

MIMETAS

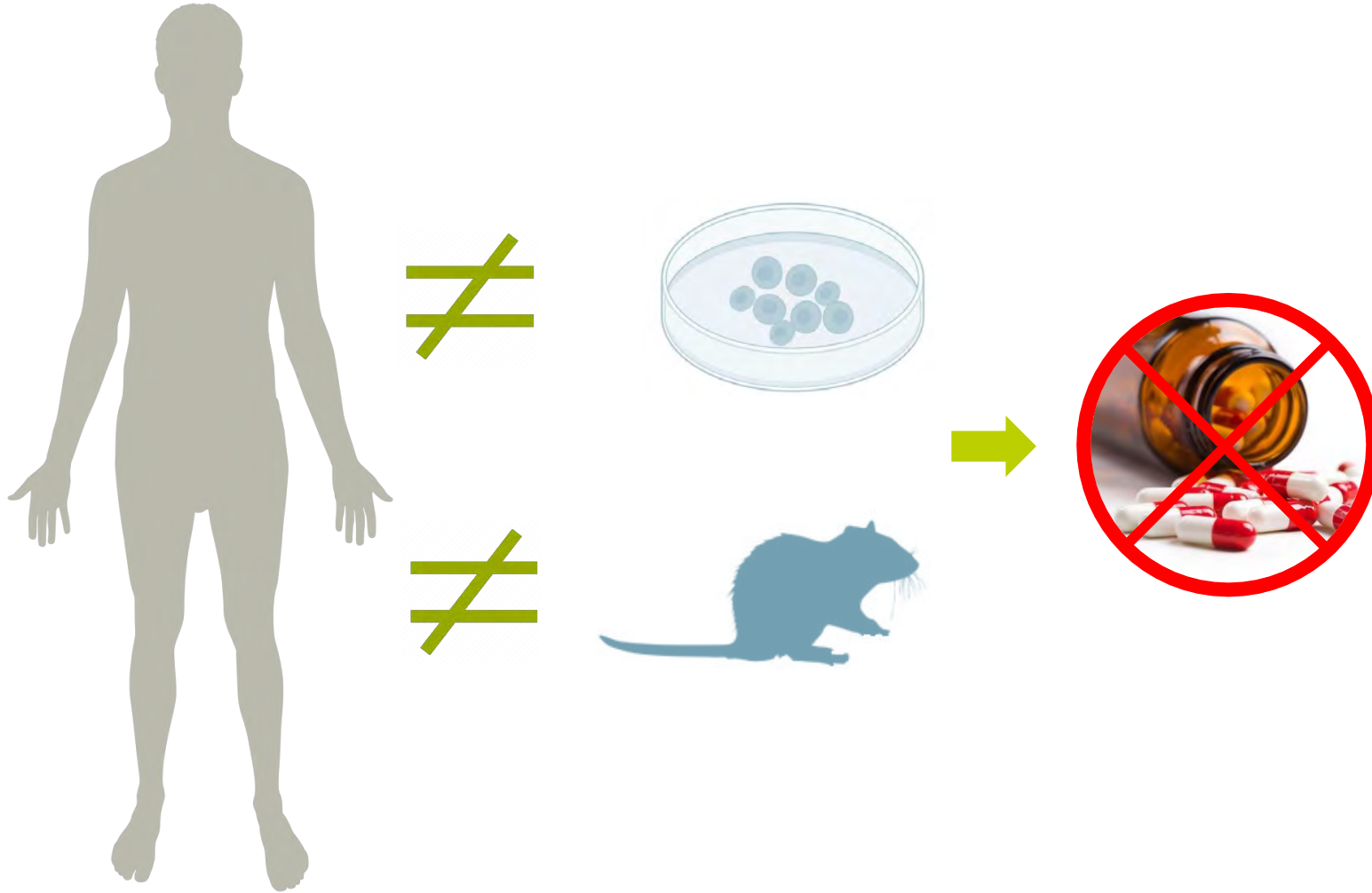
TECAN.

**Coupling organ-on-chip and multimode plate readers: Real-time assessment of snake venom toxicity on a three-dimensional blood vessel model**

Sebastian J. Trietsch

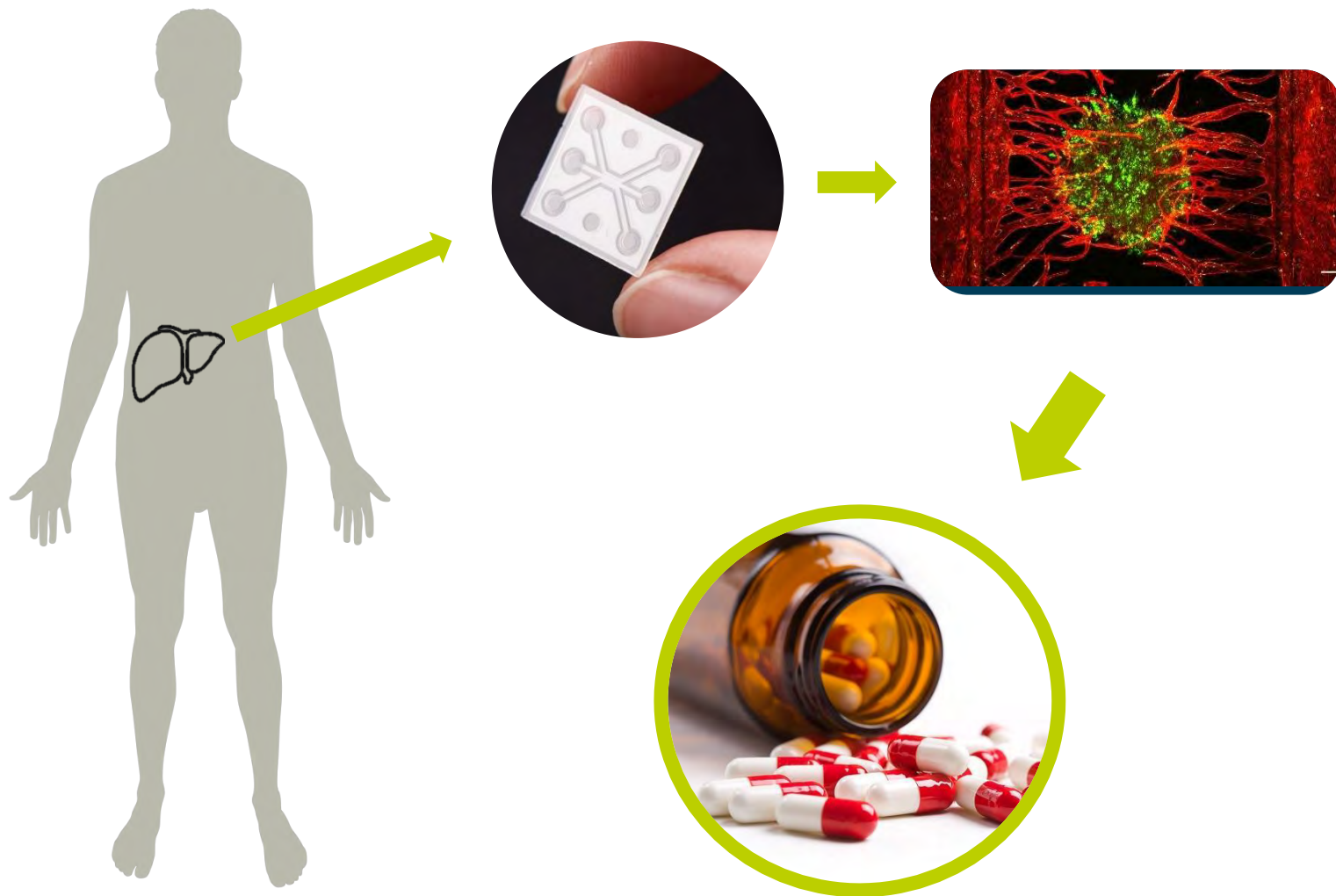
Grow. Learn. Discover.

