

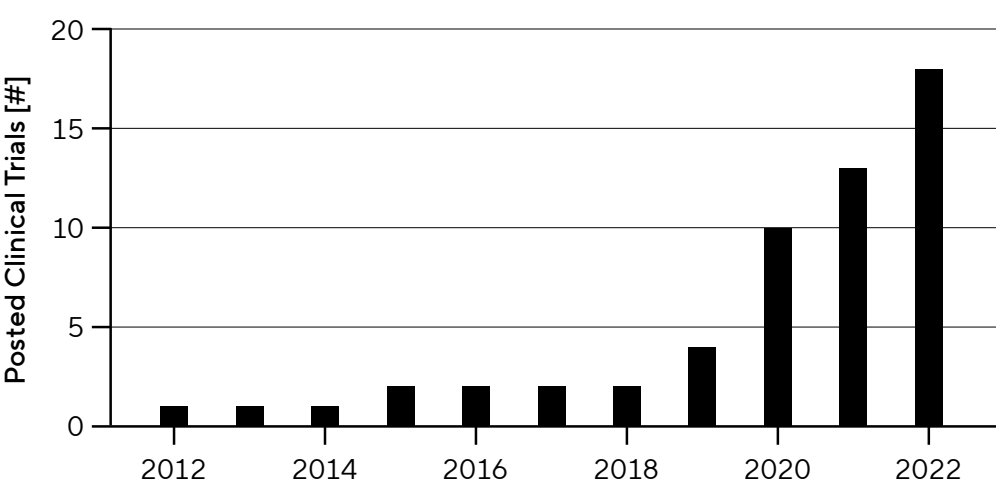
Successful Development of a Scalable and Robust Process for hMSC-EV Production

Grishma Patel, Sanket Jadhav, Rukmini Ladi, Priyanka Gupta, Stephen Lenzini, Jae Jung, Michael Boychyn, Jon Rowley, Elie Zakhem

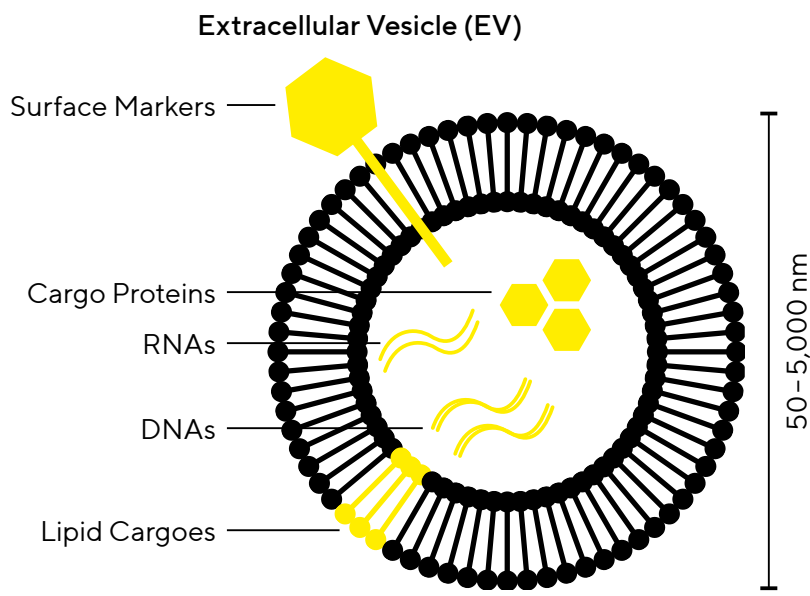
Background & Objective

- Extracellular vesicles (EVs) are nano-sized particles composed of a lipid bilayer.
- EVs are released from many different types of cells, and they contain proteins, nucleic acids (RNA and DNA), and lipid cargoes.
- EVs released from Mesenchymal Stromal | Stem Cells (hMSC-EVs) are rapidly gaining momentum in clinical trials.
- One of the challenges in EV production is scaling up the manufacturing processes using the optimal cell source and collection media.
- RoosterBio® offers reproducible and scalable processes to manufacture high-quality, xeno-free (XF) hMSCs and media for EV production.

