

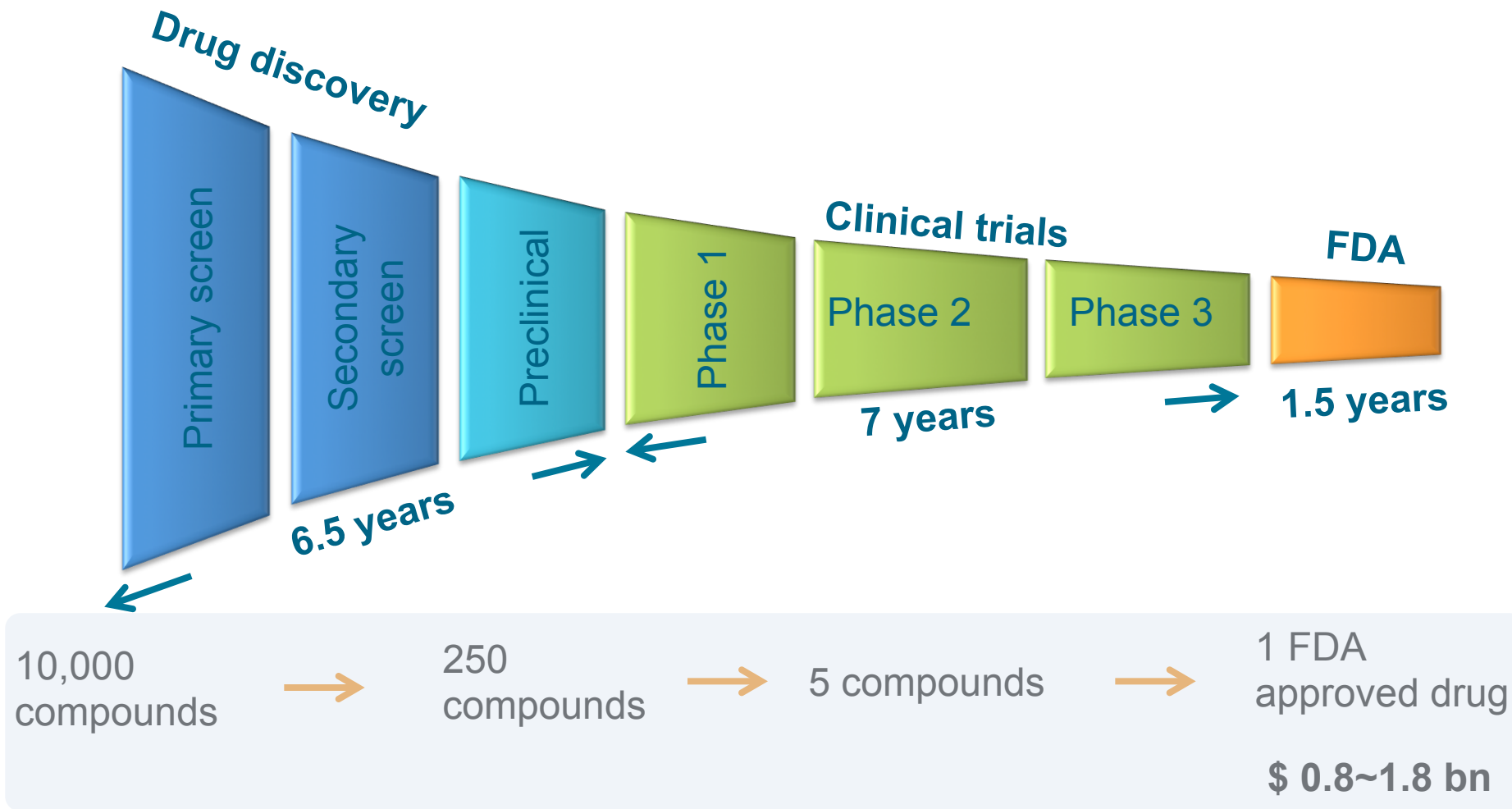
MIMETAS

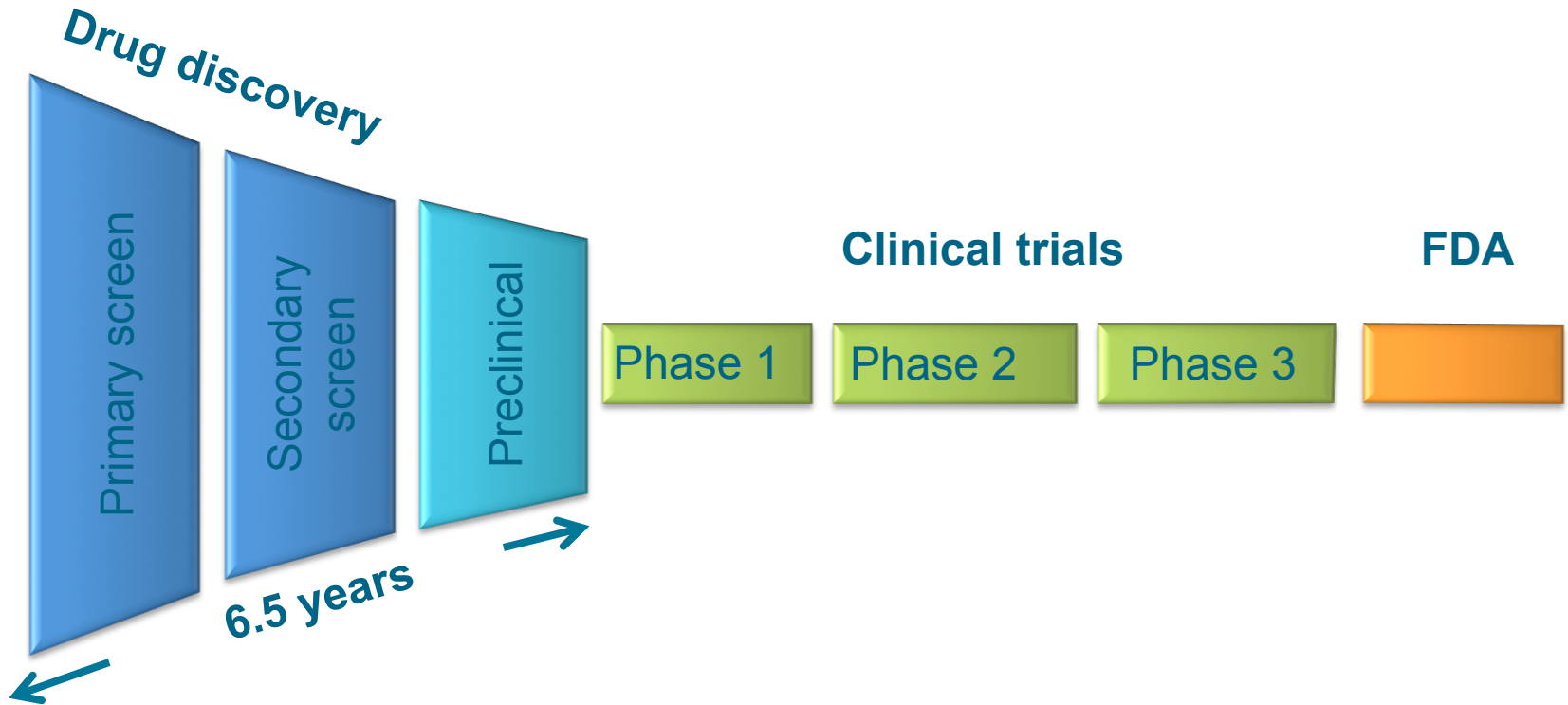
the organ-on-a-chip company

Karlijn Wilschut

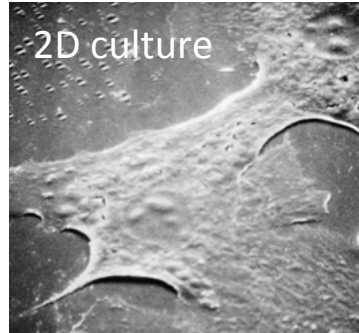
Promega Discover Glo

March 10, 2015- AMC Amsterdam

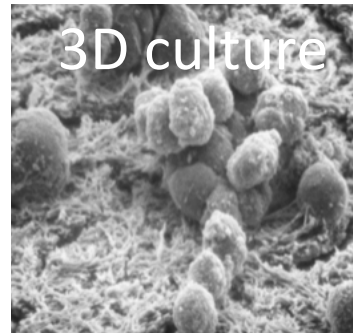
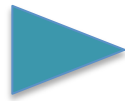
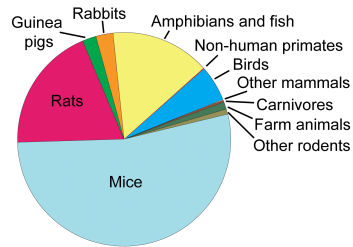




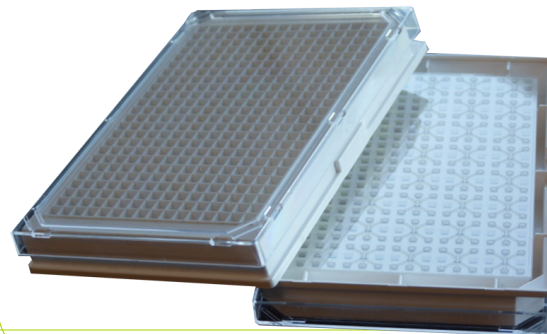
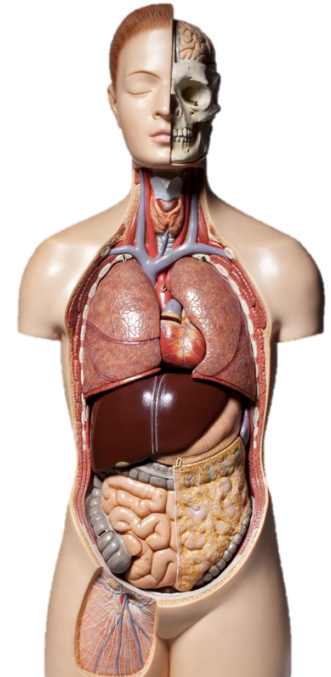
In need of better predictive models

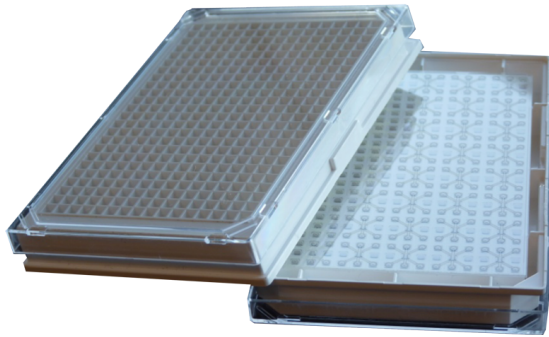


Animal testing



Human tissues
Layered co-culture
Tube formation
Boundary tissues

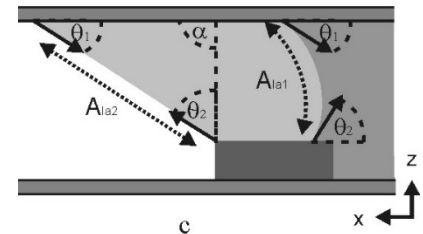
