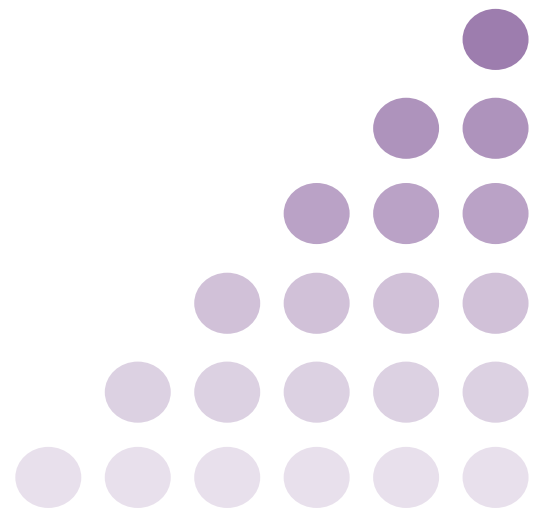

In Vivo Pharmacology

Orthotopic Mouse Tumor Models	3
Subcutaneous and SubQperior™ Mouse Tumor Models	4
Metastasis Mouse Tumor Models	6
In Vivo Hollow Fiber Model	7
Mouse-derived Isograft Tumor Models (MDI)	8
Immuno-Oncology Platform	9

Reaction Biology has established a large panel of subcutaneous, subQperior™, orthotopic and metastasis tumor models (including bioluminescence read-out) for testing the efficacy of novel therapeutic agents on primary tumors and metastases in animals. Moreover, standard and proprietary syngeneic tumor models are available, which enables efficacy testing of test compounds in immune competent animals and the investigation of the immune-modulatory effect (e.g. by flow cytometry). Tumor angiogenesis can be investigated in In vivo Angiogenesis Models. Reaction Biology's In Vivo Hollow Fiber Model allows to screen the most suitable cell line or the most promising antitumoral effective lead test compound in a cost and time reduced manner in mice. The service portfolio is supplemented by other study types (e.g. tolerability, PK, PD). In addition, new models can be established upon request.

Let's discover together.



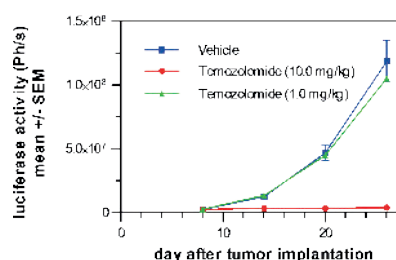
Orthotopic Mouse Tumor Models

Overview

In orthotopic tumor models, tumor cells are inoculated into the organ of their origin. This implantation site allows organo-typical interaction between tumor cells and surrounding stroma affecting growth, differentiation, and drug sensitivity of tumor cells. For human tumor cell lines immune-compromised mice are used with the advantage to study classical antitumoral test compounds. In contrast, murine tumor cell lines can be grown in immune-competent mice (syngeneic). These tumor models combine the advantage of tumor stroma interaction with a functional developed immune system to assess novel immunotherapeutic approaches.

Our Service

- Customer Cell Lines** Orthotopic tumor models can be established with cell lines provided by our customers
- Analysis of Tumor Size** Via bioluminescence imaging (BLI) using luciferase-transduced cell lines or calipering of skin and breast tumors
- Reference Cpd** 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 via BLI (1x/week) / calipering (2x/week); (v) treatment (1x/day; 5x/week); (vi) treatment period depending on model (vi) protocol & report
- Optional Services** (i) Blood sampling; (ii) paraffin embedding of tumor tissue or organs; (iii) histology & pathology (iv) MRT (v) flow cytometry



Study Example

Effect of Temozolomide (high dose = red, low dose = green) on orthotopic tumor growth of U-87 MG cells (vehicle control = blue) determined by in vivo bioluminescence imaging

