

# Upgrade your monoclonal antibody purification workflow with a new source of Protein A resin

## Introducing Thermo Scientific MabCaptureC affinity matrix

**ThermoFisher**  
SCIENTIFIC

Hendrik Adams<sup>1</sup>, Wannie Horrevoets<sup>1</sup>, Elina Kleijs<sup>1</sup>, Jenny England<sup>2</sup>, Alejandro Becerra<sup>2</sup>, Jessica de Rooij<sup>1,2</sup>, Pim Hermans<sup>1</sup>, Marien Vis<sup>1</sup>, Frank Detmers<sup>1</sup>

1. Thermo Fisher Scientific, Leiden, the Netherlands, 2. Thermo Fisher Scientific, Bedford, MA

Bioprocessing

### INTRODUCTION INTO A NEW SOURCE OF PROTEIN A

Since decades Protein A chromatography has been the method of choice for the purification of monoclonal antibodies. Modifications to the first generation of Protein A ligands has given way to more robust and efficient purification resins allowing the design of antibody purification processes with greater flexibility. With the introduction of Thermo Scientific<sup>™</sup> MabCaptureC<sup>™</sup> affinity matrix, a new Protein A resin is now available. Featuring high capacity and a highly cross-linked agarose backbone, this Protein A resin is specifically designed to help improve the productivity and efficiency of your antibody purification process.

#### MabCaptureC affinity matrix – features and benefits

Our MabCaptureC<sup>™</sup> resin is based on a new engineered and in-house produced protein A ligand, recombinantly expressed in yeast. It features:

- **High binding capacity:** >50 g/L IgG at 4.8 min residence time
- **Alkali stable ligand:** >100 cycles at 0.2M NaOH
- **Highly cross-linked agarose backbone (Praesto<sup>™</sup> jetted technology)**
- **Uniform bead size (75 μm +/- 10 μm) delivering improved performance characteristics**
- **Excellent scalability and free of animal components – allowing use in commercial manufacturing**