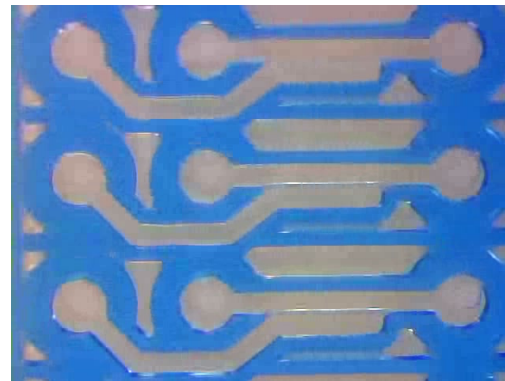




- Organ and tissue functionality
- Perfusion
- High throughput
- Low reagent consumption
- Compatible format

PhaseGuides

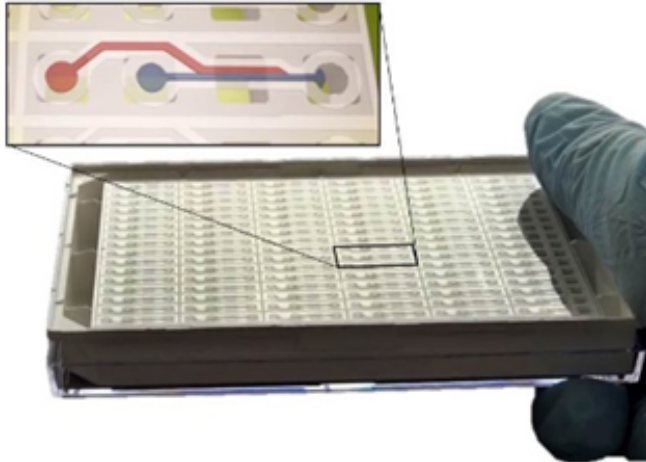
- Fully passive liquid control
- Easy filling of any shape
- Liquid & gel patterning
- Gradient formation
- Tissue layering



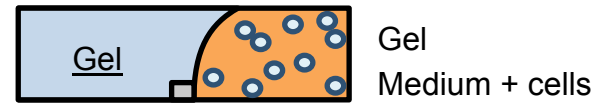
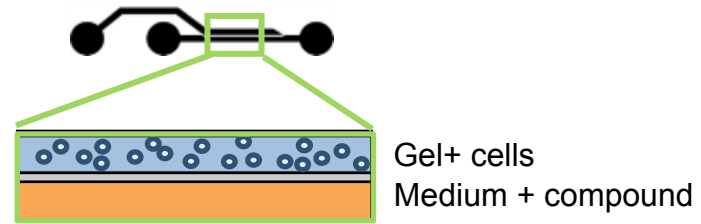


Movie: http://youtu.be/L_VEJAZ5J6U

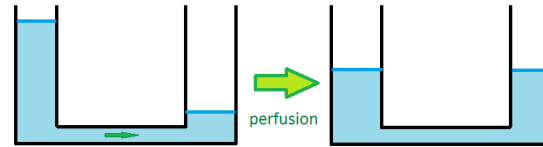




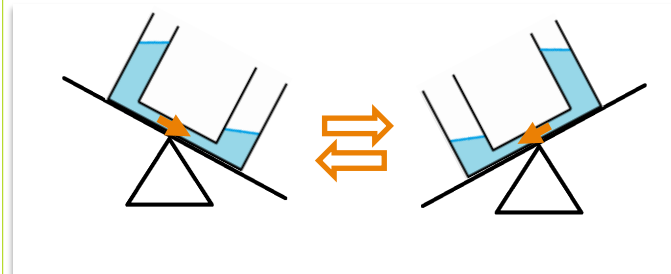
MIMETAS OrganoPlate™ 96x
2-lane chip: gel lane (red) and
perfusion lane (blue)



