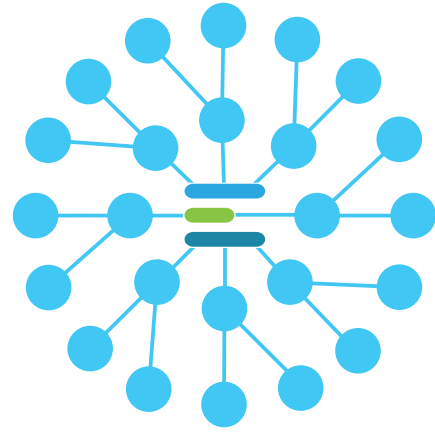


MIMETAS

the organ-on-a-chip company

Organ-on-a-Chip Workshop



In this 2-day workshop, you will learn about 3D cell culture and model development in microfluidic platforms, directly from our experienced scientists. The workshop involves both theory and practice and you will spend a full day in the lab to seed and analyze your cultures. One expert trainer per 3 participants makes this an intensive learning experience. By attending this workshop, you will learn how to design and carry out Organ-on-a-Chip experiments using the OrganoPlate®, and introduce the technology in your lab.

After the training, you will receive 2 OrganoPlates® and protocols to get you started in your own lab.

Program

Hands-on laboratory training in:

- Handling of cells and extracellular matrix gels for microfluidic devices
- Establishing 3D cultures in the OrganoPlate® (in-gel and perfused tubules)
- Fixation, staining and imaging of 3D cultures
- Assays and data analysis

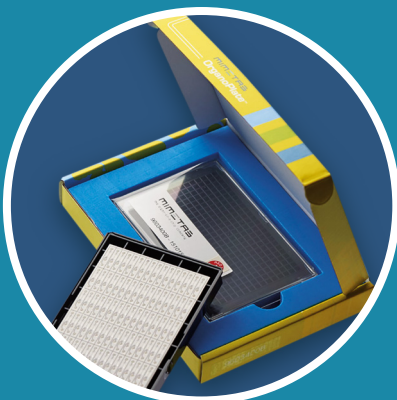
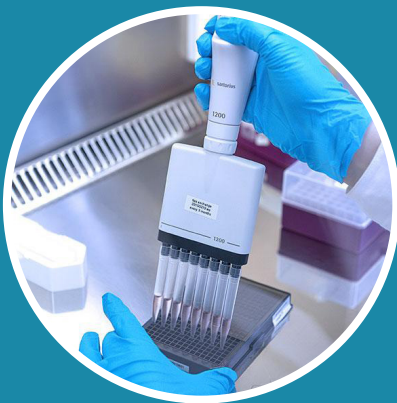
In-depth lectures on the following topics:

- Introduction to Organ-on-a-Chip and the OrganoPlate®
- Key parameters for advanced cell culture models
- Assays and (high-content) imaging of 3D cultures
- OrganoPlate® platform and applications

Materials included in the workshop

2 new OrganoPlates® to take home and start immediately in your lab

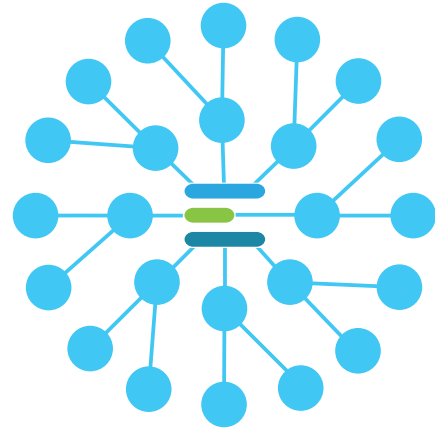
- Seminar materials and handouts
- OrganoPlates®, cells and reagents
- Use of equipment, such as the Molecular Devices ImageXpress® Confocal high-content imaging system
- Lunch on both training days
- A workshop certificate



MIMETAS

the organ-on-a-chip company

Organ-on-a-Chip Workshop



Registration information

- Info:** The workshop is co-organized with Molecular Devices Japan
- Dates:**
- May 21 - 22, 2019
 - May 23 - 24, 2019
 - August 27 - 28, 2019
 - August 29 - 30, 2019
- Location:** Molecular Devices Japan
7th Floor, Ichigo Nihonbashi East Bldg. 2-7-8
Nihonbashi Bakuro-cho, Chuo-ku,
Tokyo 103-0002, Japan
www.MolecularDevices.co.jp
www.MolecularDevices.com
- Register:** Please register per e-mail via info-jp@mimetas.com

MIMETAS – the Organ-on-a-Chip Company

Enabling you to study complex 3D tissue biology in a simple device, that's our goal. With perfused vessels, co-culture and optimized microenvironments. So easy to use that you forget you're working with a highly advanced 3D culture platform. With the OrganoPlate®, we believe we've reached our goal. Say hello to the future of 3D tissue models.