

SARTORIUS

Simplifying Progress

Sartobind® Lab
The Future of Lab Purification

April 2023



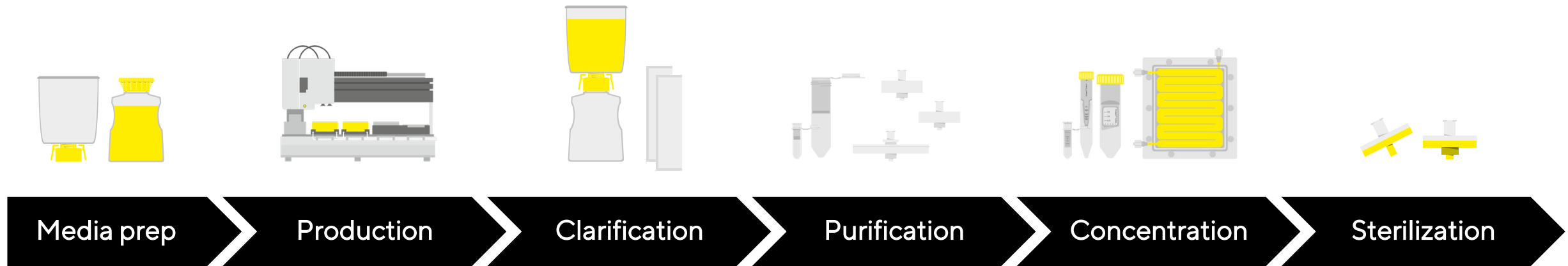
Sartobind® Lab

The Future of Lab Purification

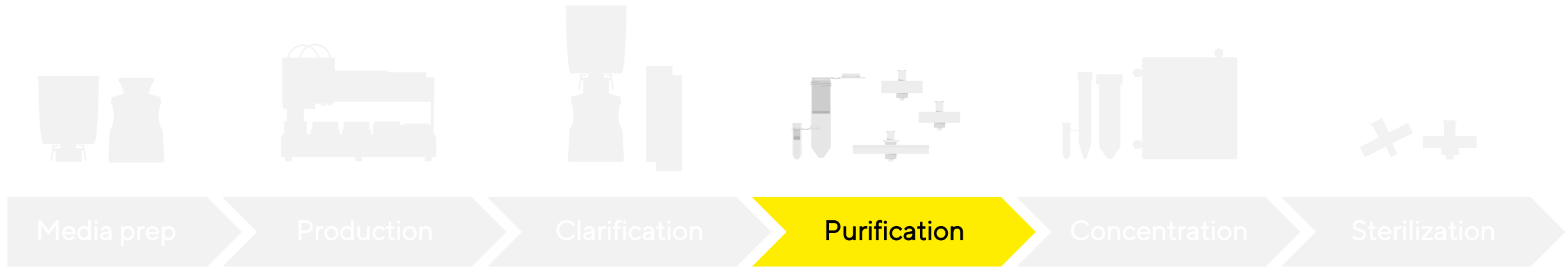
- 1 Purification Workflows
- 2 Sartobind® Lab
- 3 Chromatography
- 4 Membrane Adsorbers
- 5 Competitor Landscape
- 6 Applications and Lead Qualification



Purification in Research and Development



Purification in Research and Development

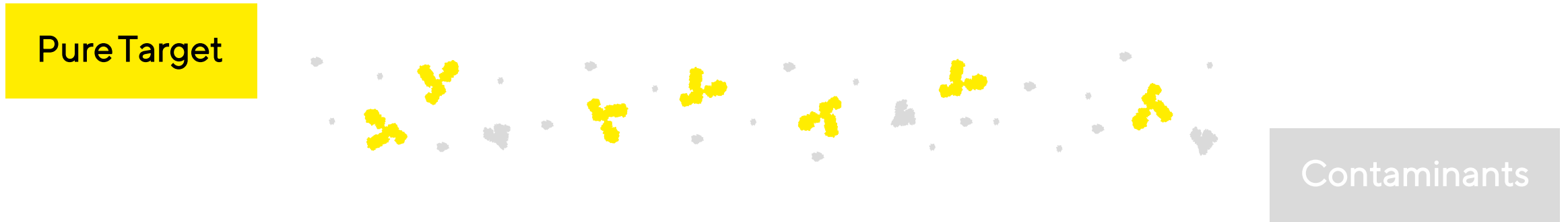


Sartobind® Lab: Purification As Simple As Filtration

- **Ready-to-use** membrane adsorber units
- Syringe filter format for **reliable, equipment-free** purification
- Choice of ligands for **ion exchange** and **affinity** chromatography
- **Faster** with **higher yields** than resin-based columns