Objective: The goal of this study was to develop a robust and scalable upstream process for hMSC-EV production in a stirred tank bioreactor system.

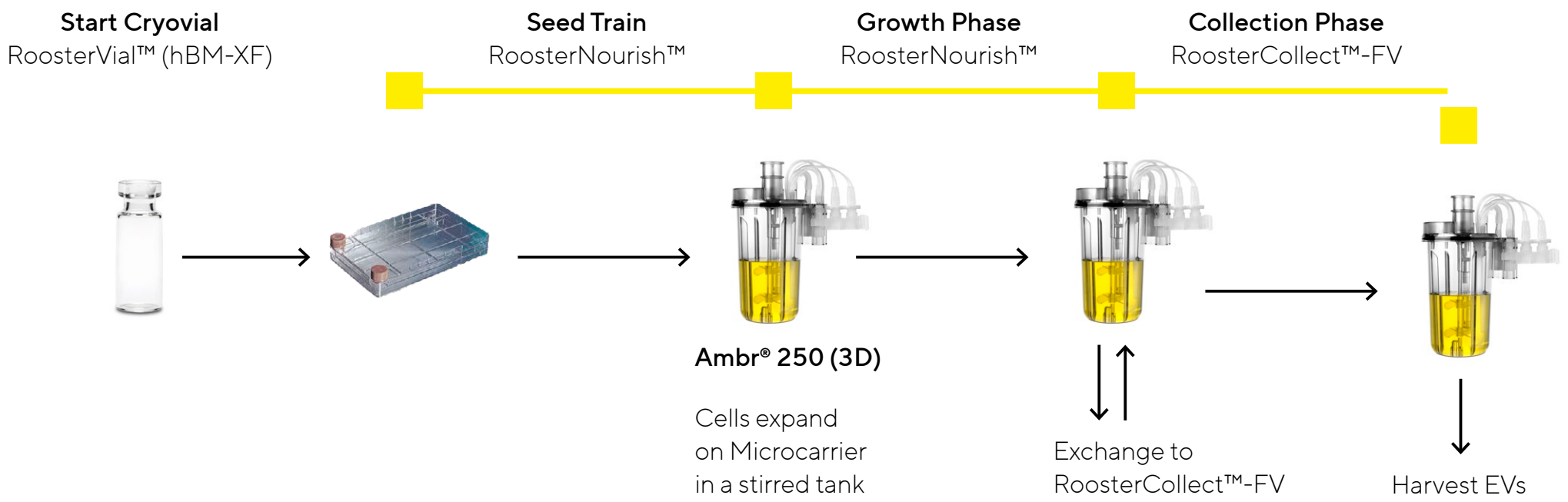


The majority of clinical trials involving the secretome and | or extracellular vesicles use MSCs as producers. (Source: Clinicaltrials.gov)



Process Development Approach

Process Flow Diagram Showing the Adapted Process for EV Manufacturing in Stirred Tank Bioreactors.

