



A scalable AAV production process from cell culture to purified bulk

Webinar about AAV workflow and production

Junichi Inagawa, Application sales, BioProcess
Cytiva-Japan





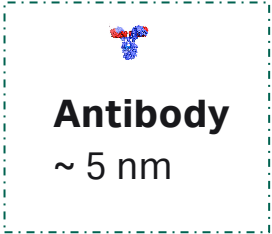
1. Introduction
2. Upstream cell culture and virus production
3. Downstream purification
4. Conclusions and next steps



1

Introduction

Sizes of common viruses



Preventive vaccines

- JE** (40–60 nm)
- Yellow Fever** (40–60 nm)
- HAV** (30 nm)
- VLP** (40–60 nm)
- Polio** (30 nm)
- Hep B** (42 nm)

Rota
(80 nm)

Flu
(90–120 nm)

Corona
(125 nm)

Mumps (200 nm)

Measles (100 × 300 nm)

Rabies (75 × 180 nm)

Recombinant virus vectors for cell and gene therapy

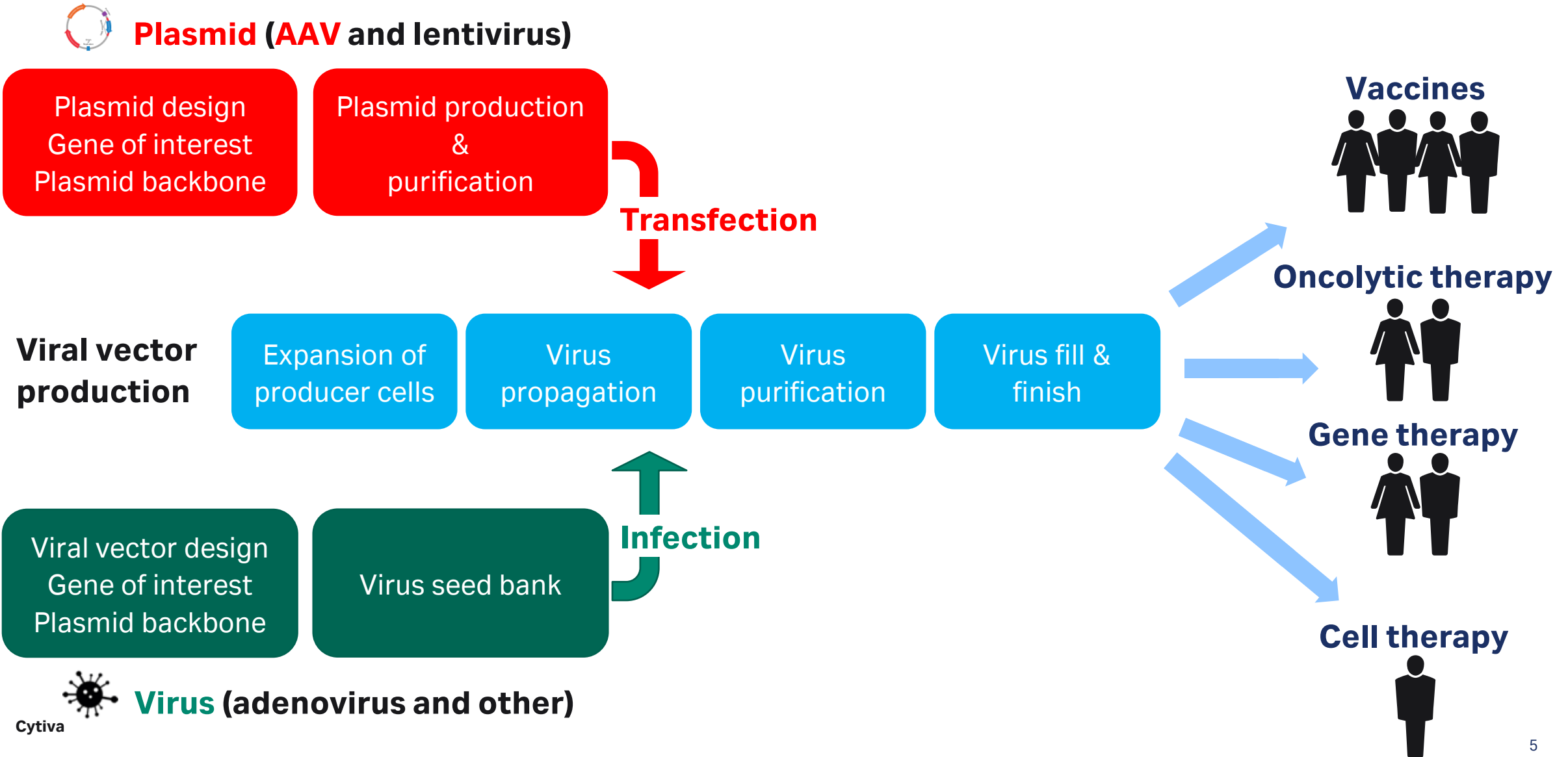
AAV
(25 nm)

Adeno
(70–90 nm)

Lenti
(80–120)

POX
(200–300)

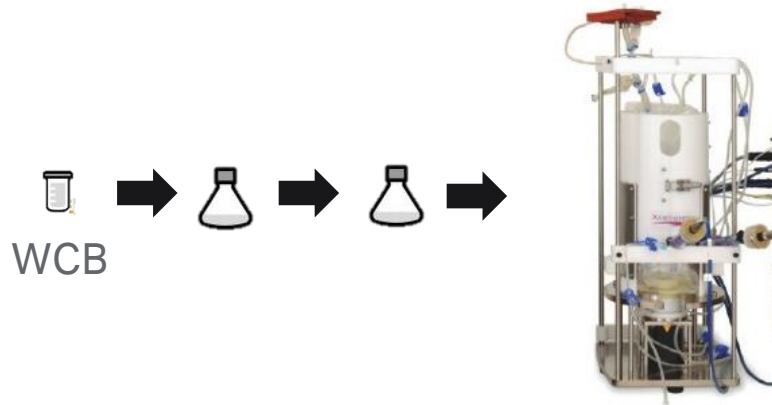
Viral vector production and clinical use