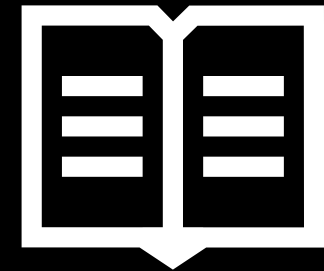
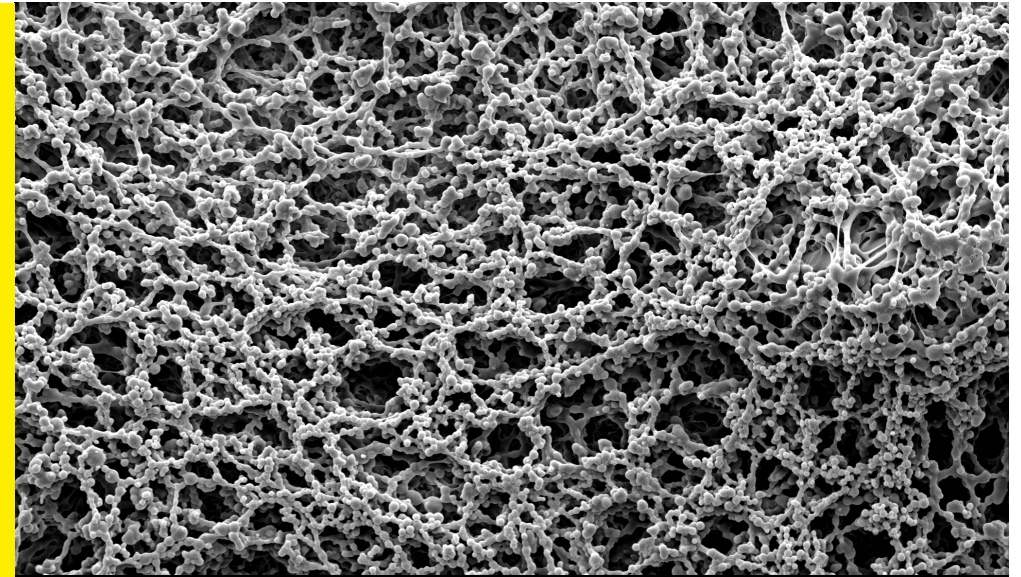


How Are Samples Purified?



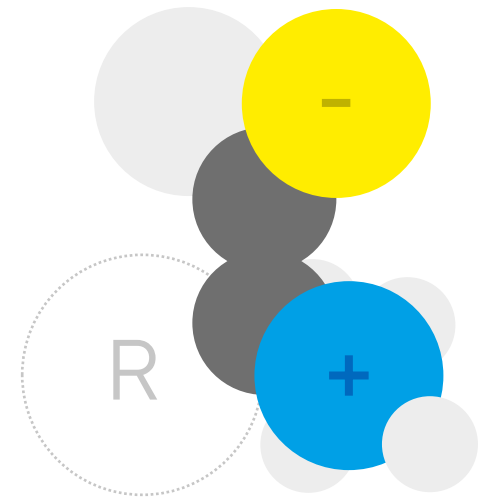
Chromatography

A technique for **separating macromolecules** by passing them through a stationary phase matrix.