# ≡ Wrong model means wrong conclusions



**Simple cell cultures & animal models fail to capture the complexity of human disease**

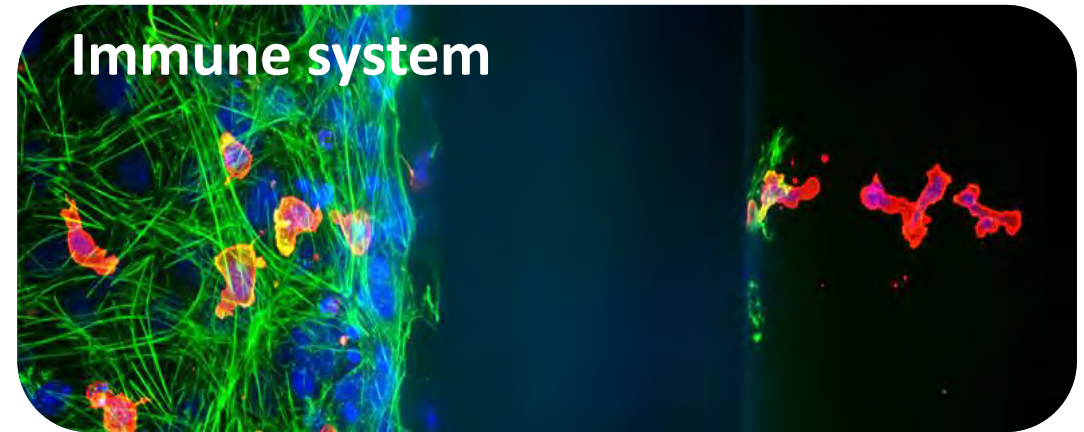
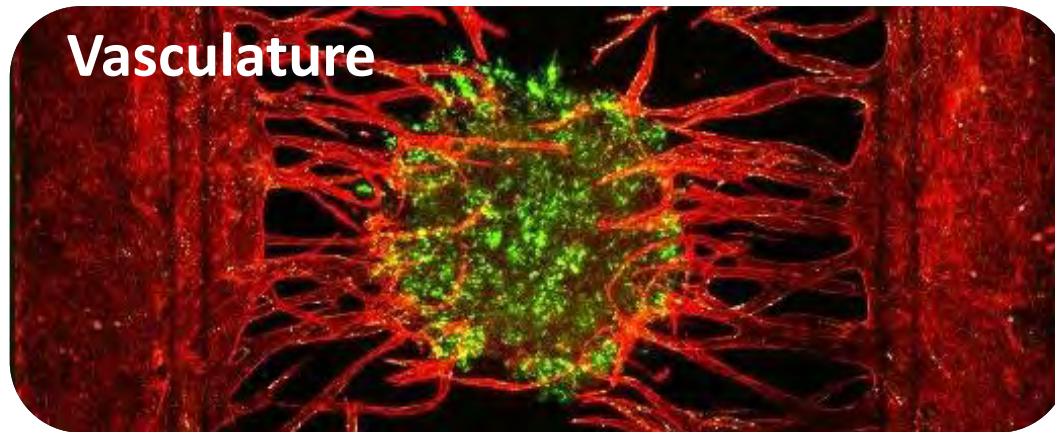
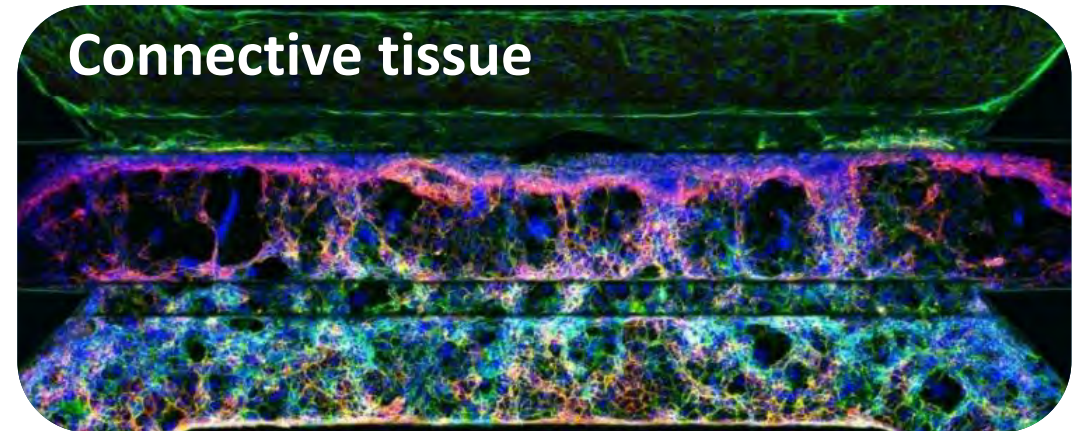
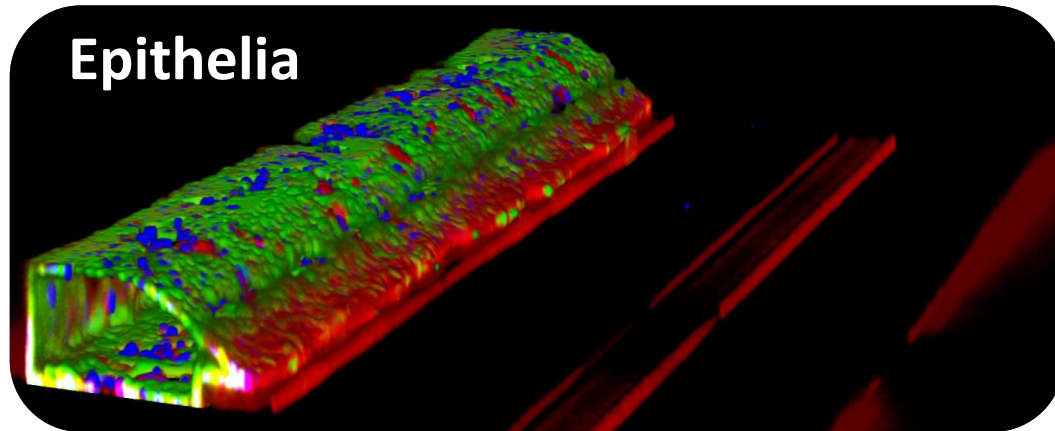
# ≡ MIMETAS provides the right model



**Human Organ-on-a-Chip technology recapitulates the human organ niche and captures the complexity of disease**



≡ The right model looks like this...

