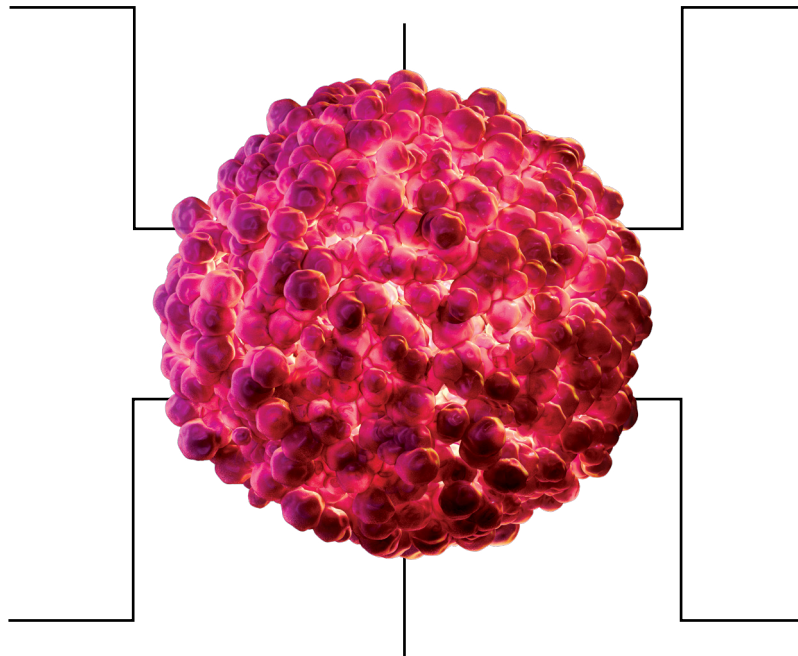


## Advanced Cell Models for Biomedical Research and Discovery

Advanced, three-dimensional (3D) cell models, such as patient-specific and iPSC-derived organoids and spheroids can uncover mechanisms of action with greater precision and detail, thereby accelerating the development of more effective treatments.

Sartorius provides solutions for automated, standardized generation, acquisition, multiplex screening and real-time, AI-driven characterization of advanced cell models.



### Accelerate

Market leading speed with cell analysis instruments and reagents



### Translate

In-depth, predictive insights that can translate into better clinical outcomes



### Simplify Workflows

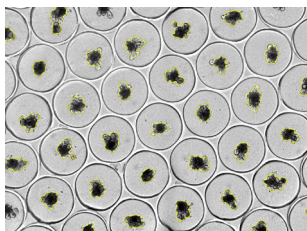
Easy assay preparation and intuitive software guides the user through each step



### Reduce Costs

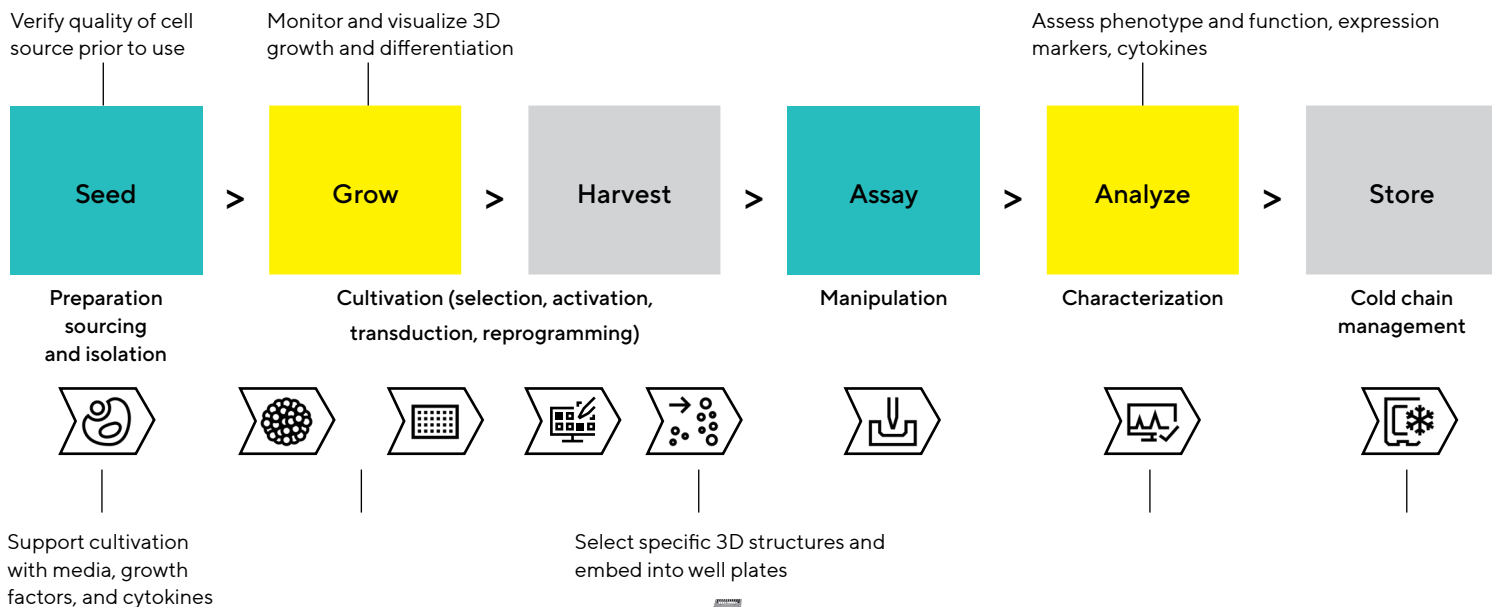
Capture unwanted effects earlier and reduce failure rates later

# 3D Cell Culture Workflow, Applications and Solutions



**Incucyte® Live-Cell Analysis System with Organoid Analysis Software** automatically locates and analyzes organoids, providing powerful insight into morphology, size, and count.

**iQue® 3 High-Throughput Cytometry Platform.** The fastest way to generate high-content phenotype and function data from small and precious 3D samples.



**CellSelector Automated Cell Selection and Retrieval Platform**

## Resources



**Poster**  
Quantify T Cell Response in 3D Tumor Spheroids Using Advanced Flow Cytometry and Live-Cell Analysis



**eBook**  
Organoid Analysis Guide



**Technical Note**  
Automated Workflows for the High-Throughput Selection and Picking of Complex 3D Structures



**Webinar**  
Maximizing the Success of 3D Cell Models for Clinical Research

### Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

### USA

Sartorius Corporation  
3874 Research Park Dr.  
Ann Arbor, MI 48108  
Phone +1 734 769 1600



For further information, visit  
[www.sartorius.com](http://www.sartorius.com)