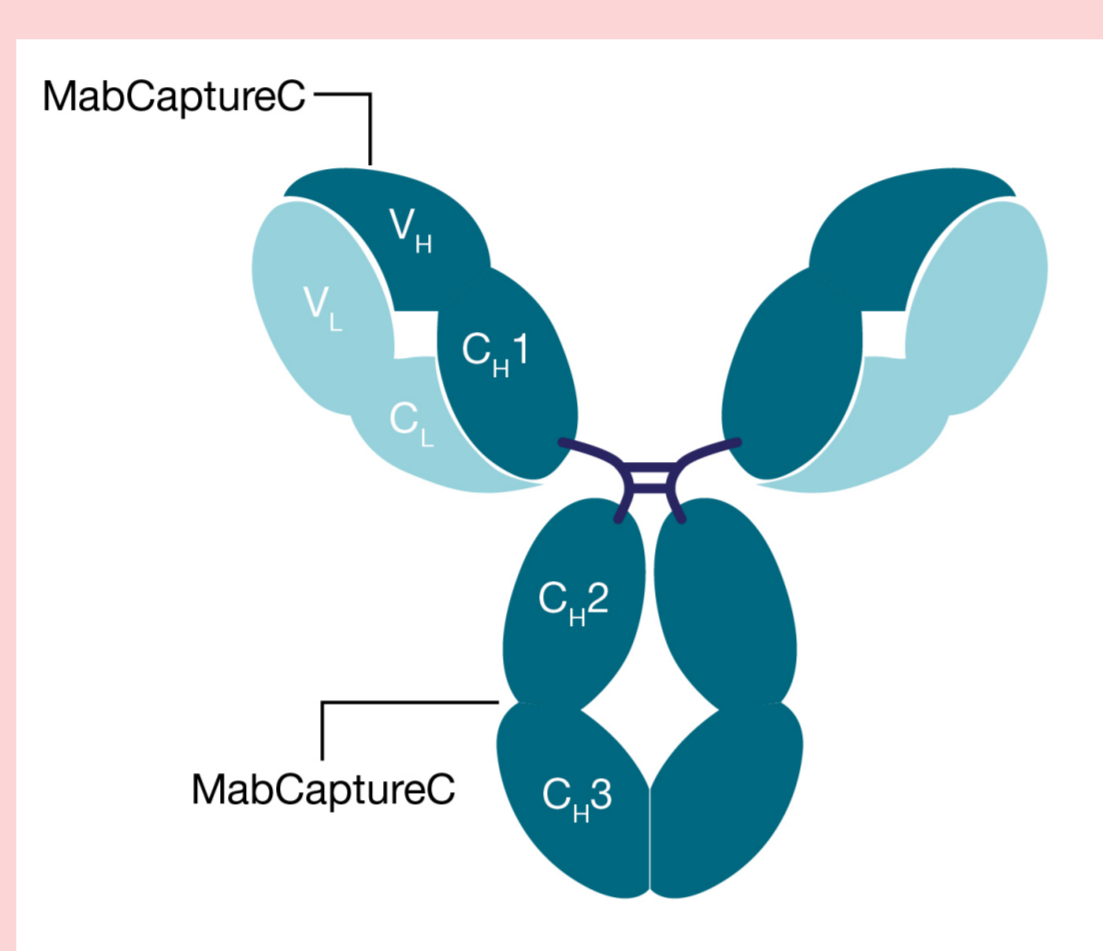


Fig 1. The MabcaptureC resin binds at the C<sub>H</sub>2-C<sub>H</sub>3 interface and V<sub>H</sub>3 region of IgG

