August 21-23 | Boston, USA

PREDIOT:3DModels

Enhancing the Predictive Confidence of Models-Based Discovery & Development

38 Expert Speakers Including:



Silva Krause Senior Scientist - Translational Medicine Momenta Pharmaceuticals



Brian Berridge Associate Director National Institute of Environmental Health Sciences – (NIEHS)



William Mattes Director, Division of Systems Biology, National Center for Toxicological Research U.S. Food and Drug Administration



Kristin Fabre Microphysiological Systems Lead, IMED Biotech Unit AstraZeneca



Terry R. Van Vleet Head of Molecular and Computational Toxicology AbbVie

EXPERTISE PARTNERS:

DRAPER









INNOVATION PARTNERS:





EXHIBITION PARTNERS:







Tel: +1 212 537 5898Mail: info@hansonwade.comIn 3D Tissue Models: Drug Discovery & Development

ABOUT

Exploring The Utility Of 3D Models Across Discovery And Development

The 3rd PREDiCT: 3D Models (Previously "3D Tissue Models") remains committed to providing a forum that unites multidisciplinary stakeholders in overcoming the remaining translational, organizational and developmental barriers to 3D modeluse.

Driven by case studies from biopharma, PREDiCT: 3D Models looks beyond the science of tissue engineering. It will reveal how 3D models can and are being successfully deployed to expedite discovery and translation to clinical research.

So whether you've already established 3D models into your efforts or are just about to, tap into a ready-made network of peers who continue to redefine this area, solving the biggest barriers to model use.

I really do hope you can join a truly diverse group of professionals from biopharmaceutical, technology and academic institutions in redefiing the use of early stage and preclinical in-vitro models.

Adam Cohen Senior Conference Director **PREDiCT Event Series**

2

Key Case Studies Not to Miss



Expert

Speakers

1 The two previous 3D Model meetings have been highly successful and key contributions to the growth of this field and community of interest. The result has been a growing alignment on opportunities, challenges and strategies to building internal confidence in the applicability of these new capabilities

Brian R. Berridge, Associate Director, National Toxicology Program, Scientific Director, NIEHS

The 2017 Predict:3D Models conference was excellent. I came away feeling that I had been brought up to date on the latest technologies, by an impressive group of leaders in the pharmaceutical, academic and technology development industries. 3D models are developing rapidly, so I am excited to attend this year's conference, to learn of the progress that has been made over the last year and share my own experiences with these models as a speaker. Highly recommended!

James Morelli, Research Investigator, Sanofi

1 appreciated the very focused topic and the format of the meeting, which allowed for significant one-one discussions with other delegates and speakers John Westwick CEO. Resonant

I thought the conference was very focused. It was time well spent. I learned a lot and had productive networking sessions Hicham Alaoui, Senior Vice President, Discovery Biology, Symic Biomedical

The upcoming meeting really does feel like one of those 'can't miss' opportunities. Once again, it will bring together champions and users across industry, academia and government to define the applicability, and explore successes in utilizing more physiologically relevant systems to better identify therapeutic targets and predict drug toxicities

Yvonne Will, Senior Director, WRD Technology Strategy and Innovation, Pfizer Inc

3

🎒 RESONA

3rd PREDiCT: 3D Models August 21-23 | Boston, USA

WHY PAST ATTENDEES ARE RETURNING











.....

.....

.....





Adrian Roth Head of Mechanistic Safety & Non-**Clinical Safety** Roche



Ajamete Kaykas Senior Investigator Novartis



Alejandro Amador Scientific Leader GlaxoSmithKline

YOUR EXPERT SPEAKERS



John Wikswo Gordon A. Cain Professor & Director VIIBRE





Andreas R. Baudy Associate Principal Scientist, Safety Assessment Merck & Co.



Anita Seshire Head of Laboratory Cellular Pharmacology & Translational Innovation Platform Oncology **Merck KGgA**



Anna M Sitarski Post Doctoral Scientist **Cyteir Therapeutics**

Matthew Wagoner Associate Director Investigative Toxicology Takeda





B. Prabhakar Pandian Chief Technology Officer SynVivo Inc.

Brian Wamhoff

Co-founder and

Associate Director -

Preclinical Hepatic

AstraZeneca

Fabian Zanella

Director of R&D

StemoniX

Safety

HemoShear



Bhushan Mahadik Assistant Director NIH/NIBIB Center for Engineering Complex Tissues (CECT) **University of** Maryland



Brian R. Berridge Associate Director, National Toxicology Program, Scientific Director NIEHS

Dik C. van Gent Associate Professor **Erasmus MC**

Rhiannon Hardwick Research Scientist - Nonclinical Safety Assessment Theravance **Biopharma**





Swati Gupta Director, Immunology Non-Clinical and Translational Sciences Allergan



