

Multi-organ and Immune-Competence on top of our Best Spheroid Microplate

Discover Akura™ Twin Microplate

Multi-organ **Immune** Response Made **Scalable**

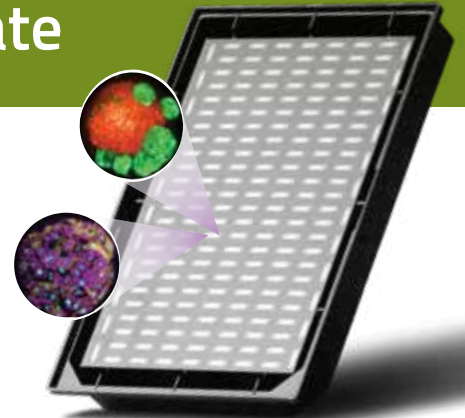
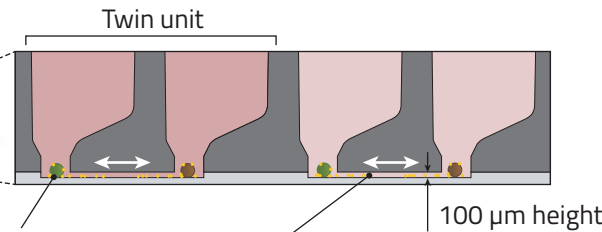
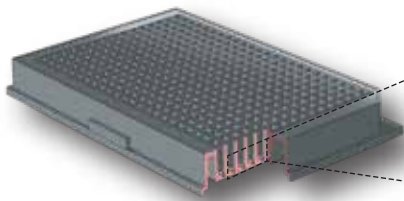


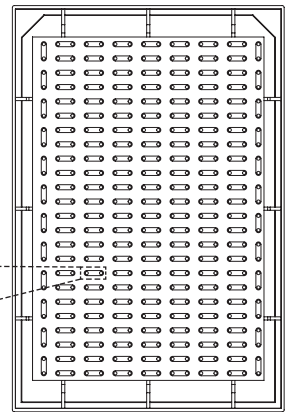
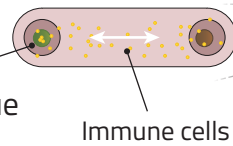
Plate Layout

- 384-well format engineered for spheroids (based on Akura™ 384)

- High-density array with **192-conditions** per including no-flow control

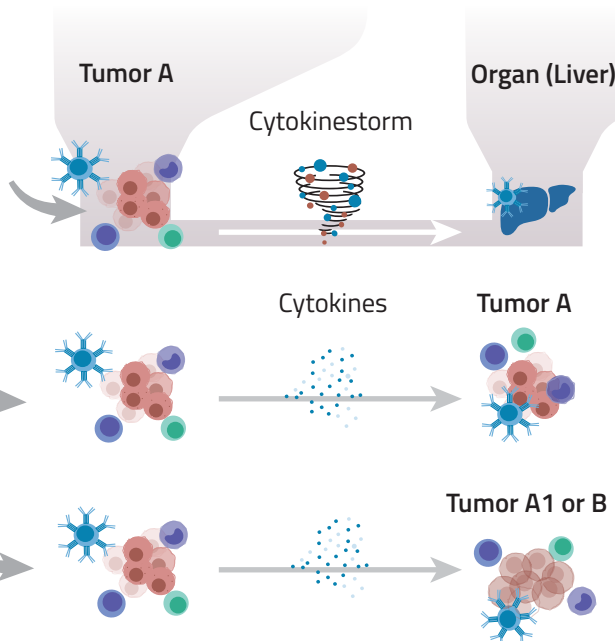


- One spheroid/organoid per well
- Two adjacent wells are interconnected by a microfluidic channel enabling active tissue-tissue cross-talk and immune cell migration



Pharmacological treatment

- Engineered cells
- Small molecules
- Antibodies



Applications

Immune Safety: iDILI Cytokine release syndrome

Organ of same tumor origin, different organ (healthy or diseased)

Test the Apscopal effect

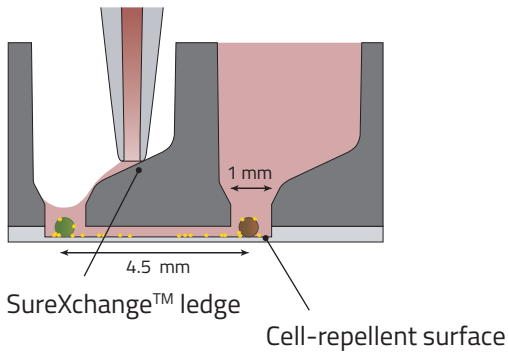
Test the ability of the treatment modalities to induce a systemic immune response allowing tumor regression in the non-treated area

Test cancer heterogeneity

Different tumor cell line, resistive tumor, tumor of other organ



Unique Features



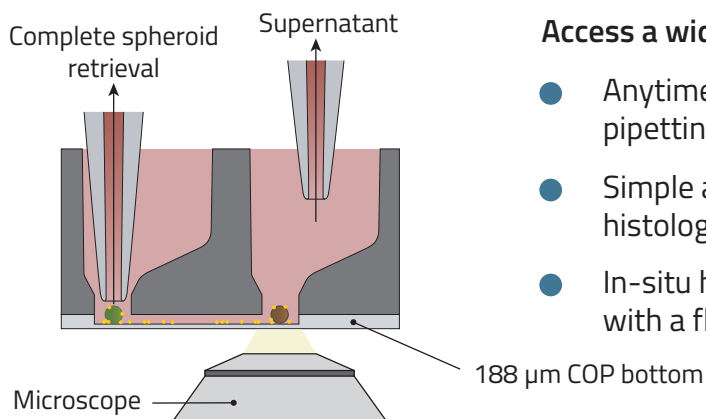
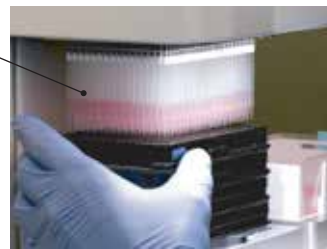
Experience an intuitive plate and well architecture

- Conical well design with SureXchange™ ledge protecting the spheroid/organoid cells from aspiration
- Near-complete medium exchange for precise dosing and low experimental variation
- Stable ULA coating in the well and microchannel preserves spheroid/organoid morphology and prevent immune cell adhesion in long-term culture

Rely on universal plate standards

- ANSI/SLAS standard format quick implementation widely used lab tools
- Full compatibility and with automation and robotic liquid handlers

Multi-channel systems



Access a wide range of read of methods

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

Workflow