### MabCaptureC resin Binding Capacity and Elution Properties

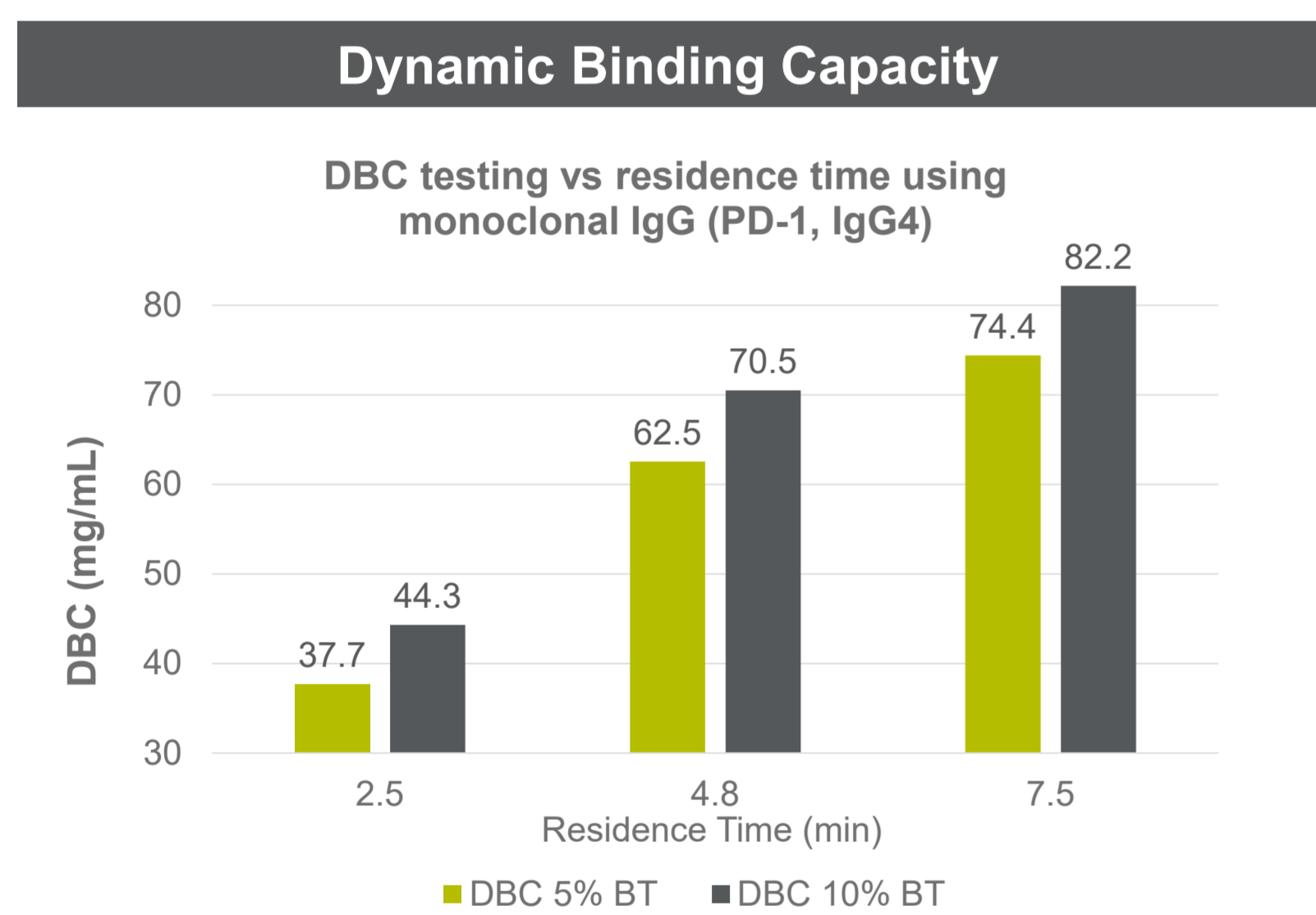


Fig 2. MabCaptureC resin dynamic binding capacity measured at 5 & 10% breakthrough at 2.5-, 4.8- and 7.5-min residence time using monoclonal IgG (load density 70mg/mL)

- ✓ **The MabCaptureC resin demonstrates high dynamic binding capacity**
- ✓ **The resin shows efficient elution (>98%) at pH 3.0-3.5**