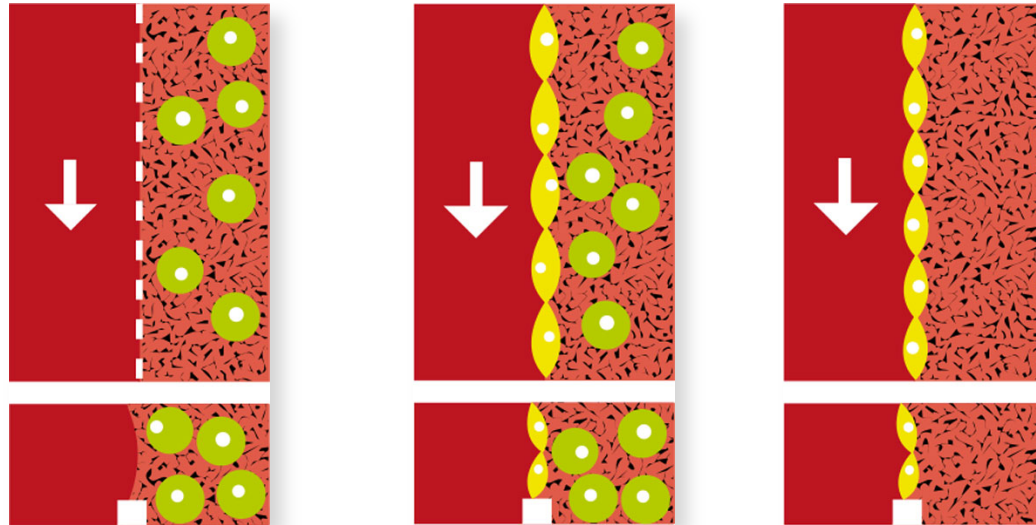
Flow created by perfusion-by-levelling



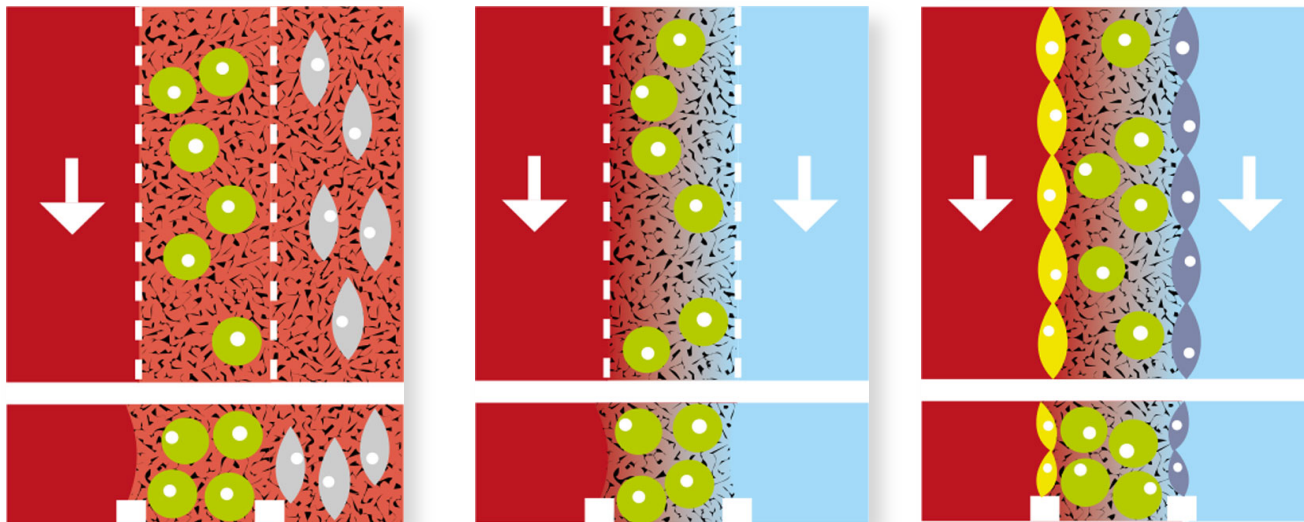
Perfusion-by-levelling on a rocker platform



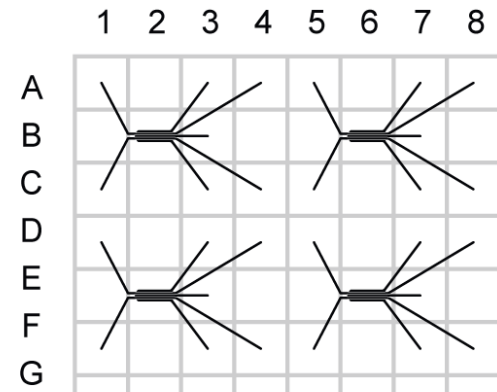
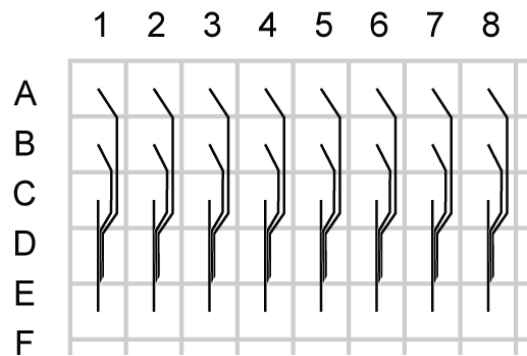
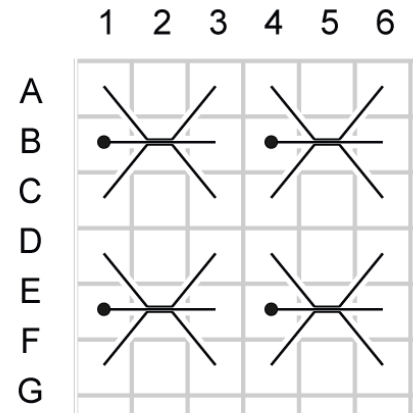
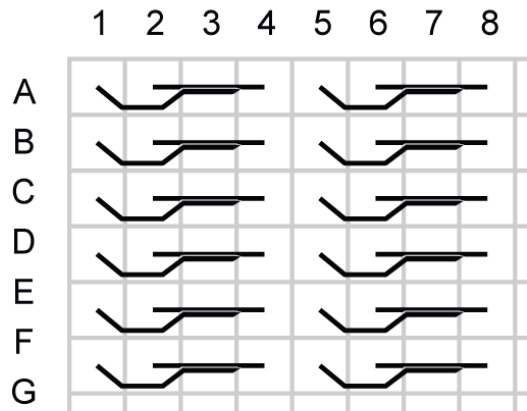
2 - lane

