

## Sartorius Stedim Biotech launches new services for mammalian cell bank manufacturing

### Integrated package of new and established services saves time and minimizes risks

Goettingen, GER | Glasgow, UK - August 8, 2019: Sartorius Stedim Biotech (SSB), a leading partner of the biopharmaceutical industry, today announced the launch of new services for mammalian cell bank manufacture under GMP conditions. These services are offered by its subsidiary, Sartorius Stedim BioOutsource, an experienced contract testing organization (CTO) based in Glasgow, UK, and in Cambridge, Massachusetts, USA.

Based on these new services, Sartorius Stedim Biotech now offers the manufacture of GMP master and working cell banks (MCB/WCB) for mammalian suspension cells. This activity will be conducted in a custom-designed 260 m<sup>2</sup> GMP cleanroom facility that has been audited and approved in 2018 by the Medicine and Healthcare Products Regulatory Agency (MHRA) according to the latest guidelines. Exclusively dedicated to mammalian suspension cell lines, the facility uses selected equipment that enables closed-system manufacture of GMP-compliant cell banks, from vial thaw to automated filling. For maximum process reliability and assurance of sterility, this entire manufacturing service is offered as a fully qualified broth technology platform.

The GMP manufacture of master cell banks is a critical step during the development of biopharmaceutical drugs where time is always key. To ensure the best possible timelines, Sartorius Stedim Biotech cell bank manufacturing is offered in a package along with cell bank characterization services that the company has been providing for more than ten years. The combination of its new manufacturing services and its established characterization service for mammalian cell banks allows clients to work with SSB as a single-source provider from vial thaw to released cell banks.

Up to 500 cell bank vials are produced during a manufacturing run, followed by full characterization, genetic stability assessment in compliance with EU and FDA guidelines, and release by the respective qualified person. The cell bank vials produced are then shipped to the client or transferred to a long-term storage facility. In addition, these cell bank manufacturing and characterization services are a perfect match for the cell line development services based on Sartorius Stedim Cellca's CHO expression platform. With all three services combined – cell line development, cell bank manufacturing, and cell bank characterization – Sartorius Stedim Biotech supports clients as a total solutions provider from DNA to released GMP cell bank within a 10-month timeline.

"We're really proud that we now offer even more comprehensive services that can be easily combined with those for CHO cell line development and biosafety. As a result, we provide our clients with an integrated package that saves considerable time and minimizes risks. Furthermore, they will benefit from client-focused customer service with a single point of contact throughout their entire development and manufacturing projects," stated Lucia Rieger, Product Manager for Protein Expression Systems.

### Image file:



Sartorius Stedim Biotech now even offers new services for mammalian cell bank manufacturing

**Download link:** [Mammalian cell bank manufacturing](#)

To find out more about SSB's new services, please click:  
<https://www.biooutsource.com/cell-bank-manufacturing>

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



**A profile of Sartorius Stedim Biotech**

Sartorius Stedim Biotech is a leading international partner of the biopharmaceutical industry. As a total solutions provider, the company helps its customers to manufacture biotech medications safely, rapidly and economically. Headquartered in Aubagne, France, Sartorius Stedim Biotech is quoted on the Eurolist of Euronext Paris. With its own manufacturing and R&D sites in Europe, North America and Asia and an international network of sales companies, Sartorius Stedim Biotech has a global reach. The Group has been annually growing by double digits on average and has been regularly expanding its portfolio by acquisitions of complementary technologies. In 2018, the company earned sales revenue of €1,212.2 million, and currently employs some 5,800 people.

---

**Contact:** Dominic Grone, Trade Press Manager, Sartorius Stedim Biotech GmbH,  
Phone: +49(0)551.308.3324, [dominic.grone@sartorius-stedim.com](mailto:dominic.grone@sartorius-stedim.com), [www.sartorius-stedim.com](http://www.sartorius-stedim.com)