

Introducing Incucyte® Proliferation Assays for Live-Cell Analysis

With Incucyte® Proliferation Assays and AI Label-Free Cell Segmentation, You Can:

- Identify and segment cells reliably and label-free
- Conduct real-time, robust image analysis for high-throughput compound testing
- Quantify proliferation and multiplex with other cell health dyes—in mono- or co-cultures

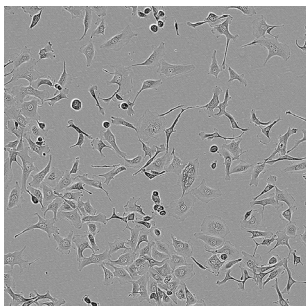
Measure Cell Proliferation Robustly with Integrated and Automated Analysis

The Incucyte® Base Analysis Software employs both Artificial Intelligence (AI) and Classic Confluence Analysis with fully automated acquisition and identification of non-adherent or adherent cells to measure confluence in a physiologically relevant environment.

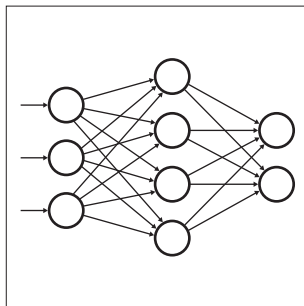
An AI-driven confluence analysis workflow offers accurate, objective analysis of cells and is trained to adapt to a wide range of cell types and morphologies while requiring less user input.