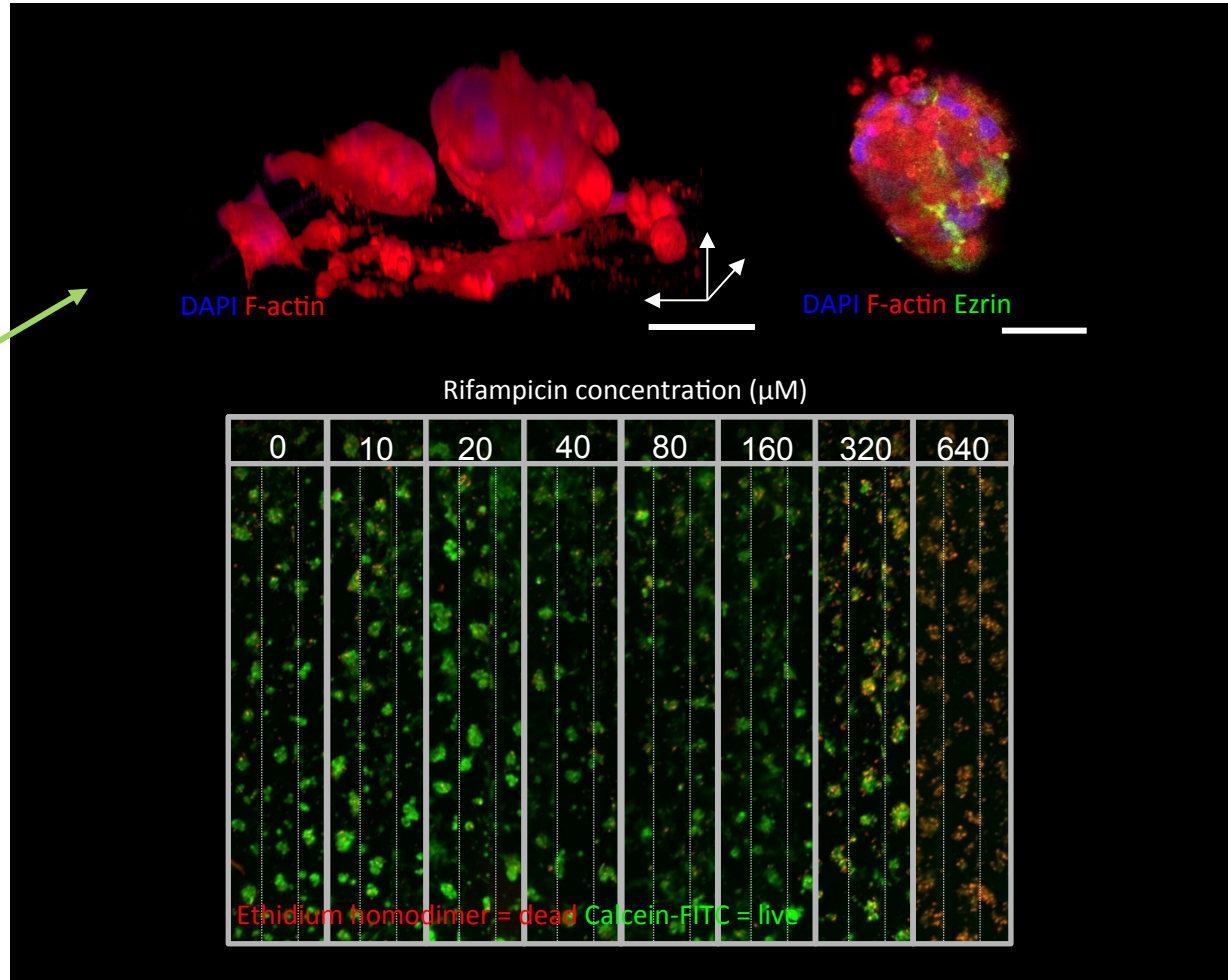
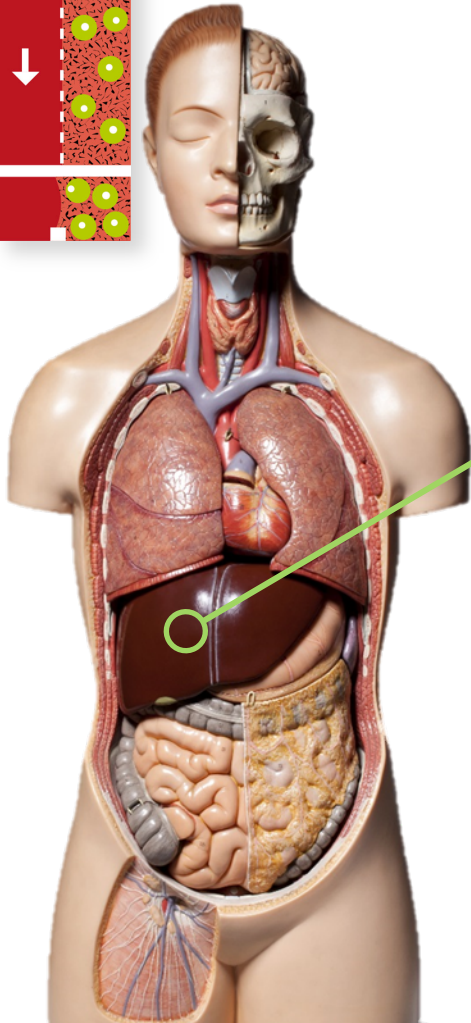
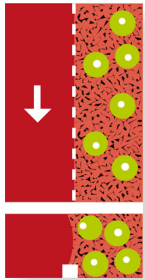


