

# Optimizing Human Hepatic Organoid Development: Leveraging iPSC Differentiation and Advanced Analytical Platforms with RUO Growth Factors and Cytokines

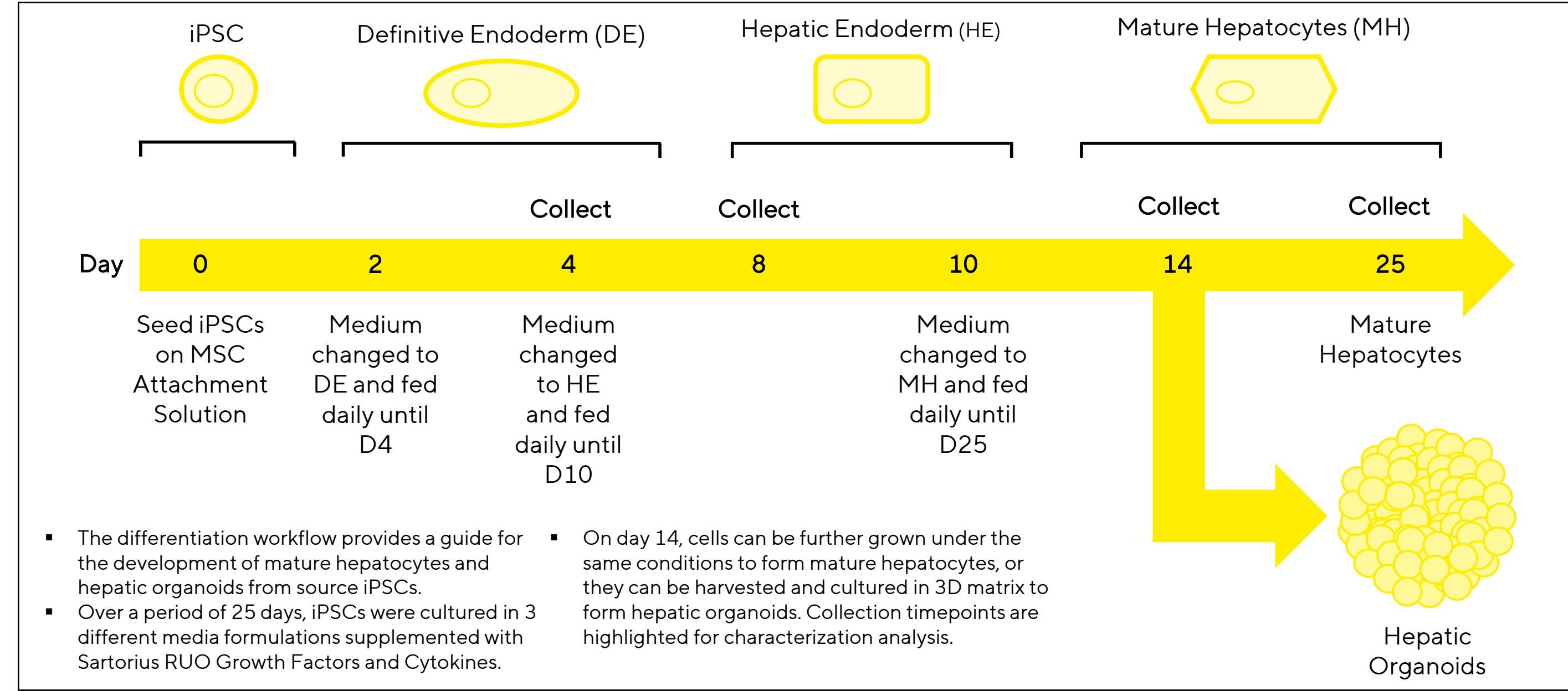


N. Lewis\*, D. Cole, A. Ward, J. Trigg and N. Bevan  
BioAnalytics, Sartorius (Essen BioScience), Royston, Hertfordshire, UK;  
\*Corresponding author: Natasha.Lewis@sartorius.com

