

Advantage of antibody based selectivity in the purification of biologics



Pim Hermans¹, Frank Detmers¹, Hendrik Adams¹, Orjana Terova² and Jessica de Rooij¹
1. Thermo Fisher Scientific, Leiden, the Netherlands
2. Thermo Fisher Scientific, Bedford, MA

Bioprocessing

INTRODUCTION

Advances in bio-therapeutics are generating an increasing range of complex molecules that present unique and often complex purification challenges. By taking advantage of antibody based selectivity, Camelid heavy-chain antibody fragments (V_HHs) have proven to be a reliable immuno-affinity chromatography (IAC) solution in the downstream process of biologics. Thermo Scientific™ CaptureSelect™ affinity products and analytical tools are developed for the discovery and manufacturing of even the most demanding bio-therapeutics. The affinity resins provide high target purity in a single step, independent of feedstock.

CAPTURESELECT TECHNOLOGY – UNIQUE AFFINITY PURIFICATION SOLUTION

- Affinity through antibody selectivity: technology based on Camelid single domain [V_HH] antibody fragments
- Unique V_HH screening technology to determine final resin properties such as target specificity, mild elution & ligand stability
- Animal origin free production process (*Saccharomyces Cerevisiae*)
- Technology used in commercial purification processes

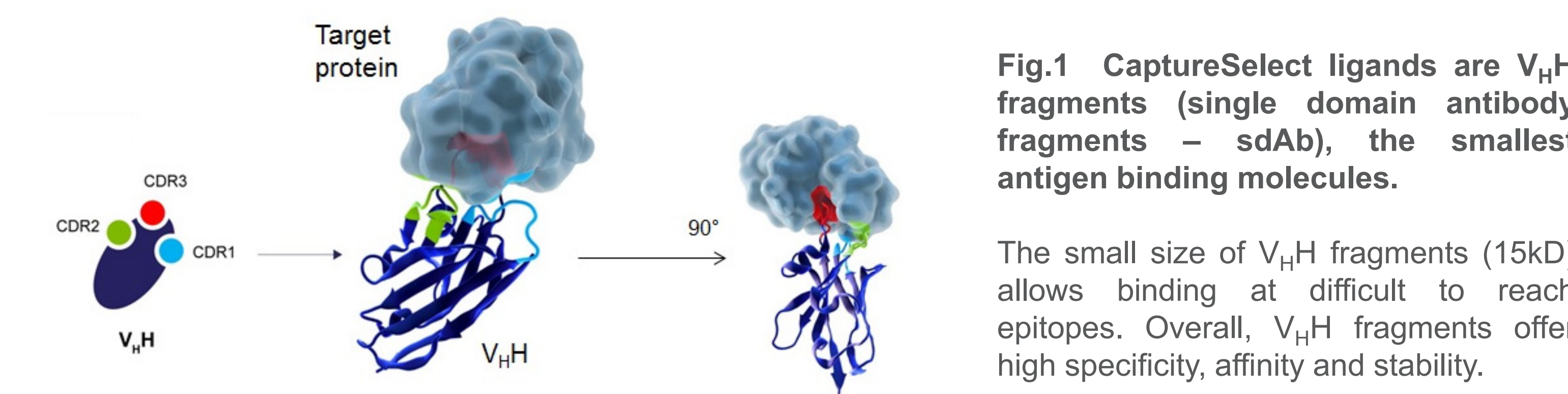


Fig.1 CaptureSelect ligands are V_HH fragments (single domain antibody fragments – sdAb), the smallest antigen binding molecules. The small size of V_HH fragments (15kD) allows binding at difficult to reach epitopes. Overall, V_HH fragments offer high specificity, affinity and stability.

