

Monitor and characterize iPSC culture and differentiation using Advanced Flow Cytometry and Live-Cell Analysis

D. Cole*, K. McBain, N. Bevan and T. Dale

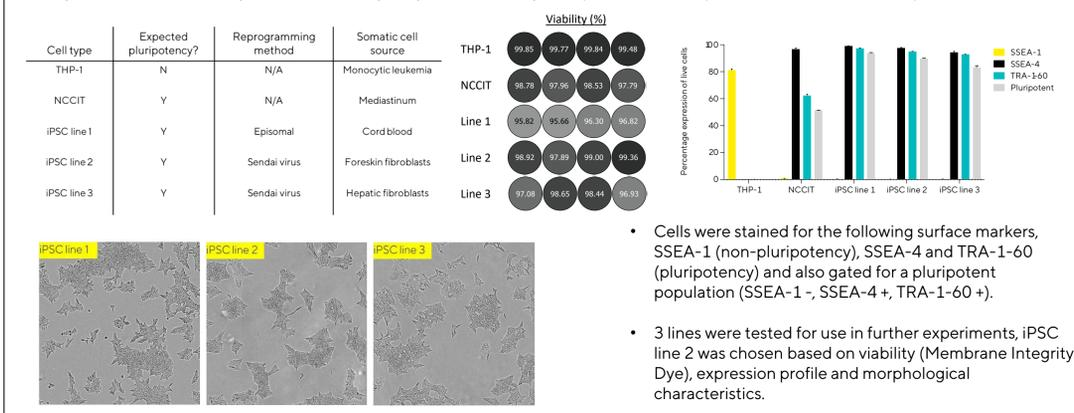
BioAnalytics, Sartorius (Essen BioScience), Royston, Hertfordshire, UK;
*Corresponding author: Daryl.Cole@sartorius.com

Introduction

- iPSCs are intrinsically valuable due to their unique characteristics and the control they afford to enable researchers to investigate early stages in cellular development.
- The major benefits of iPSCs are the variety of cell types they can be differentiated into and their capacity for infinite expansion.
- This flexibility provides opportunities for development of cell and tissue models in both 2D and 3D for pharmacological testing, cancer research, organoid modelling of tissues and neurodevelopmental biology.
- iPSCs are increasingly used in translational applications, targeting eventual clinical use via autologous cell therapies and individualized medicine approaches.
- Despite their broad application potential, iPSCs are high maintenance, expensive, and require constant monitoring to ensure they maintain pluripotency, viability, and homogeneity.
- Here we demonstrate how you can evaluate and monitor the pluripotency of iPSCs both in 2D and 3D culture systems by cell surface marker evaluation using the iQue® Advanced Flow Cytometer and morphological analysis using the Incucyte® Live-cell analysis platform.
- Long-term hepatic iPSC differentiation experiments can also be successfully monitored and validated throughout the time-course, providing a streamlined solution for differentiation quality control experiments.
- Maintenance of pluripotent phenotype can be achieved without daily media changes using Sartorius® Research Use Only (RUO) growth factors and cytokines.

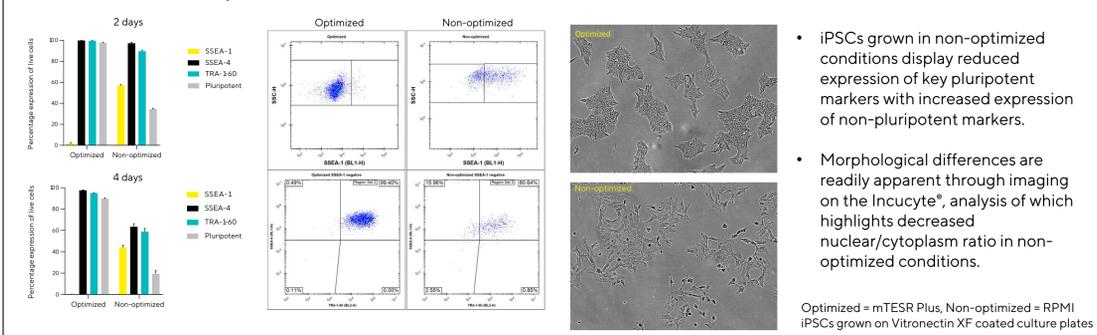
Test pluripotency

Using a combination approach, cell surface markers and morphological indicators of pluripotency can be analyzed using the iQue® Advanced Flow Cytometer and the Incucyte® Live-Cell Analysis System to identify the optimal line for expansion and downstream experiments.