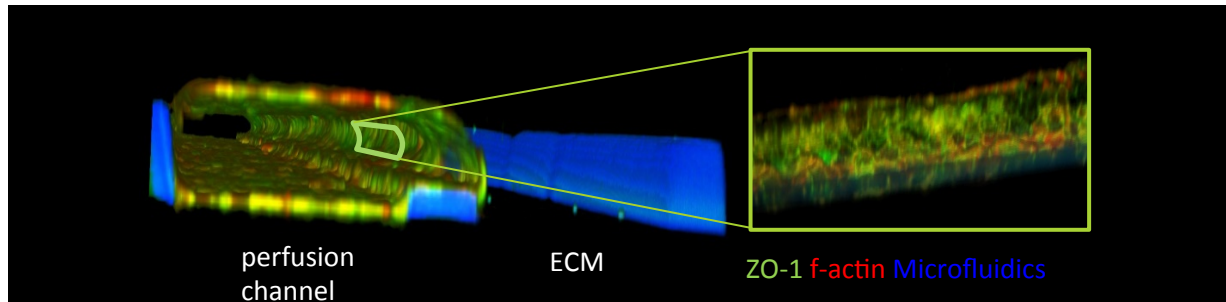
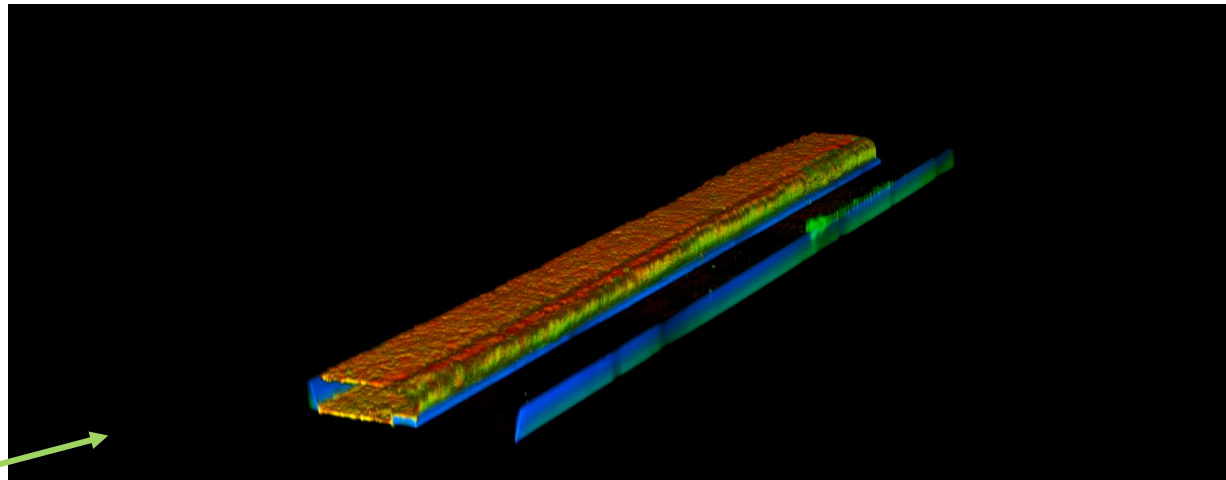
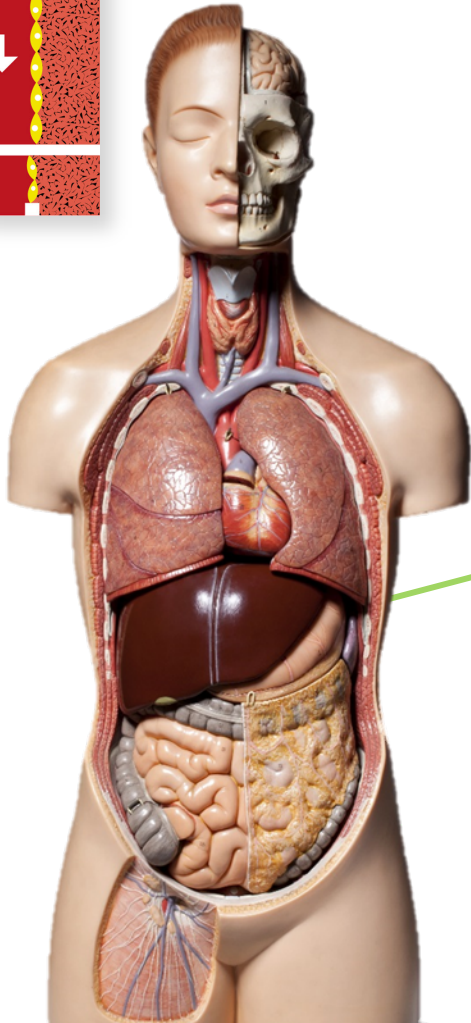
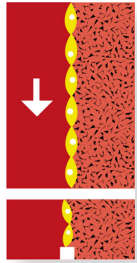
3 - lane



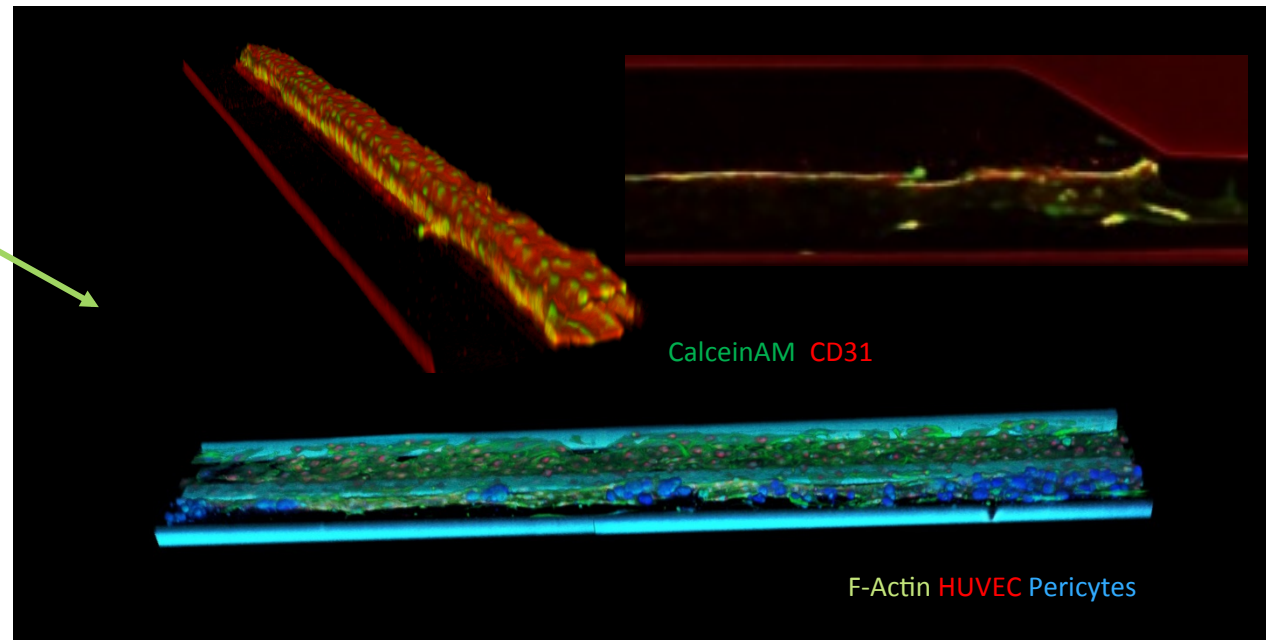
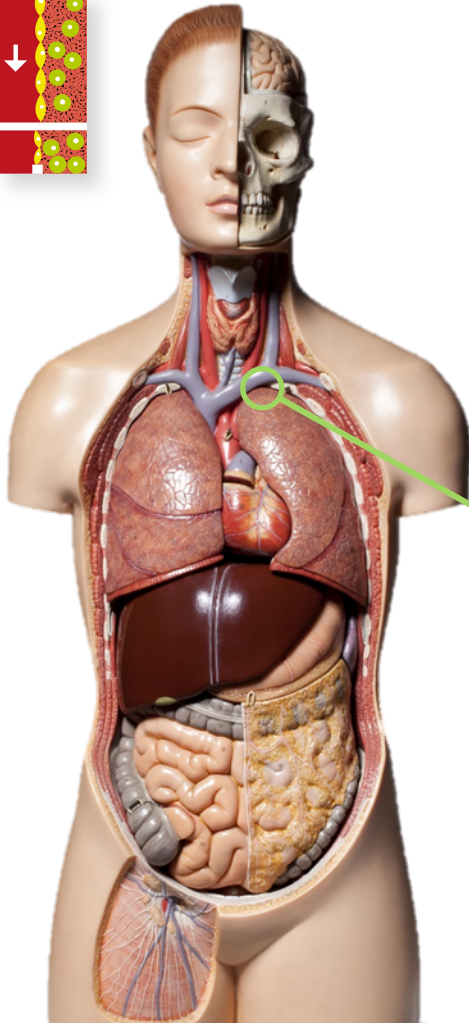
...and 4 and 5-lane setups