LARGE SCALE PURIFICATION OF AAV VECTORS

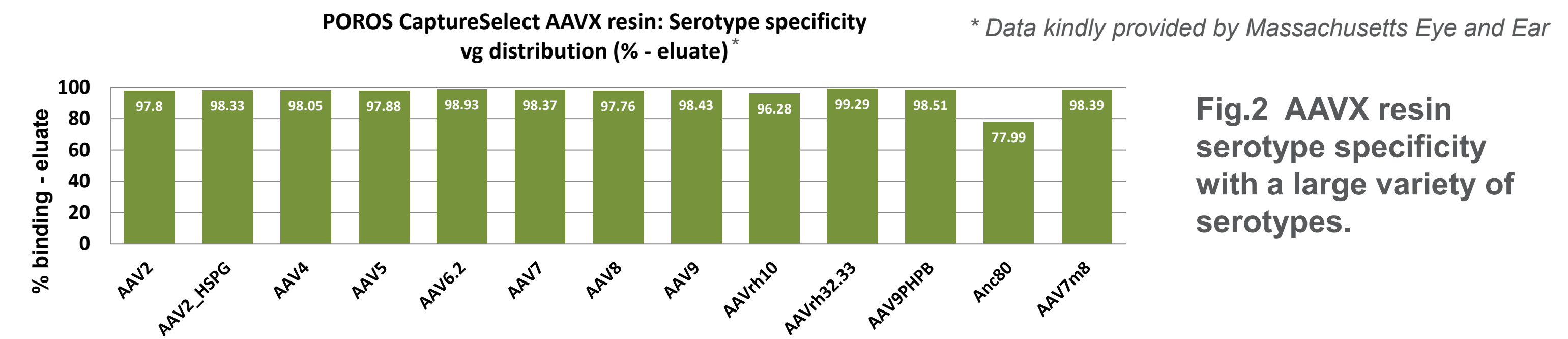
- Increased yields through the reduction of chromatography steps in the viral vector purification process
- High purity and recovery in a single purification step
- Scalable purification of multiple AAV serotypes

Thermo Scientific™ resin	Binding Capacity (vg/mL)	Serotype Affinity
POROS™ CaptureSelect™ AAV8	>10 ¹³	AAV8
POROS™ CaptureSelect™ AAV9	>10 ¹⁴	AAV9
POROS™ CaptureSelect™ AAVX	>10 ¹⁴ *	AAV1, AAV2, AAV3, AAV4, AAV5, AAV6, AAV7, AAV8, AAV9, recombinant & chimeric vectors

* viral genomes per millilitre (vg)/mL, binding capacity will vary based on serotype, feed stream, additives, and mutations to parent serotypes

POROS CAPTURESELECT AAVX: A PLATFORM FOR AAV PURIFICATION

- Broad selectivity to both natural and synthetic capsids
- High dynamic binding capacity
- High elution recovery at different flow rates
- Robust, with less process optimization steps