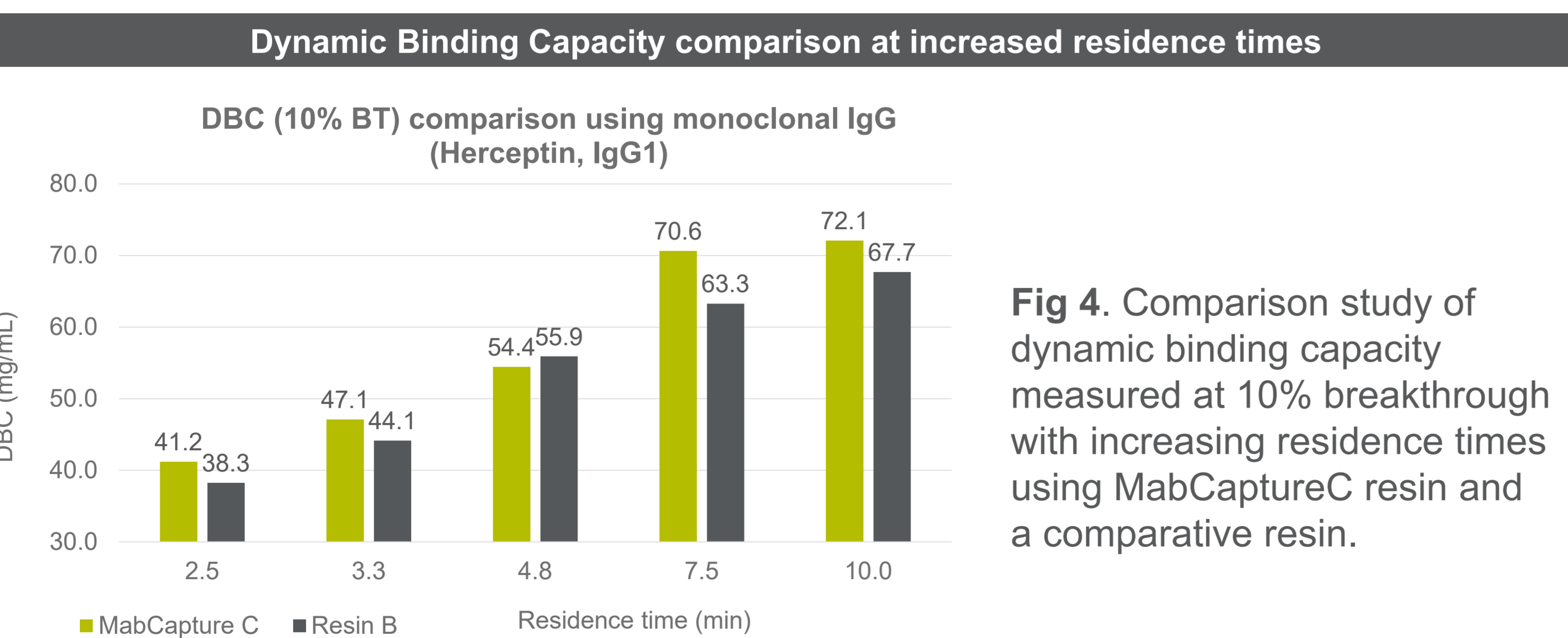


Fig 4. Comparison study of dynamic binding capacity measured at 10% breakthrough with increasing residence times using MabCaptureC resin and a comparative resin.

- ✓ **The MabCaptureC resin shows excellent binding capacity compared to comparative resins**
- ✓ **High dynamic binding capacity is shown at increased residence times, allowing processing of high-titer feedstocks**

#### TRADEMARKS/LICENSING

Caution: For research use or further manufacturing. Not for diagnostic use or direct administration into humans or animals. © 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Praesto is a registered trademark of Puroilite. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

### MabCaptureC resin cleaning, reusability and impurity removal

#### Advised cleaning strategy

- ✓ **Acid strip after every run:**  
Up to 0.5 M Citric acid or Acetic acid
- ✓ **Additional cleaning/sanitization steps**  
Process optimization mainly depending on type of feed
  - Cleaning after every run with 0.2 M NaOH
  - If needed, cleaning with 0.4 M NaOH after every 5<sup>th</sup> or 10<sup>th</sup> run