Adeno-associated virus (AAV) production process

Upstream

- Triple plasmid transfection
- HEK293T suspension
- AAV2-GFP



Xcellerex™ XDR-10
Scaleable to 2000 L

Seed train

WCB = working cell bank

Cytiva

Virus production



Downstream

Cell lysis
DNA fragmentation

Clarification

Concentration &
buffer exchange

Capture

Polishing

Concentration &
buffer exchange

Sterile filtration

Analysis

Virus infectious titer

Transduction assay: flow cytometry

Virus titer

Viral genomes: qPCR

Viral capsids: ELISA, SPR (Biacore™)

Full-empty ratio: qPCR/ELISA,
Analytic IEX, TEM

Host cell impurities

Total protein: BCA assay

Total DNA: Picogreen™ Assay

HC DNA: qPCR

HCP: ELISA

Characterization

SDS-PAGE, Western blotting

TEM SEC and IEX HPLC

2

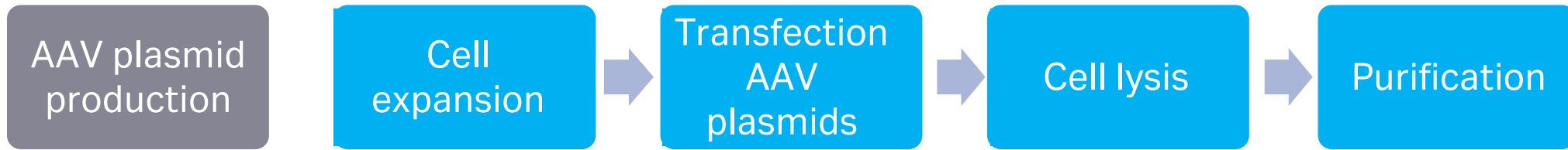
Upstream cell culture and virus propagation

AAV upstream strategy

- Adaptation to serum free suspension cell culture
- Cell culture medium evaluation
- Cell density optimization
- Optimization of AAV transfection
- Optimization of culture conditions post transfection



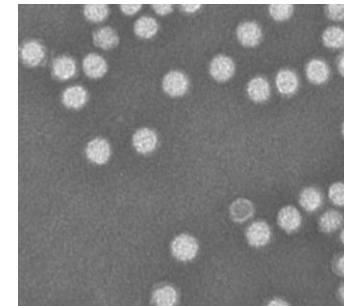
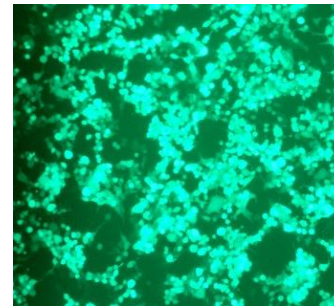
HyClone™
HyCell
TransFect-H



Cell

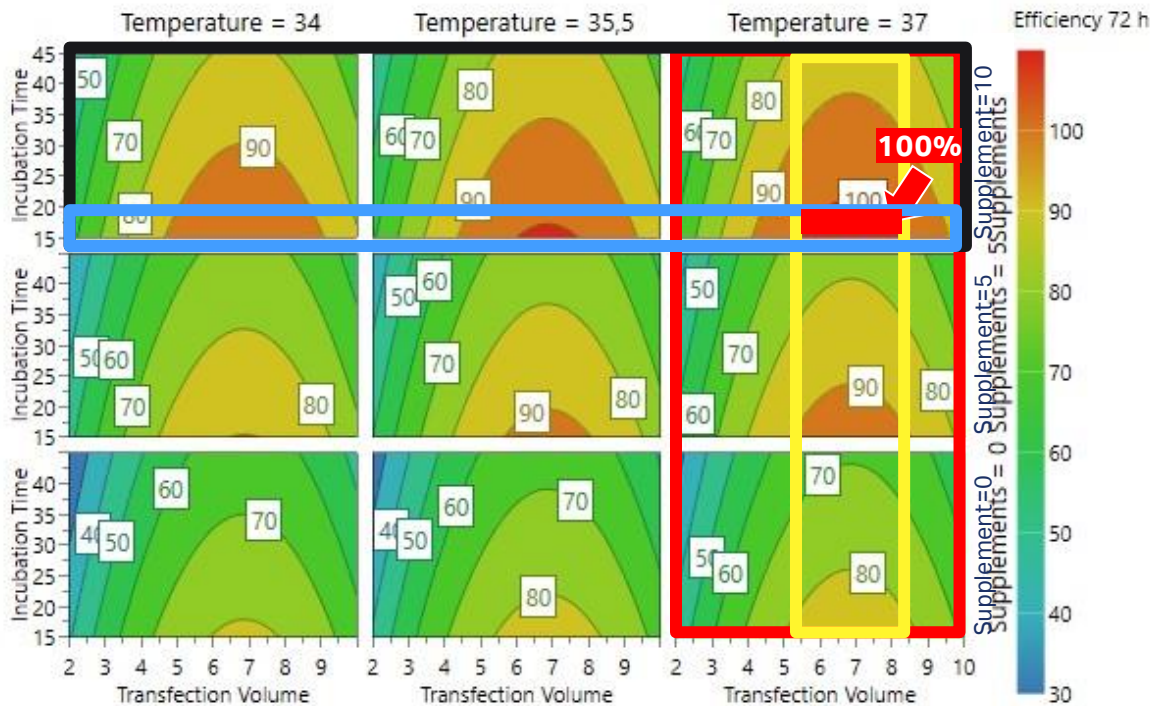


XDR bioreactor
(scalable up to 2000 L)



DoE results

Transfection efficiency



Transfection efficiency: % GFP expressing cells

Criteria:

Titer (ELISA) > 10^9 capsids/mL

Optimized transfection protocol

VCD: 1×10^6 /mL

DNA ($\mu\text{g}/\mu\text{L}$): 1

PEI/DNA ratio: 2

Transfection volume (% of total): 5

Incubation time: 20 minutes

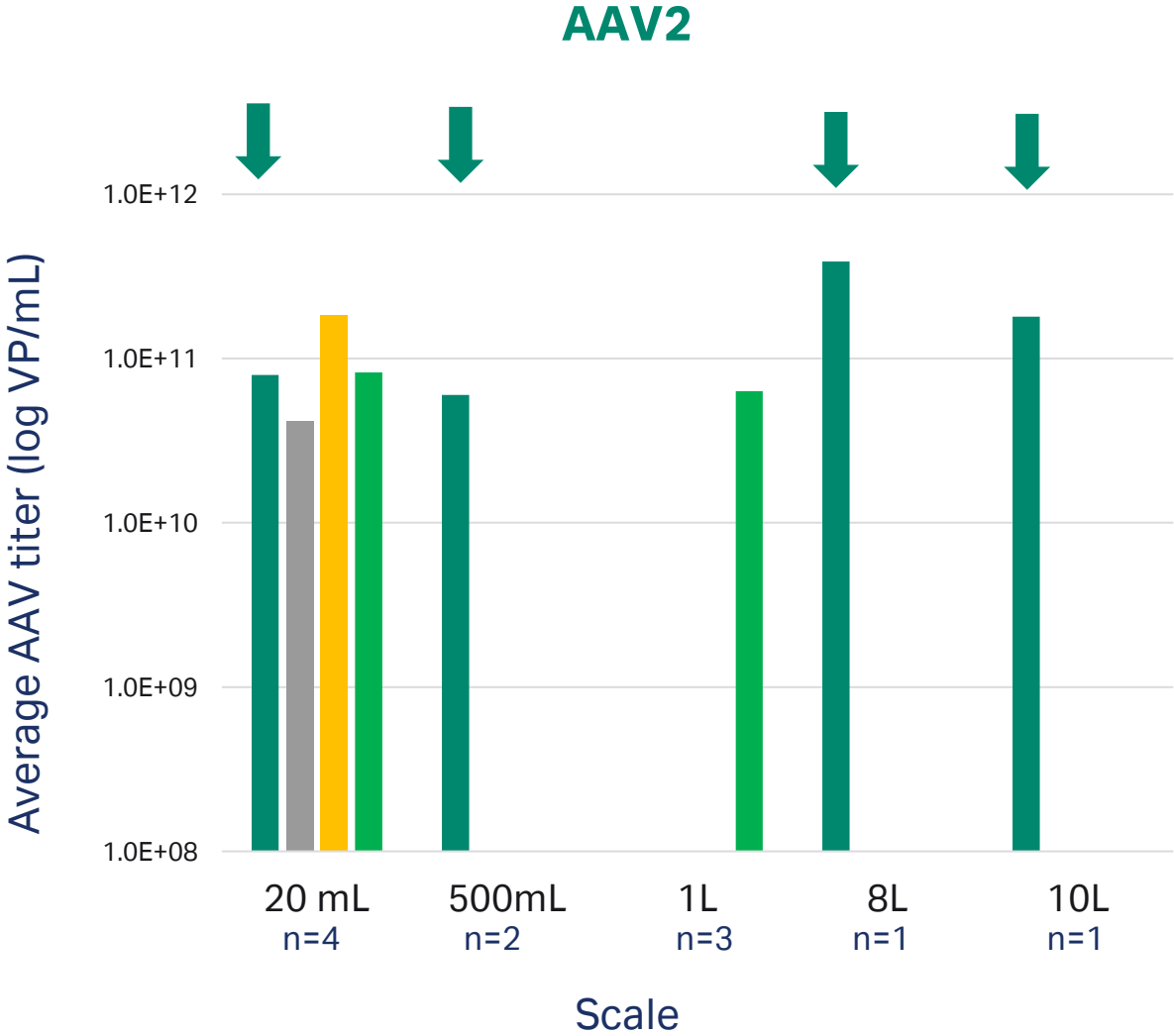
Temperature: 37°C

DNA ratio: 1:1:2

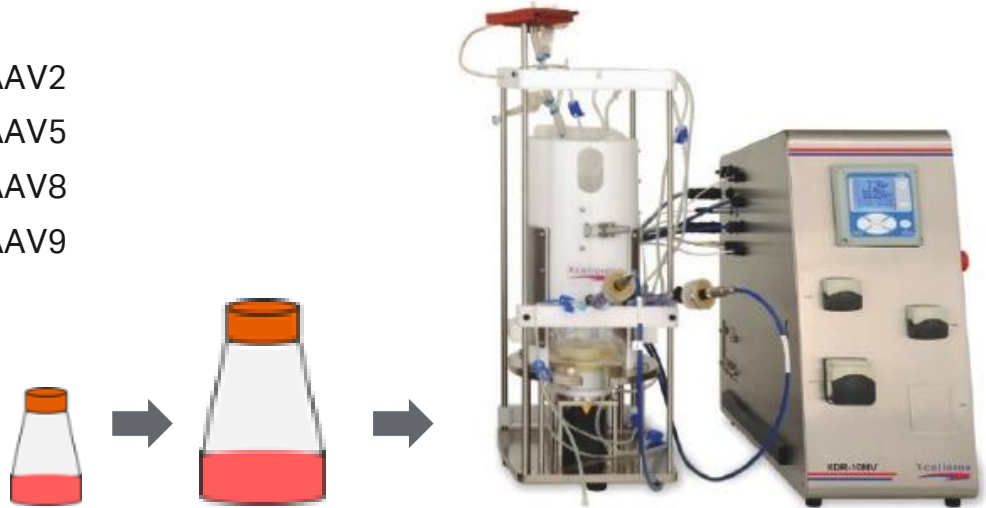
(Rep/cap: helper: transgene plasmid)

Sodium Butyrate: 5 mM

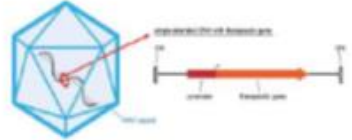
AAV productivity in different scales



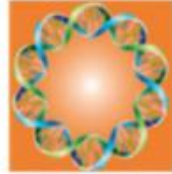
- AAV2
- AAV5
- AAV8
- AAV9



Homology's Upstream HEK293 Transfection Production Platform Shows Linear Scalability Up to 500 L



