

Cell types counted by Cellometer™ automated cell counters

Have We Counted Your Cell Type?

We have been counting cells since 2003. Browse through our list of over 1,800 cell types to see if we have counted yours.

For research use only. Not for use in diagnostic procedures.

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
697	Blood	B-lymphocytes	Suspension
721.211	unknown	unknown	unknown
5637	Bladder	Bladder epithelial cells	Adherent
A3/Kawakami	Lymphatic system	B-lymphocyte	unknown
B-lymphocyte	Blood	B-lymphocyte	Suspension
1205Lu	Skin	Melanocytes	unknown
1321N1	Brain	Astrocyte	Either
143B	Bone	Osteocyte	Adherent
143B pML BK TK	Bone	Osteocyte	Either
145M-Twi-hGALC-G270D	Brain	Glial	Adherent
16-HBE	Bronchi	Bronchial epithelial	Adherent
184B5	Mammary gland	Breast epithelial	Adherent
1A6	Bladder	Bladder epithelial cells	unknown
2008C	unknown	unknown	unknown
201T	Lung	Lung cells	unknown
22RV1	Prostate	Prostate cells	Adherent
23132/87	Stomach	Stomach cells	Adherent
273T	Lung	Lung cells	unknown
293T	Kidney	Debated	Suspension
2B8	Lymphatic system	Lymphocyte	Suspension
3A-KO	unknown	unknown	unknown
3F1	unknown	unknown	unknown
3T3	N/a	Fibroblast	Adherent
3T3-L1	N/a	Fibroblast (Adipocyte-like phenotype)	Adherent
4244135 GCT	Skin	Lung cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
42-MG-BA	Brain	Astrocytes	unknown
621-101	Kidney	Kidney cells	unknown
639-V	Urinary Tract	ureter transitional cell	Adherent
647-V	Urinary Tract	urothelial Bladder carcinoma	Adherent
66cl4	Mammary gland	Mammary gland epithelial	Adherent
67NR	Mammary gland	Mammary gland epithelial	Adherent
722-MP14	Pancreas	unknown	unknown
769-P	Kidney	Kidney epithelial cell	Adherent
786-O	Kidney	renal epithelial	Adherent
8305C	Thyroid	Thyroid cells	Adherent
84-31	unknown	unknown	unknown
8505C	Thyroid	Thyroid cells	Adherent
8-MG-BA	Brain	Astrocyte	Adherent
950MP11	Pancreas	unknown	unknown
950-MP5	Pancreas	unknown	unknown
9L	Brain	Astrocyte	Adherent
A101D	Skin	Melanocytes	Adherent
A13A	Pancreas	unknown	unknown
A172	Brain	Astrocyte	Adherent
A172-BM	Brain	Astrocyte	Adherent
A172-P	Brain	Astrocyte	Adherent
A172-sh3MSH6-BM	Brain	Astrocyte	Adherent
A172-sh5MSH6-BM	Brain	Astrocyte	Adherent
A172-shMSH6-BM	Brain	Astrocyte	Adherent
A172-shMSH6-P	unknown	unknown	unknown
A172-TR3-BM	unknown	unknown	unknown
A2.1	Pancreas	Pancreas cell	unknown
A-204	Muscle	Striated muscle cell	Adherent
A2058	Skin	Skin cell	Adherent
A-253	Salivary gland	Salivary gland epithelial	Adherent
A2780	Ovary	Ovarian epithelial cell	Adherent
A2780ADR	Ovary	Ovarian epithelial cell	Adherent
A2780cis	Ovary	Ovarian epithelial cell	Adherent
A3	Blood	T-lymphocyte	Suspension
A373-C6	Skin	Skin cell	Adherent
A-375	Skin	Skin cell	Adherent
A375.S2	Skin	Skin cell	Adherent
A-388	Skin	cutaneous squamous cell	Adherent
A-427	Lung	Lung cells	Adherent
A431	Skin	Skin cell	Adherent
A498	Kidney	Kidney cells	Adherent
A4-Fuk	Blood	B-lymphocyte	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
A549	Lung	"adenocarcinomic alveolar basal epithelial cells"	Adherent
A549cisR	Lung	"adenocarcinomic alveolar basal epithelial cells"	Adherent
A673	Muscle	Ewing sarcoma cells	Adherent
A6L	Pancreas	Pancreas cell	unknown
A704	Kidney	Kidney cells	Adherent
ABC-1	Lung	Lung cells	unknown
ABCG2-HEK293	unknown	unknown	unknown
ACC 8-2	Head & Neck	unknown	unknown
ACC2	Cervix	Cervical cell	Adherent
ACC3	Cervix	Cervical cell	Adherent
ACCS	Bladder	Bladder epithelial cells	unknown
Ace1	Prostate	Prostate cells	unknown
ACHN	Kidney	Kidney cells	Adherent
ACN	Nervous system	Neuroblast	Adherent
ACS	unknown	unknown	unknown
AdMSC cells	unknown	unknown	unknown
AESF-1	unknown	unknown	unknown
AGS	Stomach	Stomach cells	Adherent
AKR-2B	Embryo	Embryonic cells (mouse)	Adherent
ALL-PO	Blood	B-lymphocytes	Suspension
ALL-SIL	Blood	T-lymphocytes	Suspension
AM-38	Brain	Astrocytes	unknown
AML12 cells	Liver	Hepatocyte	Adherent
AML-193	Peripheral blood	Myeloid cells	Suspension
AMO-1	Peripheral blood	Plasma cell	unknown
AN3CA	Uterus	Uterine endometrial cancer cells	Adherent
ARH-77	Blood	B-lymphoblastoid cell	Suspension
ASH-3	Thyroid	Thyroid cells	unknown
ASLNCaP-FGC	unknown	unknown	unknown
AsPC-1	Pancreas	Pancreas cell	Adherent
Aspergillus niger	unknown	Fungus	unknown
ATCC	unknown	unknown	unknown
ATN-1	Blood	T-lymphocytes	unknown
AtT20	Pituitary gland	Pituitary gland cells	Adherent
AU565	Breast	Breast epithelial	Adherent
AX3	unknown	unknown	unknown
AZ-521	Duodenum	Duodenal cells	Adherent
AZD1480	unknown	unknown	unknown
B. afzelii	unknown	Bacteria	unknown
B. burgdorferi	unknown	Bacteria	unknown
B. garinii	unknown	Bacteria	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
B. thetaiotaomicron	unknown	Bacteria	unknown
B1 B cells	Blood	Blood	unknown
B10	Blood	Blood	unknown
B16-F10	Skin	Skin cell	Adherent
B2-17	Brain	Astrocyte	unknown
Ba/F3	Lymphatic system	Lymphocyte	Suspension
Ba/F3	Blood	B-lymphocytes	Suspension
Bacillus anthracis	unknown	Bacteria	unknown
Bacillus atrophaeus	unknown	Bacteria	unknown
Bacillus cereus	unknown	Bacteria	unknown
Bacteroides thetaiotaomicron	unknown	Bacteria	unknown
BAL	Bronchi	Bronchoalveolar cells	unknown
BALL-1	Blood	B-lymphocytes	Suspension
Barramundi Muscle	Muscle	Striated muscle cell	Adherent
BB30-HNC	Head & Neck	Squamous cells	unknown
BB49-HNC	Head & Neck	Squamous cells	unknown
BB65-RCC	Kidney	Kidney cells	unknown
BC-1	Lymphatic system	Lymphocyte	Suspension
BC-2	Lymphatic system	Lymphocyte	Suspension
BCIRL-AtE-CLG15	unknown	unknown	unknown
BCIRL-Cc-AM	unknown	unknown	unknown
BCIRL-Cc-JG	unknown	unknown	unknown
B-CPAP	Thyroid	Thyroid cells	unknown
BDCM	Blood	B-lymphocyte	Suspension
BE(2)-C/ADR(5)	Nervous system	Neuroblast	Adherent
BE-13	Blood	T-lymphocyte	Suspension
BEAS-2B	Bronchi	Bronchial epithelial	Adherent
Beauveria bassiana	unknown	unknown	unknown
Becker	Brain	Astrocyte	unknown
BEN	Lung	Lung cells	Adherent
BEP2D	Bronchi	Bronchial epithelial	unknown
BeWo	Placenta	Placental cells	Adherent
BFTC-905	Urinary Tract	Bladder epithelial cells	Adherent
BFTC-909	Kidney	Kidney cells	Adherent
BHK21	Kidney	Kidney cells	Adherent
BHT-101	Thyroid	Thyroid cells	Adherent
BHY	Mouth	Squamous cells	unknown
BIC-1	Colon	Colorectal epithelial cells	Adherent
BICR 10	Mouth	Squamous cells	Adherent
BICR 22	Tongue	Squamous cells	Adherent
BICR 31	Tongue	Squamous cells	Adherent
BICR 78	Oral Alveolus	Squamous cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
BL-41	Lymphatic system	B-lymphocyte	Suspension
BL-70	Lymphatic system	B-lymphocyte	Suspension
BON-1	Pancreas	Pancreas cell	unknown
BONNA-12	Leukemia*	B-lymphocyte	Suspension
bovine neutrophils	Blood	Neutrophil	unknown
BPH-1	Prostate	Prostate epithelial cell	Adherent
BSL2KA	Lymphoma	B-lymphocyte	unknown
BT-1008	Turbinate	Bovine turbinate cells	Adherent
BT-20	Breast	Breast epithelial	Adherent
BT-474	Breast	Breast epithelial	Adherent
BT-483	Breast	Breast epithelial	Adherent
BT549	Breast	Breast epithelial	Adherent
BT-B	Cervix	Cervical cell	Adherent
BU-ML	Skin	unknown	unknown
BV-173	Blood	pre-B cell	Suspension
BV-2	Brain	MicroAstrocytes	Suspension
BxPC-3	Pancreas	Pancreas cell	Adherent
C. albicans	unknown	unknown	Neither
C. elegans GERM CELLS	Gonad	Germ cell	Neither
C. glabrata	unknown	unknown	Neither
C. infirmominiatum	unknown	unknown	Neither
C. reinhardtii	unknown	unknown	Neither
C170	Colon	Colorectal epithelial cells	unknown
C2BBe1	Colon	Colorectal epithelial cells	Adherent
C2C12	Muscle	Myeloblast	Adherent
C32	Colon	Colorectal epithelial cells	Adherent
C-33 A	Cervix	Cervical cell	Adherent
C3H	Embryo	Fibroblasts	Adherent
C4	Liver	Liver cell	Adherent
C-4 I	Cervix	Cervical cell	Adherent
C-4 II	Cervix	Cervical cell	Adherent
C42	Prostate	Prostate epithelial cell	Adherent
C5/MJ	Umbilical cord	T-lymphocyte	Suspension
C-81	Prostate	Prostate epithelial cell	Adherent
C8166	Leukemia	T-lymphocyte	Suspension
Ca Ski	Cervix	Cervical epithelial cells	Adherent
CA46	Lymphoma	B-lymphocyte	Suspension
Ca9-22	Head & Neck	Squamous cells	Adherent
Caco-2	Colon	Colorectal epithelial cells	Adherent
CAKI	Kidney	Kidney	Adherent
Caki-1	Kidney	Kidney	Adherent
CAL 27	Mouth	Squamous cells	Adherent
CAL-120	Breast	Mammary epithelial cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
CAL-12T	Lung	Lung cells	Adherent
CAL-148	Breast	Mammary epithelial cells	Suspension
CAL-29	Bladder	Bladder epithelial cells	Adherent
CAL-33	Tounge	Squamous cells	Adherent
CAL-39	Vulva	Squamous cells	Adherent
CAL-51	Breast	Breast epithelial	Adherent
CAL-54	Kidney	Kidney cells	Adherent
CAL-62	Thyroid	Thyroid cells	Adherent
CAL-72	Bone	Osteocyte	Adherent
CAL-78	Muscle	Muscle cell	Adherent
CAL-85-1	Breast	Breast epithelial	Adherent
Calu-1	Lung	Lung cells	Adherent
Calu-3	Lung	Lung cells	Adherent
Calu-6	Lung	Lung cells	Adherent
CAMA-1	Breast	Breast epithelial	Adherent
Caov-3	Ovary	Ovarian carcinoma	Adherent
Caov-4	Ovary	Ovarian carcinoma	Adherent
Capan-1	Pancreas	Pancreas cell	Adherent
Capan-2	Pancreas	Pancreas cell	Adherent
CaR-1	Colon	Colorectal epithelial cells	unknown
CAS-1	Brain	Astrocyte	Adherent
Castration resistant prostate cancer cells	Prostate	Prostate epithelial cell	unknown
CC-1690	unknown	unknown	Neither
CC-3269	unknown	unknown	Neither
CC-425	unknown	unknown	Neither
CCD 1105 KIDTr	Fetus	Fetal epithelial	Adherent
CCD 841 CoN	Colon	Colorectal epithelial cells	Adherent
CCD-14Br	Bronchi	Bronchial epithelial	unknown
CCD-18Co	Colon	Colon fibroblast	Adherent
CCF-STTG1	Brain	Astrocyte	Adherent
CCK-81	Colon	Colorectal epithelial cells	unknown
CCL-2	unknown	unknown	unknown
CCL-34	Kidney	Kidney epithelial cell	Adherent
CCL-57	Skin	Skin fibroblast	Adherent
CCL-88	unknown	Epithelial	Adherent
CC-LP-1	Biliary Tract	Biliary cell	unknown
CCRF-CEM	Peripheral blood	T-lymphoblast	Suspension
CCRF-HSB-2	Peripheral blood	T-lymphoblast	Suspension
CCRF-SB	Peripheral blood	B-lymphoblast	Suspension
CCRT	unknown	unknown	unknown
CCSWI	unknown	unknown	unknown
CD-1	unknown	unknown	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
CD133+	unknown	unknown	unknown
CD14+ Monocytes	unknown	unknown	unknown
CD24	unknown	unknown	unknown
CD3	unknown	unknown	unknown
CD3+ B	unknown	unknown	unknown
CD3+ T	unknown	unknown	unknown
CD34 + cellsPrimary	unknown	unknown	unknown
CD36	unknown	unknown	unknown
CD4+ T cells	unknown	unknown	unknown
CD4+ Th2	unknown	unknown	unknown
CD44	unknown	unknown	unknown
CD51+	unknown	unknown	unknown
CD8a	unknown	unknown	unknown
CDC 1185	unknown	unknown	unknown
CEM T	Blood	T-lymphoblast	Suspension
CEM/C1	Blood	T-lymphoblast	Suspension
CEM/C2	Blood	T-lymphoblast	Suspension
CEM-ON	Blood	T-lymphoblast	Suspension
CESS	Blood	Lymphoblast	Suspension
CFPAC-1	Pancreas	Pancreas cell	Adherent
CFs rat	unknown	unknown	unknown
CG7	unknown	unknown	unknown
CGTH-W-1	Thyroid	Thyroid cells	Adherent
Ch 1 Es	unknown	Fibroblast	Adherent
Ch8	Bone	Osteocyte	unknown
CHA hESCs-H1	Blastocyst	Embryonic stem cell (human)	unknown
ChaGo-K-1	Lung	Bronchial epithelial	Adherent
CHL-1	Skin	Melanocytes	Adherent
Chlamydomonas reinhardtii	unknown	unknown	Neither
Chlorella vulgaris	unknown	unknown	Neither
CHO-EBNA GS	Ovary	Ovary cell	Adherent
CHO-K1	Ovary	Ovary cell	Adherent
CHO-K1-BH4	Ovary	Ovary cell	Adherent
Chondrocytes	Cartilage	Chondrocyte	unknown
CHP-126	Nervous system	Neuroblast	Adherent
CHP-134	Nervous system	Neuroblast	Adherent
CHP-212	Nervous system	Neuroblast	Adherent
CHSA 0011	Carilage	Chondrocyte	unknown
CHSA 0108	Carilage	Chondrocyte	unknown
CHSA8926	Carilage	Chondrocyte	unknown
CI-1	Lymphatic system	B-lymphocyte	Suspension
CIK cells	Kidney	Kidney cells	unknown
CL-1	Blood	Leukocyte	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