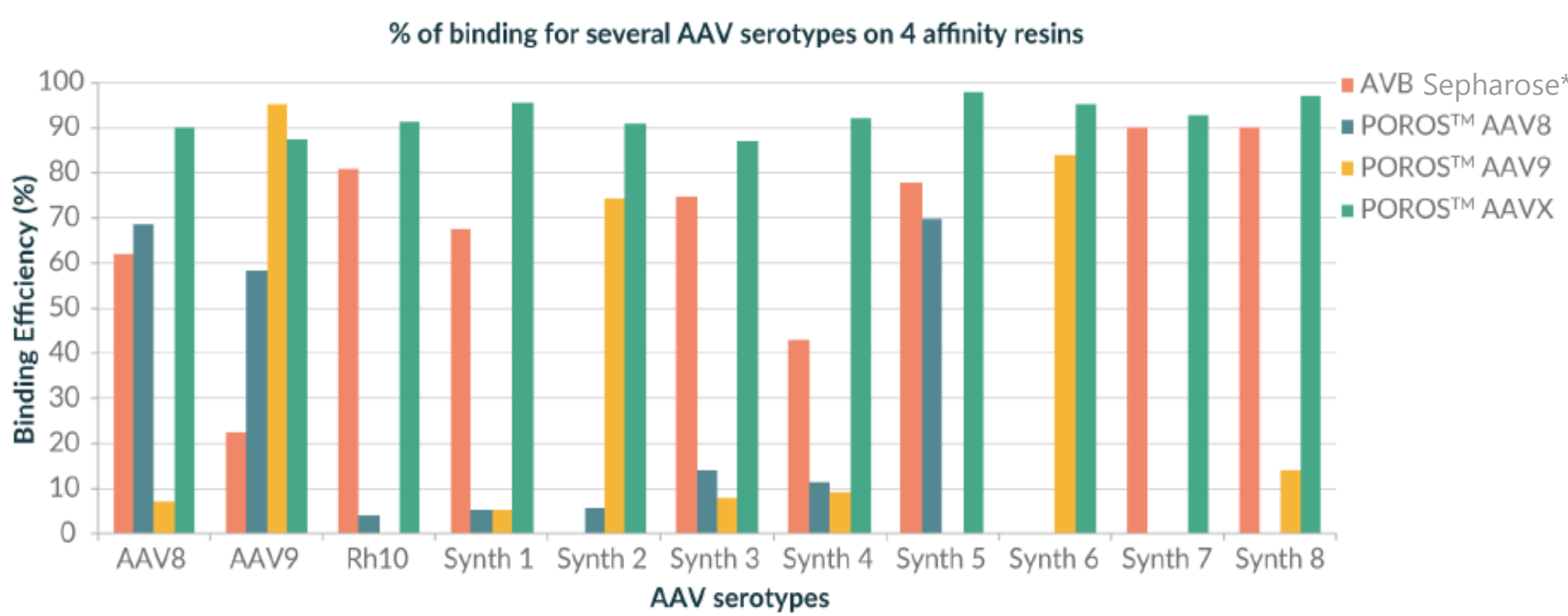
* Data kindly provided by Massachusetts Eye and Ear

Fig.2 AAVX resin serotype specificity with a large variety of serotypes.

Experimental settings. Static binding mode experiment: resin was mixed with AAV serotype in tube – no flow properties performed. VG was determined by qPCR