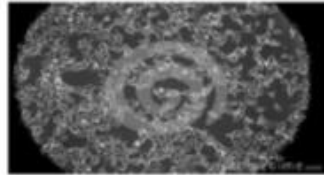
1. Construct Selection



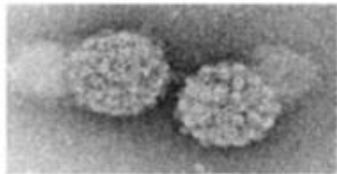
2. Plasmid Engineering & fine tuning



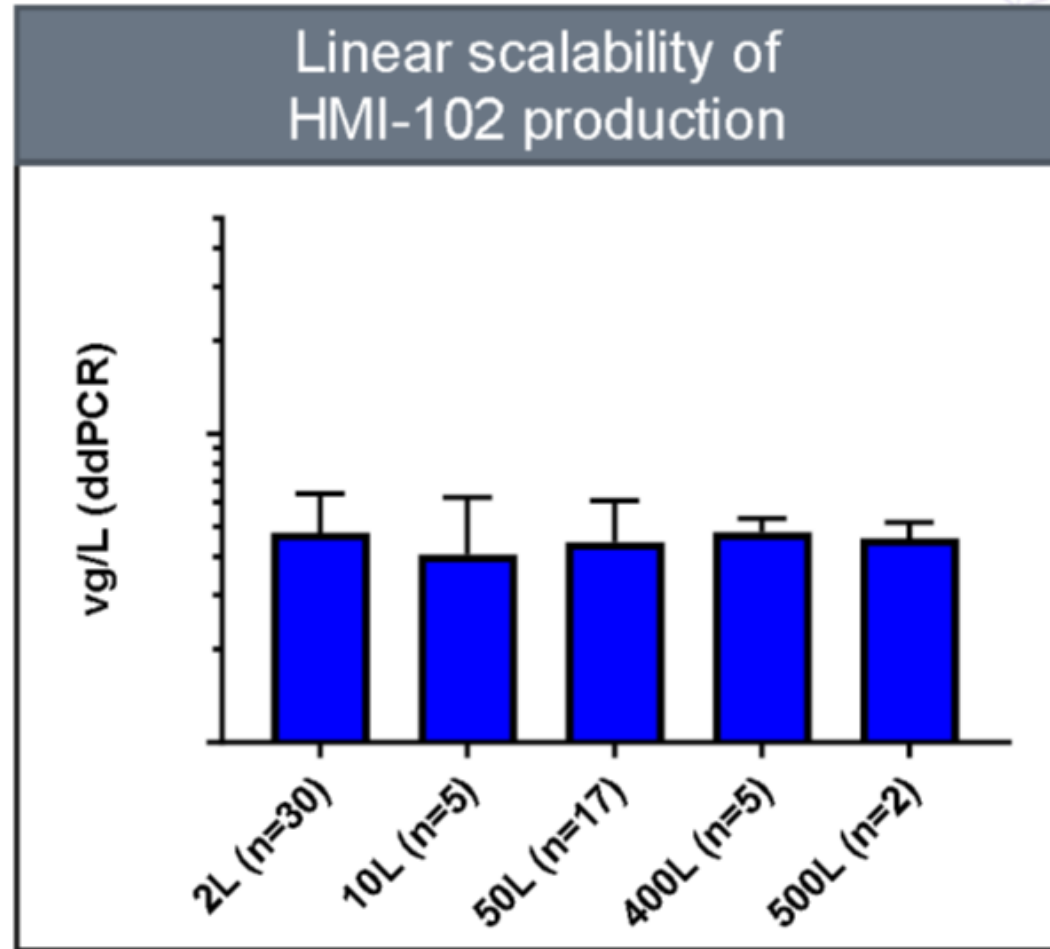
3. Cell growth & propagation



4. Transfection



5. Harvest product frozen & stored



3

Clarification Downstream purification

Harvest: Cell lysis, DNA fragmentation and Clarification



**Cell lysis
DNA fragmentation**

Clarification

Conc. and buffer exchange

Capture

Polishing

Conc. and buffer exchange

Sterile filtration

Harvest:

0.5 % Tween™ 20
300 mM NaCl
1 mM MgCl₂
40 U/mL Denarase™ (DNA nuclease)
Incubation in Bioreactor at 37°C with mixing for 4 hours

Normal flow filtration:

ULTA™ capsules 5 μm + 2 μm + 0.6/0.2 μm HC
Flow 30 to 50 LMH
Recovery 74 to 80 %



Concentration and buffer exchange: tangential flow filtration hollow fiber with 300 kDa NMWCO



Downstream

Cell lysis
DNA fragmentation

Clarification

Conc. and buffer exchange

Capture

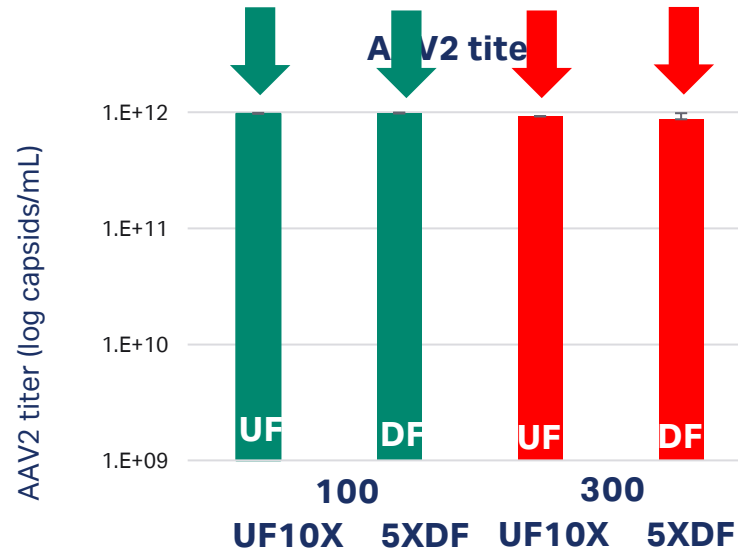
Polishing

Conc. and buffer exchange

Sterile filtration

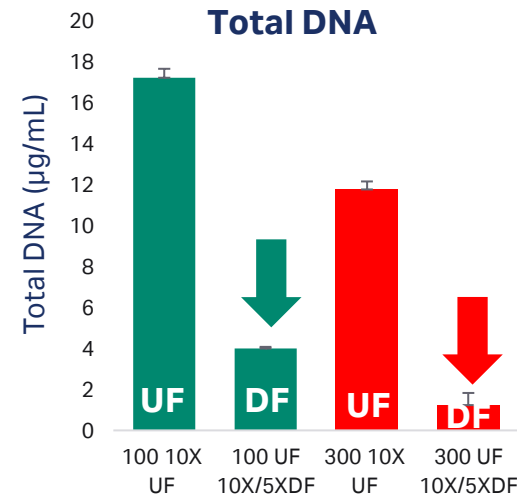
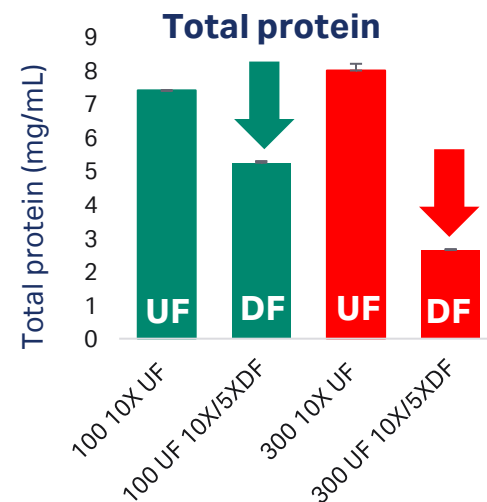
UF= Ultrafiltration
DF= Diafiltration
AIEX = Anion exchange chromatography

