

Babraham, UK | Göttingen, Germany | April 7, 2025

Sartorius partnership with bio-incubator LiveLabs empowers early-stage startups with cutting-edge solutions

- Sartorius partners with LiveLabs to provide early-stage life science companies with advanced lab solutions
- Agreement includes the Incucyte® SX5 Platform and leading essential lab technologies
- Partnership streamlines workflows and delivers high-quality data, accelerating growth and market readiness of new ventures

The life science group Sartorius announces a partnership with the bio-incubator LiveLabs at Babraham Research Campus to provide early-stage life science companies with cutting-edge lab solutions that accelerate their ideas to market. This is part of Sartorius' ongoing mission to support scientists by simplifying lab workflows and helping bring life-changing therapeutics to patients faster.

LiveLabs currently offers 2,400 sq ft of fully equipped R&D space for new ventures to nurture and grow their bioscience research, with plans for future expansion. As part of this new partnership, Sartorius is providing LiveLabs with a fully equipped Incucyte® SX5 Platform for high-throughput live-cell analysis, including all software modules. The comprehensive agreement also includes Sartorius Lab Essential products, featuring the leading Picus® 2 Electronic Pipette and Cubis® II Balance.

"Limited funding during the startup phase often leads to bottlenecks at a time when there is immense pressure to generate high quality data and achieve milestones," says Debbie Brown, Head of Life Science Sales EMEA at Sartorius. "Our partnership with LiveLabs empowers companies at this crucial stage by providing state-of-the-art solutions that simplify workflows and deliver reproducible, physiologically relevant data, accelerating the journey from bench to market."

The Incucyte® platform is a powerful asset in cell culture workflows involving sensitive or expensive cell types due to its non-invasive imaging and kinetic monitoring of live cells. Its high-throughput capability is unique in the market and valuable for startup labs aiming to maximize their data generation potential. Combined with Sartorius' lab essentials, this portfolio supports and fosters the growth of new companies in LiveLabs.

"It's a pleasure to work with the Sartorius team to support the LiveLabs community. Sartorius equipment has significantly improved LiveLabs' capabilities, plus the technical support and training the team offers are fantastic," says Ruth Campbell, LiveLabs Manager.

“We are thrilled to welcome Sartorius as a new sponsor of LiveLabs, our shared lab space.” adds Jenny Hirst, Senior Lead in Science and Entrepreneurship at Babraham Research Campus. “Their support comes at a pivotal moment in the expansion of our shared lab space, enabling us to grow the number of life science startup companies that we support. Together, we are fostering innovation and driving forward the future of scientific discovery.”

Sartorius is fostering innovation and entrepreneurship in the life sciences sector through a variety of initiatives. The primary goal is to support startups and researchers transitioning from academic settings to entrepreneurial ventures by offering resources such as mentoring programs, networking opportunities, and access to equipment. Among other activities, Sartorius is an initiator of the Life Science Factory, collaborates with the BioLabs network, and is a partner of Young Entrepreneurs in Science (YES).

More information about LiveLabs:

<https://www.babraham.com/livelabs/>

A profile of Sartorius

Sartorius is a leading international partner of life sciences research and the biopharmaceutical industry. With innovative laboratory instruments and consumables, the Group’s Lab Products & Services Division focuses on laboratories performing research and quality control at pharmaceutical and biopharmaceutical companies as well as academic research institutes. The Bioprocess Solutions Division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications, vaccines, and cell and gene therapies safely, rapidly, and sustainably. The company, based in Göttingen, Germany, has a strong global reach with around 60 production and sales sites worldwide. Sartorius regularly expands its portfolio through the acquisition of complementary technologies. In 2024, the company generated sales revenue of around 3.4 billion euros. Around 13,500 employees are working for customers around the globe.

A profile of Babraham Research Campus

Babraham Research Campus is distinct in its co-location of 60 bioscience companies with the Babraham Institute, a world-renowned research organization which receives strategic funding from the Biotechnology and Biological Sciences Research Council (BBSRC). The aim of the Campus is to support UK bioscience through academic research, but also with facilities and capabilities for early-stage and growing commercial organizations. The Campus provides companies laboratory and office space, networking and collaboration opportunities, together with access to outstanding scientific facilities in an ideal geographical location at the core of the Cambridge cluster.

Visit our [Newsroom](#) or follow us on [LinkedIn](#).

Contact

Lisa Dabney

Head of Marketing Programs

Lab Products & Services Division

+1 734-474-7117

lisa.dabney@sartorius.com