

## Intensification of mAb Processes Leveraging Sartobind® Rapid A and Full Connected Membrane-Based DSP

Fabien Rousset

Sartorius Stedim FMT S.A.S., Zone Industrielle les Paluds, 300 Avenue de la Fleuride, 13400 Aubagne, France  
Contact: Fabien.Rousset@sartorius.com

### Introduction

In the field of mAb purification, high performances chromatography membranes that are ready-to-use, for “one-batch-one device” manufacturing strategy, take off. The newly Protein A capture technology “Sartobind® Rapid A”, used in rapid cycling conditions, brings a 10-fold higher productivity (203 g/L vs 14 g/L with traditional resins), has similar performances for DBC, yield and HCP | hcDNA removal. This allows new-generation full membrane-based purification platforms.



The first milestones to fully membrane-based process is achieved by implementing a competitive double-flow-through polishing process with connected Sartobind® Q and Sartobind® S. Comparable purity and yield are obtained (> 98% for each flowthrough steps) with a strong footprint reduction of the purification process. The second step to a full membrane process is combining the Resolute® MCC multicolumn technology with protein A, AEX and CEX membranes in parallel batch mode. This results in increasing even more the productivity (> 400 g/L/h) compared to a resin-based multi-column chromatography process (< 200 g/L/h).

This innovative Sartobind® Rapid A combined with process intensification solutions demonstrates that alternative mAb purification platforms are safer and highly competitive against classic resin-based approaches.



### Sartobind® Rapid A as Convectdiff Membrane vs Purely Convective | Diffusive Materials

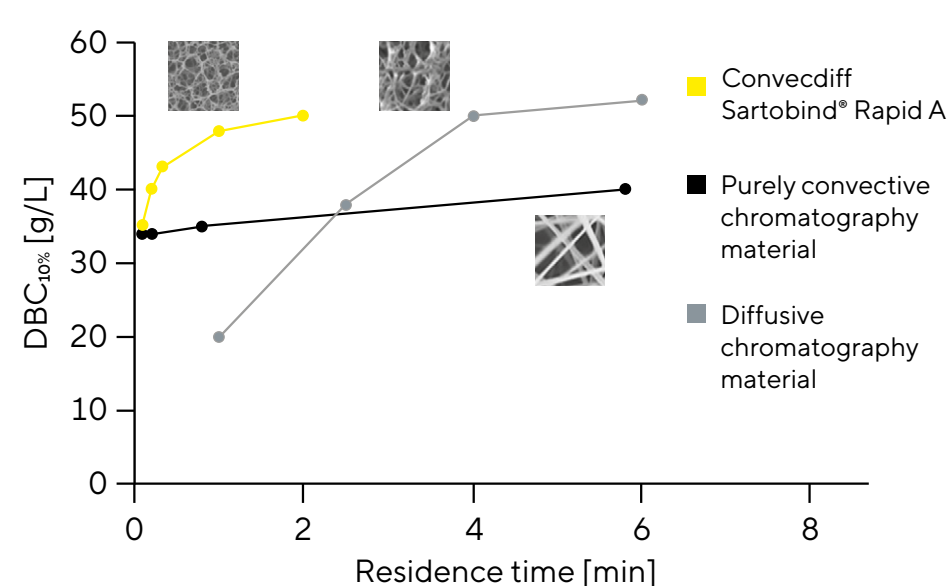
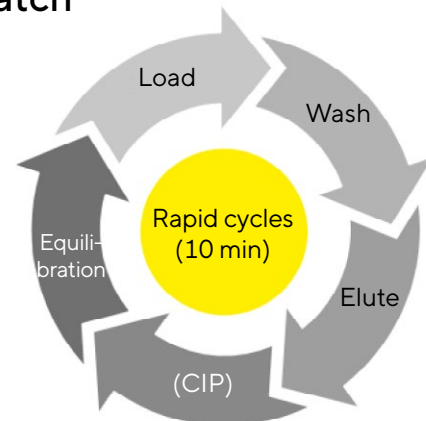


Figure 1: SDBC<sub>10%</sub> As a Function of Residence Time for Commercially Available Materials and the Convectdiff Sartobind® Rapid A.

