

Göttingen, May 1, 2020

Letter to our new customers

Dear Valued Customer:

We are pleased to announce today that we closed our transaction with the Danaher Group to acquire selected Life Science platform businesses. This acquisition covers the following product segments that now belong to Sartorius:

- Label-free biomolecular characterization business (FortéBio)
- Stainless steel and single-use chromatography skids
- BioSMB chromatography skids
- Chromatography resins
- Conventional chromatography columns
- Single-use technology tangential flow filtration (SUT TFF) skids and stainless steel hollow fibre tangential flow filtration (SS HF TFF) skids
- Microcarriers and particle validation standards (SoloHill)

Though you may already be familiar with Sartorius, please allow us to briefly introduce ourselves: Sartorius is a global partner for life science research and of the biopharmaceutical industry. Together with close to 10,000 staff members, we empower scientists and engineers to simplify and accelerate progress in drug discovery and manufacture. Our innovative tools and technologies cover key processes in drug development and production – from molecule and cell line development to cell growth, harvesting, and purification and filling steps. Over the past two decades, we have continuously added new tools to our portfolio to grow as an integrated solution provider – also the reason behind this acquisition.

What are the next steps?

Under our agreement, all customer contracts for the above portfolios will be transferred from the Danaher Group to us, while contracts that also include other products that remain with Danaher will be jointly managed by both Danaher and us to ensure a smooth transition. Any further order execution such as delivery and invoicing will be realized by the relevant Sartorius entity. Starting May 4th we will initiate the transfer of pending orders. Upon completion a new order confirmation will be provided including relevant company details such as sales entity and bank details. If your company already provided approval for such transfer, there is no need for further actions. In case this did not happen yet and your company objects to such transfer due to any limitations, please contact us within the shortest possible delay at DHR.BPS.TaskForce@Sartorius.com latest by

May 11th 2020. In such event we will reach out to you to identify a suitable solution for all parties. For any new orders our Customer Service organization will support you with any enquiries you may have.

Who am I to contact?

Please note that the FortéBio products will be supported by our Lab Products & Services (LPS) division. We kindly invite you to contact your local LPS representative for any questions related to these solutions. In case you need support to identify the right contact, please visit our website at <https://www.sartorius.com/en/general-contact/local-contact>.

All other acquired technologies will be supported through our BioProcess Solutions (BPS) division. For these technologies please find here the following contact details:

Local Customer Service contact: Please find your local contact at <https://www.sartorius.com/en/general-contact/local-contact>

Service and support assistance:

- for Americas region: BPS.Service.NA@Sartorius.com
- for EMEA region: Technical-Service-Bioprocess@Sartorius.com
- for APAC region: Service.cn@sartorius.com

The entire Sartorius team would like to welcome you as a new customer and we look forward to serving your needs and working with you as a partner. We will contact you soon to discuss our collaboration in more depth. In the meanwhile, please visit us at www.sartorius.com to learn more about our company.

Sincerely,
Sartorius Stedim Biotech GmbH



Bettina Berendsen
Head of Sales and Services BPS