

Efficient monoclonal antibody aggregate removal by Hydrophobic Interaction Chromatography (HIC)

John J. Li, Jessica De Rooij, Alejandro Becerra, and Orjana Terova, Thermo Fisher Scientific, 35 Wiggins Ave, Bedford MA, 01730 USA

INTRODUCTION

This case study shows the optimization of POROS™ HIC resin-use in high-throughput screening (HTS) and subsequently upscaling in both Bind-Elute (B/E) and Flow-Through (FT) mode. Our study shows that a well-designed process together with a robust resin are key to a successful and efficient Mab polishing process.

GOAL OF THE STUDY

Design a more efficient, robust and cost-effective polish step utilizing POROS™ HIC resins as an alternative to the mixed-mode step in the original purification process of a clinical Mab containing >10% aggregates.



POROS HIC RESINS

POROS Ethyl

Non-POROS HIC resins

POROS Benzyl

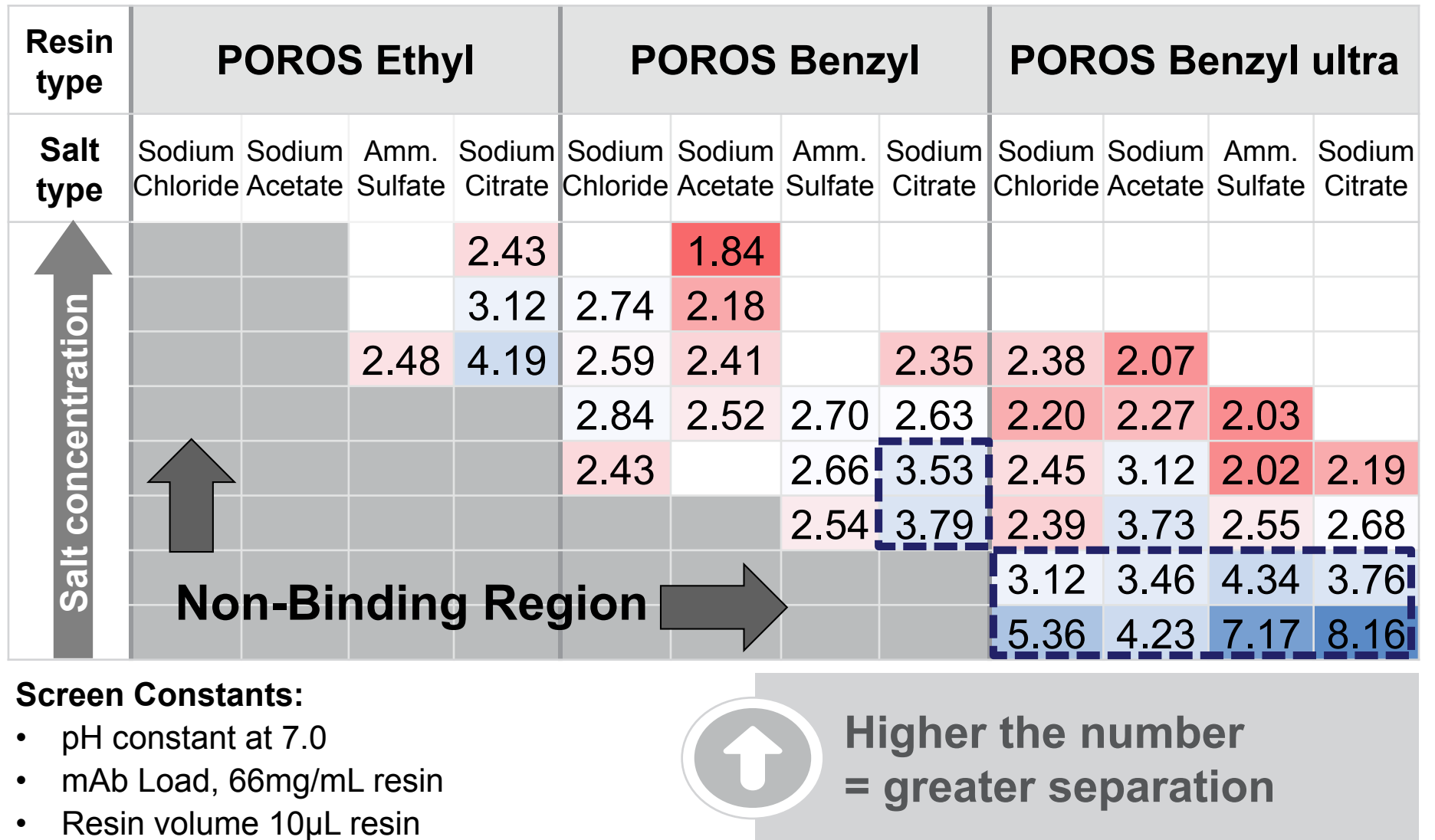
POROS Benzyl Ultra

Increasing hydrophobicity

- ✓ Designed for use with lower salt concentrations
- ✓ Differentiated selectivity and ligand chemistry
- ✓ Novel 50 µm base bead
- ✓ Improved recovery, resolution and capacity
- ✓ Superior pressure-flow characteristics
- ✓ Consistent lot-to-lot performance
- ✓ Robust chemical and base stability

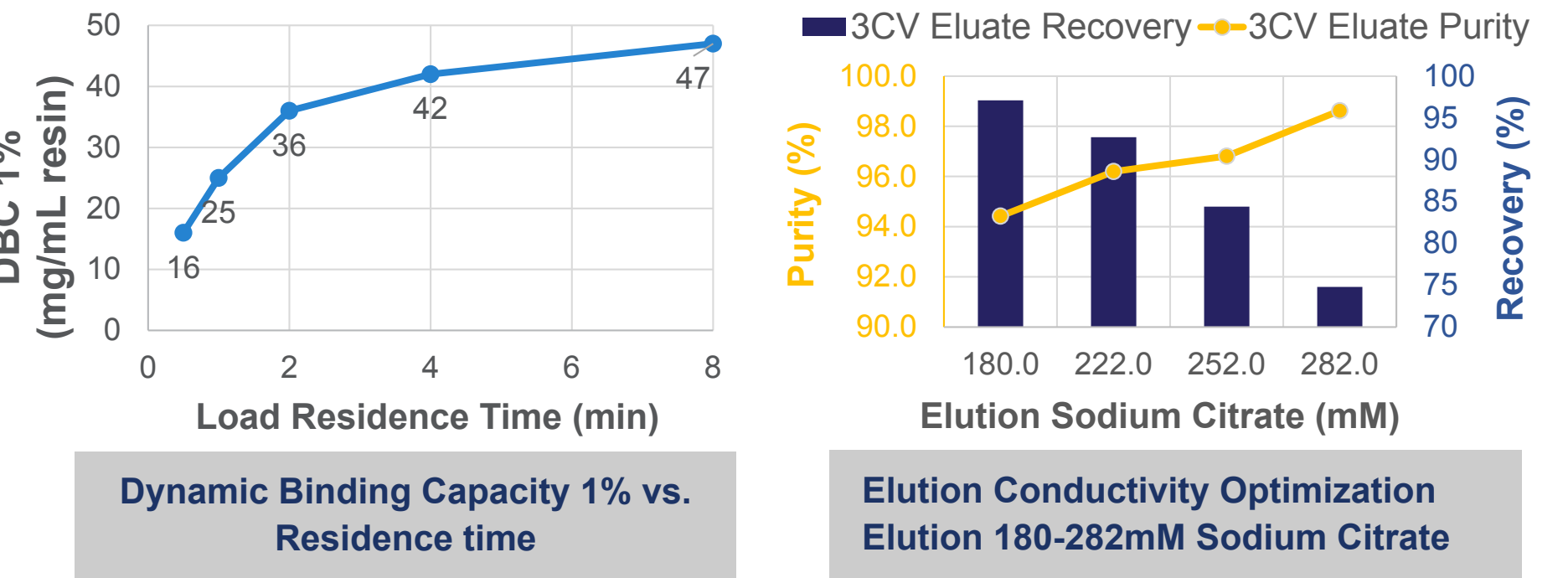
HTS: Resin Selection for Scale-Down

Screening variables used to predict conditions for scale down model: resin type, salt type and salt concentration. Partition selectivity ratio is used to determine to level of separation.



