

No.	Cell Line Name	Cancer Type	Tissue Origin	
1	A-673	Bone Osteosarcoma	Bone	
2	HOS			
3	MG-63			Ewing's Sarcoma
4	U2OS			Bone Carcinoma
5	BT474	Ductal Carcinoma	Breast	
6	MCF-7	Adenocarcinoma		
7	MDA-MB-231	Adenocarcinoma, TNBC		
8	MDA-MB-361	Breast Cancer		
9	MDA-MB-468	Breast Cancer, TNBC		
10	SK-BR-3	Adenocarcinoma		
11	T47D	Ductal Carcinoma		
12	ZR-75-1	Breast Cancer	Cervix	
13	HeLa	Adenocarcinoma		
14	C-33A	Carcinoma		
15	Ca Ski	Epidermoid Carcinoma		
16	SiHa	Squamous Cell Carcinoma		
17	HT 1080	Connective Tissue Carcinoma		
18	93T449	Retroperitoneal Liposarcoma		
19	RD	Muscle Rhabdomyosarcoma	GI Tissue	
20	5637	Bladder Cancer		
21	COLO 205	Colorectal Adenocarcinoma		
22	DLD-1	Colon Cancer		
23	HCT 116	Colorectal Carcinoma		
24	HT-29	Colon Cancer		
25	MIA PaCa-2	Pancreatic Carcinoma		
26	NCI-N87	Gastric Carcinoma		
27	PANC-1	Pancreatic Carcinoma		
28	SNU-16	Stomach Cancer		
29	SW480	Colorectal Adenocarcinoma		
30	T-24	Transitional Cell Carcinoma		
31	T84	Colorectal Carcinoma	Head and Neck	
32	A172	Glioblastoma		
33	CAL 27	Squamous Cell Carcinoma		
34	Detroit 562	Pharyngeal Carcinoma		
35	IMR-32	Brain Neuroblastoma		
36	SCC-25	Squamous Cell Carcinoma		
37	SCC-9			
38	T98G	Brain Glioblastoma		
39	U87MG			
40	Hep G2	Hepatocellular Carcinoma	Liver	
41	Hep 3B			
42	PLC/PRF/5			Hepatoma
43	SK-HEP-1			Liver Cancer
44	SNU-182			
45	A-549	Lung Cancer	Lung	
46	HCC78	Adenocarcinoma		
47	MSTO-211H	Mesothelioma		
48	NCI-H441	Adenocarcinoma		
49	NCI-H460	Carcinoma		
50	NCI-H1975	Adenocarcinoma		

51	BV-173	B cell lymphoma	Blood		
52	CCRF-CEM	Acute lymphoblastic leukemia			
53	H-9	T cell lymphoma (derivative of HuT-78)			
54	HH	T cell lymphoma			
55	HL-60	acute promyelocytic leukemia			
56	HuT 102	T cell lymphoma			
57	HuT 78				
58	Jurkat	acute T cell leukemia			
59	K-562	Leukemia (CML)			
60	KASUMI-2	B cell lymphoma			
61	MJ (G11)	Lymphoma			
62	MOLT-4	acute lymphoblastic leukemia			
63	MV-4-11	B myelomonocytic leukemia			
64	NOMO-1	Acute Myeloid Leukemia			
65	Raji	Burkitt's lymphoma			
66	REH	B cell lymphoma			
67	SU-DHL-1	ALCL lymphoma			
68	SU-DHL-6	B cell lymphoma			
69	SUP-T1	T cell lymphoma			
70	SZ-4	Helper T cell lymphoma			
71	KMS-11	Multiple myeloma			
72	MM.1S	Immunoglobulin A lambda myeloma			
73	RPMI 8226	Plasmacytoma, myeloma			
74	U266	Myeloma, plasmacytoma			
75	A2780	Carcinoma		Ovary	
76	ES-2				Clear Cell Carcinoma
77	SK-OV-3				Adenocarcinoma
78	OVCAR-3				
79	22Rv1	Prostate Cancer	Prostate		
80	DU 145				
81	LNCap				
82	PC-3				
83	VCaP	Melanoma	Skin		
84	A2058				
85	A-375			Malignant melanoma	
86	A375-P				
87	FaDu			Squamous cell carcinoma	
88	SK-MEL-2			Malignant melanoma	
89	TE354.T			Basal cell carcinoma	
90	DU4475			Skin carcinoma	
91	UACC-62	Malignant melanoma			
92	A-431	Epidermoid carcinoma	Kidney		
93	769-P	Renal Cell Adenocarcinoma			
94	786-O				
95	A-498	Kidney carcinoma			
96	ACHN	Renal Cell Adenocarcinoma	Uterus		
97	AN3 CA	Adenocarcinoma			
98	HEC-1-A				
99	KLE				
100	RL95-2	Carcinoma			