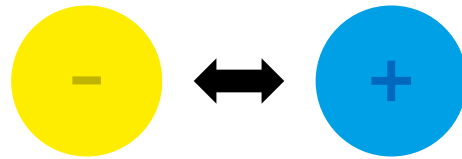


Separation Methods

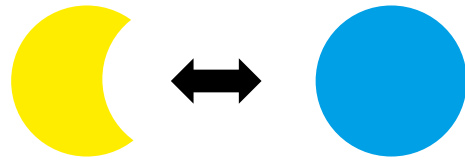
- Every macromolecule is unique
- **Charged molecules** will interact with a ligand of **opposite charge**
- Other molecules may bind to **specific ligands**
- **Ion exchange (IEX)** uses charged ligands for separation of **many molecules**
- **Affinity chromatography (AC)** is used to purify **tagged molecules** or antibodies








Separation Methods



Separation Methods



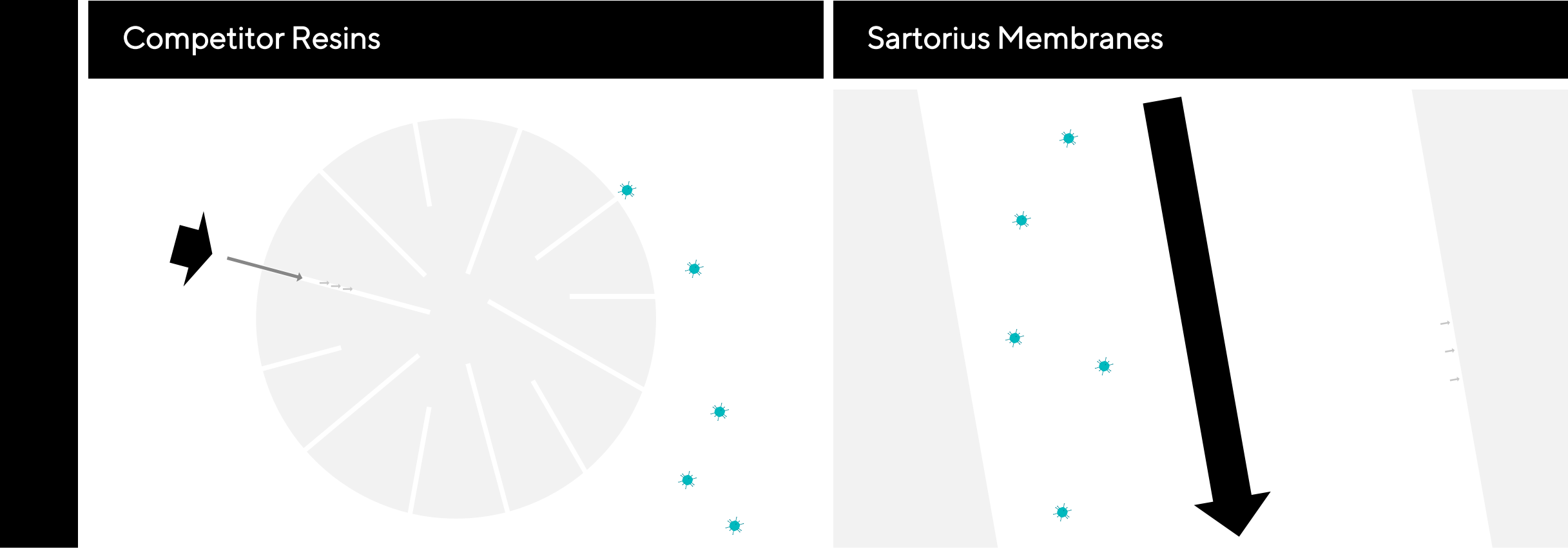
Matrix Types

	Competitor Resins	Sartorius Membranes
	Synthetic beads	Regenerated cellulose membranes
	Microporous, 15 - 40 nm	Macroporous, 3 – 5 µm
	Setup required	Plug and play
	Slow, diffusive flow	Fast, convective flow
	Capture and polishing of proteins up to 200 kDa	Capture and polishing for most macromolecules

Matrix Types

	Competitor Resins	Sartorius Membranes
		

Matrix Types



Purification Challenges



High Cost

Requires an FPLC system.
Unreliable results with reduced equipment.



Slow Setup

Labour and time intensive preparative steps.
Column packing, buffer degassing, system setup, preservative removal.