Cytiva



No virus losses using 300 kDa NMWCO
Recovery was 75 to 80 %

100 NMWCO UF10X/5XDF
300 NMWCO UF10X/5XDF



Better impurity removal using 300 kDa NMWCO

Capture: Affinity chromatography with **Capto™ AVB**



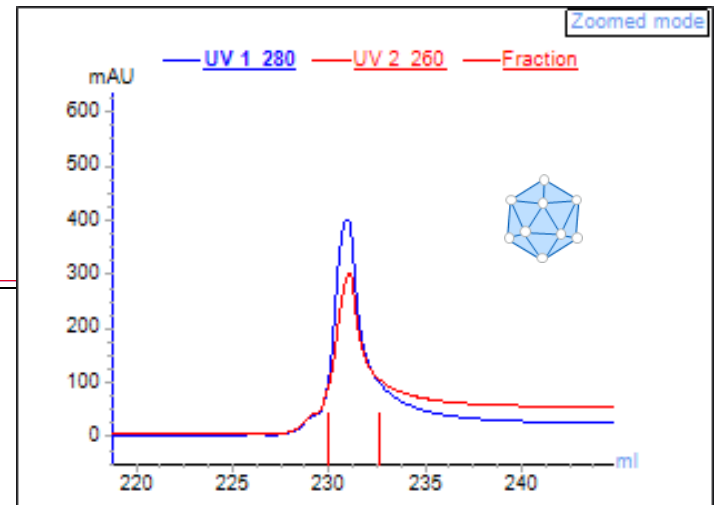
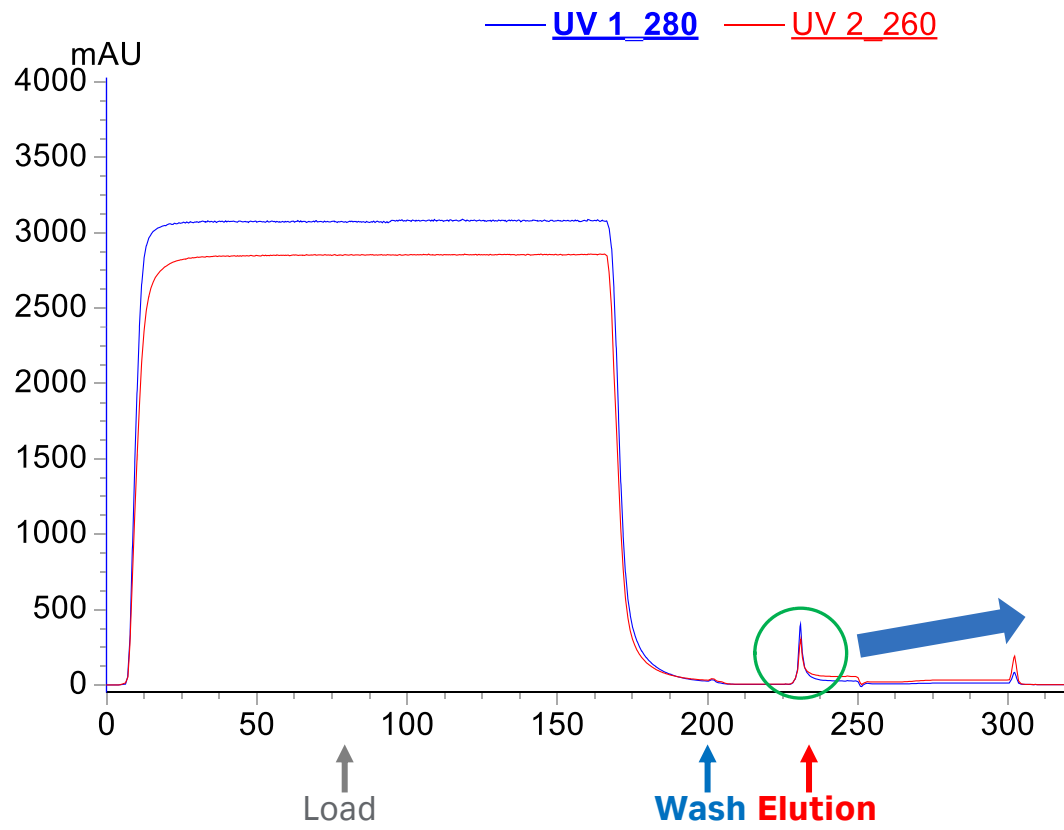
ÄKTA™ pure 25
HiTrap™ Capto AVB 5mL

Load flow: 5 mL/ min (1 min RT)
Eluate flow: 2.5 mL/min (2 min RT)

Eq buffer: 20 Mm Tris, pH 8.0 + 500 mM NaCl
Washing buf. ①50 Mm Tris, pH 8.0 + 300 mM NaCl
②20 Mm Tris pH 8.0
Elution buf: 50 mM Citrate pH 3.5 500mM NaCl
500 mM Arginine

Binding capacity: ~ 1 x 10¹⁴ capsids/mL resin
Concentration factor: ~ 100×
Recovery range: 60 to 80 %

- Cell lysis
DNA fragmentation
- Clarification
- Conc. and buffer exchange
- Capture**
- Polishing
- Conc. and buffer exchange
- Sterile filtration