Available Models

#	Tumor Cell Line	Tissue Origin	Species
1	MB-49	Bladder	Mouse
2	A172	Brain	Human
3	GL261		Mouse
4	LN-229		Human
5	U-87 MG		Human
6	4T1	Breast	Mouse
7	BT-474		Human
8	EMT6		Mouse
9	EO771		Mouse
10	HCC-1569		Human
11	JIMT-1		Human
12	MCF7		Human
13	MDA-MB-231 (Z1)*		Human
14	MDA-MB-468	Human	
15	CT26 wt	Colon	Mouse
16	MC38-CEA		Mouse
17	HCT-116		Human
18	SW620		Human

#	Tumor Cell Line	Tissue Origin	Species
19	RENCA	Kidney	Mouse
20	786-O		Human
21	Hep 3B2.1-7	Liver	Human
22	HepG2		Human
23	Hepa1-6		Mouse
24	PLC/PRF/5		Human
25	LL/2	Lung	Mouse
26	A2780	Ovary	Human
27	OVCAR-3		Human
28	SKOV-3		Human
29	AsPC1	Pancreas	Human
30	MiA-PaCa2		Human
31	Pan02		Mouse
32	PANC-1		Human
33	PC-3	Prostate	Human
34	B16-F10	Skin	Mouse
35	CB161		Human
36	CB161.9		Human

* in vivo selected subpopulation of the corresponding parental cell line

Subcutaneous and SubQperior™ Mouse Tumor Models

Overview

Subcutaneous mouse tumor models are widely used in preclinical drug development for measurement of antitumoral efficacy in a standardized and cost-effective manner.

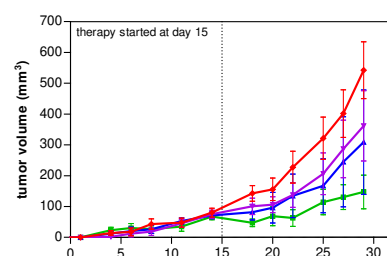
SubQperior™ tumor models are an innovative alternative overcoming many of the drawbacks of tumor models based on subcutaneous tumor cell implantation. Standard implantation via subcutaneous injection causes frequent ulceration resulting in early abrogation of studies, leaving researchers with too short treatment windows and high heterogeneity with poor statistical value of study results. SubQperior™ tumor models are based on tumor cell implantation into the mammary fat pad.

Advantages of SubQperior™ tumor models:

- Implantation results in larger tumors and longer treatment windows
- Homogeneous and reliable growth yields study outcomes with outstanding statistical value
- Tumors are measured via caliper, making the handling as easy and inexpensive as for subcutaneous models

Our Service

Established Cell Lines	Xenograft models (human cells); syngeneic models (murine cells)
Customer Cell Lines	Studies can be performed with cell lines provided by our customers
Reference Cpd	Available for most established cell lines
Standard Study	Comprises among other things: <ul style="list-style-type: none"> (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	<ul style="list-style-type: none"> (i) blood sampling; (ii) paraffin embedding of tumor tissue or organs; (iii) histology & pathology; (iv) MRT; (v) Flow cytometry