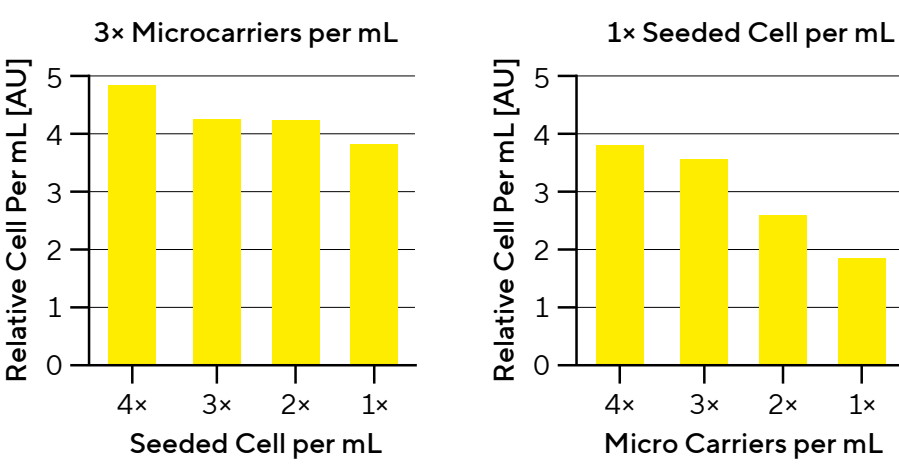
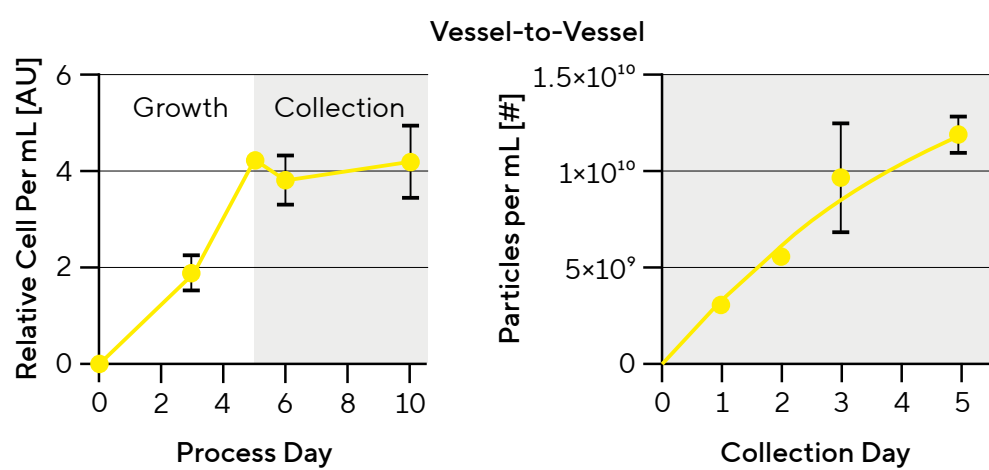
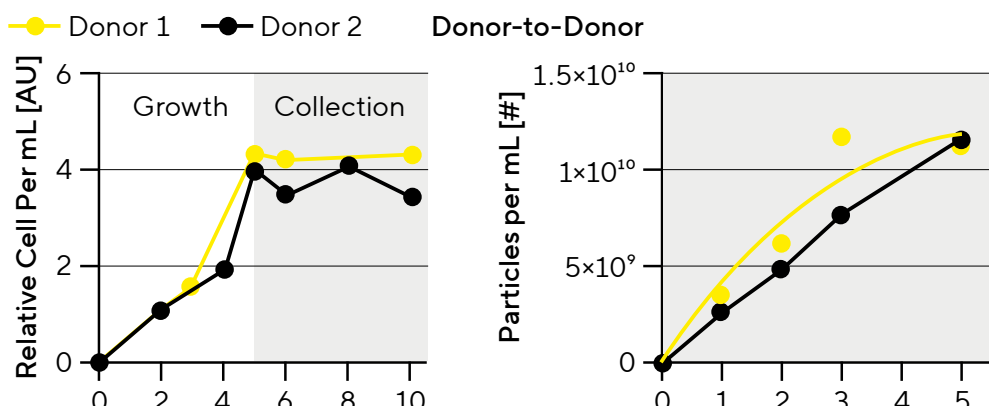
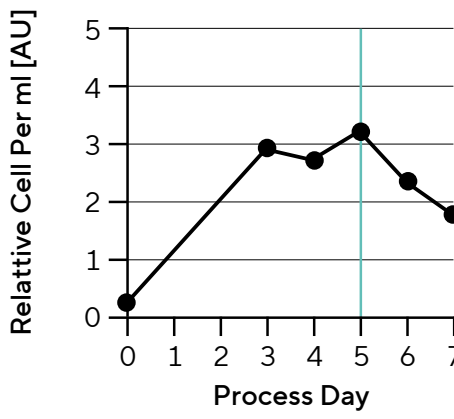
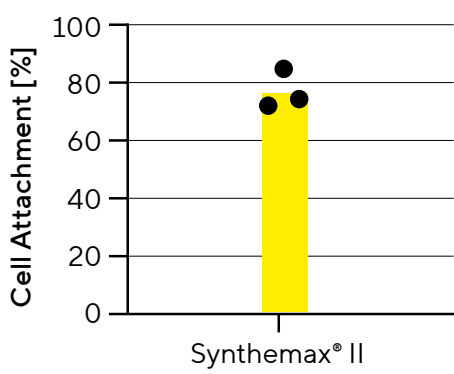


