

Label-free, real-time live cell assays for 3D Organoids embedded in Matrigel®.

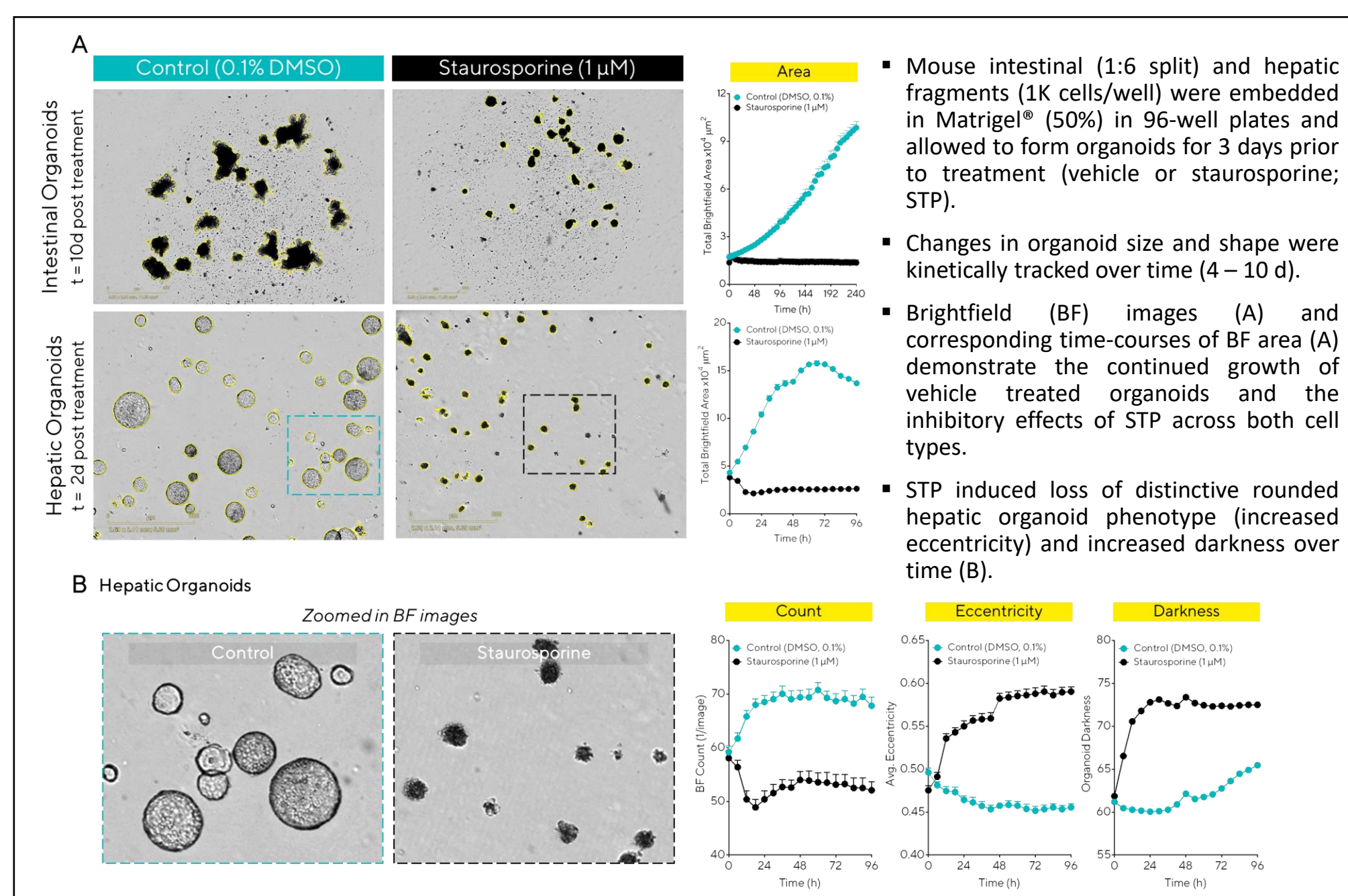
T. Jackson¹, M. Oliver¹, D. Appledorn², T. Dale¹, K. Barnes¹