#### Reusability: Life cycle study with CIP 0.2M NaOH for 100 cycles

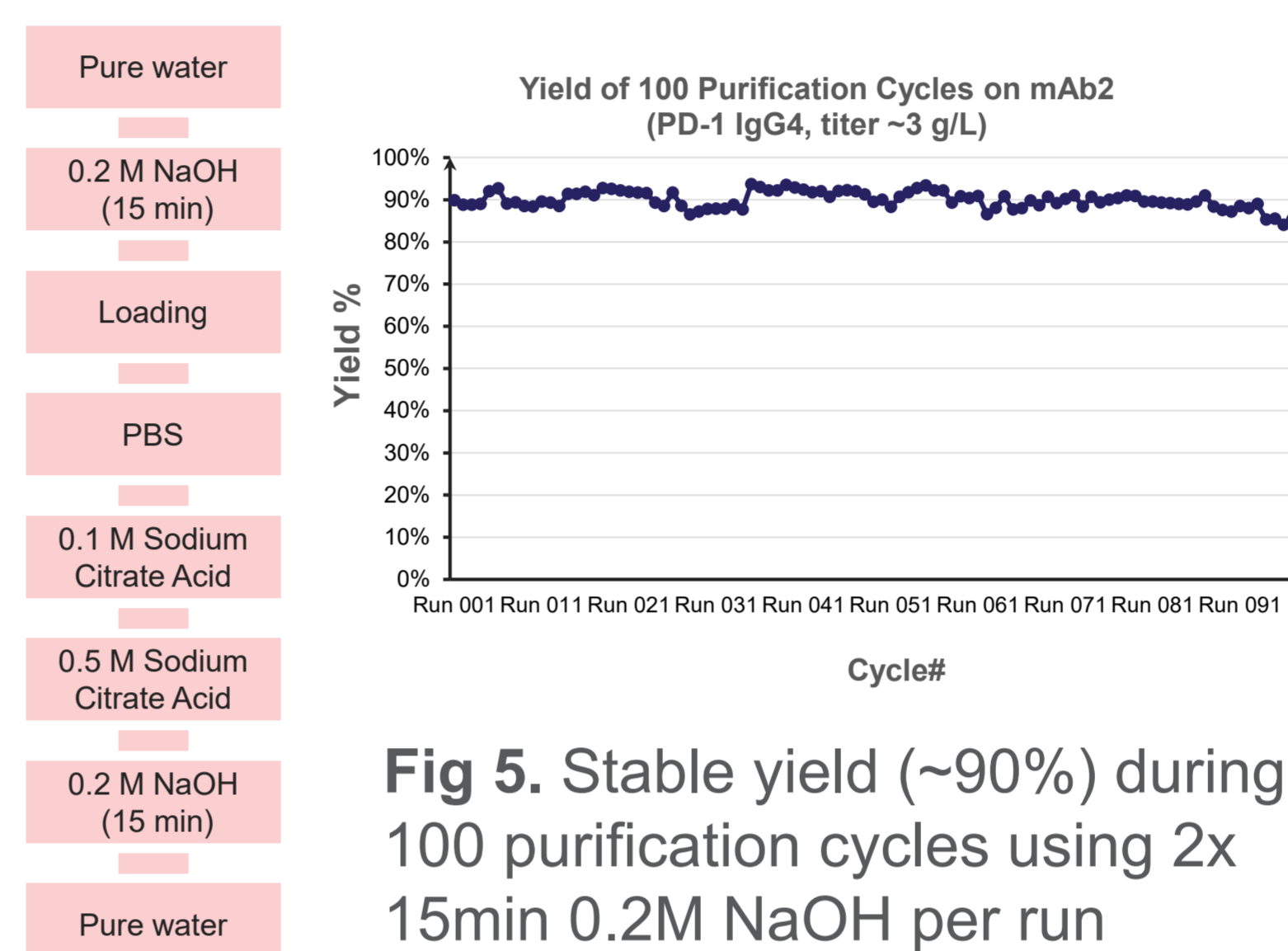


Fig 5. Stable yield (~90%) during 100 purification cycles using 2x 15min 0.2M NaOH per run