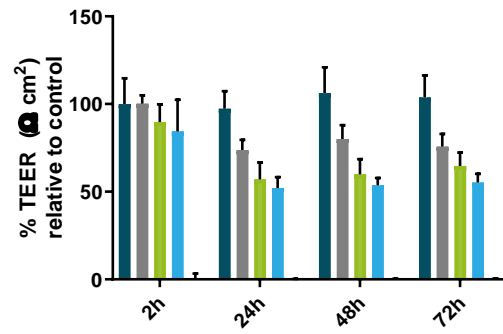


... yields robust and quantitative data

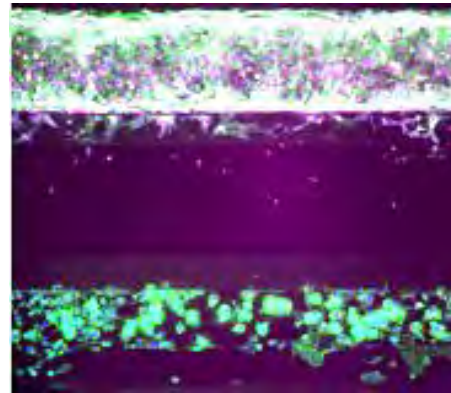
## TEER measurement



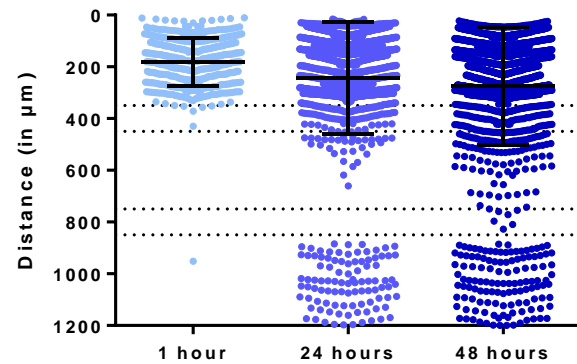
2CC Apical exposure



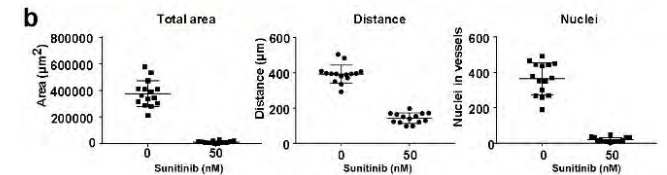
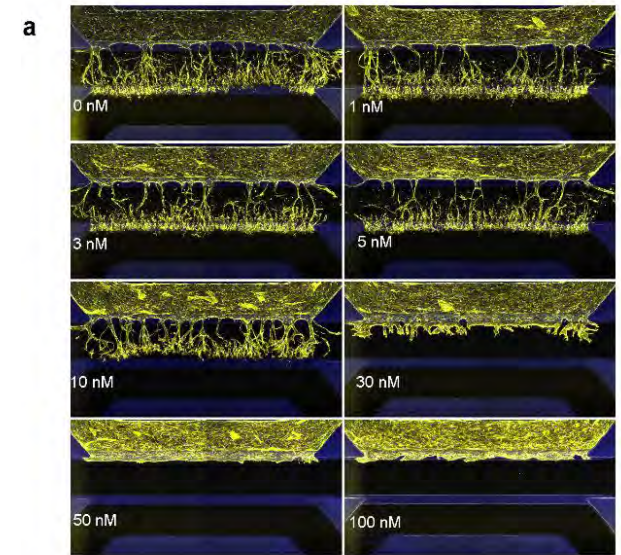
## Immune cell Migration



Actin CD45 E-selectin/CD62E Nuclei

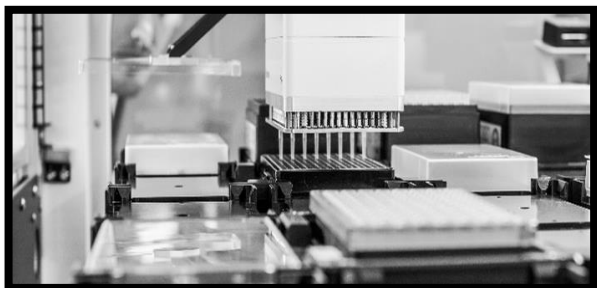


## Angiogenesis





≡ ... and can be repeated thousands of times



**Automated workflow**



**Screenable platform**



**High content screening systems and multimode readers**



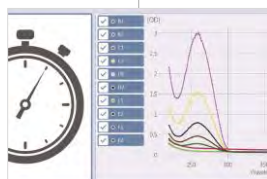
**Thousands of chips**

# Spark<sup>®</sup> Cyto – live cell multimode plate reader with real-time image cytometry

A POWERFUL TOOL FOR MULTIPLEXING



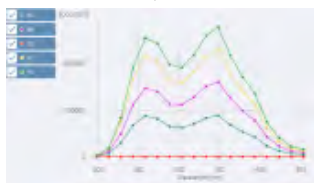
**Absorbance with high-speed monochromators**



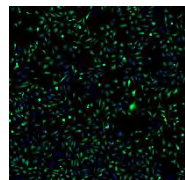
**Fluorescence with FUSION Optics**



**Luminescence with BRET & NanoBRET™**



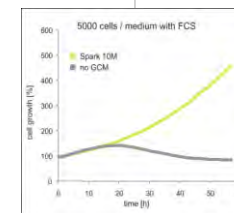
**Brightfield and Fluorescence Imaging**



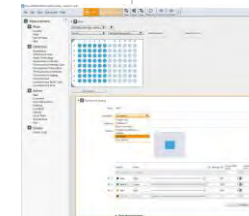
**Automate your workflow with Lid Lifter™**



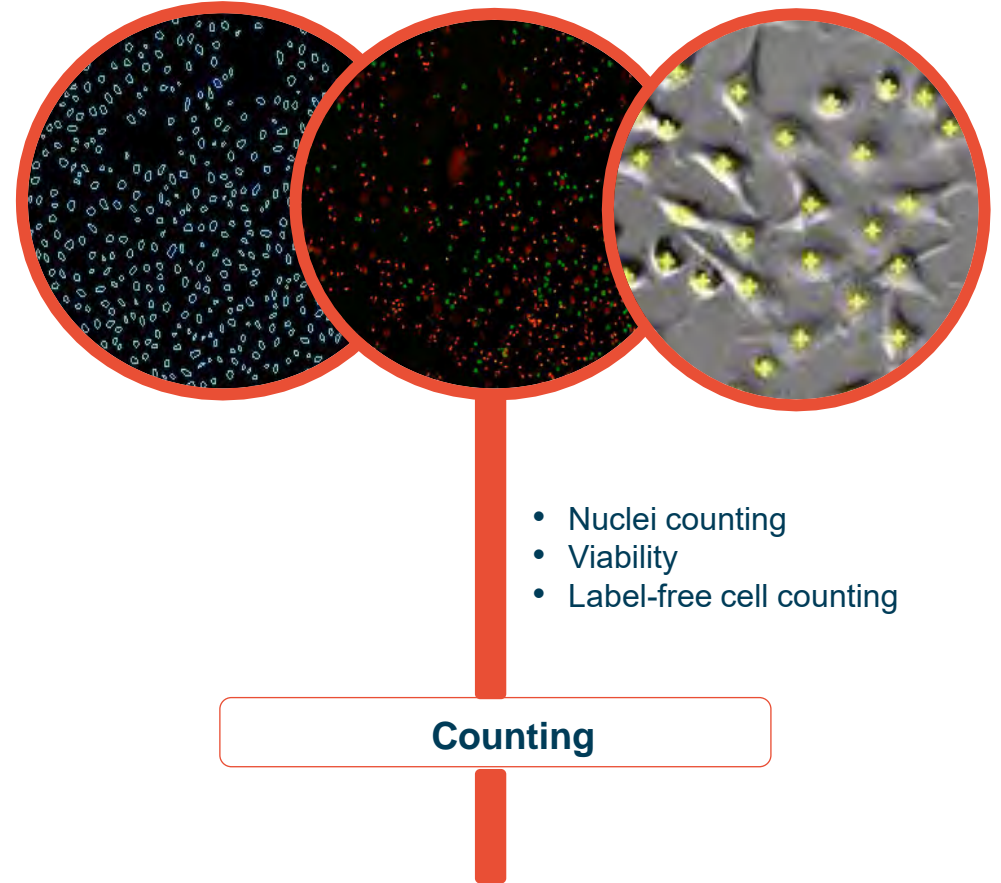
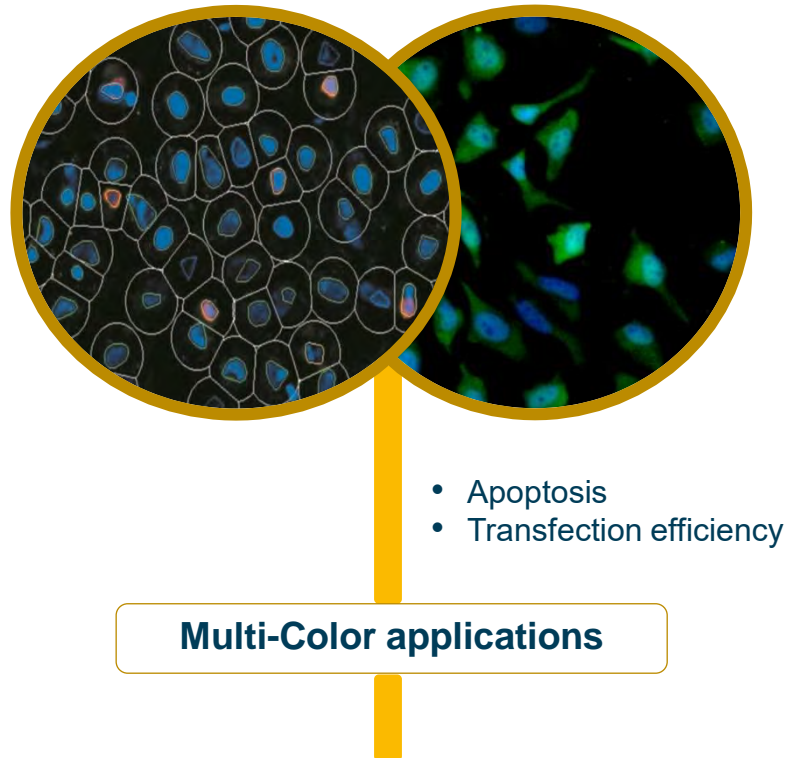
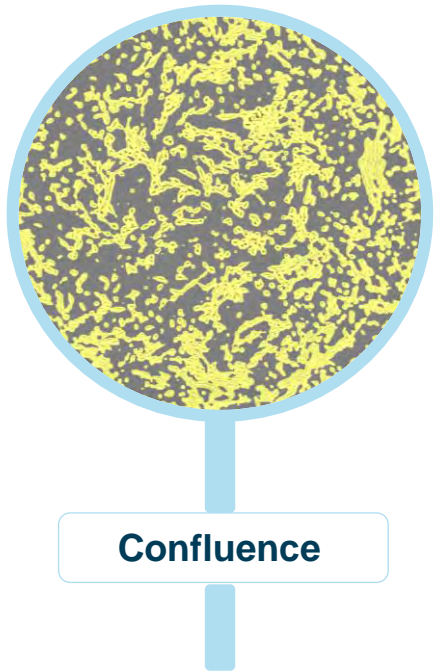
**Cell incubation with temperature/ gas control and humidity cassette**