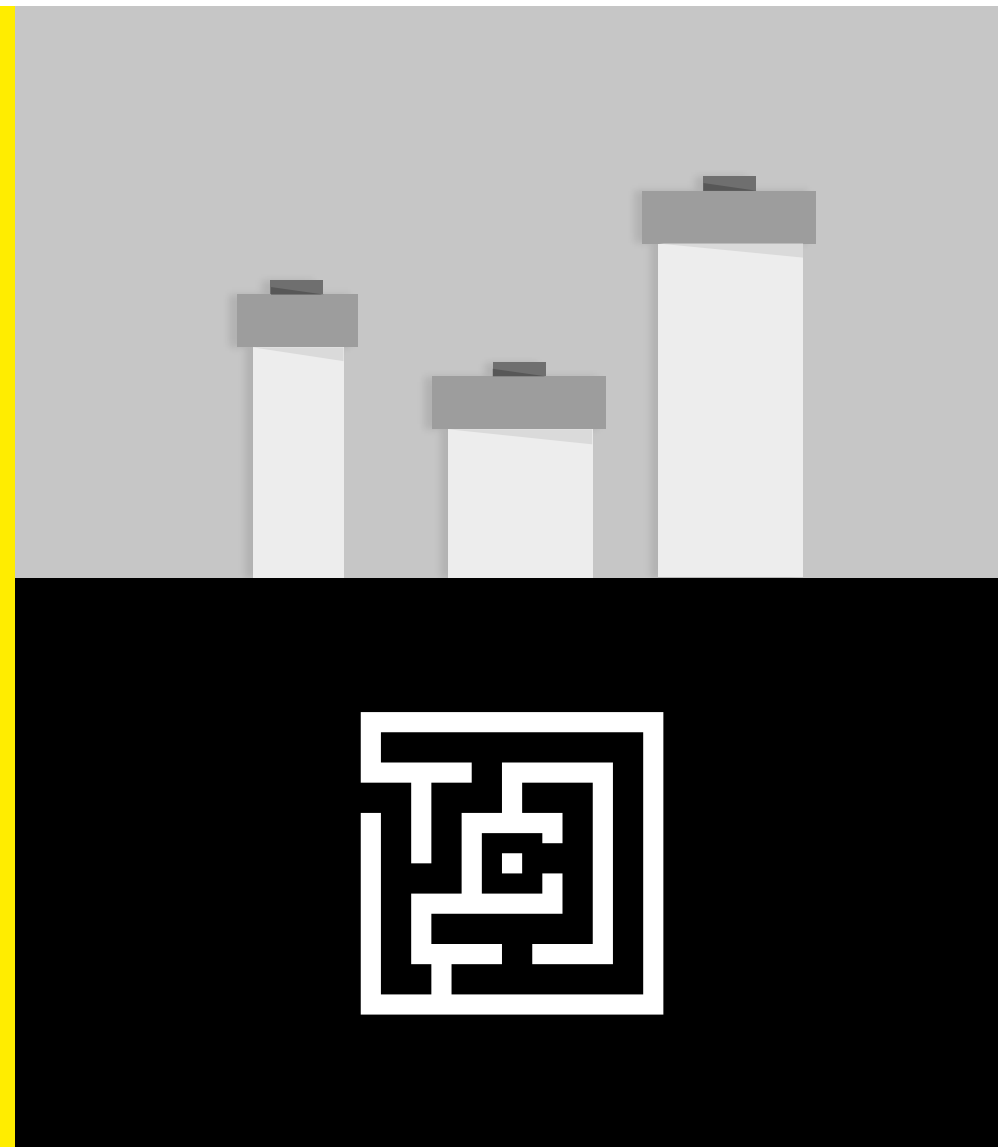


Low Yields






Diffusive flow and size exclusion effects.
Longer cycle times with reduced yields and limited productivity.

Resin Columns

Lab purification is **expensive**,
complicated and **slow**.



