

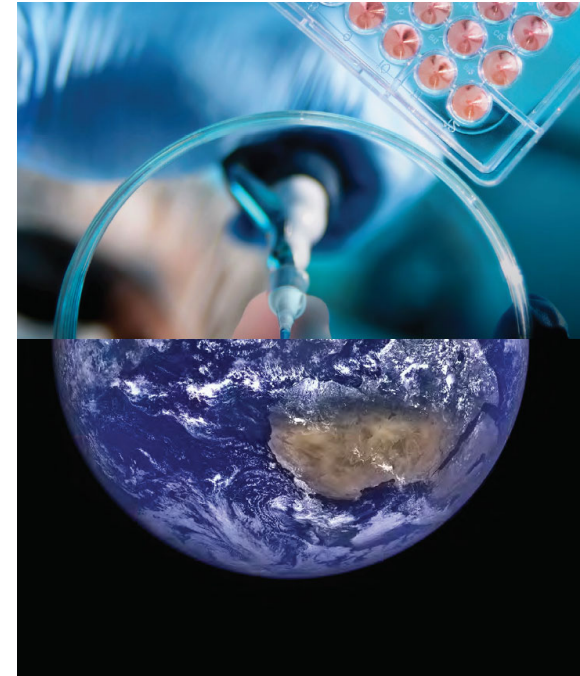
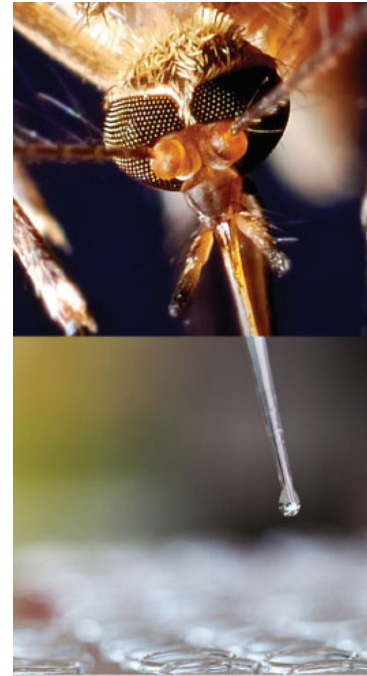
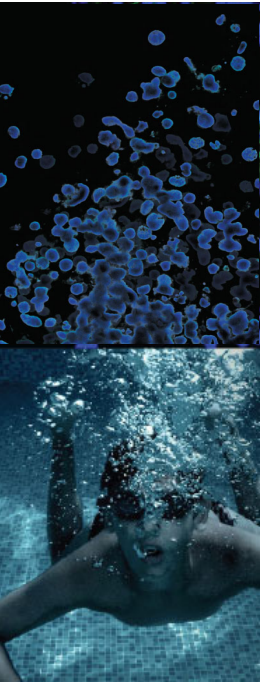


Immortalized Primary Cells: The Best of Both Worlds

Jonathan Sagal, BS
Product Manager, ATCC

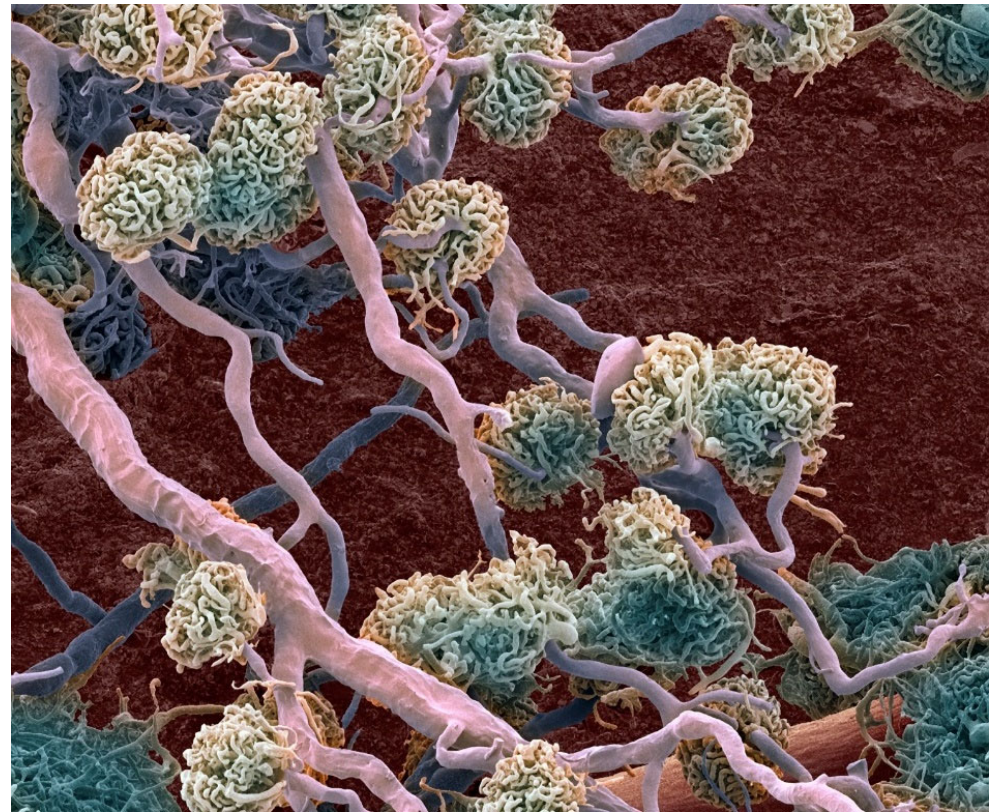
Sujoy Lahiri, PhD
Lead Scientist, ATCC

Credible Leads to Incredible™



Agenda

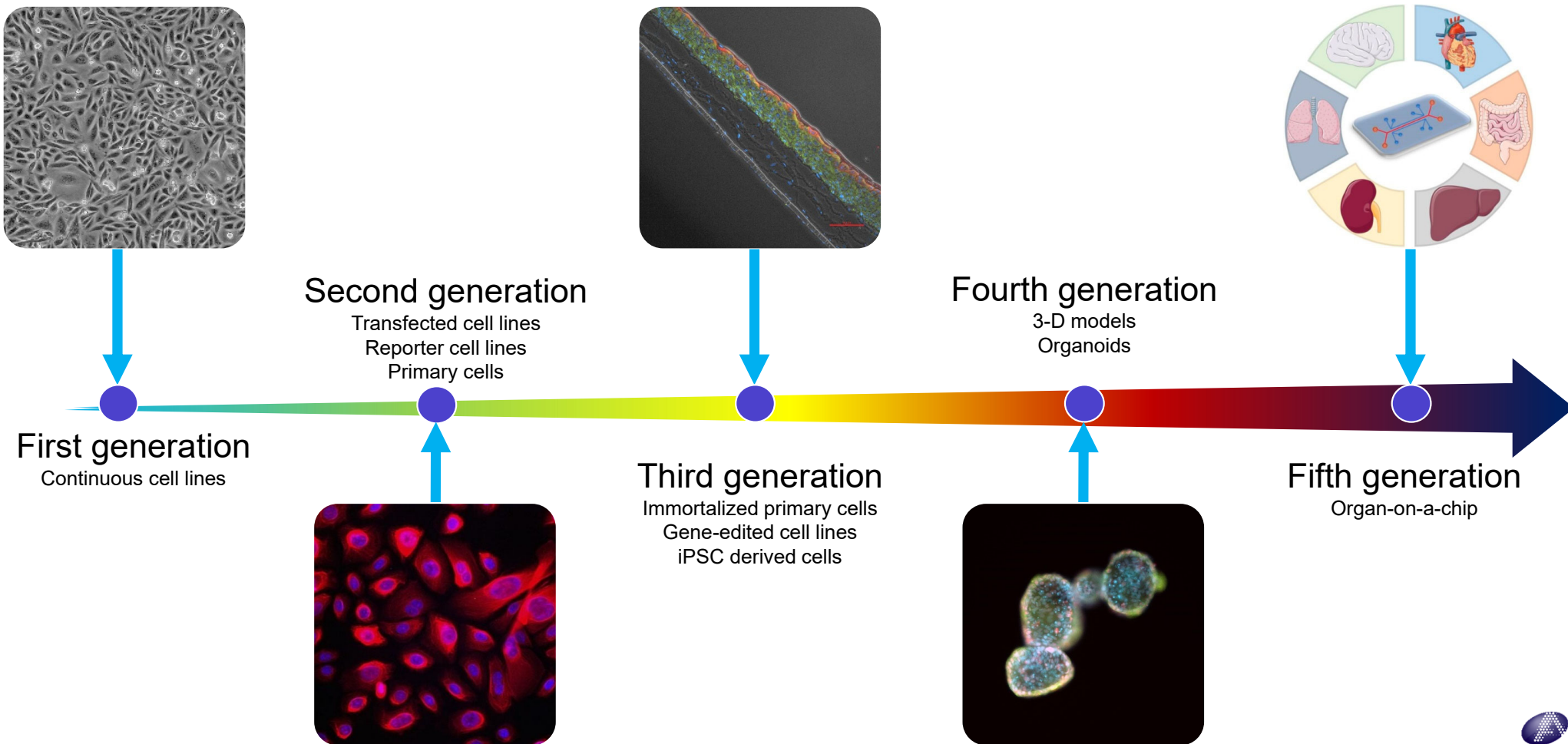
- ATCC® mission and future direction
- ATCC® immortalized primary cell portfolio
- Case studies
 - Kidney models
 - Cardiovascular models



About ATCC®

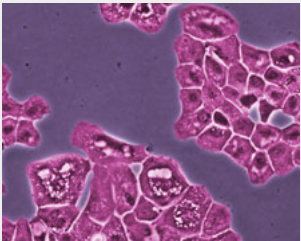
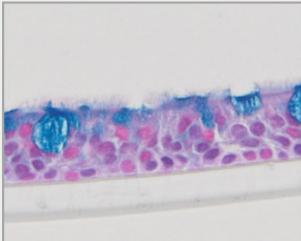
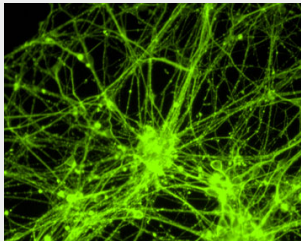
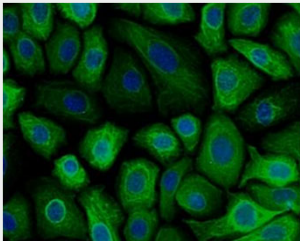
- Founded in 1925, ATCC is a non-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's largest, most diverse biological materials and information resource for cell culture – the “*gold standard*”
- Innovative R&D company featuring gene editing, differentiated stem cells, advanced models
- cGMP biorepository
- Partner with government, industry, and academia
- Global supplier of authenticated cell lines and viral and microbial standards
- Sales and distribution in 150 countries, 19 international distributors
- Talented team of 450+ employees, over one-third with advanced degrees