**Intuitive and user-friendly SparkControl™ software**



# ≡ Spark Cyto imaging - key applications at a glance





# ≡ Synergy between OrganoPlate and Tecan Spark Cyto



**OrganoPlate**



Ease of use

Scalable and throughput capable

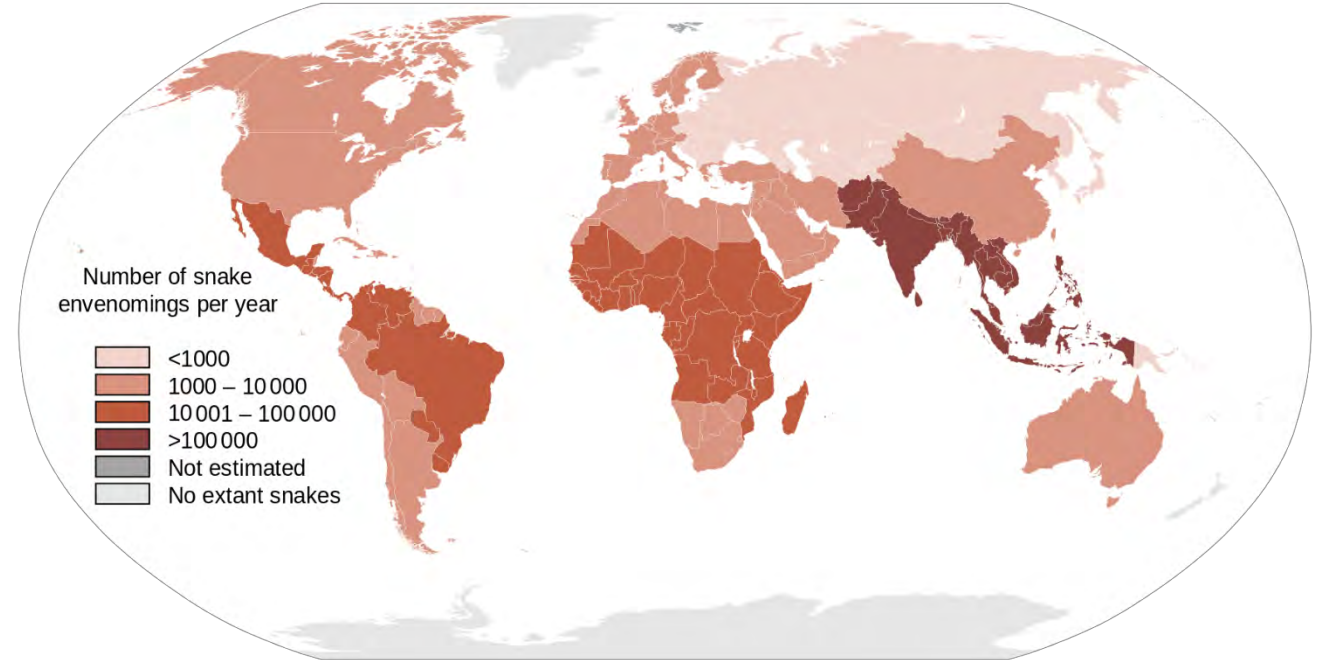
Imaging capable



**Tecan Spark Cyto**

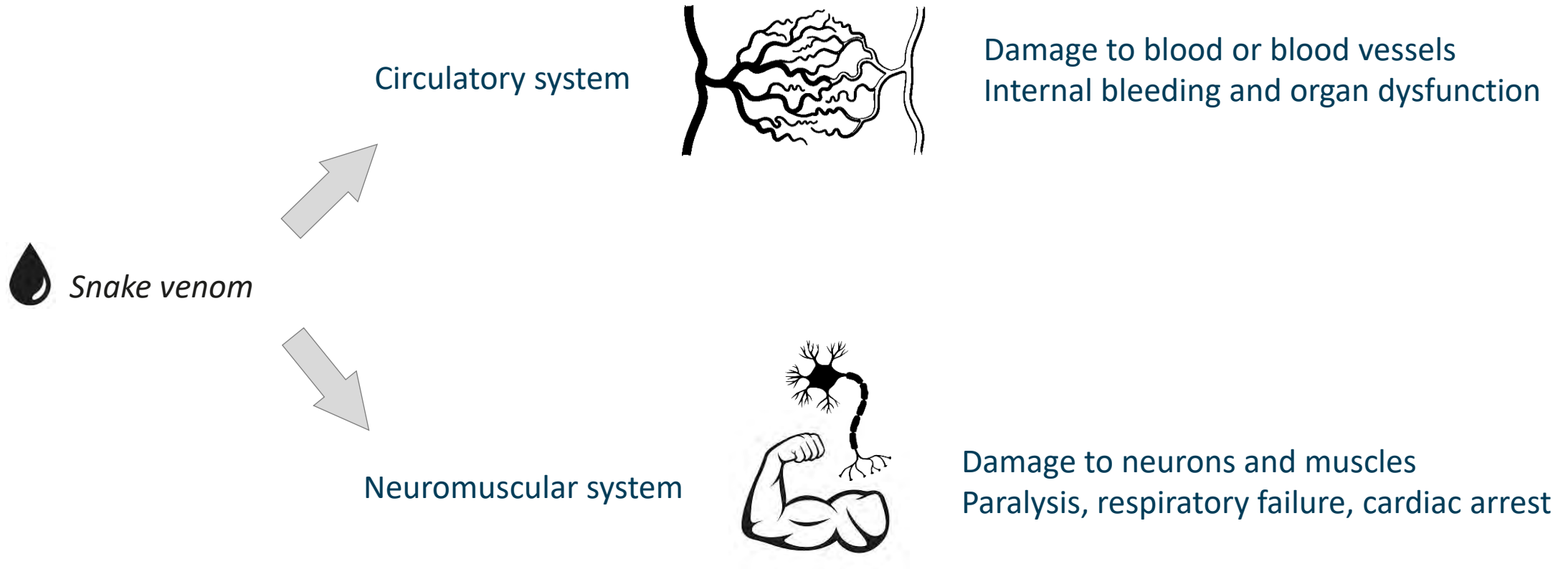


# ☰ Venomous snake bites affects millions



Snakebite seriously injure up to **2.7 million people** and claim some **125,000 lives.**  
(World Health Organization)

