627.1 | SK-BR-3-Luc | human | mammary | Luciferase stably expressing cell line (SK-BR-3 : breast c luciferase-expressing cancer cells | |
| JCRB1629 | hTf/CHO-S #17/sf | hamster, C | ovary | pREP4-hTf transf This cell line produ prod cell freezed by 10% general cells | |
| JCRB1629.01 | hTf/CHO-S #17/sf | hamster, C | ovary | pREP4-hTf transf This cell line produ prod cell freezed by Cry general cells | |
| JCRB1630 | DR-EcoScreen | mouse | liver, gallb | DR-EcoScreen is for evaluation of a rapid in vitro aryl hydri genetically-modified cells | |
| JCRB1631 | hTf/CHO-S #29/sf | hamster, C | ovary | pREP4-hTf transf This cell line produ prod cell freezed by 10% general cells | |
| JCRB1631.01 | hTf/CHO-S #29/sf | hamster, C | ovary | pREP4-hTf transf This cell line produ cell freezed by Cry general cells | |
| JCRB1632 | VV-1H831 | mouse x n | hybridoma | Mouse hybridoma producing mAb (IgG1) to hemagglutinin general cells | |

| | | | | | | | | |
|-------------|------------------------|------------|--------------------------------|---|---|---|----------------|---------------|
| JCRB1633 | HuH-6-Luc | human | liver, gallb | Luciferase stably expressing cell line (HuH-6; Human Hep luciferase-expressing cancer cells) | | | | |
| JCRB1634 | VV-B2D10 | mouse x n | hybridoma | Mouse hybridoma producing mAb to hemagglutinine (HA) general cells | | | | |
| JCRB1636 | pREP4 c/CHO-S #5/sf | chinese hz | ovary | pREP4-cont transfected CHO-S cells (cell freezed by 10% general cells | | | | |
| JCRB1636.01 | pREP4 c/CHO-S #5/sf | chinese hz | ovary | pREP4-cont transfected CHO-S cells (cell freezed by Cry: general cells | | | | |
| JCRB1638 | JHH-5-Luc | human | liver, gallb | Luciferase stably expressing cell line (JHH-5; Human hep: luciferase-expressing cancer cells) | | | | |
| JCRB1642 | KMS-11/BTZ | human | pleural effi | bortezomib-resistant maultiple myeloma cell line. | general cells | | | |
| JCRB1647 | 28SC-ES | human | peripheral endotoxin sensitive | human peripheral blood cell line (Belie | general cells | | | |
| JCRB1648 | ICH-ERMS-1 | human | pelvic tum | Human cell derived from anaplastic embryonal rhabdomyic | general cells | | | |
| JCRB1651 | VV-B6F4 | mouse x n | hybridoma | hybridoma | Mouse hybridoma producing mAb to v: general cells | | | |
| JCRB1657 | HuCCA-1 | human | intrahepatic bile duct | | Human cell line derived from cholangic | general cells | | |
| JCRB1658 | vdr2-4 | mouse | embryo | Control cell line used for the establishment of mouse hom mousehomozygous mutant ES cells | | | | |
| JCRB1670 | HT-1080/CMV-Luc | human | sarcoma | The cells contain activated N-ras oncogene. pLVSIN-luc tr | luciferase-expressing cancer cells | | | |
| JCRB1674 | HOS/CMV-Luc#2(c-1) | human | sarcoma | pLVSIN-luc transfected osteosarcoma | luciferase-expressing cancer cells | | | |
| JCRB1675 | B16-F10/CMV-Luc#2 | mouse | melanoma | pLVSIN-luc transfected B16-F10 cells. Luciferase stably e | luciferase-expressing cancer cells | | | |
| JCRB1679 | HeLa/CMV-Luc | human | uterine cei | pLVSIN-luc transfe first establish cell line in vitro from hum | luciferase-expressing cancer cells | | | |
| JCRB1680 | B16-F10/EF1a-Luc#1 | Mouse | Skin | pLVSIN-luc transfe Luciferase stably e MTA is required. | luciferase-expressing cancer cells | | | |
| JCRB1681 | MIA PaCa-2/CMV-Luc | human | pancrea | pLVSIN-luc transfected MIA PaCa-2 cells. Luciferase stab | luciferase-expressing cancer cells | | | |
