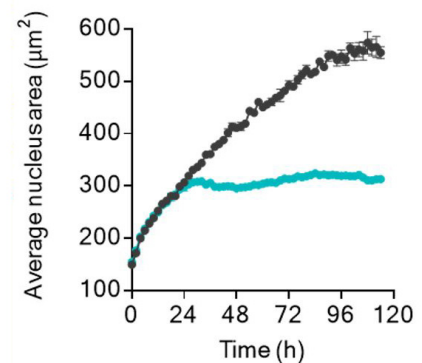
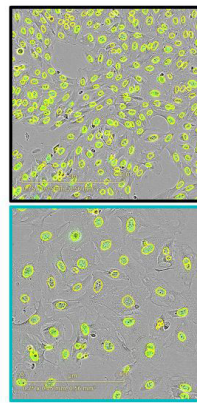


Introducing the Incucyte® AI Nuclei Detection Analysis Software Module for Live-Cell Analysis

With Incucyte® AI Nuclei Detection Analysis Software Module You Can:

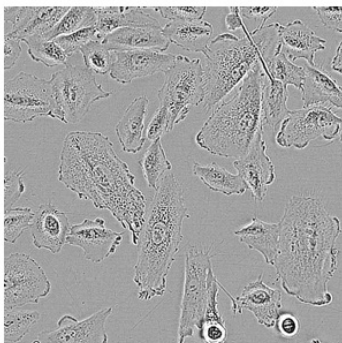
- Identify and segment live cell nuclei in HD phase contrast images—no labels or dyes required
- Extract detailed nuclear metrics such as area, eccentricity, and texture for advanced phenotypic profiling
- Enable robust, automated image analysis across diverse tumor cell lines without time-consuming parameter tuning
- Perform high-throughput, label-free analysis of cell proliferation to high confluence



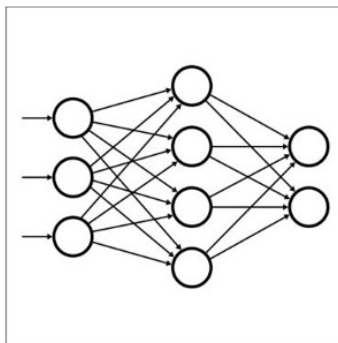
Quantify Nuclei with Integrated, Automated, and Label-Free Analysis That Is Accurate and Repeatable

The Incucyte® AI Nuclei Detection Analysis Software Module is an advanced, all-in-one image analysis tool that segments individual nuclei in HD phase-contrast images—without the need for fluorescent labels. It delivers high-accuracy nuclear metrics such as area, eccentricity, and texture, enabling objective, reproducible results with minimal user input and reduced analysis time.