BioAnalytics, Sartorius (Essen BioScience), Royston, Hertfordshire, UK¹ & Ann Arbor, MI, 48108, USA²

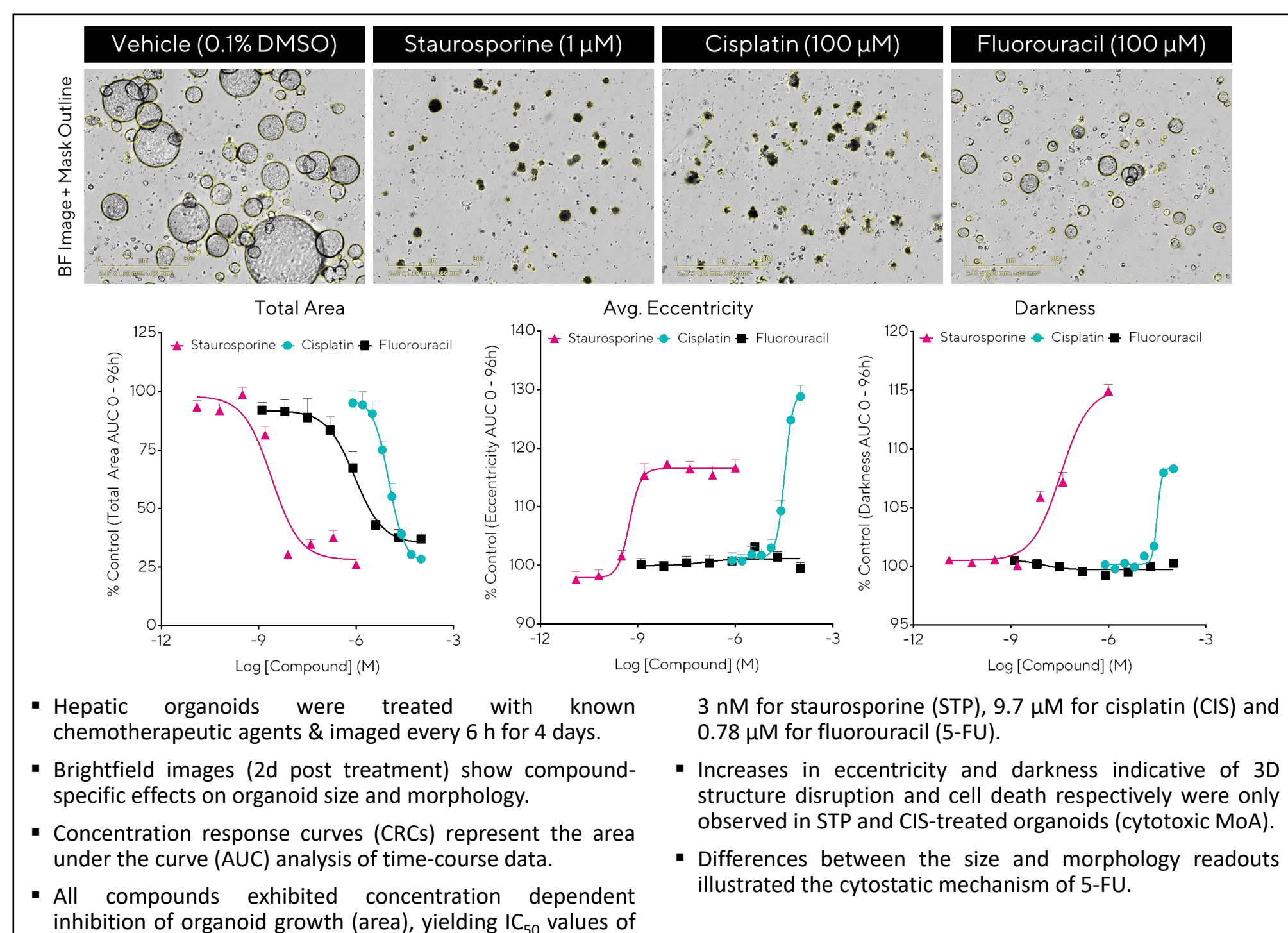
Summary & Impact

- As organoids exhibit structural, morphogenetic, and functional properties that recapitulate *in vivo* pathophysiology, they are increasingly being used *in vitro*.
- To successfully use these models across a variety of research disciplines and applications, technology pipelines to image & quantify these complex structures are key.
- Here, we demonstrate simple, robust workflows for monitoring and quantifying organoid growth, death and morphology.