Convectdiff materials ally high DBC & high flowrate to enable short cycle time

Sartobind® Rapid A enabling lifetime capacity utilization in one single-batch

- Fast cycles: 10 - 15 min
- ~30 - 150 cycles/batch



### CPP and CQA of IgG Purified With Sartobind® Rapid A and Standard Resin<sup>1</sup>

In this comparison, we show how Sartobind® Rapid A compares to standard protein A resin. Both materials were tested with the same feed material. The analyzed data show a very good comparability of Sartobind® Rapid A with the protein A resin. The membrane showed superior performance in DNA reduction and protein A leaching, with a 14.5-fold increase in productivity.

	Sartobind® Rapid A	Protein A Resin
DBC10% [g/L]	42.9±0.8	30.4±0.5
Residence time [min]	0.2	4.0
Yield [%]	94.7±0.2	96.4±0.4
HCP reduction [LRV]	2.2±0.2	2.3±0.1
hcDNA reduction [LRV]	2.9±0.2	2.3±0.1
Protein A leached [ppm]	2.7±0.7	6.7±0.3
av. Productivity [g/L*h]	203.6	14.1

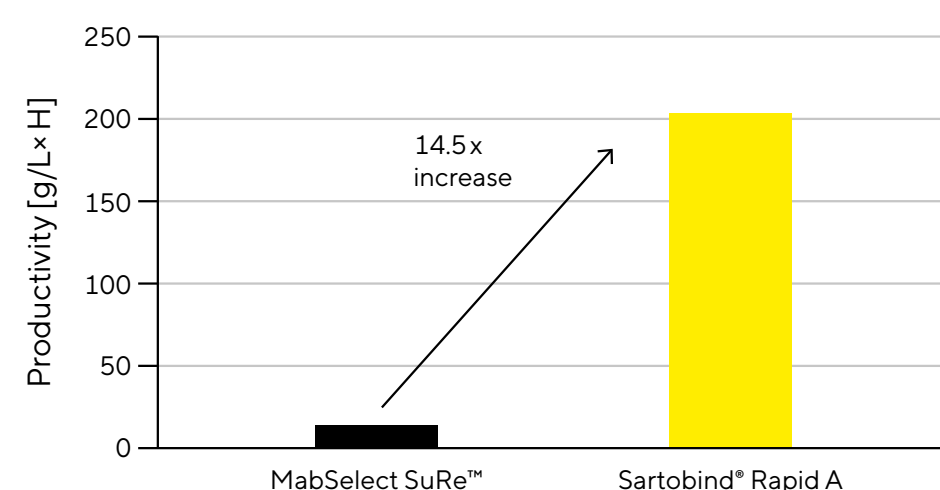


Figure 2: Productivity Comparison of HiTrap® MabSelect SuRe™ to Sartobind® Rapid A



