SARTURIUS

Success Story

Resolute® BioSMB for Increased Productivity at the mAb Capture Step

For customers processing mAbs on protein A resin in a traditional single column batch mode, Resolute® BioSMB provides a single-use multicolumn alternative for the Capture step. After an easy conversion from a single column to a multicolumn format customers reduced protein A consumption by 80% allowing them to enjoy reductions in overall cost and upfront investment in their consumables with no change to the quality of the resulting product.

Customer Challenge

- High Protein A resin cost in mAb capture
- Need to intensify single-use chromatography

Provided Solution

- Introduction to intensified processing at a small scale with the Resolute[®] BioSMB PD system
- Demonstrated scalability of the process with Resolute® BioSMB 80 | 350 systems
- Single-use Flow Path to reduce bioburden risk and cleaning validation
- Application specialist support to assist with process development and technology enablement

Case Profile

Company Type: Large Biopharma, large CMO | CDMO

Related Molecule: mAb

Related Application: mAb capture process at clinical scale using multi-column chromatography



Single Column Batch:

- Column diameter: 40 cm
- Volume of resins: 20 L
- Specific productivity: 16 g/L/h



Multi Column Chromatography:

- Column diameter: 14 cm
- Volume of resins: 3.85 L
- Specific productivity: 56 g/L/h

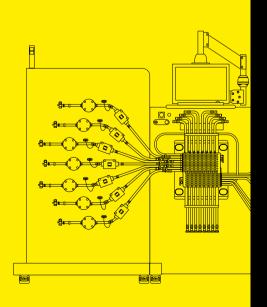
80% cost savings by reducing Protein A demands

3-5-fold increase in productivity

Scale up to sinlge-use solution

Easy integration

into existing batch process



Contact a Sartorius
Specialist or Visit
www.sartorius.com