Available Models

Human

Tumor Cell Line	Tissue Origin	Route of Application	Tumor Cell Line	Tissue Origin	Route of Application
Ramos	Ascites	metastasis (i.v.)	A549		subcutaneous; metastasis (i.v.)
UM-UC-3	Bladder	orthotopic	Calu-3		subcutaneous
HG-3		subcutaneous	Calu-6		subcutaneous
HL-60		subcutaneous	EBC1		subcutaneous
KARPAS 299		subcutaneous	LK-2		subcutaneous
LP-1		orthotopic	NCI-H1373	Lung	subcutaneous
MOLM-13	Blood	subcutaneous; metastasis (i.v.)	NCI-H1437		subcutaneous
MOLT-4		subcutaneous	NCI-H292		subcutaneous
MV4-11		subcutaneous; metastasis (i.v.)	NCI-H441		subcutaneous/subQperior
RPML-8226		orthotopic	NCI-H460		subcutaneous; metastasis (i.c.)
SU-DHL-6		subcutaneous	PC-9		subcutaneous
THP-1		subcutaneous	CAL-27	Oral	subcutaneous
SJSA-1	Bone	subcutaneous	A2780		subcutaneous; orthotopic
NCI-H929		subcutaneous; orthotopic	Hey		subcutaneous
Raji	Bone Marrow	subcutaneous; metastasis (i.v.)	OVCAR-3	Ovary	subcutaneous; orthotopic
A172		orthotopic	SKOV-3		subcutaneous; orthotopic; orthotopic ascites
LN-229		subcutaneous; orthotopic	AsPC1		subcutaneous; subcutaneous/subQperior; orthotopic
U-118 MG	Brain	subcutaneous	BxPC-3		subcutaneous
U-87 MG		subcutaneous; orthotopic	HuPT4	Pancreas	subcutaneous
BT-474		subcutaneous; orthotopic	MiAPaCo2		subcutaneous; orthotopic
HCC1187		subcutaneous	PANC-1		subcutaneous; orthotopic
HCC1569		orthotopic	Stew		subcutaneous
JIMT1		orthotopic; metastasis (i.c.)	Detroit 562	Pharynx	subcutaneous
MCF7	Breast	subcutaneous; orthotopic	PC-3	Prostate	subcutaneous; orthotopic
MDA-MB-231		subcutaneous; orthotopic; metastasis (i.c.); metastasis (i.v.)	PC3-M		subcutaneous
MDA-MB-468		orthotopic	A2058		subcutaneous
COLO 201		subcutaneous	A375		subcutaneous
COLO 205		subcutaneous	A431		subcutaneous; subcutaneous/subQperior
COLO 320DM		subcutaneous	C32		subcutaneous
DLD-1		subcutaneous	C8161	Skin	subcutaneous; orthotopic
GP2d		subcutaneous	C8161.9		orthotopic
HCT 116		subcutaneous; orthotopic; metastasis	G361		orthotopic
HT29	Colon	subcutaneous	HT144		subcutaneous
LoVo		metastasis	UACC-257		subcutaneous
LS 174T		subcutaneous	Hs746T		subcutaneous
RKO		subcutaneous	MKN-1	Stomach	subcutaneous
SW480		subcutaneous	MKN-45		subcutaneous
SW620		subcutaneous/subQperior; orthotopic	HeLa	Uterus, Cervix	subcutaneous
WiDr		subcutaneous	SK-LMS-1	Vulva	subcutaneous
HT1080	Connective Tissue	subQperior			
HuTu80	Duodenum	subcutaneous			
786-O		subcutaneous; orthotopic			
ACHN	Kidney	subcutaneous			
Hep3B2.1-7		subcutaneous; orthotopic			
HepG2		orthotopic			
Huh-7	Liver	subcutaneous; metastasis (i.v.); metastasis (i.c.)			
PLC/PRF/5		subcutaneous; orthotopic			
SNU-182		subcutaneous			

Mouse

Tumor Cell Line	Tissue Origin	Route of Application
MB-49	Bladder	subcutaneous; orthotopic; metastasis (i.c.); metastasis (i.v.)
MBT-2		subcutaneous
A20		subcutaneous; subQperior
C1498	Blood	subcutaneous; metastasis (i.v.)
E.G7-OVA		subcutaneous
EL4		subcutaneous
GL261	Brain	orthotopic; subQperior
4T1		subcutaneous; orthotopic; subQperior_orthotopic; metastasis (orthotopic); metastasis (i.v.)
EMT6	Breast	subcutaneous; orthotopic; subQperior_orthotopic; metastasis
EO771		orthotopic
HC11-NeuT		orthotopic
C26		subcutaneous
CT26wt	Colon	subcutaneous; subQperior; orthotopic
MC38		subcutaneous
MC38-CEA#		subcutaneous; subQperior; orthotopic
RENCA	Kidney	subcutaneous; subQperior; orthotopic
Hepa1-6	Liver	subcutaneous; subQperior; orthotopic
3LL		subcutaneous
AB12		subcutaneous; subQperior
LL/2 (LLC1)	Lung	subcutaneous, subQperior; orthotopic; metastasis
M109		subcutaneous
TC-1		subcutaneous
KPC-2838		subcutaneous
KPC-6419	Pancreas	subcutaneous
Pan02		subcutaneous, subQperior; orthotopic
B16-F0		subcutaneous; subQperior; orthotopic
B16-F10	Skin	subcutaneous; subQperior; orthotopic; metastasis (i.c.); metastasis (i.v.)
Clone M3/Cloudman S91		subcutaneous, subQperior