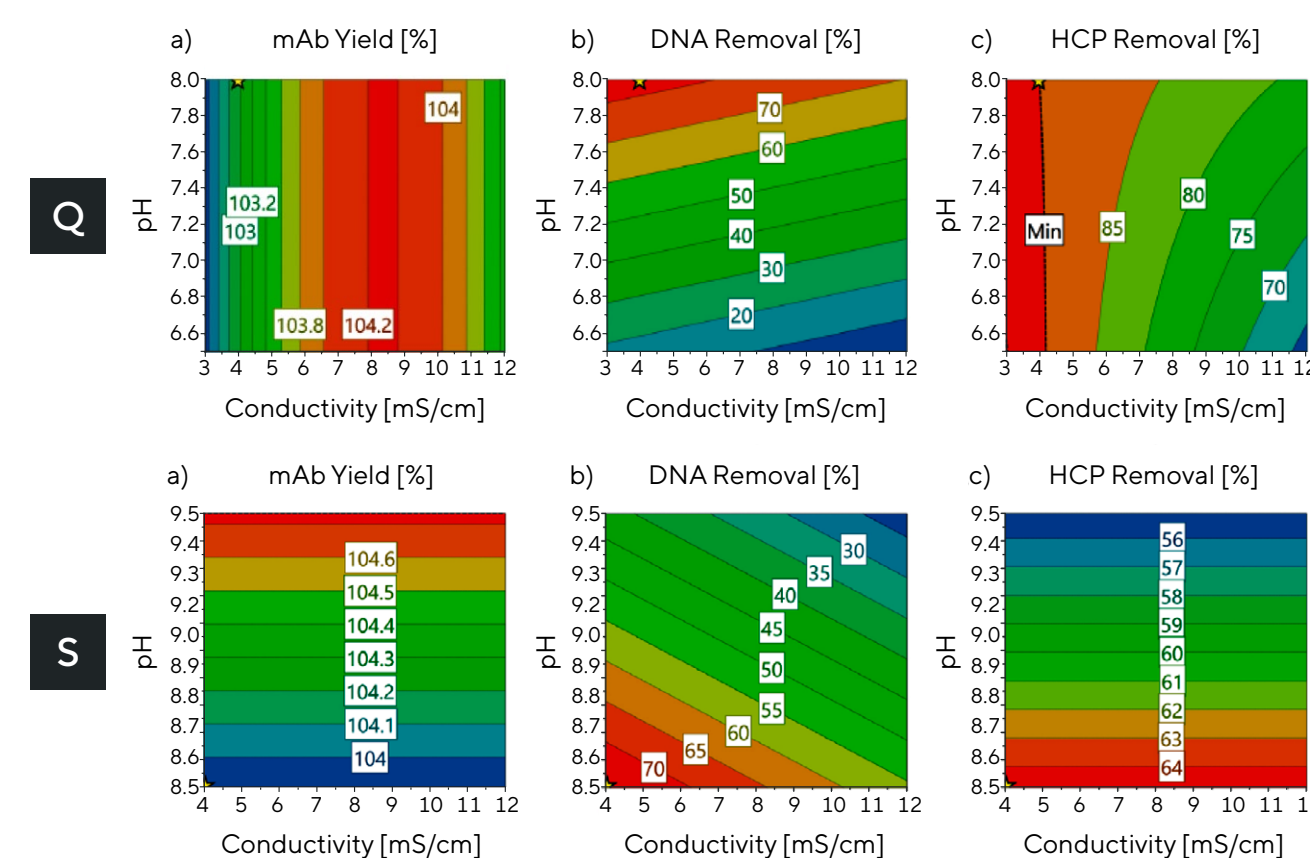
### Levels of Intensification for Downstream Processing

Level 0	Level 1	Level 2	Level 3	Level 3.1
Standard batch, standalone UO	Intensified, standalone UO increases the individual step productivity; higher cycling (RCC or MCC), improved buffer management (ILD, ILC), high DBC resins; pooling tanks	Connected process: at least 2 UO; subsequent steps started before first step is finished; could be staggered batch; may have pooling tanks; software orchestration is beneficial; might be called clustered or linked process	Continuous process: more integrated case of a connected process; steady state flow; only small intermediate (surge) tanks; software orchestration is a must have; long run times; closed processing;	Flow-through continuous process: further integrated case with complete steady state flow. All bind and elute steps are replaced with flow through mode. Molecule does not stop – ideally no intermediate (surge) tanks

