



Universal extracellular matrix supports toxicology and drug screening in 2D and 3D cell culture models from iPSC-derived cells, primary cells, engineered lines, and cancer lines

James Clinton, Ph.D., Penney McWilliams-Koeppen, M.S., Leelamma M Panicker, Ph.D., Siddhartha Paul, Ph.D., Chaozhong Zou, Ph.D., Joy Wells, B.S., and Dezhong Yin, Ph.D.
 ATCC Cell Systems, ATCC, 217 Perry Pkwy, Gaithersburg, MD, 20877 USA

ASCB Poster #: P877

Abstract

In vitro cell culture models are becoming increasingly complex in an effort to better mimic *in vivo* physiology and enhance their predictive power and relevance in areas including absorption, distribution, metabolism, excretion, and toxicity assays (ADMET), drug discovery, phenotypic screening, developmental biology, and basic research. One major advance in model development has been the recognition of the critical role of the cell microenvironment including the presence of extracellular biomolecules, cell-to-cell interactions, and physical properties of the substrate. These elements can influence a variety of cellular functions including cellular metabolism, differentiation potential, gene/protein expression, drug susceptibility, proliferation, and survival, typically with extracellular matrix (ECM)- and cell type-specificity. Herein we report the development of a "universal" Cell Basement Membrane ECM that provides a suitable microenvironment to support a wide variety of cell biology applications. Our results show that with optimized protocols the ECM permitted the long-term 3D culture of primary- and iPSC (induced pluripotent stem cells)- derived gastric and intestinal (GI) organoids from human and mouse cells and significantly enhanced spheroid forming efficiency in multiple continuous cancer lines. ATCC Cell Basement Membrane also supported angiogenic vessel formation in engineered, immortalized and primary human endothelial cells under co-culture conditions; increased cytochrome P450 enzyme activity and multidrug resistance-associated protein 2 (MRP2) expression in "sandwich" culture of primary human hepatocytes; and allowed the routine culture of iPSCs and NPCs (neural progenitor cells) under serum- and feeder-free conditions while maintaining and promoting the capacity for terminal differentiation into neurons and beating cardiomyocytes. These results demonstrate that ATCC Cell Basement Membrane provides a suitable biological matrix for advanced, physiologically relevant *in vitro* cell-based models and assays.

Methods

Extracellular matrix: ATCC Cell Basement Membrane was stored at -80°C prior to use. Protein concentration ranged from 12 to 18 mg/mL. Please refer to the certificate of analysis for the lot-specific concentration. To thaw, stock vial was placed at 4°C overnight and aliquoted and stored at -80°C. Immediately prior to use, Cell Basement Membrane aliquots were thawed and kept on ice.

Description	ATCC® No.	Size
Cell Basement Membrane; Growth factor reduced; LDEV-free	ACS-3035	5 mL

Hepatocyte culture: Cryopreserved primary human hepatocytes were thawed and plated on collagen coated 24-well plates in primary human hepatocyte media (Thermo Fisher Scientific). Six hours after plating the media was removed and an overlay of ATCC Cell Basement Membrane was added at a concentration of 250 µg/mL. Media was changed daily. Images were captured via phase contrast and immunofluorescent microscopy with a 10X objective.

iPSC and NPC culture: ATCC iPSCs were cultured according to the ATCC Stem Cell Culture Guide¹. NPCs were handled and differentiated as previously described². Both iPSCs and NPCs were routinely maintained on Cell Basement Membrane coated dishes and plates at a concentration of 150 µg/mL. ATCC iPSCs were differentiated to beating cardiomyocytes using published protocols³. Images were captured via phase contrast and immunofluorescent microscopy with a 10X or 20X objective.

Description	ATCC® No.	Size
Human Induced Pluripotent Stem Cells	Various	>1x10 ⁶ cells
Neural Progenitor Cells; Normal, Human	ACS-5003	>1x10 ⁶ cells
Neural Progenitor Cells; Normal, Human; MAP2-NanoLuc®- HaloTag®	ACS-5007	>1x10 ⁶ cells
Neural Progenitor Cell Expansion Growth Kit	ACS-3003	500 mL
Neural Progenitor Cell Dopaminergic Differentiation Kit	ACS-3004	250 mL

Organoid culture: Human iPSC-derived organoids were generated and cultured as described elsewhere⁴ and maintained in 100% Cell Basement Membrane in 3D. Primary murine small intestine-derived organoids were derived from LRG5+ stem cells cultured in 100% Cell Basement Membrane in 3D according to published protocols⁵. Images were captured via bright field and immunofluorescent microscopy or confocal microscopy with a 4X, 10X or 20X objective.