CL-11	Colon	Colorectal epithelial cells	Adherent
CL-14	Colon	Colorectal epithelial cells	Adherent
CL-34	Colon	Colorectal epithelial cells	Adherent
CLL	Blood	B-lymphocytes	Suspension
Closterium ehrenbergii	unknown	unknown	Neither
CMK	Blood	Leukocyte	Suspension
CML-T1	Peripheral blood	T-lymphocyte	Suspension
CoCM-1	Colon	Colorectal epithelial cells	unknown
Collagen type I	unknown	unknown	unknown
COLO 201	Colon	Colorectal epithelial cells	Suspension
COLO 205	Colon	Colorectal epithelial cells	Suspension
COLO 206F	Colon	Colorectal epithelial cells	Adherent and suspension
COLO 320DM	Colon	Colorectal epithelial cells	Adherent and suspension
COLO 684	Uterus	Uterine cells	Suspension
COLO 685	Uterus	Uterine cells	Suspension
COLO 741	Skin	Melanocytes	Adherent
COLO 792	Skin	Melanocytes	Adherent
COLO 853	Skin	Melanocytes	Adherent
COLO 857	Skin	Melanocytes	Adherent
COLO 858	Skin	Melanocytes	Adherent
COLO-320	Colon	Colorectal epithelial cells	Adherent and suspension
COLO-320-HSR	Colon	Colorectal epithelial cells	Adherent and suspension
COLO-668	Lung	Lung oat cells	Adherent and suspension
COLO-678	Colon	Colorectal epithelial cells	Adherent
COLO-679	Skin	Melanocytes	Adherent
COLO-680N	Esophagus	Esophagus epithelial cells	Adherent
COLO-699	Lung	Lung cells	Adherent
COLO-704	Ovary	Ovarian cells	Adherent
COLO-783	Skin	Melanocytes	Adherent
COLO-800	Skin	Melanocytes	Adherent
COLO-818	Skin	Melanocytes	Adherent
COLO-824	Breast	Mammary epithelial cells	Adherent and suspension
COLO-829	Skin	Melanocytes	Adherent
COLO-849	Skin	Melanocytes	Adherent
Conidia	unknown	unknown	Neither
COR-L 105	Lung	Lung cells	Adherent
COR-L 23/CPR	Lung	Lung cells	Adherent
COR-L23	Lung	Lung cells	Adherent
COR-L279	Lung	Lung cells	Suspension
COR-L47	Lung	Lung cells	Suspension
COR-L51	Lung	Lung cells	Suspension
COR-L88	Lung	Lung cells	Adherent and suspension
COR-L95	Lung	Lung cells	Suspension

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Corneal Endothelial Cells (HCEC)	Cornea	Corneal endothelial cells	unknown
Corneal epithelial stem cells (CESCs)	Cornea	Corneal epithelial stem cells	unknown
Cortical Neuron Culture	Brain	Cortical neurons	unknown
COS-1	Kidney	Kidney fibroblast	Adherent
COS-7	Kidney	Kidney fibroblast	Adherent
CP50-MEL-B	Skin	Melanocytes	unknown
CP66-MEL	Skin	Melanocytes	unknown
CP67-MEL	Skin	Melanocytes	unknown
CPC-N	Skin	Melanocytes	Suspension
CRL11372	Bone	Osteoblast	Adherent
CRO-AP2	Lymphatic system	B-lymphocytes	Suspension
CRO-AP3	Lymphatic system	B-lymphocytes	Suspension
CRO-AP5	Lymphatic system	B-lymphocytes	Suspension
CS1	Bone	Chondrocyte	unknown
CS1R	Bone	unknown	unknown
CS2	Bone	unknown	unknown
CS3	Bone	unknown	unknown
CT-26	Colon	Mouse Colorectal epithelial cells	unknown
CTB-1	Lymphatic system	B-lymphocyte	unknown
CTV-1	Blood	T-lymphocyte	Suspension
CW-2	Colon	Colorectal epithelial cells	Adherent
CWR-R1	Prostate	Prostate epithelial cell	Adherent
D. hansenii	unknown	unknown	Neither
D1-1	unknown	Induced pluripotent stem cell	unknown
D-245MG	Brain	Astrocyte	unknown
D-247MG	Brain	Astrocyte	unknown
D-263MG	Brain	Astrocyte	unknown
D283 Med	Brain (medulla)	Brain cells	Suspension
D-336MG	Brain	Astrocyte	unknown
D341 Med	Brain (medulla)	Brain cells	Suspension
D-392MG	Brain	Astrocyte	unknown
D-423MG	Brain	Astrocyte	unknown
D-502MG	Brain	Astrocyte	unknown
D-542MG	Brain	Astrocyte	unknown
D-566MG	Brain	Astrocyte	unknown
DAN-G	Pancreas	Pancreatic epithelial cells	Adherent
Daoy	Brain	Brain cells	Adherent
Daudi	Lymphatic system	B-lymphoblast	Suspension
DB	Lymphatic system	B-lymphoblast	Suspension
DBTRG-05MG	Brain	Astrocyte	Adherent
DEL	Lymphatic system	Lymphocytes	Suspension
Dendritic cells	Immune system	Dendritic cell	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
DERL-2	Lymphatic system	T-lymphocyte	Suspension
DERL-7	Lymphatic system	T-lymphocyte	Suspension
Detroit 562	Pharynx	Squamous cells	Adherent
DF	Central nervous system	Astrocyte	unknown
DG-75	Lymphatic system	B lymphocyte	Suspension
DH82	unknown	Macrophage	Adherent
DI TNC1 astrocytes	Brain	Fibroblast	Adherent
Dictyostelium	unknown	unknown	Neither
Dictyostelium AX3	unknown	unknown	Neither
DIFI	Colon	Colorectal epithelial cells	unknown
DIFI R10 pool 1	Intestine	unknown	unknown
DIFI R10 pool 2	Intestine	unknown	unknown
DJM-1	Skin	Squamous cells	Adherent
DK-MG	Brain	Astrocyte	Adherent
DKO	Colon	Colorectal epithelial cells	unknown
DLD-1	Colon	Colorectal epithelial cells	Adherent
DMS 114	Lung	Lung cells	Adherent
DMS 153	Lung	Lung cells	Adherent
DMS 273	Lung	Lung cells	Adherent
DMS 53	Lung	Lung cells	Adherent
DMS 79	Lung	Lung cells	Suspension
DMS454	Lung	Lung cells	Adherent
DND-41	Blood	T-lymphocytes	Suspension
DOHH-2	Lymphoma	B-lymphocytes	Suspension
DOK	Head & Neck	Keratinocyte	Adherent
DoTc2 4510	Cervix	Cervical epithelial cells	Adherent
DRG Dorsal root ganglia neurons	Central nervous system	Dorsal root ganglion	unknown
DSH1	Urinary Tract	Bladder epithelial cells	unknown
DU145	Prostate	Prostate epithelial cell	Adherent
DU175	unknown	unknown	Adherent
DU4475	Breast	Mammary epithelial cells	Suspension
Dunaliella salina	unknown	Algae	Neither
Dunaliella sp.	unknown	Algae	Neither
Dunaliella tertiolecta	unknown	Algae	Neither
DV-90	Lung	Lung cells	Adherent
E. coli	unknown	Bacterial cells	Neither
E. faecalis	unknown	Bacterial cells	Neither
E. histolytica	unknown	Unicellular protist	Neither
E006AA	Prostate	Prostate epithelial cell	Adherent
E006AA-HT	Prostate	Prostate epithelial cell	Adherent
EB	Blood	Lymphoblast	Suspension
EB1	Lymphatic system	Lymphoblast	Suspension

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
EB2	Lymphatic system	B lymphocyte	Suspension
EB-3	Lymphatic system	B lymphocyte	Suspension
EBC-1	Lung	Squamous cell (lung)	Adherent
EBS	unknown	unknown	unknown
ECC1	Uterus	Uterine endometrial cells	unknown
ECC10	Stomach	Stomach cells	unknown
ECC12	Stomach	Stomach cells	unknown
ECC4	unknown	unknown	unknown
EC-GI-10	Esophagus	Esophagus epithelial cells	Adherent
Ede	unknown	unknown	unknown
EFC-1	Embryo	Embryonic stem cell	unknown
EFE-184	Uterus	Uterine endometrial cells	Adherent
EFM-19	Mammary gland	Mammary epithelial cells	Adherent
EFM-192A	Mammary gland	Mammary epithelial cells	Adherent
EFM-192B	Mammary gland	Mammary epithelial cells	Adherent
EFM-192C	Mammary gland	Mammary epithelial cells	Adherent
EFO-21	Ovary	Ovarian cells	Adherent
EFO-27	Ovary	Ovarian cells	Adherent
EGI-1	Biliary Tract	Bile ductal epithelial cells	Adherent
EHEB	Blood	B-lymphocyte	Suspension
EJ138	Urinary Tract	Bladder epithelial cells	Adherent
EJ-30 clone B3	Urinary Tract	Bladder epithelial cells	Adherent
EJM	Peripheral blood	Plasma cell	Adherent and suspension
EKVX	Lung	Lung cells	Adherent
EL16.7	unknown	unknown	unknown
EL-4	Thymus	T-lymphocyte	Suspension
ELT3	Uterus (rat)	Uterine fibroid cells	unknown
EM-2	Peripheral blood	Leukocyte	Suspension
embryonic stem cells mouse	Embryo	Embryonic stem cell	unknown
embryonic stem cells human	Embryo	Embryonic stem cell	unknown
EMSC	unknown	unknown	unknown
EN	Uterus	Uterine endometrial cells	Adherent
ENDOTHELIAL CELLS HUMAN PRIMARY	unknown	Endothelial cells	unknown
Enterococcus faecium	unknown	Bacterial cells	Neither
EoL-1-cell	Blood	Leukocyte	Suspension
Eos-HL-60	Blood	Myeloid cells	Suspension
Eosinophil cells	Blood	Eosinophil	unknown
EPITHELIAL CANCER CELLS	unknown	unknown	unknown
EPLC-272H	Lung	Lung squamous cells	Adherent
Erythrocytes	Blood	Erythrocytes (red blood cells)	unknown
Erythroid Progenitor Cells	Peripheral blood	Erythroid Progenitor Cells	unknown
ES1	Bone	Ewing sarcoma cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
ES-2	Ovary	ovarian clear cell	Adherent
ES3	Bone	Ewing sarcoma cells	unknown
ES4	Bone	Ewing sarcoma cells	unknown
ES5	Bone	Ewing sarcoma cells	unknown
ES6	Bone	Ewing sarcoma cells	unknown
ES7	Bone	Ewing sarcoma cells	unknown
ES8	Bone	Ewing sarcoma cells	unknown
ESC	unknown	unknown	unknown
ESCCTE9	Esophagus	Esophageal squamous cells	Adherent
esophageal epithelial cells primary	Esophagus	Esophageal epithelial cells	unknown
ESS-1	Uterus	Uterine endometrial cells	Adherent
ETK-1	Liver	Liver cells	unknown
EW 7476	Bone	Osteocytes	unknown
EW-1	Bone	Ewing sarcoma cells	unknown
EW-11	Bone	Ewing sarcoma cells	unknown
EW-12	Bone	Ewing sarcoma cells	unknown
EW-13	Bone	Ewing sarcoma cells	unknown
EW-16	Bone	Ewing sarcoma cells	unknown
EW-18	Bone	Ewing sarcoma cells	unknown
EW-22	Bone	Ewing sarcoma cells	unknown
EW-24	Bone	Ewing sarcoma cells	unknown
EW-3	Bone	Ewing sarcoma cells	unknown
EW-7	Bone	Ewing sarcoma cells	unknown
Exophiala dermatitidis	unknown	Fungus	unknown
Expi293	unknown	unknown	unknown
Expi293F	unknown	unknown	unknown
F480+ macrophages	unknown	unknown	unknown
F98	Brain	Astrocyte	Adherent
Farage	Lymphatic system	B-lymphocyte	Suspension
Fetal Osteoblasts	Bone (fetus)	Osteoblast	unknown
FHs 74 Int	Intestine small	Intestinal epithelial cell fetus	Adherent
fibroblasts Primary	unknown	unknown	unknown
FIE	unknown	unknown	unknown
FLPLN293	Kidney (fetus)	Fetal kidney cell	Adherent
FLYA13	unknown	Fibrosarcoma cell	Adherent
FMC FELINE Mammary carcinoma	Mammary gland	Mammary epithelial cells	unknown
Foal Leukoctyes	Blood	Leukocyte	unknown
FoxD1-derived cells	unknown	unknown	unknown
fR2	Breast	Mammary epithelial cells	Adherent
FTC-133	Thyroid	Thyroid cells	Adherent
FTC-238	Thyroid	Thyroid cells	Adherent
FU97	Stomach	Stomach cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
FU-OV-1	Ovary	Ovary cell	Adherent
FXR-UAS-bla HEK 293T	Kidney (fetus)	Fetal kidney cell	unknown
G-292 Clone A141B1	Bone	Osteocytes	Adherent
G-361	Skin	Melanocytes	Adherent
G-401	Kidney	Kidney cells	Adherent
G-402	Kidney	Kidney cells	Adherent
G-415	Biliary Tract	Biliary cell	Adherent
G-418	unknown	unknown	unknown
GA-10	Lymphatic system	B-lymphocyte	Suspension
GA-10 (Clone 20)	Lymphatic system	B-lymphocyte	Suspension
GA-10 (Clone 4)	Lymphatic system	B-lymphocyte	Suspension
GAK	Skin	Melanocytes	Adherent
GAMG	Brain	Astrocyte	Adherent
Gastric mucosal small endocrine cells	Stomach	Gastric mucosal small endocrine cells	unknown
GB-1	Brain	Astrocyte	Adherent
GBD1	Biliary Tract	Biliary cell	unknown
GBM	Brain	Astrocyte	unknown
GCIY	Stomach	Stomach cells	Adherent
GCT	Miscellaneous	Lung	Adherent
GDM-1	Blood	monoblast	Suspension
Geobacillus	unknown	Bacterial cells	Neither
GI-1	Brain	Glial cell	Adherent
GI-ME-N	Kidney	Adrenal cell	Adherent
GL-261	Brain	Astrocyte	Adherent
Gli60I-BM	Brain	Astrocyte	Adherent
Gli60I-P	Brain	Astrocyte	Adherent
Gli60-MSH6-BM	Brain	Astrocyte	Adherent
Gli60-Vect-BM	Brain	Astrocyte	Adherent
Gli60-Vect-P	Brain	Astrocyte	Adherent
Glioblastoma stem-like cells	Brain	Astrocyte	Adherent
GMCs rat	unknown	unknown	unknown
G-MEL	Skin	Melanocytes	unknown
GMS-10	Brain	Astrocyte	Adherent
GOS-3	Brain	Astrocyte	Adherent
GOTO	Nervous system	Neuroblast	unknown
GOTO.P3	Kidney	Adrenal cell	Suspension
Gp2D	Colon	Colorectal epithelial cells	Adherent
GP5d	Colon	Colorectal epithelial cells	Adherent
GPLM	unknown	unknown	unknown
GRANTA-519	Lymphatic system	Lymph cells	Suspension
Granulocytes	Blood	Granulocytes	Suspension
GRPr	unknown	unknown	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
GR-ST	Blood	Lymphocyte	Suspension
GT3TKB	Lung	Lung squamous cells	unknown
GTL-16	Stomach	Stomach cells	Adherent
H1299	Lung	Lung cells	Adherent
H1650	Lung	Lung cells	Adherent
H1975	Lung	Lung cells	Adherent
H2052	Pleura	Mesothelial cells	Adherent
H226	Pleura	Mesothelial cells	Adherent
H2369	Pleura	Mesothelial cells	unknown
H2373	Pleura	Mesothelial cells	unknown
H2461	Pleura	Mesothelial cells	unknown
H2591	Pleura	Mesothelial cells	unknown
H2595	Pleura	Mesothelial cells	unknown
H2596	Pleura	Mesothelial cells	unknown
