

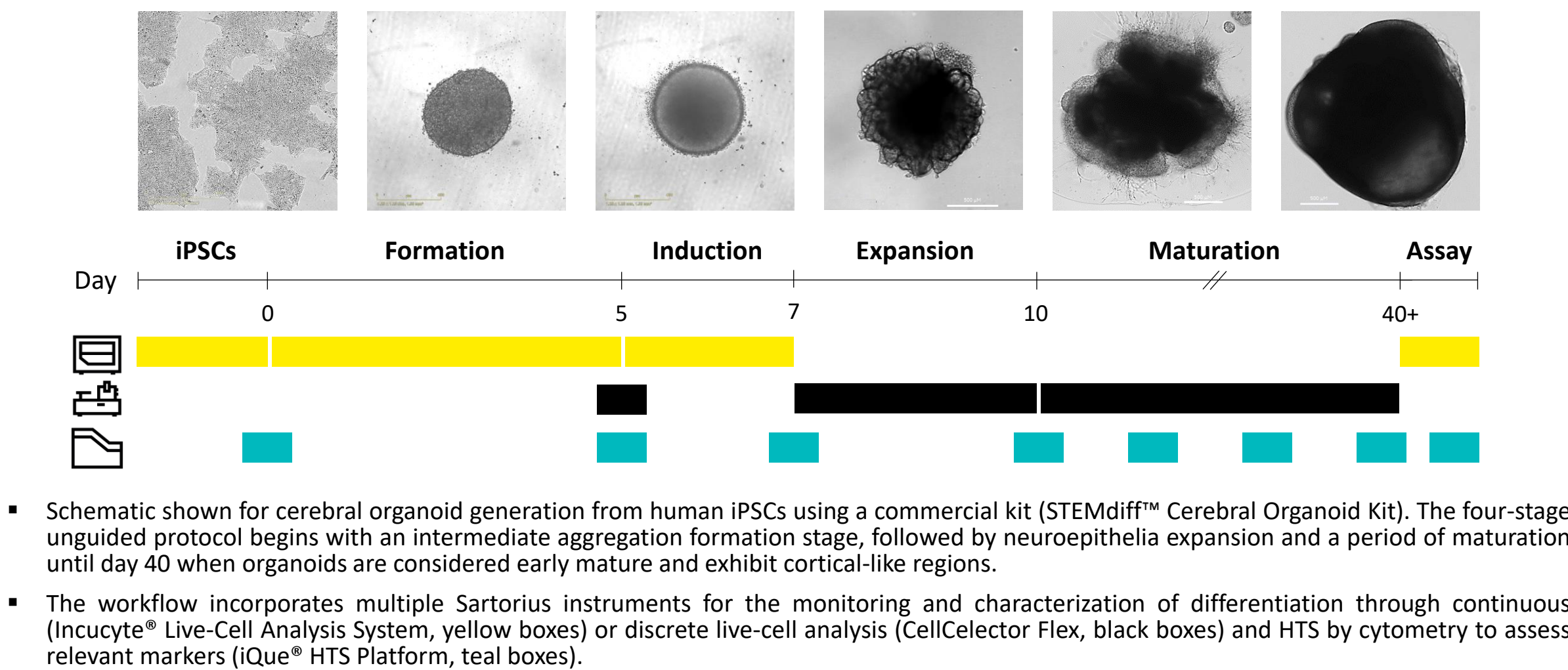
Monitoring Brain Organoid Differentiation and Development Leveraging Live-Cell Imaging and Cytometry