This study was carried out using xeno-free human bone marrow MSCs (RoosterVial™-hBM, RoosterBio®) expanded in RoosterNourish™ (RoosterBio®) in Ambr® 250 (Sartorius). EVs were collected in low-particle media, RoosterCollect™-EV (RoosterBio®).

- Critical process parameters (CPPs)** were evaluated and optimized to increase final cell yield and EV productivity:
 - Microcarrier type
 - Cell growth phase length (pre-collection)
 - Microcarrier density
 - Initial cell seeding density
 - EV collection time
- Critical quality attributes (CQAs)** were evaluated for particle count and size, identity (tetraspanins by western blot, percent lipidated particles, RNA content) and potency (in vitro wound healing assay).

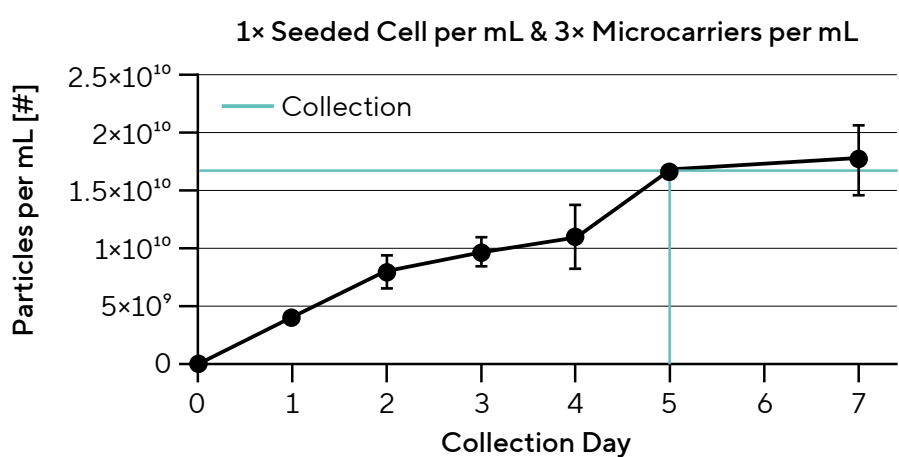
Small Scale (Ambr® 250) Process Development Results

(Right) Synthemax® II (Corning) showed high MSC attachment (~80%) and was selected as the optimal microcarrier. MSC growth reached its peak at day 5 of culture and then plateaus afterwards. Therefore, 5 days was selected as the optimal growth phase length.