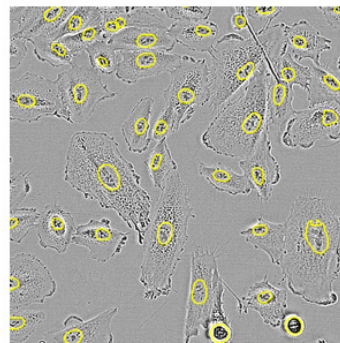
Image Acquisition



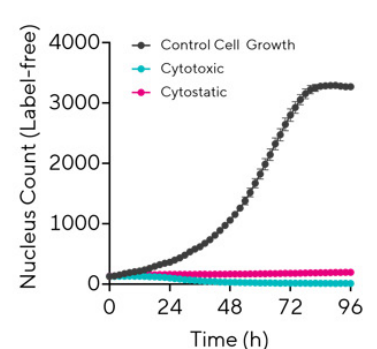
AI Nuclei Segmentation



Nuclei Mask

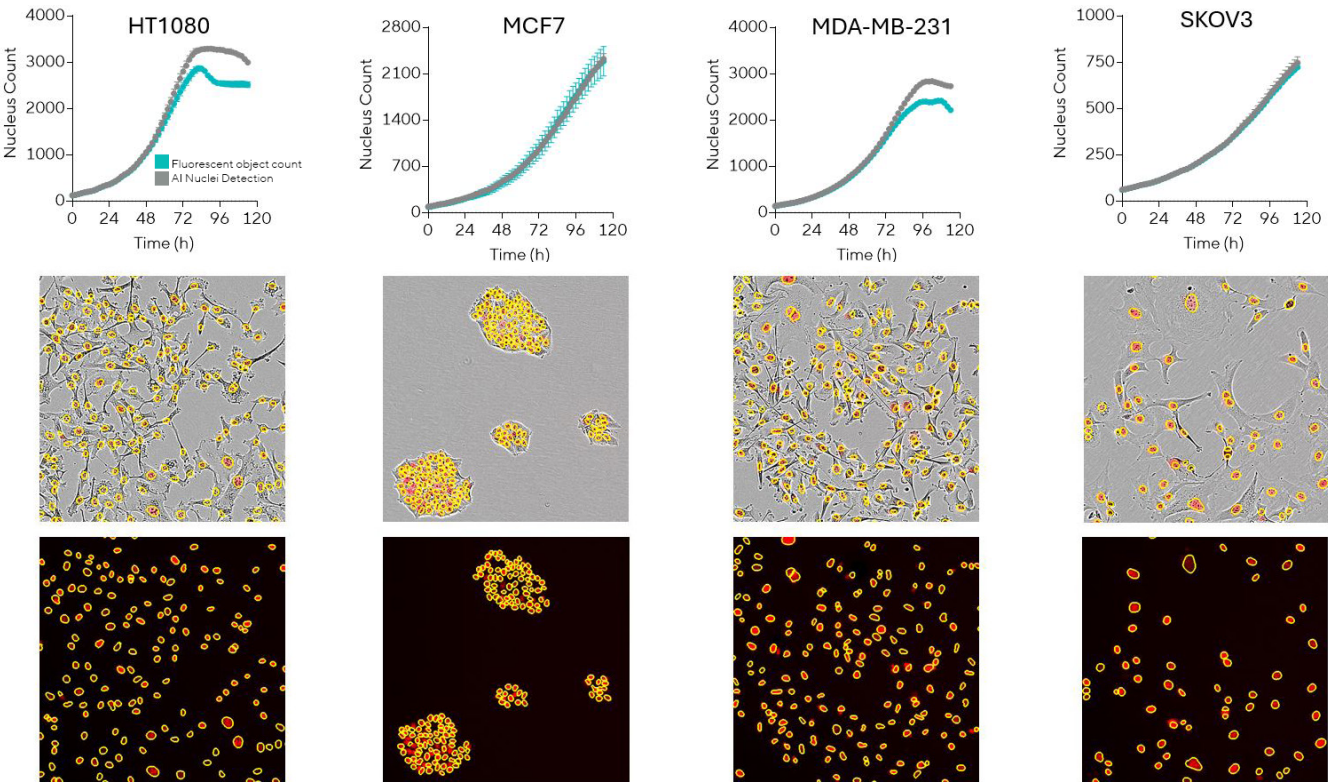


Data Visualization



AI Nuclei Detection uses a pre-trained neural network to segment the nuclei of live, adherent cells in Phase Contrast images. Nuclei masks are displayed and provide robust quantification of nuclei count and shape over time.

Monitor and Conduct Label-Free, Unbiased Analysis of Cell Proliferation



Label-free analysis of cell proliferation. Adherent cell lines HT1080, MCF7, MDA-MB-231 and SKOV3 expressing NucLight Red were seeded at low density and proliferation was measured over 5 days using either AI Nuclei Detection (grey) or fluorescent object count in Basic Analyzer (teal). Nucleus count is similar with AI Nuclei Detection more accurately identifying cells at high confluence. Phase contrast images highlight diverse morphology across cell lines; fluorescence images with AI Nuclei segmentation mask (yellow outline) indicate accurate label-free identification of the nuclei.

The Incucyte® AI Nuclei Detection Analysis Software Module is available to purchase for Incucyte® CX3, SX5, S3, and SX1 systems and requires software version 2025B or higher.

Ordering Information

Product	Description	Cat. No.	Instrument Compatibility
Incucyte® AI Nuclei Detection Analysis Software Module	Enables label-free, AI-driven detection and counting of live cell nuclei	BA-04897	Incucyte® CX3, SX5, S3 and SX1 Live-Cell Analysis System


Germany

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

Germany

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