

Automating 3D culture model systems: Delivering robust, reproducible solutions to complex tissue model culture

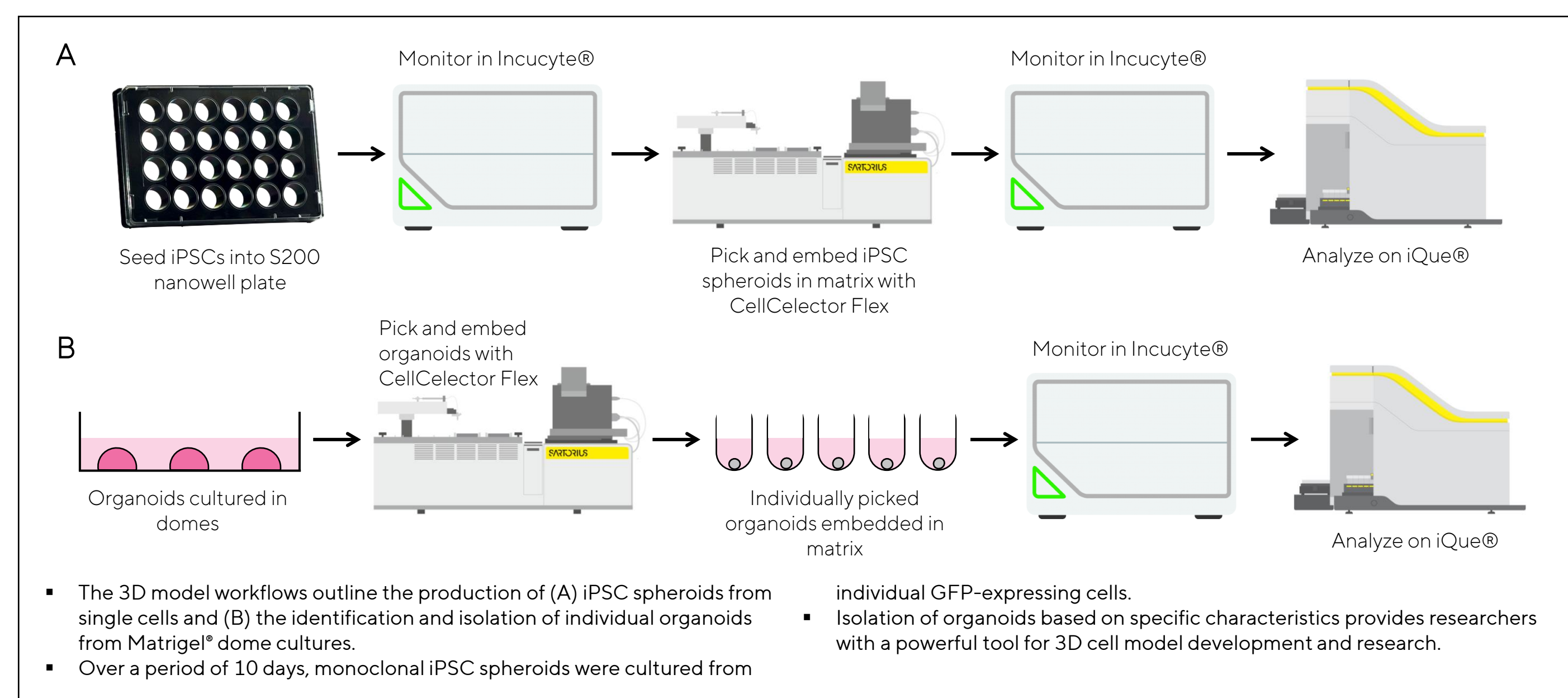
D. Cole* and N. Bevan

BioAnalytics, Sartorius (Essen BioScience), Royston, Hertfordshire, UK;
*Corresponding author: Daryl.Cole@sartorius.com