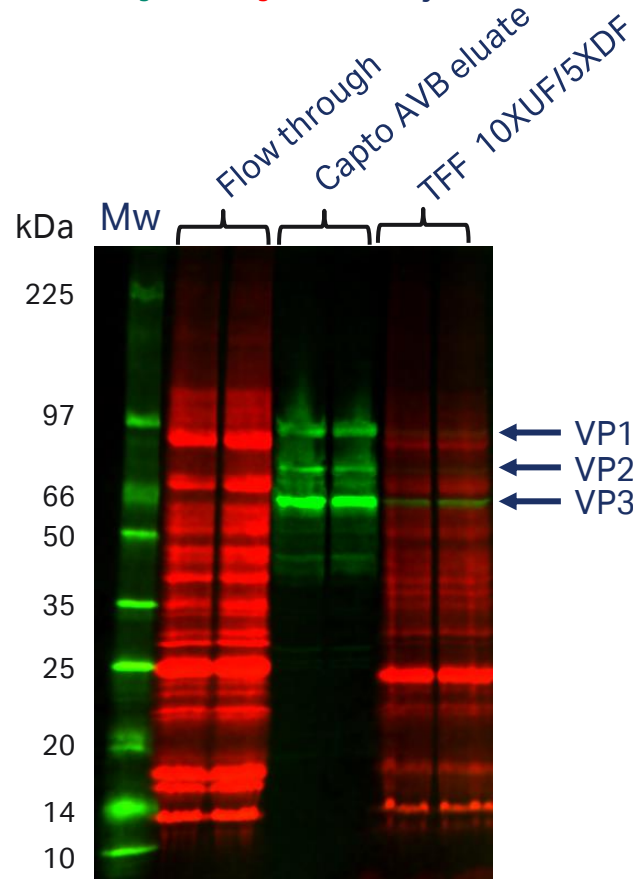
Capture: Affinity chromatography with **Capto™ AVB**



Downstream

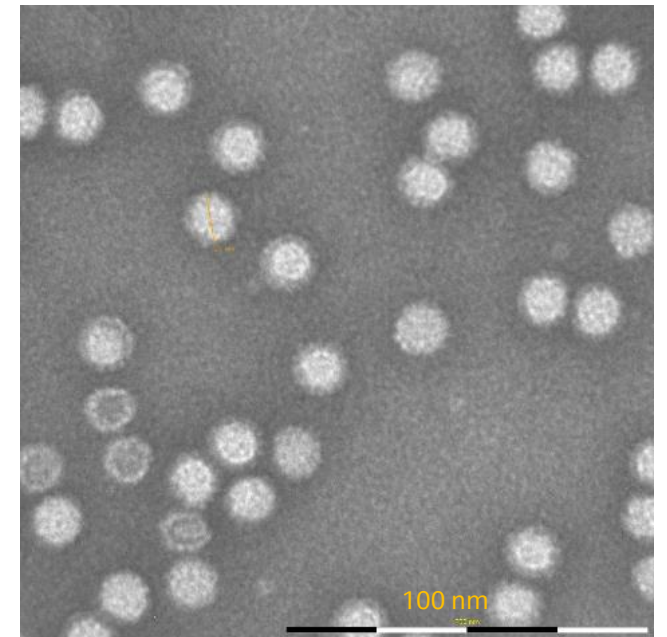
- Cell lysis
DNA fragmentation
- Clarification
- Conc. and buffer exchange
- Capture**
- Polishing
- Conc. and buffer exchange
- Sterile filtration

Membrane image
Cy™3/Cy5 overlay



Red = HCP
Green = Viral proteins

Capto AVB eluate

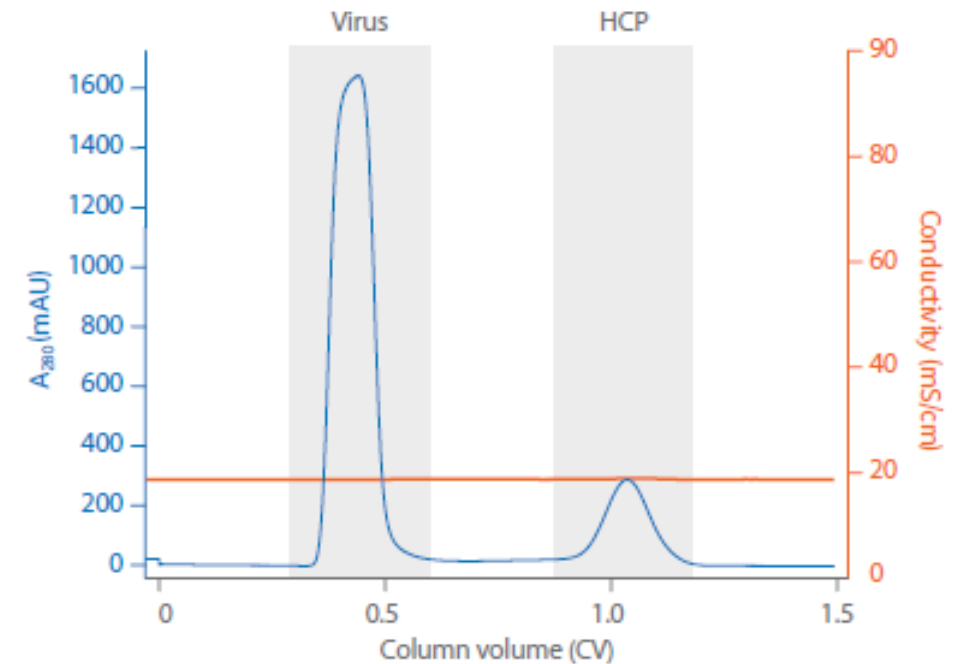
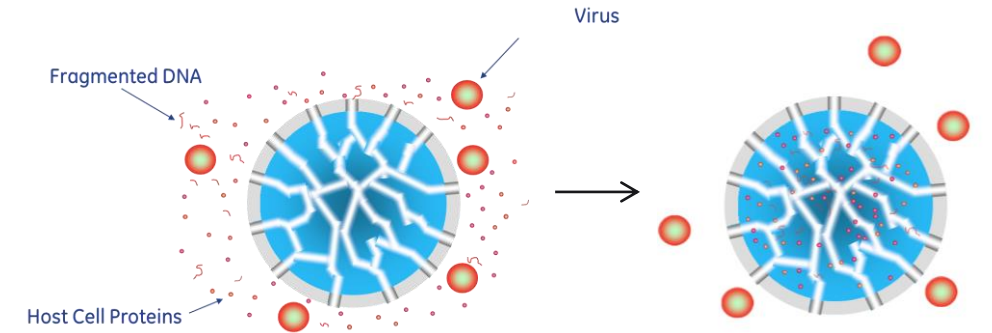


Transmission electron microscopy in collaboration with Vironova AB using MiniTEM™ system, Stockholm, Sweden

Polishing: Using **Core Technology** for reduction of impurities.