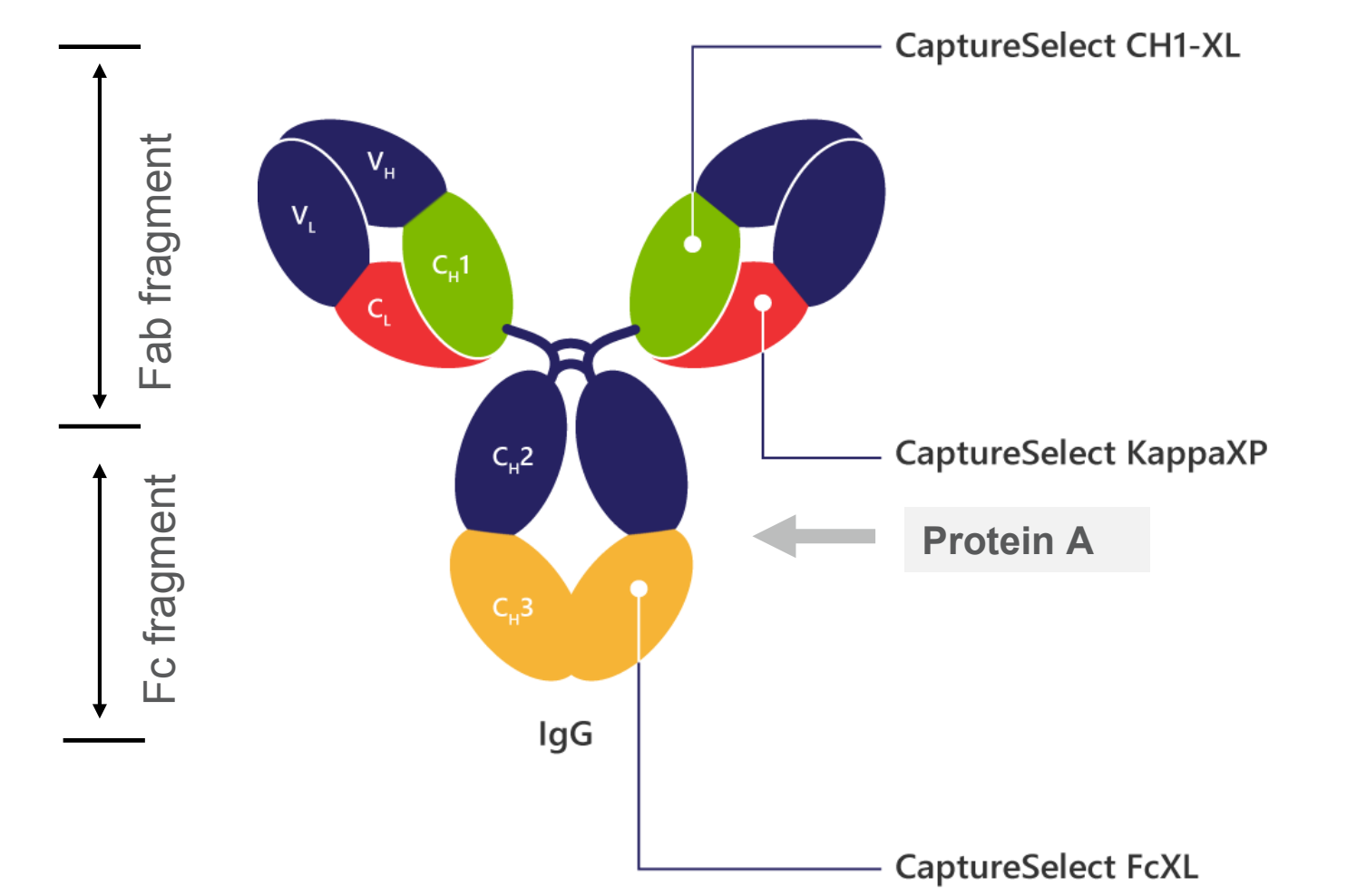
CENETHON The following data has been kindly provided by Genethon and published in Cell Gene Therapy Insights 2018; 4(7), 637-645.

Fig.3 Comparison of POROS CaptureSelect AAV resins with an alternative resin. 3 native AAV vectors (AAV8-10) and 8 synthetic vectors were tested in static binding mode with the 4 different resins.



- ✓ The use of AAV affinity resins simplifies processes and can reduce costs by a factor of 6 (customer testimonial)

PURIFICATION OF ANTIBODY THERAPEUTICS



A unique set of CaptureSelect affinity ligands has been developed (fig 4.), directed against a variety of antibody subdomains, providing tools for researchers and manufacturers to help facilitate purification of a broad range of antibody formats.

Fig.4 CaptureSelect resins: Antibody Binding Selectivity Binding regions of CaptureSelect resins for affinity purification of antibodies and antibody fragments.

ANTIBODY PURIFICATION – CAPTURESELECT CH1-XL

- CH1 binding domain ligand
- No co-purification of free light chains (only correct assembled Fabs)
- Efficient elution at milder pH (4 – 4,5)