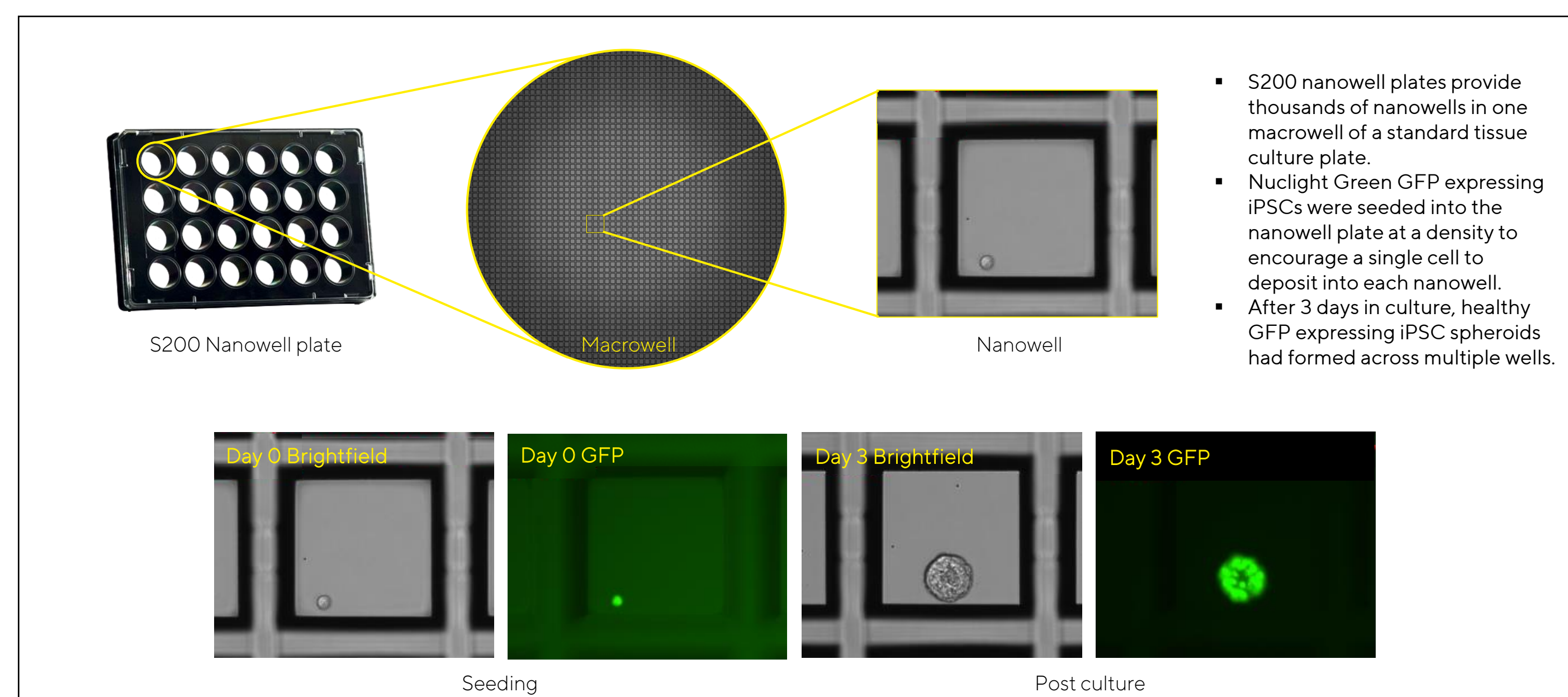
Introduction

- 3D models such as spheroids and organoids are becoming increasingly relevant for the study of disease and development.
- Generation of these models can prove to be highly complex, expensive and resource intensive, necessitating efficient and reliable methods for their development, monitoring and characterization.
- Further complexity is inherent in the production of disease models that contain populations of cells with specific markers or genetic mutations associated with disease.
- Development of such models requires accurate manipulation of cells to successfully create a physiologically relevant system.
- In this study, spheroids were cultured from individual iPSCs to produce clonal iPSC spheroids derived from a single cell, while human iPSC derived hepatic organoids were selected from Matrigel® domes based on their attributes and transferred for further culture and analysis.
- Here, we outline a simple, standardized, and robust workflow using CellCelector Flex to identify and pick 3D cell models for further development based upon key growth and morphological attributes.
- During the workflow, the Incucyte® Live-Cell Analysis System and the iQue® HTS Cytometer were used to monitor growth and morphological phenotype over time and phenotypic characterization through marker expression analysis.
- This approach simplifies the culture and generation of organoids and monoclonal spheroids from iPSCs for drug discovery, development, and toxicity studies.