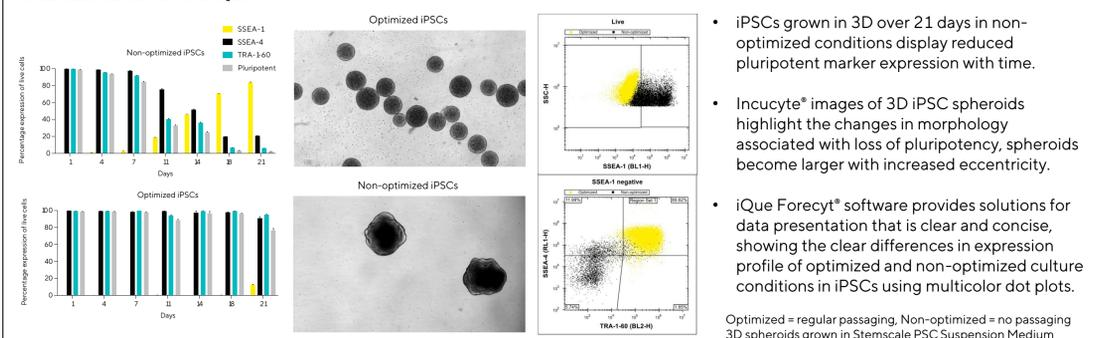
Monitor 2D growth

Culture iPSC cells in 2D and test the effects on pluripotency and morphology of different media formulations - image on the Incucyte® and stain for surface markers on the iQue®.



Monitor 3D growth

Culture iPSC cells in 3D long term in shake flasks, analyzing the effects of different culture conditions over time - image on the Incucyte® and stain for surface markers on the iQue®.

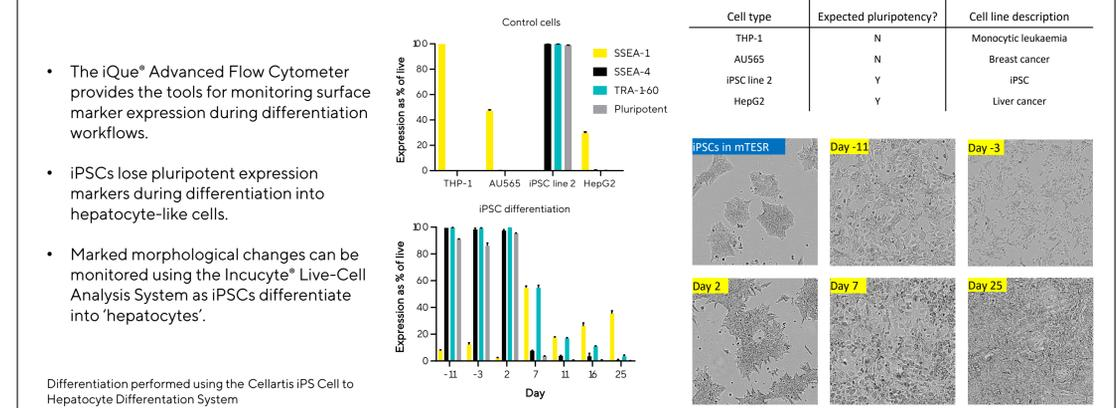


Incucyte® & iQue® Systems



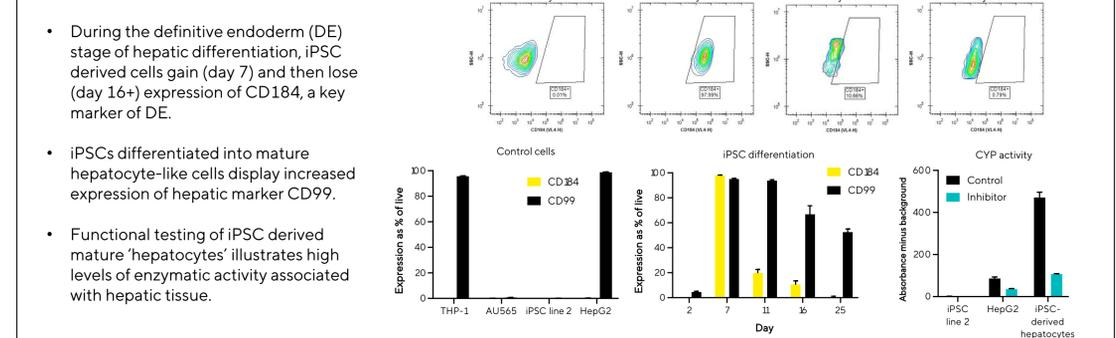
Monitor differentiation

Follow differentiation (e.g. to iPSC-derived hepatocytes) by monitoring pluripotency and target cell type marker expression using the iQue® and morphology over time using the Incucyte®.



Assess derived cell function

Use functional assays to determine whether cell has been differentiated to the desired state (e.g. luminescence based Cytochrome P450 (CYP) inhibition assay to assess hepatocyte function).



RUO growth factors and cytokines

Maintain iPSC pluripotency without the requirement for daily feeding.

