

End-user evaluation of 30 L and 300 L HyPerforma Single-Use Fermentors and scale-down model

Summary

In moving processes from bench to pilot and ultimately production fermentors, it is critical to have a reliable scaledown model. The purpose of a scale-down model is to demonstrate at the benchtop level what can be achieved in production-scale fermentors. Scale-down models can vary depending on the media, strain, and process conditions. In this case, a customer wanted a scale-down model of the Thermo Scientific™ HyPerforma™ Single-Use Fermentor (S.U.F.) in order to establish what fermentation processes could be scaled into the S.U.F., in particular for the production of plasmid DNA. The 300 L S.U.F. exceeded the customer's expectations. The customer anticipated that the test process using *E. coli* with high oxygen consumption would reach a final OD of ~80 and the HyPerforma S.U.F. exceeded expectations by reaching an OD of 147.

This application also was a successful demonstration of media preparation through culture harvest, using a jacketed 50 L Thermo Scientific™ HyPerforma™ Single-Use Mixer (S.U.M.) (60 L working volume and heated to 67°C), 200 L Thermo Scientific™ HyPerforma™ DS 300 Single-Use Mixer (S.U.M.), Thermo Scientific™ Powdertainer™ BioProcess Container (BPC) system, 2 L–50 L Thermo Scientific™ 2D Labtainer™ BPC systems, pre-irradiated filters, sterile funnel BPC, seed BPC, 30 L and 300 L HyPerforma S.U.F.s, and Thermo Scientific™ CentriPAK™ harvest BPCs.

After evaluating additional leading single-use fermentors from two other suppliers, the customer purchased the Thermo Scientific 300 L HyPerforma S.U.F..

Purpose

To evaluate the 300 L HyPerforma S.U.F. for GMP plasmid production process using a customer's standard evaluation procedure and test strain.

Procedure

The *E. coli* culture was maintained at 37°C ± 0.1, pH 7 ± 0.05, maximum pressure of 0.55 psi, and a minimum dissolved oxygen (DO) of 30%. DO control cascaded at 200–375 rpm, 200–500 slpm air, and 0–400 slpm of oxygen as needed to maintain the minimum set point of 30% DO.

After scaling up the model procedure to the 300 L S.U.F., the power-to-volume input, $k_L a$, and gassing rates were scaled down into the 1 L glass fermentors. The same original procedure was then followed but using the scaled-down power-to-volume input, $k_L a$, and gassing rates for S.U.F. comparison. The scale-down model could then be used for estimating performance with all production strains for this customer.

The procedure was then repeated with the 30 L S.U.F. and showed scalability within the HyPerforma product line.

E. coli cells were collected with the Sorvall BIOS 16 centrifuge using CentriPAK BPC singles with quick connect at 5,373 x g for 15 minutes in 15 L batches. The 30 L S.U.F. was harvested as closed system in two batches using CentriPAK manifolds. The 300 L S.U.F. was harvested within 3 hours using two BIOS 16 centrifuges and CentriPAK singles filled to about 1.87 L each.

During this feasibility study, the off-gas analyses of the cultures were compared. It was noticed that the off-gas trend was comparable. The off-gas analyzer sample line to each S.U.F. can be placed inside the exhaust filter exit cavity or connected to a small sterile filter connected to the S.U.F.'s filter chamber.

Equipment and supplies

Strain:

- HMS174(DE3)

Chemicals:

- Base
- Acid
- Antifoam C8840 (New London Chemicals)
- 60 L nitrogen source, 30% w/v feed solution
- 50 L carbon source, 50% w/v feed solution
- Batch medium

Bioreactor setup

The S.U.F. systems were set up according to the user guide and controlled by Thermo Scientific™ TruBio™ Software, powered by the DeltaV™ Distributed Control Platform from Emerson, utilizing the Thermo Scientific™ TruFluor DO single-use sensors and Hamilton or Mettler Toledo single-use pH sensors. Operating parameters are listed in Table 2.

Bioreactor inoculum, cultivation, and scale-up

The inoculum was cultured in an incubator at 37°C, 250 rpm, 1 inch arc, and for 16 hr.

Table 1. Equipment and materials.

Description	Cat. No.
30 L S.U.F. Hardware	S.U.F.0030.AAA.BAAABB0C00
30 L S.U.F. BPC	SH3B11722.01
300 L S.U.F. Hardware	S.U.F.0300.AAA.DAAABB0C00
300 L S.U.F. BPC	SH3B11861.01
100 L S.U.M., jacketed with touchscreen console	SUM0100.9002
200 L plastic drum	SH30959.03
Drum dolly	SH30958.01
Nalgene™ polyethylene 5 gal tank liner	11100-0005
Nalgene cylindrical 5 gal tank	43050-0005
1.5 L funnel	SH3B14865.01
PowderFill or funnel stand	129752
HQ incubator shaker	11-676-235
Seed BPC (3 L working volume)	SH3B9830.01
Seed BPC clip	122554
0.5, 1, 2, 5, 10, and 50 L Labtainer BioProcess Container (BPCs)	SH30712.01-.02 and SH30963.01-.03
Powdertainer BPCs	SH30737.01 and SH30737.02
HyPerforma™ G3 Controllers	NA
Prima™ BT off-gas analyzer	NC1256292
Bios 16 Sorvall Bioprocessing Centrifuge	2 x L85007685
CentriPAK BPC Adapter 2 L bucket liner	4 x 75003873
CentriPAK BPC 6 x 1.7 L harvest manifold	2 x 75003880
CentriPAK BPC Single with Quick connect	14 x 75003891

Table 2. 30 L and 300 L S.U.F. operating conditions.