AbbVie

Terry R. Van Vleet Head of Molecular and Computational Toxicology



Yvonne Will Senior Director, WRD Technology Strategy and Innovation **Pfizer Inc**



1 really enjoyed the conference, and I learned a lot, both from presentations and informal interactions. Presentations were well balanced between science and technical innovation, and speakers focused on highlighting pros and cons of each application, which is helpful for someone just starting to explore the field of 3D models

Monica Gostissa, 3D Tissue Models: Oncology 2018 Attendee

Head of Innovation **Therapeutics Dominic Williams**

Darrell Kotton Seldin Professor of Medicine & Director, Center for Regenerative Medicine (CReM) **Boston University** and Medical Center



Eric McDuffie Scientific **Director & Head** of Investigative & Mechanistic Toxicology Janssen

Hicham Alaoui Senior Vice President, **Discovery Biology Symic Biomedical**





4

James Morelli Research Investigator Sanofi



Scientist Squibb

Jinping Gan Senior Principal **Bristol-Myers**

Graham Marsh

Scientist I -

Sciences

Biogen

Translational

Tel: +1 212 537 5898 Mail: info@hansonwade.com 3d-tissuemodels.com in 3D Tissue Models: Drug Discovery & Development

















John K. Westwick

President and CEO

3rd PREDiCT: 3D Models August 21-23 | Boston, USA

Joseph Charest Program Manager **Draper Laborator**



Kristin Fabre Microphysiological Systems Lead, IMED Biotech Unit AstraZeneca

Misti Ushio CEO **TARA Biosystems**



Piyush Bajaj Scientist II. Drug Safety Research and Evaluation Takeda

Sarah Hoyle Senior Research Scientist **Epistem Ltd**



Silva Krause Senior Scientist - Translational Medicine Momenta **Pharmaceuticals**

Seyoum Ayehunie Vice President of Immunological Systems MatTek Corporation



Szczepan Baran Global Head of Animal Welfare Compliance Training Novartis

Thomas Knudsen Researcher **US Environmental Protection Agency**



William Mattes Director, Division of Systems Biology, National Center for Toxicological Research **U.S. Food and Drug** Administration

Zhen Ma Assistant Professor **Syracuse University**



AGENDA AT A GLANCE

7

8am			Conference Day August 22nd, 8.20a	т 1 m-7pm		Conferen August 23rd	
			3D Systems Informing Disease Biology			Guidi &	
10am	August 21st, 9am-4	pm				Informin	
	Organs-on-Chips: A Developer's	3D Printing & Bioprinting for	Accelerating Target & Phenotypic Discovery	Enhancing the Confidence of Safety		Early S Discov	
12pm	Masterclass	Engineering	Discovery	Assessments		Towards	
2pm	3D Models for	Stem Cell Oraanoid Models		Predicting		Outco	
	ADME Tox Studies	for Drug	Optimising Lead Selection	Safety Liabilities In		Engi	
4pm		Screening		Vitro		12 Presen	
	Presenting Organizations Include:					Janssen 🕇	
	VANDERBILIT		Overcoming Core Challenges with 3D Models			Bristol-Myers Squibb	
6pm	MARYLAND	SUUS SUUS SUUS SUUS SUUS SUUS SUUS SUU	23 Presenting Orgo	anizations Include:			
			AstraZeneca	sk Roche			
8pm			SANOFI HERAP	SHEAR Takeda			

6

rence Day 2

23rd, 8.30am-3.30pm uiding Indication Selection & Biomarker Discovery

ming Your ly Stage scovery

Key Design Criteria for Safety Testing

rds In-Vivo tcomes

Advancing **DMPK** Profiling

Engineering Future Success

esenting Organizations Include:

UNOVARTIS

Biogen



Pre-Conference Workshop Day | August 21st

Workshop A

Organs-on-Chips: A Developer's Masterclass

9.00am - 12.00pm

Following the success and overwhelming feedback from the session in the last two event iterations. This workshop will be back once more to provide an in depth analysis of how to approach and develop microfluidic devices with a focus on interconnected systems and their use within industry.

This workshop will help you realize the unlimited potential of engineered tissue models for replicating organ system signaling and dynamics.

Workshop Leader

John Wikswo Gordon A. Cain Professor & Director VIIBRE

Leave this workshop with:

- A greater understanding of the key challenges facing the development of microfluidic systems and how to overcome them
- An in depth insight into how to model, interconnect, test, and control these biological systems consistent with allometric, biochemical and/or functional scaling
- Knowledge of how best to achieve noninvasive readouts with the latest advances in imaging, sensor, and multi-omic technologies

I really enjoyed the conference, and I learned a lot, both from presentations and informal interactions. Presentations were well balanced between science and technical innovation, and speakers focused on highlighting pros and cons of each application, which is helpful for someone just starting to explore the field of 3D models

Monica Gostissa, 3D Tissue Models: Oncology 2018 Attendee



Workshop B

3D Printing & Bioprinting for Complex Tissue Engineering

9.00am - 12.00pm

Bioprinted tissue constructs have great potential as platforms to improve our understanding of cell and tissue-level biology for various applications. However, numerous challenges exist in adopting and utilizing the variety of technologies and approaches within development workflows. The objective of this workshop is to introduce tools and techniques related to 3D Printing and Bioprinting in order to fabricate complex tissue architecture that can serve as in vitro models capable of recapitulating native biology.