J. Trigg*, D. Cole, K. Barnes, and N. Bevan
Sartorius, Royston, UK; *Corresponding author: Jasmine.Trigg@Sartorius.com

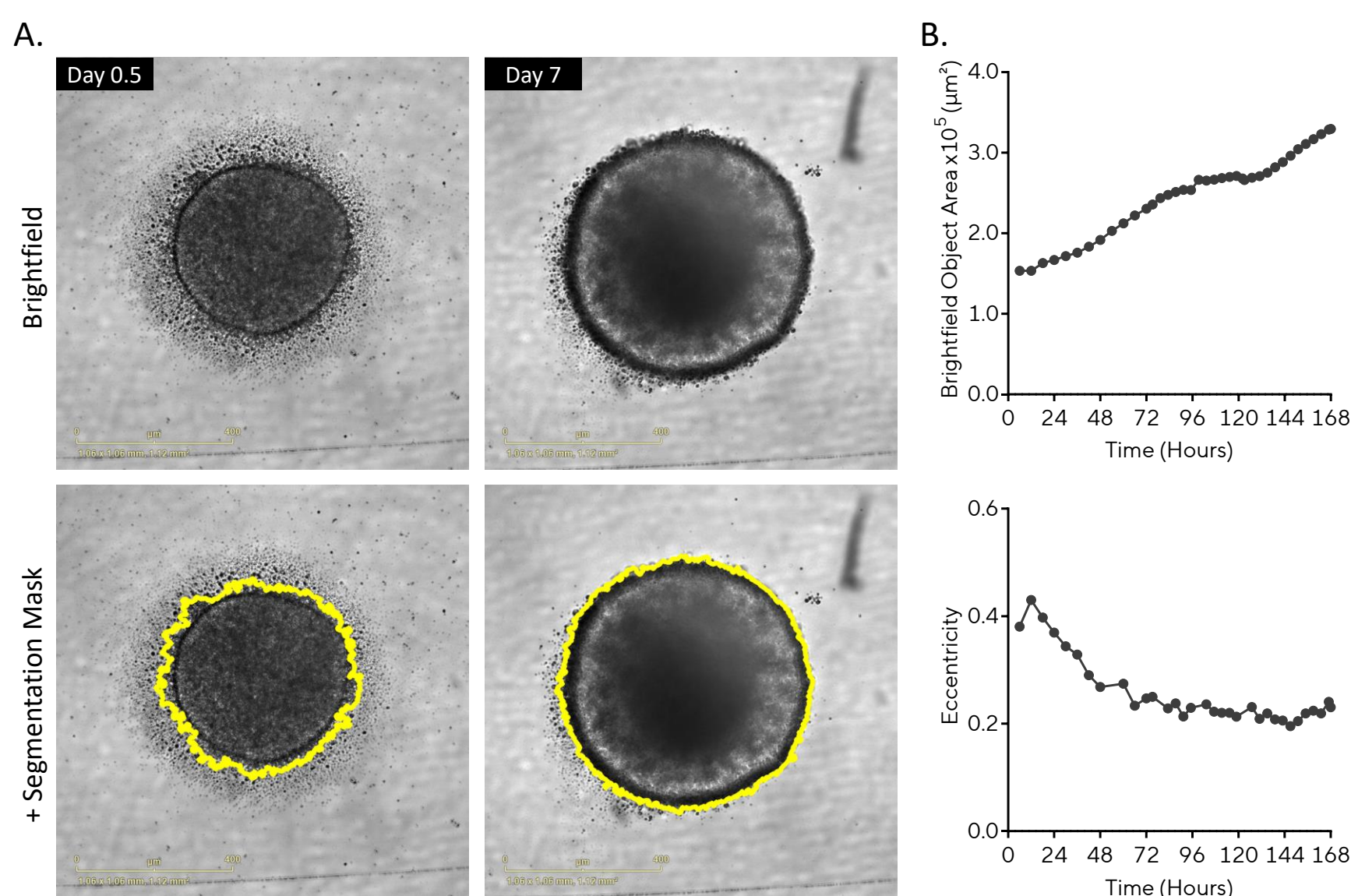
Introduction

- hPSC-derived brain organoids are physiologically relevant 3D *in vitro* models that mimic human brain development and cellular organization.
- They are increasingly employed to study neurodevelopment, neuropathology, and drug screening but face challenges due to long protocols, sensitivity, large size, heterogeneity, and limited temporal methods for objective monitoring.
- This workflow incorporates live-cell analysis and high-throughput screening (HTS) by cytometry to monitor cerebral organoid differentiation.
- The data exemplifies how this approach enables objective assessment of cerebral organoid differentiation by temporally monitoring growth, morphology, and marker expression, to support data-driven QC decisions and neurodevelopmental or toxicity studies.