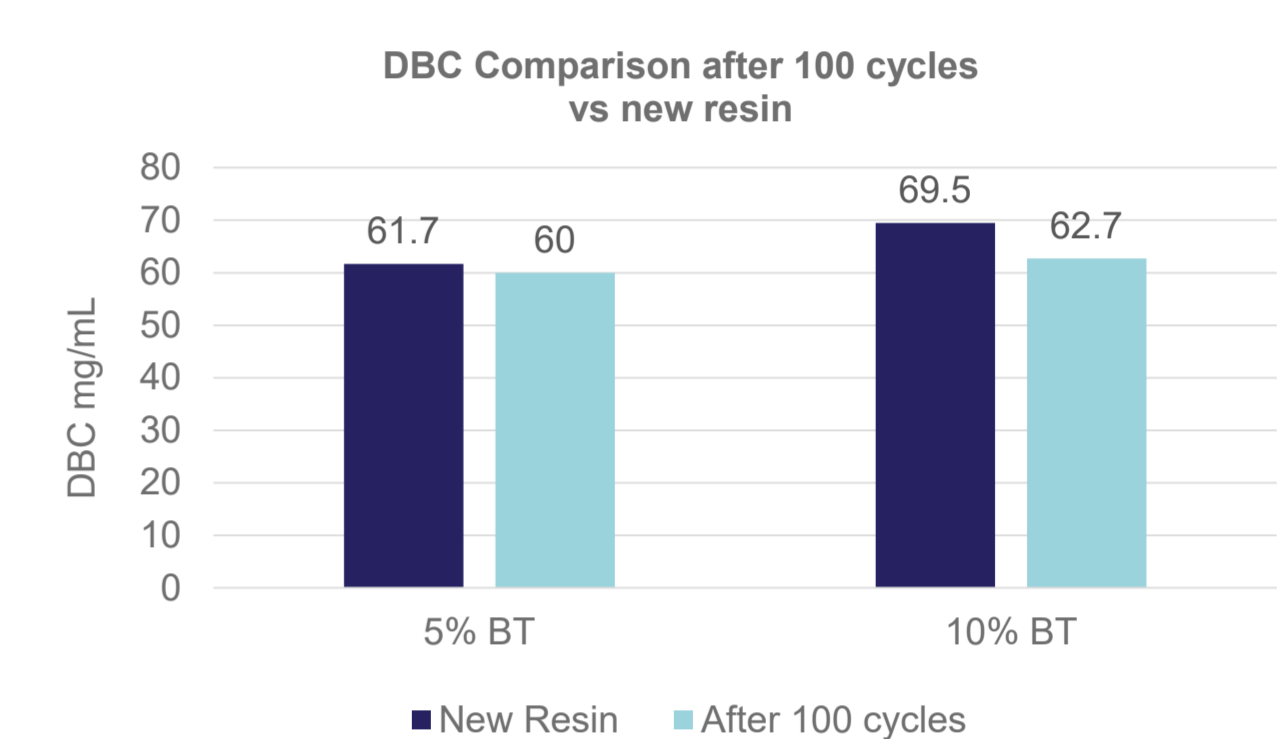


Fig 6. Excellent reusability: less than 10% capacity loss after 100 cycles

- ✓ **Excellent alkaline stability – No decline in capacity is observed after cleaning with 0.2M NaOH (2x 15 min/cycle) for 100 cycles**