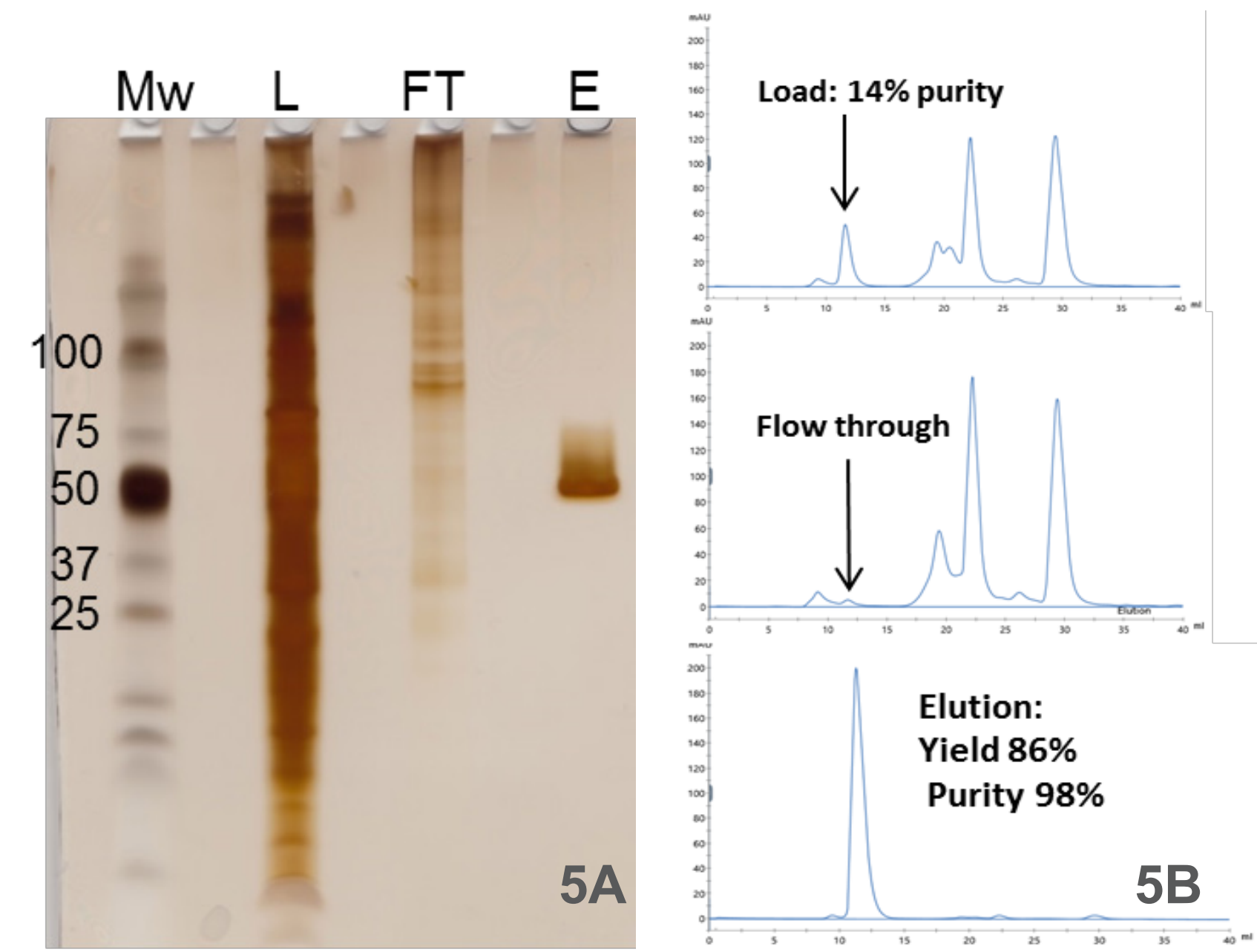


Fig.5 Ranibuzimab feed from HEK293 cells. Analysis of the fractions after purification with CaptureSelect CH1-XL resin shows high yield and purity in a single step.

5A: SDS-PAGE silver staining of the load (L), flow through (FT) and elution (E) fractions, showing no presence of light chains in the elution pool.