Metastasis Mouse Tumor Models

Overview

In metastatic tumor models, tumor spreading originates either from a primary tumor or is artificially induced by intravenous or intracardial tumor cell injection. The metastatic pattern is dependent on the tumor cell line with a preference for lung metastasis in the case of intravenous injection. For human tumor cell lines immune-compromised mice are used with the advantage to study classical antitumoral test compounds. In contrast, murine tumor cell lines can be grown in immune-competent mice (syngeneic), providing a functional immune system to assess novel immunotherapeutic approaches.

Our Service

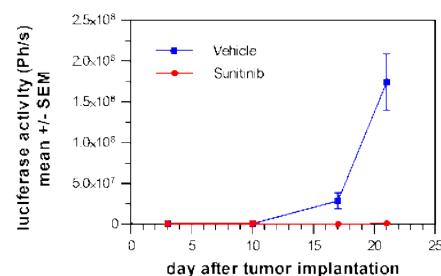
Customer Cell Lines Metastatic tumor models can be established with cell lines provided by our customers

Analysis of Metastasis Via bioluminescence imaging (BLI) using luciferase-transduced cell lines

Reference Cpd 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 burden via BLI (1x/week); (v) treatment (1x/day; 5x/week); treatment period depending on model; (vi) protocol & report

Optional Services (i) Blood sampling; (ii) paraffin embedding of organs; (iii) histology & pathology (iv) MRT (v) flow cytometry (vi) ex vivo analysis of organs



Available Models

#	Tumor Cell Line	Tissue Origin	Route of application	Species
1	Ramos	Ascites	i.v.	Human
2	MB-49	Bladder	i.c., i.v.	Mouse
3	C1498		i.v.	Mouse
4	MOLM-13	Blood / Leukemia	i.v.	Human
5	MV4-11		i.v.	Human
6	Raji	Bone Marrow	i.v.	Human
7	4T1 (M3)*		orthotopic	Mouse
8	4T1		i.v.	Mouse
9	EMT6	Breast	i.v.	Mouse
10	JIMT-1		i.c.	Human
11	MDA-MB-231 (Z1)*		i.c., i.v.	Human
12	HCT 116		orthotopic	Human
13	LoVo	Colon	orthotopic	Human
14	Huh-7	Liver	i.c., i.v.	Human
15	A549		i.v.	Human
16	LL/2	Lung	intra-splenic	Mouse
17	NCI-H460		i.c.	Human
18	B16-F10	Skin	i.c., i.v.	Mouse

* in vivo selected subpopulation of the corresponding parental cell line

In Vivo Hollow Fiber Model

Overview

The In Vivo Hollow Fiber Model is a fast and economical in vivo screening approach which can be used for two different purposes:

- Selection of a compound with the best in vivo activity against a tumor cell line of interest (“compound screening”)
- Identification of the most suitable cell line for an in vivo efficacy study with a selected test compound (“tumor model screening”)

The method allows simultaneous evaluation of test items against up to three different tumor cell lines in two different compartments (s.c. & i.p. implantation) within the same mouse. Due to the low variability of the assay, studies can be performed with three mice per group. These characteristics predestine the Hollow Fiber Model for drug or tumor model screening in vivo.

Our Service

Typical Examples of Custom-Tailored Projects:

Objective	Short Description of a Standard Study
In vivo compound screening	Testing of up to 14 compounds in one study. Cell culturing; encapsulation of tumor cell line of interest in hollow fibers; subcutaneous & intraperitoneal implantation of hollow fibers; in vivo study with treatment (study duration of 14 days); hollow fiber harvesting; measurement of tumor cell viability by CellTiter-Glo; protocol & report
In vivo tumor model screening	Testing of up to 12 tumor cell lines in one study. Cell culturing; encapsulation of three different cell types in hollow fibers; subcutaneous & intraperitoneal implantation of hollow fibers; in vivo study with treatment (study duration of 14 days); hollow fiber harvesting; measurement of tumor cell viability by CellTiter-Glo; protocol & report