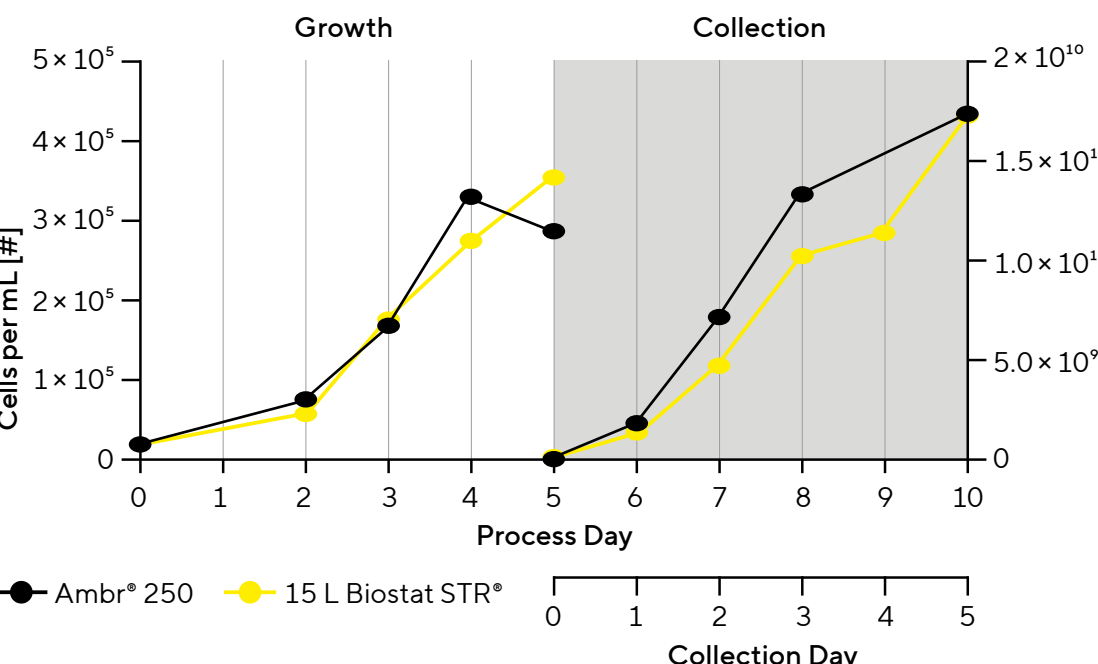
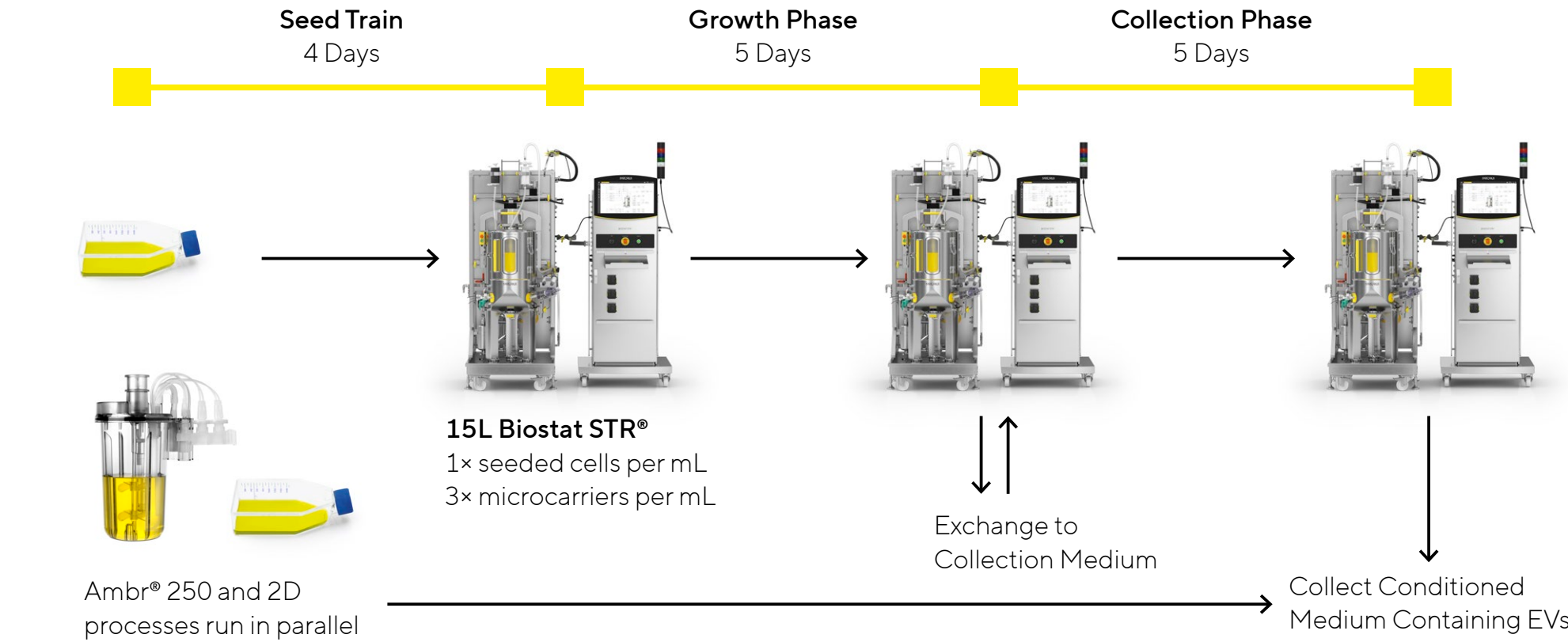
(Above) Increasing microcarrier density up to 3x significantly increases cell yield by day 5, and 1x cell seeding density was sufficient to reach maximal cell yield without cell | microcarrier aggregation.