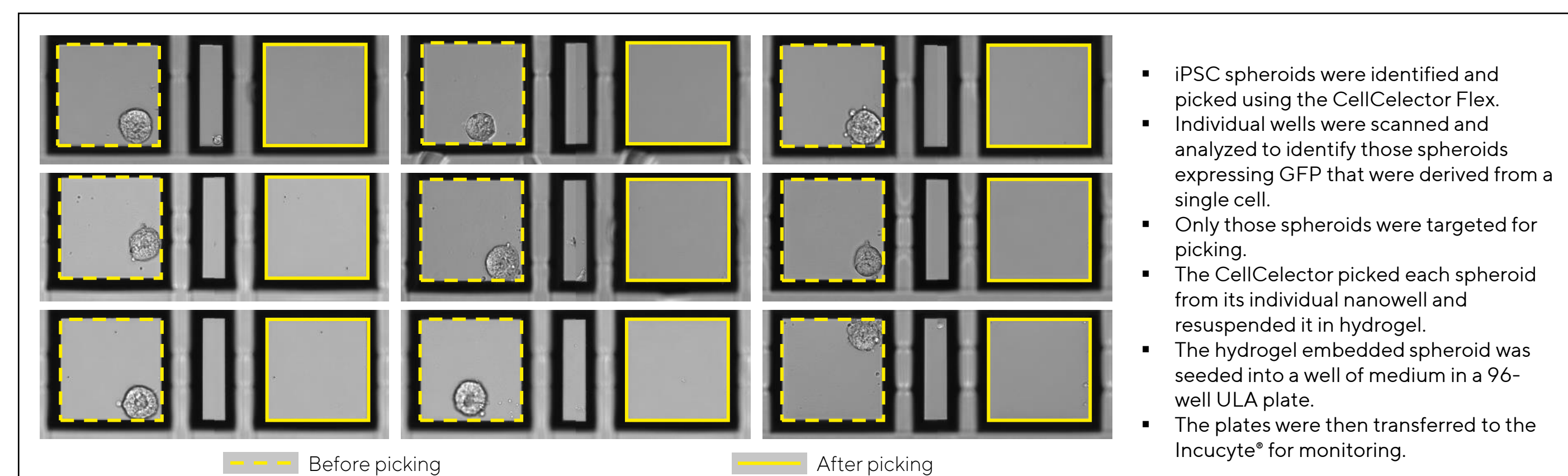
3D model workflows



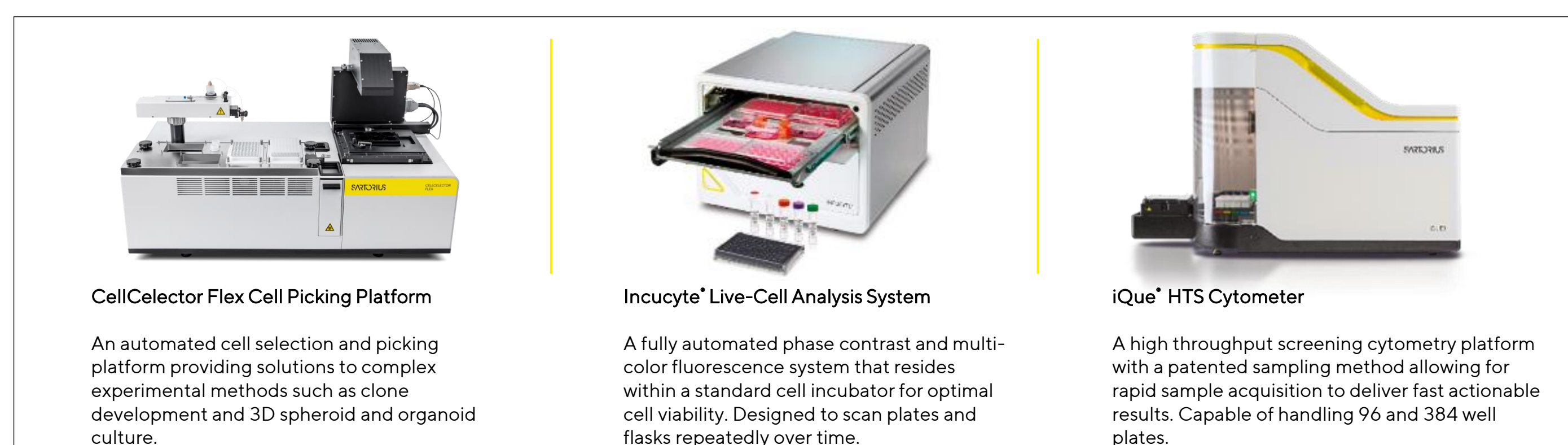
Monoclonal iPSC spheroid generation