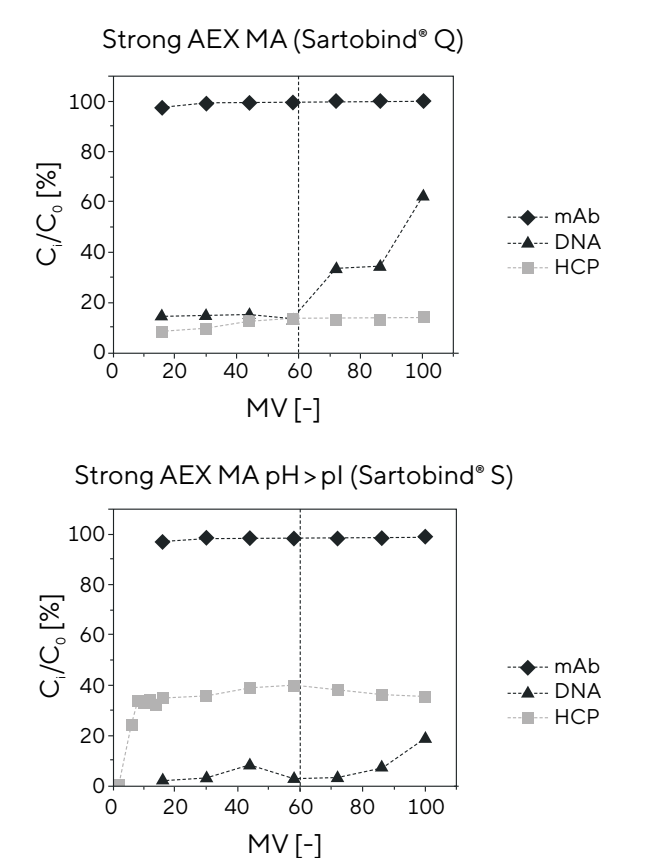


### Double Flow-Through With Membrane Adsorbers<sup>2</sup>

#### Step 1: DOE to Define Buffer Conditions



#### Step 2: Breakthrough Curves

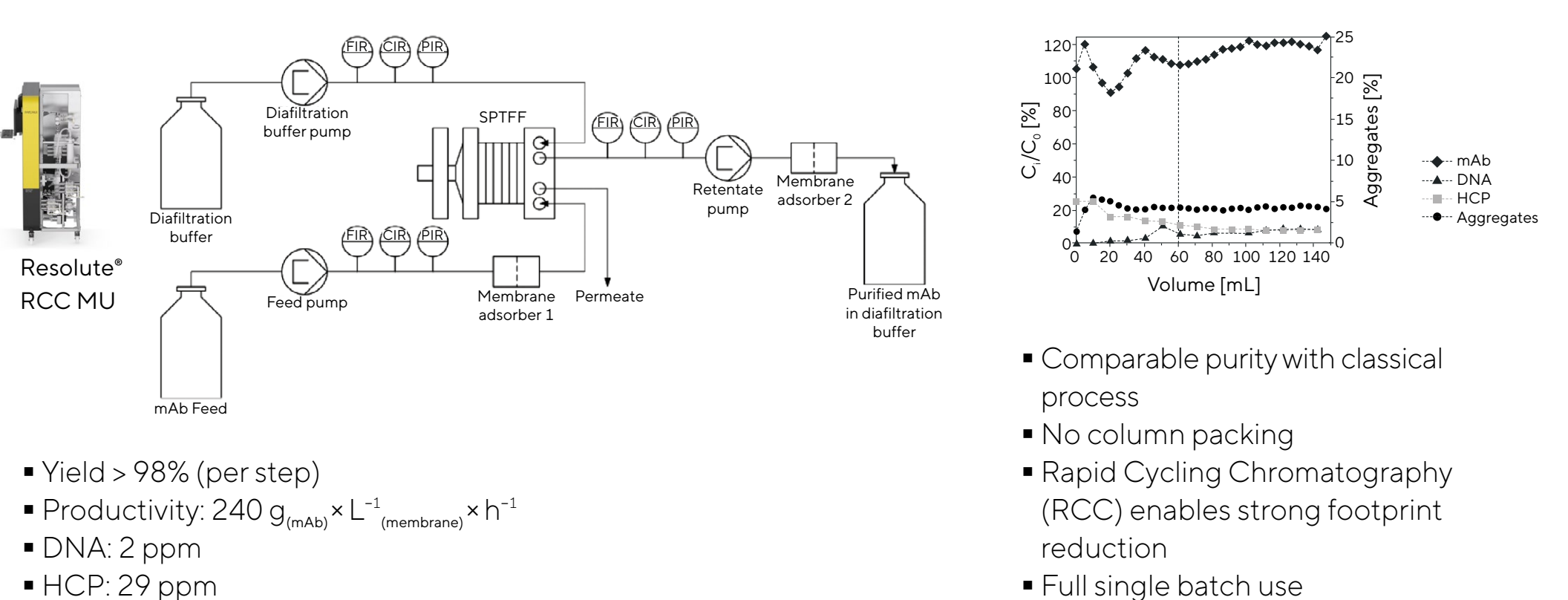


Note: Sartobind Rapid A - Beta Test Opportunity, DOE done with MODDE<sup>®</sup> 13