- Incucyte®s Organoid Analysis Software Module enables the ability to kinetically visualize and quantify distinct organoid morphologies embedded in Matrigel®.
- These validation data demonstrate the ability to characterize the differentiation and maturation of organoid cultures in 24-well plates and assess treatment effects on organoid growth in 96-well microplates.
- Integrated, label-free size and morphology metrics enabled real-time elucidation of compound mechanisms of action and assessment of CFTR function *in vitro*.
- These data exemplify the amenability of this approach for real-time compound profiling across a range of disease areas.

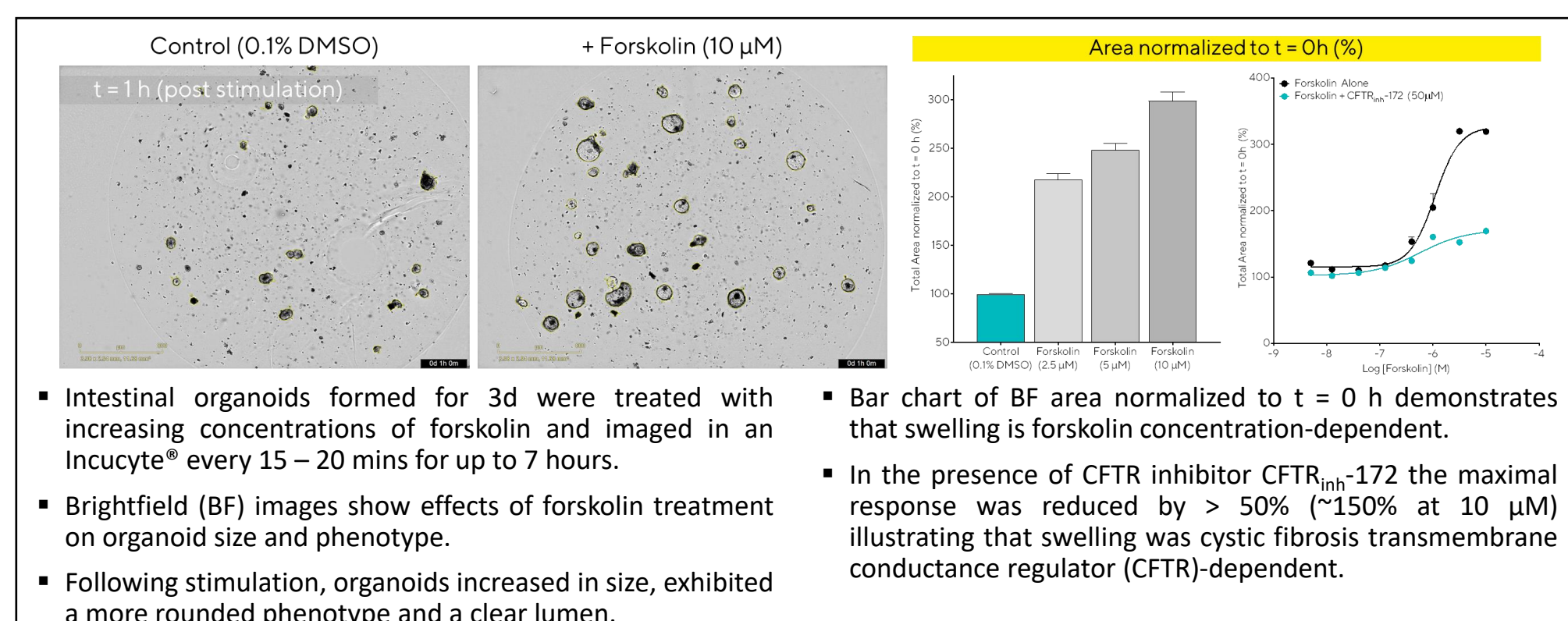
Label free quantification of organoid growth and death



