

Simplify Upstream Process Intensification From PD to Manufacturing

Prity Bengani-Lutz, Priyanka Gupta

Sartorius Stedim North America Inc., 565 Johnson Avenue, Bohemia, NY 11716

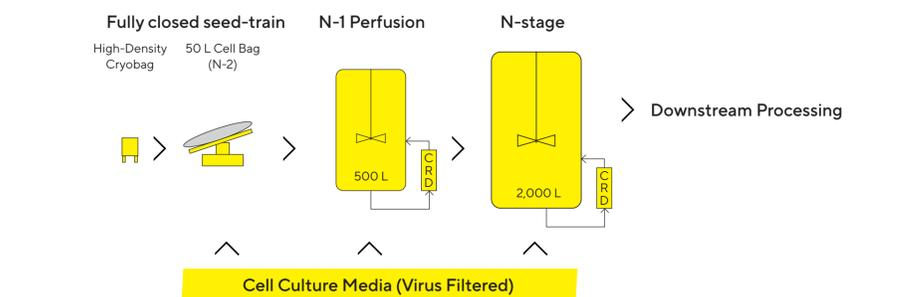
1. Process Intensification Satisfies Key Business Drivers

Low Capex Investment	< 50%	Process intensification is a holistic framework to maximize overall productivity of the unit operation(s), the manufacturing process and/or the facility output for biomanufacturing.
Faster Buildout Time	< 2 years	
Flexible Smaller Footprint	50 - 70%	Benefits of Process Intensification: <ul style="list-style-type: none"> Enables faster drug development Increases the efficiency and productivity in GMP manufacturing Applicable to any process irrespective of molecule Can be step-wise (per unit operation) or end-to-end for maximum impact Other terms used in the same context: "continuous", "connected" and "integrated"
Higher Productivity	2 - 3x	
Lower COGS	> 30%	

Source: Data derived from Biosolve modelling

2. Upstream Process Intensification Strategies

Requirements include robust cell line and media suitable for long culture durations & high VCDs of perfusion, scalable performance from clone screening to manufacturing along with recipes & PAT sensors for process control

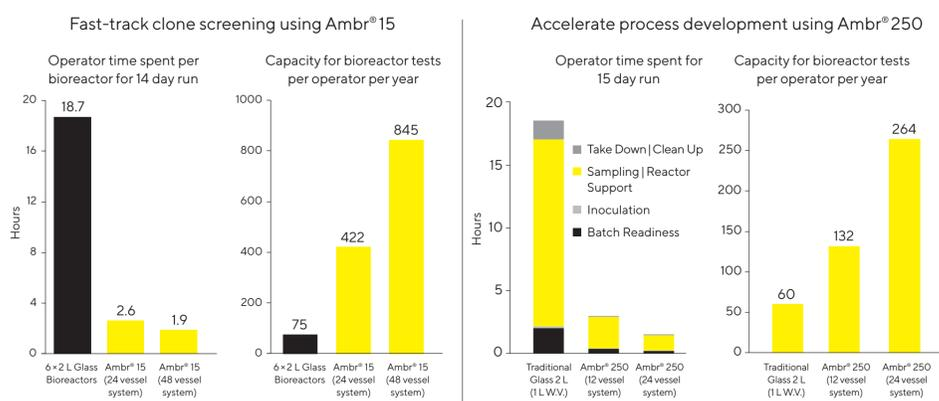


Seed Train Intensification	Main Bioreactor Intensification	Clarification Intensification
<ul style="list-style-type: none"> HCD cell banks direct thaw in Rocking Motion (RM) RM Perfusion (before N-3) to skip N-2 (Fed-Batch) N-1 Perfusion using RM or stirred tank reactor (STR) 	<ul style="list-style-type: none"> High inoculation FB from N-1 Perfusion Intensified Fed-Batch: <ol style="list-style-type: none"> Concentrated FB = Product in bioreactor (Dynamic) Perfusion = Product in permeate 	<ol style="list-style-type: none"> Single use centrifuge (e.g., Sartorius Ksep®) Multi column chromatography (e.g., Sartorius BioSMB)

Ref: BPOG Technology Roadmap 1st edition, 2017; 3. Process Technologies Bio® BioPhorum Operations Group, Ltd.