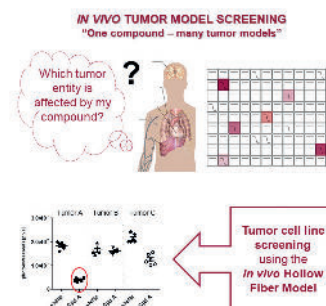
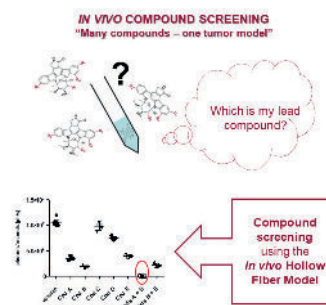
The Test Model
The key element of the assay are semi-permeable fibers which allow access of test items (<500 kDa: small molecules, antibodies etc.) to the encapsulated tumor cells

First Description in the Literature
Hollingshead et al (1995), Life Sciences 57, pp. 131 - 141

Available Cell Lines
Cell lines can be selected from Reaction Biology’s cell lines or from our in vivo panel of established subcutaneous or orthotopic tumor models

Customer Cell Lines
Studies can also be performed with cell lines provided by our customers

Readout
Level of tumor cells in the hollow fibers are quantified by CellTiter-Glo



Mouse-derived Isograft Tumor Models (MDI)

Overview

Today, cancer research tends to focus on the development of novel cancer immunotherapies using syngeneic mouse models for the development of such compounds. While classical syngeneic mouse models are based on the implantation of cultured tumor cell lines, for Reaction Biology's mouse-derived isograft (MDI) tumor models are propagated subcutaneously in mice only. Implanted tumor pieces of low in vivo passage are originally derived from spontaneous- or carcinogen-induced mouse tumors. Hence, the major advantage of these novel and unique MDI tumor models is the preservation of primary tumor phenotype and intratumoral immune cell populations.

Our Service

Characterization

- All models were histologically characterized
- Growth curves were established for all models
- All models were tested with reference compounds
- RNA sequence data available for selected models

Flow Cytometric Analysis

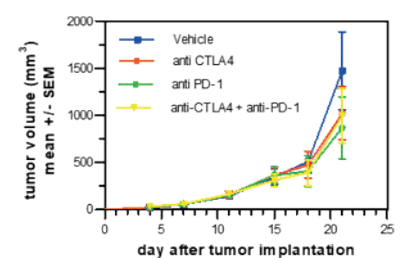
- Analysis of tumor infiltrating leukocytes and cells isolated from spleen and/or lymph nodes
- Multicolor 17 marker panel: T cells, MDSCs, Macrophages, NK cells, B cells, DCs
- Customized staining procedures are possible

Standard Study

- Comprises among other things:
- Subcutaneous implantation of tumor pieces
 - Measurement of animal weight (3x/week)
 - Determination of tumor size

Optional Services

- Blood sampling
- Characterization of abundance and relative distribution of different immune cell subsets in the tumor and lymphatic tissues by flow cytometry
- Paraffin embedding of tumor tissue
- Histological & pathological analysis
- Cytokine determination
- Provision of tumor tissue for target validation
- MRT



Available Models

MDI tumor models:

#	Model	Origin	Creation	RNAseq	anti-PD1	anti-CTLA-4	Combination	Therapeutic window	Relevant immune cells
1	JA-0009	adeno-carcinoma	spontaneous	yes	low	low	low	2 weeks	M2 macrophages
2	JA-0017		spontaneous	ongoing	nd	nd	nd	>6 weeks	nd
3	JA-0032		spontaneous	ongoing	nd	nd	nd	4 weeks	nd
4	JA-2011	sarcoma	carcinogen	yes	low	low	nd	2 weeks	neutrophils
5	JA-2019		carcinogen	ongoing	high	high	nd	2 weeks	MDSCs/Tcells
6	JA-2041		carcinogen	ongoing	moderate	moderate	nd	3 weeks	MDSCs/Tcells
7	JA-2042		carcinogen	yes	moderate	moderate	high	2 weeks	Treg cells

nd = not determined

Immuno-Oncology Platform

Overview

Our immuno-oncology platform supports the discovery of immune-modulatory drugs and the investigation of immune effects from conventional therapies. From functional in vitro assays through immunocompetent mouse models with comprehensive biomarker analysis, we connect mechanistic insights to in vivo outcomes

