

# >> A novel non-invasive workflow for differentiation and characterization of iPSC-derived hepatic organoids

Nathalie Opdam – van de Laar<sup>1,✉</sup>, Inge Thijssen – van Loosdregt<sup>1</sup>, Denise Sullivan<sup>2</sup>

<sup>1</sup>Axion BioSystems, Eindhoven, the Netherlands; <sup>2</sup>Axion BioSystems, GA, USA ✉ Email: nathalie.opdam@axionbio.com

## Omni: Kinetic cell tracking

### Automated whole-vessel imaging

*In vitro* models are essential for studying diseases and development. While traditional 2D cell culture models have provided valuable insights, they often fail to replicate *in vivo* complexity. This has led to increased interest in 3D models such as spheroids and organoids, which better mimic *in vivo* conditions.

Live-cell provides a powerful technique for studying these 3D models, enabling real-time visualization and analysis at defined time intervals.

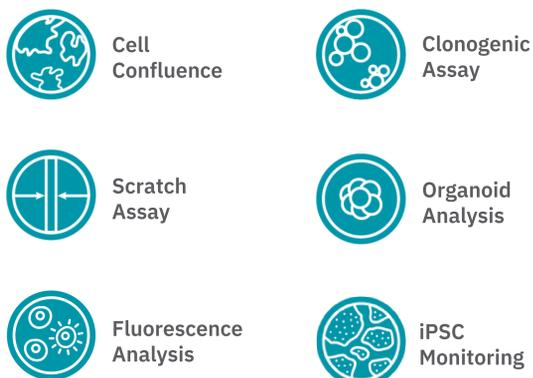


### The Omni product family

- >> Assay your cells in brightfield and fluorescence
- >> Track every moment, straight from your incubator
- >> See every cell by movement of the camera
- >> Monitor and analyze your cells remotely
- >> Get started quickly

### AI-Driven imaging software for powerful, yet simple analysis

The Omni platform software modules simplify assay setup, offer real-time cellular visualization, and enable fast analysis.



## Real-time Monitoring of iPSCs

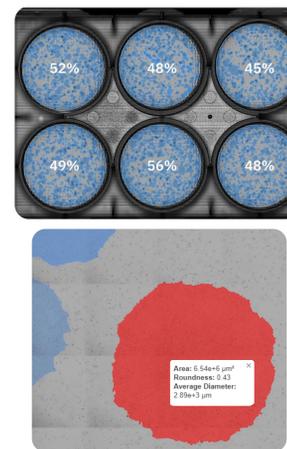
### Background

The liver plays vital roles in detoxification, protein synthesis, metabolism, and hormone regulation. While it regenerates efficiently *in vivo*, expanding hepatocytes *in vitro* is difficult. Induced pluripotent stem cells (iPSCs) provide a versatile source for generating hepatic cells and organoids that mimic liver structure and function. This study presents a novel non-invasive workflow for monitoring iPSC-derived hepatic organoid development using the Omni.

### iPSC Module tracks iPSC growth

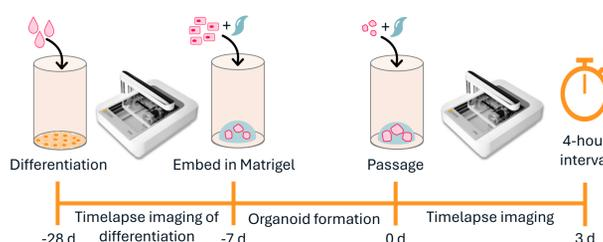
The Omni is a live-cell analysis platform is capable of continuous multi-well imaging directly from the incubator.

Example whole-well brightfield images of iPSC colonies acquired by the Omni in a 6-well plate with the confluency map overlay and a close up of a single colony with the metrics provided by the iPSC Module including area, diameter, and roundness.



### Methods

iPSCs were differentiated into hepatic progenitor cells (HPCs) and hepatocyte-like cells (HLCs) and subsequently formed into organoids. Organoid area, diameter, and roundness were determined every 4 h for 3 days.



## iPSC-derived organoid workflow

### iPSC differentiation

Distinct morphological changes occurred during differentiation: iPSCs were compact, definitive endoderm cells smaller, HPCs larger with a cobblestone appearance, and HLCs the largest with a similar arrangement.

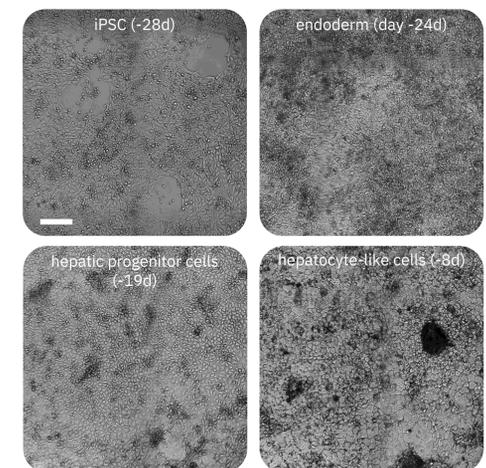


Figure 1: Morphological changes during differentiation of iPSC towards hepatocyte-like cells. Scalebar is 200 µm and accounts for all images.

### Hepatic organoid formation

HPC- and HLC-derived organoids were spherical with lumens and had similar roundness (0.85). Initially smaller, HPC-derived organoids grew 3.6-fold over 72 hours, compared to 3.0-fold for HLC-derived organoids, reflecting differences that may be related to cell maturity.

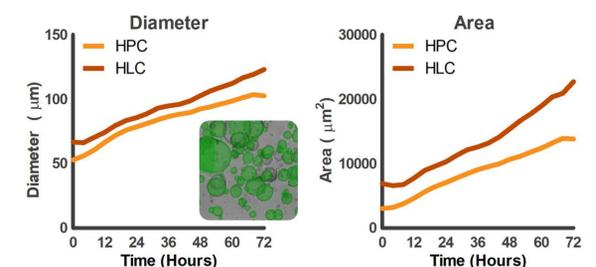


Figure 2: Change in organoid diameter and area over time for the hepatic progenitor cell derived organoids (HPC) and the hepatocyte-like cell derived organoids (HLC), including an example of organoid detection (green overlay) at 72h by the Omni.

### Conclusion

This workflow demonstrates the power of live-cell imaging for real-time monitoring of iPSC differentiation and hepatic organoid formation, enhancing liver research and applications in disease modelling and therapy.