Evolution of in vitro cell models



Sources of cell models

Cells are one of the most important resources for predicative in vitro models

Cell lines	Primary cells	iPS-derived cells	Immortalized primary cells
<ul style="list-style-type: none"> Highly proliferative Easy to culture and transfect Differ genetically and phenotypically from their tissue origin Accepted for specific functions, i.e., Caco-2 is well accepted for drug transporters studies; however, it lacks physiological CYP activities 	<ul style="list-style-type: none"> Most accepted cell types; physiological relevance Maintain many of the important markers and functions Higher predictability Donor-donor variability Finite lifespan and limited expansion capacity Difficult to source or preserve some tissue types (alveolus, cardiomyocytes, neurons) 	<ul style="list-style-type: none"> Advantages of precision medicine It might maintain many of the important markers and functions Multiple cell types from the same donor Need experienced users to culture and maintain Cryopreservation is an issue Poor differentiation/ maturation 	<ul style="list-style-type: none"> Maintain physiological feature of primary cells Allows for extended cultivation Improved supply and reproducibility Amenable to genetic editing Need experienced users to generate and characterize
 <p>Chordoma Cell Line</p>	 <p>Bronchial Epithelial Cells</p>	 <p>Neural Progenitor Cells</p>	 <p>hTERT Bronchial Epithelial Cells</p>

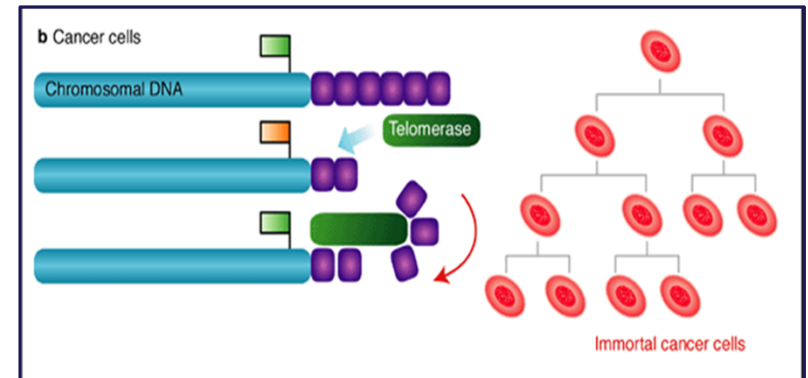
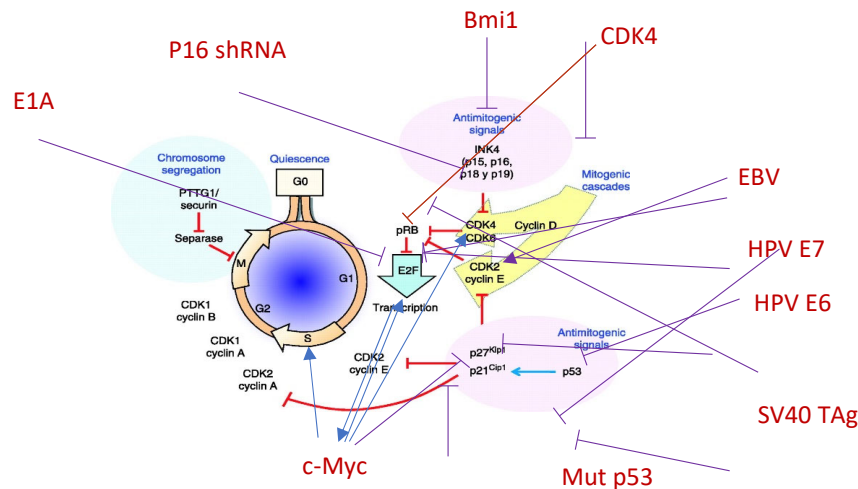
Immortalization - Cell cycle regulation

- Primary cells can only replicate a handful of times, or not at all. Cellular senescence, donor availability, and donor-to-donor variability limit the applications

Immortalizing genes:

- hTERT**
- Viral oncogenes: SV40T Antigen, EBV, HPV E6/E7, E1A
- Other cellular genes: CDK4, **Bmi1**, P16, c-Myc, mutant p53

Cell cycle regulation by immortalizing genes:

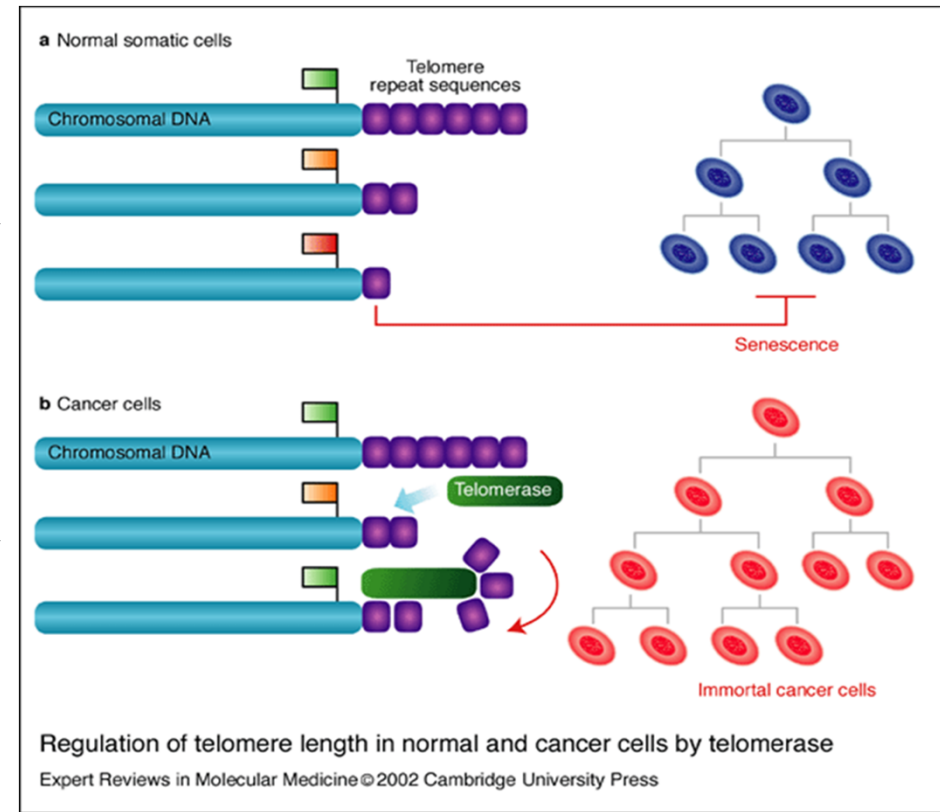
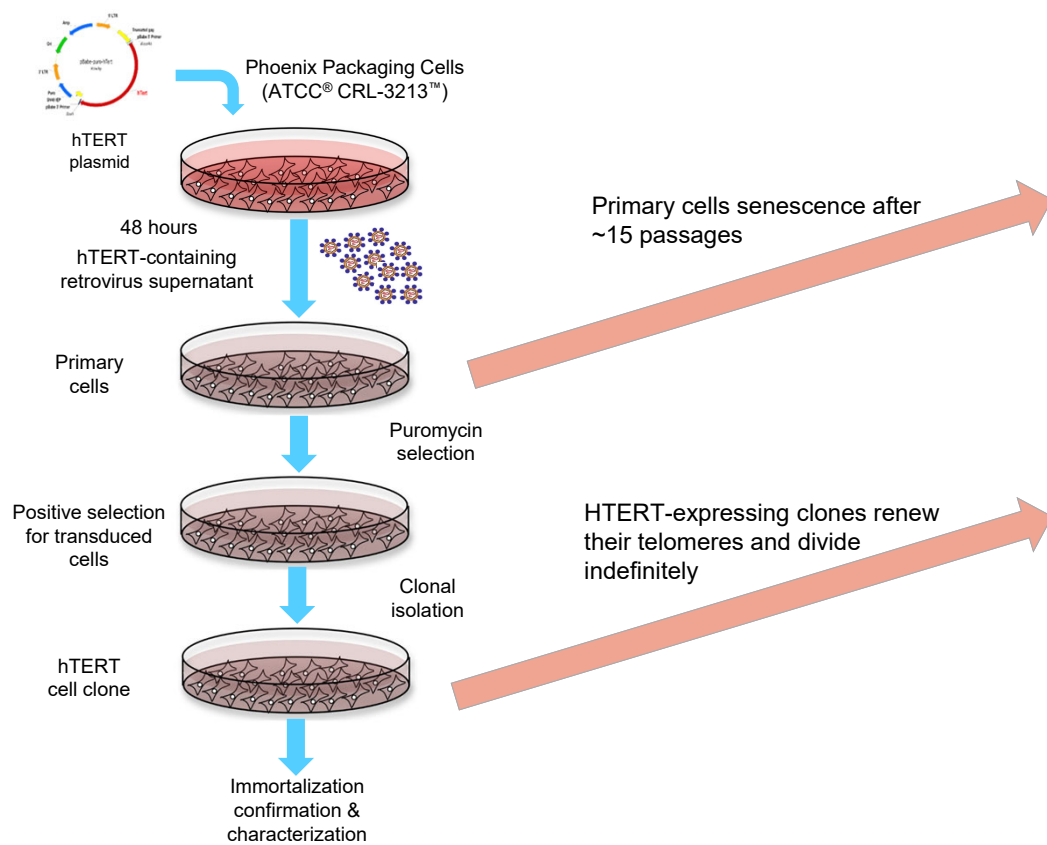


Modified from Victor Q, Malumbres M. J Mol Endocrinol 42(2):75-86, 2009.

Keith W, et al. Expert Rev Mol Med 22;4(10):1-25, 2002.

Cell immortalization process - hTERT alone

Retroviral Transduction of telomerase (hTERT) gene and Clone Selection



Keith W, et al. Expert Rev Mol Med 22;4(10):1-25, 2002.



