

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

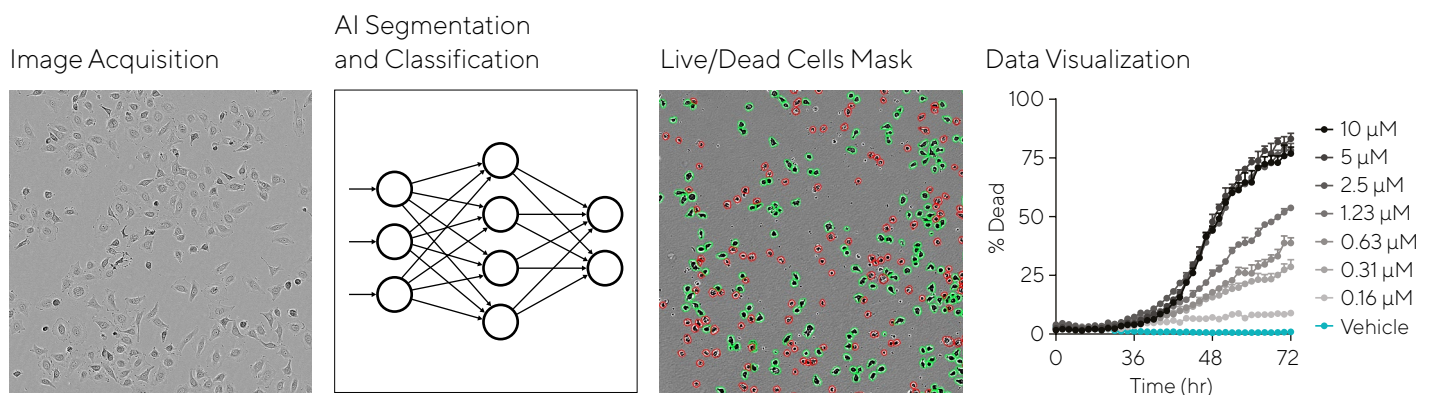
With Incucyte® AI Cell Health Analysis Software Module and AI Label-Free Viability Assay, You Can:

- Identify and segment cells to quantify live vs dead without the need for labels
- Combine with additional fluorescent indicators of cell biology for deeper insight
- Conduct real-time, robust image analysis for unbiased high-throughput compound testing
- Maximize sophisticated AI image processing algorithms with a powerful GPU co-processor update to your Incucyte® controller



Measure Cell Viability Accurately With Integrated and Automated Analysis That Is Objective and Replicable

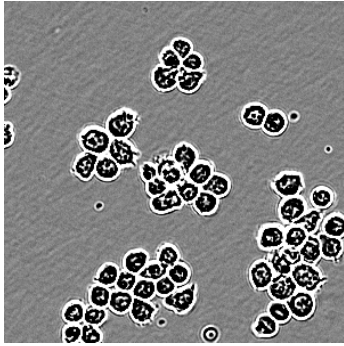
The Incucyte® AI Cell Health Analysis Software Module is an all-in-one analysis tool that segments cells and classifies them as live or dead. It allows users to perform cell health analyses without the need for fluorescent labels, generating accurate and objective data with reduced time requirements and cost.



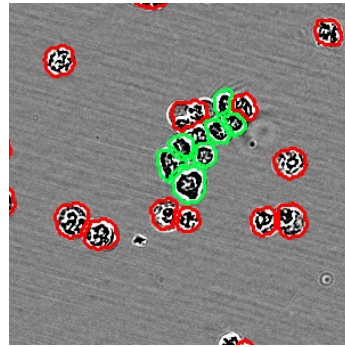
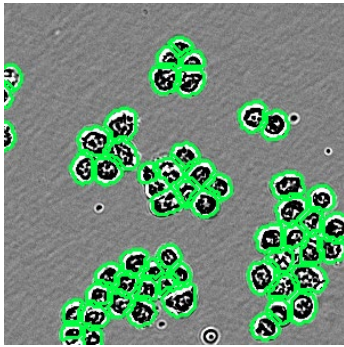
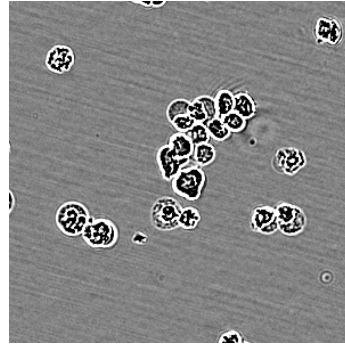
Images are captured at specified intervals, then processed using our purpose-built, validated AI algorithms to segment and classify cells, producing AI cell segmentation and live/dead classification masks. The resulting data can be visualized over a time course and EC_{50} values calculated.

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

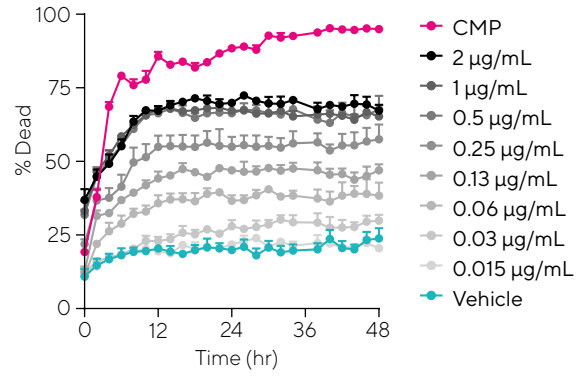
Ramos Untreated
(48 hr)



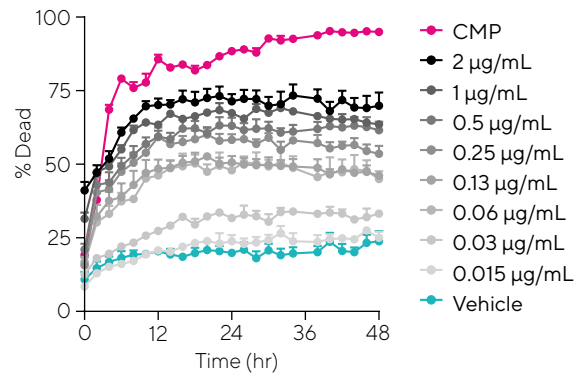
Ramos + Rituximab
(2 µg/mL, 48 hr)



Rituximab



Truxima®



Label-free analysis of cell viability. Ramos B-cell Lymphoma cell line was treated with increasing concentrations of anti-CD20 antibody Rituximab and the biosimilar, Truxima® and cell death was quantified using Incucyte® AI Cell Health Analysis. Biosimilar mAbs Rituximab and Truxima® induce death in Ramos cells with similar efficacy. Untreated cells are healthy and classed as Live (green segmentation) while treated cells show partial cell death (red classification).

The Incucyte® AI Cell Health Analysis Software Module is available to purchase for all S-series instruments and requires a GPU co-processor installed as a drop-in hardware upgrade to the controller and 2022B software version.

Ordering Information

Product	Description	Cat. No.	Instrument Compatibility
Incucyte® AI Cell Health Analysis Software Module	Enables label-free, AI-driven classification of live dead cells and cells counts of adherent and non-adherent cells	BA-04871	Incucyte® SX5, S3 and SX1 Live-Cell Analysis System
Incucyte® GPU (Graphics Processing Unit)	Provides enhanced processing speed, software sold separately	BA-04870	Incucyte® SX5, S3 and SX1 Live-Cell Analysis System

Germany:

Sartorius Lab Instruments
Phone +49 551 308 0

USA:

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