Immuno-Oncology Models

Syngeneic Mouse Models: 28 cell line-derived models with validated checkpoint inhibitor response data. Placement options include subcutaneous, SubQperior™, orthotopic, and metastasis models. SubQperior™ implantation into the mammary fat pad eliminates ulceration issues, extending treatment windows for reliable immune-modulating drug evaluation.

Humanized Models: Immunodeficient mice reconstituted with human immune cells for testing human-specific antibodies and cell therapies. Five characterized PBMC donors available with flow cytometry immune profiling.

CART Platform: End-to-end workflow from T cell engineering (lentivirus or transfection) through cytotoxicity assays, cytokine profiling, and in vivo validation in hematological or solid tumor models.

Immuno-Oncology Assays

T Cell Killing Assays: Co-culture tumor cells with T cells, PBMCs, or CAR-T cells to measure cytotoxicity. Primary readout via xCELLigence real-time impedance monitoring; alternative formats include high-throughput luciferase-based (384-well), flow cytometry-based, and 3D tumor spheroid penetration assays.

T Cell Functional Assays: Proliferation, activation, and cytokine determination assays reveal immune response triggers and mechanism of action, highlighting resistance mechanisms for immunotherapies.

NK Cell & Macrophage Assays: NK cell killing and ADCC assays for antibody-dependent therapies; macrophage differentiation and phagocytosis assays for tumor-associated macrophage biology and CD47/SIRPα pathway evaluation.

Immunophenotyping Services

Flow Cytometry: Quantify immune cell populations with up to 20-marker multiplexing using our Sony ID7000 spectral platform. Off-the-shelf panels include NK/T Cell, T Cell Activation Marker, DC, and All-in-One panels covering T cells, NK cells, B cells, MDSCs, DCs, macrophages, and neutrophils. Detection of 27 activation and exhaustion markers including PD-1, TIM-3, LAG-3, and functional markers.

Multiplex Immune Assays: Simultaneous measurement of 10+ cytokines and chemokines from plasma, serum, or tissue lysates via MSD platform—revealing systemic immune responses and tumor microenvironment signaling.

IHC & Tissue Microarrays: Microscopy-based detection of immune cells and checkpoint proteins directly in tumor tissue. Tissue cores from all syngeneic models available in tumor tissue microarrays (TMA) for up-front characterization of target immune cell presence and spatial distribution.

Notes

Notes

LET'S DISCOVER TOGETHER.

Recombinant Proteins

- Kinase proteins
- Epigenetic proteins
- Substrates
- Custom-tailored protein production



Biochemical Assays

- Kinases, Epigenetic Enzymes
- Protein: Protein Interaction assays
- Metabolic & Pathway
- Receptors & Channels



Cell-Based Assays

- 2D and 3D proliferation assays
- Drug combination screening
- Migration assays
- Angiogenesis assay



Biophysical Assays

- Surface Plasmon Resonance
- Thermal Shift Assay
- Isothermal Titration Calorimetry
- Microscale Thermophoresis



In Vivo Pharmacology

- In Vivo Hollow Fiber Model
- Xenograft models
- Orthotopic models
- Metastasis models



Safety & Toxicology

- In Vitro Safety Panel
- Cardiac Safety Panel
- Cytochrome P450
- Maximum-Tolerated Dose
- GLP Regulatory Toxicology



Integrated Solutions

- Antibody Drug Conjugates
- RAS Drug Discovery
- Targeted Protein Degradation



Biomarker Discovery

- Genomic biomarkers
- Protein biomarkers
- Immunophenotyping



Immuno-Oncology

- In Vitro Killing Assays
- Syngeneic Mouse Models
- Proprietary Tumor Models
- Immunophenotyping