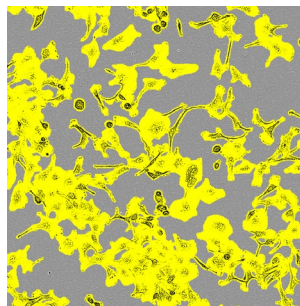
Image Acquisition



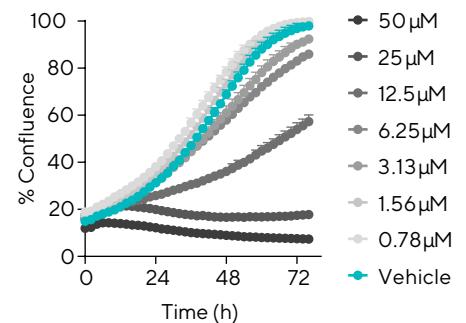
AI-Neural Network



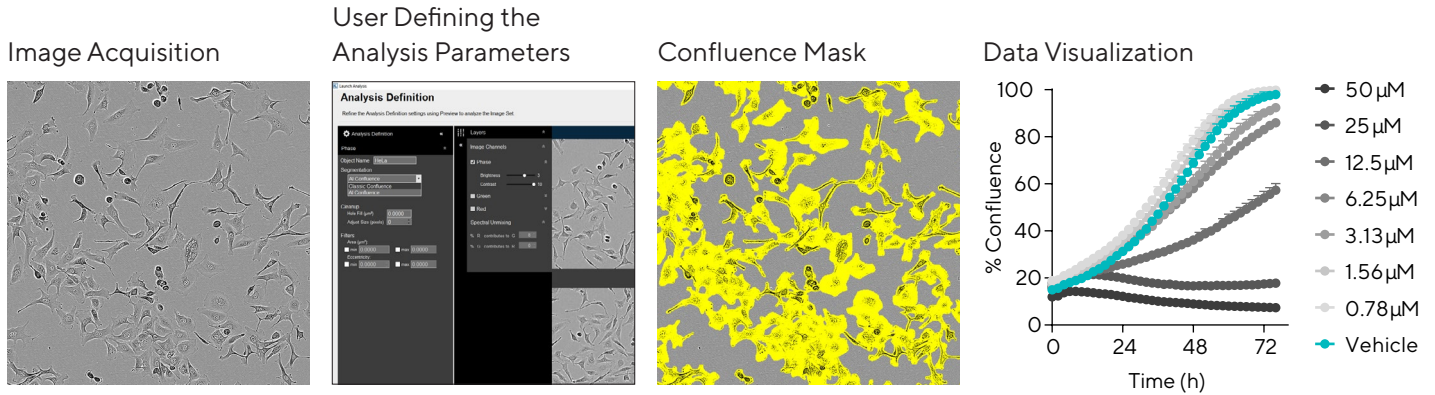
Confluence Mask



Data Visualization

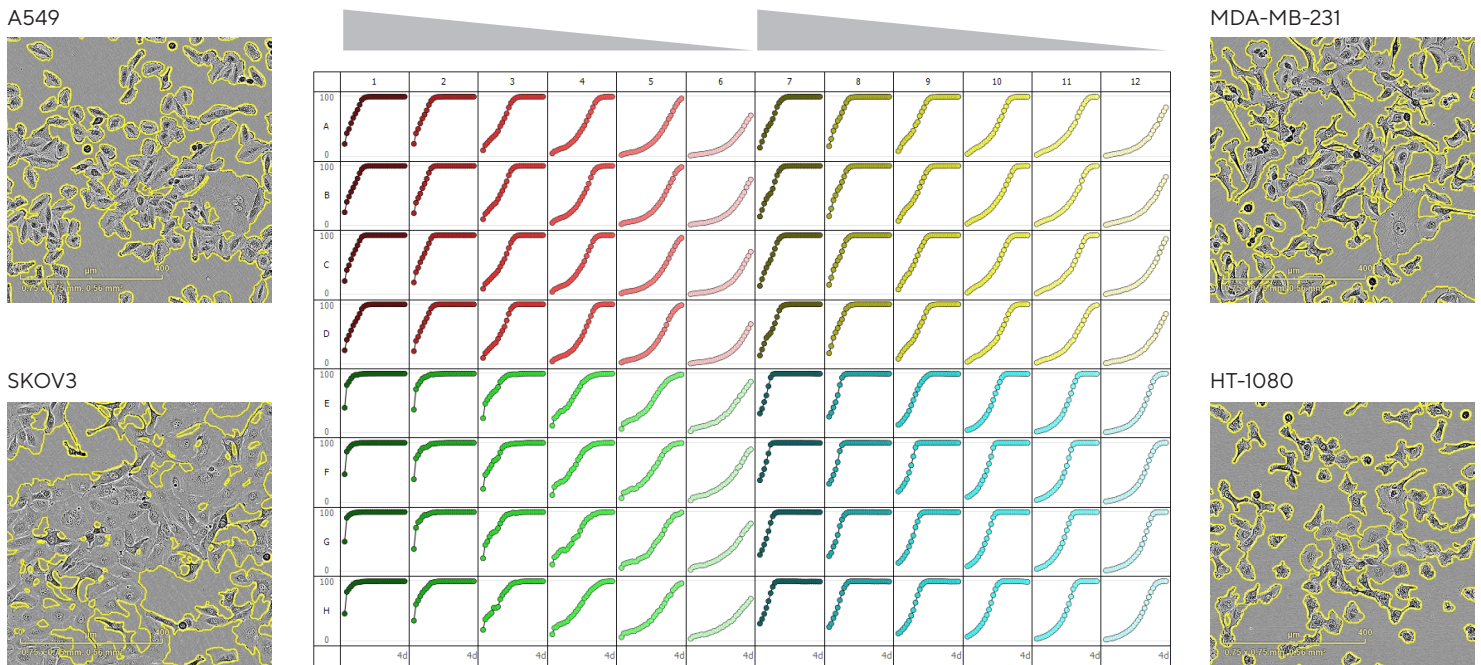


A classic confluence analysis workflow incorporates user-defined segmentation based on experimental conditions. Fluorescence image acquisition and analysis of cell health or viability assays are also available with Incucyte® Base Analysis Software, using classic analysis definitions.



Kinetically Monitor and Conduct Robust Analysis of Cell Confluence with High-Throughput

Robust Segmentation of Cell Confluence Using AI Confluence Analysis



Note. A549, MDA-MB-231, SKOV3 and HT-1080 cells were seeded into a 96-well plate at a range of densities (0.5–10 K cells/well). High definition (HD) phase-contrast images were acquired every 2 hours and cell proliferation was measured over 4 days. Yellow outline indicates AI Confluence segmentation and plate view reports % Confluence over 4 days.

Germany:
Sartorius Lab Instruments
Phone +49 551 308 0

USA:
Sartorius Corporation
Phone +1 631 254 4249
Toll-free +1 800 635 2906

Learn more: www.sartorius.com/incucyte

For questions, email: orders.US07@sartorius.com