AAV

	Capto Core 400	Capto Core 700
Matrix	Highly cross-linked agarose	
Average particle size (d_{50v})	90 μm	85 μm
Ligand	Octylamine	
Binding capacity*	22 mg ovalbumin/mL resin	13 mg ovalbumin/mL resin
Average molecular weight cutoff	M_r 400 000	M_r 700 000
Maximum flow velocity	700 cm/h in column with 20 cm bed height at < 2 bar (0.2 MPa)	500 cm/h in column with 20 cm bed height at < 2 bar (0.2 MPa)
pH stability Operational [†] CIP [‡]	3 to 13 3 to 14	
Chemical stability	All commonly used aqueous buffers, 1 M sodium hydroxide (NaOH) [§] , 6 M guanidine hydrochloride, 30% isopropanol, and 70% ethanol.	
Avoid	Oxidizing agents, citrate buffers	
Storage	20% ethanol at 4°C to 30°C	

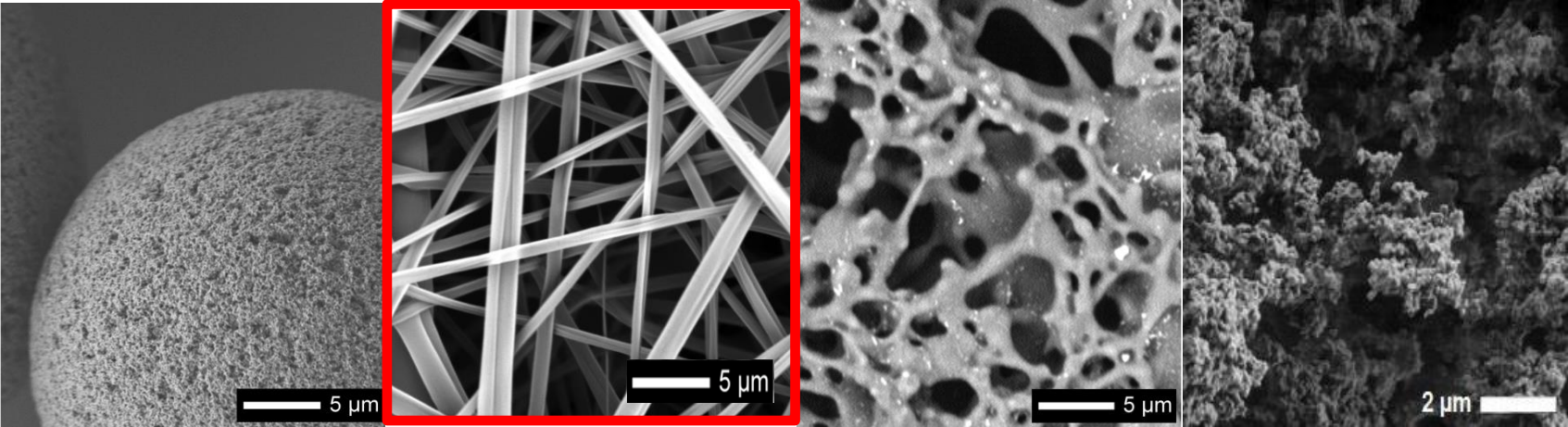


4

Technology formats

Evolution in formats :

Fibro – well-suited for viral vector chromatography



Bead resin

Fibro

Membrane adsorber

Monolith

Parameter	Bead resin	Fibro	Membrane	Monolith
Pore size	15–40 nm	0.2–2.0 μm	3–5 μm	2 μm
Surface area	~ 40 m ² /g	~ 10 m ² /g	< 2 m ² /g	< 7 m ² /g

Fibro has high surface area for high binding capacity and macroporosity needed for viruses

Fibro prototype for AAV5 purification

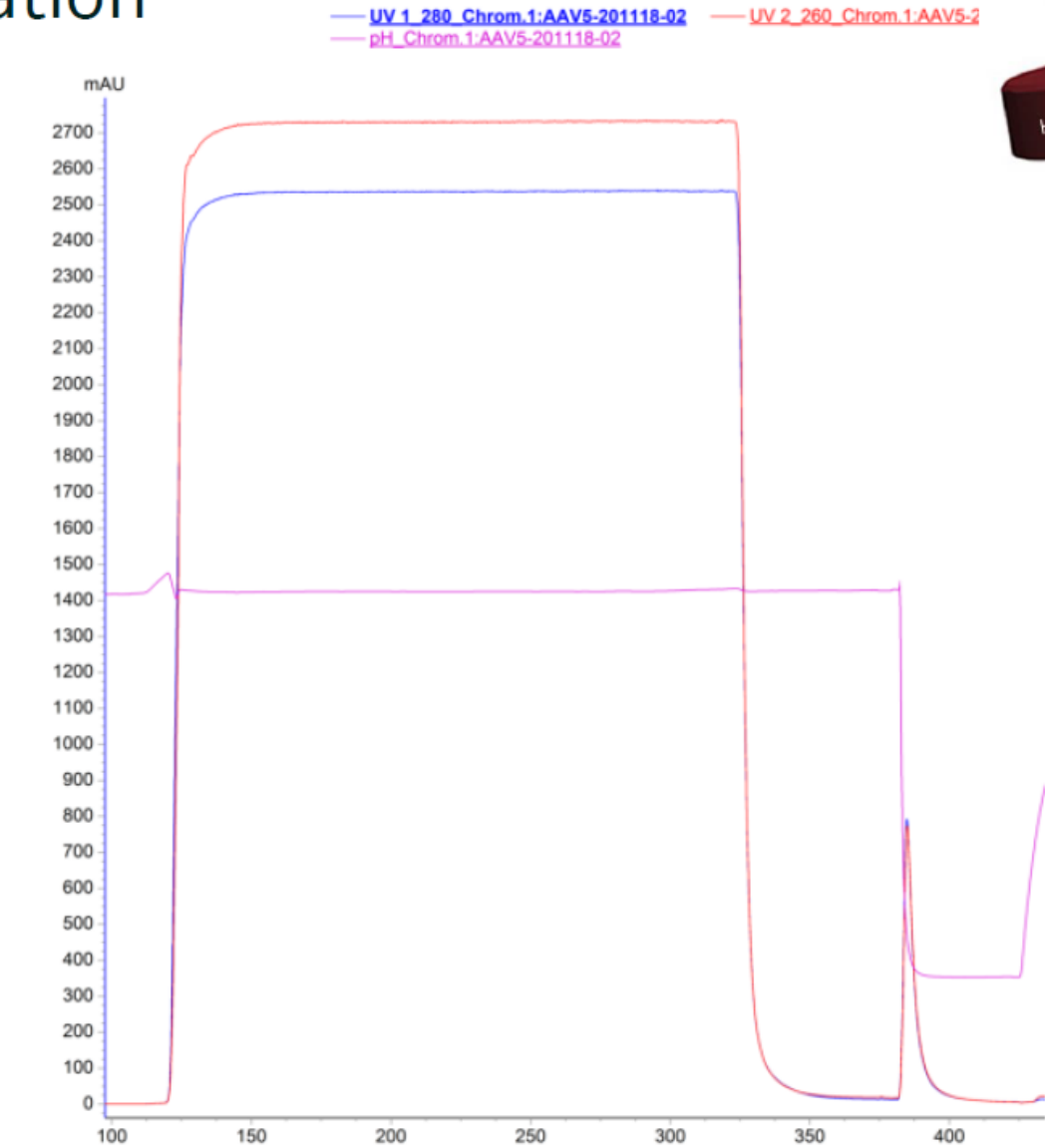
Biomanufacturer case study: clarified sample