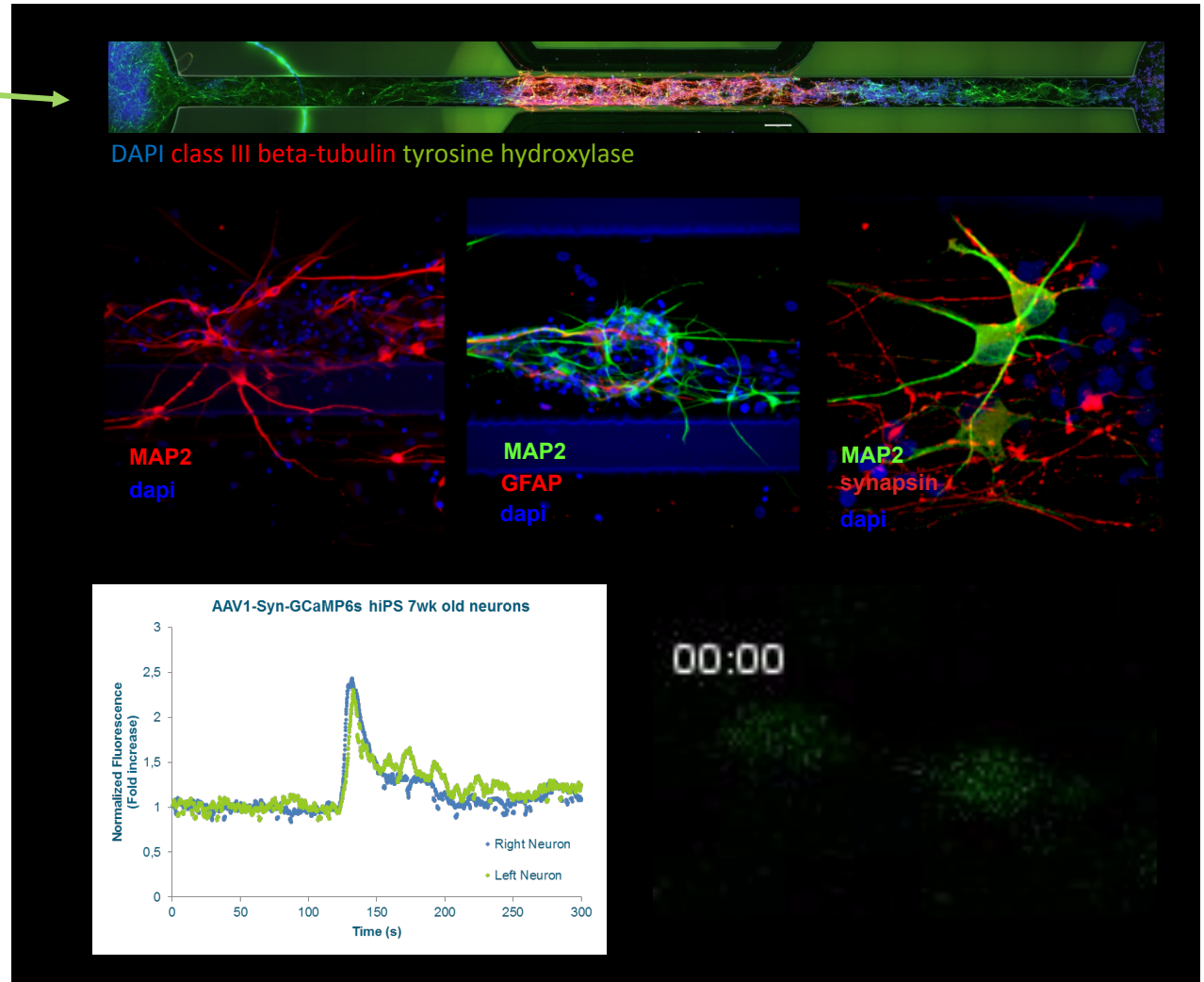
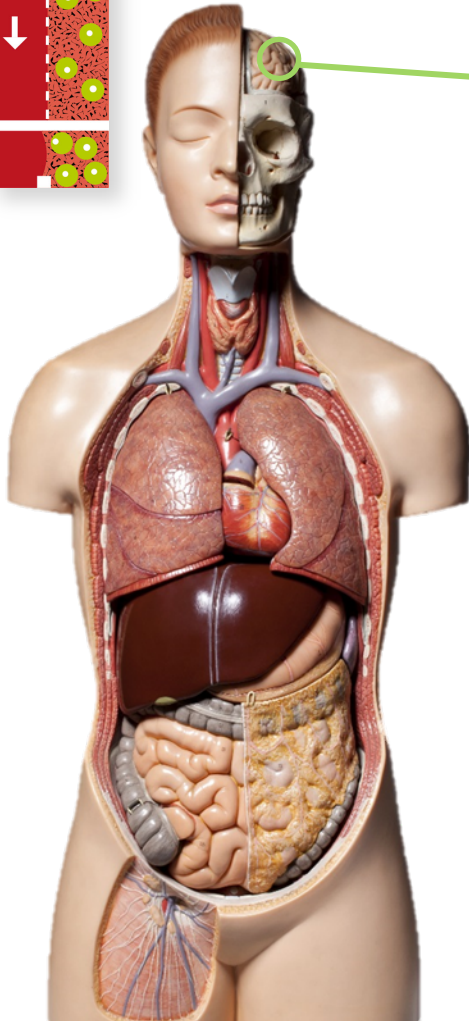
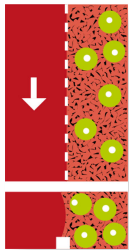


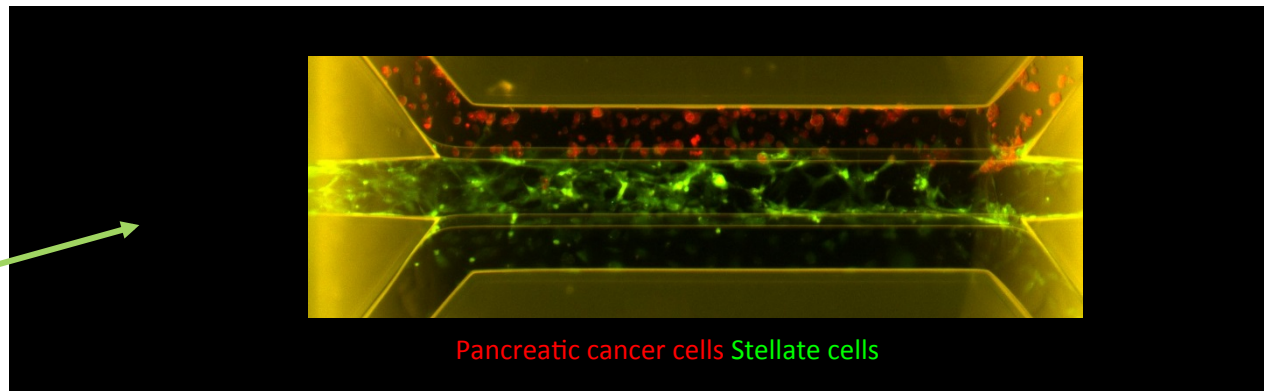
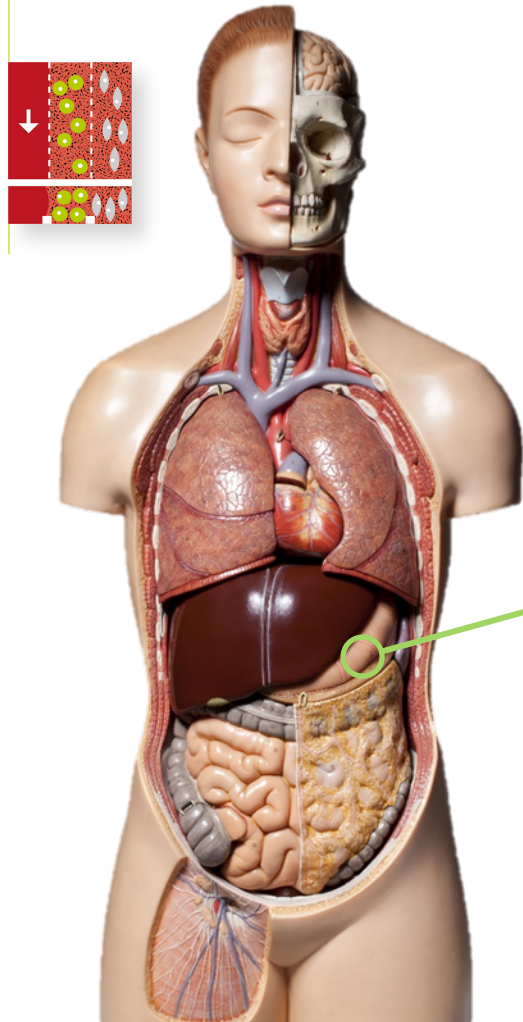