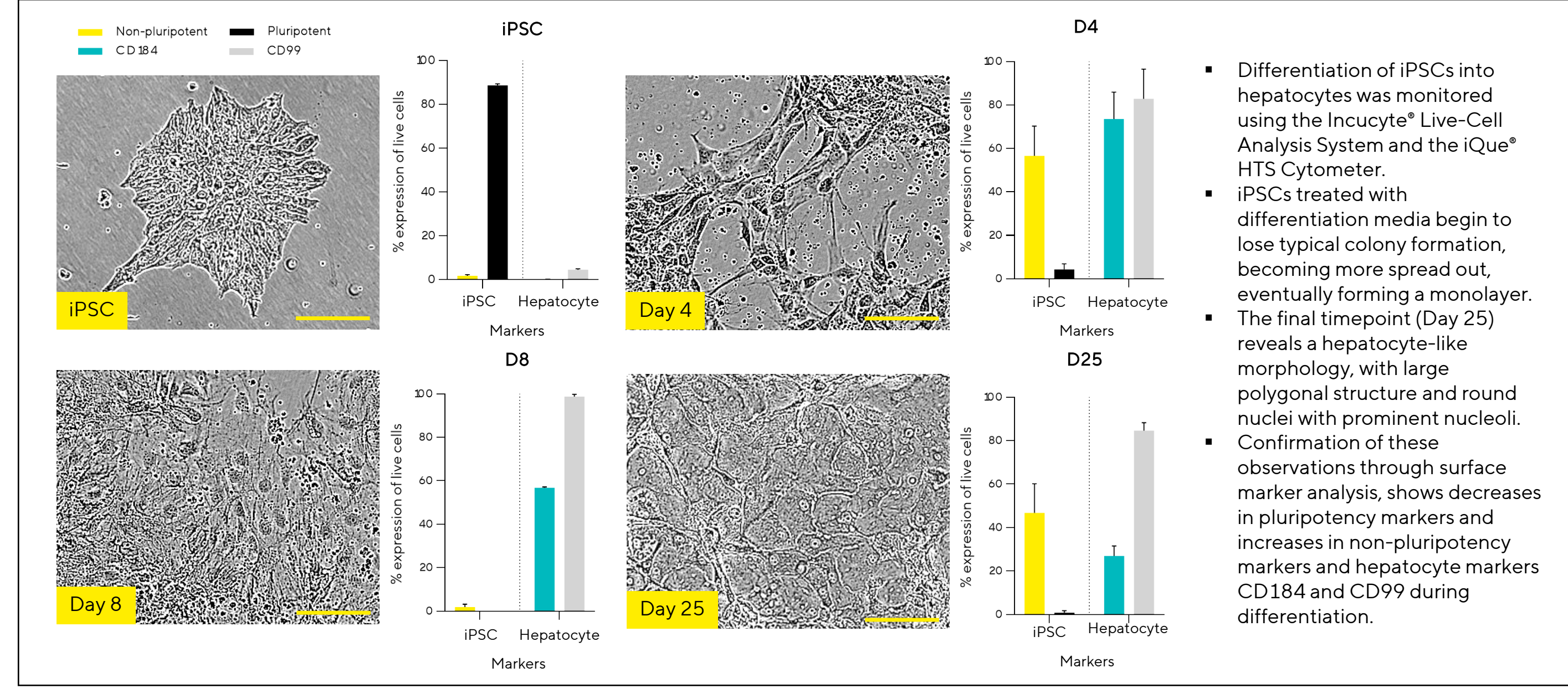
## Introduction

- Increasing requirement for the generation of species-specific, physiologically relevant systems for the study of disease and development has led to the expansion of 3D models, such as organoids, used in research.
- These biologically accurate tissue models can be expensive and time-consuming to produce, necessitating efficient and reliable methods for their development, monitoring and characterization.
- The utility of stem cells in laboratory and clinical research has been clearly demonstrated through an extensive number of studies exemplifying stem cells as a model for a multitude of diseases and therapeutic development.
- Differentiation of induced pluripotent stem cells (iPSCs) into somatic tissues requires specific conditions dependent on the desired tissue type.
- In this workflow, human iPSCs were differentiated into hepatocytes using three distinct media types with their own cytokine and growth factor supplementation at key stages of differentiation, providing a relatively simple method for producing mature hepatic cells.
- This enabled the production of human hepatic organoids from these source cells, demonstrating a robust method for unlimited organoid development.
- Here, we outline a simple, standardized, and robust workflow using Sartorius RUO Growth Factors and Cytokines in combination with the Incucyte® Live-Cell Analysis System and the iQue® HTS Cytometer for phenotypic characterization and monitoring of growth.
- This approach simplifies the generation of functional hepatic cells and organoids from iPSCs for drug discovery, development, and toxicity studies.