Immortalized primary cells – Key characteristics

ATCC® offers primary and immortalized cell solutions that are authenticated via rigorous QC and tested for common biomarker expression and cell performance

■ Growth

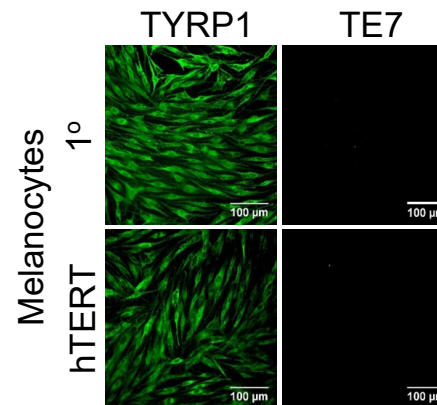
- Cells retain replicative capacity (“immortalized”)
- Population doubling rate is comparable to primary cells

■ Characterization

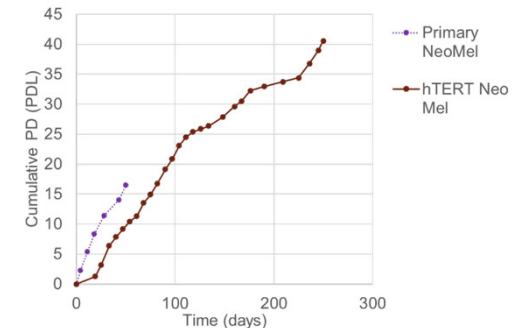
- Morphology and marker expression similar to primary cells
- Sterility testing
- STR profile

■ Functional responses

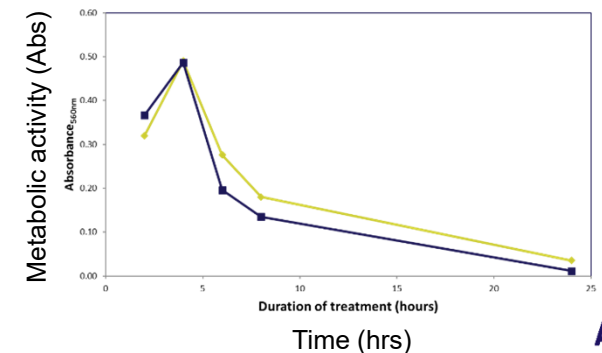
- Within expected range
- Analogous to primary cells

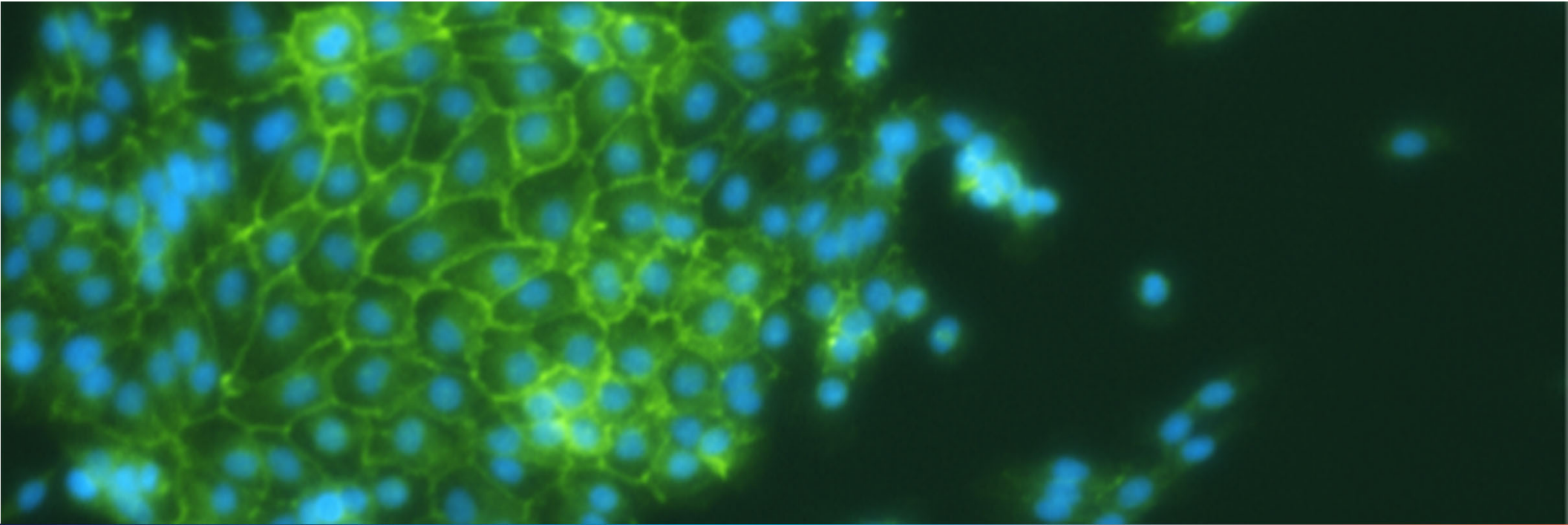


Growth of ATCC® CRL-4064™ Neonatal Dermal Melanocytes



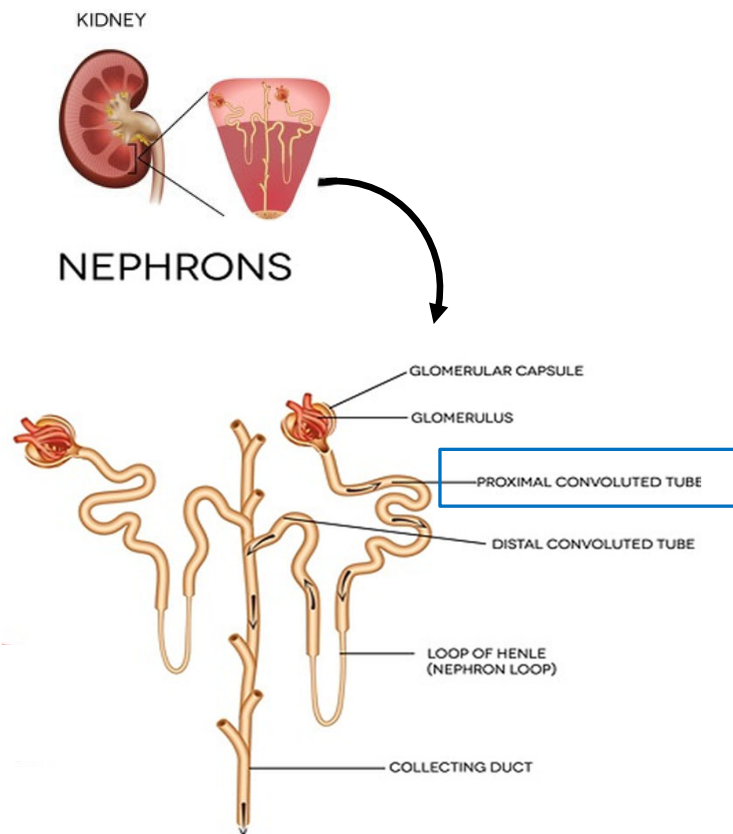
Metabolic reduction by 3-D organotypic skin culture in Triton-X





Kidney Models and Functionality

Kidney drug toxicity and transporters



Nephron, the functional unit of the kidney. Credit: Modified image by TefiM

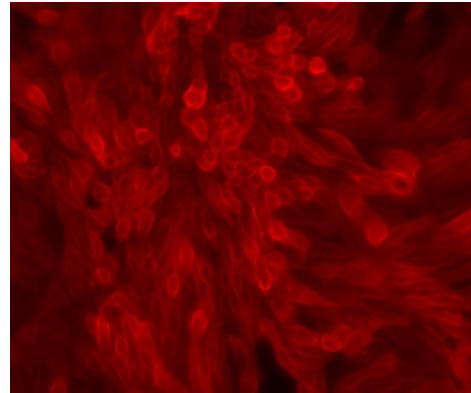
- **The kidney is one of the major target organs for drug-induced toxicity**
 - Large functional reserve of the kidney
 - Nephrotoxic effects become obvious only after regulatory approval
- **Nephrotoxic potential**
 - Often underestimated when new drugs are available
 - Leads to clinical complications such as COX2 inhibitors
- **Renal proximal tubule (PT, blue box) is a major target for drug-induced toxicity due to its role in:**
 - Glomerular filtrate concentration
 - Transport of drugs and organic compounds

Kidney models

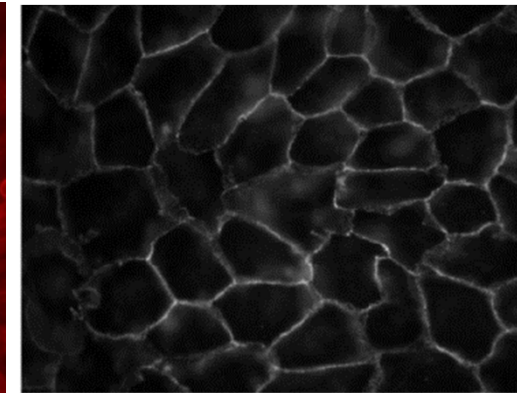
Renal proximal tubule epithelial cells

- hTERT-RPTEC–immortalized renal proximal tubule epithelial cells
- Key characteristics:
 - Uniform expression of E-cadherin and CD13 (aminopeptidase N)
 - Formation of dome-like structures
 - Stabilized transepithelial electrical resistance (TEER)

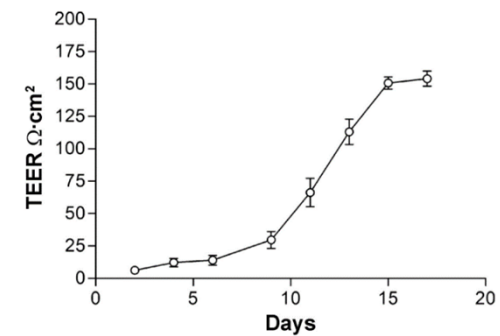
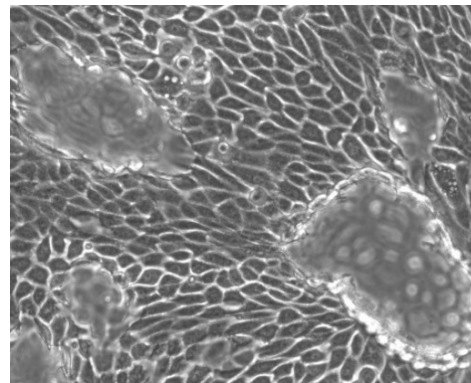
RPTEC/TERT1: CD13



