

Field of Application

Subcutaneous mouse tumor models are widely used in preclinical drug development for measurement of antitumoral efficacy in a standardized and cost-effective manner. The use of immune-compromised mice offers the advantage to study effects on human tumor cell lines (xenografts, CDXs). In contrast, murine tumor cell lines can be grown in immune-competent mice (syngeneic models). These syngeneic tumor models are advantageous providing a functional immune system to assess novel immunotherapeutic approaches.

Our Service

➤ Established Cell Lines

#	Tumor Cell Line	Tissue Origin	Species	#	Tumor Cell Line	Tissue Origin	Species	
1	HG3**	Blood/ Leukemia	Human	24	A549	Lung	Human	
2	HL-60		Human	25	Calu-3**		Human	
3	KARPAS 299		Human	26	Calu-6		Human	
4	MOLM-13		Human	27	H460		Human	
5	MOLT-4**		Human	28	LL/2		Human	
6	MV4-11		Human	29	NCI-H1437**		Human	
7	SJSA-1	Bone	Human	30	NCI-H292**		Human	
8	LN-229	Brain	Human	31	NCI-H441**		Human	
9	U-87 MG		Human	32	SKOV-3		Ovary	Human
10	4T1	Breast	Mouse	33	AsPC1		Pancreas	Human
11	MCF7		Human	34	BxPC-3	Human		
12	MDA-MB 231		Human	35	MiA-PaCa2	Human		
13	HeLa	Cervix	Human	36	LNCaP (Z2)*	Prostate	Human	
14	Colo-205	Colon	Mouse	37	PC-3		Human	
15	COLO-320DM**		Human	38	A2058	Skin	Human	
16	CT26 wt		Mouse	39	A375		Human	
17	HCT-116		Human	40	B16-F10		Mouse	
18	HT29		Mouse	41	C8161**		Human	
19	LS 174T		Human	42	Clone M3		Mouse	
20	MC38-CEA		Mouse	43	HT-144**		Human	
21	SW620		Human	44	Hs746T		Stomach	Human
22	HuTu80		Duodenum	45	MKN-1			Human
23	RENCA		Kidney	46	MKN-45			Human

* in vivo selected subpopulation of corresponding parental cell line

** model under development

Models with Genetically Engineered Cells

#	Tumor Cell Line	Target	Species
1	MEF-IGF1-Rrep*	IGF1 Receptor	Mouse
2	NIH3T3-ErbB2-Rrep*	ErbB2 Recepto	Mouse

➤ Customer Cell Lines

Studies can be performed with cell lines provided by our customers.

➤ Reference Compound

Available for most established cell lines.

➤ Standard Study

Comprises among other things: (i) cell culturing & cell implantation; (ii) measurement of animal weight (up to 3x/week); (iii) determination of tumor size (2x/week); (iv) treatment (1x/day; 5x/week); treatment period depending on model; (v) protocol & report.

➤ Optional Services

(i) Blood sampling; (ii) paraffin embedding of tumor tissue or organs; (iii) histology & pathology; (iv) Magnetic resonance tomography and (v) PET/CT.