Competitors

Company	Products	Pricing ^[1]	Technology	Processing	Capacity ^[2]
	<ul style="list-style-type: none"> Vivapure, Sartobind IEX and AC 	<ul style="list-style-type: none"> Value-based € 0.02 per mg 	<ul style="list-style-type: none"> Membrane Adsorber Syringe, FPLC 	<ul style="list-style-type: none"> 5 - 30 MV/min Up to 112 cycles/hr 	<ul style="list-style-type: none"> 60 - 80 mg BSA 7.5 mg his-protein
	<ul style="list-style-type: none"> HiTrap, SpinTrap IEX and AC 	<ul style="list-style-type: none"> Competitive € 0.03 per mg 	<ul style="list-style-type: none"> Resin FPLC (syringe) 	<ul style="list-style-type: none"> 1 - 4 CV/min Up to 12 cycles/hr 	<ul style="list-style-type: none"> 55 - 160 mg BSA 10 - 40 mg his-protein
	<ul style="list-style-type: none"> EconoFit IEX and AC 	<ul style="list-style-type: none"> Competitive € 0.07 per mg 	<ul style="list-style-type: none"> Resin FPLC (syringe) 	<ul style="list-style-type: none"> 0.3 - 6 CV/min Up to 13 cycles/hr 	<ul style="list-style-type: none"> 37 - 180 mg BSA 15 - 40 mg his-protein
	<ul style="list-style-type: none"> MiniChrom, POROS ^[4] IEX and AC 	<ul style="list-style-type: none"> Cost-plus € 0.81 per mg 	<ul style="list-style-type: none"> Resin FPLC 	<ul style="list-style-type: none"> 0.33 CV/min Up to 1 cycle/hr 	<ul style="list-style-type: none"> 120 - 190 mg BSA
	<ul style="list-style-type: none"> Acrodisc ^[5] IEX 	<ul style="list-style-type: none"> Cost-plus € 1.36 per mg 	<ul style="list-style-type: none"> Membrane Adsorber Syringe (FPLC) 	<ul style="list-style-type: none"> 5 MV/min Single use 	<ul style="list-style-type: none"> 10 - 60 mg BSA

^[1] Pricing for Sartobind Lab 75 | 100 and competitor equivalents collected from mfr. websites November 2022. Price/mg based on 1 hour production at shortest cycle times.

^[2] Binding capacities for Sartobind Lab 75 | 100 and competitor 1 mL devices collected from mfr. datasheets and user manuals November 2022.

^[3] Cycle times based on manufacturer recommended min. volumes and max. flow rates, where specified. One cycle = equilibrate + wash + elute + regenerate.

^[4] PL Repligen OPUS MiniChrom Column platform available with all listed ligands from Merck, or all excluding C from Thermo Fisher Scientific.

^[5] Acrodisc Mustang | Mustang XT offer direct scale up in SU units, with smaller, 0.65 - 0.8 µm pores; S sold separately only as XT version (kit with Q, S and E for "standard").

Sartorius' Solution



Out-of-the-Box Flexibility

Connect directly to a syringe or FPLC system.
Reliable, equipment-free purification.