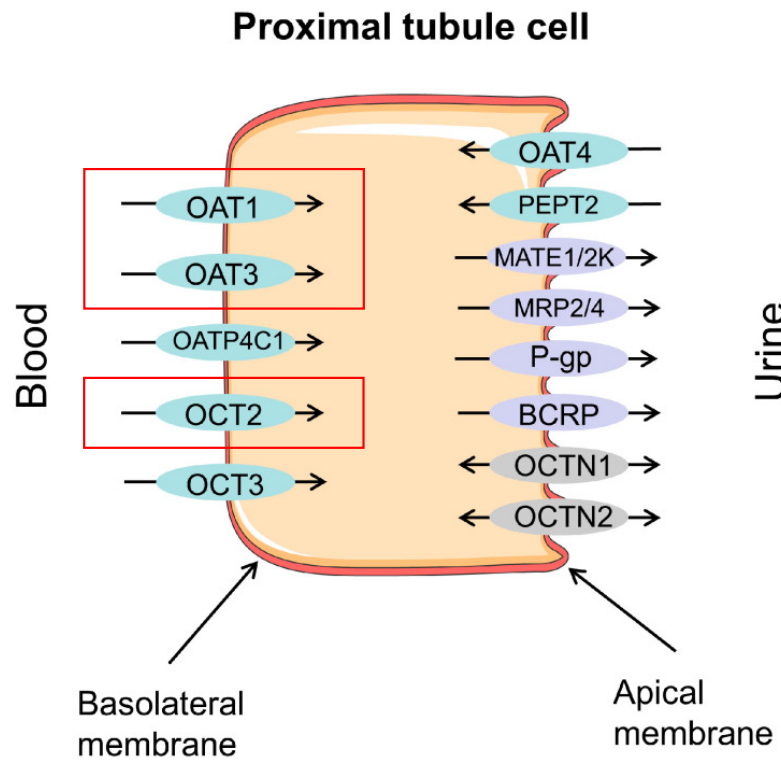
RPTEC/TERT1: E-cadherin



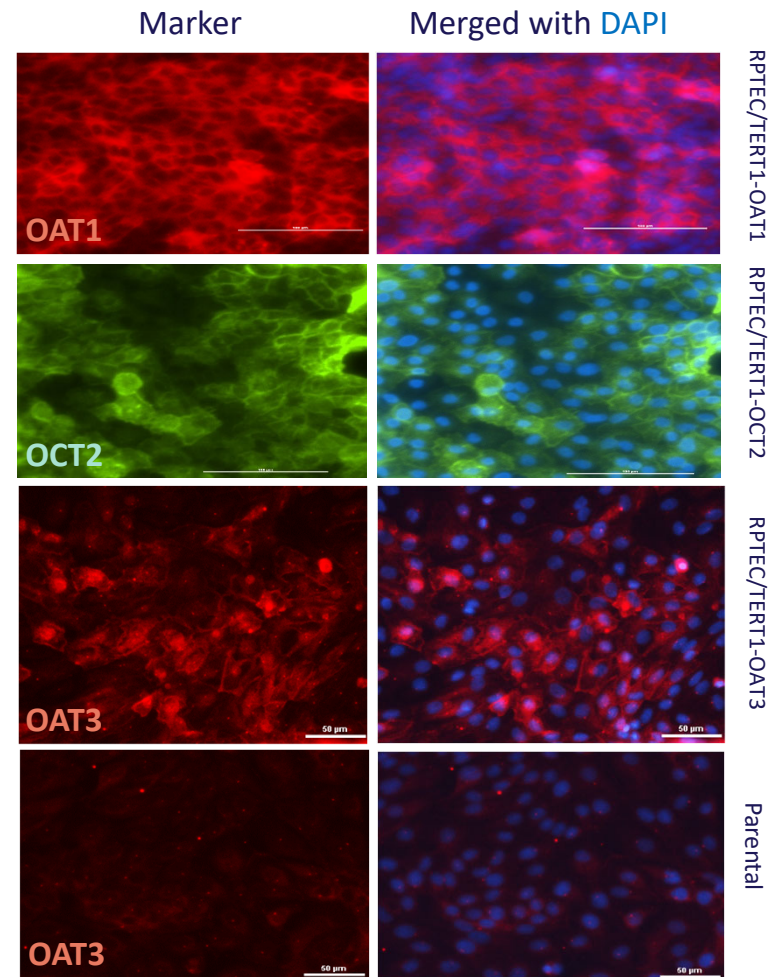
Dome formation



Kidney cells – Modeling solute transporters

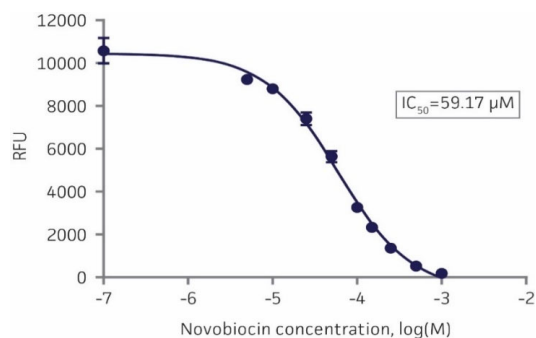
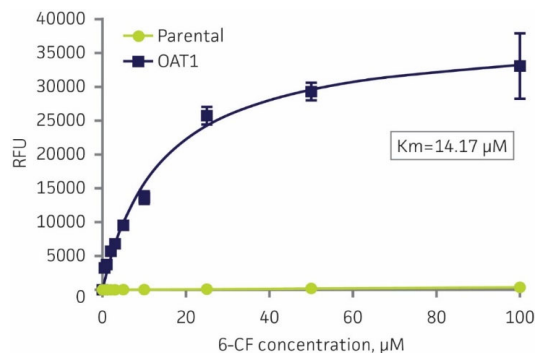


Lin K, et al. *Molecules* 28(13): 5252, 2023.

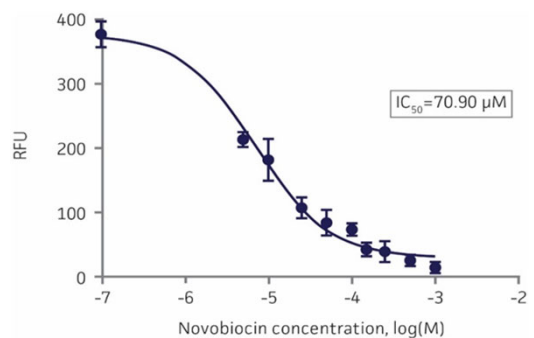
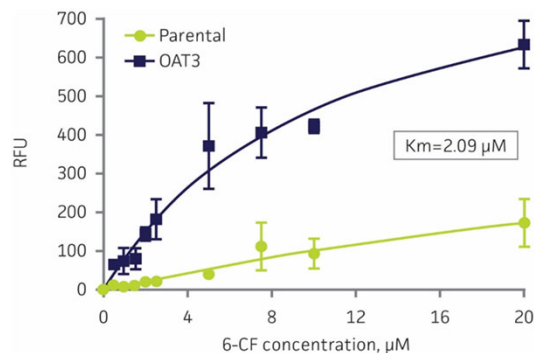


Functionality – Drug uptake & inhibition assay

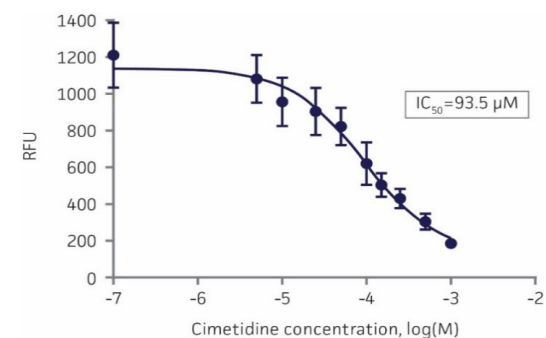
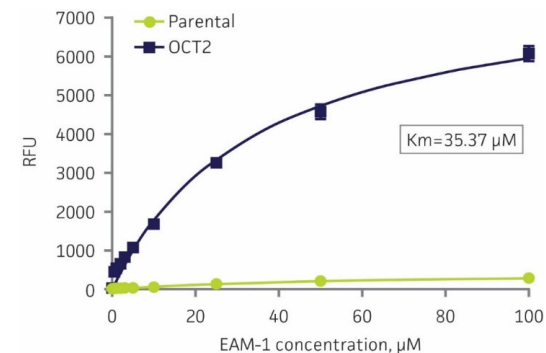
6-CF uptake inhibition in HTERT/RPTEC OAT-1



6-CF uptake inhibition in HTERT/RPTEC OAT-3



EAM-1 uptake inhibition in HTERT/RPTEC OCT-2



ASSAY PROTOCOL

- Equal numbers of both parental and transporter cells were seeded into 96-well plates in triplicate for 24 hours
- Increasing concentrations of 6-CF or EAM1 with or without inhibitors were added, and cells were incubated for 20 minutes at 37°C
- After washing with cold HBSS 4 times, cells were lysed and uptake intensity was measured

RPTEC-OCT2 – Drug-drug interactions (DDI) application

The problem

- A multitude of disease- and therapy-related factors drive the frequent development of renal disorders in cancer patients
- Many cancer patients have comorbidities such as urinary tract infections, tuberculosis, and diabetes
- Commonly prescribed medications such as levofloxacin (TEA) or metformin can interact with chemotherapeutics
- These common medications can block the renal uptake of the candidate via organic cation SLC transporters



The solution: Use SLC transporter cells to identify DDI

- Incubate candidate drugs with radiolabeled known SLC substrate drugs in RPTEC-OCT2 cultures then monitor uptake
- If uptake of radiolabeled compounds is inhibited, then DDI is indicated