H2691	Pleura	Mesothelial cells	unknown
H2722	Pleura	Mesothelial cells	unknown
H2731	Pleura	Mesothelial cells	unknown
H2795	Pleura	Mesothelial cells	unknown
H28	Pleura	Mesothelial cells	Adherent
H2803	Pleura	Mesothelial cells	unknown
H2804	Pleura	Mesothelial cells	unknown
H2810	Pleura	Mesothelial cells	unknown
H2818	Pleura	Mesothelial cells	unknown
H2869	Pleura	Mesothelial cells	unknown
H290	Pleura	Mesothelial cells	unknown
H292	Lung	Lung cells	Adherent
H3118	Head & Neck	Parotid gland cells	unknown
H3255	Lung	Lung cell	unknown
H33HJ-JA1	Blood	unknown	Suspension
H4	Brain	Astrocyte	Adherent
H460	unknown	Hybridoma	unknown
H513	Lung	Lung squamous cells	unknown
H69	Lung	Lung cells	Suspension
H69AR	Lung	Lung cells	Adherent
H69V	Lung	Lung cells	Adherent
H9	Lymphatic system	T-lymphocyte (cutaneous)	Suspension
H929	Lymphatic system	B-lymphocyte	Suspension
H9c2	Myocardium	Myeloblast	Adherent
HA7-RCC	Kidney	Kidney cells	unknown
HaCaT	Skin	Skin cell	Adherent
HAL-01	Leukemia	B-lymphoblasts	Suspension
HARA	Lung	Lung squamous cells	unknown
hASCs	Adipose tissue	Adipocytes	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
HBE135-E6E7	Lung	Lung cells	Adherent
HBE4-E6/E7	Lung	Lung cells	Adherent
HBE4-E6/E7-C1	Lung	Lung cells	Adherent
HBG3	Lung	unknown	unknown
HBL100	Breast	Mammary epithelial cells	Adherent
Hbmob	Peripheral blood	Human Peripheral blood osteoblasts	Adherent
HBR cells	unknown	unknown	unknown
HC-1	Blood	Leukocyte	Suspension
HC11	Mammary gland	Mammary epithelial cells	Adherent
HCC 2998	Colon	Colorectal epithelial cells	unknown
HCC 827	Colon	Colorectal epithelial cells	Adherent
HCC1008	Mammary gland	Mammary epithelial cells	Adherent
HCC1143	Mammary gland	Mammary epithelial cells	Adherent
HCC1187	Mammary gland	Mammary epithelial cells	Adherent and suspension
HCC1187 BL	Blood	B-lymphoblast	Suspension
HCC1395	Mammary gland	Mammary epithelial cells	Adherent
HCC1419	Mammary gland	Mammary epithelial cells	Adherent
HCC1428	Mammary gland	Mammary epithelial cells	Adherent
HCC-15	Lung	Lung cells	Adherent
HCC1500	Mammary gland	Mammary epithelial cells	Adherent
HCC1569	Mammary gland	Mammary epithelial cells	Adherent
HCC1599 BL	Blood	B-lymphocytes	Suspension
HCC1806	Mammary gland	Mammary epithelial cells	Adherent
HCC1937	Mammary gland	Mammary epithelial cells	Adherent
HCC1954	Mammary gland	Mammary epithelial cells	Adherent
HCC1954 BL	Blood	B-lymphoblast	Suspension
HCC202	Mammary gland	Mammary epithelial cells	Adherent
HCC2157	Mammary gland	Mammary epithelial cells	Adherent
HCC2157 BL	Blood	B-lymphoblast	Suspension
HCC2218	Mammary gland	Mammary epithelial cells	Suspension
HCC2218 BL	Blood	B-lymphoblast	Suspension
HCC2279	Lung	Lung cells	unknown
HCC2998	Colon	Colorectal epithelial cells	unknown
HCC-33	Lung	Lung cells	Suspension
HCC-366	Lung	Lung cells	Adherent
HCC38	Mammary gland	Mammary epithelial cells	Adherent
HCC4006	Lung	Lung cells	Adherent
HCC-44	Lung	Lung cells	Adherent
HCC-56	Colon	Colorectal epithelial cells	unknown
HCC70	Mammary gland	Mammary epithelial cells	Adherent
HCC-78	Lung	Lung cells	Adherent
HCC-827	Lung	Lung cells	Adherent
HCE7	Esophagus	Esophagus epithelial cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
HCE-T	Cornea	Corneal epithelial cells	Adherent
HCI-1173	unknown	unknown	unknown
HCI-128	unknown	unknown	unknown
HCI-1299	unknown	unknown	unknown
HCI-1373	unknown	unknown	unknown
HCI-1725	unknown	unknown	unknown
HCI-1944	unknown	unknown	unknown
HCI2077	unknown	unknown	unknown
HCI-2087	unknown	unknown	unknown
HCI-211	unknown	unknown	unknown
HCI-2122	unknown	unknown	unknown
HCI-23	unknown	unknown	unknown
HCI-3122	unknown	unknown	unknown
HCI-322	unknown	unknown	unknown
HCI-358	unknown	unknown	unknown
HCI-460	unknown	unknown	unknown
HCI-592	unknown	unknown	unknown
HCI-620	unknown	unknown	unknown
HCI-678	unknown	unknown	unknown
HCI-69	unknown	unknown	unknown
HCI-720	unknown	unknown	unknown
HCI-77	unknown	unknown	unknown
HCI-82	unknown	unknown	unknown
HCI-820	unknown	unknown	unknown
HCI-854	unknown	unknown	unknown
HCT116	Colon	Colorectal epithelial cells	Adherent
HCT-15	Colon	Colorectal epithelial cells	Adherent
HCT-8	Colon	Colorectal epithelial cells	Adherent
HDFs	Skin	Dermal fibroblasts	unknown
HDLM-2	Lymphatic system	Reed-sternberg cells	Suspension
HD-MY-Z	Lymphatic system	Reed-sternberg cells	Adherent and suspension
HDQ-P1	Mammary gland	Mammary epithelial cells	Adherent
HE8935	unknown	unknown	unknown
HEART-ATTAC transgenic primary cardiomyocytes	unknown	unknown	unknown
HEC-1	Uterus	Uterine endometrial cells	unknown
HECV	Umbilical cord	Umbilical cord cells	Adherent
HEIOC1	unknown	unknown	unknown
HEI-OC1 murine-derived auditory cells	unknown	unknown	unknown
HEK293	Kidney	Debated	Adherent
HEK293-A	Kidney	Debated	Adherent
HEK293-F	Kidney	Debated	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
HEK293FT	Kidney	Debated	Adherent
HEK293T	Kidney	Debated	Adherent
HEL	Lymphatic system	Lymphocyte	Suspension
HEL 92.1.7	Lymphatic system	Lymphocyte	Suspension
HeLa	Cervix	Cervical epithelial cells	Adherent
HeLa 229	Cervix	Cervical epithelial cells	Adherent and suspension
HeLa AG	Cervix	Cervical epithelial cells	unknown
HeLa Bu25	Cervix	Cervical epithelial cells	Adherent
HeLa K11	Cervix	Cervical epithelial cells	unknown
HeLa S3	Cervix	Cervical epithelial cells	Adherent
HeLa S3 (sc)	Cervix	Cervical epithelial cells	unknown
HeLa TG	Cervix	Cervical epithelial cells	unknown
HeLa TG Cap	Cervix	Cervical epithelial cells	unknown
HeLa.P3(S)	Cervix	Cervical epithelial cells	unknown
Hela-USU	Cervix	unknown	unknown
HEMANGIOSARCOMA CELLS	Cervix	unknown	unknown
Hematopoietic stem cells	unknown	unknown	unknown
H-EMC-SS	Cartilage	Cartilage cell	Adherent
hemocytes	Blood	Blood cell	unknown
Hep 3B2.1-7	Liver	Hepatocyte	Adherent
Hep G2	Liver	Hepatocyte	Adherent
Hep-2 laryngeal carcinoma	Cervix	Cervical epithelial cell	Adherent
Hepa1-6	Liver	Hepatocytes	Adherent
hepatic progenitor cells	Liver	Hepatic progenitor cells	unknown
Hepatocytes	Liver	Hepatocytes	unknown
Hep-G2	Liver	Hepatocytes	Adherent
Hep-G2/C3A	Liver	Hepatocytes	Adherent
hES	Cervix	Cervical cells	unknown
hESC HUES6 cells	unknown	unknown	unknown
hESCs-H7	Embryo	Embryonic stem cell (human)	Adherent
HEY	Ovary	Ovarian cells	unknown
HFL-1	Lung (fetus)	Lung cells (fetus)	Adherent
hFOB	Bone (fetus)	Osteoblast (fetus)	Adherent
HGC-27	Stomach	Stomach cells	Adherent
HH	Peripheral blood	T lymphocyte	Suspension
HIC-1466	unknown	unknown	unknown
HISTIOCYTIC CELLS	Peripheral blood	Histiocytic cells	unknown
HK-2	Kidney	Kidney cells	Adherent
HL	Lung	Lung fibroblasts	Adherent
HL1-hT1	Heart	Cardiomyocytes	Adherent
HL-60	Peripheral blood	Promyeoblast	Suspension
HL-60 clone 15	Kidney	Lymphoblast	Suspension
HL60(S)	Peripheral blood	Promyeoblast	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
HL-60/MX1	Peripheral blood	Promyeoblast	Suspension
HL-60/MX2	Peripheral blood	Promyeoblast	Suspension
HL60RG	Peripheral blood	Promyeoblast	Suspension
HLCL-1	Peripheral blood	Lymphocyte	Suspension
HLE	Liver	Hepatocyte	Adherent
HLF	Liver	Hepatocyte	Adherent
HMCB	Skin	Melanocyte	Adherent
HMEC-1	Skin	dermal microvascular endothelium	Adherent
hMSC	unknown	Mesenchymal stem cells human	unknown
HMVEC-Lly	Lung	Lung Microvascular Endothelial Cells	unknown
HMVII	Skin (vaginal epithelium)	Vaginal epithelial (skin) cells	Adherent
HN	Mouth	Oral squamous cells	Adherent
hNP (Neural Progenitor derived from hES)	Central nervous system	Shwann cells	unknown
HO-1-N-1	Mouth	Oral squamous cells	Adherent
HO-1-u-1	Mouth	Oral squamous cells	Adherent
HO8910	Ovary	Ovarian cells	unknown
HOP-62	Lung	Lung cells	unknown
HOP-92	Lung	Lung non-small cells	unknown
HOS	Bone	Osteoblast	Adherent
HPAC	Pancreas	Pancreatic cells	Adherent
HPAF-II	Pancreas	Pancreatic ductal cells	Adherent
Hpsc	unknown	Pluripotent stem cells human	unknown
HRT-18	Colon	Colorectal epithelial cells	Adherent
Hs 1.Int	Colon (fetus)	Colon fetal fibroblasts	Adherent
Hs 14.T	Bone	Osteoblast	Adherent
Hs 172.T	Bladder	Bladder epithelial cells	unknown
Hs 195.T	Bladder	Bladder epithelial cells	unknown
Hs 219.T	Rectum	Rectal epithelial cells	unknown
Hs 228.T	Bladder	Bladder fibroblasts	Adherent
Hs 257.T	Colon	Colorectal epithelial cells	Adherent
Hs 313.T	Lymphatic system	Lymphocyte	Adherent
Hs 387.T	Bone	Osteoblast	Adherent and suspension
Hs 417.Lu	? (fetus)	unknown	unknown
Hs 445	Lymphatic system	B-lymphocyte	Suspension
Hs 505.T	unknown	Lymphocyte	Adherent
Hs 518.T	unknown	Lymphocyte	unknown
Hs 564(E).Mg	Breast	Mammary epithelial cells	Adherent
Hs 573.T	Lung	Lung epithelial cells	Adherent
Hs 578Bst	Breast	Mammary epithelial cells	Adherent
HS 578T	Breast	Mammary epithelial cells	Adherent
Hs 579.Mg	Breast	Mammary epithelial cells	Adherent
Hs 587.Int	Colon	Colorectal epithelial cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Hs 588.T	Cervix	Cervical epithelial cells	Adherent
Hs 600.T	Skin	Melanocyte	Adherent
Hs 602	Lymphatic system	Lymphocyte	Suspension
Hs 604.T	Lymphatic system	Lymphocyte	Adherent
Hs 611.T	Lymphatic system	Lymphocyte	Adherent and suspension
Hs 616.T	Lymphatic system	Lymphocyte	Adherent
Hs 633T	unknown	Fibroblast	Adherent
Hs 683	Brain	Oligodendrocyte	Adherent
Hs 704.T	Bone	Osteoblast	Adherent
Hs 707(A).T	Bone	Osteoblast	Adherent
Hs 735.T	Bone	Osteoblast	unknown
Hs 737.T	Bone	Osteoblast	Adherent
Hs 738.St/Int	Stomach and intestine (fetus)	Fibroblast (fetus)	Adherent
Hs 746T	Stomach	Stomach epithelial cells	Adherent
Hs 751.T	Lymphatic system	Lymphocyte	unknown
Hs 755(B).T	Bone	Osteoblast	Adherent
Hs 766T	Pancreas	Pancreatic cells	Adherent
Hs 777.T	Lymphatic system	Lymphocyte	Suspension
Hs 781.T	Bone	Osteoblast	Adherent
Hs 792(B).T	Bone	Osteoblast	unknown
Hs 805.T	Bone	Osteoblast	unknown
Hs 811.T	Bone	Osteoblast	unknown
Hs 819.T	Cartilage	Cartilage cell	Adherent
Hs 834.T	Skin	Melanocyte	Adherent
Hs 839.T	Skin	Melanocyte	Adherent
Hs 845.T	Bone	Osteoblast	unknown
Hs 852.T	Skin	Melanocyte	Adherent
Hs 860.T	Bone	Osteoblast	Adherent
Hs 863.T	Bone	Osteoblast	Adherent
Hs 865.Sk	Skin	Fibroblast	Adherent
Hs 866.T	Bone	Osteoblast	unknown
Hs 870.T	Bone	Osteoblast	Adherent
Hs 888.T	Bone	Osteoblast	Adherent
Hs 889.T	Bone	Osteoblast	unknown
Hs 890.T	Bone	Osteoblast	Adherent
Hs 891.T	Kidney	Renal cell	Adherent
Hs 894(A).T	Bone	Osteoblast	Adherent
Hs 894(B).T	Bone	Osteoblast	unknown
Hs 894(C).T	Bone	Osteoblast	unknown
Hs 894(D).T	Bone	Osteoblast	unknown
Hs 894(E).Lu	unknown	unknown	unknown
Hs 895.T	Skin	Melanocyte	Adherent
Hs 899(A).T	Bone	Osteoblast	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Hs 899(B).T	Bone	Osteoblast	unknown
Hs 899(C).T	Bone	Osteoblast	Adherent
Hs 899(D).T	Bone	Osteoblast	Adherent
Hs 906(A).T	Skin	Melanocyte	Adherent
Hs 908.T	Skin	Melanocyte	unknown
Hs 926.T	Kidney	Renal rhabdomyosarcoma cell	Adherent
Hs 936.T	Skin	Melanocyte	unknown
Hs 939.T	Skin	Melanocyte	unknown
Hs 940.T	Skin	Melanocyte	Adherent
Hs 944.T	Skin	Melanocyte	Adherent
HSC-1	Skin	Squamous skin cell	Adherent
HSC-2	Mouth	Squamous cell oral	Adherent
HSC-3	Tongue	Squamous cell tongue	Adherent and suspension
HSC-39	Stomach	Signet ring cell	unknown
HSC-4	Tongue	Squamous cell tongue	Suspension
HSFs	unknown	unknown	unknown
HSPCs	Blood	Hematopoietic stem cell	unknown
HS-Sultan	Lymphatic system	B-lymphocyte	Suspension
HT	Lymphatic system	B-lymphoblast	Suspension
HT 1080	Connective tissue	Fibroblast	Adherent
HT 1080	Connective tissue	Fibroblast	Adherent
HT 1376	Bladder	Bladder epithelial cells	Adherent
HT 1417	Lymphatic system	Lymphocyte	unknown
HT 728.T	Bone	Osteoblast	unknown
HT-1197	Bladder	Bladder epithelial cells	Adherent
HT-144	Skin	Melanocyte	Adherent