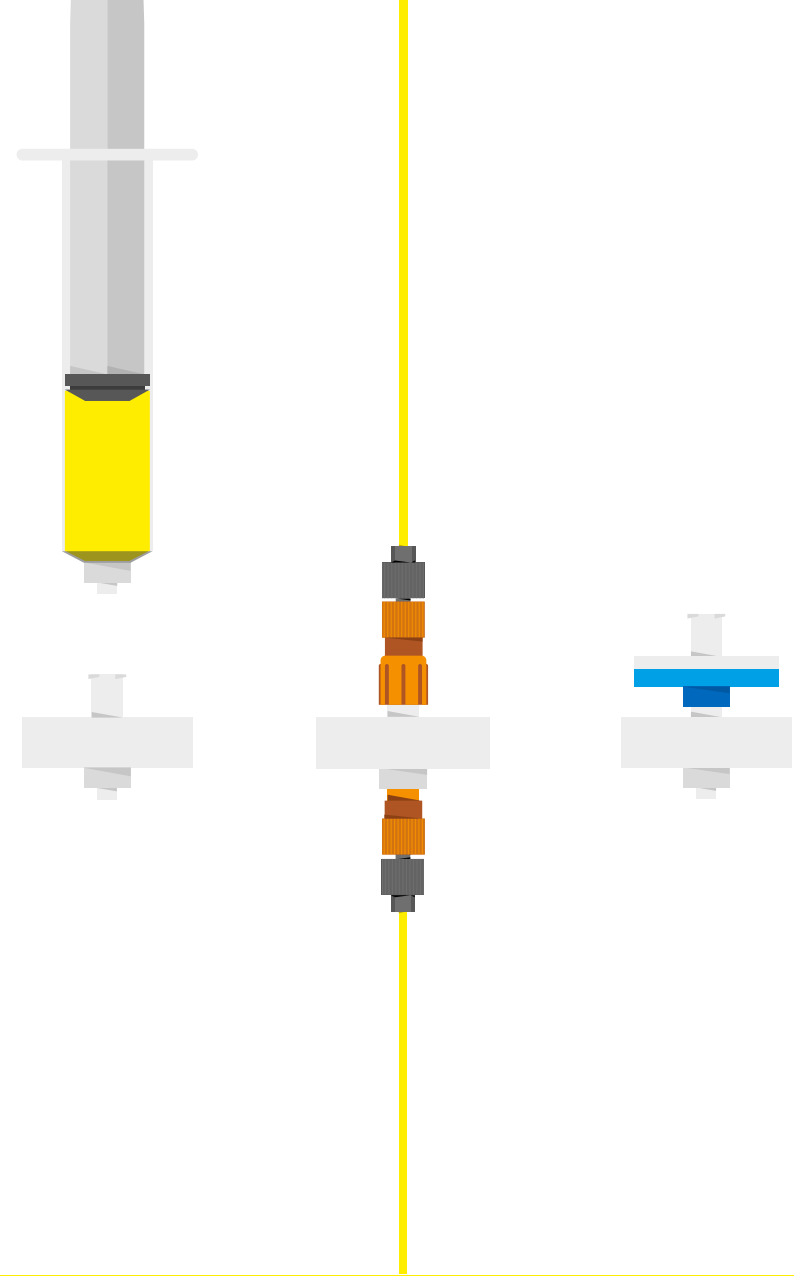
Minimum Setup Time

Ready-to-use, syringe filter format.
Avoid and even combine preparative steps.



High Yields in Seconds

Rapid, convective flow.
Faster flow rates, shorter cycle times and higher productivity^[1].



^[1] Productivity (g/hr) calculated from manufacturer recommended maximum flow rates and DBCs, and min. equilibration, wash, elution and regeneration volumes for standard reference proteins (BSA or lysozyme for anion or cation exchangers, respectively). Dotted bars indicate data from alternative reference proteins.

Sartorius' Solution



Out-of-the-Box Flexibility

Connect directly to a syringe or FPLC system.
Reliable, equipment-free purification.