Spheroid formation: Continuous cancer cell lines were thawed and cultured according to the respective ATCC product sheet. In ultra-low attachment round bottom plates, 1000 cells per well were seeded in complete growth medium containing 2.5% Cell Basement Membrane. Spheroids appeared within 24 hours and were monitored for 96 hours. Images were captured via phase contrast microscopy with a 10X objective.

Description	ATCC® No.
MCF7; Mammary epithelial adenocarcinoma	HTB-22
A549; Lung epithelial carcinoma	CCL-185
MDA-MB-231; Mammary epithelial adenocarcinoma	HTB-26
Hep G2; Hepatocellular carcinoma	HB-8065

Angiogenesis: Primary or immortalized endothelial cells, some constitutively expressing GFP, were seeded on Cell Basement Membrane coated (10 mg/mL) multiwell plates as previously described [6]. Tubules formed within 6 hours and could persist for 10+ days in culture. Images were captured via fluorescent microscopy with a 4X objective.

Description	ATCC® No.	Size
Primary Umbilical Vein Endothelial Cells; Normal, Human	PCS-100-010	>0.5x10 ⁶ cells
Primary Aortic Endothelial Cells; Normal, Human	PCS-100-011	>0.5x10 ⁶ cells
TeloHAEC-GFP; immortalized human airway epithelial cells expressing GFP	CRL-4054	>1x10 ⁶ cells
TIME-GFP; immortalized human dermal microvascular endothelial cells expressing GFP	CRL-4045	>1x10 ⁶ cells

© 2016 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. NanoLuc®, HaloTag® and Promega are registered trademarks or trademarks of the Promega Corporation. Thermo Fisher is a trademark of Thermo Fisher Scientific, Inc.

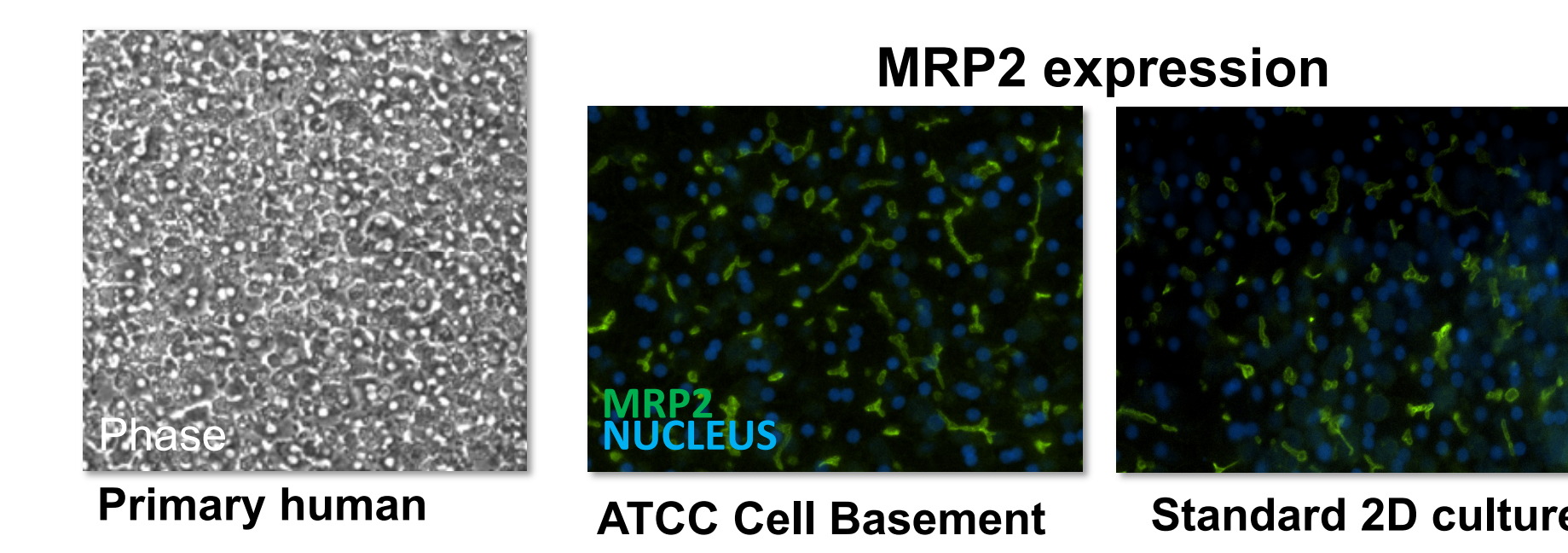
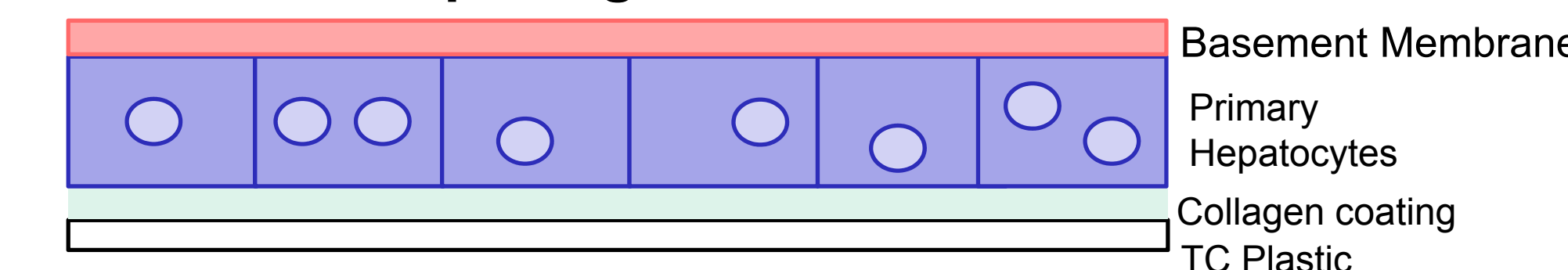
Results