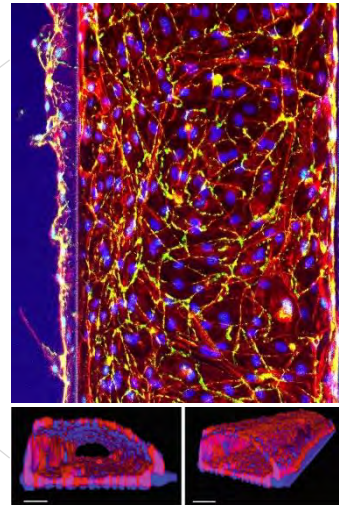
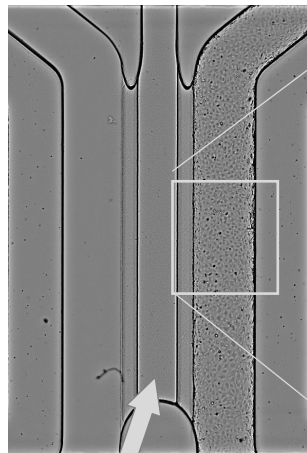
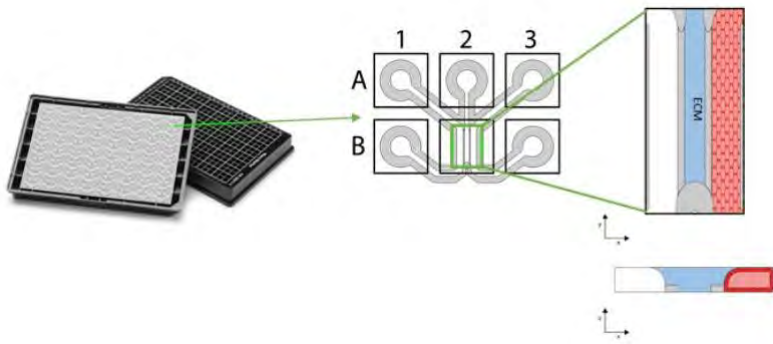
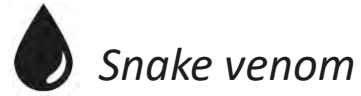
# ☰ Toxicity of snake venoms



*Snake venoms are comprised of a complex cocktail of bioactive substances. They can have multiple effects, including affecting the circulatory or neuromuscular system. The effect of many venoms on the human body is unknown*



# ☰ Venom-induced vascular toxicity model



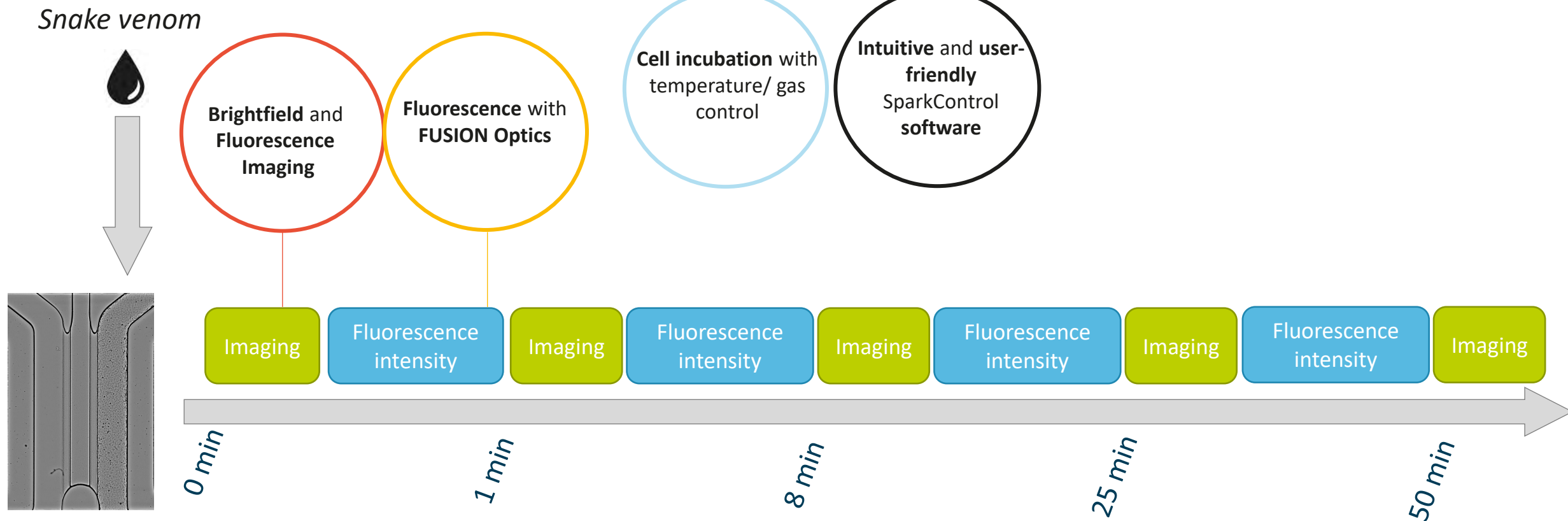
Collagen 1 HUVECs

Vascular leakage  
Morphology  
Viability  
Toxicity

# Integrated multiplex workflow



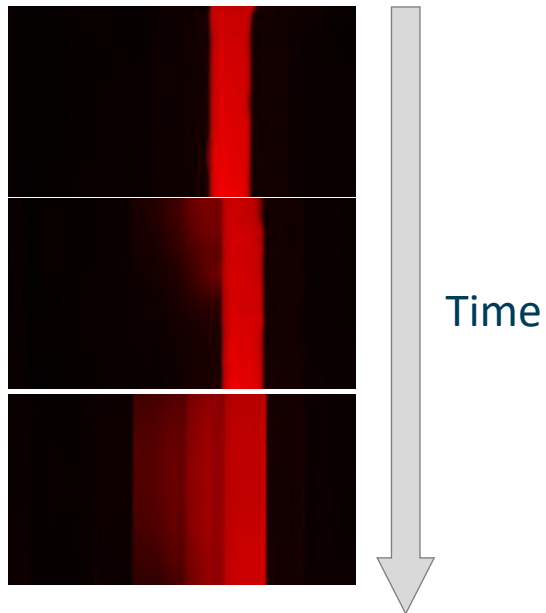
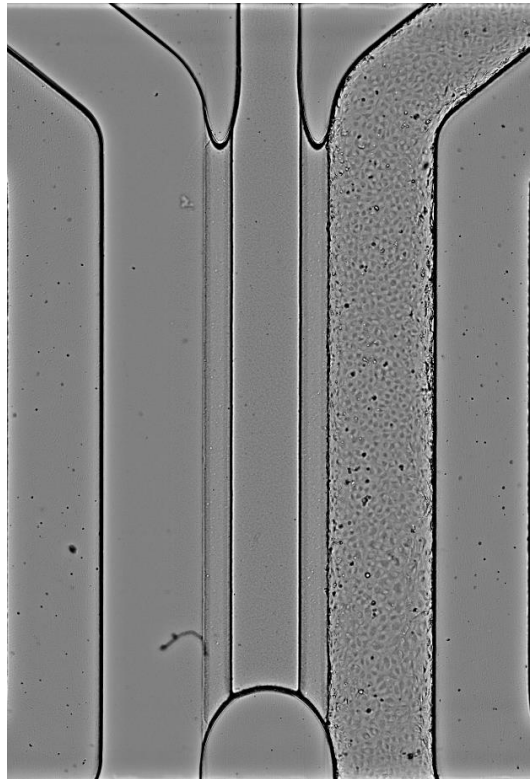
- One integrated workflow protocol
- Multiplexed fluorescence intensity measurements and imaging
- Temperature controlled environment



