

Most Scalable Multi-organ System

Discover Akura™ Flow 384

Flexible Pick-and-Place **Organ-Organ** Crosstalk

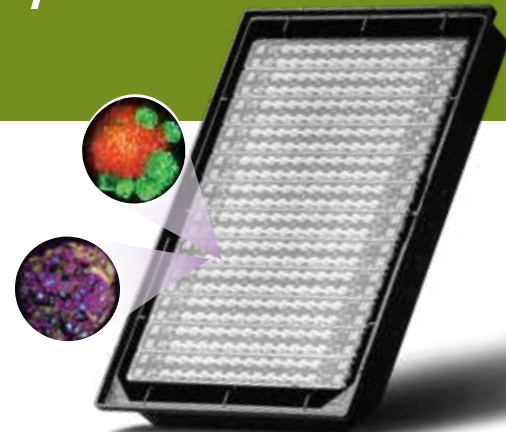
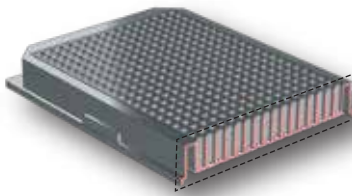
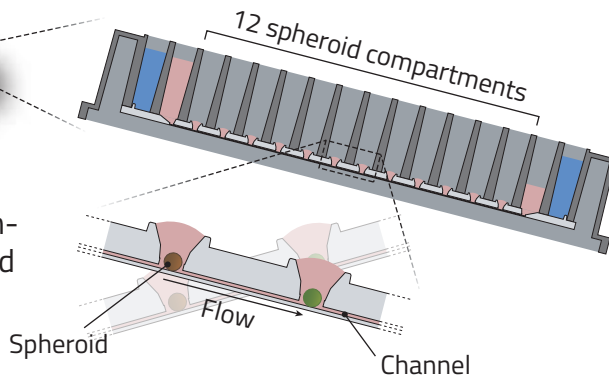


Plate Layout

- 384-well format engineered for spheroids

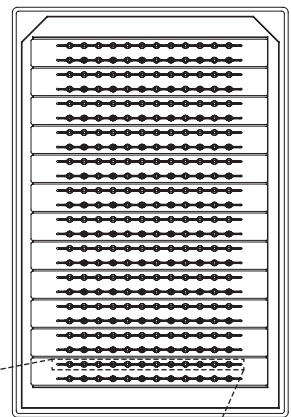


- Gravity-based perfusion-control through repeated back and forth tilting

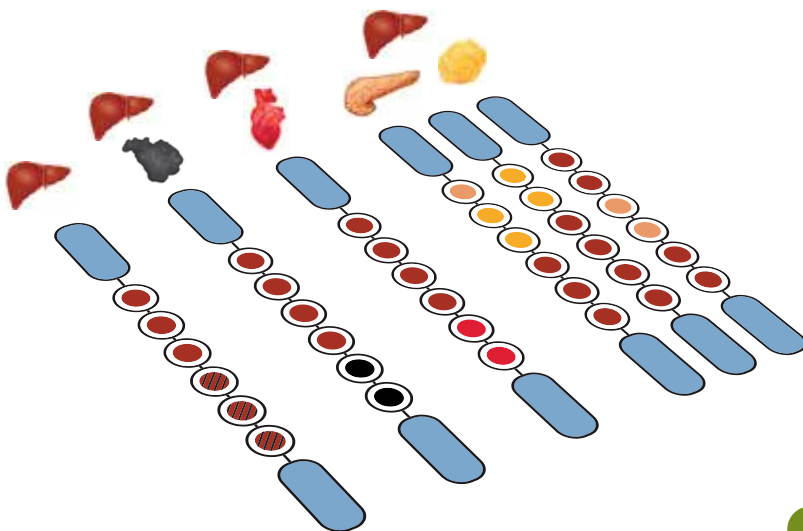


- Interconnecting up to 12 spheroids by a microfluidic channel

- High-density arrangement of **24x 12-tissue** channels



Applications



Signaling between Healthy and Diseased Tissue

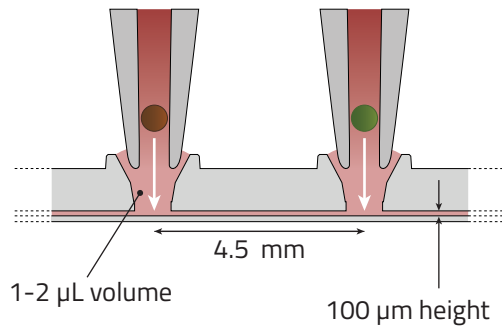
Pro-drug Activation and Drug-drug Interaction

Metabolite Toxicity on On-target and Off-target Tissues

Metabolite Disease Modeling



Unique Features



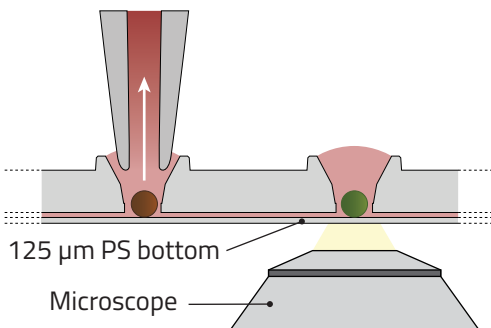
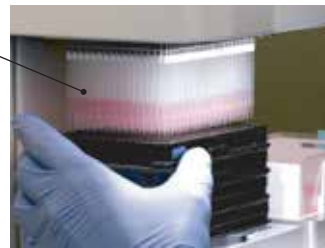
Experience a highly flexible multi-organ platform

- Conical well design with SureXchange™ ledge for easy gravity-based loading and access of spheroids
- Quick plug-and-play assembly of different microtissue types and seamless integration of additional models in the future
- Spheroid compartment with minimal dead volume enabling very precise medium exchange and drug treatment and low experimental variation

Rely on universal plate standards and low operational complexity

- ANSI/SLAS standard format quick implementation widely used lab tools
- Full compatibility and with automation and robotic liquid handlers
- Gravity-based and pumpless flow control for simple plate handling

Multi-channel systems



Access a wide range of readout methods

- Anytime access to supernatants using manual or automated pipetting
- Parallel retrieval of spheroids for lytic assays, histology or RNA sequencing
- In-situ high-content imaging with minimal optical aberration with a flat and transparent PS bottom

Workflow