Parameter	30 L	300 L
Initial volume	24 L	240 L
Final volume	~30 L	~300 L
Temperature	37°C	37°C
pH	7.0 ± 0.05	7.0 ± 0.05
Agitation	300–600 RPM	200–375 RPM
DO setpoint	30%	30%
DO cascade	Cascade RPM, air, then supplement oxygen	Cascade RPM, air, then supplement oxygen
Gas flow	3–60 standard liter per min (slpm)	30–500 slpm
Antifoam	3 mL, more as foam detected	30 mL initially, more as foam detected

Results

The standard procedure for expression in *E. coli* model was used for technical transfer to the S.U.F. The culture conditions in the 300 L S.U.F. are seen in Figure 1, showing feasibility of plasmid production in the 300 L S.U.F. as per the customer’s standard procedure. The customer stopped the feed in 300 L S.U.F. to be able to harvest right away, where in the 30 L the feed ran out hence the 30 L reached a higher density.

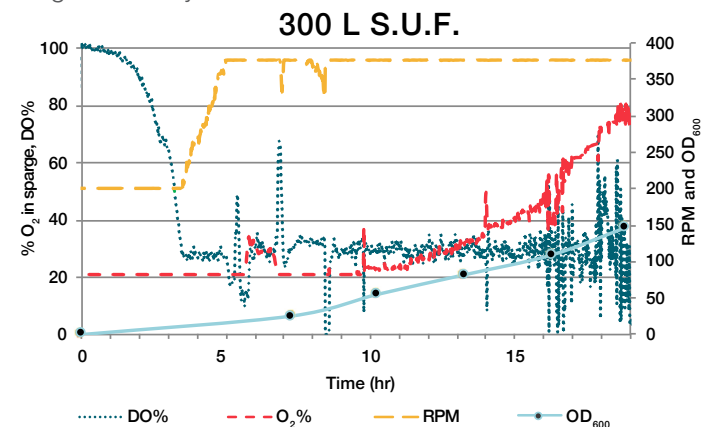


Figure 1. Data from 300 L S.U.F. (n = 1) after one trial scale-up following customer’s procedure. The process in the HyPerforma S.U.F. reached an OD of 147 by 19 hr. Samples showed a wet cell weight (WCW) of 167 g/L was achieved.

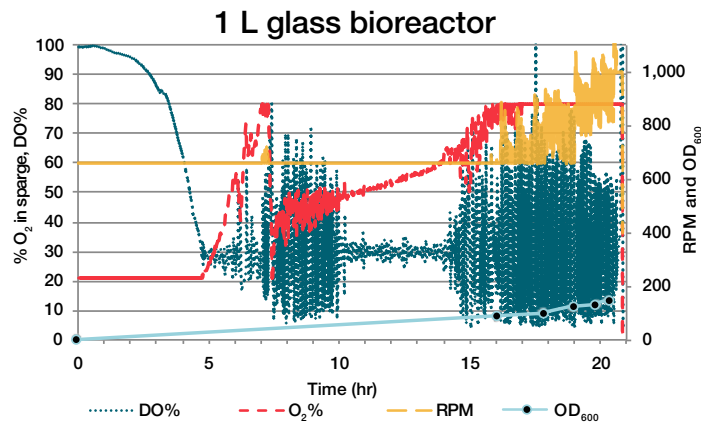


Figure 2. Data from 1 L glass vessel (n = 1) used as a working-volume model for the scale-up procedure. The WCW was 170 g/L for the final samples, same as that achieved in the 300 L S.U.F.

The power-to-volume inputs was calculated for reproducing the conditions from a 300 L S.U.F. to the customer’s 1 L benchtop glass fermentors. The same original procedure was then followed, but using the scaled-down power-to-volume input, $k_L a$, and gassing rates. Culture conditions of one of the two scale-down 1 L fermentors are shown in Figure 2. The procedure was then scaled to the 30 L S.U.F. (Figure 3).

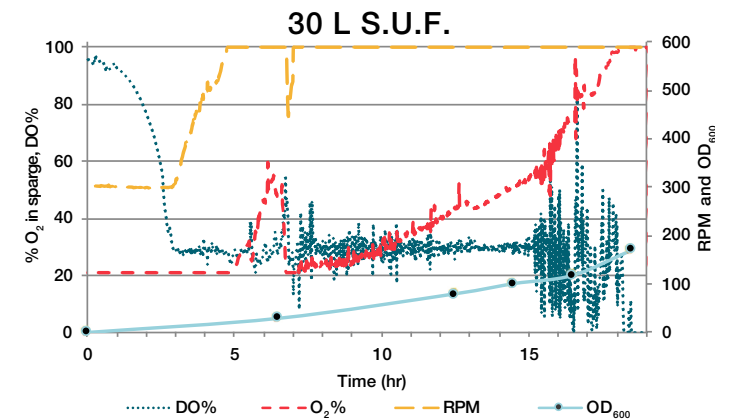


Figure 3. Data from 30 L S.U.F. (n = 1)—second scale-up following customer’s procedure. In the 30 L S.U.F., an OD of 170 was reached in our offline measurements. The WCW was 221 g/L.

Conclusion

The HyPerforma 300 L S.U.F. exceeded the customer’s expectations and reached a WCW of 167 g/L before feed was stopped. It was determined that the 300 L and 30 L S.U.F.s are adequate for production following the procedures for most of the customer’s strains.

Authors

1. Jason Brown, Process Development Engineer, Research and Development, Thermo Fisher Scientific
2. Nephi Jones, Senior Manager, Research and Development, Thermo Fisher Scientific

Find out more at thermofisher.com/suf