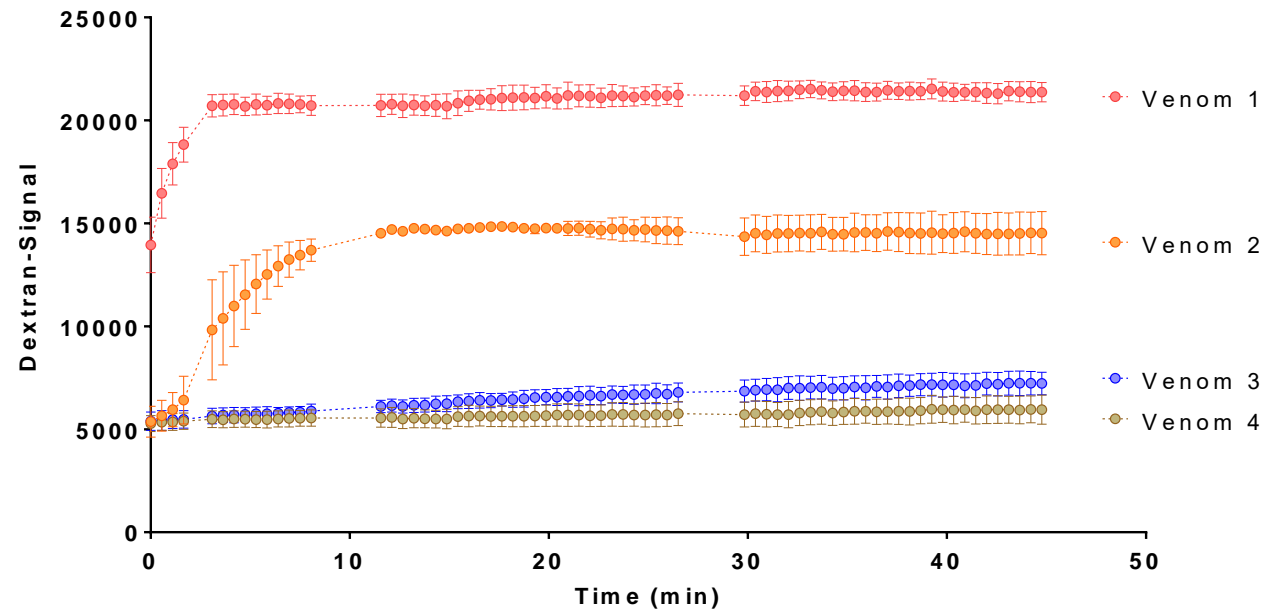
# Vascular leakage assessment

TRITC-Dextran  
Snake venom

Dextran is retained in the blood vessel lumen but permeates in the adjacent channel if the barrier is disrupted



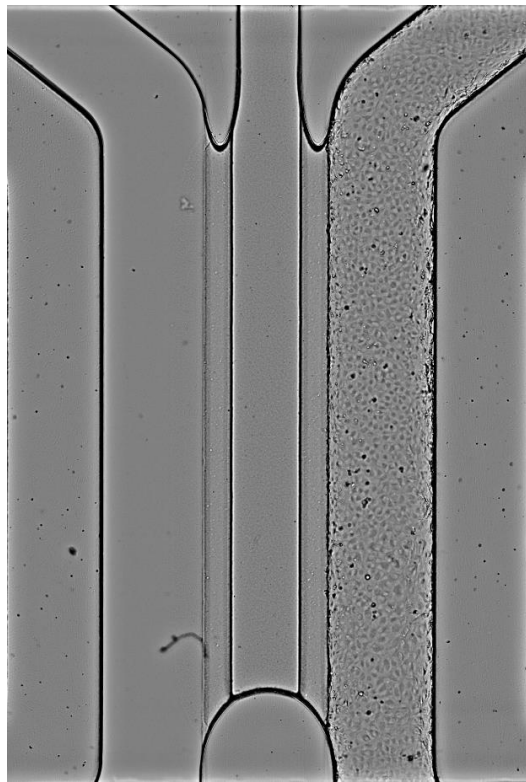
Fluorescence intensity measurement      Vascular leakage



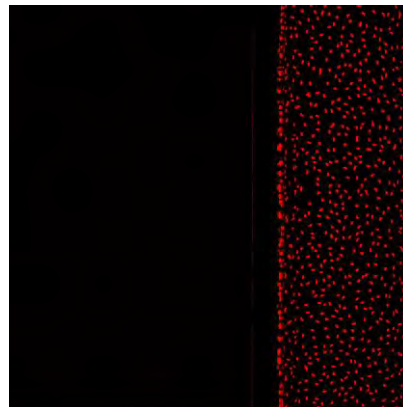
*Venom 1 and Venom 2 appear to induce massive leakage seconds or minutes after addition. Venom 3 shows slight leakage over time while Venom 4 did not induce vascular leakage in this timeframe.*



# Dead nuclei count

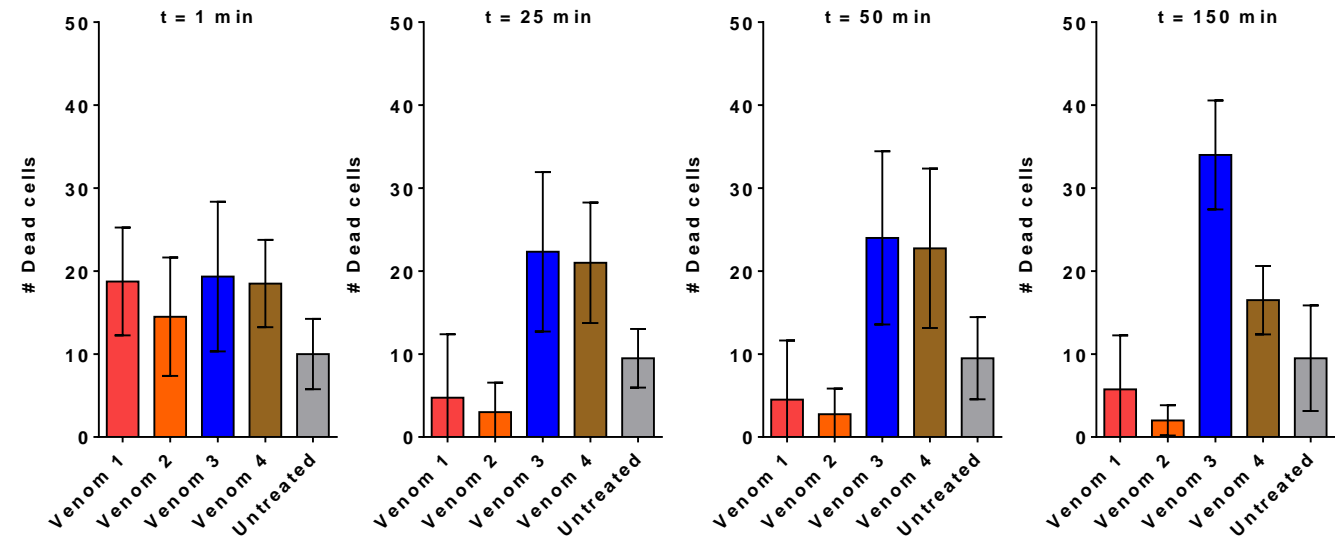


1% Triton control (100% dead cells)