Workshop Leader Bhushan Mahadik

Assistant Director NIH/NIBIB Center for Engineering Complex Tissues (CECT) **University of Maryland**

Leave this workshop with:

- An in-depth insight into a variety of tools and techniques in additive manufacturing for generating tissue models
- Exploration into issues and strategies to mitigate material selection and printability
- A greater understanding of 3D printing and Bioprinting principles for getting started with complex in vitro models developed

Workshop C

Leveraging 3D In-vitro Models for ADME-Tox Studies

1.00pm - 4.00pm

Active drug metabolites and drug-drug interactions are attributed for a large percentage of drug failures. As such, representative ADME studies are essential to the drug discovery process and yet current models fall short on their pharmacokinetic profile, largely due to their over simplicity and lack of physiological relevance. Better recapitulating the physiological and metabolic properties of the in vivo environment, 3D models represent significant advantages for detection of drug induced toxicities.



interest 🗾

Brian R. Berridge, Associate Director, National Toxicology Program, Scientific Director, NIEHS, PREDiCT:3D Models 2016, 2017 & 2018 Meeting Chair



Workshop D

Development of Stem Cell Organoid Models for Drug Screening Applications

1.00pm - 4.00pm

Human in vitro cardiac models predictive of human drug response would be a significant advancement for understanding and developing new drugs. Recent technical advances in human iPSC-derived 3D organoids have presented huge opportunities to bridge the in vitro and in vivo gap. Providing insights, beyond basic descriptors at the cellular level, these models are offering new insights into complex organ development and have vast applications within pharmaceutical research. With specific reference to 3D hiSPC cardiac organoids, this session will look to explore the technical advances in designing these models, applications within the basic and pharmaceutical setting as well as strategies for overcoming remaining issues to use.



Zhen Ma Assistant Professor **Syracuse University** This workshop will review current advances in 3D model development and use for ADME-tox with reference to single and coupled 3D microfluidic systems

Leave this workshop with:

- Novel approaches utilizing chip based models for metabolite identification and enzyme mapping
- Strategies utilizing 3d microtissues in the clearance prediction of low turnover compounds

1 This was a great scientific meeting and networking opportunity. It brought together key players in the field to discuss important strategic issues of mutual



Leave this workshop having:

- Envisioned the integration of hiPSC biology, tissue engineering and gene editing technologies to develop next-gen 3D tissue models and organ-onchip systems
- Reviewed current biofabrication technologies for creating 3D tissue models for high-throughput and high-content preclinical drug evaluation
- Highlights on the integration of in vitro model with in silicon model for multifaceted data analysis and drug toxicity prediction.





Conference Day One | August 22nd 2018

	Brian R. Berridge Associate Director, National Toxicology Program, Scientific Director NIEHS	8.20	Chairs Opening Remarks
	Terry R. Van Vleet Head of Molecular and Computational Toxicology AbbVie	8.30	 IQ Consortium Microphysiological Systems Working Group: History & Updates Describing the consortium and key past contributions Reviewing current changes and insights into ongoing activities of the consortium Highlighting future direction & objectives
ŀ	Adding Human Re	levanc	y into the Pipeline – 3D Systems Informing Disease Biology
F	Joseph Charest Program Manager Draper Laboratory	9.00	 Leveraging Microfabrication for Tissue Models with a Path to Validation Highlighting how microfabricated models can influence and quantify tissue or function Illustrating two examples of microfabricated models: A complex in vitro model to culture and evaluate tissues or functions with barrier function A system to quantify and evaluate interactions between patient-derived tissues and immune components
	Darrell Kotton Seldin Professor of Medicine & Director, Center for Regenerative Medicine (CReM) Boston University and Medical Center	9.30	 Pluripotent Stem Cell Models of Lung Disease in 3D Review the state-of-the-art methods in using patient specific iPSC's for lung epithelial differentiation in 3D Application of these models for understanding airway and alveolar lung diseases Developing drug screens and precision medicine approaches using iPSC lung cell models
	Fabian Zanella Director of R&D StemoniX	10.00	 Adding Structure to HTP hiPSC-Derived Cardiac and Neuronal Models for Improved Physiological Relevance and Performance. microBrain 3D and microHeart are iPSC-based screening platforms that have been structurally engineered with greater physiological relevance aimed to elevate performance in drug discovery and development applications Both platforms are tested and validated for performance in screening and safety applications Case studies are provided for the successful implementation of these platforms in screening and safety applications
10.30	Speed Networking		