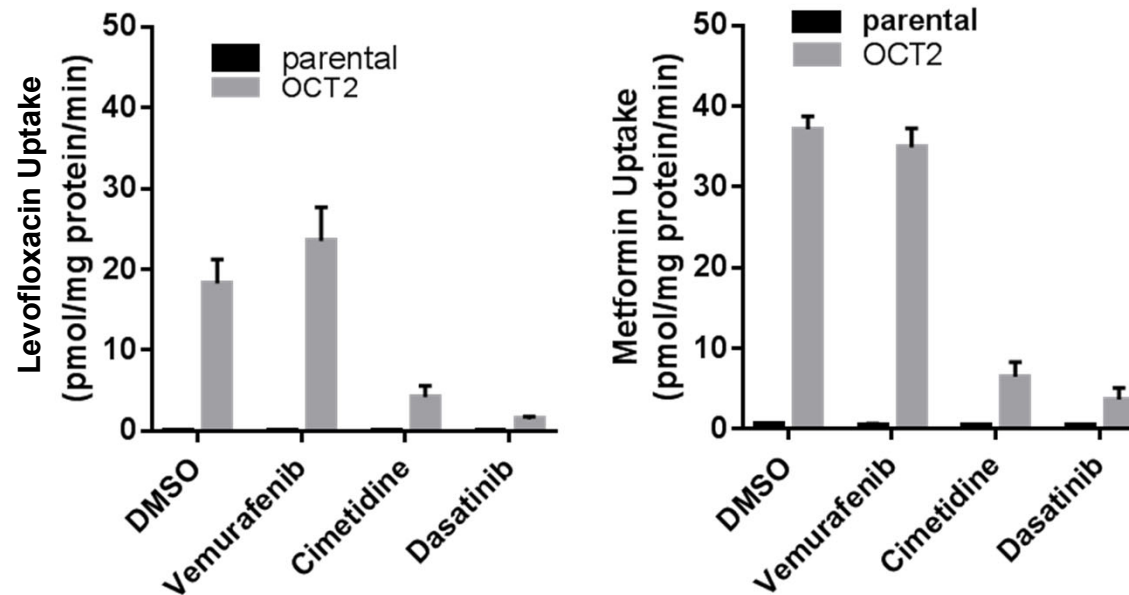
Uptake inhibition assay protocol

- Aspirate growth media and wash once with warm 1X PBS; remove PBS and add 250 μ L of cold inhibitors (prepared serum-free DMEM, 0.5 μ M) and incubate for 15 minutes
- Remove inhibitors and add 250 μ L of radio-labeled TEA or metformin (prepared serum-free DMEM, 4.5 μ M) and incubate for 15 minutes
- Remove drug and wash 3 times with cold PBS; lyse the cells and count



RPTEC-OCT2 – Drug-drug interactions (DDI) application

Drug-drug interactions



Data kindly provided by:

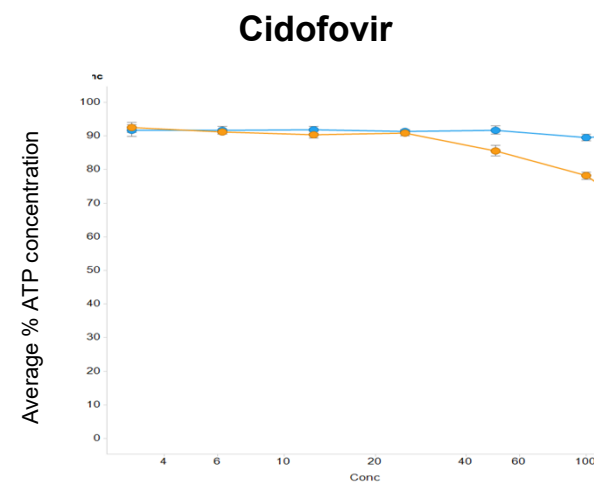
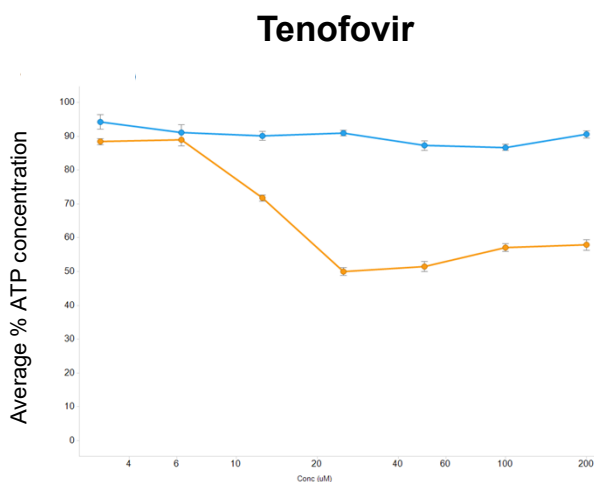
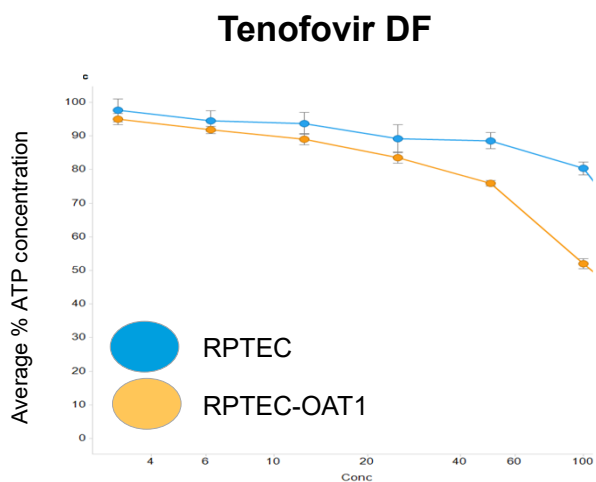
Kevin Huang, *Graduate Research Associate*, Ohio State University, College of Pharmacy
Alice Gibson, Ph.D., *Senior Research Specialist*, Ohio State University, College of Pharmacy

Nephron toxicity testing

Data kindly provided by: Merck & Co., Inc.

Use SLC transporter cells to identify nephron toxicity

- Targeted therapeutics can cause renal dysfunction through on- and off-target mechanisms
- Incubate candidate molecules with RPTEC-OAT1 or parental RPTEC cells, then monitor toxicity to identify mechanism of action



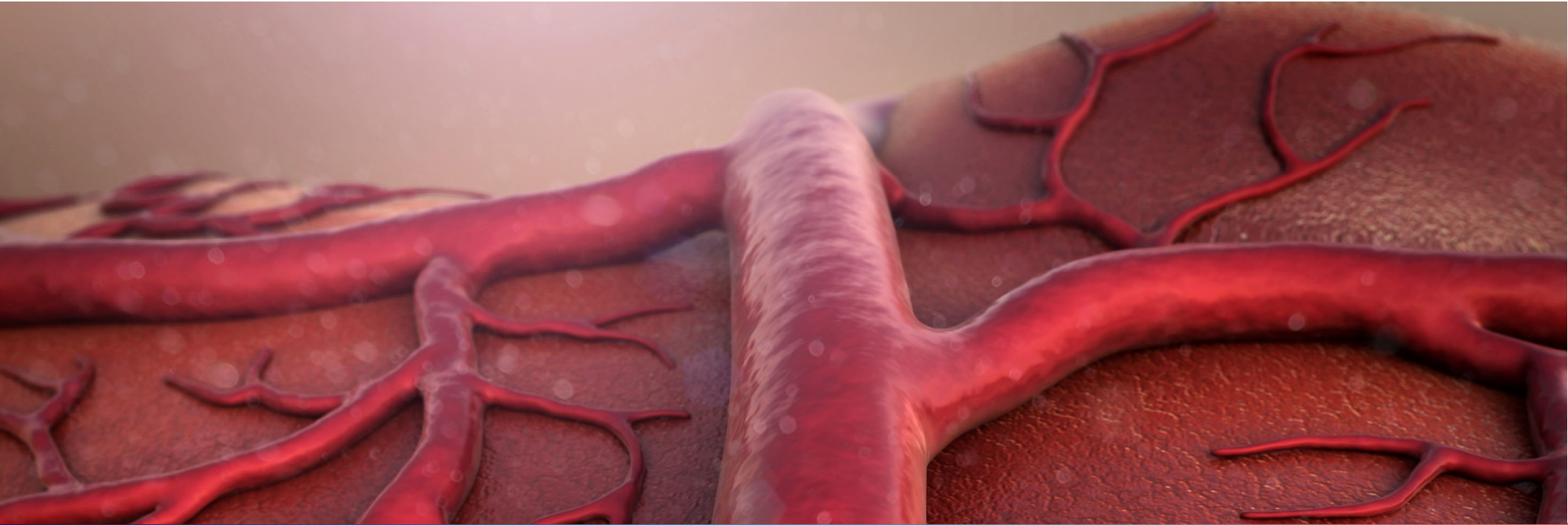
ASSAY PROTOCOL

- About 35000 cells were seeded per well in triplicate in a 96-well plate and incubated overnight
- Cells were incubated with a series of compounds at various concentrations for 3 days
- Cell viability was determined using a cell viability assay per the manufacturer's instructions

Summary of kidney models

- ATCC® primary and hTERT-immortalized RPTECs display many key in vivo characteristics
- We enhanced hTERT-immortalized RPTEC with organic anion/cation transporter proteins
- The hTERT-immortalized RPTEC models have been evaluated for
 - Uptake of specific fluorescent substrates
 - Selective substrate drug uptake effects by known transporter protein inhibitors
 - Drug-drug interactions (DDI) and nephron toxicity applications



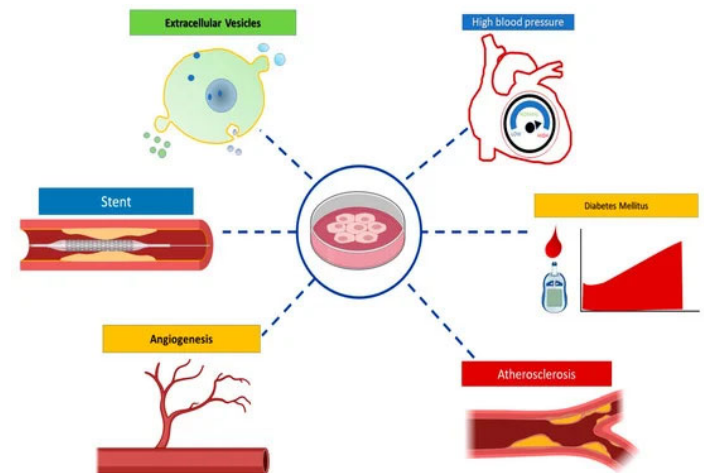
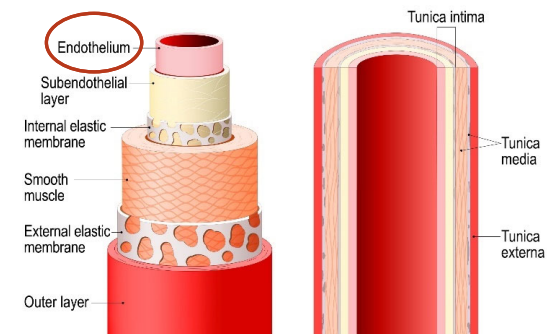


Cardiovascular Models and Functionality

Cardiovascular research

- Cardiovascular disease causes one-third of deaths worldwide and represents an urgent threat to global health
- Efforts to treat and cure cardiovascular disease depend upon advances in our understanding of the etiology and molecular mechanisms affecting the disease
- Areas of active research include:
 - Basic cardiovascular development and angiogenesis
 - Metabolism
 - Regenerative medicine, including tissue bioengineering
 - Cardiovascular and blood diseases
 - Cardiovascular effects of inflammation
 - Vascular and lymphatic disorders
 - Hypertension & atherosclerosis
- As a key cell in vascular tissue, endothelium cells are central to this research activity

Anatomy of arteries



Medina-Leyte DJ, et al. Appl Sci 10(3), 938, 2020.

