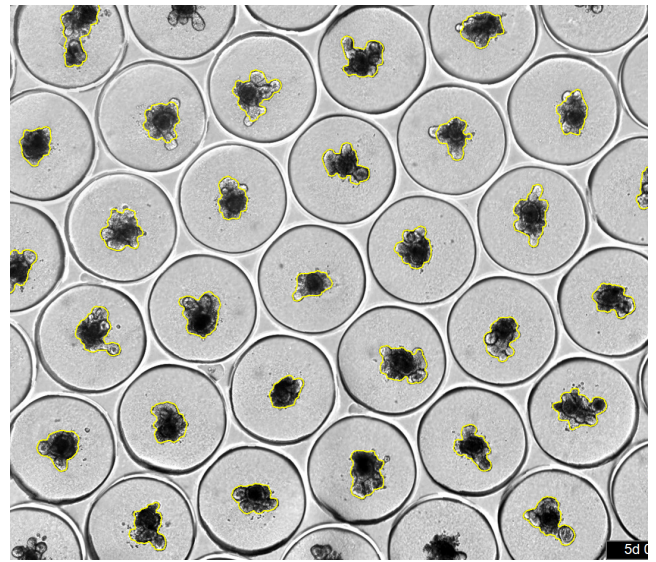


## Incucyte® Organoid Analysis Software Module

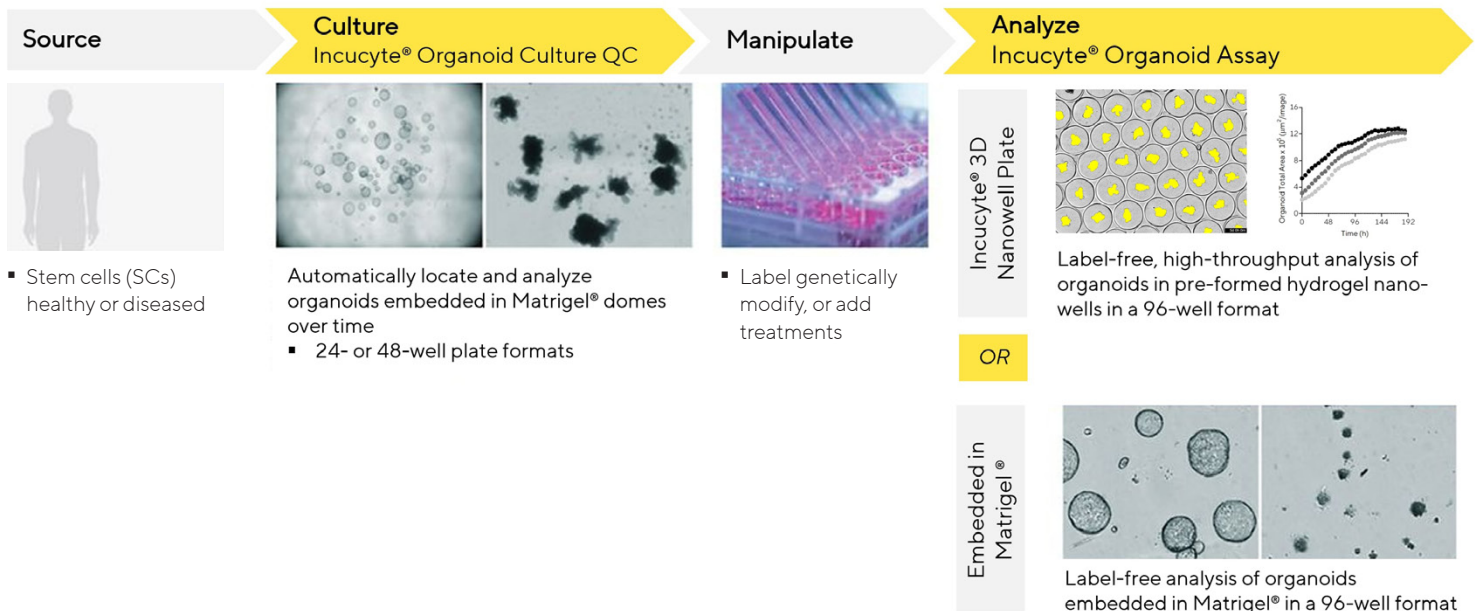
Make Informed Decisions About Complex Organoid Cultures

With the Incucyte® Organoid Analysis Software Module, You Can:

- **NEW!** Uniformly form organoids using Incucyte® 3D Nanowell Plates for high-throughput compound screening
- Automatically locate and analyze organoids embedded in Matrigel®
- Visualize differentiation, maturation (24-, 48-well plates) or death (96-well plates) of organoids
- Perform continuous, label-free analysis in psychologically relevant conditions
- Use quantitative data to support and document passaging decisions
- Access treatment effects on organoid growth in 96-well plates



Standardize your entire organoid workflow, from generation, maintenance and passaging to the final assay analysis, with image-based, label-free measurement of organoid count, size and morphology in a relevant environment using Incucyte's automated image acquisition and integrated data analysis.