- Residence time in seconds rather than minutes
- Loading time reduced from hours to minutes
- Recovery > 80% (based on vg)
- Purification performance similar to affinity resins
- Simple elution

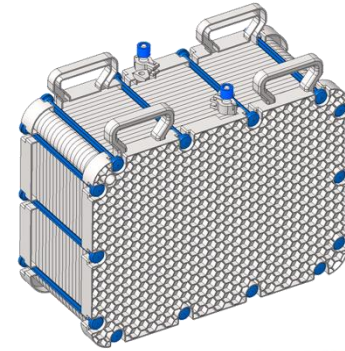
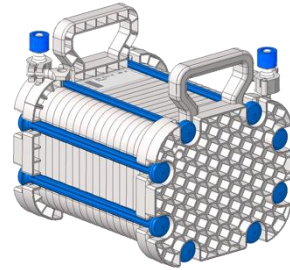
Clarification:	NFF + uncharged DF
Load:	9×10^{13} vg clarified AAV5
Load volume:	500 Fibro volumes ($500 \times 400 \mu\text{L} = 200 \text{ mL}$)
Load rate:	25 Fibro volumes/min
Residence time:	2.4 s
Load time:	20 min
Elution:	pH 2.5

Cytiva

Data from collaboration with [Belief BioMed](#)



Fibro formats and ÄKTA™ systems



Lab-scale Fibro unit

- Screening and HTPD tool
- 0.4 mL adsorbent volume
- ~ 1 L AAV feed

Small Fibro unit

- PD tool – scale-down mimic
- 4 mL adsorbent volume
- ~ 10 L AAV feed

Medium Fibro unit

- GMP compatible
- Up to 160 mL adsorbent volume
- Up to 500 L AAV feed

Large Fibro process unit

- GMP compatible
- Up to 2400 mL adsorbent volume

6

Conclusions

Conclusions

- ✓ Scalable suspension cell culture process with chemically defined medium
- ✓ High AAV titers with optimized transfection protocol ($\sim 10^{11}$ capsids/mL)
- ✓ TFF with 300 NMWCO hollow fiber
- ✓ Efficient affinity capture chromatography with Capto™ AVB
- ✓ Fibro technology improves AAV capture

Thank you



www.cytivalifesciences.co.jp

Cytiva (サイティバ)

法人名： グローバルライフサイエンステクノロジーズジャパン株式会社

バイオダイレクトライン

TEL: 03-5331-9336 / FAX: 03-5331-9370

e-mail: Tech-JP@cytiva.com

本資料の使用については、お客様施設内での使用に限ります。他社への転送、譲渡等は禁じます。本資料の著作権その他の知的財産権は、グローバルライフサイエンステクノロジーズジャパン株式会社に帰属します。無断転載、無断コピー、改ざん、二次利用を禁じます。掲載されている内容は2021年5月現在のもので、掲載されている製品は試験研究用以外には使用しないでください。掲載されている内容は予告なく変更される場合がありますのであらかじめご了承ください。掲載されている社名や製品名は、各社の商標または登録商標です。お問合せに際してお客さまよりいただいた情報は、お客さまへの回答、弊社サービスの向上、弊社からのご連絡のために利用させていただく場合があります。弊社は、資料の掲載内容の正確性を記すべく、情報を随時更新しておりますが全ての情報が最新であることを保証するものではありません。したがって、当資料上の掲載内容に誤りがあった場合でも弊社は責任を負いかねます。



Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Xcellerex, Biacore, HyClone, ULTA, ÄKTA, HiTrap, Cy and Capto are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Denarase is a trademark of c-Lecta GmbH.

MiniTEM is a trademark of Vironova AB. PicoGreen is a trademark of Thermo Fisher Scientific. Tween is a trademark of Croda International PLC. All other third party trademarks are the property of their respective owner.

Cy5 is covered under US patent number 6,828,116 and equivalent patents and patent applications in other countries in the name of GE Healthcare UK Limited. The purchase of CyDye products includes a limited license to use the CyDye products for internal research and development but not for any commercial purposes. A license to use the Cy and CyDye trademarks for commercial purposes is subject to a separate license agreement with GE Healthcare. Commercial use shall include:

1. Sale, lease, license or other transfer of the material or any material derived or produced from it.
2. Sale, lease, license or other grant of rights to use this material or any material derived or produced from it.
3. Use of this material to perform services for a fee for third parties, including contract research and drug screening.

If you require a commercial license to use the Cy and CyDye trademarks, please contact LSlicensing@ge.com.

© 2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)