(Below) After exchange into EV collection media, particle yield increased and peaked at day 5 of the collection phase.



Specifications subject to change without notice.
©2025 Copyright Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Göttingen, Germany
Status: 01 | 2025

15 L Scale-Up in Biostat STR®

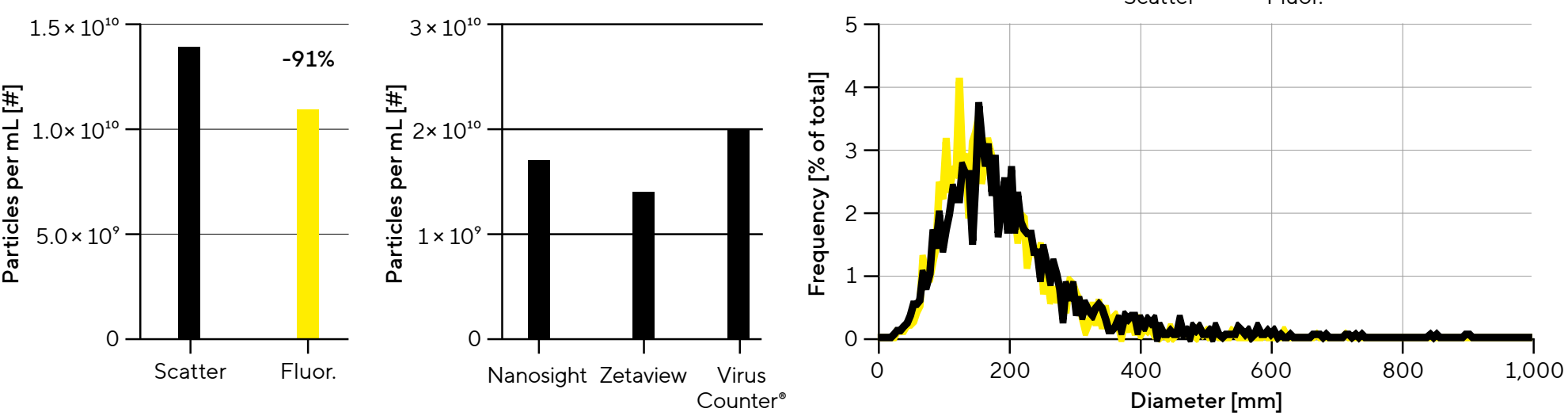


Critical Process Parameters identified at the small scale were verified at the large scale (15 L) in the Sartorius Biostat STR® bioreactor.