ATCC Cell Basement Membrane is purified from the murine Engelbreth Holm Swarm sarcoma. Major components include laminin, collagen IV, entactin, and heparin sulfate proteoglycan. Cell Basement Membrane was found to support or enhance a broad range of *in vitro* 2D and 3D cell culture models for use in drug discovery, toxicology, and basic research by providing a more *in vivo*-like microenvironment for improved physiological functionality, response, and relevance.

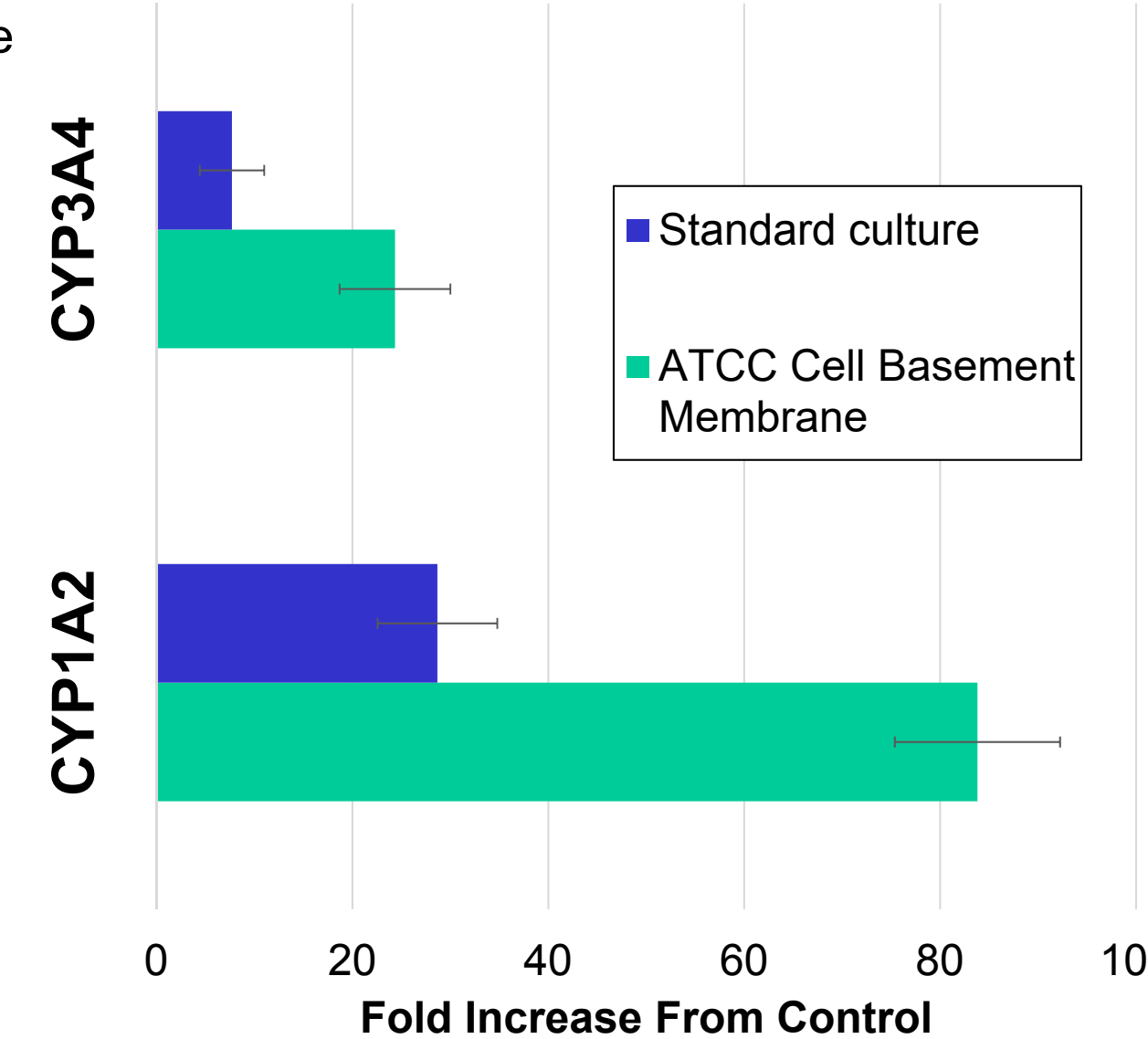
Hepatocytes

Sandwich culture of primary human hepatocytes using Cell Basement Membrane overlay provides an *in vivo*-like microenvironment that results in enhanced *in vitro* functionality.

Schematic depicting sandwich culture conditions



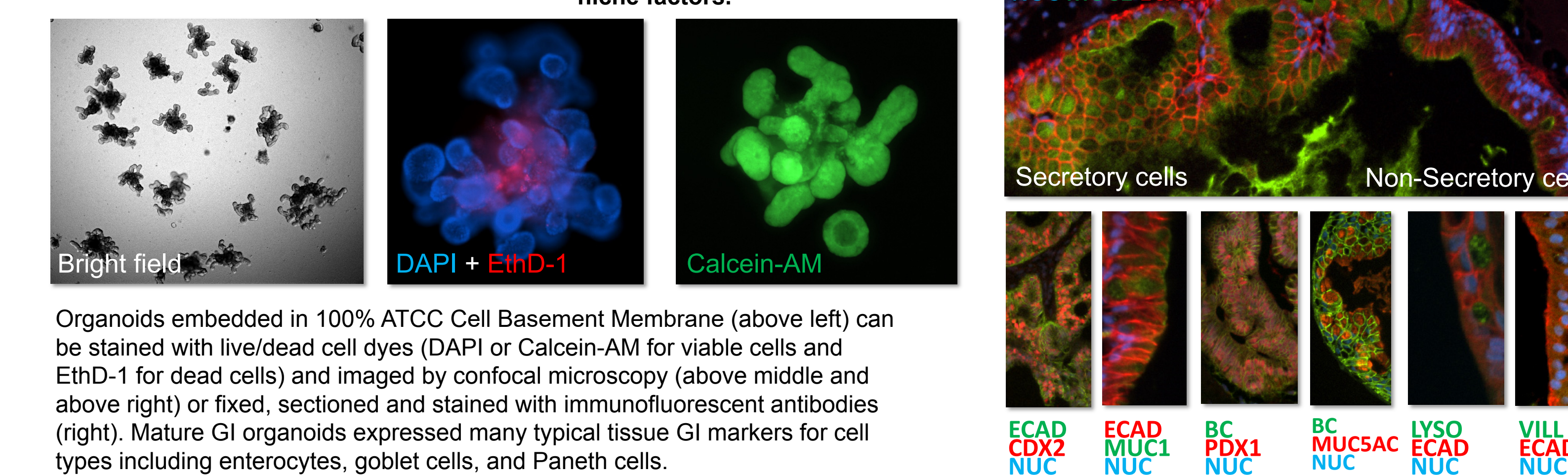
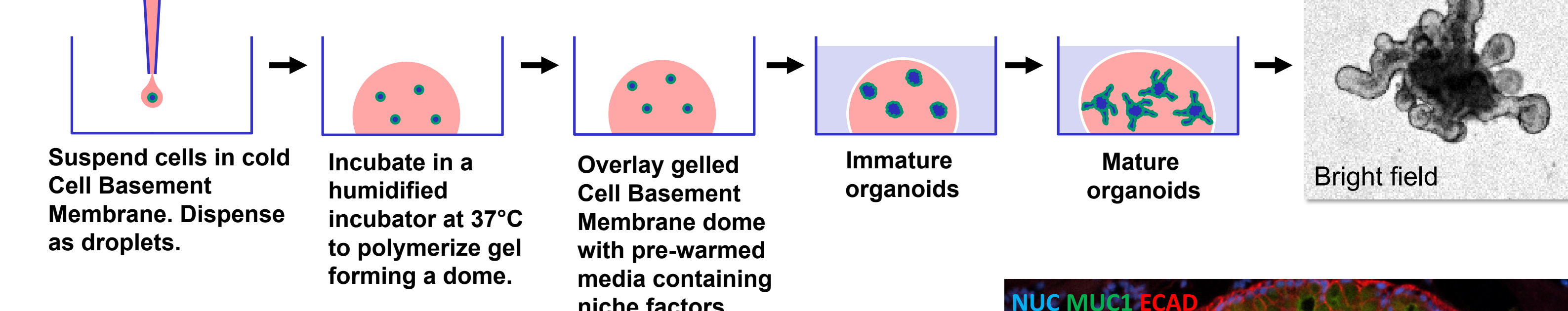
Cytochrome P450 Enzyme Induction