| JCRB1683 | HARA CMV-Luc | human | lung cancer | | This cell line expre | luciferase-expressing cancer cells | | |
| JCRB1685 | HSC-3 CMV-Luc | human | oral | pLVSIN-luc transfe squamous carcinoma, tongue, age 64, | luciferase-expressing cancer cells | | | |
| JCRB1686 | 4T1 CMV-Luc #6 | mouse | mammary | pMSCV-luc transfected. Luciferase stably expressing cell | luciferase-expressing cancer cells | | | |
| JCRB1687 | G-361/CMV-Luc#2(c-2.5) | human | melanoma | malignant melanoma, melanin production, pLVSIN-luc tr | luciferase-expressing cancer cells | | | |
| JCRB1691 | RERF-Luc-Sq1CMV-Luc | human | lung squamous carcinoma | Established from human lung squamu | luciferase-expressing cancer cells | | | |
| JCRB1692 | HuO9 CMV-Luc | human | osteosarc | Liver/bone/kidney-1 Secretes osteocalcin & responds to vit | luciferase-expressing cancer cells | | | |
| JCRB1716 | Ex-3LL/CMV-Luc#1 | mouse | lung | pLVSIN-luc transfe Lewis lung carcinoma derived cell line | luciferase-expressing cancer cells | | | |
| JCRB1718 | H1718/CMV-Luc#6 | human | lung | pLVSIN-luc transfected H1718 cells. Luciferase stably exq | luciferase-expressing cancer cells | | | |
| JCRB1725 | KIK4 | Human | skin | intron 12, c.1845+2 Human fibroblast cell line with mutatig | general cells | | | |
| JCRB1727 | YS2 | human | skin | LDL receptor muta | LDL receptor mutation, Intron 12, c.18- | general cells | | |
| JCRB1729 | RMG-V | human | ovary | | Human ovarian clear cell adenocarcin | general cells | | |
| JCRB1731 | 293_hEcad/hSyn2 | human | kidney | G418 resistant (0.5 | μM cell line into whic | genetically-modified cells Only for users in non-profit institute | | |
| JCRB1733 | 293_hEcad | Human | kidney | Promycin resistor | 293 cell line into wt Approval required. | genetically-modified cells | | |
| JCRB1735 | 293_hSyn4 | human | kidney | G418 (0.5mg/mL) | 293 cell line into wt genetically-modified cells | Only for users in non-profit institute | | |
| JCRB1736 | 293_Syn1 | Human | kidney | G418 resistant (0.5 | 293 cell line into wt Approval required. | genetically-modified cells | | |
| JCRB1738 | XYFMGG01A | human | blood | 46, XY | EBV transformed E XY female, Androg | general cells | | |
| JCRB1742 | Vn-324 | Syrian har | brain | | Choroid plexus papilloma cell line of S: | general cells | | |
| JCRB1743 | XYFMGG18B | human | blood | 46, XY | EBV transformed E XY female | general cells | | |
| JCRB1744 | XYFMGG03D | human | blood | 46, XY | EBV transformed E XY female | general cells | | |
| JCRB1745 | XYFMGG23G | human | blood | 46, XY | EBV transformed E XY female | general cells | | |
| JCRB1746 | XYFMGG24K | Human | blood | 46, XY | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1747 | XYFMGC50H | human | blood | 46, XY | EBV transformed E XY female | general cells | | |
| JCRB1748 | XYFMGG06K | Human | blood | 46, XY | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1750 | XYFMGC07K | human | blood | 46, XY, r(9) | EBV transformed E XY female | general cells | | |
| JCRB1751 | XYFMGG19M | Human | blood | 46, XY | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1752 | XYFMGG16M | Human | blood | 46, XY | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1753 | XYFMGC44W | Human | blood | 46, X, Y fragment | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1754 | XYFMGC10S | Human | blood | 46, X, idic(Yp) | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1756 | XYFMGG04DM | Human | blood | | Parent of XYFMGC | EBV transformed B-cell lymphoblastoic | general cells | |