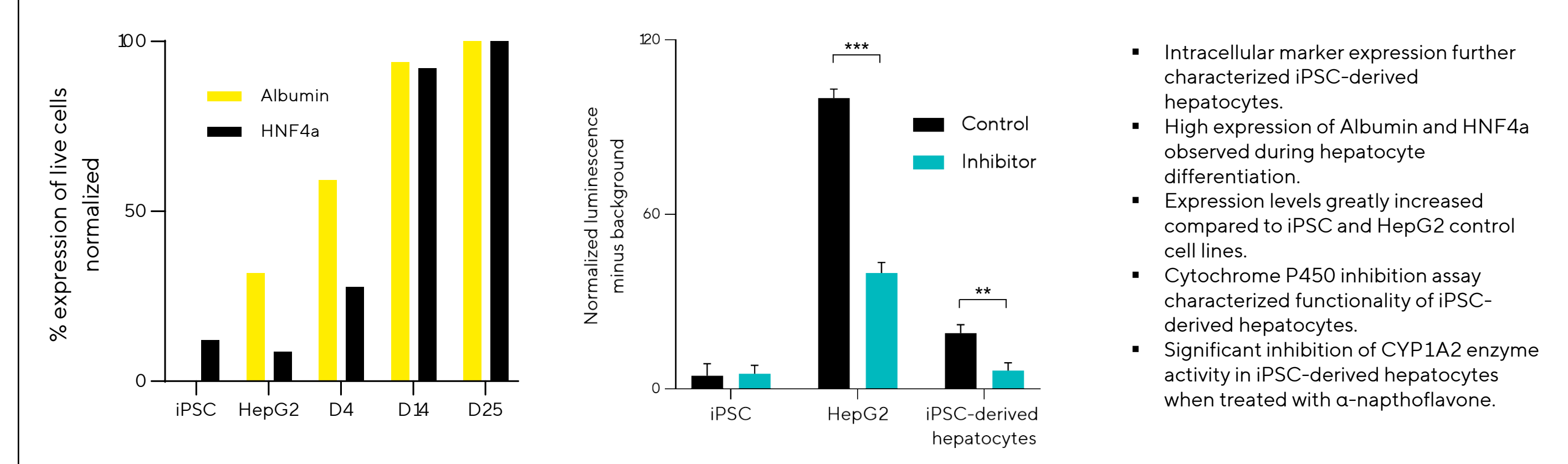
## Differentiation Workflow



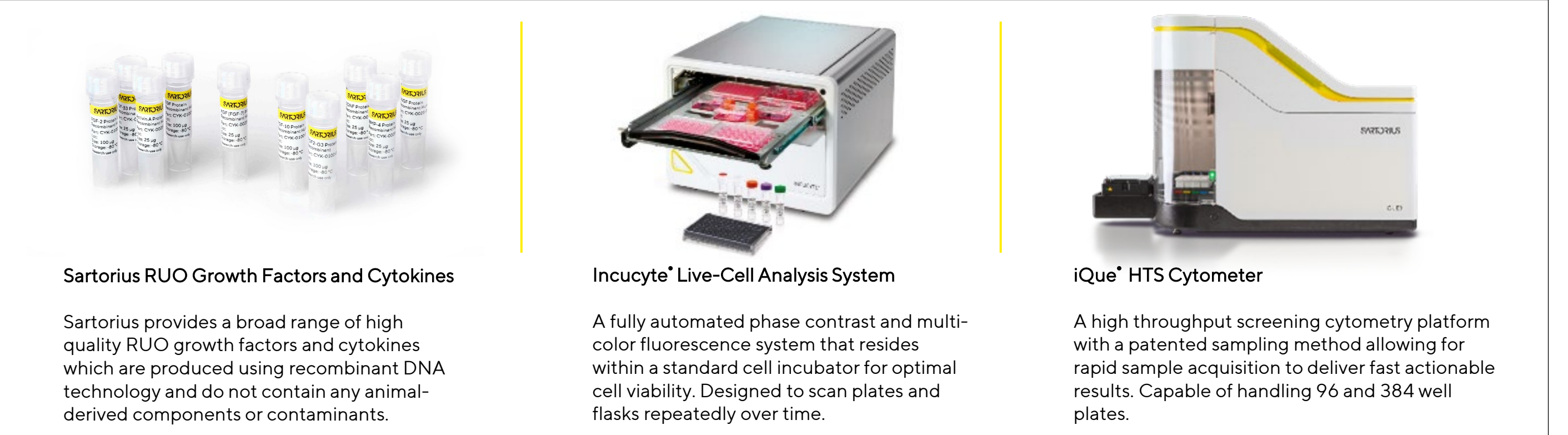
## Monitoring iPSC to hepatocyte