Organoids

GI organoids from primary tissue and multiple human ATCC iPSCs exhibited gastric- and intestine-like features after culture in serum free, defined media and embedded in Cell Basement Membrane.

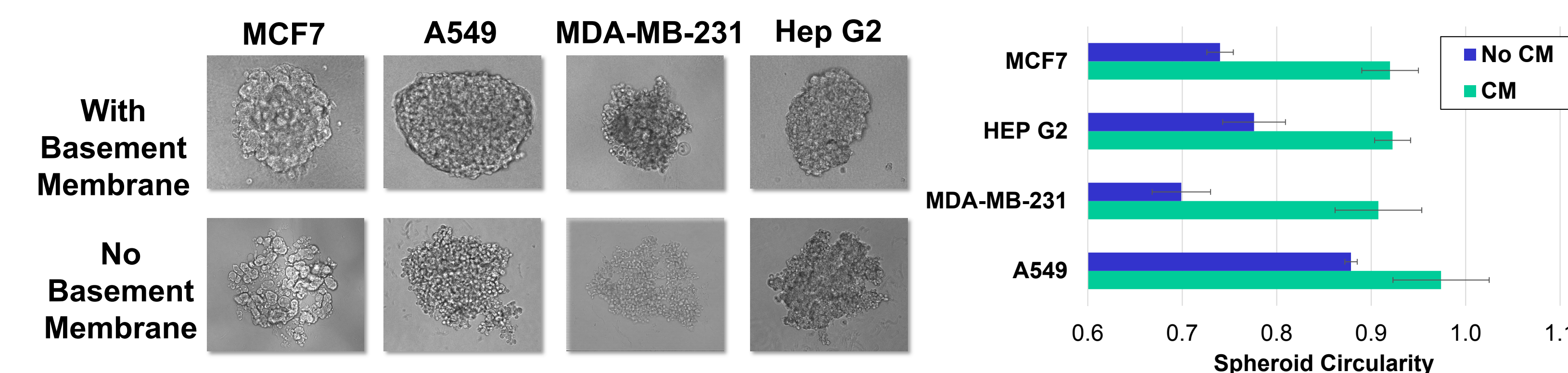
Embedded 3D culture of organoids in 100% ATCC Cell Basement Membrane



Organoids embedded in 100% ATCC Cell Basement Membrane (above left) can be stained with live/dead cell dyes (DAPI or Calcein-AM for viable cells and EthD-1 for dead cells) and imaged by confocal microscopy (above middle and above right) or fixed, sectioned and stained with immunofluorescent antibodies (right). Mature GI organoids expressed many typical tissue GI markers for cell types including enterocytes, goblet cells, and Paneth cells.

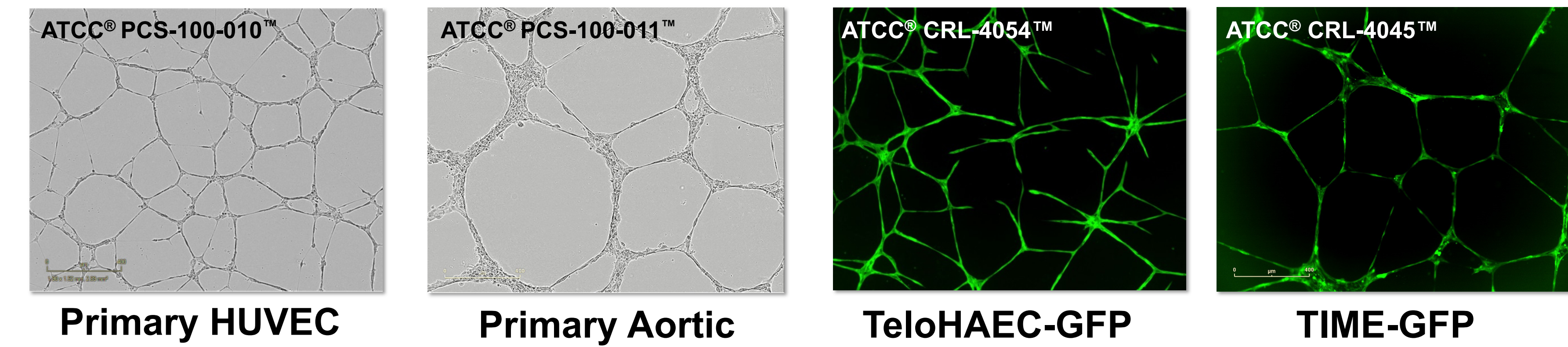
Spheroids

Spheroid formation is enhanced by the addition of Cell Basement Membrane in multiple cancer cell lines. Spheroids formed with Cell Basement Membrane are more spherical (below) and develop faster (data not shown) than when cultured alone.