HT-29	Colon	Colorectal epithelial cells	Adherent
HT-3	Cervix	Cervical epithelial cells	Adherent
HT55	Colon	Colorectal epithelial cells	Adherent
HTC-C3	Thyroid	Thyroid epithelial cells	Adherent
hTCEpi	Cornea	Corneal epithelial cells	Adherent
Htert- hpne	Pancreas	Pancreas intermediary cells formed during acinar-to-ductal metaplasia	Adherent
hTSC	unknown	unknown	unknown
hUC	Bladder	Human Urothelial Cells	Adherent
HuCCT-1	Liver bile duct	Liver bile duct epithelial cells	Adherent
huH-1	Liver	Hepatocytes	unknown
HuH28	Liver bile duct	Liver bile duct epithelial cells	Adherent
HUH-6 Clone 5	Liver/gallbladder	Hepatoblasts	Adherent
HuH-7	Liver	Hepatocytes	Adherent
Human Astrocytes Primary	Brain	Astrocytes	unknown
Human Melanoma	Skin	Melanocyte	unknown
human monocytes	Blood	Monocytes	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
human T lymphocytes	Blood	T-lymphocytes	unknown
HuO-3N1	Bone	Osteoblast	Adherent
HuO9	Bone	Osteoblast	Adherent
HuO9N2	Bone	Osteoblast	Adherent
HUP-T3	Pancreas	Pancreas epithelial cells	Adherent
HUP-T4	Pancreas	Pancreas epithelial cells	Adherent
HuT 102	Lymphatic system	T-lymphocytes	Suspension
HuT 78	Lymphatic system	T-lymphocytes	Suspension
HUTU-80	Duodenum	Duodenal epithelial cells	Adherent
HUVEC	Umbilical vein/ vascular endothelium	Umbilical vascular epithelial cells	Adherent
I 2.1	Blood	T-lymphocytes	Suspension
I 9.2	Blood	T-lymphocytes	Suspension
IA-LM	Lung	Lung large-cells	Adherent
IEC-6	Small intestine	Intestinal epithelial cells	Adherent
IFT27RNAi Knockdown cells	unknown	unknown	unknown
IGR-1	Skin	Melanocyte	Adherent
IGR-37	Skin	Melanocyte	Adherent
IGR-39	Skin	Melanocyte	Adherent
IGROV-1	Ovary	Ovarian cells	Adherent
IHH-4	Thyroid	Thyroid epithelial cells	Adherent
IL-2	unknown	unknown	unknown
IM-9	Peripheral blood	B-lymphocyte	Suspension
IM-95	Stomach	Gastric epithelial cells	Adherent
IM-95m	Stomach	Gastric epithelial cells	Adherent
IMR-32	Brain	Neuroblast	Adherent
IMR-5	Nervous system	Neuroblast	Adherent
IMR-90	Lung (fetus)	Fibroblast	Adherent
IMR-90-40	Lung (fetus)	Fibroblast	Adherent
inflammatory breast cancer (IBC) spheroids	Breast	Mammary epithelial cells	unknown
INS113	unknown	unknown	unknown
INS1832	Pancreas	Pancreatic beta cells	unknown
INS-1E	Pancreas	Pancreatic beta cells	Adherent
IP-1	unknown	unknown	unknown
IPC-298	Skin	Melanocyte	Adherent
IPTP/98	unknown	unknown	unknown
iRin	Pancreas	Pancreatic beta cells	Adherent
Ishikawa	Uterus	Uterine endometrial cells	Adherent
IST-MES1	Mesothelium	Squamous epithelial cells	Adherent
IST-SL1	Lung	Lung small-cells	Adherent
IST-SL2	Lung	Lung small-cells	Suspension
ITO-II	Testicle	Testicular germ cells	unknown
ITSC	unknown	Induced totipotent stem cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
J.21.WT	unknown	unknown	unknown
J.RT3-T3.5	Blood	T-lymphocytes	Suspension
J1	Blastocyst	Embryonic stem cell	Colonies
J45.01	Peripheral blood	T-lymphocytes	Suspension
J82	Bladder	Bladder epithelial cells	Adherent
JAK2 V617F	unknown	unknown	unknown
JAR	Placenta	Trophoblasts	Adherent
JEG-3	Placenta	Trophoblasts	Adherent
JEKO-1	Peripheral blood	Lymphoblast	Suspension
JH10	unknown	Soil-dwelling ameoba	unknown
JHH-1	Liver	Hepatocytes	Adherent
JHH-2	Liver	Hepatocytes	Adherent
JHH-4	Liver	Hepatocytes	Adherent
JHH-6	Liver	Hepatocytes	Adherent
JHH-7	Liver	Hepatocytes	Adherent
JHU-011	Larynx	Laryngeal epithelial cells	unknown
JHU-012	Mouth	Squamous cells	unknown
JHU-013	Pharynx	Squamous cells	unknown
JHU-019	Prostate	Prostate cells	unknown
JHU-022	Mouth	Squamous cells	unknown
JHU-028	Lung	Lung cells	unknown
JHU-029	Larynx	Laryngeal epithelial cells	unknown
JIMT-1	Breast	Mammary epithelial cells	Adherent
Jiyoye	Lymphatic system	B-lymphocytes	Suspension
JJN-3	Peripheral blood	Plasma cells	Suspension
JKT-beta-del	Peripheral blood	T-lymphocytes	Suspension
JM1	Peripheral blood	B-lymphocytes	Suspension
JSC-1	Peripheral blood	Lymphocyte	unknown
JU547	unknown	Freshwater cyanobacteria	Free-floating
JURKAT	Blood	T-lymphocyte	Suspension
Jurkat E6	Blood	T-lymphocyte	Suspension
Jurkat E6-1	Blood	T-lymphocyte	Suspension
JURL-MK1	Bone marrow	Myeloid cells	Suspension
JURL-MK2	Bone marrow	Myeloid cells	Suspension
JVM-2	Lymphatic system	Lymphoblast	Suspension
JVM-3	Lymphatic system	B-lymphocytes	Suspension
K. flaccidum	unknown	Algae	unknown
K. lactis	unknown	Yeast	unknown
K. waltii	unknown	Yeast	unknown
K1	Thyroid	Thyroid epithelial cells	Adherent
K19	Skin	Melanocyte	unknown
K2	Skin	Melanocyte	unknown
K23	unknown	unknown	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
K25	unknown	Induced pluripotent stem cell	unknown
K4	Skin	Melanocyte	unknown
K5	Thyroid	Thyroid epithelial cells	unknown
K562	Peripheral blood	Lymphoblast	Suspension
K562/ADM	Peripheral blood	Lymphoblast	Suspension
K562-derived T cells	Peripheral blood	T-Lymphoblast	Suspension
K8	unknown	unknown	unknown
KALS-1	Brain	Glial cell	Adherent
KARPAS 299	Peripheral blood	T-lymphoblast	Suspension
KARPAS-1106P	Lymphatic system	B-lymphocyte	Suspension
KARPAS-231	Lymphatic system	B-lymphocyte	Suspension
KARPAS-299	Lymphatic system	T-lymphoblast	Suspension
KARPAS-422	Lymphatic system	B-lymphocyte	Suspension
KARPAS-45	Bone marrow	T-lymphocytes	Suspension
KARPAS-620	Peripheral blood	Plasma cells	Suspension
Kasumi-1	Peripheral blood	Myeloblast	Suspension
Kasumi-3	Peripheral blood	Myeloblast	Suspension
Kasumi-4	Peripheral blood	Myeloblast	Suspension
Kasumi-6	Peripheral blood	Myeloblast	Suspension
KATO II	Stomach	Signet ring cells	Adherent and suspension
KATO III	Stomach	Signet ring cells	Adherent and suspension
KB-3-1	Cervix	Cervical epithelial cells	Adherent
KB-ChR-8-5-11	Cervix	Cervical epithelial cells	Adherent
KCL-22	Peripheral blood	Myeloblast	Suspension
KE-37	Peripheral blood	T-lymphocyte	Suspension
KELLY	Nervous system	Neuroblast	Adherent
KG-1	Bone marrow	Myeloblast	Suspension
KG-1a	Bone marrow	Myeloblast	Suspension
KG-1-C	Brain	Mixed glioma	Adherent
KGN	Ovary	Ovarian granulocyte	Adherent
KHM-3S	Lung	Lung non-small cells	Suspension
KHOS/NP	Bone	Osteoblast	Adherent
KHOS-240S	Bone	Osteoblast	Adherent
KHOS-312H	Bone	Osteoblast	Adherent
Kidney	Kidney	Renal cells	unknown
KIDNEY EPITHELIAL CELLS FELINE	Kidney	Renal epithelial cells	unknown
KINGS-1	Brain	Astrocytes	Adherent
KIT-transduced IC2 cells	unknown	unknown	Adherent
KLE	Uterus	Uterine endometrial cells	Adherent
KM12	Colon	Colon epithelial cells	unknown
KMCH	unknown	unknown	unknown
KMCH-1	Liver and bile duct	Liver cells; bile duct epithelial cells	Adherent
KMH-2	Thyroid	Thyroid epithelial cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
KM-H2	Lymphatic system	Reed-sternberg cells	Suspension
KMOE-2	Blood	Erythroblasts	Suspension
KMRC-1	Kidney	Renal cells	Adherent
KMRC-2	Kidney	Renal cells	Adherent
KMRC-20	Kidney	Renal cells	Adherent
KMRC-3	Kidney	Renal cells	Adherent
KMRM-M1	Kidney	Renal cells	Adherent
KMS-12-BM	Peripheral blood	Myeloblast	Suspension
KMS-12-PE	Peripheral blood	Myeloblast	Suspension
KNS-42	Brain	Glial cell	Adherent
KNS-62	Bronchus	Bronchial squamous cell	Adherent
KNS-81-FD	Brain	Glial cell	Adherent
KO52	Peripheral blood	Myeloblast	Suspension
KON	Mouth	Oral squamous cells	Adherent
KOPN-8	Peripheral blood	B-lymphoblast	Suspension
KOSC-2	Head and neck	Squamous cells	Adherent
KP-1N	Pancreas	Pancreatic ductal cells	Adherent
KP-1NL	Pancreas	Pancreatic ductal cells	Adherent
KP-2	Pancreas	Pancreatic ductal cells	Adherent
KP-3	Pancreas	Pancreatic ductal cells	Adherent
KP-3L	Pancreas	Pancreatic ductal cells	Adherent
KP-4	Pancreas	Pancreatic ductal cells	Adherent
KPL-1	Breast	Mammary ductal epithelial cells	Adherent
KP-N-RT-BM-1	Nervous system	Neuroblast	Adherent
KP-N-YN	Nervous system	Neuroblast	Adherent
KP-N-YS	Nervous system	Neuroblast	Adherent
KRJ-I	Digestive system	unknown	Adherent
KS-1	unknown	unknown	unknown
KU-19-19	Bladder	Bladder epithelial cells	Adherent
KU812	Peripheral blood	Basophil	Suspension
KU812E	Peripheral blood	Myeloblast	Suspension
KURAMOCHI	Ovary	Ovarian serous epithelial cell	Adherent
KY821	Peripheral blood	Myeloblast	Suspension
KY821A3	Peripheral blood	Myeloblast	Suspension
KYM-1	Neck	Somatic cell (not otherwise specified)	Adherent
KYSE-140	Esophogus	Esophageal squamous cells	Adherent
KYSE-150	Esophogus	Esophageal squamous cells	Adherent
KYSE-180	Esophogus	Esophageal squamous cells	Adherent
KYSE-220	Esophogus	Esophageal squamous cells	Adherent
KYSE-270	Esophogus	Esophageal squamous cells	Adherent
KYSE-30	Esophogus	Esophageal squamous cells	Adherent
KYSE-410	Esophogus	Esophageal squamous cells	Adherent
KYSE-450	Esophogus	Esophageal squamous cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
KYSE-50	Esophagus	Esophageal squamous cells	Adherent
KYSE-510	Esophagus	Esophageal squamous cells	Adherent
KYSE-520	Esophagus	Esophageal squamous cells	Adherent
KYSE-70	Esophagus	Esophageal squamous cells	Adherent
L-1236	Lymphatic system	Reed-sternberg cells	Adherent and suspension
L-363	Peripheral blood	Myeloblast	Suspension
L-428	Lymphatic system	Reed-sternberg cells	Suspension
L-540	Lymphatic system	Reed-sternberg cells	Suspension
L929	Connective tissue	Fibroblasts	Adherent
LAMA-84	Peripheral blood	Myeloblast	Suspension
LAMA-87	Peripheral blood	Myeloblast	Suspension
LAN-5	Nervous system	Neuroblast	Adherent
LAN-6	Nervous system	Neuroblast	Adherent
LB1047-RCC	Kidney	Renal cells	Adherent
LB2241-RCC	Kidney	Renal cells	Adherent
LB2518-MEL	Skin	Melanocyte	unknown
LB373-MEL-D	Skin	Melanocyte	unknown
LB647-SCLC	Lung	Lung small cells	unknown
LB771-HNC	Head and neck	Head and neck squamous cells	unknown
LB831-BLC	Bladder	Bladder epithelial cells	unknown
LB996-RCC	Kidney	Kidney cells	unknown
LBCL11832	Lymphatic system	B-lymphoblast	unknown
LC-1 sq	Lung	Lung squamous cells	Adherent
LC-1F	Lung	Lung squamous cells	Suspension
LC-2-ad	Lung	Lung cells	Adherent
LC4-1	Peripheral blood	B-lymphoblast	Suspension
LCL	Peripheral blood	Lymphoblast	unknown
LCLC-103H	Lung	Lung large cells	Adherent
LCLC-97TM1	Lung	Lung large cells	Adherent
LE/6	Liver	Precursor epithelial (oval) cells from liver	unknown
Leukocytes	Blood	Leukocyte	Suspension
LI90	Blood vessels	Blood vessel endothelial cell	Adherent
Liposarc3269	unknown	unknown	unknown
LK-2	Lung	Lung squamous cells	Adherent
LLCPK1	Kidney	Kidney epithelial cell	Adherent
LM7	Bone	Osteoblast	Adherent
LN	Brain	Glial cell	Adherent
LN-18	Brain	Glial cell	Adherent
LN-229	Brain	Glial cell	Adherent
LN-405	Brain	Astrocyte	Adherent
LNCaP	Prostate	Prostate cells	Adherent
LNCaP clone FGC	Prostate	Prostate cells	Adherent
LNCaP-FGC	Prostate	Prostate cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
LNZTA3WT11	Brain	Astrocyte	Adherent
LNZTA3WT4	Brain	Astrocyte	Adherent
Loucy	Peripheral blood	T-lymphocyte	Suspension
LOU-NH91	Lung	Lung squamous cells	Adherent
LoVo	Colon	Colon epithelial cells	Adherent
LOX-IMV1	Skin	Melanocyte	Adherent
LOXIMVI	Skin	Melanocyte	Adherent
LP-1	Peripheral blood	Plasma cell	Suspension
LPs	unknown	unknown	unknown
LS-1034	Cecum	Cecum epithelial cells	Adherent
LS-123	Colon	Colon epithelial cells	Adherent
LS174T	Colon	Colon epithelial cells	Adherent
LS180	Colon	Colon epithelial cells	Adherent
LS-411N	Colon	Colon epithelial cells	Adherent
LS-513	Cecum	Cecum epithelial cells	Adherent
Lu1205	Skin	Melanocyte	Adherent
LU-134-A	Lung	Lung small cells	Adherent
Lu-134-A-H	Lung	Lung small cells	Adherent
Lu-134-B	Lung	Lung small cells	Suspension
Lu-135	Lung	Lung small cells	Suspension
LU-139	Lung	Lung small cells	Suspension
LU-165	Lung	Lung small cells	Suspension
LU65	Lung	Lung giant cells	Adherent
LU65A	Lung	Lung giant cells	Adherent
LU65B	Lung	Lung giant cells	Adherent
LU65C	Lung	Lung giant cells	Adherent