Cardiovascular research - Angiogenesis

- Angiogenesis, or blood vessel formation, is the sprouting of new capillaries from pre-existing vessels
- Plays a central role in numerous human diseases, including cancer and age-related macular degeneration
- Numerous anti-angiogenesis drugs have been developed, and several have received FDA approval

In vitro angiogenesis models can be used for

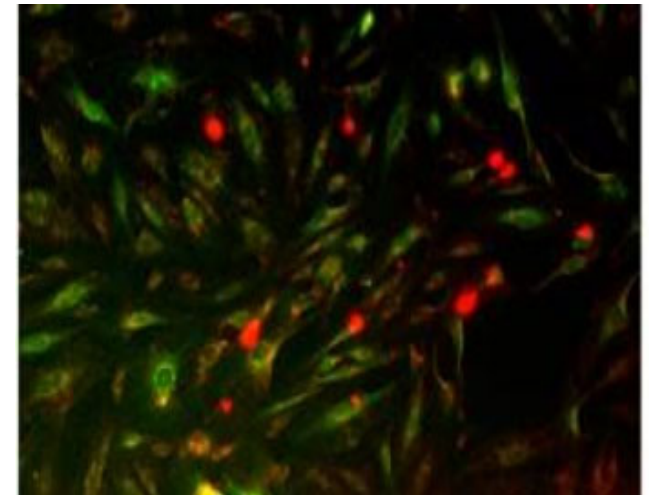
- Screening of activation and inhibition compounds
- Vessel formation in wound healing study and regenerative compound screening
- Angiogenesis under hypoxia situation study and compounds screening



Immortalized primary endothelial cells

Broad and growing collection of cell models

- HUVEC/TERT2 (ATCC® CRL-4053™)
- TIME (Microvascular endothelial cells; ATCC® CRL-4025™)
- TIME-GFP (GFP-expressing microvascular endothelial cells; ATCC® CRL-4045™)
- NFκB-TIME (ATCC® CRL-4049™)
- TeloHAEC (ATCC® CRL-4052™)
- TeloHAEC-GFP (Aortic endothelial cells; ATCC® CRL-4054™)
- HPAEC-BMI1 (ATCC® CRL-4065™) **New!**



TIME-GFP stained for Ac-LDL (red) and CD31

Pulmonary hypertension and hPAEC

- Pulmonary arterial hypertension (PAH) is a medical condition with a 1% global incidence rate
 - 100,000 patients in US/year
- High blood pressure in the lungs occurs due to narrowed and stiffened pulmonary arteries.
- Human Pulmonary Artery Endothelial Cells (hPAEC) play a key role in the etiology of the disease

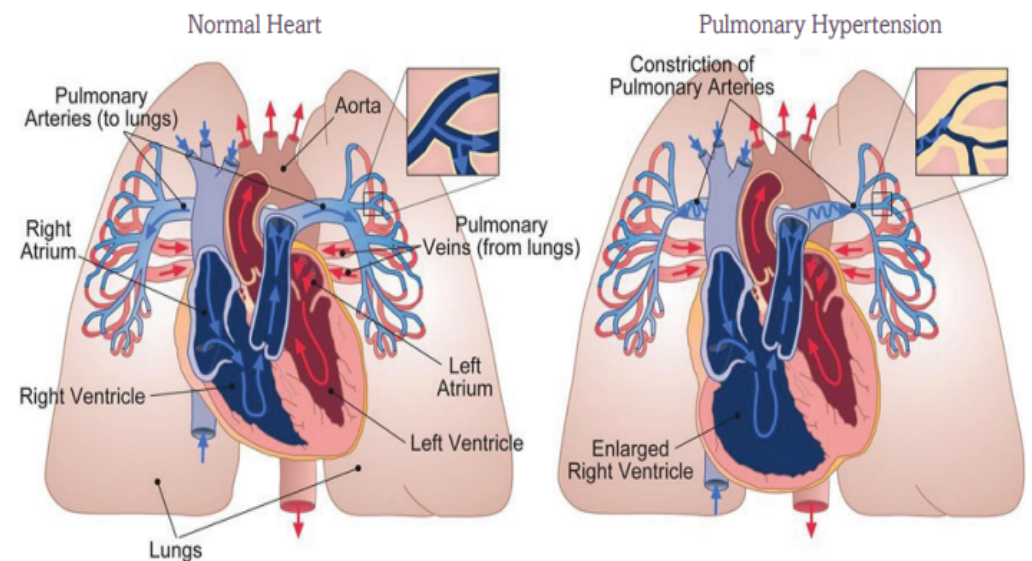


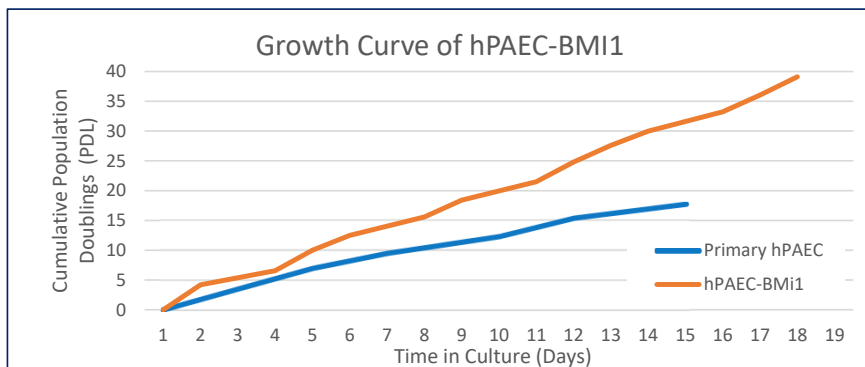
Image courtesy of the Columbus Ohio Adult Congenital Heart Disease Program at Nationwide Children's Hospital Heart Center, Columbus, Ohio

HPAEC-BMI1– Key characteristics

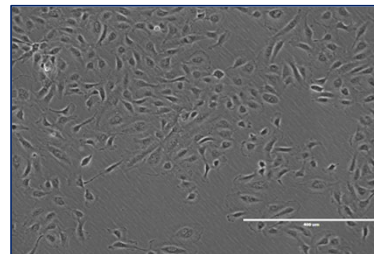
ATCC® offers primary and immortalized cell solutions that are authenticated with our rigorous QC and tested for common biomarker expression and cell performance

Growth & Morphology

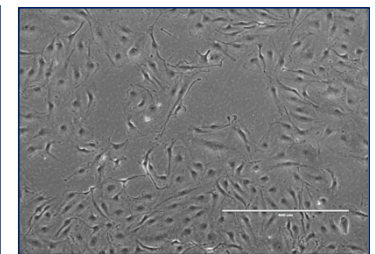
- Cells retain replicative capacity (“immortalized”)
 - Morphology similar to primary cells
- Key biological functions similar to primary cells
 - Formation of capillary-like tubes
 - 82% of cell uptake acLDL



Primary hPAEC

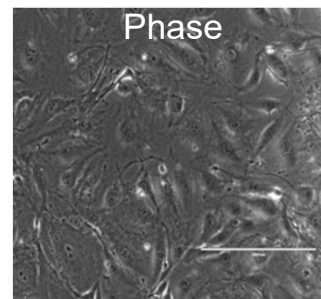


hPAEC BMI1

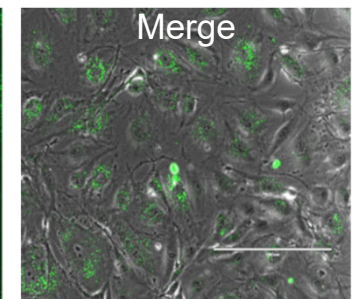
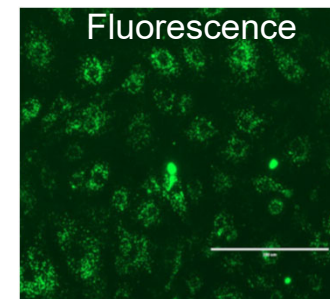


10x magnification

acLDL Uptake



20x magnification



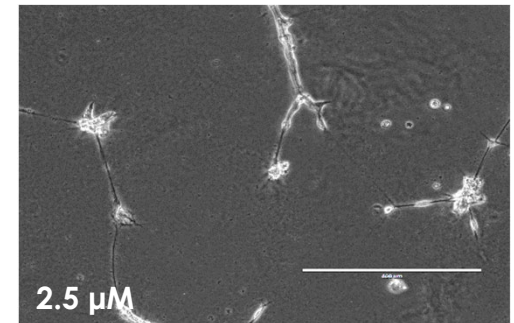
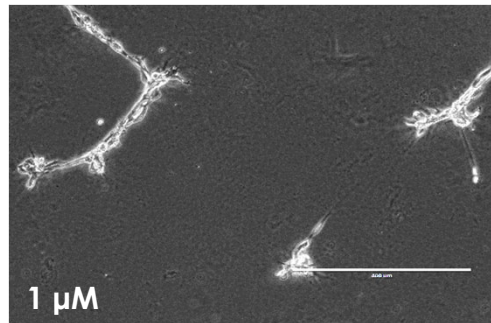
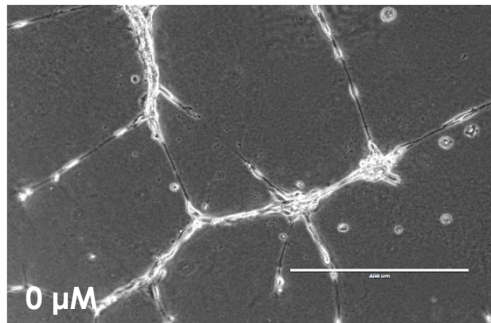
HPAEC-BMI1– Key characteristics