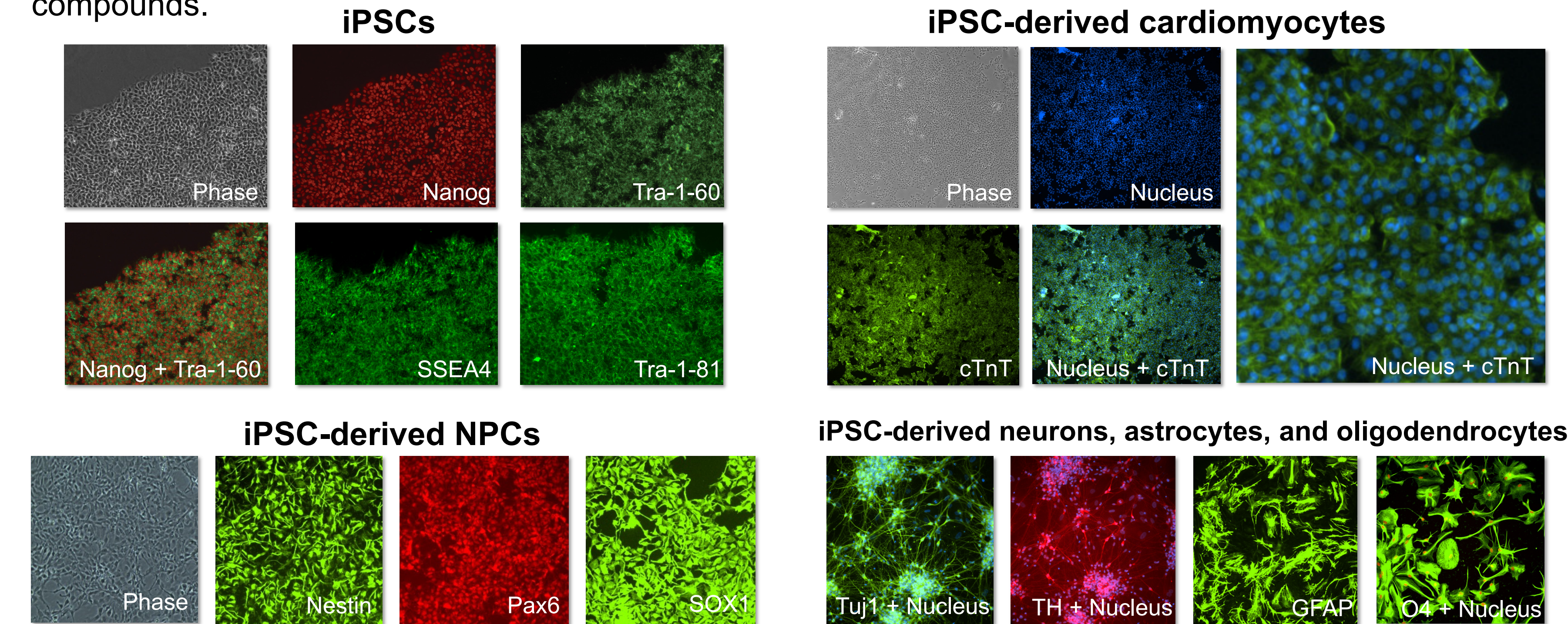
Angiogenesis

Primary, immortalized, and engineered endothelial cells constitutively expressing GFP form networks of tubules when seeded on Cell Basement Membrane. Angiogenesis can be stimulated or inhibited with known regulators such as VEGF and suramin (data not shown) and monitored in real-time.