LU99A	Lung	Lung giant cells	Adherent
LU99B	Lung	Lung giant cells	unknown
LU99C	Lung	Lung giant cells	unknown
LUDLU-1	Lung	Lung squamous cells	Adherent
Luise	unknown	unknown	unknown
Lung Leukocytes	Lung	Leukocyte	unknown
Lung mononuclear cells	Lung	Lung mononuclear cells	unknown
LX-2	Liver	Liver cells	unknown
LXF-289	Lung	Lung cells	Adherent
Lymphoblast	Lymphatic system	Lymphoblast	unknown
Lymphocytes	Lymphatic system	Lymphocyte	unknown
M. fructicola	unknown	Fruit-dwelling fungus	unknown
M. tuberculosis	unknown	Pathogenic bacteria	unknown
M059J	Brain	Glial cell	Adherent
M059K	Brain	Glial cell	Adherent
M-07e	Peripheral blood	Myeloblast	Suspension
M-14	Skin	Melanocyte	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Macrophage	Blood	Macrophage	unknown
Macrophages bone derived	unknown	unknown	unknown
MALME-3M	Skin	Melanocyte	Adherent and suspension
MAMMARY TUMOR MOUSE TRANSGENIC	Mammary glands	Mammary ductal epithelial cells	unknown
Mast Cells	Blood	Mast cell	unknown
MAT-Ly-Lu (CRL2376)	unknown	unknown	unknown
MBDK	Kidney	Kidney cells	Adherent
MBSa1	Connective tissue	unknown	unknown
MC/CAR	Peripheral blood	B-lymphocyte	Suspension
MC116	Lymphatic system	B-lymphocyte	Suspension
MC38	Colon	Colon epithelial cells	Adherent
MC38-CEA	Colon	Colon epithelial cells	Adherent
MC3T3E1	Bone	Osteoblast	Adherent
MCAS	Ovary	Ovarian cells	Adherent
MCC13	Skin	Merkel cell	Adherent
MCC26	Skin	Merkel cell	Adherent
MCEC	Heart	Cardiac endothelial cell	unknown
MCF 10A	Mammary glands	Mammary ductal epithelial cells	Adherent
MCF 10F	Mammary glands	Mammary ductal epithelial cells	Adherent
MCF-12A	Mammary glands	Mammary ductal epithelial cells	Adherent
MCF7	Mammary glands	Mammary ductal epithelial cells	Adherent
MCF-7 MX20	Breast	Mammary ductal epithelial cells	Adherent
MCF7LUC	Mammary glands	Mammary ductal epithelial cells	Adherent
MC-IXC	Nervous system	Neuroblast	Adherent
MCL-AANAT	Bile ducts	Murine cholangiocytes	unknown
MDA PCa 2b	Prostate	Prostate cells	Adherent
MDA-2A	unknown	unknown	unknown
MDA-H2774	Ovary	Ovarian endometrial cells	Adherent
MDA-MB-134	Breast	Mammary lobular cells	Adherent
MDA-MB-134-VI	Breast	Mammary lobular cells	Adherent
MDA-MB-157	Mammary glands	Mammary ductal epithelial cells	Adherent
MDA-MB-157	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-175	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-175-VII	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-231	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-330	Breast	Mammary lobular cells	Adherent
MDA-MB-361	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-415	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-435	Skin	Melanocyte	unknown
MDA-MB-435S	Skin	Melanocyte	Adherent
MDA-MB-436	Breast	Mammary ductal epithelial cells	Adherent
MDA-MB-453	Breast	Mammary ductal epithelial cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
MDA-MB-468	Breast	Mammary ductal epithelial cells	Adherent
MDA-MET 28	Breast	Mammary ductal epithelial cells	Adherent
MDA-PCa-2b	Prostate	Prostate cells	Adherent
MDBK	Kidney	Kidney epithelial cell	Adherent
MDCK	Kidney	Kidney epithelial cell	Adherent
MDDC	Blood	Monocyte-derived dendritic cells	unknown
MDSC myeloid-derived suppressor cells	Blood	myeloid-derived suppressor cells	unknown
MDST8	Colon	Colon epithelial cells	Adherent
ME-1	unknown	unknown	unknown
ME-180	Cervix	Cervical squamous cells	Adherent
MEC-1	Peripheral blood	B-lymphocyte	Suspension
MEC-2	Peripheral blood	B-lymphocyte	Suspension
Medullary thymic epithelial cells	Thymus	Medullary thymic epithelial cells	unknown
MEF	Connective tissue	Embryonic fibroblasts	unknown
MEG-01	Bone marrow	Megakaryoblast	Suspension
Melan-a	unknown	unknown	unknown
Melan-c	unknown	unknown	unknown
MEL-F-NEO	Skin	Neonatal melanocyte	unknown
MEL-HO	Skin	Melanocyte	Adherent
MEL-JUSO	Skin	Melanocyte	Adherent
Mesenchymal cells	Connective tissue	Mesenchymal cells	unknown
Mesenchymal Stem cells	Connective tissue	Mesenchymal stem cells	unknown
mesenchymal stem cells umbilical cord blood-derived	Umbilicall cord	Mesenchymal stem cells	unknown
Mesenteric adipose tissue adipocytes	Adipose tissue	Mesenteric adipose tissue adipocytes	unknown
MES-SA	Uterus	Uterine connective tissue cells	Adherent
MES-SA/Dx-5	Uterus	Uterine connective tissue cells	Adherent
Metschnikowia fructicola	unknown	Yeast	unknown
MeWo	Skin	Melanocyte	Adherent
MFE-280	Uterus	Uterine endometrial cells	Adherent
MFE-296	Uterus	Uterine endometrial cells	Adherent
MFE-319	Uterus	Uterine endometrial cells	Adherent
MFH-ino	Uterus	Uterine non-endometrial cells	Adherent
MFM-223	Breast	Mammary ductal epithelial cells	Adherent
MG-63	Bone	Osteoblast	Adherent
MG63.3	Bone	Osteoblast	Adherent
MGH-BA-1	Skin	Melanocyte	unknown
MGH-BO-1	Skin	Melanocyte	unknown
MGH-MC-1	Skin	Melanocyte	unknown
MGH-MCC-1	Skin	Melanocyte	unknown
MGH-PO-1	unknown	unknown	unknown
MGH-QU-1	Skin	Melanocyte	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
MGH-ST-1	Skin	Melanocyte	unknown
MGH-SW-1	Skin	Melanocyte	unknown
MGH-TH-1	unknown	unknown	unknown
MH-1	Lymphatic system	Lymphocytes	unknown
MH-2	Skin	Melanocyte	unknown
MHH-CALL-2	Blood	B-precursor cells	Suspension
MHH-ES-1	Muscle	Ewing sarcoma cells	Adherent
MHH-NB-11	Nervous system	Neuroblasts	Adherent
MHH-PREB-1	Lymphatic system	B-lymphocytes	Suspension
MIA	Pancreas	Pancreatic cells	Adherent
MIN6	Pancreas	Pancreatic cells	Adherent and suspension
MIN6-4N	Pancreas	Pancreatic Beta cells	unknown
MJ	Peripheral blood	T-lymphocyte	Suspension
MK2206	unknown	unknown	unknown
MKL-1	Skin	Merkel cell	Suspension
MKL-1 subclone 2	Skin	Merkel cell	Suspension
MKL-2	Skin	Merkel cell	Suspension
MKN1	Stomach	Gastric epithelial cells	Adherent
MKN28	Stomach	Gastric cells	Adherent
MKN45	Stomach	Gastric cells	Adherent and suspension
MKN7	Stomach	Gastric cells	Adherent
MKN74	Stomach	Gastric cells	Adherent
ML-1	unknown	unknown	unknown
ML-2	Peripheral blood	Myeloid cells	Suspension
MLLAF9	unknown	unknown	unknown
MLL-ENL	unknown	unknown	unknown
MLMA	Blood	Leukocyte	Suspension
MM455	Skin	Melanocyte	unknown
MM608	Skin	Melanocyte	unknown
MMAC-SF	Skin	Melanocyte	unknown
MMP12	unknown	unknown	unknown
MN-60	Peripheral blood	B-lymphocyte	Suspension
MNNG/HOS Cl #5	Bone	Osteoblast	Adherent
Mo	Blood	B-lymphoblast	Suspension
Mo-B	Blood	B-lymphoblast	Suspension
MOG-G-CCM	Brain	Astrocyte	Adherent
MOG-G-UVW	Brain	Astrocyte	Adherent
MOLM-13	Peripheral blood	Myeloid cells	Suspension
MOLM-16	Peripheral blood	Myeloid cells	Suspension
MOLM-20	Blood	Neutrophil	Suspension
MOLP-8	Peripheral blood	B-lymphocyte	Suspension
MOLT-13	Peripheral blood	T-lymphocyte	Suspension
MOLT-14	Peripheral blood	T-lymphocyte	Suspension

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
MOLT-16	Peripheral blood	T-lymphocyte	Suspension
MOLT-17	Peripheral blood	T-lymphocyte	Suspension
MOLT-3	Peripheral blood	Lymphocyte	Suspension
MOLT-4	Peripheral blood	Lymphocyte	Suspension
MOLT-4F	Peripheral blood	T-lymphocyte	Suspension
MONOCYTES BLOOD	Blood	Monocyte	unknown
MONO-MAC-1	Blood	Monocyte	Suspension
MONO-MAC-6	Blood	Monocyte	Suspension
Mononuclear cells	Peripheral blood	Mononuclear cells	unknown
Mo-T	Tongue	Squamous cells	unknown
MP2	unknown	unknown	unknown
MPM04/BAR	unknown	unknown	unknown
MPM10/CORO	unknown	unknown	unknown
MPM11/DEL	unknown	unknown	unknown
MPM17/ GAG	unknown	unknown	unknown
MPM24/MART	unknown	unknown	unknown
MPM25/MAY	unknown	unknown	unknown
MPM28/MLD	unknown	unknown	unknown
MPM59/LLA	unknown	unknown	unknown
MPM60/MASS	unknown	unknown	unknown
MPP-89	Pleura	Mesothelial cells	Adherent
MRC5	Lung	Lung cells	Adherent
MRC-9	Lung	Lung fibroblast	Adherent
MRK-nu-1	Breast	Mammary ductal epithelial cells	Adherent
MS-1-L	Lung	Lung cells	Adherent
MS751	Cervix	Cervical squamous cells	Adherent
MSC-1	unknown	unknown	unknown
MSTO-211H	Pleura	Mesothelial cells	Adherent
MT-3	unknown	unknown	unknown
mucor indicus	unknown	Fungal cell	unknown
multinucleated osteoclastic cells	Bone	Osteoclast	unknown
MUTZ-1	Lymphatic system	B-lymphocytes	Adherent and suspension
MUTZ-2	Peripheral blood	Myeloid cells	Suspension
MUTZ-3	Peripheral blood	Myeloid cells	Suspension
MV-4-11	Peripheral blood	Macrophage	Suspension
MVEC	Blood vessels	Microvascular endothelial cells	unknown
Mvt-1	Breast	Mammary ductal epithelial cells	Adherent
MY	Peripheral blood	Mononuclear cells	Suspension
MY-M12	Peripheral blood	Mononuclear cells	Suspension
MY-M13	Peripheral blood	Mononuclear cells	Suspension
MZ1-PC	Pancreas	Pancreas cell	unknown
MZ2-MEL	Skin	Melanocyte	unknown
MZ7-mel	Skin	Melanocyte	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
N. gaditana	unknown	Algae	unknown
N.crassa	unknown	Fungal cell	unknown
N488D	unknown	unknown	unknown
NA6	unknown	unknown	unknown
NAE	unknown	unknown	unknown
NALM-1	Peripheral blood	Lymphoblast	Suspension
NALM-19	Peripheral blood	B-lymphocyte	Suspension
NALM-6	Peripheral blood	B-lymphocyte	Suspension
NAMALWA	Lymphatic system	B-lymphoblast	Suspension
NAMALWA.CSN/70	Lymphatic system	B-lymphoblast	Suspension
NAMALWA.IPN/45	Lymphatic system	B-lymphoblast	Suspension
NAMALWA.KN2	Lymphatic system	B-lymphoblast	Suspension
NAMALWA.PNT	Lymphatic system	B-lymphoblast	Suspension
natural human dermal fibroblast (NHDF)	Skin	Fibroblast	unknown
NB(TU)1-10	Nervous system	Neuroblast	Adherent
NB-1	Mouth (inner cheek)	Squamous cells	unknown
NB14	Nervous system	Neuroblast	unknown
NB16	Nervous system	Neuroblast	unknown
NB17	Nervous system	Neuroblast	unknown
NB-4	Peripheral blood	Promelocyte	Suspension
NB5-B2	Nervous system	Neuroblast	unknown
NB69	Nervous system	Neuroblast	Adherent
NBL1	unknown	unknown	unknown
NBsusSR	Nervous system	Neuroblast	Adherent
NC-37	Lymphatic system	B-lymphoblast	Suspension
NCC16-P11	Cervix	Endo-cervical columner cells	Adherent
NCC-IT-A3	Testicle	Testicular germ cells	Semi-adherent
NCE 16IIA	Cervix	Endo-cervical columner cells	Adherent
NCI/ADR-RES	Ovary	Ovary cell	unknown
NCI/ADR-RES-EGFP	Ovary	Ovary cell	unknown
NCI-2347	Lung	Lung cells	Adherent
NCI-ADR-RES	Ovary	Ovary cell	unknown
NCI-BL1607	Peripheral blood	B-lymphoblast	Suspension
NCI-H1048	Lung	Lung small cells	Adherent
NCI-H1092	Lung	Lung small cells	Suspension
NCI-H1105	Lung	Lung small cells	Suspension
NCI-H1155	Lung	Lung non-small cells	Suspension
NCI-H1184	Lung	Lung small cells	Suspension
NCI-H1238	Lung	Lung small cells	Suspension
NCI-H128	Lung	Lung small cells	Suspension
NCI-H1299	Lung	Lung large cells	Adherent
NCI-H1304	Lung	Lung small cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
NCI-H1341	Lung	Lung small cells	Suspension
NCI-H1355	Lung	Lung cells	Suspension
NCI-H1385	Lung	Lung non-small cells	Suspension
NCI-H1395	Lung	Lung cells	Adherent
NCI-H1404	Lung	Lung cells	Suspension
NCI-H1417	Lung	Lung small cells	Suspension
NCI-H1435	Lung	Lung non-small cells	Adherent
NCI-H1436	Lung	Lung small cells	Suspension
NCI-H1437	Lung	Lung non-small cells	Adherent
NCI-H146	Lung	Lung non-small cells	Suspension
NCI-H1522	Lung	Lung small cells	Adherent and suspension
NCI-H1563	Lung	Lung non-small cells	Adherent
NCI-H1568	Lung	Lung non-small cells	unknown
NCI-H157	Lung	Lung squamous cells	unknown
NCI-H1581	Lung	Lung large cells	Adherent
NCI-H1618	Lung	Lung small cells	Suspension
NCI-H1623	Lung	Lung non-small cells	Adherent
NCI-H1648	Lung	Lung cells	Adherent
NCI-H1650	Lung	Lung cells	Adherent
NCI-H1651	Lung	Lung non-small cells	Adherent
NCI-H1666	Lung	Lung cells	Adherent
NCI-H1672	Lung	Lung small cells	Suspension
NCI-H1688	Lung	Lung small cells	Adherent
NCI-H1693	Lung	Lung cells	unknown
NCI-H1694	Lung	Lung small cells	Suspension
NCI-H1703	Lung	Lung non-small cells	Adherent