#### Protein A ligand leakage

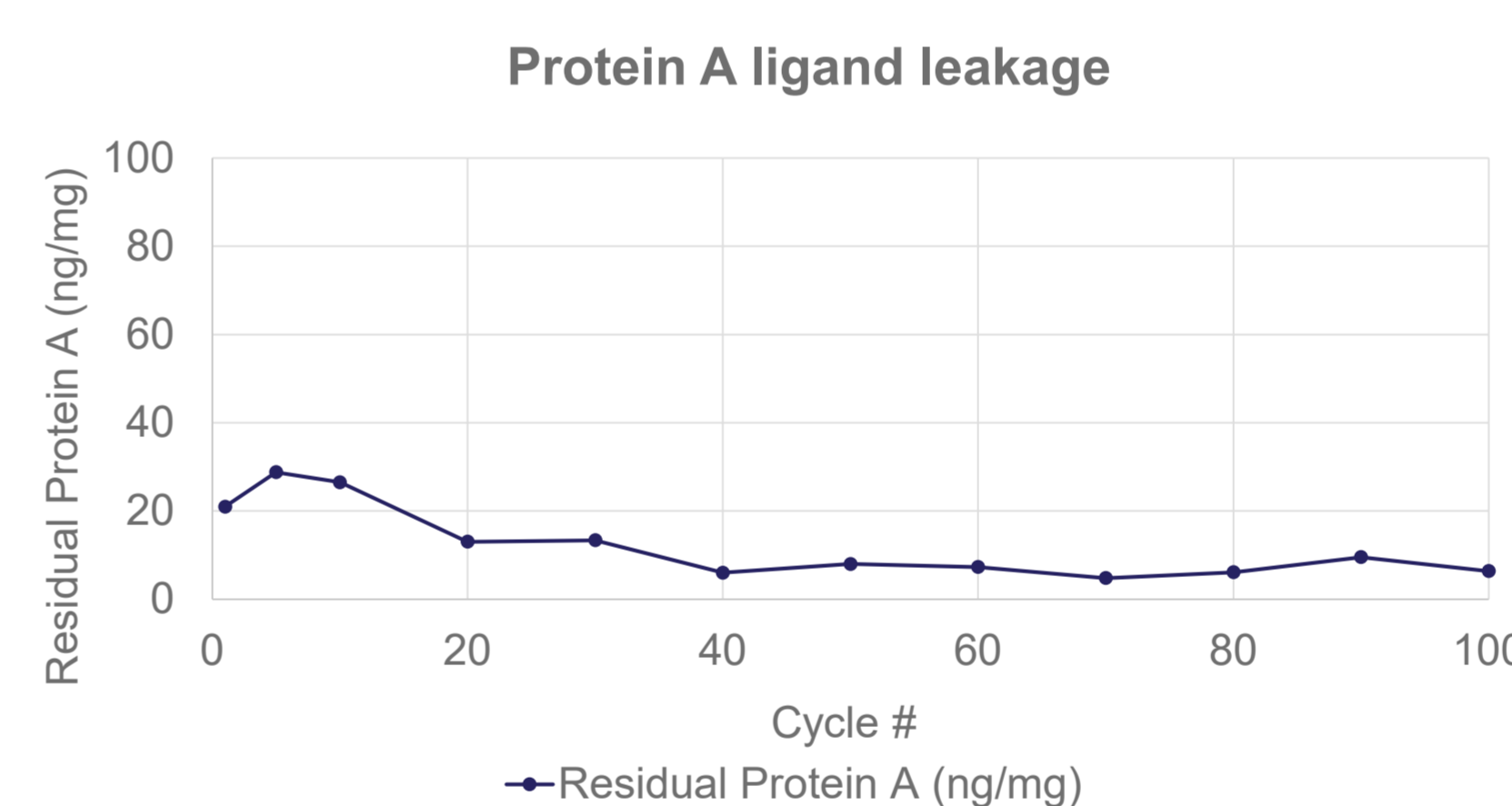


Fig 7. Ligand leakage (ppm) was measured over a 100 cycles using the Cygnus rec. Protein A ELISA kit demonstrating leakage on average of 12ppm over 100 cycles.

