

List of Cell Lines for Cell Proliferation Assays



No.	Cell specification	Entity of origin	No.	Cell specification	Entity of origin	No.	Cell specification	Entity of origin	No.	Cell specification	Entity of origin		
1	J82	Bladder	37	HCC 1569	Breast	73	Hep3B2.1-7	Liver	109	OVCAR-3	Ovary		
2	HL-60	Blood	38	HCC38		74	HepG2		Lung	110		OVK18	
3	K562		39	JIMT-1		75	A427	111		RL95-2			
4	KARPAS 299		40	MCF-7		76	A549	112		SK-OV3			
5	KG1		41	MDA MB 231		77	BEN	113		SNU840			
6	KG1a		42	MDA MB 468		78	Calu-6	114		AsPC-1	Pancreas		
7	KMS-12-BM		43	SK.BR-3		79	COR-L279	115		BxPC3			
8	LP-1		44	T47D		80	DV90	116		Mia PaCA2			
9	M07e		45	ZR-75-1		46	HCC202	81		EPLC-272H	117	PANC-1	Prostate
10	Molm-13		47	C33a		82	H1299	83		H460	118	DU-145	
11	MOLT4		48	HeLa	49	SiHa	84	HCC827		119	LnCap		
12	MV4-11		49	SiHa	50	Colo 205	85	LOU-NH91	120	PC3	Skin		
13	NALM-6	51	CX-1	52	DLD-1	86	N417	121	A2058				
14	OCI-AML5	53	HCT-116	54	HCT-15	87	NCI-H292	122	A375				
15	OPM-2	55	HT-29	56	LOVO	88	NCI-H441	123	A431				
16	P31/FUJ	57	LS174T	58	RKO	89	NCI-H838	124	MDA MB 435	Stomach			
17	RPMI 8226	59	SW480	60	SW620	90	NCI-H1048	125	SK-MEL-3				
18	RS4 11	61	SW948	62	T-84	91	NCI-H1437	126	Hs 746T				
19	SU-DHL5	63	Caco-2	64	Colo-320DM	92	NCI-H1563	127	MKN-1				
20	U-266	65	Hutu 80	66	HEC-1-A	93	NCI-H1573	128	MKN-45	Stomach			
21	U-937	66	HEC-1-A	67	HEC-1-B	94	NCI-H1581	129	NCI-N87				
22	WSU-NHL	68	Ichikawa	69	HT-1080	95	NCI-H1703	130	SCH				
23	Saos-2 EC	69	HT-1080	70	A498	96	NCI-H1838	131	SNU-1				
24	SJSA-1	70	A498	71	Caki-1	97	NCI-H2009			Lung			
25	SK-ES-1	71	Caki-1	72	Caki-2	98	NCI-H2110						
26	U2OS	72	Caki-2			99	NCI-H2286						
27	A172					100	RERF-LC-Ad2						
28	H4					101	RERF-LC-MS						
29	LN229					102	SCLC-21H						
30	SK-N-FI					103	SK-LU-1						
31	SK-N-MC					104	COV434						
32	SK-N-SH					105	EFO-27						
33	U118MG					106	MCAS						
34	U87MG					107	NCI-ADR						
35	BT-20					108	OV56						
36	EFM-192A												

Cell Lines