| JCRB1757 | XYFMGG04DP | Human | blood | | Parent of XYFMGC | EBV transformed B-cell lymphoblastoic | general cells | |
| JCRB1760 | XYFMGG04D | Human | blood | 46, X, i(Yq) | EBV transformed B-cell lymphoblastoic | general cells | | |
| JCRB1761 | AR-EcoScreen GR KO M1 | Chinese h | ovary | This cell is that stat AR-EcoScreen is f1 | MTA is required. | genetically-modified cells | | |
| JCRB1771 | KKK-D049 | human | liver, gallbladder | | Human well differentiated tubular aden | general cells | | |
| JCRB1772 | KKU-452 | human | bile duct | | Human poorly differentiated cholangio | general cells | | |
| JCRB1775 | KKK-D068 | human | liver, gallbladder | | Human well differentiated tubular aden | general cells | | |
| JCRB1776 | OCUG-1-Luc#5 | human | liver, gallb | Chromosome numl Luciferase stably e | Cells are tumoriger | luciferase-expressing cancer cells | | |
| JCRB1777 | KKK-D131 | human | liver, galbladder | | Human well differentiated tubular aden | general cells | | |
| JCRB1778 | KKU-023 | human | bile duct | | Human well differen | The depositor's apj | general cells | |
| JCRB1779 | KKK-D138 | human | liver, gallbladder | | Human adenosquamous carcinoma | ce | general cells | |
| JCRB1780 | BLACK-93A | Human | Pleural effusion | | Burkitt lymphoma cell line | | general cells | |
| JCRB1781 | TREE-92 | Human | Pleural effusion | | Burkitt lymphoma cell line | | general cells | |
| JCRB1782 | BALM-7 | human | blood | B cell leukemia | cell line | | general cells | |
| JCRB1783 | BALM-6 | human | bone marr | Karyotype informat | B cell leukemia | cell line | general cells | |
| JCRB1784 | BALM-8 | human | blood | B cell leukemia | cell line | | general cells | |
| JCRB1785 | BALM-9 | human | blood | B cell leukemia | cell line | | general cells | |
| JCRB1787 | BALM-9K | human | blood | B cell leukemia | cell Immnoglobulin phenotype kappa+,lamb | general cells | | |
| JCRB1788 | BALM-9KL | human | blood | B cell leukemia | cell Immnoglobulin phenotype kappa+,lamb | general cells | | |
| JCRB1790 | BALM-9N | human | blood | B cell leukemia | cell Immnoglobulin phenotype kappa-,lamb | general cells | | |
| JCRB1791 | BALM-13 | human | bone marr | Karyotype informat | B cell leukemia | cell line | general cells | |
| JCRB1792 | KU-Lu-MPPt3 | human | lung | EGFR exon19 del | Human lung adenocarcinoma | cell line | general cells | |
| JCRB1794 | BALM-9L | human | peripheral | Karyotype informat | B cell leukemia | cell line | general cells | |
| JCRB1795 | BALM-16 | human | blood | B cell leukemia | cell lacking expression of immunoglobulin | i | general cells | |
| JCRB1796 | BALM-18 | human | peripheral | Karyotype informat | B cell leukemia | cell line | Ig negative, A | general cells |
| JCRB1797 | BALM-20 | human | blood | B cell leukemia | cell Immnoglobulin phenotype and cell surf | general cells | | general cells |
| JCRB1803 | BALM-26 | human | bone marrow | | B cell leukemia | cell line, ALL-L2 cell lin | general cells | |
