

List of Cell Lines for Cell Migration 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									
2	SJSA-1	Bone									
3	LN299	Brain									
4	SK-N-SH										
5	U118MG										
6	U87MG										
7	MCF-7	Breast									
8	MDA-MB 231										
9	HeLa	Cervix									
10	DLD-1	Colon									
11	HCT116										
12	RKO										
13	HT-1080	Fibrosarcoma									
14	A498	Kidney									
15	Caki-2	Kidney									
16	A427	Lung									
17	A549										
18	Calu-6										
19	H1299										
20	HCC827										
21	MCAS	Ovary									
22	SK-OV3										
23	DU-145	Prostate									
24	PC3										
25	A375	Skin									
26	A431	Stomach									
27	Hs746T										
28	MKN-1										

 Cell Lines