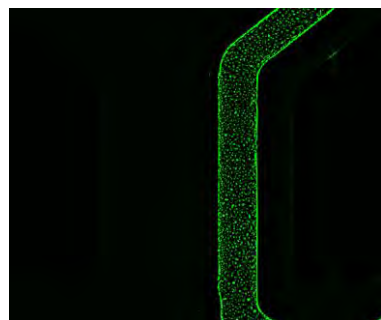
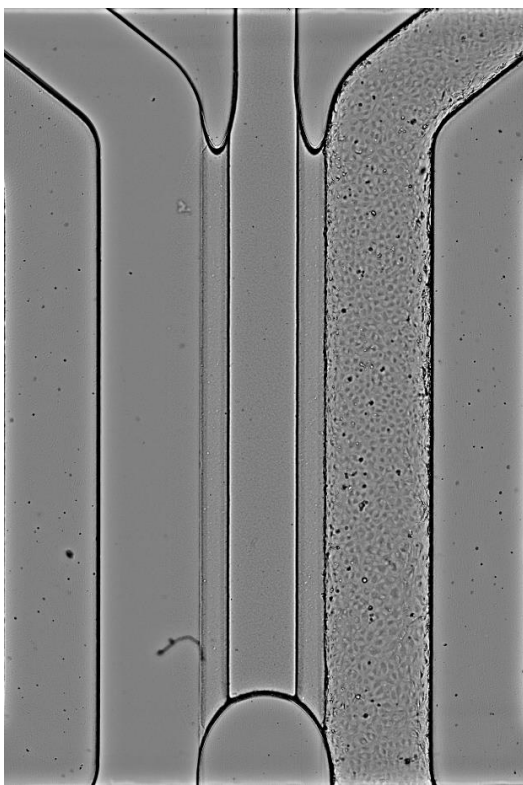
Imaging

Dead nuclei count



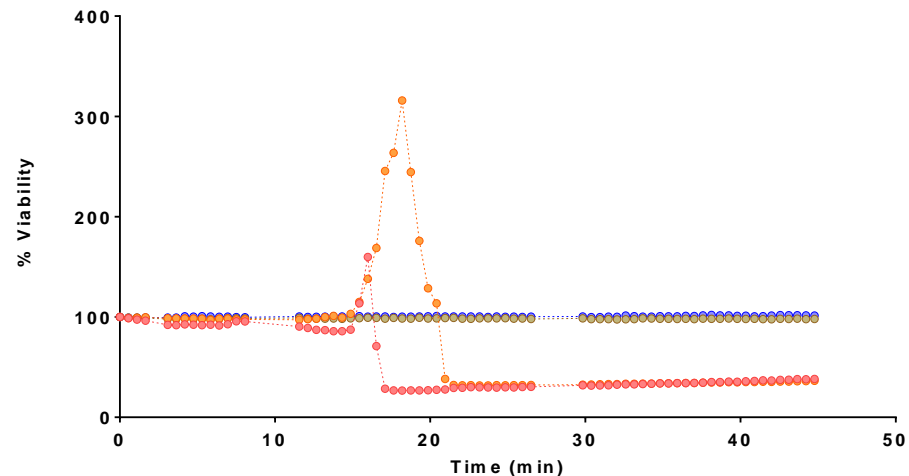
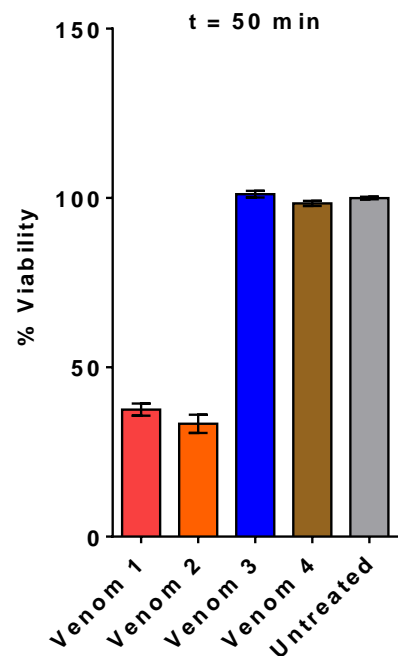
Although **Venom 1** and **Venom 2** induce massive leakage, no increase in dead cells was observed. **Venom 3** and (to a lesser extent) **Venom 4** appear to be causing cell death over time

# Viability assessment



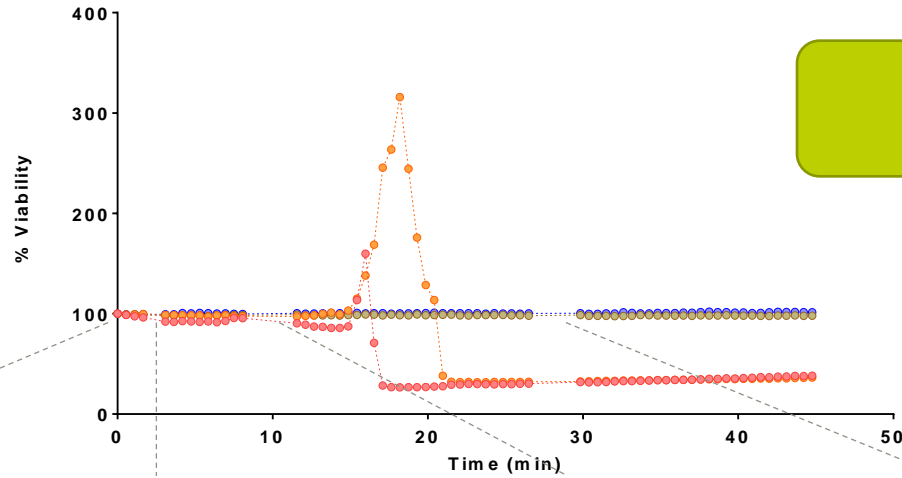
Fluorescence intensity measurement

Viability / morphology



After 50 min incubation, Calcein-AM signal was significantly reduced for **Venom 1** and **Venom 2**. Over-time assessment indicates the emergence of distinct “spikes” followed by a substantial drop in signal. **Venom 3** and **4** do not seem to significantly affect viability.