| JCRB2201 | A9(Neo1) | human chi | human mc | 93.6% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2202 | A9(Neo2) | human chi | human mc | 82.8% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2203 | A9(Neo3) | human chi | human mc | 88.6% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2204 | A9(Neo4) | human chi | human mc | 93.8% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2205 | A9(Neo5) | human chi | human mc | 90.4% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2206 | A9(Neo6) | human chi | human mc | 76.3% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2207 | A9(Neo7) | human chi | human mc | 93.2% cells include skin fibroblast, AG- | Human chromoson | general cells | | |
| JCRB2208 | A9(Neo8) | human chi | human mc | 98.7% cells include skin fibroblast, AG- | Human chromoson | general cells | | |

| | | | | |
|----------|----------------------|---|--|---|
| JCRB2209 | CHO(His9) | human chi human mc 70.2% cells include Cells include huma Human chromoson general cells | | |
| JCRB2210 | A9(Bsr10) | human chi human mc 93.3% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2211 | A9(neo11) | human chi human mc 91.5% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2212 | A9(Neo12) | human chi human mc 85.4% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2213 | A9(Hygro13) | human chi human mc 94.7% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2214 | A9(Hygro14) | human chi human mc 90.4% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2215 | A9(Neo15) | human chi human mc 96.3% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2216 | A9(Neo16) | human chi human mc 89.3% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2217 | A9(Neo17) | human chi human mc 42.7% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2218 | A9(Neo18) | human chi human mc 100% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2219 | A9(Neo19) | human chi human mc 90.0% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2220 | A9(Neo20) | human chi human mc 79.0% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2221 | A9(Hygro21) | human chi human mc 93.3% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2222 | A9(Hygro22) | human chi human mc 91.4% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB2223 | A9(Bsr\X) | human chi human mc 28.0% cells include skin fibroblast, AG- Human chromoson general cells | | |
| JCRB9002 | BW5147.G.1.4.OUA-R.1 | mouse hemo-lymp T cell lymphoma | Prepared from CRI general cells | |
| JCRB9003 | NCTC Clone 929 | mouse skin connective tissue, connective tissue, | Prepared from CCI general cells | |
| JCRB9004 | HeLa | human uterine cervix | first establish cell li EV=1.000 against i general cells | |
| JCRB9005 | BALB/3T3 clone A31 | mouse embryo | contact inhibited | Prepared from CCI general cells |
| JCRB9008 | MRC-5 | human fetus lung embryonal lung, djj fetal lung, normal d | Prepared from CCI general cells | N/A via Xenotech |
| JCRB9009 | A-431 | human epidermoi epidermoid carcino epidermoid carcino | Prepared from CRI general cells | |
| JCRB9010 | HeLa S3 | human uterine cei epitheloid carcinon adaptable to a susj | Prepared from CCI general cells | |
| JCRB9012 | RAJI | human hemo-lymp Burkitt lymphoma | Burkitt lymphoma | Prepared from CCI general cells |
| JCRB9014 | 3T3 L1 | mouse embryo | differentiate to adip The 3T3 L1 is a clonal subline of mous general cells | |
| JCRB9018 | CHO-K1 | hamster, C ovary | used as a chromos | Prepared from CCI general cells |