ATCC® offers primary and immortalized cell solutions that are authenticated with our rigorous QC and tested for common biomarker expression and cell performance

Application Data

- Capillary-like tube formation assay is responsive to well know stimulatory and inhibitory agents

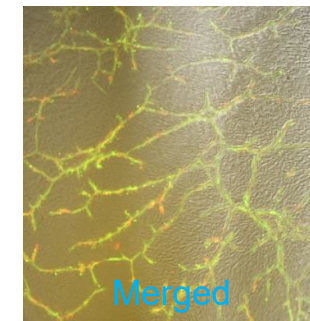
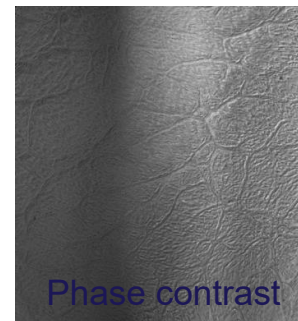
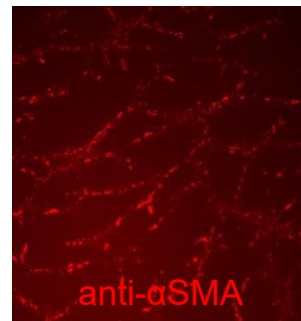
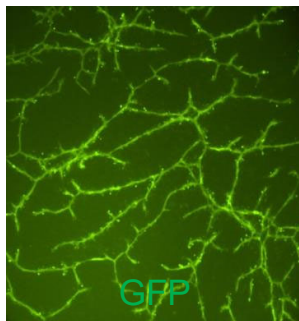
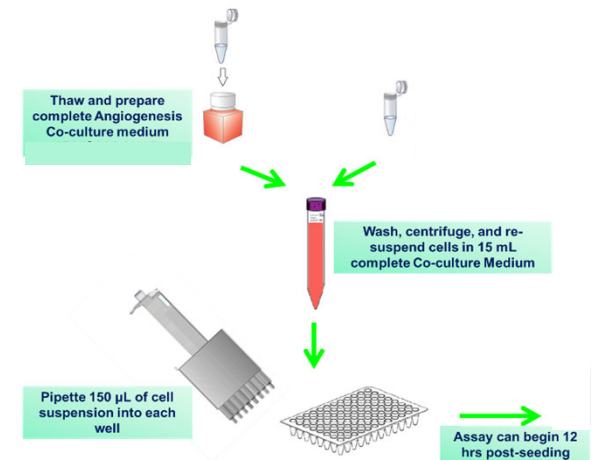
Sunitinib Inhibition



10x magnification

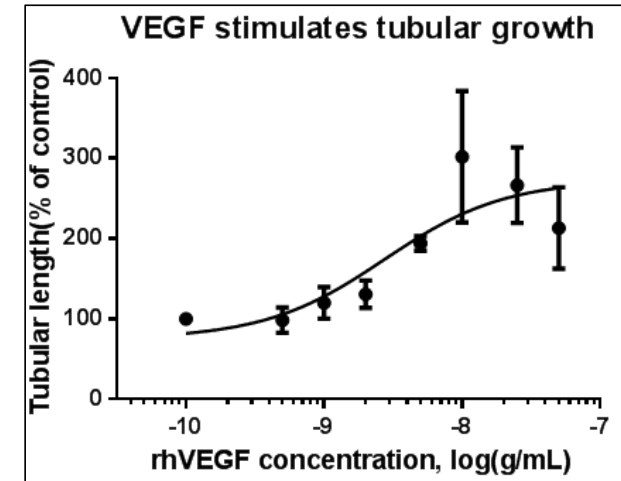
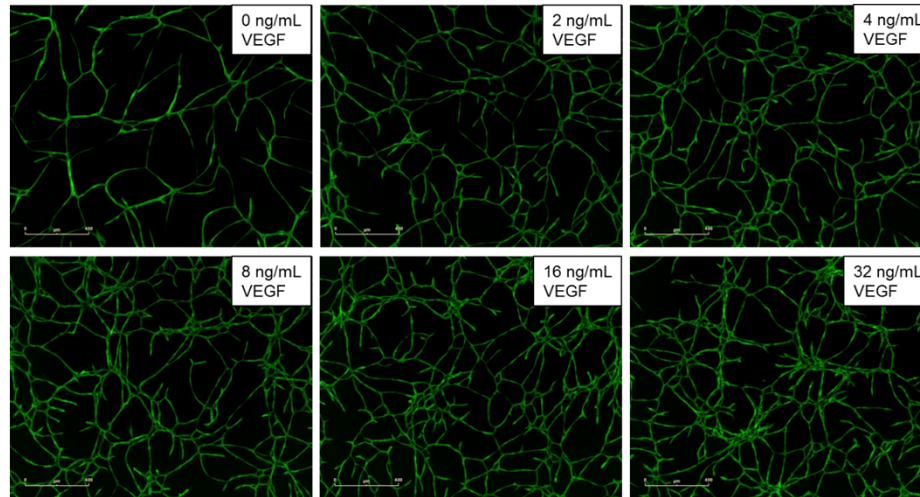
An advanced co-culture system

- Co-culture of endothelial cells with fibroblasts allows for
 - Formation of more heterogeneous tubules
 - More defined culture conditions, no need for ECM coating
- Scientists at ATCC® optimized a co-culture system using:
 - hTERT-immortalized mesenchymal stem cells ASC52telo
 - GFP-labeled TeloHAEC, an immortalized aortic endothelial cell
- Can form tubular structures in < 7 days instead of 14 days compared with standard co-culture methods
- Immunofluorescence shows MSCs surrounding the microvascular structures differentiate into smooth muscle

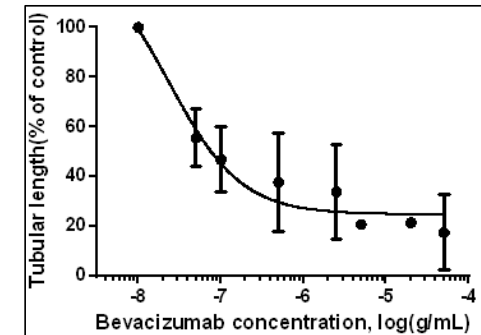
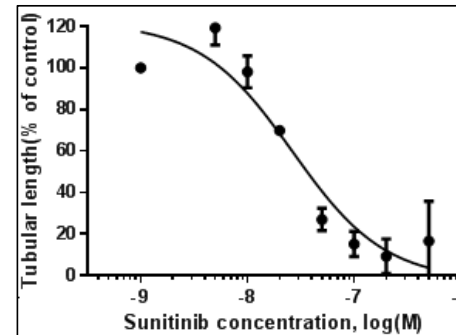
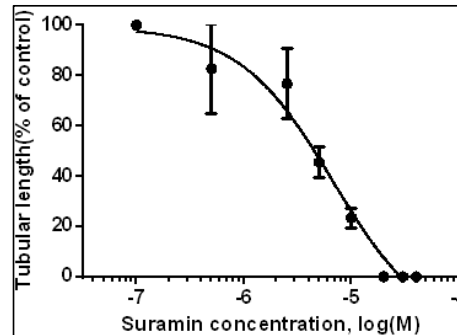


An advanced co-culture system – Response to stimuli

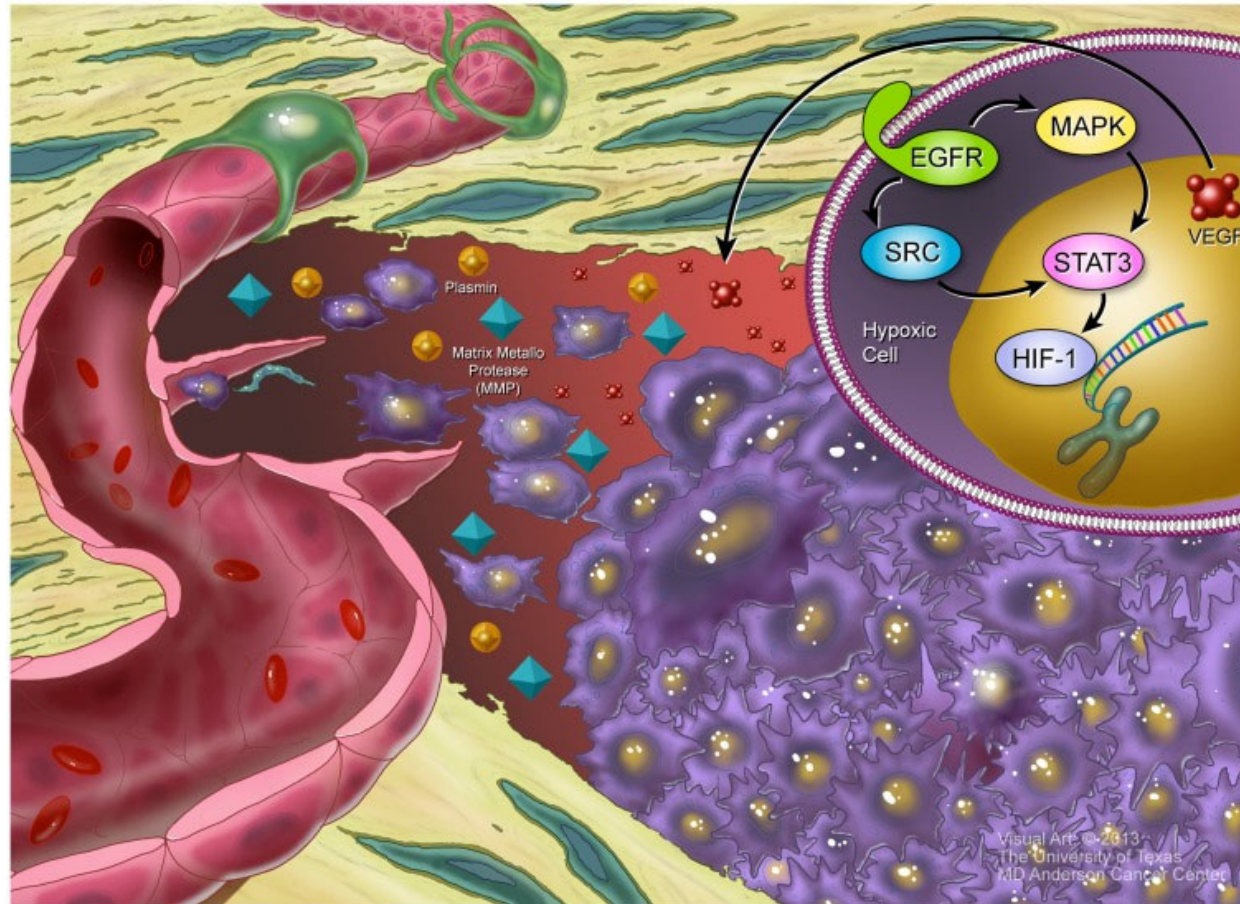
Positive response to VEGF stimulation



Response to angiogenesis inhibitors



Screening tumor angiogenesis inhibitors



Zhu L, et al. *Front Oncol* 3: 230, 2013.

