

Incucyte® Live-Cell Analysis System All-In-One Warranty



Ensure Peace of Mind With Our All-In-One-Warranty

Your Incucyte® Live-Cell Analysis System is continuously powered on and operates in a challenging incubator environment. Get the protection you need for your Incucyte® by purchasing an additional 1, 2 or 3 years all-in-one warranty.

What's Included:

Priority Response Time Over Non-Warranty Customers

Ensure rapid response from our global

support team in the event of any issue. Our all-in-one warranty covers costs including remote diagnostics and troubleshooting, all travel costs and labor, on-site labor and any required spare parts for preventative maintenance and or break fix service.

Maintenance and Inspection

Yearly preventative maintenance and inspection visits can be arranged on request to keep your Incucyte® live-cell analysis system operating optimally. Our 17-point preventative maintenance plan includes:

- A comprehensive system health check to identify and resolve issues before they impact instrument performance.

- Thorough inspection of the camera assembly, pulleys and belts.
- Review of image storage capacity.
- Full ethanol wipe down of interior components and inspection for mold.

Free Software Updates

Receive free base software updates for your Incucyte® live-cell analysis system as they become available, giving you access to new instrument capabilities, system features and the latest vessel databases.

The warranty product can be applied to the Incucyte® SX5, S3, SX1 and ZOOM systems.

Sales and Service Contacts

Save 30% off your all-in-one warranty when you purchase with an instrument.

To purchase an all-in-one warranty, simply contact your local sales representative, or call our offices at

North America

Phone: +1 734 769 1600 Ext. 3

Europe

Phone +44 (0) 1763 227400

Or e-mail us at:

incucyte.support@sartorius.com