| JCRB9019 | 3T3-Swiss albino | mouse embryo | embryo fibroblasts embryo fibroblasts, Prepared from CCI general cells | |
| JCRB9020 | BHK-21(C-13) | hamster, E kidney | kidney kidney | Prepare from CCL general cells |
| JCRB9021 | U937 | human hemo-lymp express many mon Receptors for Fc & | Prepared from CRI general cells | |
| JCRB9023 | CCRF-CEM | human hemo-lymp peripheral blood, a peripheral blood, a A | maximum of 1-2 general cells | |
| JCRB9026 | L1210 | mouse hemo-lymp lymphocytic leuken | lymphocytic leuken | Prepared from CCI general cells |
| JCRB9027 | KB | human HeLa-cont | epidermoid carcino epidermoid carcino | Prepared from CCI general cells |
| JCRB9029 | MDCK (NBL-2) | dog kidney | kidney, epithelial lik | kidney, epithelial lik Prepared from CCI general cells |
| JCRB9030 | P3X63-Ag8.653 | mouse hemo-lymp non-secreting myel | non-secreting myel | Prepared from CRI general cells |
| JCRB9031 | MOLT-4 | human hemo-lymp peripheral blood, | a peripheral blood, a A | maximum of 1-2 general cells |
| JCRB9032 | SKJ-237-71 | mouse antibody-p hybridoma | cell line producing monoclo | Prepared from CRI general cells |
| JCRB9033 | STK1 | mouse antibody-p hybridoma, | monocl hybridoma, monocl | Prepared from CR general cells |
| JCRB9034 | LLC-RK1 | rabbit kidney | an aneuploid cell li susceptible to virus | Prepare from CCL general cells |
| JCRB9036 | 104C1 | guinea pig fetal body | fetal carcass, benzo(a)pyrene transfor | Prepared from CRI general cells |
| JCRB9039 | Pt K1 | marsupial kidney | Kidney | Kidney Prepared from CCI general cells |
| JCRB9040 | PK(15) | pig kidney | Kidney | Kidney Prepared from CCI general cells |
| JCRB9041 | T98G | human neural | G1 arrested when anchorage-indeper | Prepared from CRI general cells |
| JCRB9042 | WI-26 VA4 | human lung | lung, SV-40 virus tr | lung, SV-40 virus tr Prepared from CCI general cells |
| JCRB9044 | Detroit 532 | human skin | Down's syndrome | (Prepared from CCI general cells |
| JCRB9045 | Detroit 539 | human skin | Skin, Down's syndr | Skin, Down's syndr Prepared from CCI general cells |
| JCRB9046 | GH1 | rat pituitary | Pituitary tumor | Pituitary tumor, sor |
| JCRB9047 | GH3 | rat pituitary | Pituitary tumor, Mt | Pituitary tumor, Mt Prepared from CCI general cells |
| JCRB9048 | MOLT-3 | human hemo-lymp | Peripheral blood, a Peripheral blood, | Peripheral blood, a Prepared from CRI general cells |
| JCRB9050 | IMR-32 | human neural | N-myc amplification | transplantable to n |
| JCRB9051 | KG-1 | human hemo-lymp | bone marrow, acut bone marrow, acut | Prepared from CCI general cells |
| JCRB9052 | Pt K2 (NBL-5) | marsupial kidney | kidney | kidney Prepared from CCI general cells |
| JCRB9053 | L-2 | rat lung | adult, lung, female | adult, lung, female Prepared from CCI general cells |
| JCRB9054 | IMR-90 | human lung | fetal lung, diploid, f | fetal lung, diploid, f Prepared from CCI general cells |
| JCRB9055 | R2C | rat testis | Leydig cell testicul | Leydig cell testicul Prepared from CCI general cells |
| JCRB9056 | Y-1 | mouse adrenal cc | adrenal tumor | adrenal tumor Prepared from CCI general cells |
| JCRB9057 | WI-38 VA13 sub 2 RA | human lung | lung, SV40 virus tr | lung, SV40 virus tr; Prepared from CCI general cells |
| JCRB9058 | RPMI 2650 | human nasal sept | quasi-diploid | squat quasi-diploid squat |
| JCRB9059 | 3T6-Swiss albino | mouse embryo | embryo, collagen-s | embryo, collagen-s Prepared from CCI general cells |