11.00 Morning Refreshments

10

Discovery Track

Accelerating Target & Phenotypic Based Discovery

- 11.30 Creating Physiological Disease Models for Drug Discovery
- Utilizing a novel disease relevant 3D model to interrogate disease pathways and identify new targets for rare metabolic diseases
- Highlighting the use of parallel in-silico models to generate rationally driven hypotheses and drive the discovery of more relevant data
- Exploring concepts and methods applied for model validation
- Brian Wamhoff, Co-founder and Head of Innovation, HemoShear Therapeutics

Development Track

Enhancing the Predictive Confidence of Preclinical Safety Assessments

- 11.30 Application of 3D Hepatic Cultures for Risk Assessment of Drug Induced Liver Injury
- Recapitulating healthy and disease liver statesengineering models to identify patient specific toxicities
- Exploring the capabilities of model-incorporated cell types and their impact on therapeutic and DILI assessment
- Demonstrating cross validated platforms utilized within AstraZeneca's pipeline and its potential implications for DILI assessment of novel therapeutic modalities **Dominic Williams,** Associate Director - Preclinical Hepatic Safety, **AstraZeneca**

12.00 Advancing Drug Discovery Through the Technological Convergence of Organs-On-Chips & Mass Spectrometric Multi-Omics

- Exploring how microfluidics and untargeted multiomic reconstruction of drug MoA, through advanced mass spectrometry and machine learning, are set to revolutionise discovery
- Highlighting its potential impact in revealing problematic human haplotypes and drug-drug interactions for small molecules
- Demonstrating its utility in the discovery of novel mechanisms of human diseases, identification of novel compounds, and the discovery of on- and off-target of drug candidate MoAs

John Wikswo, Gordon A. Cain Professor & Director, VIIBRE

12.30 Active-Learning Strategies for High Content Screening of 3D Tumor Cell Models

- Introducing novel 3D tumor in vitro models, high content imaging and cell-based screening technologies improving clinical relevance in the testing of oncology drugs
- Showcasing the design and build of an intelligent, automated platform to accelerate the identification of high value small molecule leads, and the delivery of quality candidates into clinical development
- Discussing computational active-learning approaches used in high content screenings of a large number of 3D tumor spheroid images
- Alejandro Amador, Scientific Leader, GlaxoSmithKline

1.00 Lunch

Novel Screening Applications to Optimise Lead Selection

2.00 Bringing 3D Cell Culture into Cancer Drug Development & High Throughput Drug Screening

- High throughput assays in 3D drug screening
- Spheroids as a next step from 2D to 3D in assay development in chemotherapeutics
- Scaffolds (biomaterials) into the high throughput world
- Challenges in characterizing a system, measurements, reproducibility and expense

Anna M Sitarski, Post Doctoral Scientist, Cyteir Therapeutics

12.00 Intestinal 3D organoids – A Predictive Model of Intestinal Toxicity

- Brief history and description of intestinal organoids and how they compare to in vivo structures
- Overview of the species from which organoids have been made
- Examples of toxicity studies in different species

Sarah Hoyle, Senior Research Scientist, Epistem Ltd

- 12.10 Exploring the Utility of Spheroids for Improving the Prediction of Hepatotoxicity Studies
- Describing an alternative method of spheroid culture that results in more spheroid micro-tissues per well, compared to hanging drop or single well methods
- Presenting data comparing rat and human hepatic spheroids, allowing for species-specific comparisons of hepatotoxicity in vitro to known outcomes in vivo
- Reviewing this methodology in comparison with others, giving a brief review of best practices, and the predictive power of 3D hepatic models over traditional models

James Morelli, Research Investigator, Sanofi

12.40 Primary Human Endothelial Cells-Based 3D Models & Assays for Detection of Innate Immune Responses

- Comparing models and assays to monitor inflammation responses following interactions with host proteins and impurities.
- Developing vascularized environments for investigation of immune cell behaviour
- Demonstrating the ability for real-time visualization and quantitation of vascular leakiness

Swati Gupta, Director, Immunology Non-Clinical and Translational Sciences, Allergan

12.50 iPSC Derived Mature Cardiac Tissue for Drug Development

- Description of BiowireTM II Platform
- Characteristics of Cardiac Tissue Maturity
- Applications for Drug Development

Misti Ushio, CEO, TARA Biosystems

Improving In-Vitro Prediction of Safety Liabilities

2.00 Adapting 3D Microtissues for The Detection & Mechanistic Evaluation of Drug-Induced Toxicity

- Exploring how to model clinically relevant tissue responses through the generation of a 3D liver model with architecture closely resembling that of in vivo human tissue
- Confirming viability and functionality over prolonged culture periods as well as recapitulation of clinical liver toxicity not detected in conventional models
- Demonstrating how 3D liver tissues can both effectively model DILI and distinguish between highly related compounds with differential profiles
- Adrian Roth, Head of Mechanistic Safety & Non-Clinical Safety, Roche