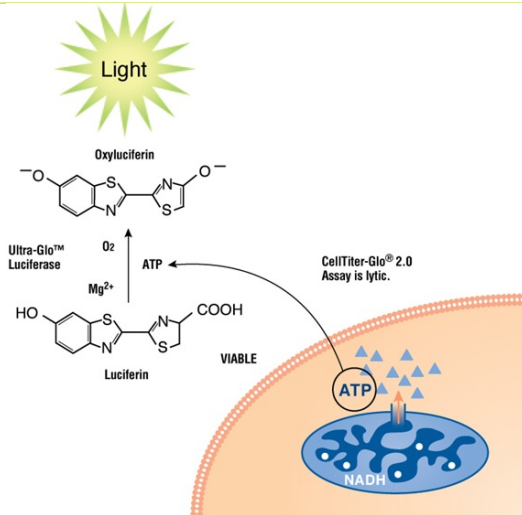


Blood vessels, capillaries

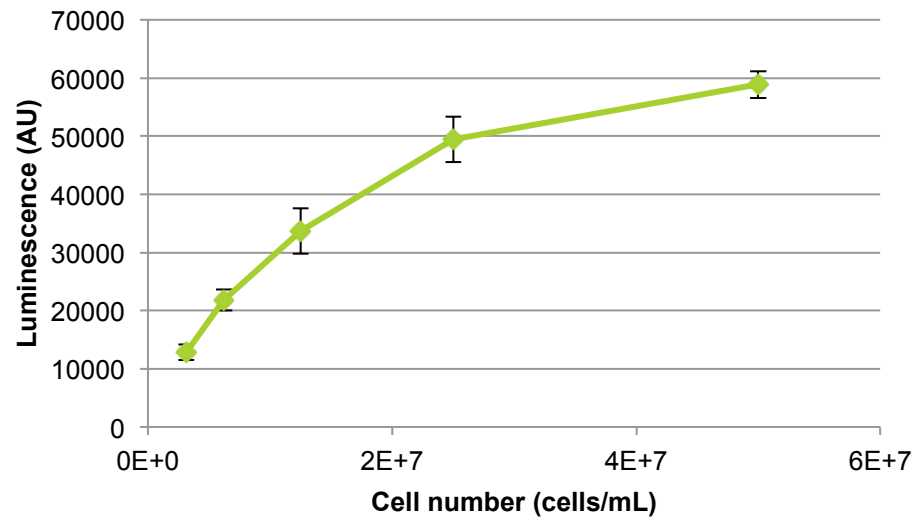
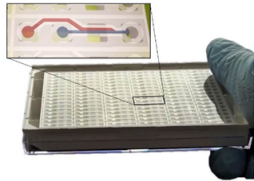






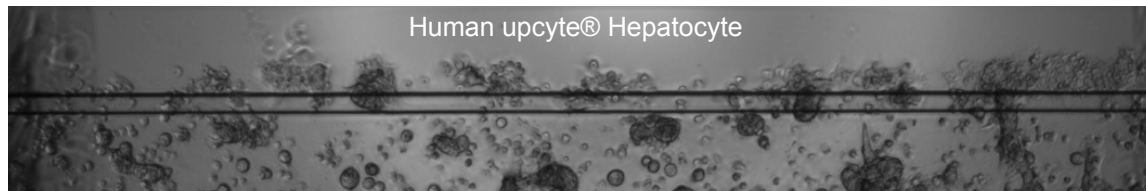


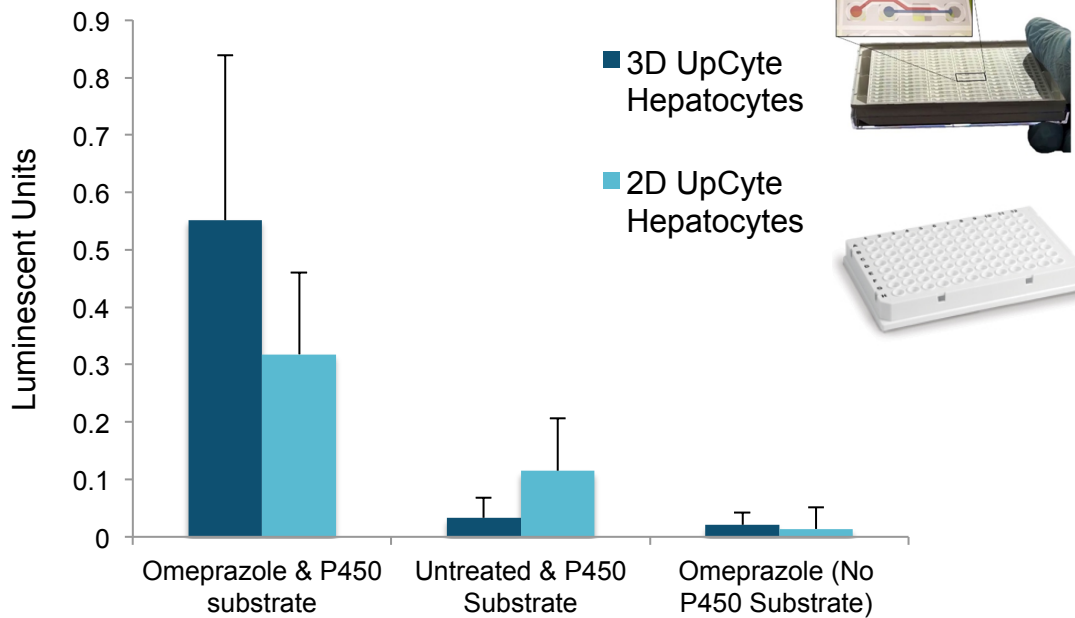
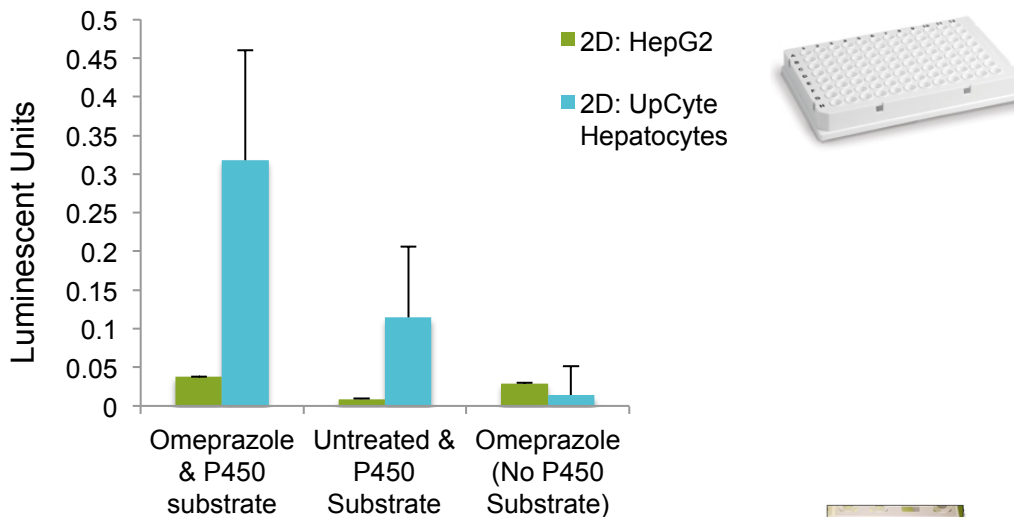
- Luminescent method for measuring for ATP in the OrganPlate™
- Different HepG2 cell densities
- Luminescence microplate reader