NCI-H1734	Lung	Lung non-small cells	Adherent
NCI-H1755	Lung	Lung non-small cells	Adherent
NCI-H1770	Lung	Lung non-small neuroendocrine cells	Adherent
NCI-H1781	Lung	Lung cells	Adherent
NCI-H1792	Lung	Lung cells	Adherent
NCI-H1793	Lung	Lung non-small cells	Adherent
NCI-H1819	Lung	Lung cells	Adherent and suspension
NCI-H1836	Lung	Lung small cells	Suspension
NCI-H1838	Lung	Lung non-small cells	Adherent
NCI-H1869	Lung	Lung squamous cells	Adherent
NCI-H187	Lung	Lung small cells	Suspension
NCI-H1876	Lung	Lung small cells	Adherent
NCI-H1882	Lung	Lung small cells	Adherent
NCI-H1915	Lung	Lung large cells	Adherent
NCI-H1926	Lung	Lung small cells	unknown
NCI-H1930	Lung	Lung small cells	Suspension
NCI-H1944	Lung	Lung non-small cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
NCI-H196	Lung	Lung small cells	Adherent
NCI-H1963	Lung	Lung small cells	Suspension
NCI-H1975	Lung	Lung cells	Adherent
NCI-H1993	Lung	Lung cells	Adherent
NCI-H1994	Lung	Lung cells	unknown
NCI-H2009	Lung	Lung cells	unknown
NCI-H2023	Lung	Lung non-small cells	unknown
NCI-H2029	Lung	Lung small cells	Adherent
NCI-H2030	Lung	Lung non-small cells	Adherent
NCI-H2059	Lung	Lung small cells	unknown
NCI-H2066	Lung	Lung squamous cells	Adherent
NCI-H2073	Lung	Lung non-small cells	Adherent
NCI-H2081	Lung	Lung small cells	Suspension
NCI-H2085	Lung	Lung non-small cells	Adherent
NCI-H2087	Lung	Lung non-small cells	Adherent
NCI-H209	Lung	Lung non-small cells	Suspension
NCI-H2106	Lung	Lung large cells	unknown
NCI-H2107	Lung	Lung small cells	Suspension
NCI-H2108	Lung	Lung small cells	Suspension
NCI-H211	Lung	Lung small cells	Suspension
NCI-H2110	Lung	Lung small cells	unknown
NCI-H2122	Lung	Lung cells	Adherent
NCI-H2126	Lung	Lung large cells	Adherent
NCI-H2135	Lung	Lung large cells	unknown
NCI-H2141	Lung	Lung small cells	Suspension
NCI-H2170	Lung	Lung squamous cells	Adherent
NCI-H2171	Lung	Lung small cells	Suspension
NCI-H2172	Lung	Lung large cells	Adherent
NCI-H2195	Lung	Lung small cells	Adherent
NCI-H2196	Lung	Lung small cells	unknown
NCI-H2198	Lung	Lung small cells	unknown
NCI-H220	Lung	Lung small cells	Suspension
NCI-H2227	Lung	Lung small cells	Adherent and suspension
NCI-H2228	Lung	Lung large cells	Adherent
NCI-H226	Pleura	Pleural squamous cells	Adherent
NCI-H2286	Lung	Lung small cells	Adherent
NCI-H2291	Lung	Lung large cells	Adherent
NCI-H23	Lung	Lung large cells	Adherent
NCI-H2330	Lung	Lung small cells	Suspension
NCI-H2342	Lung	Lung large cells	Adherent
NCI-H2347	Lung	Lung large cells	Adherent
NCI-H2405	Lung	Lung large cells	Adherent and suspension
NCI-H2444	Lung	Lung large cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
NCI-H2452	Pleura	Mesothelial cells	Adherent
NCI-H250	Lung	Lung small cells	Suspension
NCI-H292	Lung	Lung epidermal cells	Adherent
NCI-H3122	Lung	Lung cells	unknown
NCI-H322	Lung	Lung cells	unknown
NCI-H322M	Lung	Lung cells	unknown
NCI-H345	Lung	Lung small cells	Adherent and suspension
NCI-H358	Lung	Lung cells	Adherent
NCI-H378	Lung	Lung small cells	Suspension
NCI-H441	Lung	Papillary lung cells	Adherent
NCI-H446	Lung	Lung small cells	Adherent and suspension
NCI-H460	Lung	Lung large cells	Adherent
NCI-H510A	Lung	Lung small cells	Adherent and suspension
NCI-H520	Lung	Lung squamous cells	Adherent
NCI-H522	Lung	Lung large cells	Adherent
NCI-H524	Lung	Lung small cells	Suspension
NCI-H526	Lung	Lung small cells	Suspension
NCI-H596	Lung	Lung adenosquamous cells	Adherent
NCI-H630	Undefined (contaminated)	Undefined (contaminated)	unknown
NCI-H64	Lung	Lung small cells	Suspension
NCI-H647	Lung	Lung adenosquamous cells	Adherent
NCI-H650	Lung	Lung large cells	Suspension
NCI-H660	Prostate	Prostate small cell	Adherent and suspension
NCI-H661	Lung	Lung large cells	Adherent
NCI-H676B	Lung	Lung cells	Suspension
NCI-H69	Lung	Lung small cells	Suspension
NCI-H711	Lung	Lung small cells	Suspension
NCI-H716	Cecum	Cecum epithelial cells	Suspension
NCI-H719	Lung	Lung small cells	Suspension
NCI-H720	Lung	Lung cells	Suspension
NCI-H727	Lung	Lung cells	Adherent
NCI-H735	Lung	Lung small cells	Suspension
NCI-H740	Lung	Lung small cells	Suspension
NCI-H747	Cecum	Cecum epithelial cells	Adherent and suspension
NCI-H748	Lung	Lung small cells	Suspension
NCI-H774	Lung	Lung small cells	Suspension
NCI-H810	Lung	Lung large cells	Adherent
NCI-H82	Lung	Lung small cells	Suspension
NCI-H820	Lung	Papillary lung cells	Adherent
NCI-H835	Lung	Lung cells	Suspension
NCI-H838	Lung	Lung large cells	Adherent
NCI-H841	Lung	Lung small cells	Adherent and suspension
NCI-H847	Lung	Lung small cells	Suspension

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
NCI-H865	Lung	Lung small cells	Suspension
NCI-H889	Lung	Lung small cells	Suspension
NCI-H920	Lung	Lung large cells	Suspension
NCI-H929	Peripheral blood	B lymphocyte	Suspension
NCI-H969	Lung	Lung large cells	unknown
NCI-N87	Stomach	Gastric epithelial cells	Adherent
NEC14	Testicle	Testicular germ cell	Adherent
NEC8	Testicle	Testicular germ cell	Adherent
Neonatal ventricular cardiomyocytes Primary	Heart (ventricle)	Cardiomyocytes	unknown
Neural progenitor cells (NPCs)	Brain	Neural progenitor cells	unknown
Neuro-2a	Nervous system	Neuroblast	Adherent
Neuro-2a	Nervous system	Neuroblast	Adherent
Neuroblastoma	Nervous system	Neuroblast	unknown
Neuroscreen-1	Nervous system	Chromaffin cells	unknown
Neutrophil	Blood	Neutrophil	unknown
Neutrophil extracellular traps NETs	unknown	unknown	unknown
Newborn Keratinocytes	Skin	Keratinocytes	unknown
NH-12	Nervous system	Neuroblast	Adherent
NH-6	Nervous system	Neuroblast	Adherent
NHC	Bile ducts	Bile duct epithelial cells	unknown
NHDF	Skin	Dermal fibroblasts	Adherent
NIH	unknown	unknown	unknown
NK CELLS	Blood	Natural killer cells	unknown
NK-92	Peripheral blood	Natural killer cells	Suspension
NK-92MI	Peripheral blood	Natural killer cells	Suspension
NKM-1	Peripheral blood	Myeloid cells	Suspension
NL20	Lung	Bronchial epithelial cell	Adherent
NL20T-A	Lung	Bronchial epithelial cell	Adherent
NMC-G1	Brain	Astrocyte	Adherent
NOMO-1	Peripheral blood	Monocyte	Suspension
NOMO-1/ADM	Peripheral blood	Myeloid cells	Suspension
NOS-1	Bone	Osteoblast	Adherent
NrasG12D	unknown	unknown	unknown
NsAtWR11 cells	unknown	unknown	unknown
NSCs	Central nervous sysem	Neural stem cell	Adherent
NSPCs	unknown	unknown	unknown
nsTCs	unknown	unknown	unknown
NTERA-S-cl-D1	Testicle	Germ cells	Adherent
Nthy-ori 3-1	Thyroid	Thyroid follicular epithelial cells	Adherent
NU-DUL-1	Lymphatic system	B-lymphocytes	Suspension
NUGC-2	Stomach	Gastric epithelial cells	Adherent
NUGC-3	Stomach	Gastric epithelial cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
NUGC-4	Stomach	Gastric epithelial cells	Suspension
NY	Bone	Osteoblast	Adherent
O157:H7	unknown	unknown	unknown
OAW28	Ovary	Ovary cell	Adherent
OAW42	Ovary	Ovary cell	Adherent
OAW42	Ovary	Ovarian epithelial cells	Adherent
OC-314	Ovary	Ovary cell	Adherent
OC-314	Ovary	Ovarian cells	Adherent
OCI-AML2	Peripheral blood	Myeloid cells	Suspension
OCI-AML2	Peripheral blood	Myeloid cells	Suspension
OCI-AML3	Peripheral blood	Myeloid cells	Suspension
OCI-AML3	Peripheral blood	Myeloid cells	Suspension
OCI-AML5	Peripheral blood	Myeloid cells	Suspension
OCI-AML5	Peripheral blood	Myeloid cells	Suspension
OCI-Ly1	Lymphatic system	B-lymphocytes	Suspension
OCI-Ly1	Leukemia	B lymphocyte	Suspension
OCI-LY10	Lymphatic system	B-lymphocytes	unknown
OCI-Ly10	Leukemia	B lymphocyte	Suspension
OCI-LY19	Lymphatic system	B-lymphocytes	Suspension
OCI-Ly-19	Leukemia	B lymphocyte	Suspension
OCI-LY2	Lymphatic system	B-lymphocytes	unknown
OCI-Ly-2	Leukemia	B lymphocyte	unknown
OCI-LY3	Lymphatic system	B-lymphocytes	Suspension
OCI-Ly-3	Leukemia	B lymphocyte	Suspension
OCI-M1	Peripheral blood	Myeloid cells	Suspension
OCI-M1	Leukemia	B lymphocyte	Suspension
OCUB-M	Breast	Mammary epithelial cells	Adherent
OCUB-M	Breast	Mammary epithelial cells	Adherent
OCUG-1	Gallbladder	Gallbladder cells	Adherent
OCUG-1	Gallbladder	Gall bladder epithelial cells	Adherent
OE19	Esophagus	Esophageal epithelial cells	Adherent
OE19	Esophagus	Esophageal epithelial cells	Adherent
OE21	Esophagus	Esophageal squamous cells	Adherent
OE21	Esophagus	Esophageal squamous epithelial cell	Adherent
OE33	Esophagus	Esophageal epithelial cells	Adherent
OE33	Esophagus	Esophageal epithelial cells	Adherent
OMC-1	Cervix	Cervical squamous cells	Adherent
OMC-1	Cervix	Squamous epithelial cells	Adherent
ONCO-DG-1	Ovary	Ovary cell	Adherent
ONCO-DG-1	Ovary	Ovary epithelial cells	Adherent
Onda 10	Brain	Glial cell	Adherent
Onda 11	Brain	Glial cell	Adherent
ONS-76	Neural/Brain	Neuroepithelial cell	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
OPM-2	Bone marrow	Plasma cells	Suspension
OS187	Brain	Colon	unknown
OSA 4463997	unknown	unknown	unknown
OSA 4512351	Leukemia	unknown	unknown
OSA 7764	unknown	unknown	unknown
OSA 80	Bone	unknown	unknown
OSC-19	Tongue	Tongue squamous epithelial cells	unknown
OSC-20	Mouth	Squamous epithelial cells	Adherent
OS-RC-2	Kidneys	Renal epithelial cells	Adherent
OUMS-23	Colon/rectum	Colorectal epithelial cells	Adherent
OUS-11	Lungs	Fibroblast-like cells	Adherent
OV-1063	unknown	unknown	unknown
OV-90	Ovary	Ovarian epithelial cells	Adherent
OVCAR-3	Ovary	Ovarian epithelial cells	Adherent
OVCAR-4	Ovary	Ovarian epithelial cells	Adherent
OVCAR-5	Ovary	Ovarian epithelial cells	Adherent
OVCAR-8	Ovary	Ovarian epithelial cells	Adherent
OVISE	Ovary	Ovarian epithelial cells	Adherent
OVKATE	Ovary	Ovarian epithelial cells	Adherent
OVMANA	Ovary	Ovarian epithelial cells	Adherent
OVMIU	Ovary	Ovarian epithelial cells	Adherent
OVSAYO	Ovary	Ovarian epithelial cells	Adherent
OVTOKO	Ovary	Ovarian epithelial cells	Adherent
P116	Blood	T lymphocyte	Suspension
P116.cl39	Blood	T lymphocyte	Suspension
P12-ICHIKAWA	Blood	T lymphocyte	Suspension
P30/OHK	Blood	B lymphocyte	Suspension
P31/FUJ	Blood	Myeloid cells	Suspension
P32/ISH	Blood	B lymphocyte	Suspension
P3HR-1	Ascites	B lymphocyte	Suspension
PA-1	Ovary	Ovarian epithelial cells	Adherent
PaCa-2	unknown	unknown	unknown
Panc 02.03	Pancreas	Pancreatic epithelial cells	Adherent
Panc 02.13	Pancreas	Pancreatic epithelial cells	Adherent
Panc 03.27	Pancreas	Pancreatic epithelial cells	Adherent
Panc 04.03	Pancreas	Pancreatic epithelial cells	Adherent
Panc 08.13	Pancreas	Pancreatic epithelial cells	Adherent
Panc 10.05	Pancreas	Pancreatic epithelial cells	Adherent
Panc06.05	Pancreas	unknown	Adherent
Panc08.13	Pancreas	Pancreatic epithelial cells	Adherent
PANC-1	Pancreas	Pancreatic epithelial cells	Adherent
Panc10.05	Pancreas	Pancreatic epithelial cells	Adherent
PARG null	unknown	unknown	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Patellar Tendon Stem Cells (PTSC)	unknown	unknown	unknown
PA-TU-8902	Pancreas	Pancreatic epithelial cells	Adherent
PA-TU-8988S	Pancreas	Pancreatic epithelial cells	Adherent
PA-TU-8988T	Pancreas	Pancreatic epithelial cells	unknown
PBMC peripheral blood mononuclear cells	unknown	unknown	unknown
PBMCs human	unknown	unknown	unknown
PC12	Adrenal medulla	Adrenal cells	Adherent
PC-14	Lung	Lung epithelial cells	Adherent
PC-3	Prostate	Prostate epithelial cells	Adherent
PC3-EMT	Prostate	Prostate epithelial cells	Adherent
PC-9	Lung	Lung adenocarcinoma	Mixed
PC9(IR)1	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)10	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)11	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)2	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)3	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)5	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)6	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)7	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(IR)9	Lung	Lung adenocarcinoma	unknown