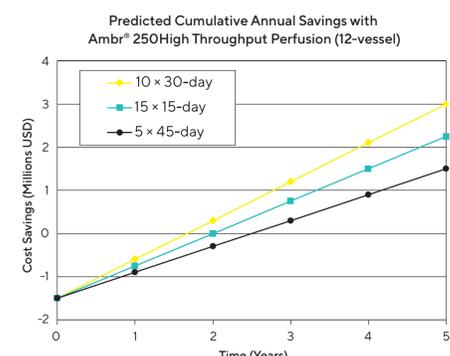
3. Clone Selection to Process Development

Save Time, Run More Experiments and Obtain Useful Data Using High Throughput, Multi-Parallel Ambr® Systems



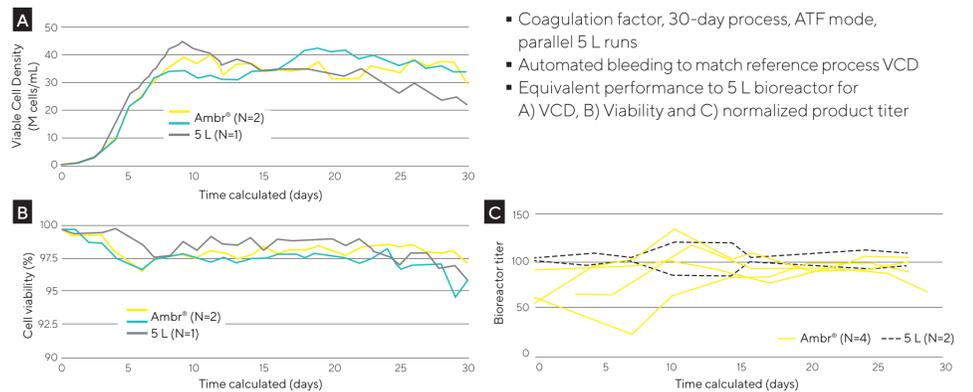
4. Process Development & Characterization

Ambr® 250 Perfusion is Efficient & Cost Effective Compared to Traditional Benchtop Perfusion



- System payback period 1-2 years
- The major source of cost savings is media cost reduction with Ambr® 250 High Throughput Perfusion
- Annual costs were calculated for three experiment capacity scenarios, for both:
 - 1x (Ambr® 250 High Throughput Perfusion - 12 vessels)
 - 2x (6x 2 L benchtop bioreactors)
- Annual cost savings calculated as:
 - cost (12x Ambr® 250) - cost (12x 2 L)
- Cumulative annual cost savings include initial CAPEX investment

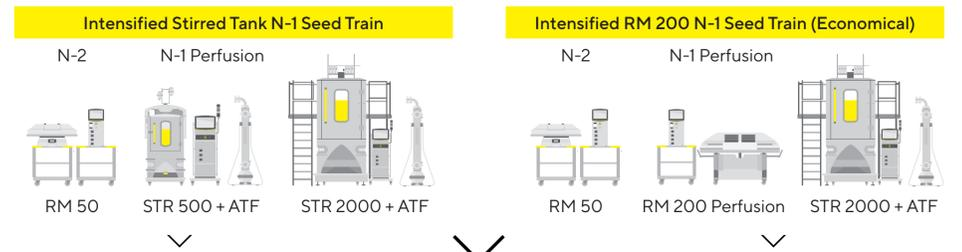
5. Ambr® 250 Perfusion: Good Match to 5 L for VCD & Titer



- Coagulation factor, 30-day process, ATF mode, parallel 5 L runs
- Automated bleeding to match reference process VCD
- Equivalent performance to 5 L bioreactor for A) VCD, B) Viability and C) normalized product titer

