

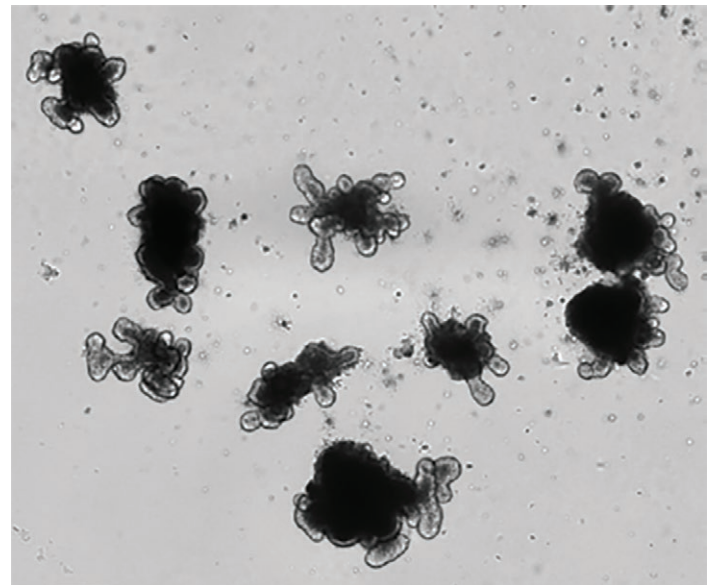
Incucyte® Organoid Analysis Software Module

Make Informed Decisions About Complex Organoid Cultures

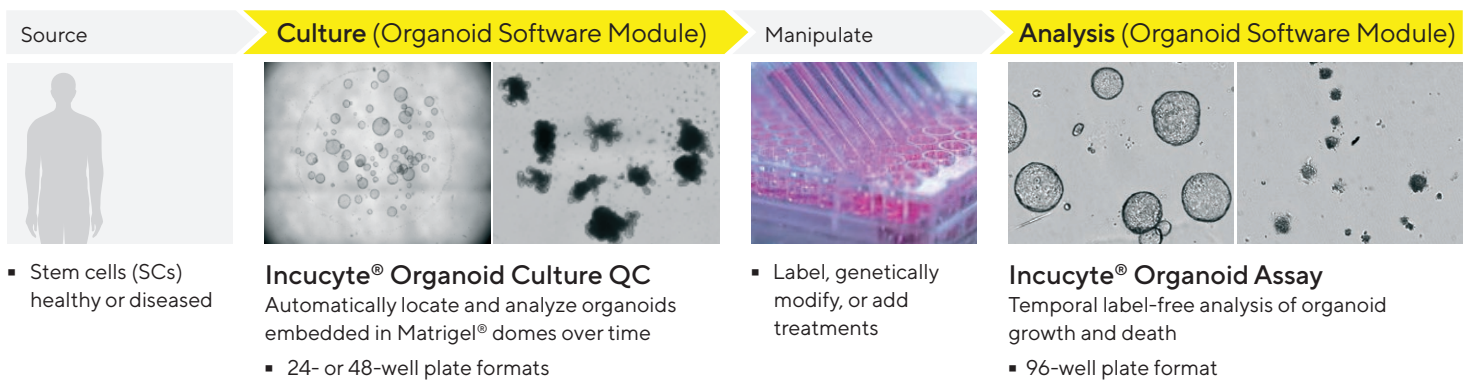
With the Incucyte® Organoid Analysis Software Module, You Can:

- Locate and analyze Matrigel® embedded organoids automatically
- Perform continuous label-free analysis in physiologically relevant conditions
- Characterize differentiation and maturation of organoids in Matrigel® domes in 24- or 48-well plates
- 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.



Incucyte® Organoid Workflow



Model complex diseases and make data driven decisions using Incucyte's automated image acquisition and integrated data analysis.

- 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

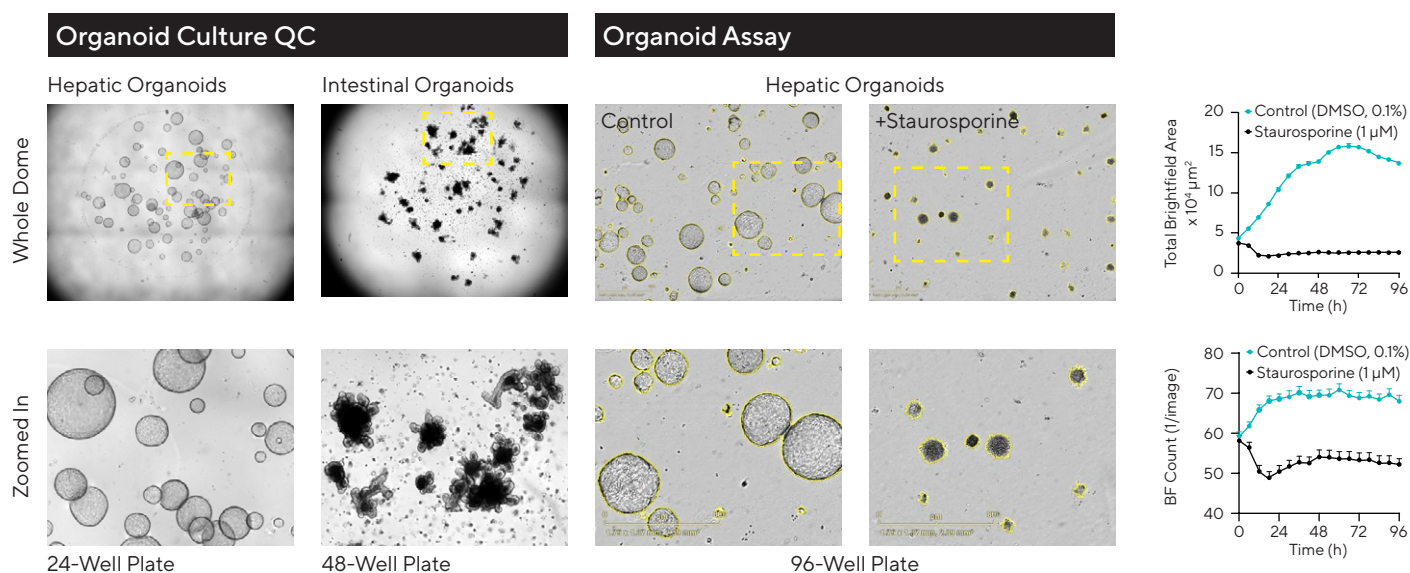
Enables label-free analysis of organoids embedded in Matrigel® in 24-, 48- or 96-well plates

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, hepatic organoids were embedded in 50% Matrigel® and treated with 1 μM Staurosporine over 4 days. Time-course data of organoid growth demonstrated that Staurosporine induced organoid cell death and reduced size of hepatic organoids.

Find out more:

www.sartorius.com/incucyte

Email orders.US07@sartorius.com

North America +1 734 769 1600, ext. 3

Europe +44 7515 947101

Asia Pacific +81 3 6478 5202

Specifications subject to change without notice. © 2021. All rights reserved. Incucyte, Essen BioScience, and all names of Essen BioScience products are registered trademarks and the property of Essen BioScience unless otherwise specified. Essen BioScience is a Sartorius Company. Publication No. 8000-0772-A00 Version 03 | 2021