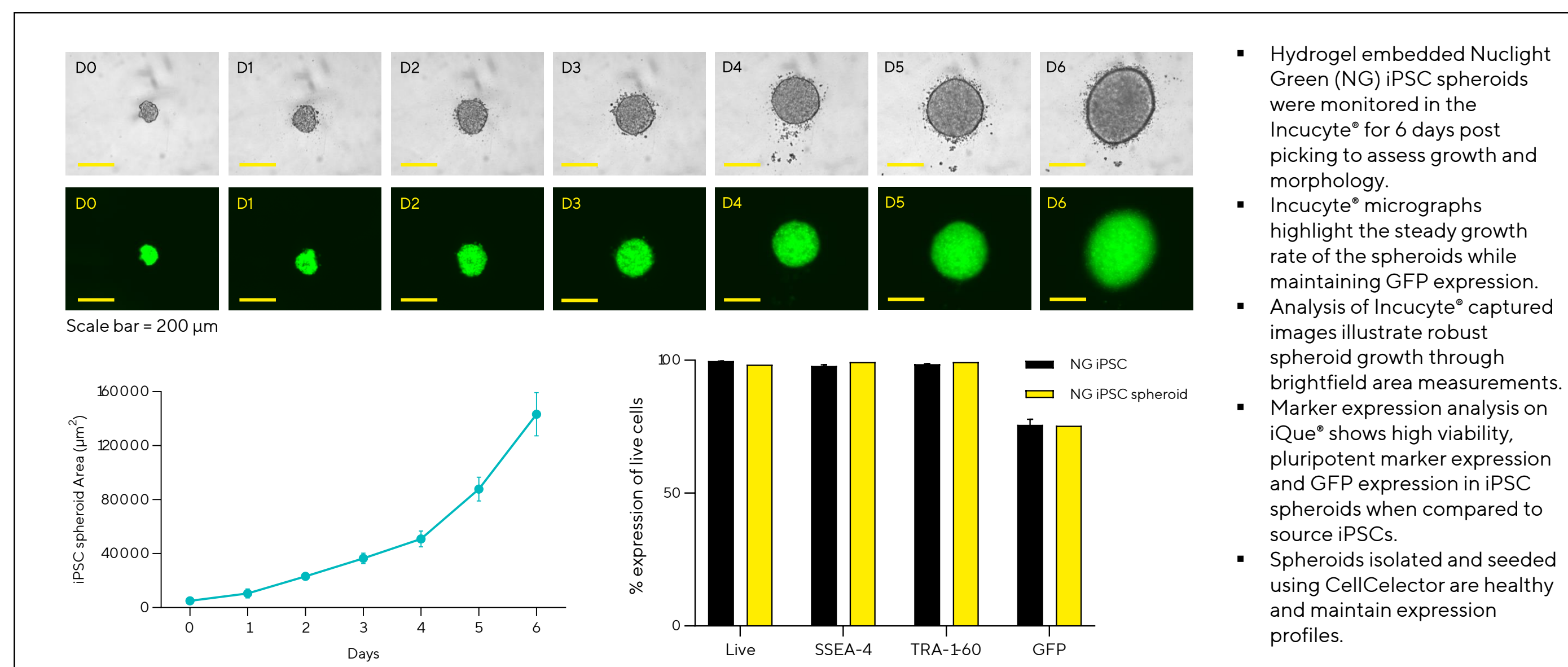
Picking and embedding iPSC spheroids



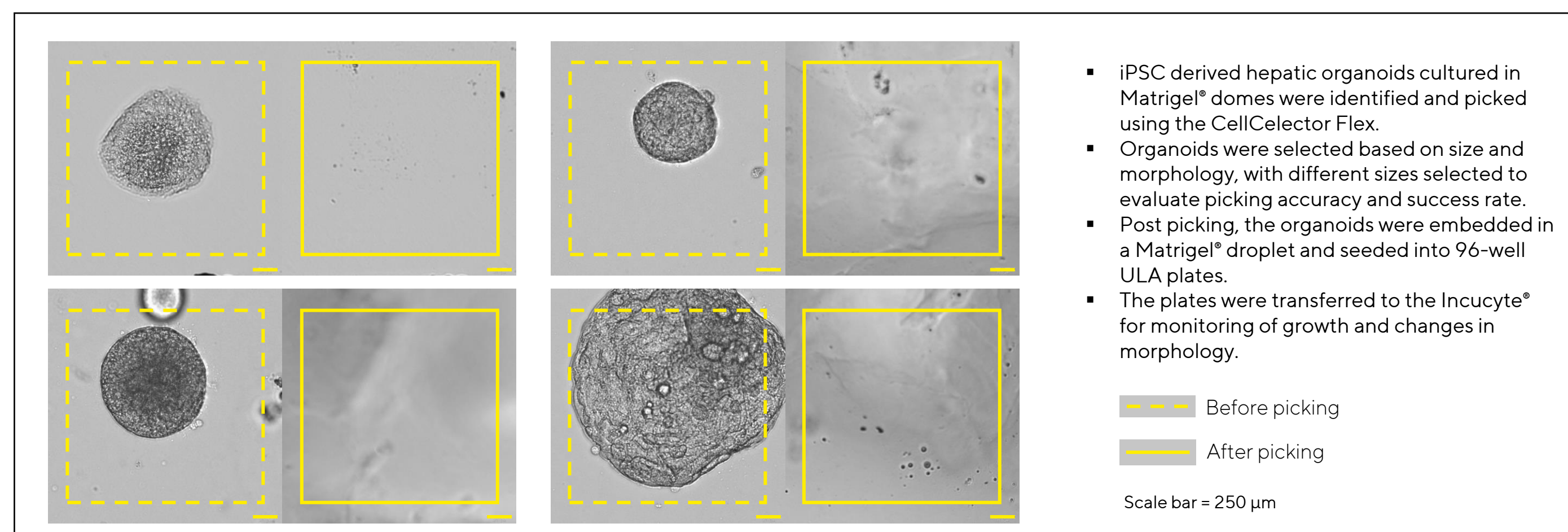
CellCelector Flex, Incucyte® & iQue® Systems



Monitoring and characterizing iPSC spheroids



Identification and isolation of organoids



Monitoring and characterizing organoid development