# Morphological assessment

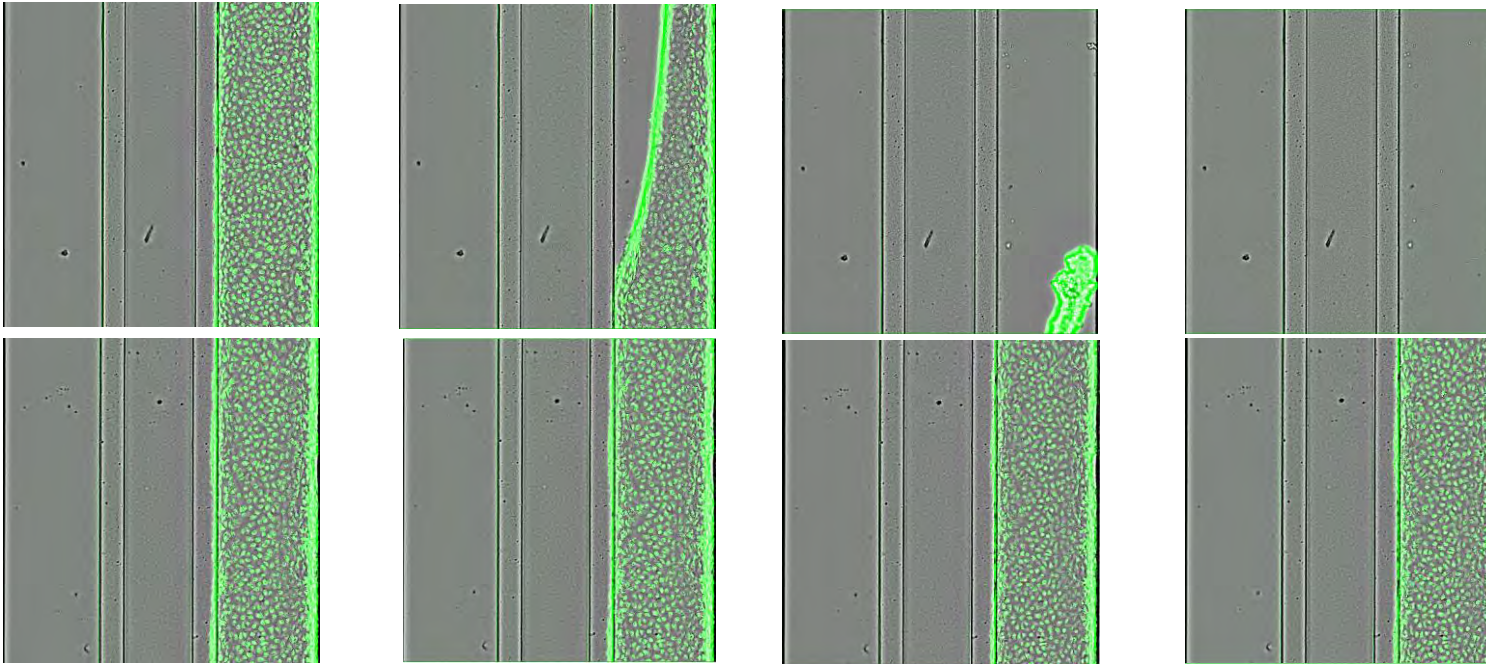


Imaging

Morphology

Venom 2

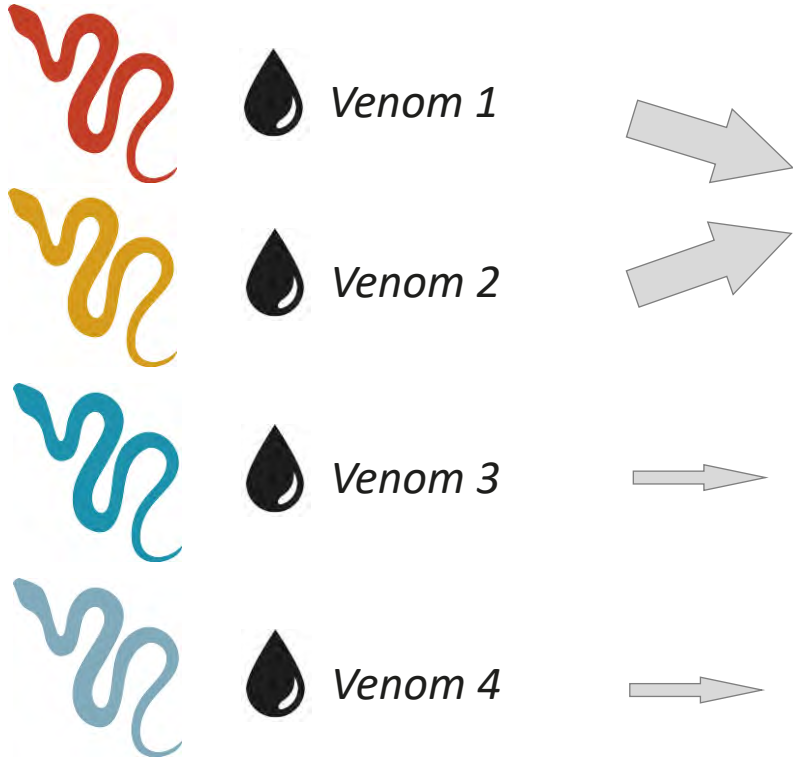
Untreated



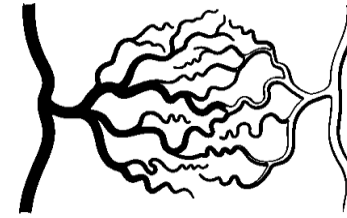
*The action of the venom does not cause the cells to die but rather to contract, hence an increase in signal is measured. Subsequently, the integrity of the whole structure is compromised, the tubule is lost and probably “flushed” out of the channel, leading to the loss of signal.*



# Summary



Circulatory system



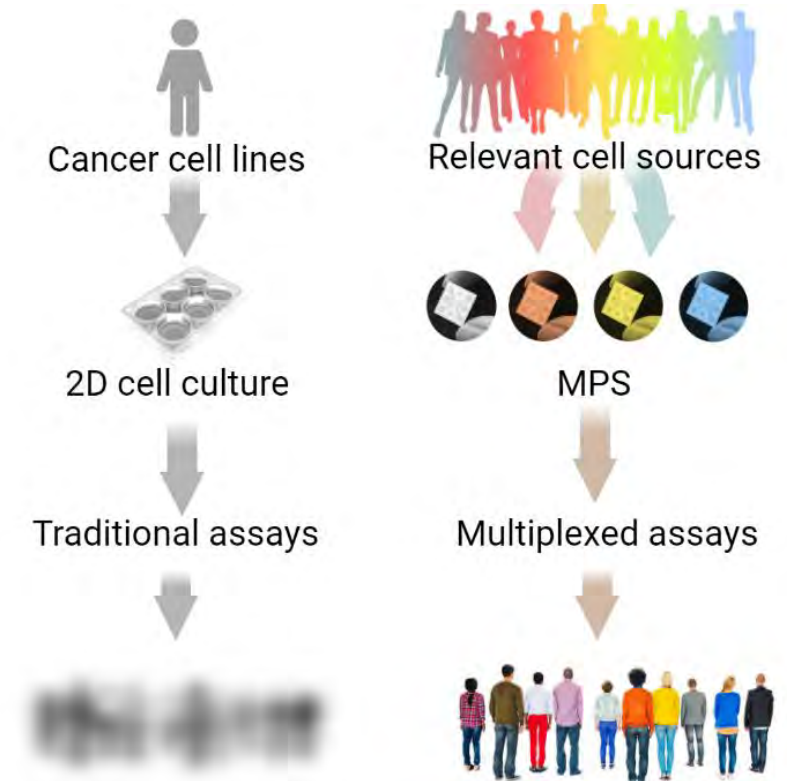
No apparent cell necrosis but massive blood vessel structure damage and high blood vessel barrier disruption

Indication of endothelial cell necrosis and elevated blood vessel leakage

Slight indication of endothelial cell necrosis but no clear sign of blood vessel leakage

# Conclusion

- Increasing importance of *in vitro* pre-clinical test platforms
- Complex organ-on-chip models are used to perform functional assessment of bioactive substances, such as snake venoms
- This requires appropriate, automation capable and relevant 3D organ-on-chip models
- This also requires appropriate, multi-mode data acquisition systems
  - Inclusion of imaging expands the understanding of biological processes
  - Multiplexing with imaging allows for high sensitivity and temporal resolution while allowing morphological assessment and signal localization



# Contributors

Flavio Bonanini<sup>1</sup>, Mátyás Bittenbinder<sup>2,3,4</sup>, Jeroen Kool<sup>3,4</sup>, Freek J. Vonk<sup>2,3,4</sup>, Sebastiaan J. Trietsch<sup>1</sup>, Dorota Kurek<sup>1</sup>

1 Mimetas, Leiden, The Netherlands

2 Naturalis Biodiversity Center, 2333 CR Leiden, The Netherlands

3 Amsterdam Institute of Molecular and Life Sciences, Division of Bioanalytical Chemistry, Department of Chemistry and Pharmaceutical Sciences, Faculty of Science, Vrije Universiteit Amsterdam, De Boelelaan 1085, 1081HV Amsterdam, The Netherlands

4 Centre for Analytical Sciences Amsterdam (CASA), 1098 XH Amsterdam, The Netherlands

 **TECAN.**

**MIMETAS**

Spark Cyto is for Research Use Only. Not for use in diagnostic procedures.