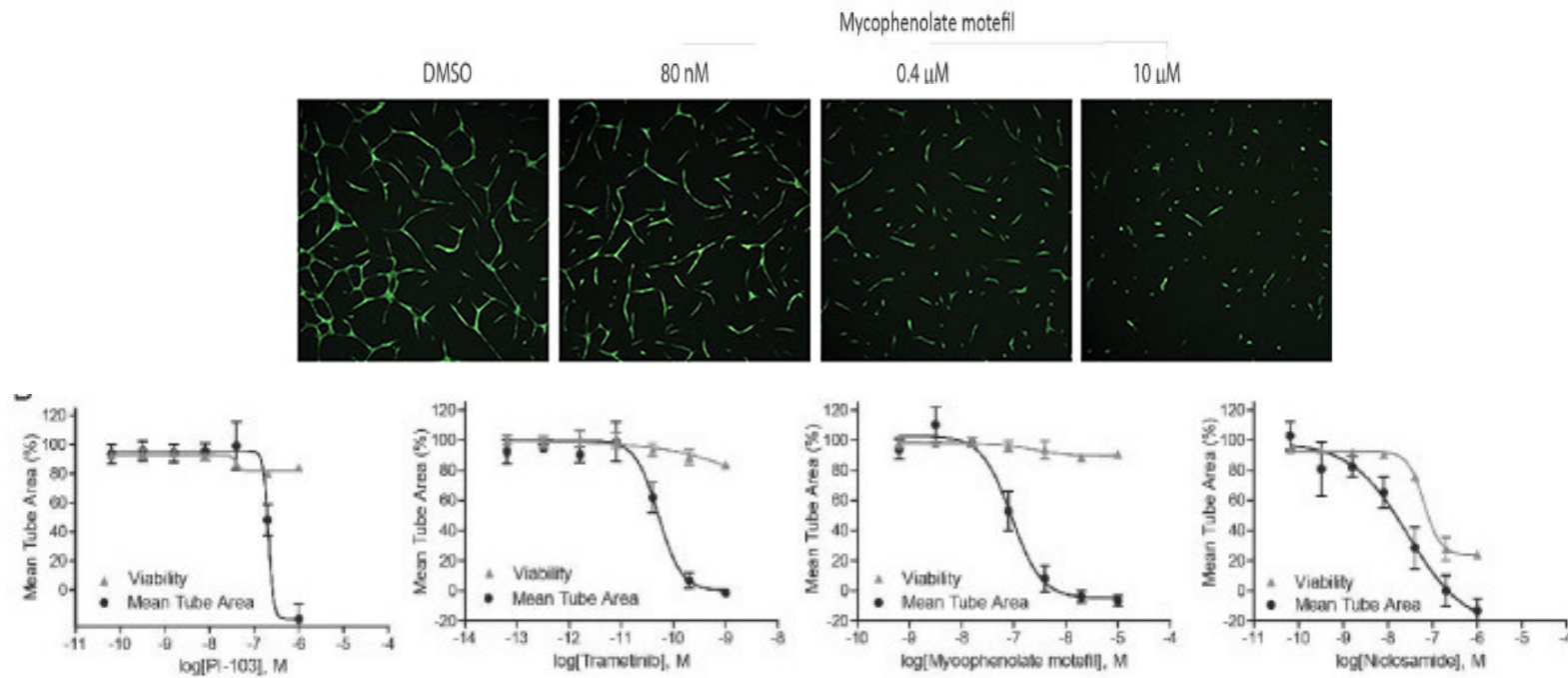
Modelling Hypoxia and tumor angiogenesis

www.impactjournals.com/oncotarget/

Oncotarget, Vol. 7, No. 7

Identification of approved and investigational drugs that inhibit hypoxia-inducible factor-1 signaling

Chia-Wen Hsu, Ruili Huang, Thai Khuc, David Shou, Joshua Bullock, Suzanne Grooby, Sue Griffin, Chaozhong Zou, Annette Little, Holly Astley, Menghang Xia



Advanced co-culture system– High-throughput screening

Identification of Angiogenesis Inhibitors Using a Co-culture Cell Model in a High-Content and High-Throughput Screening Platform

Shuaizhang Li , Chia-Wen Hsu , Srilatha Sakamuru, Chaozhong Zou, Ruili Huang, and Menghang Xia

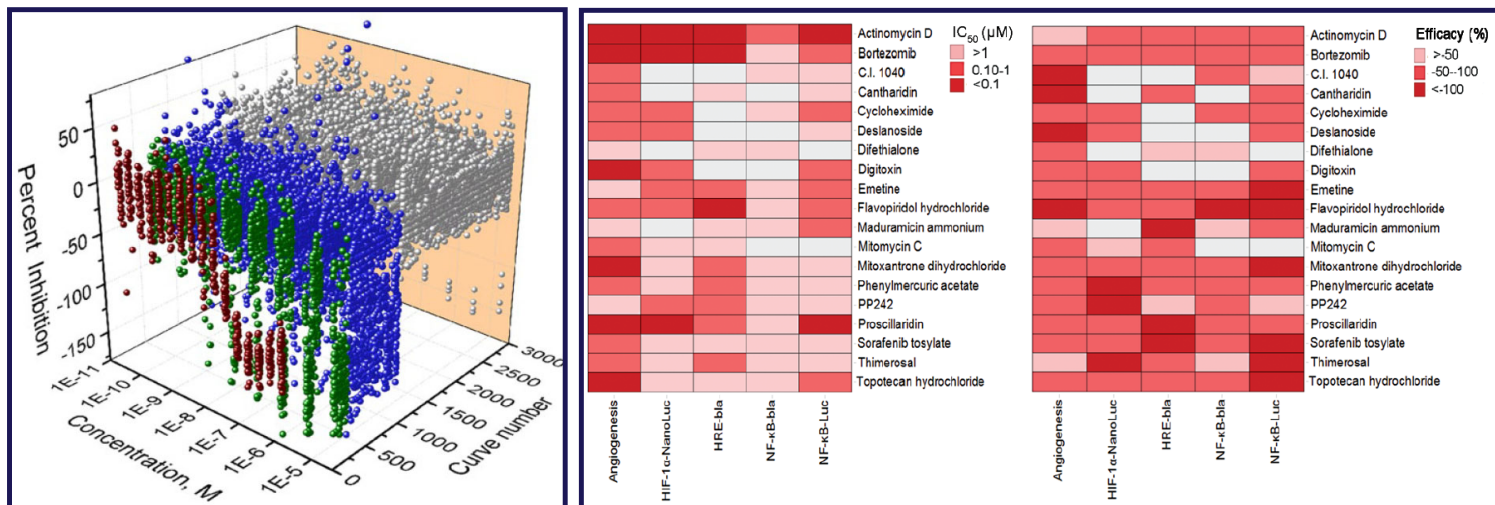
SLAS Technology
2018, Vol. 23(3) 217–225
© 2017 Society for Laboratory
Automation and Screening



DOI: 10.1177/2472630317729792
journals.sagepub.com/home/jla



Screening of 2,816 drugs on 1,536-well format using ATCC's angiogenesis co-culture model.



Summary of cardiovascular models

- ATCC® immortalized endothelial cells display vivo characteristics of primary cells
- Immortalized primary cell alone or in combination with other cells are a user-friendly solution for building reliable and predictive cell models
- In vitro co-culture models provide physiologically relevant tool to examine angiogenesis
 - The tubules formed in co-culture assays more closely resemble capillaries than those formed on a gel matrix
 - Screening of activation and inhibition compounds
 - Vessel formation in wound healing study and regenerative compound screening
 - Angiogenesis under hypoxia situation study and compounds screening



Summary and resources

- ATCC® offers a variety of immortalized cell models
 - Cell lines are authenticated for immortalization and karyotype stability
 - Including several difficult to immortalize cell types
 - Time and costs are saved in trying to immortalize these difficult cell types on one's own*
- Immortalized primary cells offer:
 - Primary cell functionality
 - Continuous cell line longevity
 - Improved supply and reproducibility for a standardized model
 - A suitable cell model for stable genetic modifications
 - Immortalization eliminates donor variability*
- Immortalized primary cells alone or in combination with other cells are a user-friendly solution for nephrotoxicity or cardiovascular studies
- Multiple resources are available at www.atcc.org/hTERT

The collage displays several ATCC research articles and guides. Key titles include:

- ATCC HUMAN BRONCHIAL/TRACHEAL EPITHELIAL CELLS: IMPROVING FUNCTIONAL STUDIES** (Chang, Cheng, and Shaper, 2010)
- ATCC® hTERT IMMORTALIZED CELL CULTURE GUIDE** (tips and techniques for culturing hTERT immortalized cells)
- IN VITRO ANGIOGENESIS ASSAY USING THE ATCC® ANGIO-READY™ SYSTEM** (Chang, Cheng, and Shaper, 2010)
- COMPREHENSIVE GENE EXPRESSION ANALYSIS AND NEURAL PROGENITOR CELLS AND NEURONS** (Larson, et al., 2010)
- PRIMARY HUMAN DERMATOLOGICAL CELLS** (ATCC, 2010)
- WELL CHARACTERIZED, HIGH PERFORMANCE PRIMARY CELLS** (ATCC, 2010)

 Each article snippet shows an abstract, methods, and figures, illustrating the variety of cell models and assays available from ATCC.