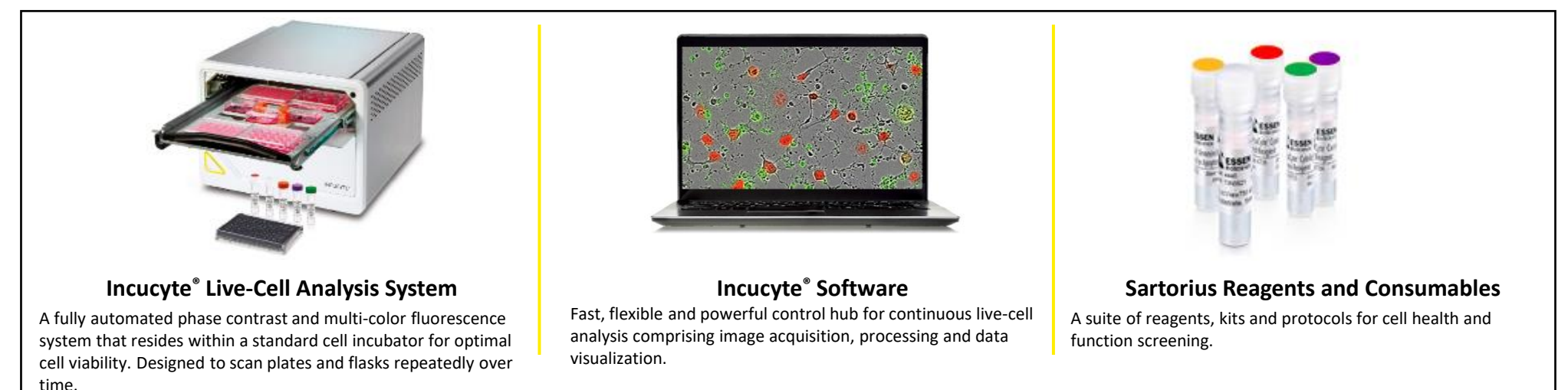
Probing mechanisms of action using morphology measurements



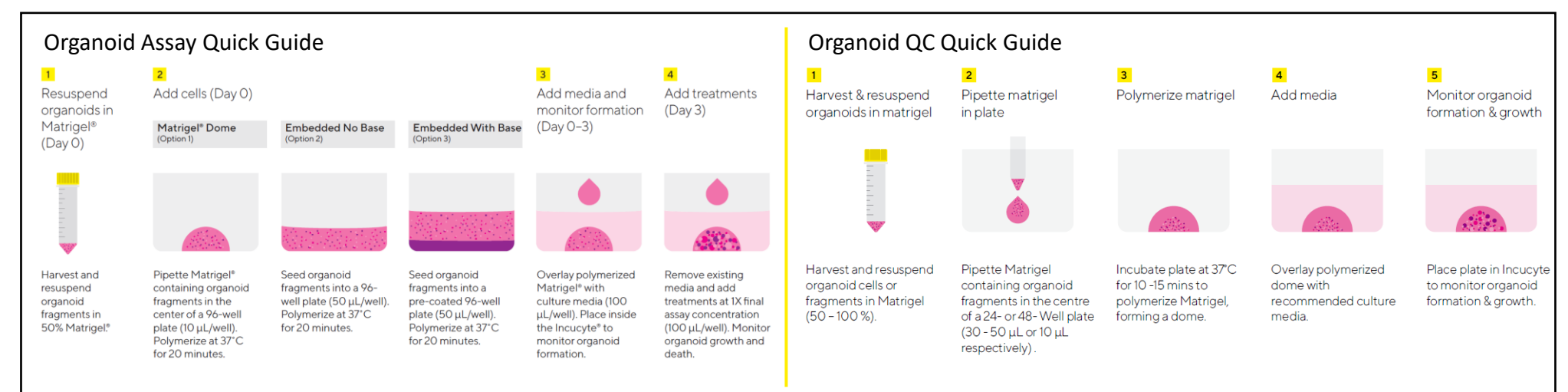
Organoid swelling in response to Forskolin stimulation