| JCRB9060 | MRC-9 | human lung | fetal lung, diploid | f fetal lung, diploid |
| JCRB9061 | PA-1 | human teratocarc | ascitic fluid cells, o | ascitic fluid cells, o May be prepared fr general cells |
| JCRB9062 | HS-Sultan | human hemo-lymphocytic | | Originally thought e general cells |
| JCRB9063 | WIL2-NS | human hemo-lymp | B lymphocyte, non- | |
| JCRB9064 | LC540 | rat testis | leidig cell testicula | leidig cell testicula Prepared from CCI general cells |
| JCRB9065 | G-401 | human kidney | Wilms' tumor (emb | Prepared from CRI general cells |
| JCRB9066 | Chang Liver | human HeLa-cont | liver, HeLa marker | This cell line was r |
| JCRB9067 | NRK-49F | rat kidney | normal kidney | fibro normal kidney fibro Prepared from CRI general cells |
| JCRB9068 | 293 | human kidney | | transformed embr |
| JCRB9069 | SW-13 | human adrenal cortex | adenocarcinoma, A | Prepared from CCI general cells |
| JCRB9070 | G-402 | human kidney | renal leiomyoblast | renal leiomyoblastic |
| JCRB9071 | Daudi | human hemo-lymp | Burkitt lymphoma | Burkitt's lymphoma Prepared from CCI general cells |
| JCRB9072 | RD | human sarcoma | rhabdomyosarcoma | Prepared from CCI general cells |
| JCRB9074 | G-361 | human melanoma | malignant melanon | malignant melanon Prepared from CRI general cells |
| JCRB9075 | NCTC Clone 1469 | mouse liver | liver | liver CCL 9.1. |
| JCRB9076 | FHs 74 Int | human intestine | fetal small intestine | fetal small intestine Prepared from CCI general cells |
| JCRB9077 | MMT 060562 | mouse mammary | spontaneous | mami spontaneous |
| JCRB9078 | De Gin | human skin | skin, osteoporosis | skin, osteoporosis Prepared from CRI general cells |
| JCRB9079 | Ran Nor | human skin | skin, osteoporosis | skin, osteoporosis Prepared from CRI general cells |
| JCRB9080 | C3H/10T1/2 clone 8 | mouse embryo | embryo, contact se | embryo, contact se Prepared from CCI general cells |
| JCRB9081 | L6 | rat muscular | : skeletal muscle | skeletal muscle my Prepared from CRI general cells |
| JCRB9083 | LoVo | human colon | colon, adenocarcin | colon, adenocarcin Prepared from CCI general cells |
| JCRB9084 | Sp2/O-Ag14 | mouse hemo-lymp | hybridoma, non-se | hybridoma, non-se Prepared from CRI general cells |

| | | | | | |
|----------|----------------------|-----------------------|--|---|---|
| JCRB9085 | P3IX63Ag8U.1 | mouse | hemo-lym ⁺ myeloma derivative | myeloma derivative | Prepared from CRL general cells |
| JCRB9086 | HeLa 229 | human | uterine cervix epithelial | carcinoma epithelial | Prepared from CCL general cells |
| JCRB9087 | RPMMI 1846 | hamster, \pm | melanoma melanotic | melanoma melanotic | Prepared from CCL general cells |
| JCRB9089 | CCRF-S-180 II | mouse | sarcoma | sarcoma 180 deriv | Prepared from CCL general cells |
| JCRB9090 | Ehrlich ascites | mouse | ascites | Ehrlich-Lettre ascit | many floating cells |
| JCRB9091 | Y3-Ag 1.2.3 | rat | hemo-lym ⁺ myeloma, kappa \pm | myeloma, kappa \pm | Prepared from CRL general cells |
| JCRB9092 | 7D4 | mouse x r; antibody-p | hybridoma, anti-mu | hybridoma, anti-mu | Prepared from CRL general cells |
| JCRB9093 | WEHI-3 | mouse | hemo-lym ⁺ myelomonocytic | lei | myelomonocytic lei |
| JCRB9094 | DLD-1 | human | colon | adenocarcin colon, adenocarcin | Prepared from CCL general cells |