2.30 High-Throughput Antibody Identification Based On the Tumor Microenvironment	2.30 Novel 3-D Primary Human Gut Model to Predict Drug Toxicity and Host-Microbe Interaction	5.00 Round Challenge Tables
 Highlighting the development of novel tumor microenvironment models, and Resonant's approach linking target and therapeutic candidate identification and functional validation Combining high-throughput fidelity phenotypic screens and machine learning strategies to accelerate the process, and result in unprecedented speed in the identification of novel, efficacious anti-tumor monoclonal antibodies Explore case studies demonstrating the process and resultant therapeutic candidates currently in development John K. Westwick, President and CEO, Resonant Therapeutics, Inc. 	 Drug toxicity results from 3D-Intestinal tissues correlate better to human responses compared to standard preclinical animal models. 3D intestinal tissues showed adhesion and invasion by symbionts and pathogenic microorganisms. Microbes or microbial components initiate innate immune responses in of 3D intestinal tissue model that mimic human responses. Seyoum Ayehunie, Vice President of Immunological Systems, MatTek Corporation 	 Building Confidence in the Utility of 3D When can studies from 3D models be considered critical path? How should we most effectively be demonstrating the internal business case? Does 3D always mean more predictive? How should we define the context of use to address this? What technolog
 3.00 Utilising Multicellular 3D Models for Preclinical Drug Discovery Developing primary 3D cell cultures from patient material Addition of monocytes or fibroblasts to identify a potential trophic role Transplantation of multicellular spheres to create tumours with microenvironments and enable the exploration of the influence of the TME on stemness Anita Seshire, Head of Laboratory Cellular Pharmacology & Translational Innovation Platform Oncology, Merck KGgA 	 3.00 Microphysiological Systems in Drug Discovery: Advances in CNS & Immune-mediated Liver Injury Sharing our ongoing efforts to validate microbrain and minibrain models with clinical data from 100 medicines. Presenting data on our novel medium throughput model of immune DILI complete with lymphocytes, Kupffer cells, endothelial cells & hepatocytes. Deficiencies in our current immunotoxicity models will also be presented in the hopes that someone in the audience has a solution to what seems to be a universal problem Matthew Wagoner, Associate Director - Investigative Toxicology, Takeda 	 How should we be working with external collaborators for PoC studies? How should we address issues of robustness and reproducibility across lab environments? Hicham Alaoui, Senior Vice President, Discovery Biology, Symic Biomedical Human Based vs Animal Models for Decision-Making: Pros And Cons Brian Wamhoff, Co-founder and Head of Innovation, HemoShear Therapeutics
5.30 Extended Q&A	 3.30 Physiologically Relevant 3D Tissue Models for Translational Studies Utilizing vascularized tumor microenvironment to evaluate cancer-stromal-immuneendothelium interactions Showcasing a blood brain barrier model for real-time visualization and quantitation of cellular interactions and barrier functionality Using an integrated assay for mechanistic understanding of vascular-immune-tissue interactions Assessing the use of layered multi-cellular architecture for complex organ physiology and toxicology B. Prabhakar Pandian, Chief Technology Officer, SynVivo Inc. 	6.00 Round Table Moderator Feedback Panel 6.15 Close of Conference Day One Poster Session After the formal presentations have finish is an informal part of the conference ager atmosphere and continue to forge new are be presented on 3D models validating inn

Thomas Knudsen Researcher **US Environmental Protection Agency**

12

Computer Modeling & Simulation to Reconstruct the Basis of 4.30 **Developmental Toxicity**

13

3rd PREDiC	T: 3D Models
August 21-23	Boston, USA

Overcoming	the	Strategies to Mitigate		
Analytical Hurdles for		Cell Sourcing and		
Testing Endpoints In		Maturity Issues		
 Testing Endpoints In Organotypic Cultures How can we best adapt the low volumes to analytical techniques such as transcriptomics and proteomics? How can we best optimise readout methods for sensitivity? What technologies can we apply? Szczepan Baran, Global Head of Animal Welfare Compliance Training, Novartis 		 Maturity Issues What are the benefits/risks of isolating key cell types within you organization versus commercial sourcing? New opportunities to acquire "hard to isolate" or unavailable cell types. How do "mature" cells need to be before incorporation into a model, and what are the opportunities for cell maturation within a system? Setting clear and manageable cell acceptance criteria to drive reproducibility. Rhiannon Hardwick, Research Scientist - Nonclinical Safety Assessment, Theravance 		
s for	What Constit	utes a Validated Model: An		
ns	Industry Pers	pective		
nnovation	How should defin	ne the standards for these		
nnovation,	 Can we develop set compound libraries to testmodels against? 			
	 Where can we establish confidence in both specificity and sensitivity to identify relevant endpoints 			
	 What key compo developing a pre 	at key components should be considered when veloping a predictive 3D model		
Yvonne Will, Ser and Innovation,		or Director, WRD Technology Strategy		

resentations have finished, the learning and networking carries on. The Poster Session of the conference agenda, allowing you to connect with your peers in a relaxed continue to forge new and existing relationships. During this session scientific posters will 5D models validating innovative targets, disease models highlighting novel mechanisms, nforming investigative toxicology and safety assessments.