differentiation



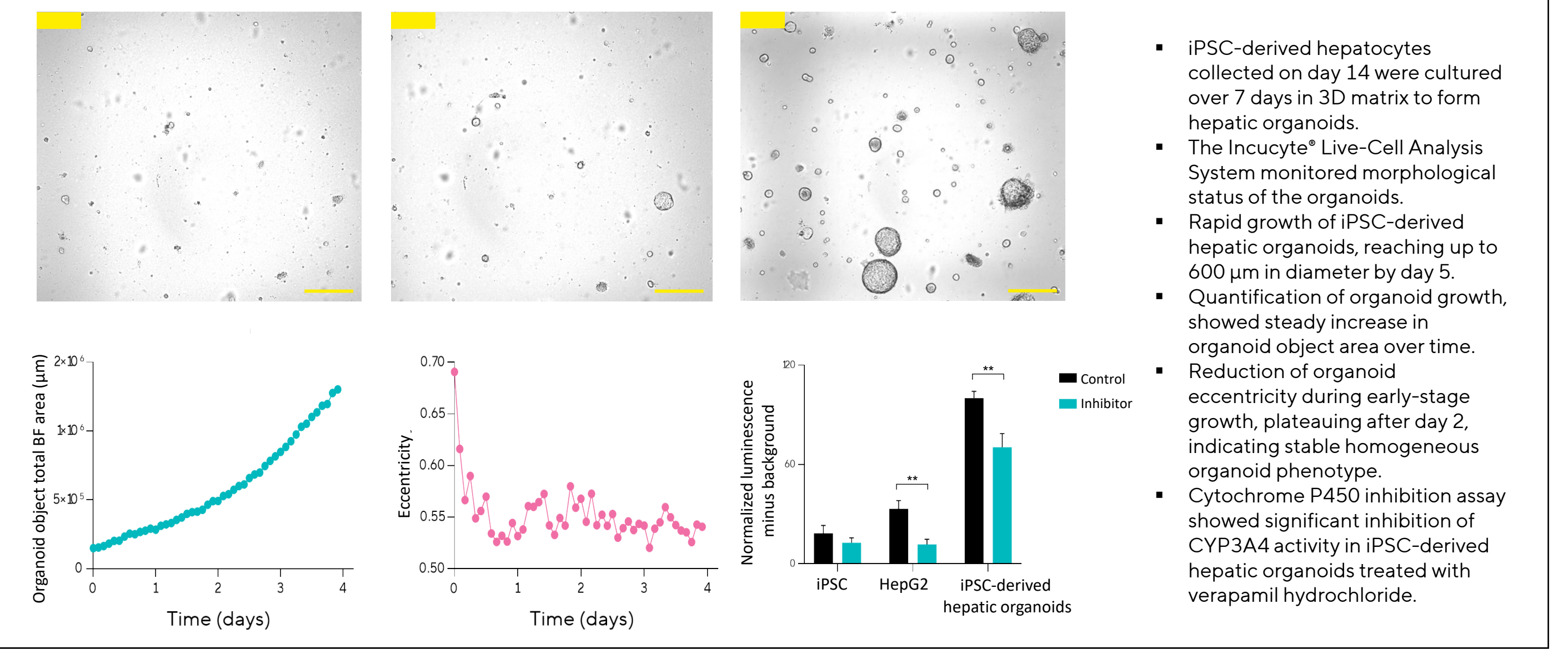
## Characterizing expression and functional activity of iPSC-derived hepatocytes