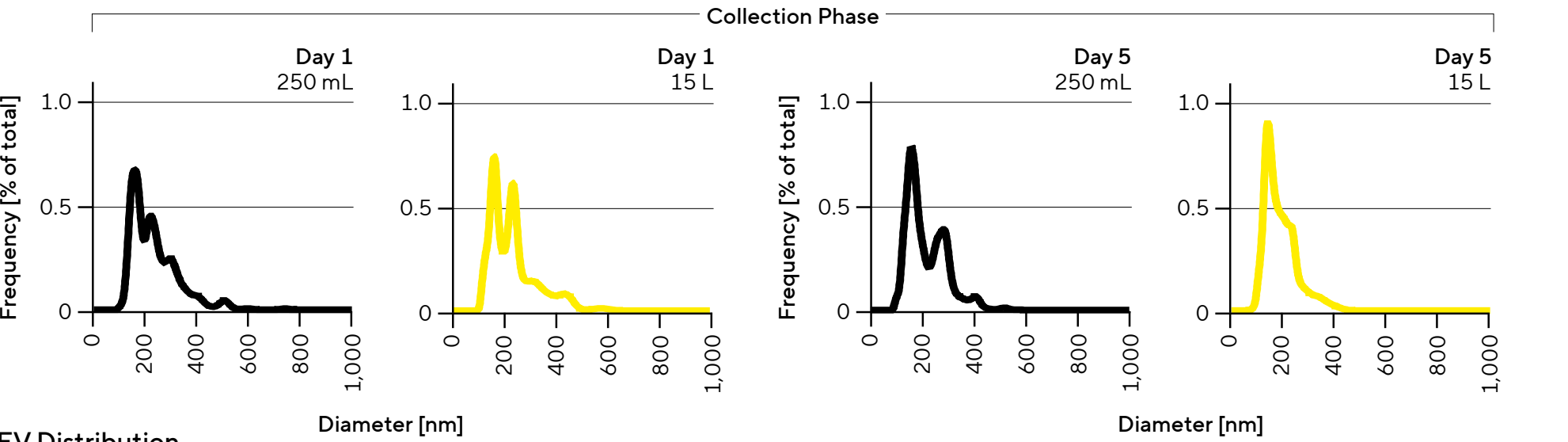
In Biostat STR®, MSCs reached up to ~400,000 cells per mL by the end of the growth phase (day 5), which was comparable to Ambr® 250. Particle production reached ~1.7 x 10¹⁰ per mL by day 5 of the EV collection phase, which was also consistent with the Ambr® 250 data. Ambr® 250 and 15 L Biostat STR® comparability demonstrate process scalability.

Critical Quality Attribute (CQA) Comparability

Nanoparticle Tracking Analysis (NTA)

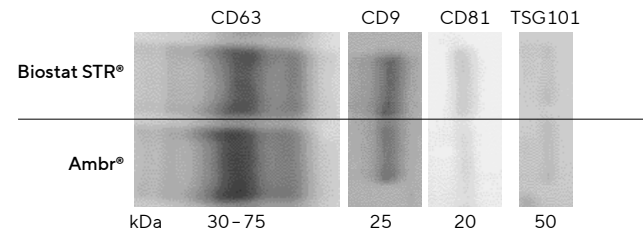


EV Count and Integrity: With NTA, the great majority (~91%) of particles stained positive for the lipid membrane dye Memglow, indicating 91% of the particles were EVs. Similar results are obtained across instruments. Virus Counter® (Sartorius) detects particles positive for lipid membrane dye CellMask Orange.

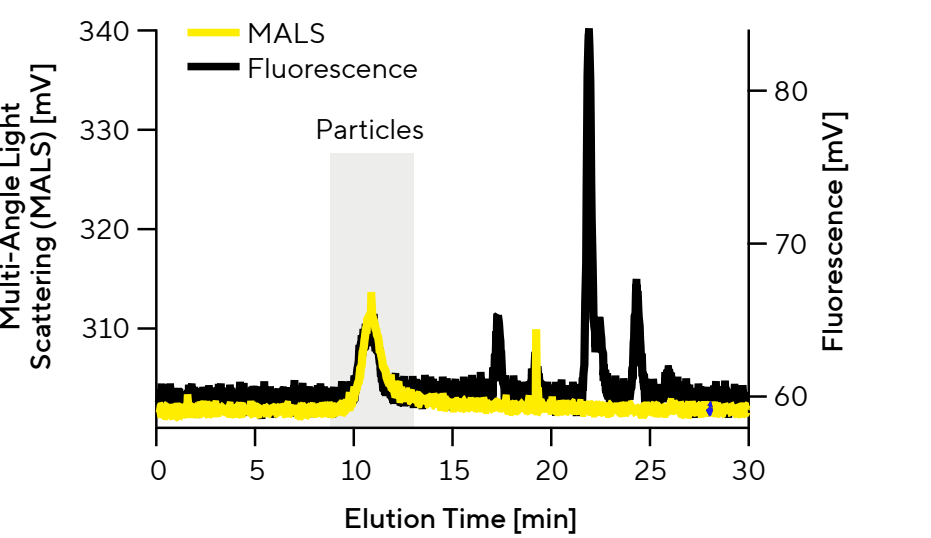


EV Distribution
Particle diameters ranged from 50 - 450 nm with median ~180 nm. Distributions are similar across platforms and collection phase duration.