# Model Complex Diseases and Make Data Driven Decisions

- **NEW!** Reproducibly form organoids and evaluate treatment effects on viability within the same plate using Incucyte® 3D Nanowell 96-well Plates, which contain a hydrogel printed array of microwells for high-throughput compound screening.
- Generate consistent and reproducible results using optimized protocols and robust image acquisition of embedded organoids
- Characterize and assess organoid maturation and morphology and probe the effects of treatments through unbiased assessment of size, count and morphology

## Ordering Information

### Product

Incucyte® Organoid Analysis Software Module

### Description

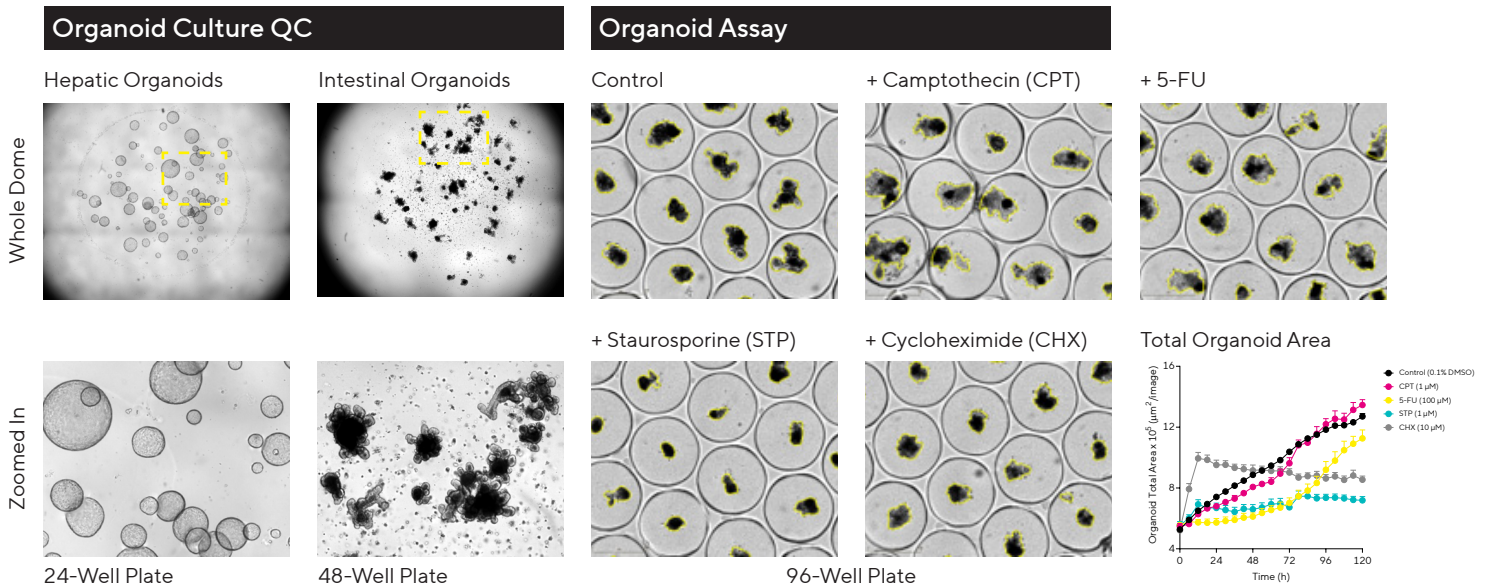
Enables label-free analysis of organoids grown and assayed in Incucyte® 3D Nanowell Plates or organoids embedded in Matrigel® in 24-, 48-, or 96-well plates

### Cat. No.

9600-0034-A00

### Instrument Compatibility

Incucyte® SX5, S3, SX1 Live-Cell Analysis Systems



Kinetically monitor and quantify organoid differentiation (Organoid Culture QC) or growth and death (Organoid Assay) undisturbed inside your incubator. Capture distinct organoid morphology, track cell death and analyze changes in size (Total Brightfield Area) using Incucyte® Organoid Analysis Software Module. Mouse hepatic and intestinal organoids were embedded in Matrigel® domes in 24- or 48-well plates respectively and imaged on the Incucyte®. In a 96-well assay format, intestinal organoids were cultured in an Incucyte® 3D Nanowell Plate and treated over 5 days with Staurosporine (1 µM, STP), Camptothecin (1 µM), Cycloheximide (10 µM), and 5-FU (100 µM). Images and time-course data of Total Organoid Area show the effects of the multiple compounds and control conditions over time.

Find out more:

[sartorius.com/incucyte-organoid](https://www.sartorius.com/incucyte-organoid)

Email [ips.opm.na@sartorius.com](mailto:ips.opm.na@sartorius.com)

North America +1 734 769 1600, ext. 3

Europe +44 1763 227400

Asia Pacific +65 6872 3966

Specifications subject to change without notice. ©2024. All rights reserved. All names of Sartorius products are registered trademarks and the property of Sartorius AG and/or one of its affiliated companies. Incucyte-Organoid-Analysis-Software-Module-1-Sheet-en-L-Sartorius. Version 07 | 2024