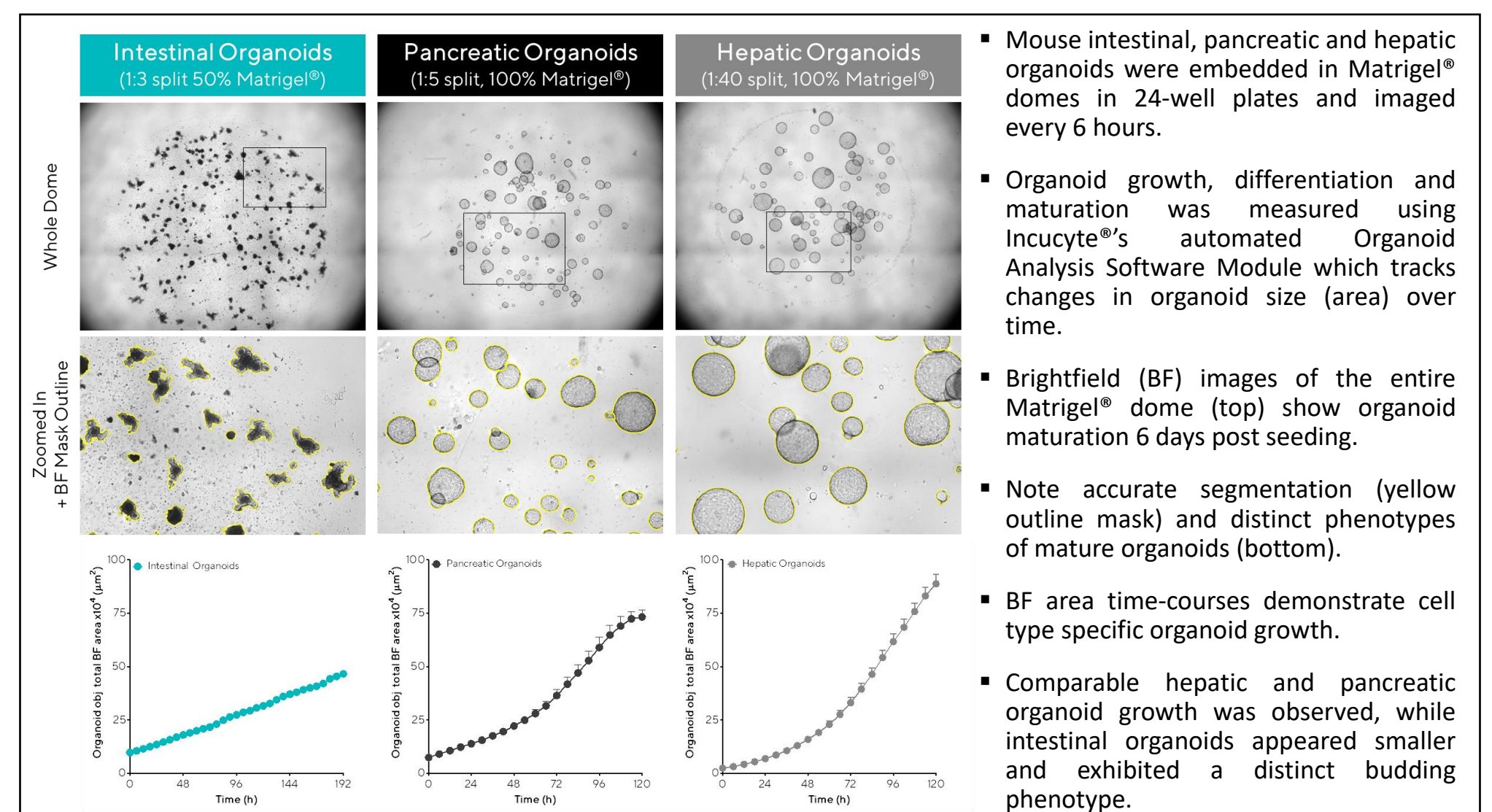
Incucyte® System for continuous live-cell analysis: Methodology