PC9(TAR)1	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)10	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)11	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)2	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)3	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)4	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)5	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)6	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)7	non-small cell carcinoma	Lung:NSCLC	unknown
PC9(TAR)8	non-small cell carcinoma	Lung:NSCLC	unknown
PC9/GR6	Lung	Lung cells	Adherent and suspension
PCI 13	Upper aerodigestive tract or mouth	Squamous epithelial cells	Adherent
PCI-15A	Upper aerodigestive tract, head, or neck	Squamous epithelial cells	Adherent
PCI-15B	Upper aerodigestive tract, head, or neck	Squamous epithelial cells	Adherent
PCI-30	Upper aerodigestive tract, head, or neck	Squamous epithelial cells	Adherent
PCI-37B	Larynx	Laryngeal epithelial cells	Adherent
PCI-38	Upper aerodigestive tract, head, or neck	Squamous epithelial cells	Adherent
PCI-4A	Larynx	Laryngeal epithelial cells	unknown
PCI-4B	Tongue	Squamous epithelial cells	Adherent
PCI-6A	Tonsils	Tonsillar squamous epithelial cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
PCI-6B	Neck lymph node	Squamous epithelial cells	Adherent
PCRP-IRF3-6C2	unknown	unknown	Adherent
PCS-200-013	Skin	Melanocytes	Adherent
PDX	Head & Neck	unknown	unknown
PE/CA-PJ15	Tongue	Tongue epithelial cells	Adherent
PEER	Blood	T lymphocytes	Suspension
Penicillium digitatum	unknown	unknown	Adherent
Penicillium expansum	unknown	unknown	unknown
Penicillium expansum	unknown	unknown	Adherent
Peripheral blood leukocyte (PBLs)	Blood	Leukocyte	Suspension
peripheral blood mononuclear cells	Blood	PBMCs	Suspension
Peripheral double negative (DNT) cells	Blood	T lymphocyte	Suspension
Peritoneal polymorphonuclear leukocytes	Blood	Leukocytes	Suspension
peritoneal cavity exudate cells (PEC)	unknown	unknown	unknown
PF-382	Blood	Leukocytes	Suspension
Pfeiffer	Blood	Lymphoblast B lymphocyte	Suspension
PFSK-1	Blood	Fibroblast cells	Adherent
PGVSMCs rat	unknown	unknown	unknown
Pichia fermentans	unknown	unknown	unknown
PL18	Lymphoma	unknown	unknown
PL-21	Blood	Myeloid cells	Suspension
PL4	Pancreas	Pancreatic epithelial cells	Adherent
PL45	Pancreas	Pancreatic epithelial cells	Adherent
Placental Tissue Cells	Placenta	unknown	Adherent
Plasma cells	Blood	B lymphocyte	Suspension
Platelet	Blood	Myeloid cells	unknown
PLC/PRF/5	Liver	Liver epithelial cells	Adherent
PNM	unknown	unknown	unknown
PNT1A	Prostate	Prostate epithelial cells	Adherent
PNT2	Prostate	Prostate epithelial cells	Adherent
primitive erythroid progenitor cells	unknown	Erythrocytes	unknown
progenitor cells (HSC)	Blood	Hematopoietic stem cell	Suspension
PSN1	Pancreas	Pancreatic epithelial cells	Adherent
PT67 retrovirus packaging	unknown	unknown	Adherent
Pulmonary epithelial lining fluid PELF	Lungs	Lung epithelial cells	unknown
Pure Cholangiocytes	Biliary tree	Epithelial cells	unknown
PWR-1E	Prostate	Prostate epithelial cells	adherent
QGP-1	Pancreas	Pancreatic epithelial cells	adherent
QIMR-WIL	Blood	Myeloid cells	Suspension
R1 ESC	unknown	unknown	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Raji	Maxilla	B Lymphocytes	Suspension
Ramos	Ascitic fluid	B Lymphocytes	Suspension
Ramos-2G6-4C10	Ascitic fluid	B Lymphocytes	Suspension
RAW264.7	Blood	Macrophage/Monocyte	Partial Adherent
RBC	unknown	unknown	unknown
RBE	unknown	unknown	unknown
rBMSC	Bone marrow	Mesenchymal stem cells	Adherent
RCC10RGB	Kidney	Renal epithelial cells	Adherent
RCH-ACV	Bone marrow	B cells	Suspension
RC-K8	Peripheral blood	B lymphocyte	Suspension
RCM-1	Rectum	Rectal epithelial cells	Adherent
RD	Muscle	Striated muscle cell	Adherent
RD-ES	Muscle	Ewing sarcoma cells	Adherent
REC-1	Lymphatic system	B-lymphocyte	Suspension
Reh	Blood	B-lymphocyte	Suspension
RERF-GC-1B	Stomach	Stomach cells	Adherent
RERF-LC-Ad1	Lung	Lung cells	Adherent
RERF-LC-Ad2	Lung	Lung cells	Adherent
RERF-LC-FM	Lung	Lung small cells	Suspension
RERF-LC-KJ	Lung	Lung cells	Adherent
RERF-LC-MA	Lung	Lung small cells	Adherent
RERF-LC-MS	Lung	Lung cells	Adherent
RERF-LC-Sq1	Lung	Lung squamous cells	Adherent
RF-48	Peripheral blood	Lymphocyte	unknown
RH-1	Muscle	Ewing sarcoma cells	Adherent
RH-18	Muscle	Striated muscle cell	Adherent
RH-30	Muscle	Striated muscle cell	Adherent
RH-41	Muscle	Striated muscle cell	Adherent
Rhabdomyosarcoma cell line (RMS)	Muscle	Striated muscle cell	unknown
RK-13	Kidney	Kidney epithelial cell	Adherent
RKN	Ovary	Ovarian cells	Adherent
RKO	Colon	Colorectal epithelial cells	Adherent
RL	Lymphatic system	B-lymphocytes	Suspension
RL95-2	Uterus	Uterine endometrial cells	Adherent
RLM	unknown	unknown	unknown
RMG-I	Ovary	Ovarian epithelial cell	Adherent
RMG-II	Ovary	Ovarian epithelial cell	Adherent
RO82-W-1	Thyroid	Thyroid follicular epithelial cells	Adherent
ROS-50	Peripheral blood	B-lymphocytes	Suspension
RPMI 1788	Peripheral blood	B-lymphocytes	Suspension
RPMI 2650	Head & Neck	Squamous cells	Adherent
RPMI 6666	Peripheral blood	B-lymphoblast	Suspension

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
RPMI 8226	Peripheral blood	B-lymphoblast	Suspension
RPMI-7951	Skin	Melanocytes	Adherent
RPMI 8226	Peripheral blood	B-lymphoblast	Suspension
RPMI-8402	Blood	T-lymphocytes	Suspension
RPMI-8866	Peripheral blood	B-lymphoblast	Suspension
RS4;11	Blood	B-lymphocyte	Suspension
RS453	unknown	unknown	unknown
RT-112	Urinary tract	Urinary transitional cell	Adherent
RT112/84	Bladder	Bladder epithelial cells	Adherent
RT4	Urinary tract	Urinary transitional cell	Adherent
RTC renal tubular cells	Kidney	Renal tubular cells	unknown
RVH-421	Skin	Melanocytes	Adherent
RWPE-1	Prostate	Prostate epithelial cells	Adherent
RWPE-2	Prostate	Prostate epithelial cells	Adherent
RWPE2-W99	Prostate	Prostate epithelial cells	Adherent
RXF 393	Kidney	Renal cells	unknown
S. bayanus	unknown	Yeast	unknown
S. bayanusuvarum	unknown	Yeast	unknown
S. castellii	unknown	Yeast	unknown
S. cerevisiae	unknown	Yeast	unknown
S. japonicus	unknown	Yeast	unknown
S. kluyveri	unknown	Yeast	unknown
S. mikatae	unknown	Yeast	unknown
S. paradoxus	unknown	Yeast	unknown
S. pombe	unknown	Yeast	unknown
S-117	Thyroid	Thyroid cells	Adherent
SACC-83	Salivary gland	Salivary gland cells	unknown
SaOS-2	Bone	Osteocyte	Adherent
Sarc9371	Bone	Osteocyte	Adherent
SAS	Tongue	Squamous cell tongue	Adherent
SAT	Mouth	Head and neck squamous cells	Adherent
Satellite cells	Muscle	Satellite cells	unknown
SBC-1	Lung	Lung small cells	Suspension
SBC-2	Lung	Lung small cells	Suspension
SBC-3	Lung	Lung small cells	Adherent
SBC-5	Lung	Lung small cells	Adherent
SC-1	Lymphatic system	B-lymphocyte	Suspension
SCaBER	Bladder	Bladder squamous cells	Adherent
SCC-15	Tongue	Squamous cell tongue	Adherent
SCC-25	Tongue	Squamous cell tongue	Adherent
SCC-3	Peripheral blood	B-lymphocyte	Suspension
SCC-4	Tongue	Squamous cell tongue	Adherent
SCC-9	Tongue	Squamous cell tongue	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
SCC90	Tongue	Squamous cell tongue	Adherent
SCCH-196	Muscle	Ewing sarcoma cells	Adherent
SCCH-26	Nervous system	Neuroblast	Adherent
SCH	Stomach	Placental cells	Adherent
Schizochytrium	unknown	Protist cell	unknown
Schneider 2	Embryo	Embryonic stem cell	unknown
Sc-1	Lymphatic system	B-lymphocyte	Suspension
SCLC-21H	Lung	Lung small cells	Suspension
SEG-1	Lung	Lung large cells	Adherent
Serratia marcescens	unknown	Bacterial cell	unknown
SET-2	Peripheral blood	unknown	Suspension
SF126	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF126-BM	Brain	Astrocyte	Adherent
SF268	Brain	Astrocyte	unknown
SF268-BM	Brain	Astrocyte	unknown
SF268-P	Brain	Astrocyte	unknown
SF268-TR10-BM	Brain	Astrocyte	unknown
SF268-TR1-BM	Brain	Astrocyte	unknown
SF268-TR2-BM	Brain	Astrocyte	unknown
SF268-TR4-BM	Brain	Astrocyte	unknown
SF268-TR6-BM	Brain	Astrocyte	unknown
SF268-TR7-BM	Brain	Astrocyte	unknown
SF295	Brain	Astrocyte	Adherent
SF295-BM	Brain	Astrocyte	Adherent
SF295-TR2-BM	Brain	Astrocyte	Adherent
SF539	Brain	Astrocyte	Adherent
SF767	Cervix	Cervical squamous cells	unknown
Sf9	Ovary	Ovary cells	Suspension
SFCC	Cervix	Cervical cells	unknown
SG231	Liver	Cholangiocyte	Adherent
SGBS	Adipose tissues	Adipocytes	Adherent
SGBS\y	unknown	unknown	unknown
SGBS\vy	unknown	unknown	unknown
SGBS\987	unknown	unknown	unknown
SGBS\uy	unknown	unknown	unknown
SH-4	Skin	Melanocytes	Adherent
SH-EP	Nervous system	Neuroblast	Adherent
SHP-77	Lung	Lung small cells	Suspension

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
SH-SY5Y	Nervous system	Neuroblast	Adherent and suspension
SIG-M5	Peripheral blood	Myeloid cells	Suspension
SiHa	Cervix	Cervical squamous cells	Adherent
SIMA	Nervous system	Neuroblast	Adherent
SISO	Cervix	Cervical squamous cells	Adherent
RH30	Muscle	Striated muscle cell	Adherent
SJNB-10	Nervous system	Neuroblast	unknown
SJNB-12	Nervous system	Neuroblast	unknown
SJNB-7	Nervous system	Neuroblast	unknown
SJSA-1	Bone	Osteocyte	Adherent
SK-BR-3	Breast	Mammary ductal epithelial cells	Adherent
SK-CO-1	Colon	Colorectal epithelial cells	Adherent
Skeletal muscle cells	Muscle	Skeletal muscle	unknown
SK-ES-1	Muscle	Ewing sarcoma cells	Adherent
SKG-IIIa	Cervix	Cervical squamous cells	Adherent
SKG-IIIb	Cervix	Cervical squamous cells	Adherent
SK-HEP-1	Liver	Hepatocyte	Adherent
SK-LMS-1	Vulva	Smooth muscle	Adherent
SK-LU-1	Lung	Lung cells	Adherent
SKM-1	Peripheral blood	Myeloid cells	Suspension
SK-MEL-1	Skin	Melanocytes	Adherent
SK-MEL-119	Skin	Melanocytes	Adherent
SK-MEL-131	Skin	Melanocytes	Adherent
SK-MEL-2	Skin	Melanocytes	Adherent
SK-MEL-24	Skin	Melanocytes	Adherent
SK-MEL-28	Skin	Melanocytes	Adherent
SK-MEL-3	Skin	Melanocytes	Adherent
SK-MEL-30	Skin	Melanocytes	Adherent
SK-MEL-37	Skin	Melanocytes	unknown
SK-MEL-39	Skin	Melanocytes	unknown
SK-MEL-5	Skin	Melanocytes	Adherent
SK-MEL-63	Skin	Melanocytes	Adherent
SK-MES-1	Lung	Squamous cell (lung)	Adherent
SK-MG-1	Brain	Astrocyte	unknown
SK-MM-2	Peripheral blood	Plasma cell	Suspension
SKN	Uterus	Uterine endometrial cells	Adherent
SKN-3	Head and neck	Head and neck squamous cells	Adherent
SK-N-AS	Nervous system	Neuroblast	Adherent
SK-N-BE	Nervous system	Neuroblast	Adherent and suspension
SK-N-BE(2)	Nervous system	Neuroblast	Adherent and suspension
SK-N-DZ	Nervous system	Neuroblast	Adherent
SK-NEP-1	Muscle (previously thought to be kidney)	Ewing sarcoma cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
SK-N-FI	Nervous system	Neuroblast	Adherent
SK-N-MC	Chest wall	Neuroepithelial cells	Adherent
SK-N-SH	Nervous system	Neuroblast	Adherent
SK-OV-3	Ovary	Ovarian cells	Adherent
SK-PN-DW	Retroperitoneal embryonal tissue	Primitive neuroectodermal cell	Adherent
SK-UT-1	Uterus	Uterine endometrial cells	Adherent
VL51	Spleen	Lymphocyte	Suspension
SMCs	Smooth muscle	Smooth muscle cells	unknown
SN12C	Kidney	Renal epithelial cells	unknown
SNB13	Nervous system	Neuroblast	unknown
SNB-19	Nervous system	Neuroblast	Adherent
SNB-75	Nervous system	Astrocyte	unknown
SNG-M	Uterus	Uterine endometrial cancer cells	Adherent
SNU-1	Stomach	Stomach epithelial cells	Suspension
SNU-16	Stomach	Stomach epithelial cells	Suspension
SNU-182	Liver	Hepatocyte	Adherent
SNU-387	Liver	Hepatocyte	Adherent
SNU-398	Liver	Hepatocyte	Adherent and suspension
SNU-423	Liver	Hepatocyte	Adherent
SNU-449	Liver	Hepatocyte	Adherent and suspension
SNU-475	Liver	Hepatocyte	Adherent
SNU-5	Stomach	Stomach epithelial cells	Suspension
SNU-C1	Colon	Colorectal epithelial cells	Adherent
SNU-C2B	Cecum	Cecum epithelial cells	Suspension
SOB-15	Liver	Liver epithelial cells	Adherent