#### Residual DNA and HCP removal

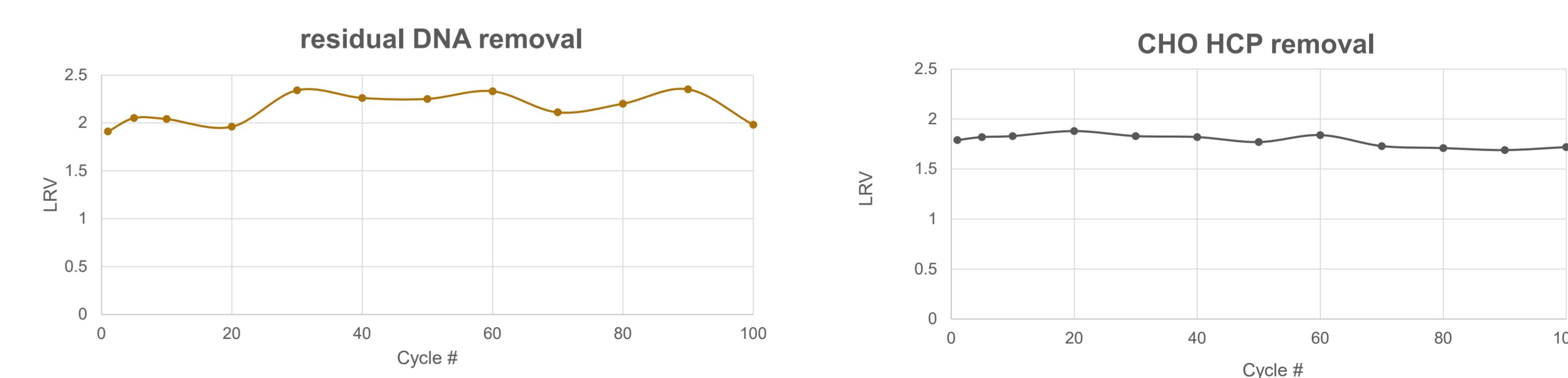


Fig 8. Consistent residual DNA and CHO HCP removal over a 100 cycles

#### Pressure-Flow characteristics

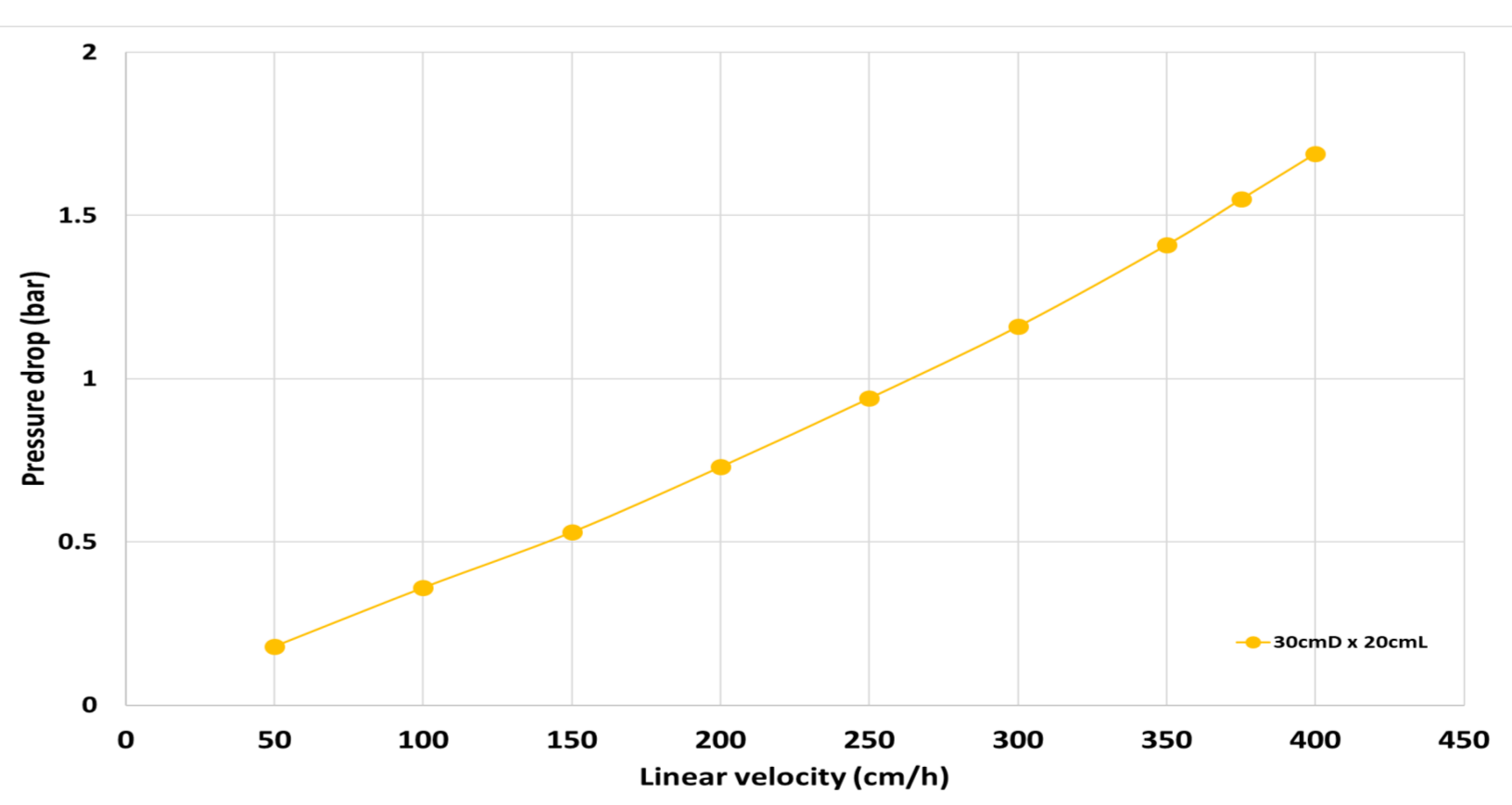


Fig 9. The MabCaptureC resin demonstrates high permeability - <2 bar at 400 cm/h in a 30 cm diameter column x 20 cm bed height

#### Available products

Cat nr.	Product
1963662250*	MabCaptureC Affinity Matrix 250ml
196366201L*	MabCaptureC Affinity Matrix 1L
196366205L*	MabCaptureC Affinity Matrix 5L
5943662001	MabCaptureC MiniChrom 1 ml
5943662005	MabCaptureC MiniChrom 5 ml
5943662200	MabCaptureC RoboColumn 200 ul
5943662600	MabCaptureC RoboColumn 600 ul

\* Products come with regulatory support (RSF)

### CONCLUSION

Featuring high capacity, excellent alkaline stability and increased productivity, the MabCaptureC affinity matrix is your protein A resin of choice in the monoclonal antibody purification workflow.