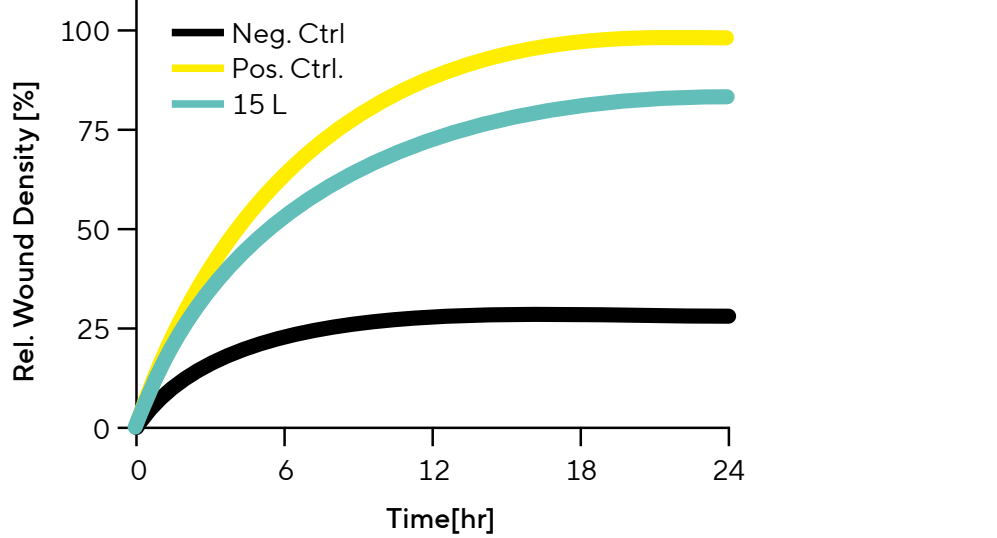
EV Identity
Harvested conditioned medium showed presence of conventional EV markers tetraspanins CD63, CD9, and CD81, as well as intraluminal marker TSG101.



PATfix® SEC-MALS Biochromatography and EV Potency Assay



PATfix® SEC-MALS: Earlier elution during size-exclusion chromatography is indicative of larger particles that did not enter the pores of the column and instead eluted in the void volume. Molecules that have entered the pores elute in order of decreasing size. Overlapping MALS and fluorescence signal (~10 - 12 min) indicates larger particles that are stained positive for lipid membrane dye CellMask Orange and are likely EVs.



EV Potency: EV conditioned medium from 15 L Biostat STR® harvest showed potency effect greater than unconditioned medium (Neg. Ctrl) and effect similar to growth medium (Pos. Ctrl) in an *in vitro* HUVEC cell wound healing assay. Cells are grown to a monolayer, a scratch wound is generated, and closure is measured over time using the IncuCyte® (Sartorius) platform. Data fit to a one-phase association model.

Conclusion and Next Steps

- We successfully developed a reproducible and scalable microcarrier-based stirred tank process for EV production using RoosterBio xeno-free hMSCs (RoosterVial™-hBM) coupled with RoosterBio® MSC expansion media (RoosterNourish™) and RoosterBio low particle EV collection media (RoosterCollect™-EV).
- The process was developed at the 250 mL scale (Ambr® 250) and its scalability was successfully demonstrated at the 15 L scale in Biostat STR®.
- The developed process consistently generates a conditioned medium containing ~1 x 10¹⁰ - 1.5 x 10¹⁰ particles per mL.
- Particles generated from this process were in the range of 50 - 450 nm diameter, stained positive with lipid membrane dyes across three independent analytical methods (NTA, Virus Counter, PATfix®), confirming that they are extracellular vesicles, expressed the EV specific tetraspanin markers (CD81, CD63 and CD9) and demonstrated biofunctionality through an *in vitro* wound healing assay.

Overall, the study demonstrated the successful development of a scalable and robust upstream process for hMSC-EV production.