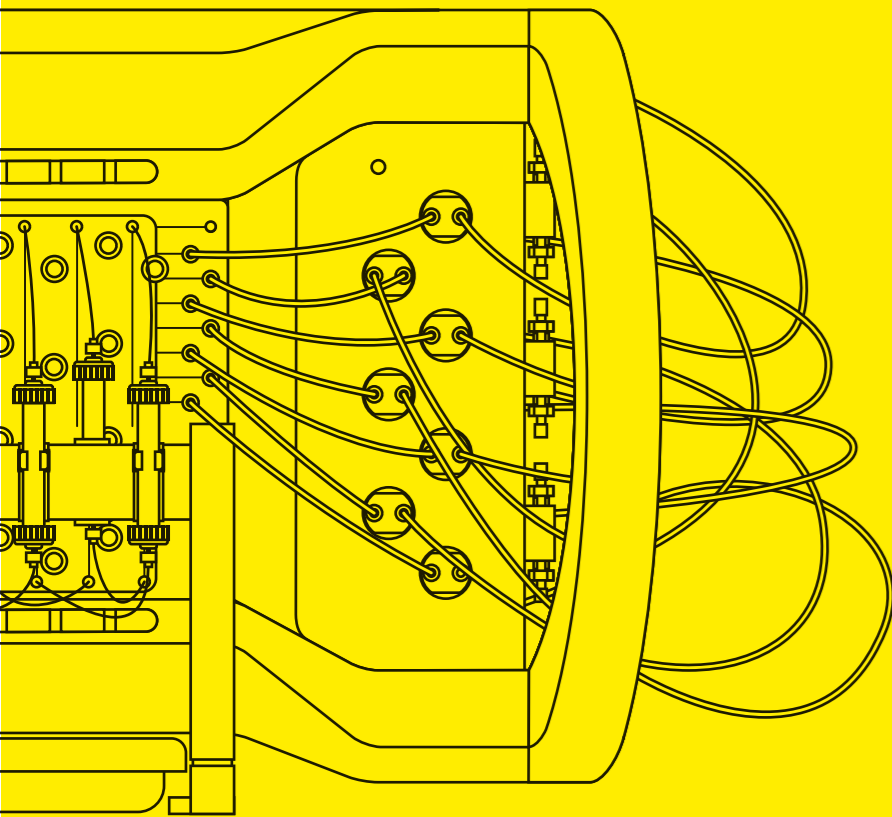




Your First Week of Resolute® BioSMB

Simplify Adoption and Let Sartorius Optimize Your Process Faster
With Step-By-Step Expert Support

See the multi-column chromatography Resolute® BioSMB PD system in action with your product and process. Sartorius can guide you through installation, setup, and implementation. Simply provide our experts with standard breakthrough curves, and they will do the rest, culminating in a complete translation of your process from batch to MCC.



Generate Standard Breakthrough Curves at Different Flow Rates

Using a breakthrough curve modeling tool, Sartorius will support the parameter generation of an optimized process on the Resolute® BioSMB. This helps you achieve a fully optimized process faster with reduced risk.

Optimize With the Resolute® BioSMB Design Tool and Phase Editor

Use the Resolute® BioSMB Design Tool and Phase Editor to enter breakthrough curve fitting outputs and other process information. This tool generates an MCC process tailored to your specific product. The Sartorius Resolute® BioSMB Phase Editor accepts the outputs of the Design Tool and quickly programs the process for automated execution.

Run the Process, See the Results

The design process is transferred to the Resolute® BioSMB system for execution. The system provides a physical readout of the process, demonstrating the benefits of MCC.

Analyze, Optimize, Repeat

Implement small changes based on results to further refine and optimize your process design. Preparation and support from Sartorius experts reduce the time needed to achieve multi-column chromatography process optimization.

Benefits of Resolute® BioSMB Expert Support



Get to an optimized process up to 4x faster



Increase confidence to work with the system



Reduce risk through expert support



Contact us for Resolute® BioSMB demo and trial:
www.sartorius.com/multi-column-chromatography