Conference Day Two | August 23rd 2018

0	Brian R. Berridge
125	Associate Director,
~	National Toxicology
	Program, Scientific
	Director
	NIEHS

8.25 Chairs Opening Remarks

3D Models Guiding Indication Selection & Biomarker Discovery

	Silva Krause Senior Scientist - Translational Medicine Momenta Pharmaceuticals	0.0
696	Dik C. van Gent	9.0

Erasmus MC

8.30 Utilizing a 3D Tissue Platform from Bench to Bedside for Mechanistic **Studies & Identifying Translational Biomarkers**

- · Demonstrating the use of a 3D model for understanding the MoA of a multi-targeted drug · Highlighting comparative similarities/differences and validation of animal models
- · Revealing the potential utility of the system in biomarker ID by referencing how protein

alterations within the system were representative of changes observed in patient plasma samples

0 Predicting Clinical Outcomes Through the Development 3D Assays on **Tumor Biopsies**

- Revealing how 3D tumor tissue slices can be cultured and treated as ex-vivo as a predictive model for therapy response
- · Clinical validation of ex-vivo assays on tumor biopsies is ongoing

· Exploring how the miniaturization and standardization in a cancer-on-chip set up may enable truly personalized medicine (testing multiple treatments, co-clinical trial)

Morning Refreshments Networking 9.30

Discovery Track

In Vitro Platforms Informing **Early Stage Discovery Efforts**

- **10.30** Interrogating Phenotypic Changes of 3D Neural **Models to Discover & Validate Biomarker** Candidates
- · Exploring contemporary microphysiological models to generate discovery compound safety profiles
- Leveraging imaging technologies and in vitro models to inform taraet assessment
- Presenting case studies: Monitoring for translational safety biomarker changes in vitro
- Eric McDuffie, Scientific Director & Head of Investigative & Mechanistic Toxicology, Johnson & Johnson

11.00 Using Stem Cells for Neuroscience Target Discovery

- Discussing the generation of a large bank of iPSC lines from schizophrenic and autistic patients
- · Highlighting strategies for the large-scale differentiation of these patient lines into cortical excitatory neurons
- · Exploring the characterization of the progenitors and neurons with omics-based technology, one which being single cell RNA-seq
- Ajamete Kaykas, Senior Investigator, Novartis

11.30 Microfluidic Models of the Blood Brain Barrier

- Exploring how researchers are building neuronal models with the goal of increasing the complexity and physiological relevance of the microphysiological models we are developing
- Reviewing novel platform technologies that are being developed to increase the throughput and scalability of our models

Graham Marsh, Scientist I - Translational Cell Sciences, Biogen

14

Development Track

Design Criteria for Optimised Safety Efficacy Testing

- 10.30 Key Criteria of Advanced 3D Human Liver Models for the Improvement of Safety Testing and Widespread Adoption
- Demonstrating how in-vivo human biological metrics must be taken into account when building an advanced in vitro model
- For liver, rates of albumin production and urea synthesis are important first considerations
- More complex is not always best. Showcasing how models that can answer key scientific questions while maintaining reasonable throughput will be at the forefront of future research

Andreas R. Baudy, Associate Principal Scientist, Safety Assessment, Merck & Co

11.00 Designing iPSC Derived Organoids for Cardiac Safety **Efficacy Testing**

- Highlighting the progress and promise human iPSCs with 3D models hold for precision medicine and companion diagnostics applications
- Demonstrating the development of a 3D cardiac organoid model mimicking early human heart development
- Showcasing its utility as a drug testing platform to evaluate the embryotoxicity of pregnancy medication to fetal development
- Engineering a fiber-based human cardiac tissue model to recapitulate more adult-like cardiac tissue structure for drug screening and disease modelling purposes
- Zhen Ma, Assistant Professor, Syracuse University

11.30 Extended Q&A

12.00 Lunch

Enhancing The Fidelity of In-Vitro Models Towards In-Vivo Outcomes

1.00 Microfluidic Technologies Impacting how AstraZeneca Optimizes Drug Scheduling

- Understanding how we are "scheduling discovery" to optimize dosing regimens for efficacy and safety
- · Exploring the use of dynamic in vitro systems can allow us to better understand efficacy and safety for drug combination studies
- Innovating capabilities to maximize output from MPS studies

