

List of Cell Lines for Soft Agar 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	HCT-15	Colon	73	NCI-H69	Lung				
2	HL-60	Blood	38	HT-29		74	NCI-H838					
3	Karpas 299		39	Hutu 80		75	NCI-H889					
4	MOLM-13		40	LOVO		76	RERF-LC-MS					
5	P31/FUJ		41	LS174T		77	SCLC-21H					
6	RPMI 8226		42	RKO		78	SK-LU-1					
7	U-937		43	SW480		79	COV434	Ovary				
8	SJSA-1		44	SW620		80	EFO-27					
9	SK-ES-1	Bone	45	SW948		81	MCAS					
10	LN229	Brain	46	T84		82	NCI-ADR					
11	SK-N-FI		47	HEC-1-A	83	OV56						
12	SK-N-MC		48	HEC-1-B	84	OVK18						
13	SK-N-SH		49	HT-1080	85	RL95-2						
14	U118MG		50	A498	86	SK-OV3						
15	U87MG	Kidney	51	Caki-2	87	AsPC-1	Pancreas					
16	CAL-148		52	HepG2	88	Mia PaCA 2						
17	EFM-19		Liver	53	A427	89	DU-145	Prostate				
18	EFM-192A			54	A549	90	LnCap					
19	HCC1569			55	BEN	91	PC3					
20	HCC38			56	Calu-6	92	A2058	Skin				
21	KPL-1			57	CORL279	93	A375					
22	HCC 1569			58	DV90	94	MDA MB 435					
23	JIMT-1			59	H1299	95	SK-MEL-3					
24	MCF-7			60	H460	96	Hs 746T		Stomach			
25	MDA MB 231	61		HCC827	97	MKN-1						
26	MDA MB 468	62		N417	98	MKN-45						
27	SK.BR-3	63	NCI-H1048	99	NCI-N87							
28	T47D	64	NCI-H1299	100	SCH							
29	ZR-75-1	65	NCI-H1437	101	SNU-1							
30	C33a	Cervix	66	NCI-H1573								
31	HeLa		67	NCI-H1581								
32	SiHa		68	NCI-H1703								
33	Colo 205	Colon	69	NCI-H2009								
34	CX-1		70	NCI-H2106								
35	DLD-1		71	NCI-H2110								
36	HCT116		72	NCI-H441								

Cell Lines