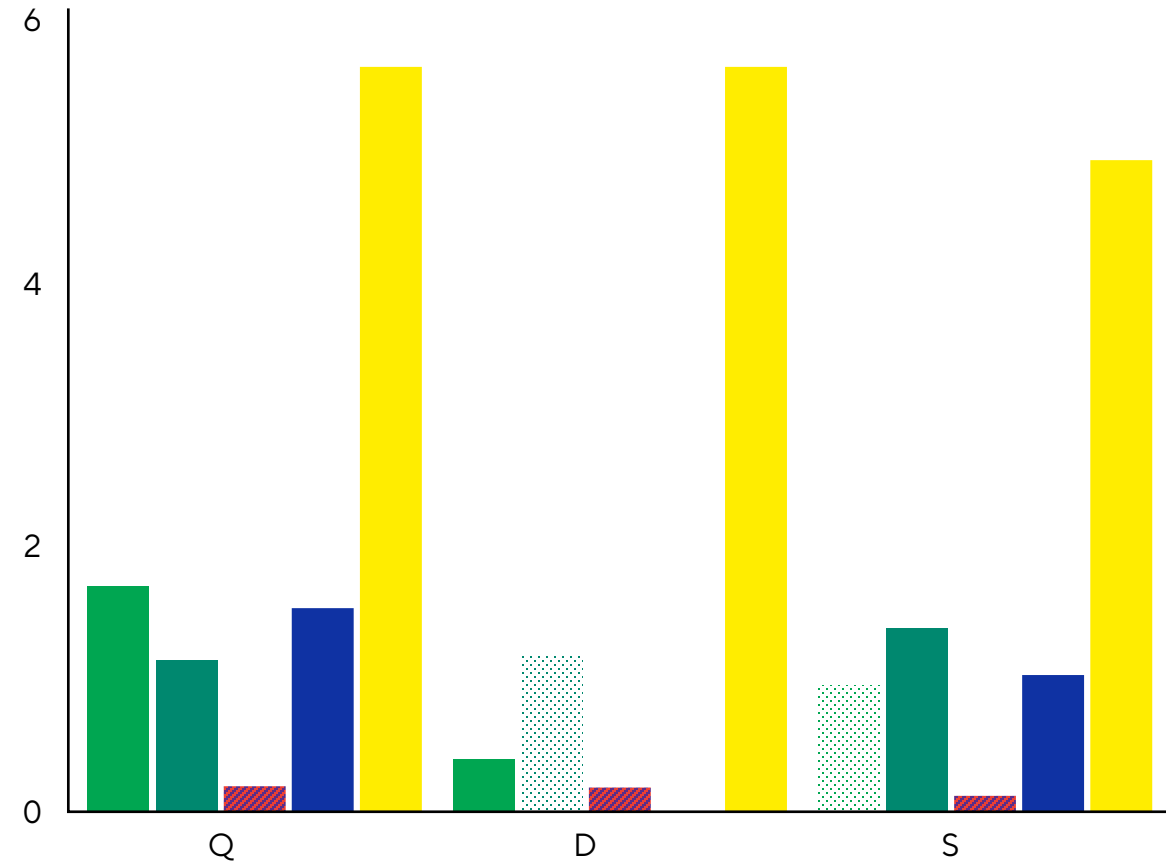
Minimum Setup Time

Ready-to-use, syringe filter format.
Avoid and even combine preparative steps.



High Yields in Seconds

Rapid, convective flow.
Faster flow rates, shorter cycle times and higher productivity^[1].



^[1] Productivity (g/hr) calculated from manufacturer recommended maximum flow rates and DBCs, and min. equilibration, wash, elution and regeneration volumes for standard reference proteins (BSA or lysozyme for anion or cation exchangers, respectively). Dotted bars indicate data from alternative reference proteins.

Use Cases, Customers and Applications

- **Screening, optimization and preparative purification in Academic and Biopharma Research Labs**
- **Capture Purification**
 - Proteins
 - Nucleic acids
 - Viruses
- **Flow Through Polishing**
 - HCPs and DNA
 - Viruses
 - Endotoxins



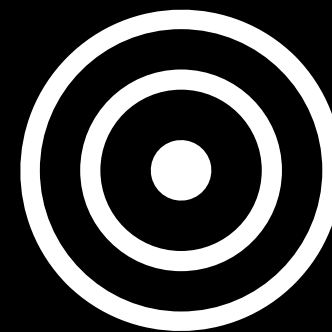
Lead Qualification

- Look out for **FPLC systems** from Cytiva, Bio-Rad or Practichem
- Listen out for terms like **cell culture**, **purification** or **chromatography**
- Ask if your customers need **pure proteins**, **nucleic acids** or **viruses**
- Find out **how they currently purify** these molecules
- Ask if it would be beneficial to **purify equipment-free**



Sartobind® Lab

Flexible, easy and fast
purification for R&D.



Thank you.

Dr. John Cashman
Product Manager – Lab Ultrafiltration
Marketing
Mobile +44 7896 664061
john.cashman@sartorius.com

The Sartorius logo is displayed in a bold, black, sans-serif font. The letters are closely spaced, and the 'S' and 'A' are particularly prominent. The logo is centered within a bright yellow rectangular area that occupies the bottom right portion of the slide.