Kristin Fabre, Microphysiological Systems Lead, IMED Biotech Unit, AstraZeneca

1.30 Advanced Predictive In-Vitro Liver & Kidney **Models for Drug Discovery**

- Opportunities for advanced predictive 3D models to address some of the challenges faced by the pharma industry
- · Impact examples of advanced liver and kidney models over conventional 2D static culture platforms Piyush Bajaj, Scientist II, Drug Safety Research and Evaluation, Takeda

2.00 **Afternoon Refreshments**

Mastermind Summary Panel: Engineering Future Success for 3D Models 2.30



Microphysiological Systems Lead, IMED

Szczepan Baran Training Novartis

3.00 Mastermind Discussion Groups - Collaborating To Advance Model Uptake

Following the conclusion of the summary panel the room will split into smaller groups to individually discuss and define the below points. The key aim of this session is to facilitate in-depth discussions between participants in an informal environment. Responses from each group will be collated and shared following the event.

- · What is the best way to establish effective communication strategies to advise model use in discovery and development?
- e.g tox to discovery?
- What is the most effective way to collaborate and advance the field?

3.30 **Chairs Closing Remarks**

Close of Conference 3.35

15

Advancing Drug Metabolism & Pharmacokinetic Profiling

- 1.00 Recent Advances in DMPK Applications of 3D **Tissue Models**
- Highlighting the real advances in the development and application of 3D models for DMPK during the last few years
- Showcasing areas of application including PK prediction of low clearance compounds, and complex metabolism and transport of drugs in multiple organs of interest
- Emerging trends for modular, higher throughput, and imaging compatible tissue models

Jinping Gan, Senior Principal Scientist, Bristol-Myers

- **1.30 ADMET Modeling Via Linked Microphysiological** Systems
- HUnderstanding the role of multi-step processes in druginduced toxicities
- Revealing insights into gene X environment interactions and inter-individual susceptibility
- Edward Kelly, Associate Professor, University of Washington

Global Head of Animal Welfare Compliance



Matthew Wagoner Associate Director -Investigative Toxicology Takeda

· How can different disciplines interact more effectively to leverage advances and lessons learned from early successes

· How should we be working external providers to facilitate model use and development within the pharma setting?



WHO ATTENDS?

PREVIOUS ATTENDEES INCLUDE



*Expected attendance figures based on actual attendance from past PREDICT: 3D Models events (2016-18)

▲ There was a diverse set of presentations, covering multiple aspects of challenges and opportunities setting up and using 3D models. A lot of attention was given to time for networking and exchanging experiences. All in all, a very efficient and useful conference

Marleen van Loenen, Director Non Clinical Development, Gadeta Bv 3D Tissue Models: Oncology Attendee

BECOME A PARTNER

16



Mo Langhi Partnerships Director **T:** +1 212 537 5898 **E:** sponsor@hansonwade.com

Tel: +1 212 537 5898Mail: info@hansonwade.com3d-tissuemodels.comin 3D Tissue Models: Drug Discovery & Development

DRAPER

Expertise Partner

Draper is an independent, not-for-profit corporation, which means its primary

commitment is to the success of clients' missions rather than to shareholders. For either government or private sector customers, Draper leverages its deep experience and innovative thinking to be an effective engineering research and development partner, designing solutions or objectively evaluating the ideas or products of others. Draper will partner with other organizations...

www.draper.com

SYNVIVO

Program Partner

SynVivo offers 3D cell-based models that enable real-time study of cell and drug interactions

and accelerates discovery by providing a biologically realistic platform that more accurately depicts in vivo reality. Complex in vivo microvascular environment including scale, morphology, hemodynamics and cellular architecture are re-created in an in vitro format. SynVivo's 3D models for blood brain barrier, tumor, inflammation and toxicology are morphologically and physiologically realistic and feature a side-by-side architecture enabling real-time visualization using standard analytical instrumentation

www.synvivobio.com

TARA

Innovation Partner

TARA Biosystems, Inc. provides predictive, in vitro human cardiac tissue models for use in drug

discovery and safety assessment. TARA Biosystems offers a high fidelity solution that is based on human stem cell-derived cardiac tissue matured to physiologically relevant adult-like levels and provides direct measures of cardiac functionality, including contractile force. The TARA solution will provide safety data earlier in the development cycle and a high fidelity discovery model. www.tarabiosystems.com

Exhibition Partner



AxoSim's Nerve-on-a-Chip technology predicts clinical results from the benchtop, helping pharmaceutical companies develop safer and more

effective drugs. Employing micro-engineering techniques and novel biomaterials, AxoSim has developed a 3D iPSC-based model which mimics living tissue in both form and function, enabling in vivo metrics from an in vitro platform. The Nerveon-a-Chip platform delivers high-content data as a more accurate, less expensive, and faster alternative to animal testing.

