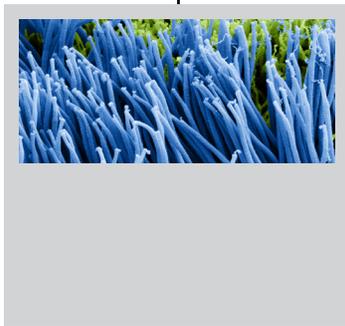
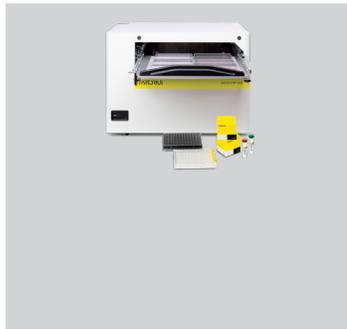


Advanced Cell Models for Biomedical Research and Discovery

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.

▼ Click any product to learn more



Accelerate

Market leading speed with cell analysis instruments and reagents



Translate

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



Simplify Workflows

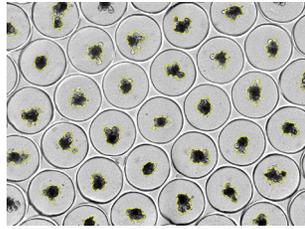
Easy assay preparation and intuitive software guides the user through each step



Reduce Costs

Capture unwanted effects earlier and reduce failure rates later

3D Cell Culture Workflow, Applications and Solutions

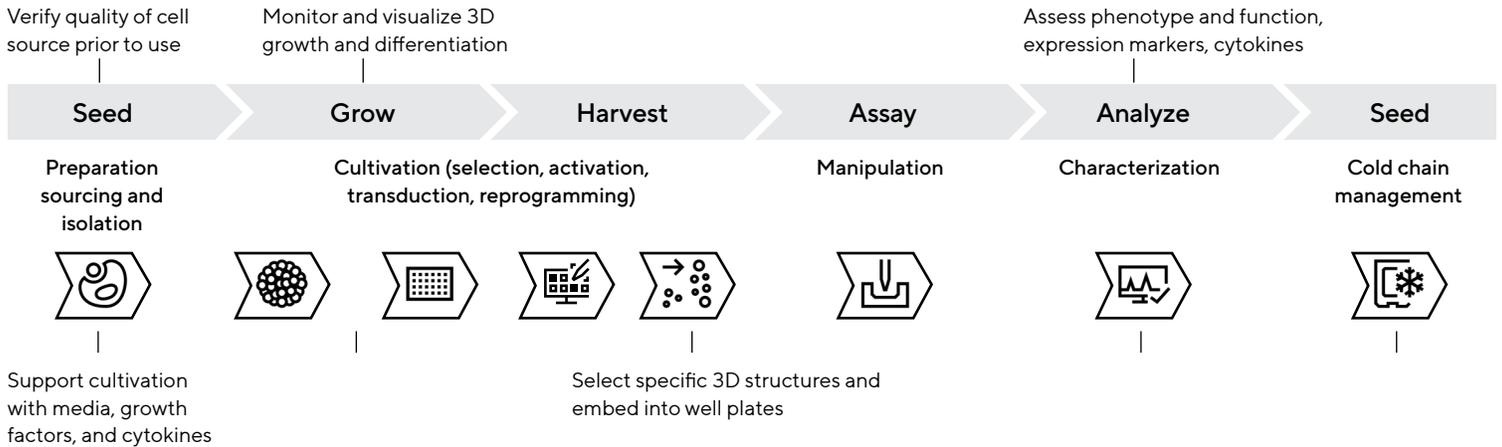


Incucyte® CX3 Live-Cell Analysis System with Organoid Analysis Software

Automatically locates and analyzes organoids, providing powerful insight into morphology, size, and count.

iQue® 5 High-Throughput Cytometry Platform

The fastest way to generate high-content phenotype and function data from small and precious 3D samples.



CellCelector Automated Cell Selection and Retrieval Platform

Isolate complex organoids from various liquid and semi-solid media.



Resources

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