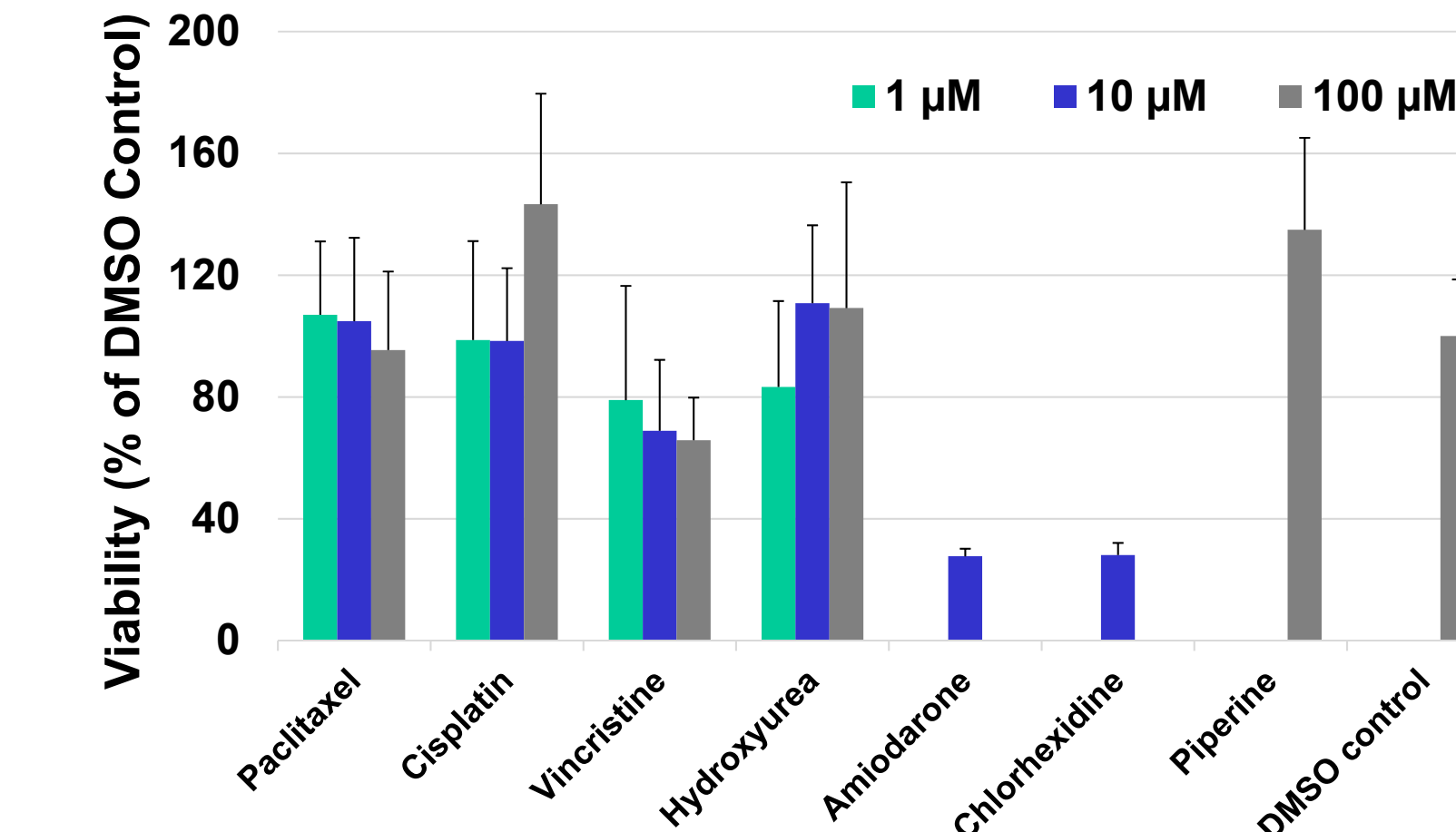


iPSC-derived NPCS, neurons, astrocytes, and cardiomyocytes

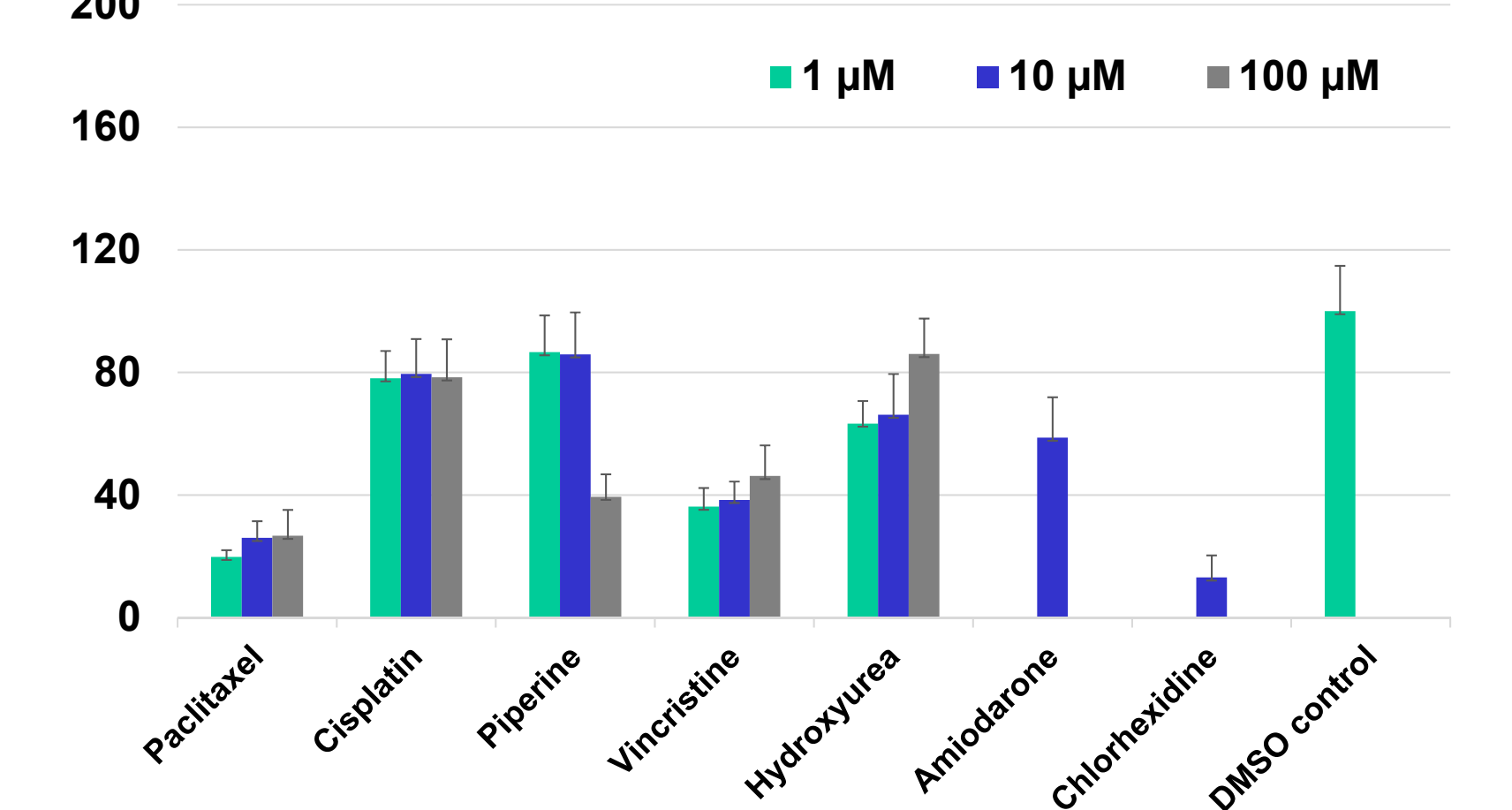
ATCC Cell Basement Membrane provided a supportive microenvironment for the xeno- and serum-free culture of iPSCs as well as iPSC-derived cell types including NPCS, neurons, and spontaneously beating cardiomyocytes. Differentiated cells (neurons vs. NPCS) show differential sensitivity to various compounds.



iPSC-derived NPC response to drugs



iPSC-derived neurons response to drugs



Summary

ATCC Cell Basement Membrane Gel supports numerous cell biology applications that are relevant for drug screening, toxicology, and basic research including hepatocyte culture, organoid generation, cancer research, 3D culture of cancer cells, and human stem cell maintenance and differentiation into neural progenitor cells, neurons, astrocytes, oligodendrocytes, and cardiomyocytes. Visit the ATCC website and our ASCB booth for additional details on the various Cell Basement Membrane applications described here.

References

- ATCC® Stem Cell Culture Guide: https://www.atcc.org/Guides/Stem_Cell_Culture_Guide.aspx
- ATCC Application Note 23, Comprehensive gene expression analysis and neurotoxicity testing of human iPSC-derived neural progenitor cells and neurons. <http://www.atcc.org>
- Burridge P W, et al. Chemically defined generation of human cardiomyocytes. Nat Methods 11(8): 855-860, 2014. PubMed: 24930130
- ATCC Application Note 26, Directed differentiation of gastrointestinal epithelial organoids using ATCC Cell Basement Membrane from multiple human ATCC iPSC lines. <http://www.atcc.org>
- Sato T, et al. Single Lgr5 stem cells build crypt villus structures *in vitro* without a mesenchymal niche. Nature 459(7244): 262-265, 2009. PubMed: 19329995
- ATCC Application Note 4, Cell Basement Membrane Gel supports *in vitro* angiogenesis assays. <http://www.atcc.org>



Scan code to download a PDF of this poster