iPSC-derived Cerebral Organoid Differentiation Workflow

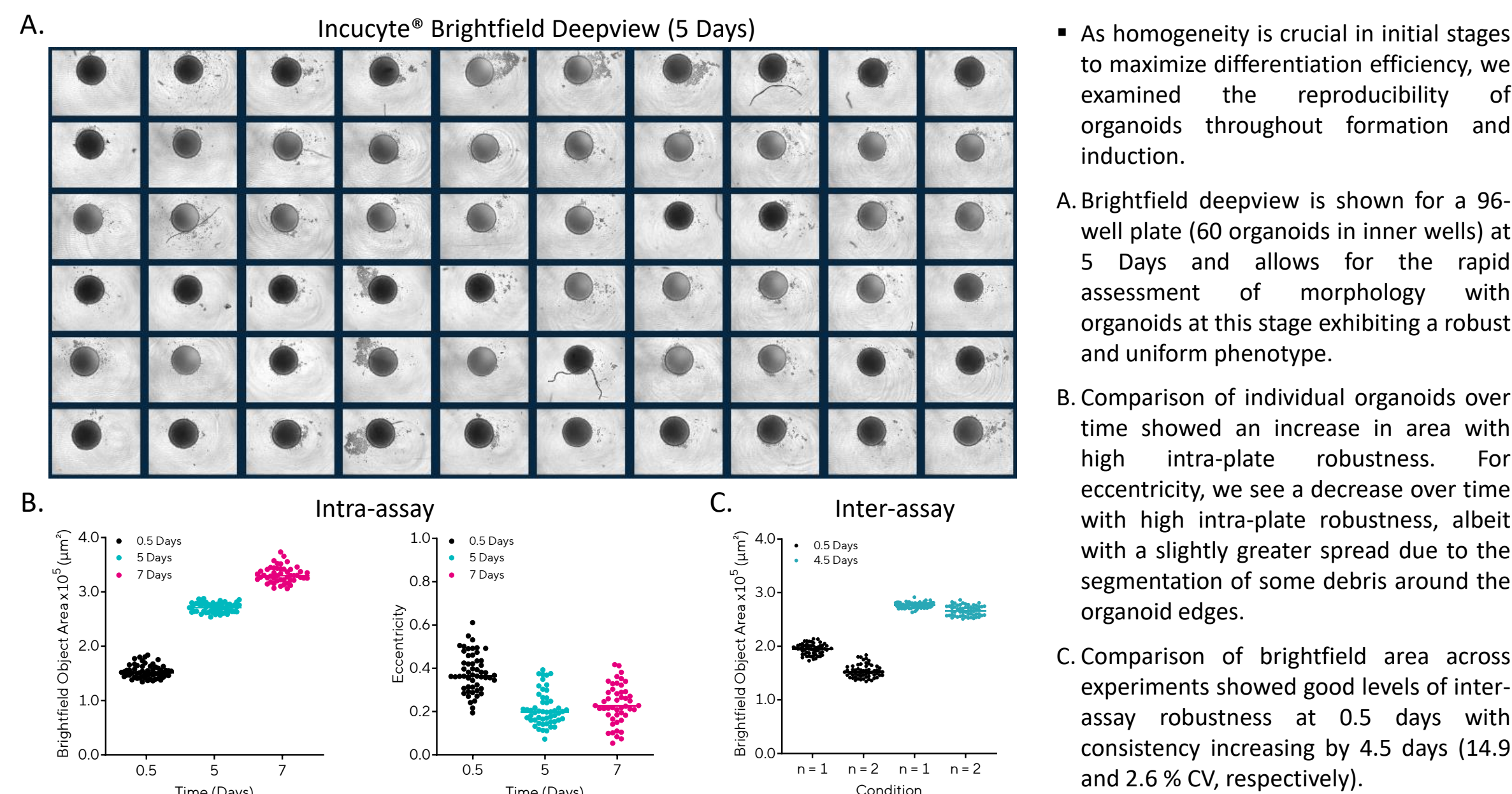


Continuous Live-Cell Analysis of Organoid Formation and Induction

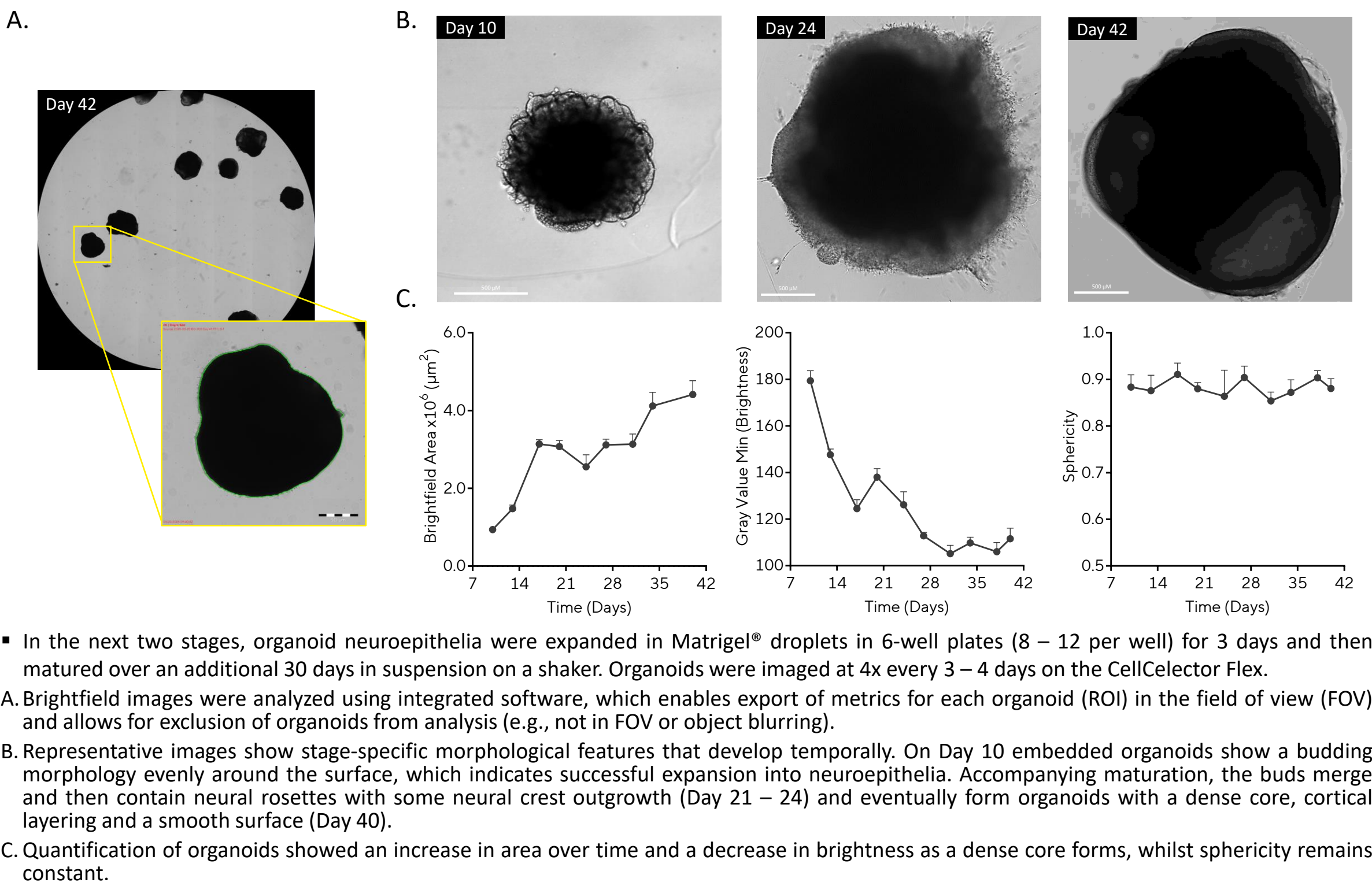


- Organoid formation and induction were performed in 96-well round-bottom ULA plates and images were acquired using the Incucyte® spheroid scan type every 6 hours over 7 days.
- Representative timelapse brightfield images show rapid aggregation of iPSCs into compact organoids with the expected morphology. Organoids were symmetrical and exhibited smooth round edges with rosette-like patterning by day 7. Integrated analysis software accurately segmented the organoids (yellow outline).
- Quantification revealed a temporal increase in organoid area (top graph) and an initial decrease in eccentricity (bottom graph) with aggregation that then remained relatively stable.