| JCRB9095 | HEPM | human | embryo | embryonic palatal r | embryonic palatal r |
| JCRB9096 | C6 | rat | neural | glial cell tumor | glial cell tumor |
| JCRB9097 | I-10 | mouse | testis | leydig cell testicula | This cell line is a cl |
| JCRB9098 | ATT-20 | mouse | pituitary | ACTH producing | Cells are subcultured |
| JCRB9099 | SV-T2 | mouse | embryo | SV40 virus-transf | SV40 virus-transf |
| JCRB9100 | Indian Muntjac | Indian Muntjac | skin | Muntjak is a primit | Prepared from CCL general cells |
| JCRB9101 | CCD-14Br | human | lung | bronchus, diploid | bronchus, diploid |
| JCRB9102 | Detroit 525 | human | skin | skin, Turner's Sync | skin, Turner's Sync |
| JCRB9103 | KD | human | oral | lip, fibroblast | lip, fibroblast |
| JCRB9104 | Detroit 510 | human | skin | skin, galactosemia | skin, galactosemia |
| JCRB9105 | P3IX63Ag8 | mouse | hemo-lym ⁺ | resistant to 10 ⁻⁴ M IgG1(gamma,kappa) | Prepared from TIB general cells |
| JCRB9106 | MPC 11 OUA | mouse | hemo-lym ⁺ myeloma | myeloma | Prepared from TIB general cells |
| JCRB9107 | P3/NS1/1-Ag4-1(NS-1) | mouse | hemo-lym ⁺ non-secreting | myel non-secreting | myel Prepared from TIB general cells |
| JCRB9108 | J774A.1 | mouse | hemo-lymphocytic | monocyte-macroph | Prepared from TIB general cells |
| JCRB9109 | P3.6.2.8.1 | mouse | hemo-lym ⁺ IgG1 secreting myt | IgG1 secreting myt | Prepared from TIB general cells |
| JCRB9110 | PC-3 | human | prostate g | adenocarcinoma | general cells |
| JCRB9111 | BeWo | human | placenta | trophoblastic cell lli | The first human enc |
| JCRB9113 | HT-1080 | human | sarcoma | The cells contain activated N-ras | oncogene. |
| JCRB9114 | BRL-3A | rat | liver | liver | liver producing MS. Additional Informat |
| JCRB9115 | SW837 | human | rectum | Rectum, adenocarcinoma | general cells |
| JCRB9117 | BCL1 Clone CW13.20 | mouse | hemo-lym ⁺ | murine B cell leuk | murine B cell leukemia, subclone 3B3 |
| JCRB9118 | SCC-4 | human | oral | tongue,squamous | tongue,squamous |
| JCRB9119 | Ramos(RA1) | human | hemo-lym ⁺ | Burkitt lymphoma (| Burkitt lymphoma (EBV negative) |
| JCRB9122 | SIRC | rabbit | cornea | susceptible to rubella virus, toxicolog: | general cells |
| JCRB9124 | M.dunn(Clon III8C) | pygmy m tail | Cloned at P14 usin | Supports the growth of all 4 classes of | general cells |
| JCRB9125 | PG-4(S+L-) | cat | neural | Transformed with | Support Type C retroviruses. |
| JCRB9128 | Mv 1 Lu | mink | lung | This cell line is useful for focus forming | general cells |

| JCRB No. | Cell Name | ANIMAL | TISSUE | GENETICS | CHARACT | COMMENT | celltype | Order type |
|------------|---------------------------|-----------------|--------------------|---|---|---|------------------------------------|--|
| JCRB0030 | CHL/IU(IVGT) | Chinese hamster | lung | This line is derivative of the WIL-2 cell line. The cells are heterozygous at the thymidine kinase (TK) locus, and can be used to quantitatively detect forward mutation at 3 loci. | Standard cell line for chromosome aberation test. | Same as the cell line in ECACC | general cells | Only for researchers in Asian countries |
| JCRB1435 | TK6(IVGT) | human | spleen | | | Same as the cell line in ECACC | general cells | Only for researchers in Asian countries |
| JCRB1771 | KKK-D049 | human | liver, gallbladder | | Human well differentiated tubular adenocarcinoma cell line established from biliary tract. | | general cells | |