Sperm	Testes	Sperm	Suspension
Sperm cells from honey bees (<i>Apis mellifera</i>)	Testes	Sperm	Suspension
Spinous cells	Skin	Spinous cells	unknown
Spleen cells mouse	Spleen	Splenocytes	unknown
Splenocytes	Spleen	Splenocytes	unknown
SR	Lymphatic system	T-cells	Suspension
H. Pylori SS1	unknown	Bacterial cell	unknown
SSP-25	Liver	Hepatocytes	unknown
ST486	Lymphatic system	B-lymphocyte	Suspension
STA-NB-6	Nervous system	Neuroblast	unknown
STC 1	Lung	Lung cells	Adherent
Streptococcus gallolyticus	unknown	Bacterial cell	unknown
Streptococcus mutans	unknown	Bacterial cell	unknown
STS 0421	Stomach	Stomach epithelial cells	unknown
SU.86.86	Pancreas	Pancreas cell	Adherent
SUBMUCOSAL GLANDULAR EPITHELIAL CELLS	Small intestine	Submucosal glandular epithelial cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
SU-DHL-1	Peripheral blood	Histiocytes	Suspension
SU-DHL-10	Peripheral blood	B-lymphocyte	Suspension
SU-DHL-16	Peripheral blood	B-lymphocyte	Suspension
SU-DHL-4	Peripheral blood	B-lymphocyte	Suspension
SU-DHL-5	Peripheral blood	B-lymphocyte	Suspension
SUDHL-6	Peripheral blood	B-lymphocyte	Suspension
SU-DHL-8	Peripheral blood	B-lymphocyte	Suspension
SUIT-2	Pancreas	Pancreas cell	Adherent
SUM102PT	Breast	Breast epithelial	unknown
SUM149PT	Breast	Breast epithelial	unknown
SUM159	Breast	Breast epithelial	unknown
SUP-B15	Peripheral blood	B-lymphocytes	Suspension
SUP-B8	Peripheral blood	B-lymphocytes	Suspension
SUP-HD1	Lymphatic system	Reed-sternberg cells	Suspension
SUP-M2	Peripheral blood	Hystiocytes	Suspension
SupT1	Peripheral blood	T-lymphocytes	Suspension
SVC Stromal vascular cells	Stromal blood vessels	Stromal vascular cells	unknown
SVF cells	Stromal vascular fraction	Stromal vascular fraction cells	unknown
SVG p12	Brain (fetus)	Glial cells (Fetus)	Adherent
SVts-8	Connective tissue	Fibroblast	Adherent
SW 1088	Brain	Astrocyte	Adherent
SW 1116	Colon	Colorectal epithelial cells	Adherent
SW 1271	Lung	Lung cells	Adherent
SW 13	Adrenal gland	Adrenal gland cortex cells	Adherent
SW 1417	Colon	Colorectal epithelial cells	Adherent
SW 1463	Rectum	Rectal epithelial cells	Adherent
SW 156	Kidney	Renal epithelial cells	Adherent
SW 1573	Lung	Alveolar epithelial cells	Adherent
SW 1783	Brain	Astrocyte	Adherent
SW 1990	Pancreas	Pancreas cell	Adherent
SW 48	Colon	Colorectal epithelial cells	Adherent
SW 626	Colon	Colorectal epithelial cells	Adherent
SW 780	Bladder	Bladder epithelial cells	Adherent
SW 900	Lung	Squamous cell (lung)	Adherent
SW 1116	Colon	Colorectal epithelial cells	Adherent
SW-1710	Bladder	Bladder epithelial cells	Adherent
SW403	Colon	Colorectal epithelial cells	Adherent
SW872	Connective tissue	Adipocytes	Adherent
SW900	Lung	Lung squamous cells	Adherent
SW-948	Colon	Colorectal epithelial cells	Adherent
SW954	Vulva	Squamous cell	Adherent
SW962	Vulva	Squamous cell	Adherent
SW982	Synovium	Synovial cell	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
Swiss-3T3	Embryo	Embryonal fibroblast	Adherent
SWRC-GRO	unknown	unknown	unknown
Symbiodinium cells	unknown	Algae	unknown
Synechococcus elongatus PCC 7942	unknown	unknown	unknown
Synechocystis sp.	unknown	Water-dwelling bacteria	unknown
Synechocystis sp. PCC 6803	unknown	Water-dwelling bacteria	unknown
Synechocystis sp. PCC 6803	unknown	unknown	unknown
Synechocystis sp. ΔglgC	unknown	Water-dwelling bacteria	unknown
T castaneum	unknown	unknown	unknown
T.T	Esophogus	Esophageal squamous cells	Adherent
T.Tn	Esophogus	Esophageal squamous cells	Adherent
T1-73	Bone	Osteocyte	Adherent
T24	Urinary tract	Urinary transitional cell	Adherent
T-47D	Breast	Mammary ductal epithelial cells	Adherent
T47D-KBLuc	Breast	Mammary ductal epithelial cells	Adherent
T47DPAC	Breast	Mammary ductal epithelial cells	Adherent
T47DPAR	Breast	Mammary ductal epithelial cells	Adherent
T84	Colon	Colorectal epithelial cells	Adherent
T95-45D	Skin	Merkel cell	Adherent
T98G	Brain	Astrocyte	Adherent
Takigawa	Stomach	Stomach cells	unknown
TALL-1	Blood	T-lymphocyte	Suspension
TALL-104	Blood	T-lymphocyte	Suspension
TANOUE	Blood	B-lymphocyte	Suspension
TASK1	Chest wall	Neuroepithelial cells	Adherent
TBV2	Embryo	Embryonic stem cells	unknown
TC-71	Muscle	Ewing sarcoma cells	Adherent
TcA-CLG1	unknown	unknown	unknown
TCCSUP	Urinary tract	Urinary transitional cell	Adherent
TCO-1	Cervix	Cervical squamous cells	unknown
TC-YIK	Cervix	Cervical cells	Suspension
TE 161.T	Lymphatic system	Lymphocyte	unknown
TE 417.T	Bone	Osteocyte	unknown
TE 418.T	Bone	Osteocyte	unknown
TE-1	Esophogus	Esophageal squamous cells	Adherent
TE-10	Esophogus	Esophageal squamous cells	Adherent
TE-11	Esophogus	Esophageal squamous cells	Adherent
TE-12	Esophogus	Esophageal squamous cells	Adherent
TE-15	Esophogus	Esophageal squamous cells	Adherent
TE-441-T	Muscle	Striated muscle cell	Adherent
TE-5	Esophogus	Esophageal squamous cells	Adherent
TE-6	Esophogus	Esophageal squamous cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
TE7	Esophogus	Esophageal squamous cells	Adherent
TE-8	Esophogus	Esophageal squamous cells	Adherent
TE-9	Esophogus	Esophageal squamous cells	Adherent
TECs	Thymus	Thymic epithelial cells	unknown
Tera-1	Testicle	Testicular germ cells	Adherent
Testis cells equine	Testicle	Testicular cells	unknown
TF-1.CN5a.1	Peripheral blood	Erythrocyte	Suspension
TF-1a	Peripheral blood	Erythrocyte	Suspension
TFK-1	Bile ducts	Bile duct epithelial cells	Adherent
TGBC11TKB	Stomach	Stomach cells	unknown
TGBC14TKB	Gallbladder	Gallbladder cells	Adherent
TGBC18TKB	Bile ducts (Ampulla of Vater)	Bile duct epithelial cells	Adherent
TGBC1TKB	Gallbladder	Gallbladder cells	Adherent
TGBC24TKB	Gallbladder	Gallbladder cells	Adherent
TGBC24TKB	Gallbladder	Gallbladder cells	Adherent
TGBC2TKB	Gallbladder	Gallbladder cells	Adherent
TGBC51TKB	Bile ducts (Ampulla of Vater)	Bile duct epithelial cells	Adherent
TGW	Nervous system	Neuroblast	Adherent
Th17-DP cells	Blood	T-helper 17 (Th17) cells	unknown
THLE-2	Liver	Liver epithelial cells	Adherent
THLE-3	Liver	Liver epithelial cells	Adherent
THP-1	Peripheral blood	Monocytes	Suspension
Thymocytes	Thymus	Thymocyte	unknown
thymus cells mouse	Thymus	Thymocyte	unknown
TI-73	Bone	Osteocyte	Adherent
TK	unknown	unknown	unknown
TK-10	Kidney	Kidney	unknown
TK-6	Peripheral blood	Myeloid cells	Suspension
TKKK	Liver	Cholangiocyte	Adherent
T-lymphoblastoid cells	Peripheral blood	T-lymphocyte	unknown
TMK-1	Stomach	Stomach cells	unknown
TNBC	unknown	unknown	unknown
TO 175.T	Lymphatic system	Reed-sternberg cells	Adherent
TO 203.T	Bone	Osteocyte	unknown
Toledo	Lymphatic system	B-lymphocyte	Suspension
TOV-112D	Ovary	Ovarian epithelial cell	Adherent
TOV-21G	Ovary	Ovarian epithelial cell	Adherent
TPC1	Thyroid	Thyroid epithelial cells	unknown
TRIP13	unknown	unknown	unknown
Trophoblasts	Embryo	Trophoblast	unknown
TS1/18.1.2.11	unknown	B-lymphocyte	Suspension
TSC	Embryo	Trophoblast stem cells	unknown
TSC1-/-MEF	Embryo	Trophoblast stem cells	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
TSC1+/-MEF	Embryo	Trophoblast stem cells	unknown
TSC2-/-MEF	Embryo	Trophoblast stem cells	unknown
TSC2+/-MEF	Embryo	Trophoblast stem cells	unknown
TSCC15 cells	unknown	unknown	unknown
TSCCD	unknown	unknown	unknown
TSCM Cells	Peripheral blood	Stem memory T cells	unknown
TS-ELT-3	unknown	Smooth muscle cells	unknown
TT	Thyroid	Thyroid cells	Adherent
TT2609-C02	Thyroid	Thyroid cells	Adherent
TUR	Peripheral blood	Histiocyte	Suspension
TYK-nu	Ovary	Ovarian cells	Adherent
U031	Kidney	Kidney cells	unknown
U-118 MG	Brain	Astrocyte	Adherent
U-138 MG	Brain	Astrocyte	Adherent
U-2 OS	Bone	Osteocyte	Adherent
U251	Brain	Astrocyte	Adherent
U-251 MG	Brain	Astrocyte	Adherent
U251-BM	unknown	unknown	unknown
U251-P	unknown	unknown	unknown
U251-sh1MSH6-BM	unknown	unknown	unknown
U251-shMSH6-BM	unknown	unknown	unknown
U251-shMSH6-P	unknown	unknown	unknown
U251T	Brain	Astrocyte	Adherent
U266	Peripheral blood	Plasma cell	unknown
U266B1	Peripheral blood	Plasma cell	Suspension
U373 MG	Brain	Astrocyte	Adherent
U-698-M	Lymphatic system	B-lymphocyte	Suspension
U-87 MG	Brain	Astrocyte	Adherent
U-937	Peripheral blood	Monocytes	Suspension
UACC-257	Skin	Melanocytes	unknown
UACC-62	Skin	Melanocytes	Adherent
UACC-812	Breast	Breast epithelial	Adherent
UACC-893	Breast	Breast epithelial	Adherent
UACC903	Skin	Melanocytes	unknown
UDSCC2	Pharynx (lower)	Pharyngeal squamous cells	Adherent
UE6E7T-2	Bone marrow	Mesenchymal stem cells	Adherent
UE6E7T-3	Bone marrow	Mesenchymal stem cells	Adherent
UE6E7T-3	Skin	Merkel cell	Adherent
UMC-11	Lung	Lung cells	Adherent
UM-UC-3	Urinary tract	Urinary transitional cell	Adherent
UO-31	Kidney	Kidney cells	unknown
UOK 257	Kidney	Kidney cells	Adherent

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
UOK262	Kidney	Kidney cells	unknown
UT-7	Peripheral blood	Myeloid cells	Suspension
UTEX 297	unknown	unknown	unknown
UtLM	unknown	unknown	unknown
UTSW-ACC112	Salivary gland	Salivary gland cells	unknown
Uveal Melanoma (UM) cells	Iris	Uveal melanocytes	unknown
V1	unknown	Plasma cell	unknown
V79	Lung	Fibroblast	Adherent
VA-ES-BJ	Skin	Mesenchymal cells	Adherent
VAL	Lymphatic system	B-lymphocyte	Suspension
Vascular Smooth Muscle Cells VSMC	Circulatory system	Vascular smooth muscle cells	unknown
VCaP	Prostate	Prostate cells	Adherent
Vero	Kidney	Kidney cells	Adherent
VK-1:NS	unknown	Ameoba	unknown
VM-CUB1	Urinary tract	Urinary transitional cell	Adherent
VMRC-LCD	Lung	Lung cells	Adherent
VMRC-LCP	Lung	Lung squamous cells	Adherent
VMRC-MELG	Skin	Melanocytes	Adherent
VMRC-RCW	Kidney	Kidney cells	Adherent
VMRC-RCZ	Kidney	Kidney cells	Adherent
VP16	unknown	unknown	unknown
Vx-2	unknown	unknown	unknown
WB-f344	Liver	Liver epithelial cells	unknown
WERI-RB-1	Retina	Immature retinal cells	Suspension
WHO VERO cell line	Kidney	Kidney cells	Adherent
Whole blood	Blood	Whole blood	unknown
WI 38	Lung (fetus)	Lung fibroblast (fetus)	Adherent
WI-26 VA4	Lung (fetus)	Lung fibroblast (fetus)	Adherent
WI-38 VA13 sub 2 RA	Lung (fetus)	Lung fibroblast (fetus)	Adherent
WI-38-40	Lung (fetus)	Lung fibroblast (fetus)	Adherent
WiDr	Colon	Colorectal epithelial cells	Adherent
WIL2 NS	Lymphatic system	B-lymphocyte	Suspension
WIL2.NS.6TG	Lymphatic system	B-lymphocyte	Suspension
WILCL	Lung	Lung squamous cells	Suspension
WM 266-4	Skin	Melanocytes	unknown
WM-115	Skin	Melanocytes	unknown
WM1158	Skin	Melanocytes	unknown
WM1552C	Skin	Melanocytes	unknown
WM164	Skin	Melanocytes	unknown
WM239A	Skin	Melanocytes	unknown
WM278	Skin	Melanocytes	unknown

Cell type	Anatomical origin (organ)	Anatomical origin (cell type)	Adherent or suspension
WM35	Skin	Melanocytes	unknown
WM793B	Skin	Melanocytes	unknown
WM852	Skin	Melanocytes	unknown
WM902B	Skin	Melanocytes	unknown
WPE1-NA22	Prostate	Prostate epithelial cells	Adherent
WPE1-NB11	Prostate	Prostate epithelial cells	Adherent
WPE1-NB14	Prostate	Prostate epithelial cells	Adherent
WPE1-NB26	Prostate	Prostate epithelial cells	Adherent
WSU-DLCL2	Lymphatic system	B-lymphocyte	Suspension
WSU-NHL	Lymphatic system	B-lymphocyte	Suspension
XPA-1	Pancreas	Pancreas cell	unknown
XPA-3	Pancreas	Pancreas cell	unknown
XPA-4	Pancreas	Pancreas cell	unknown
Y. lipolytica	unknown	Fungus	unknown
Y79	Retina	Immature retinal cells	Suspension
YAPC	Pancreas	Pancreas cell	Adherent
Yersinia pestis	unknown	Bacteria	unknown
YH-13	Brain	Astrocyte	Adherent
YKG-1	Brain	Astrocyte	Adherent
YMB-1-E	Breast	Breast epithelial	Adherent
YSCCC	Bile duct	Cholangiocyte	Adherent
YT	Peripheral blood	Natural killer cells	Suspension
Zebrafish tumor cells	unknown	unknown	unknown
ZR-75-1	Breast	Mammary ductal epithelial cells	Adherent
ZR-75-30	Breast	Mammary ductal epithelial cells	Adherent