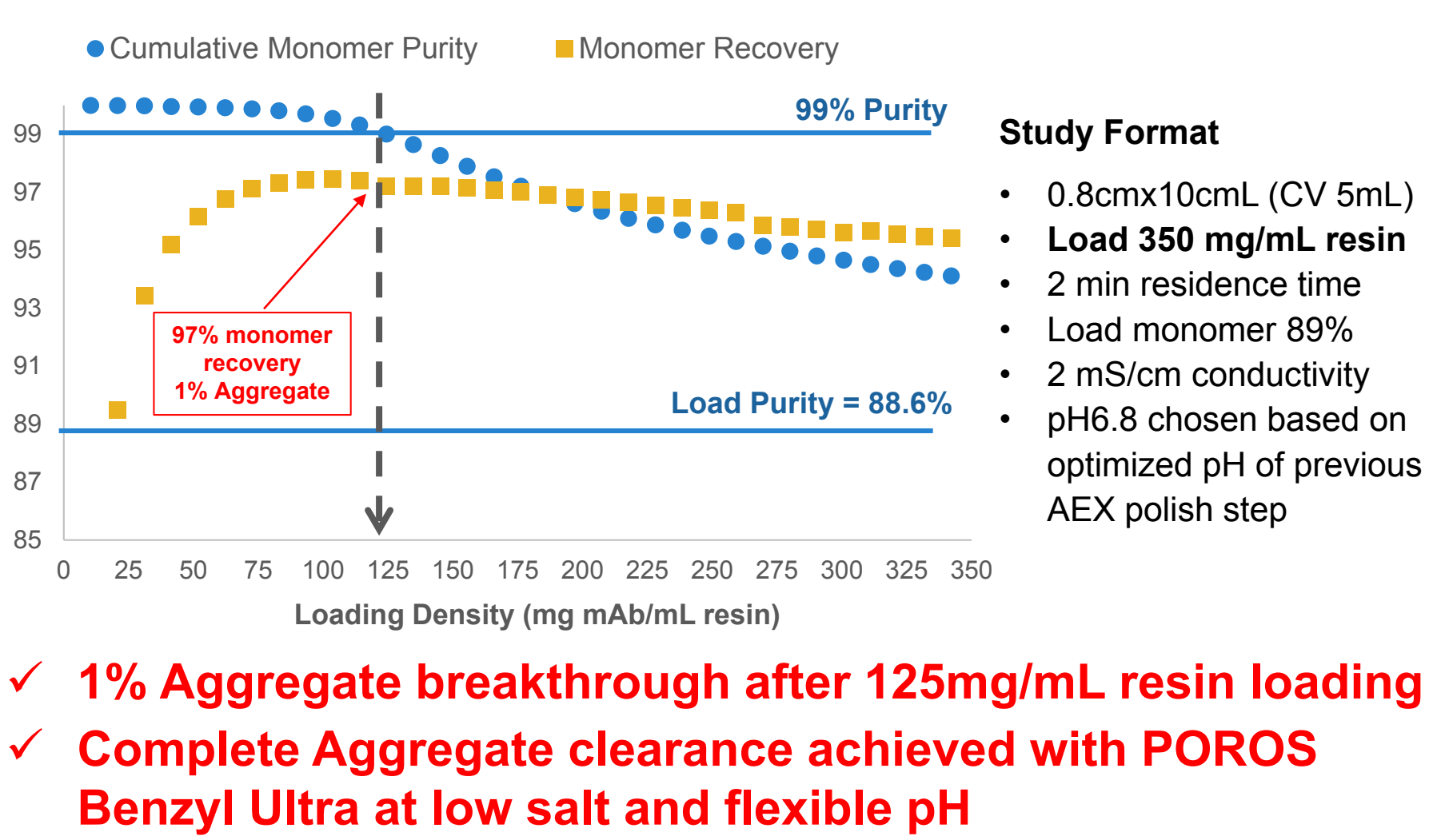
- ✓ POROS Benzyl selected for polish in bind/elute
- ✓ POROS Benzyl Ultra selected for polish in flow-through

POROS Benzyl in Bind/elute mode



- ✓ High DBC at short residence time (2min, 32mg/L resin)
- ✓ High recovery and purity in Elution (265mM Sodium Citrate)

POROS Benzyl Ultra in Flow-through mode



RESULTS SUMMARY AND CONCLUSIONS

Process Summary	Mixed-Mode (Clinical Process)	POROS Benzyl Bind-Elute Mode	POROS Benzyl Ultra Flow-through Mode
Load Monomer Purity (%)	90	89	85.5
Load Density (g/L resin)	25	32	100
Monomer Purity Pool (%)	99	99	>99
Monomer Recovery (%)	90	>99	98
HCP (ppm)	NA	120 to 12ppm	100 to 35 ppm
Residence time (min)	6	2	1.2
Pool Volume (50-50mAu)	5CV	4CV	NA
MMV Clearance	NA	2	1
XmuLV Clearance	NA	>5	>5
Productivity (g/L/hr)	7	27	89

POROS HIC resins drastically improve Mab polish step:

- ✓ Increased load density
- ✓ Improved monomer recovery
- ✓ Shorter residence time
- ✓ 4-12 times higher process step efficiency