| JCRB1772 | KKU-452 | human | bile duct | | Human poorly differentiated cholangiocarcinoma cell line established from biliary tract. | | general cells | |
| JCRB1775 | KKK-D068 | human | liver, gallbladder | | Human well differentiated tubular adenocarcinoma cell line established from biliary tract. (with micropapillary foci) | | general cells | |
| JCRB1776 | OCUG-1-Luc#5 | human | liver, gallbladder | Chromosome number distribute in a broad range from 52 to 139. | Luciferase stably expressing cell line (OCUG-1; human gall bladder carcinoma). | Cells are tumorigenic in nude mice. Poorly differentiated adenocarcinoma. | luciferase-expressing cancer cells | |
| JCRB1777 | KKK-D131 | human | liver, gallbladder | | Human well differentiated tubular adenocarcinoma cell line established from biliary tract. | | general cells | |
| JCRB1778 | KKU-023 | human | bile duct | | Human well differentiated cholangiocarcinoma cell line established from biliary tract. | The depositor's approval is required. | general cells | |
| JCRB1779 | KKK-D138 | human | liver, gallbladder | | Human adenosquamous carcinoma cell line established from biliary tract. | | general cells | |
| JCRB1782 | BALM-7 | human | blood | B cell leukemia cell line | | | general cells | |
| JCRB1783 | BALM-6 | human | bone marrow | Karyotype information included in established papers | B cell leukemia cell line | | general cells | |
| JCRB1784 | BALM-8 | human | blood | B cell leukemia cell line | | | general cells | |
| JCRB1785 | BALM-9 | human | blood | B cell leukemia cell line | | | general cells | |
| JCRB1787 | BALM-9K | human | blood | B cell leukemia cell line | Immnoglobulin phenotype kappa+,lambda- | | general cells | |
| JCRB1788 | BALM-9KL | human | blood | B cell leukemia cell line | Immnoglobulin phenotype kappa+,lambda+ | | general cells | |
| JCRB1790 | BALM-9N | human | blood | B cell leukemia cell line | Immnoglobulin phenotype kappa-,lambda- | | general cells | |
| JCRB1791 | BALM-13 | human | bone marrow | Karyotype information included in established papers | B cell leukemia cell line | | general cells | |
| JCRB1792 | KU-Lu-MPPt3 | human | lung | EGFR exon19 del E746-A750 | Human lung adenocarcinoma cell line with a micropapillary pattern. | | general cells | |
| JCRB1794 | BALM-9L | human | peripheral blood | Karyotype information included in established papers | B cell leukemia cell line | | general cells | |
| JCRB1795 | BALM-16 | human | blood | B cell leukemia cell line, Ig negative, ALL-L3 cell line, t(8;22)(q24;q11), hypercalcemia, mRNA expression of PTHrP | lacking expression of immunoglobulin chains. | | general cells | |
| JCRB1796 | BALM-18 | human | peripheral blood | Karyotype information included in established papers | B cell leukemia cell line, Ig negative, ALL-L3 cell line, t(8;14)(q24;q32) | | general cells | |
| JCRB1797 | BALM-20 | human | blood | B cell leukemia cell line | Immnoglobulin phenotype and cell surface marker information are included in established paper. | | general cells | |
| JCRB1803 | BALM-26 | human | bone marrow | | B cell leukemia cell line, ALL-L2 cell line, hypercalcemia | | general cells | |
| JCRB0709.2 | L5178Y TK+/- 3.7.2c(IVGT) | mouse | lymph node | Mouse TK+/- heterozygotic subline of the L5178Y. | | | general cells | Not available (Contact JCRB if customer need it) |