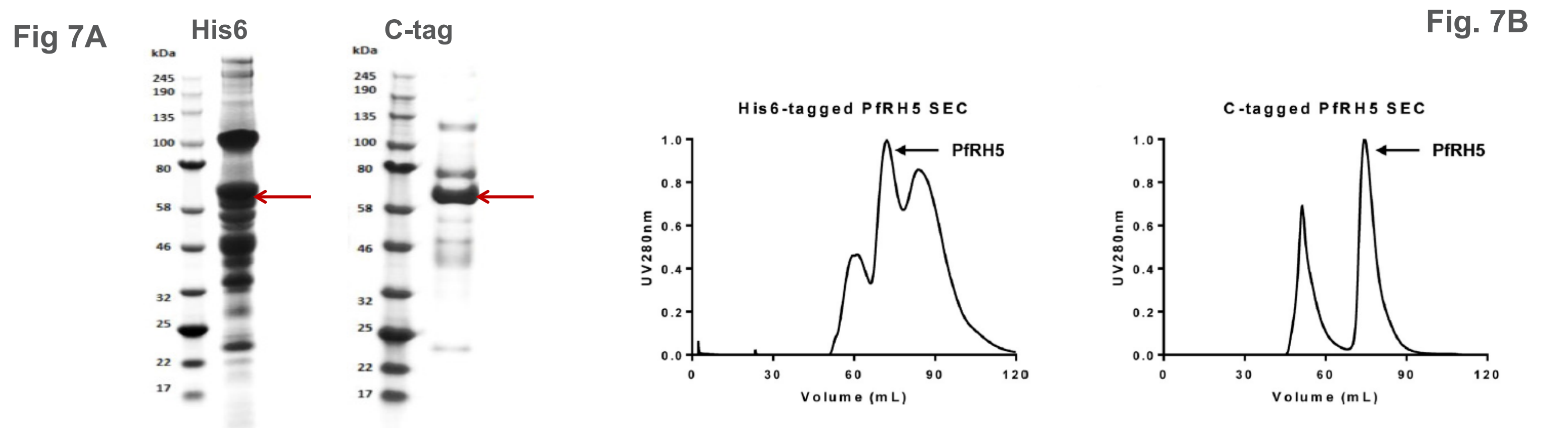
5B: Gel filtration analysis showing 98% purity of the Fab fragment in the elution fraction with a yield of 86%

- ✓ A true platform for Fab fragment purification

C-TAG – A REVOLUTIONAIRY AFFINITY TAG

- Unique affinity tag based on a 4 residue C-terminal peptide sequence: E-P-E-A
- Limited effect on protein functionality unlike larger tags 9GST, HIS6)
- Enabling high purity and yield from complex mixtures in a “one-step” process
- Mild elution, protecting the protein of interest

cGMP PURIFICATION OF THE PfRH5 PROTEIN-BASED MALARIA VACCINE



Process yield (after)	His6-tagged construct	C-tagged construct
Culture supernatant	100%	100%
Tangential Flow Filtration	82.1%	91.0%
Affinity Chromatography	52.5%	77.4%
Size Exclusion Chromatography	25.5%	43.3%
Overall purity	85-90%	>99%

Fig. 7 Purification of a recombinant malaria vaccine (PfRH5) from insect cells feedstock using a C-terminal fused His6-tag or C-tag.* A. Purity assessment: Improved purity compared to hexa-histidine tag purification B. UV280 absorbance chromatograms after Size Exclusion Chromatography (left His6, right C-tag) C. Process yields after each purification step. C-tag clearly outperforms His6-tag (table)

*Data Obtained from: Jin, J., et al., 2017. Accelerating the clinical development of protein-based vaccines for malaria by efficient purification using a four amino acid C-terminal 'C-tag'. Int J Parasitol. 47(7), 435-446

- ✓ The first affinity tag to be left on a therapeutic protein, approved to enter clinical trials

TRADEMARKS/LICENSING

Caution: For Research Use or further manufacturing, not for diagnostic use or direct administration in humans or animals. © 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.