SVISCISVAS

Incucyte[®] Label-Free Analysis Software Modules for Live-Cell Analysis

With Incucyte® Label-Free Analysis Software Modules, You Can:

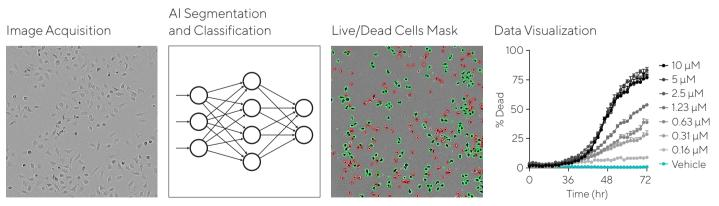
- Identify and segment adherent or non-adherent cells to quantify confluence
- Perform direct, label-free cell counts utilize fluorescent markers for deeper biological insight
- Classify cells in real-time with machine-learning image analysis for unbiased morphological assessment
- Maximize sophisticated Artificial Intelligence (AI)-driven image processing with Live/Dead classification of cells and subsequent subclassification capabilities



New Software Module Spotlight:

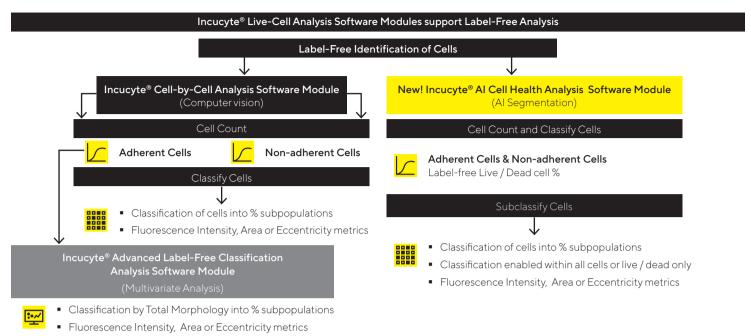
Incucyte® AI Cell Health Analysis Software Module

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.



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

Incucyte[®] Live-Cell Analysis Software Modules support Label-Free Analysis, simplifying complex analyses with a guided user interface and purpose-built analysis tools. These combine to provide a powerful workflow that supports your entire research team making the process of acquiring, viewing, analyzing, and sharing of images of living cells easier than ever before using Classic or Al-driven Analysis.



Ordering Information

Product	Description	Cat. No.	Instrument Compatibility
Incucyte® AI Cell Health Analysis Software Module	Enables label-free, Al-driven classification of live/dead cells and cells counts on adherent and non-adherent cells. Requires BA-04870 and 2022B or newer software version.	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
Incucyte® Advanced Label-Free Classification Analysis Software Module	Monitor adherent cell morphology changes to determine live/dead cell counts or classify cells based on morphology via label-free multivariate analysis. Module requires Incucyte [®] Cell-By-Cell Analysis Software Module (Cat. No. 9600-0031).	BA-04867	Incucyte® SX5, S3 and SX1 Live-Cell Analysis System
Incucyte® Cell-by-Cell Analysis Software Module	Perform label-free cell counts on adherent and non-adherent cells and subsequent cell-by-cell classification based on shape, size or fluorescence intensity to quantify dynamic changes in subpopulations within a mixed culture.	9000-0031	Incucyte [®] SX5, S3 and SX1 Live-Cell Analysis System

The Incucyte® AI Confluence Analysis included in Incucyte® Base Software with Incucyte® Software v2022A or newer software version. The Incucyte® AI Cell Health Analysis Software Module is available to purchase for all S-series instruments and requires both a GPU co-processor installed as a drop-in hardware upgrade to the controller and 2022B or newer software version.

North America Sartorius Corporation Phone +1734 769 1600 Email: Ips.opm.na@sartorius.com **Europe** Sartorius UK Ltd. Phone +44 1763 227400 Email: orderhandling.lps.ne@sartorius.com Asia Pacific Sartorius Singapore Pte. Ltd. Phone : + 65 6872 3966 Email : sartoriusap@sartorius.com

Learn more: Live Cell Analysis Software | Sartorius

For questions, email: lps.opm.na@sartorius.com