High Assay Robustness for Organoid Formation and Induction

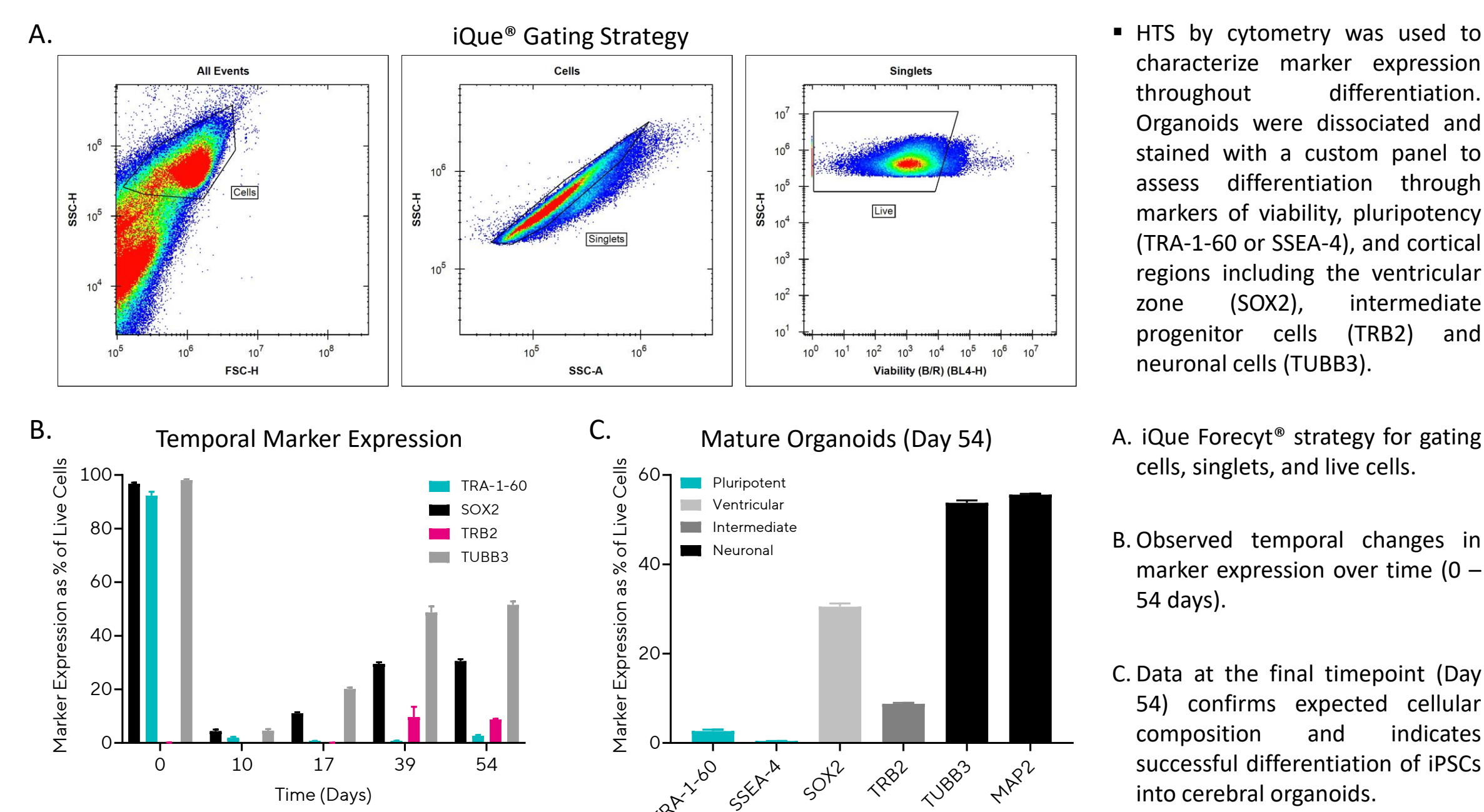


Quantification of Organoid Maturation using Discrete Live-Cell Analysis



- In the next two stages, organoid neuroepithelia were expanded in Matrigel® droplets in 6-well plates (8 – 12 per well) for 3 days and then matured over an additional 30 days in suspension on a shaker. Organoids were imaged at 4x every 3 – 4 days on the CellCelector Flex.
- Brightfield images were analyzed using integrated software, which enables export of metrics for each organoid (ROI) in the field of view (FOV) and allows for exclusion of organoids from analysis (e.g., not in FOV or object blurring).
- Representative images show stage-specific morphological features that develop temporally. On Day 10 embedded organoids show a budding morphology evenly around the surface, which indicates successful expansion into neuroepithelia. Accompanying maturation, the buds merge and then contain neural rosettes with some neural crest outgrowth (Day 21 – 24) and eventually form organoids with a dense core, cortical layering and a smooth surface (Day 40).
- Quantification of organoids showed an increase in area over time and a decrease in brightness as a dense core forms, whilst sphericity remains constant.

Characterization of Dynamic Changes in Marker Expression



Assessment of Ethanol Exposure During Brain Organoid Development