www.axosim.com

17

PROUD TO PARTNER WITH

Stemoni**X**

Expertise Partner

Our mission is to accelerate the discovery of new medicines with physiologically-relevant,

structured human cells in high density plates. We develop and manufacture human induced pluripotent stem cell-derived cardiac and neuronal platforms for high throughput drug toxicity and efficacy screening. Predictive, accurate, and consistent, our human models enable scientists to economically conduct research in a simplified workflow...

www.stemonix.com



Program Partner

MatTek Corporation is a global leader in the production of the highest quality in vitro human tissue models for safety and

efficacy testing for the chemical, cosmetic, household product, and pharmaceutical industries. For over 25 years, MatTek has established a reputation for the consistent manufacture of quality in vitro human tissue models, and the continuous development of innovative tissue engineering technologies and assays.

www.mattek.com.



Innovation Partner

Epistem provided specialist preclinical and discovery solutions for drug development. The company applies its

knowledge of epithelial tissue and adult stem cells to the area of epithelial repair, oncology, mucositis, inflammation and dermatology and has innovative platforms for both biomarkers development and molecular diagnostics.

www.epistemservices.com



Exhibition Partner

East River BioSolutions is pioneering novel tissue-specific extracellular matrix (ECM) biomaterials to meet research needs in stem cell and

cancer research, drug testing, tissue engineering, and regenerative medicine. East River Bio's TissueSpec[™] ECM products are commercially available in the form of matrix hydrogels, and scaffolds for 3D cell culture applications and media supplements, and surface coatings for 2D cell culture applications. Standard and custom-order TissueSpec[™] Matrix Kits are available for in vitro life sciences research.

www.eastriverbio.com



Exhibition Partner

OlympusOlympus Scientific Solution Group has been providing various imaging solutions adapted to customer needs. With advanced

optical technology crafted over a century, we can accurately visualize cells without losing important information in 3D samples, which have emerged as critical tools for physiologically relevant drug discovery. As a companion technology, Olympus provides true 3D cell analysis to quantitatively measure 3D samples.

www.olympus-lifescience.com/en



3rd PREDICT: 3D Models August 21-23 | Boston, USA

READY TO REGISTER?

3 EASY WAYS TO BOOK

www.3d-tissuemodels.com

Tel: +1 212 537 5898

Email: register@hansonwade.com

Identify the practical applications of 3D models in drug discovery and development

Leverage the expertise in the room to streamline model on-boarding and avoid errors in PoC studies

Experience a greater volume of case studies to enhance discovery efforts and identify early stage safety concerns

SECURE YOUR PLACE

Model & Solution

Provider Pricing

Conference + 2 Workshops

Conference + 1 Workshop

Conference Only

Workshop Only

Industry & Academic Pricing	Register & Pay Before 20th July 2018	Standard Prices
Conference + 2 Workshops	\$2,897 (Save \$400)	\$3,097
Conference + 1 Workshop	\$2,598 (Save \$300)	\$2,798
Conference Only	\$2,099 (Save \$200)	\$2,299
Workshop Only	\$59	9

Register & Pay Before

20th July 2018

\$3,797 (Save \$400)

\$3,298 (Save \$300)

\$2,799 (Save \$200)

Team Discounts*

- 10% discount 2-3 delegates
- 20% discount 4 or more delegates

Please note that discounts are only valid when three or more delegates from one company book and pay at the same time. Contact:

register@hansonwade.com

The 3D Tissue Models conference was a fantastic platform for industry and academics to come together and discuss how this field should and will move forward Moody Suliman, Nortis 3D Tissue Models 2016/2017 Attendee

 Image: Contract of the second seco

\$599

Standard Prices

\$3,997

\$3,498

\$2.999

TERMS & CONDITIONS

Full payment is due on registration. Cancellation and Substitution Policy: Cancellations must be received in writing. If the cancellation is received more than 14 days before the conference attendees will receive a full credit to a future conference. Cancellations received 14 days or less finctuding the fourteenth dayl prior to the conference will be liable for the full fee. A substitution from the same organization can be made at any time. Changes to Conference & Agenda: Hanson Wade reserves the right to postpone or cancel an event, to change the location or alter the advertised speakers. Hanson Wade is not responsible for any loss or damage or costs incurred as a result of substitution, alteration, postponement or cancellation of an event for any reason and including causes beyond its control including without limitation, acts of God, natural disasters, sabotage, accident, trade or industrial disputes, terrorism or hostilities.

Data Protection: The personal information shown and/or provided by you will be held in a database. It may be used to keep you up to date with developments in your industry. Sometimes your details may be obtained or made available to third parties for marketing purposes. If you do not wish your details to be used for this purpose, please write to: Database Manager, Hanson Wade, Suite A, 6 Honduras Street, London ECIY 0TH