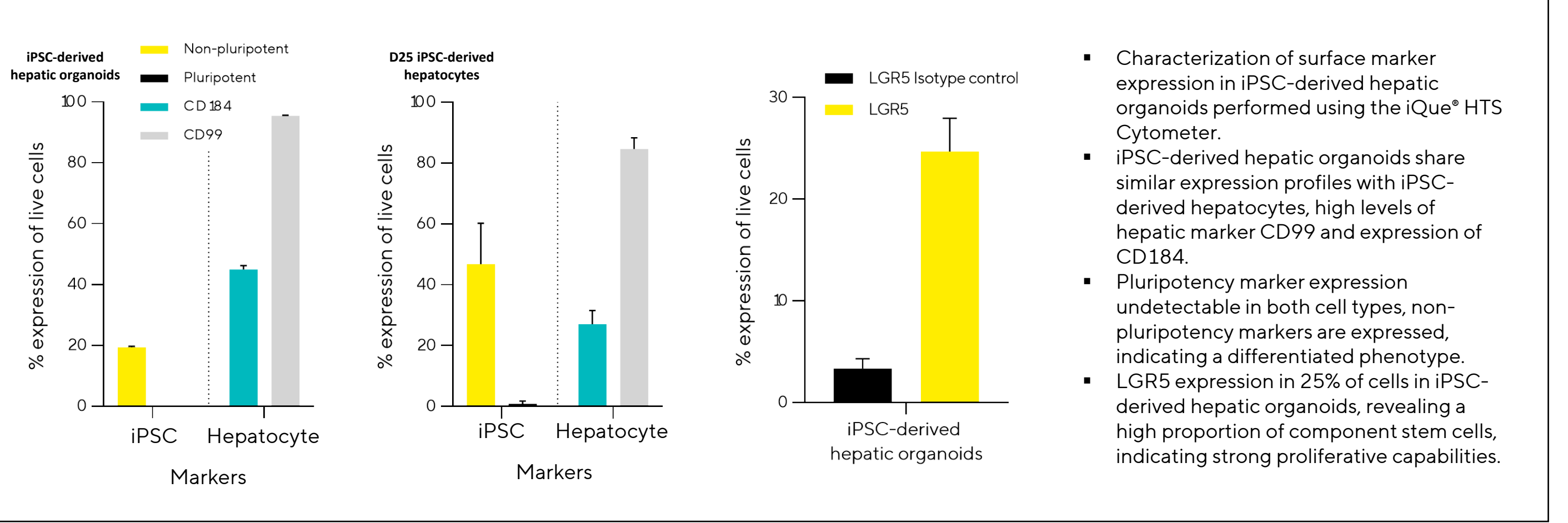
## RUO Growth Factors and Cytokines, Incucyte® & iQue® Systems



## Monitoring iPSC-derived hepatic organoid development



## Characterizing iPSC-derived hepatic organoids



## RUO growth factors and cytokines for organoid culture