6. Flexible Seed Train Options

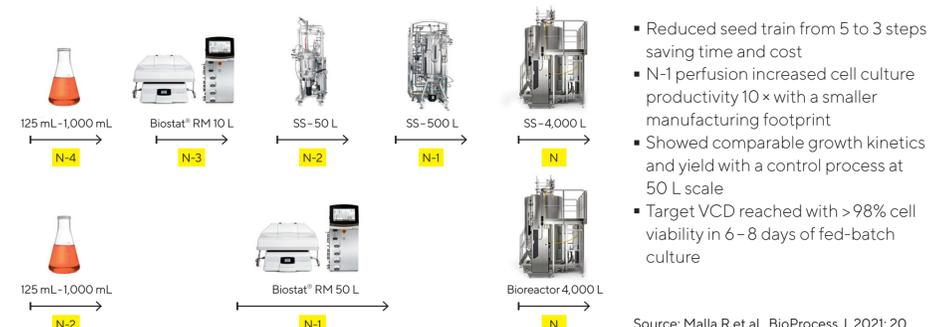
- Intensifying N-1 Seed Train is the easiest to implement with minimum change to the Fed-Batch process
- It results in upto 2x higher titer, upto 50% increase in throughput & lower COGS compared to Fed-Batch



- N-1 perfusion enables high inoculation of production bioreactor
- Reduce culture length of production bioreactor (~40%) - 17 seed train days (25-29 total USP days)
- 3 Bioreactors & 3 steps (instead of 4 each in FB)**
- 3-6 more batches/year/train**
- Upto 25% reduction in investment cost
- 50% lower consumables cost per batch
- Upto 20% reduction in production bioreactor culture length: 14 seed train days (22-26 total USP days)

7. Intensified N-1 Seed Train: 10x Productivity Increase

Traditional Batch (Top) Versus Perfusion (Below) Seed Train Cell Culture Process Workflow at Intas

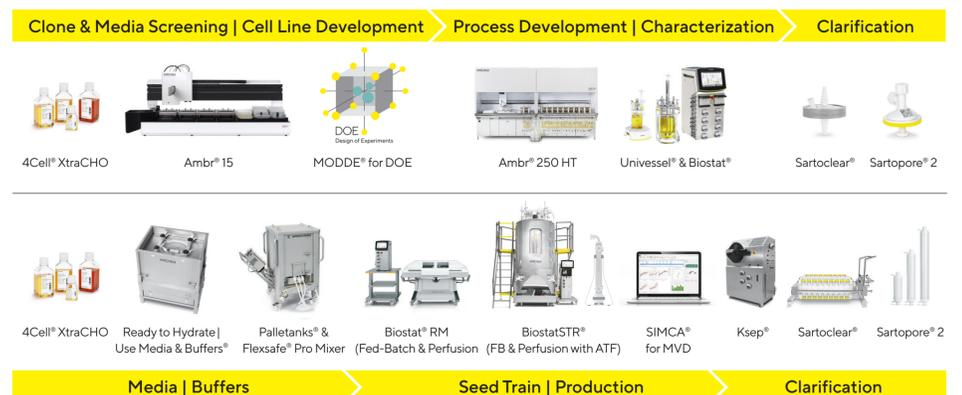


- Reduced seed train from 5 to 3 steps saving time and cost
- N-1 perfusion increased cell culture productivity 10x with a smaller manufacturing footprint
- Showed comparable growth kinetics and yield with a control process at 50 L scale
- Target VCD reached with >98% cell viability in 6-8 days of fed-batch culture

Source: Malla R et al., BioProcess J, 2021; 20.

8. Process Scale-up & Summary

Sartorius Upstream Process Intensification Platform: PD to Manufacturing



Summary

- Process Intensification maximizes productivity and flexibility and can be implemented step-wise or end-to-end
- High throughput Ambr® systems fast-track biopharm perfusion process development saving time and money
- N-1 Perfusion results in upto 2x high titer, lower COGS and is the easiest to implement in an existing facility
- PAT sensors enable culture monitoring, automated inoculation of subsequent cultures & better process control