- Luminescent method for measuring cytochrome P450 activity (CYP enzyme)
- Treatment with 50mM Omeprazole for 72hrs
- Detection of CYP 1A2 metabolic activity with P450-Glo™ Assay
- Fluoroskan Ascent™ FL Microplate Luminometer (Thermoscientific)





Pump-less perfusion

Co-culture

Boundary formation

384-well format

Glass bottom plate

Design flexibility

Low absorbance

Increased tissue longevity

Stable culture conditions

Complex, biomimetic models

Epithelial function models

Compatible with readers

Suitable for automation

Medium – high throughput

Excellent optical properties

Wide range of models

Compound compatible

▷ Immunofluorescence

▷ Luminescence assays

▷ Viability assays

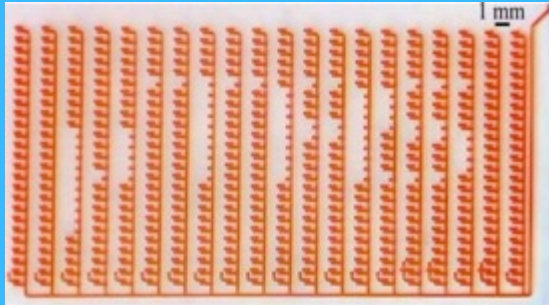
▷ Metabolic assays



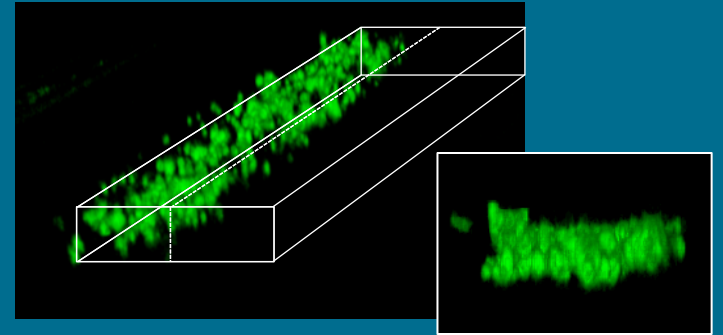
- ◆ Founded in 2011 by Paul Vulto and Jos Joore
- ◆ MIMETAS BV established in 2013 in Leiden, NL
- ◆ Microfabrication in Enschede
- ◆ MIMETAS Inc in Rockville, MD, USA



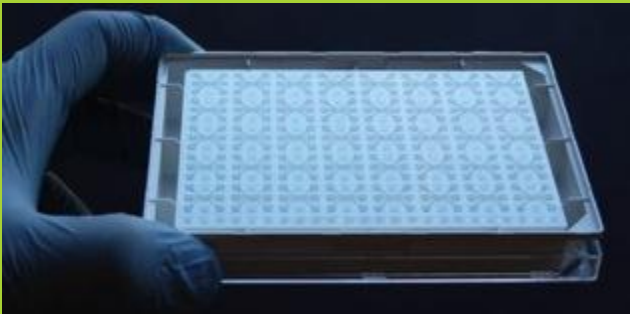
Massive parallel



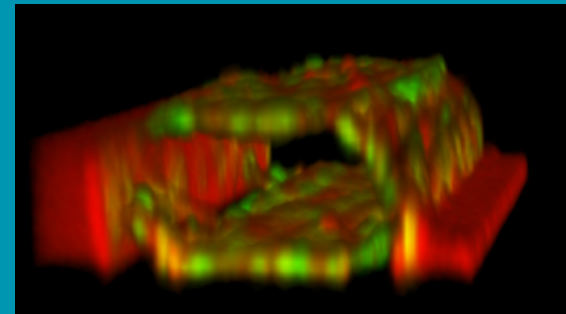
3D cell culture

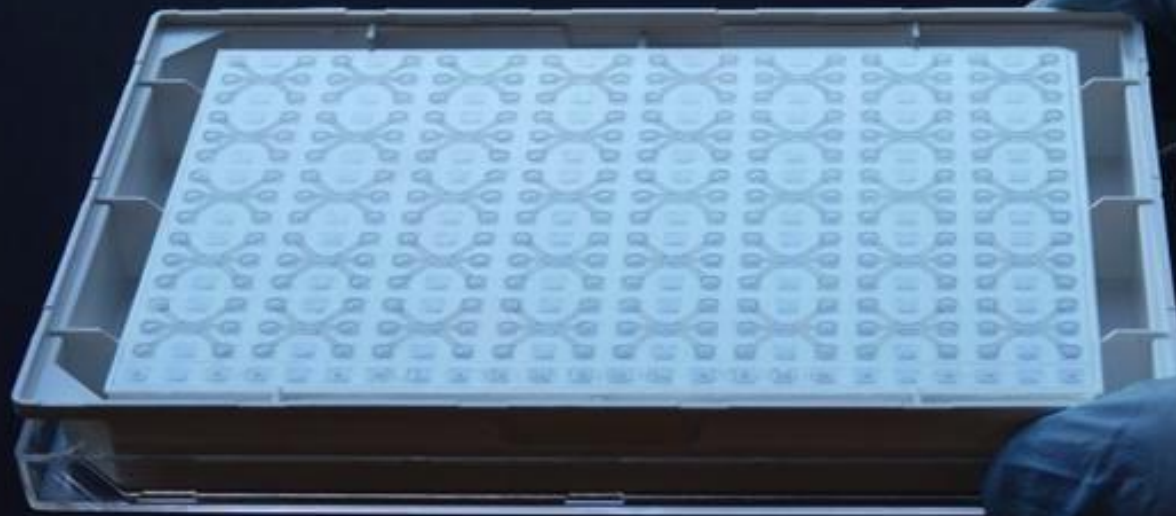


Compatible & affordable



Biomimetic





MIMETAS

the organ-on-a-chip company

k.wilschut@mimetas.com

www.mimetas.com