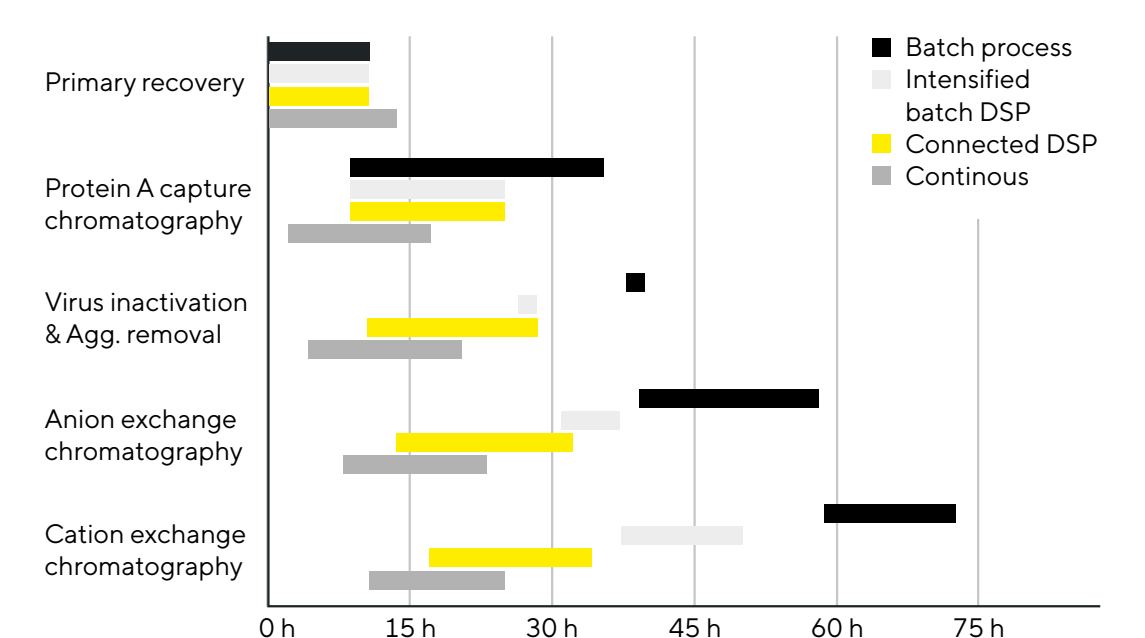
### Chromatography Membrane – Towards Connected Process<sup>2</sup>

#### Step 3: Connected Processing



### Downstream Intensification Reduces Processing Times

- Each step starts before the previous one ends
- This enables processing the sub-batches from protein A elution
- Reduction of intermediate tanks and chromatography columns' size
- Lower footprint
- Lower OPEX



### The Power of Connected Membrane Processes

Cascade	Resin (MCC Connected Process)	Sartobind®
Number of steps	3	3
Final yield [%]	83	86
DNA [ppb]	<3	<3
HCP [ppm]	3	30
HMW [%]	0.2	0.3
Total process time [h]	10.4	4.0
Av. Productivity [g/h]	179	465

- Connecting the process and using MCC or parallel batch multiplies productivity or drops costs
- Comparable purity and yield
- Lower footprint

Clarified harvest
Sartobind® Rapid A (9.6 L @ 45 g/L) 3 × 3.2 L membrane
Sartobind® Q (1.6 L @ 500 g/L)
Sartobind® S (1.6 L @ 500 g/L)



### Conclusion

Due to inherent structural characteristics, Sartobind® Rapid A offers unique possibilities in the area of full membrane-based ultra fast mAb purification process

- Ready-to-use and One batch – One device manufacturing strategy enabled thanks to Sartobind®
- DBC highly competitive
- Short cycle time (< 30 min)
- High number of cycles per batch (up to 150)
- Double Flow-Through polishing with Sartobind® Q | S
- Full membrane process with ultra high productivity – 3x compared to connected resin-based process
- Innovative agarose platform, scalable, robust for a wide variety of mAbs with limited back-pressure at large scale
- Availability of modular cassette format enables scaling to large production processes