



## Press Release

# Sartorius and the German Research Center for Artificial Intelligence launch joint research laboratory

- **The Sartorius AI Lab (SAIL) conducts research in the use of artificial intelligence in Sartorius platform solutions for the biopharmaceutical industry**

Kaiserslautern | Göttingen, December 19, 2019 - Sartorius, a leading international partner of life science research and the biopharmaceutical industry, and the German Research Center for Artificial Intelligence (DFKI) have established the Sartorius AI Lab (SAIL) research laboratory. On the DFKI campus in Kaiserslautern, the use of artificial intelligence (AI) in Sartorius products and platform solutions is being experimentally tested and further developed at SAIL.

"So far, advanced methods of data analysis have been used only to a very limited extent in the biopharmaceutical industry, in drug research and in production. This is one of the reasons that development timelines and costs for medical drugs are ever increasing. Better use of data through AI is a particularly promising approach for making significant progress in this area. Many customers are already using the powerful software we have been offering for two and a half years through our Data Analytics unit for specific applications in biopharmaceutical production. We aim to strongly expand this business activity and are therefore glad to have found DFKI, one of the most important international centers of excellence for AI and deep learning with more than 1,000 scientists, as strong partners to join us in developing such solutions," said CEO Dr. Joachim Kreuzburg.

SAIL is assigned to the DFKI research department of Smart Data & Knowledge Services headed by Professor Andreas Dengel and uses DFKI's deep learning hardware and expertise, which is unique throughout Europe. As a protected data room and independent data laboratory, SAIL is also open to Sartorius partners and customers as part of cooperation agreements, and the first projects have already been initiated. In addition, Sartorius and DFKI will use SAIL for training purposes, and DFKI employees will also be able to further their education in life science applications at Sartorius research sites.

Sartorius and DFKI intend to develop and employ machine learning and image and pattern recognition processes, among other technologies, for life science applications. For example, their researchers are working on new deep learning algorithms and methods for image recognition of cells and organoids, analysis and modeling of biological systems and for simulation and optimization of biopharmaceutical production processes. Sartorius provides extensive data for this purpose and sends specialists from its Corporate Research and Product Development units to SAIL.



## Press Release

Prof. Oscar-Werner Reif, Chief Technology Officer at Sartorius, emphasized: "Within the framework of this cooperation, we are creating the foundations for developing digital tools and methods based on AI. The results obtained will be incorporated into the further development of Sartorius products so that our application solutions become more and more adaptive and powerful. Our aim is to use better methods of data analysis and the increasing computer capacities to map and simulate the development and production of biopharmaceuticals in computers in the future. As a result, development times and costs for new therapies will improve dramatically by an accelerated timeline from idea to patient. In addition, we are pleased to offer Sartorius employees and DFKI students and scientists the opportunity to exchange ideas and gain further qualifications in AI and life science applications."

Prof. Andreas Dengel, DFKI site manager in Kaiserslautern, said: "The DFKI transfer labs are developing new service modules with artificial intelligence for various markets and target groups. For two years, we have been using our AI hardware equipment, which is unique in Europe, for application-oriented research we conduct together with renowned companies such as Allianz, Continental, Hitachi and IAV. I am very pleased that we are entering the life science tools market as a partner with the global player Sartorius."

Follow Sartorius on [Twitter](#) @Sartorius\_Group and on [LinkedIn](#).

### A profile of Sartorius

The Sartorius Group is a leading international partner of biopharmaceutical research and the industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services Division concentrates on serving the needs of laboratories performing research and quality control at pharma and biopharma companies and those of academic research institutes. The Bioprocess Solutions Division with its broad product portfolio focusing on single-use solutions helps customers to manufacture biotech medications and vaccines safely and efficiently. The Group has been annually growing by double digits on average and has been regularly expanding its portfolio by acquisitions of complementary technologies. In fiscal 2018, Sartorius earned sales revenue of around 1.57 billion euros. Currently, more than 8,900 people work at the Group's approximately 60 manufacturing and sales sites, serving customers around the globe.

### A profile of the DFKI

The German Research Center for Artificial Intelligence (DFKI) was founded in 1988 as a non-profit Public Private Partnership (PPP). It has locations in Kaiserslautern, Saarbrücken and Bremen, a project office in Berlin, a laboratory in Lower Saxony and a branch office in St. Wendel. DFKI is Germany's leading business-oriented research institute in the field of innovative software technologies based on artificial intelligence methods. In the international scientific community, DFKI is one of the most important "Centers of Excellence". Currently about 1,000 employees from about 60 nations are researching innovative software solutions. Their success: more than 130 professors from our own ranks and more than 90 spin-off companies with approx. 2,500 highly qualified jobs.



## Press Release

---

### Press contacts

Petra Kirchhoff Corporate Communications Sartorius Email: <a href="mailto:petra.kirchhoff@sartorius.com">petra.kirchhoff@sartorius.com</a> Phone: +49 551 308 1686	Christian Heyer Corporate Communications DFKI GmbH Email: <a href="mailto:uk-kl@dfki.de">uk-kl@dfki.de</a> Phone: +49 631 20575 1710
---	---

### Scientific contacts

Prof. Dr. Oscar-Werner Reif Head of Corporate Research Chief Technology Officer Sartorius Email: <a href="mailto:oscar.reif@sartorius.com">oscar.reif@sartorius.com</a> Phone: +49 551 308 2401	Prof. Dr. Prof. h.c. Andreas Dengel Head of the Research Department Smart Data & Knowledge Services DFKI GmbH Email: <a href="mailto:andreas.dengel@dfki.de">andreas.dengel@dfki.de</a> Phone: +49 631 20575 1010
--	--