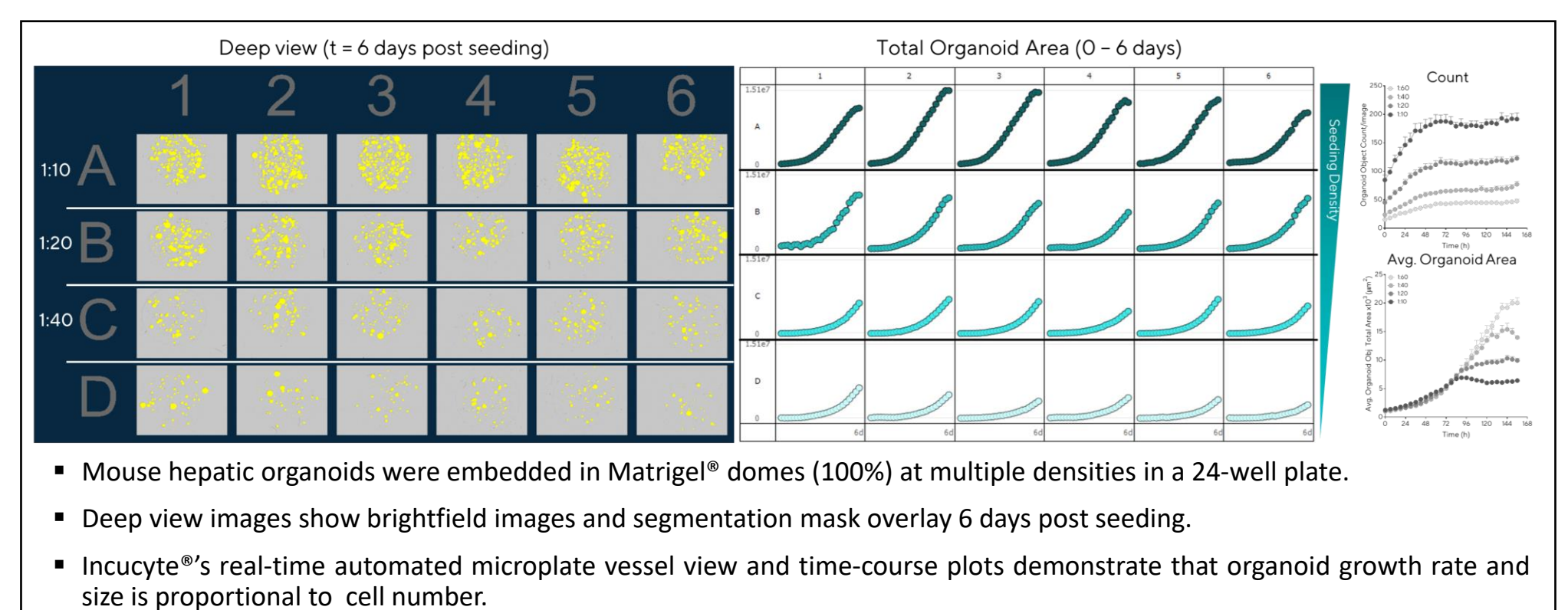
Assay Workflows



Monitoring and quantifying organoid growth in Matrigel® domes



Real-time visualisation & quantification of culture conditions



Growth & differentiation efficiency across multiple passages

