

November, 2017

Keywords or phrases:

Univessel® SU, single-use bioreactors, scalability, mesenchymal stem cells, microcarrier, advanced therapies

Microcarrier-Based Expansion of Human Mesenchymal Stem Cells in the Univessel® SU

Anita Ott¹, Valentin Jossen¹, Ann Siehoff², Christian van den Bos², Dieter Eibl¹, Regine Eibl¹

¹ Zurich University of Applied Sciences, School of Life Sciences and Facility Management, Institute of Biotechnology, Wädenswil, Switzerland

² Lonza Cologne GmbH, Cologne, Germany

Correspondence

Email: bioprocess.support@sartorius.com

Abstract

This note describes the culture conditions and setpoints for expansion of human adipose-derived mesenchymal stem cells (hAD-MSC) on microcarriers in the Sartorius Biostat® 2 L single-use Univessel. Typically, such culture conditions can be scaled to 50 L and larger Biostat STR® reactors. Reliable, scalable expansion of MSCs is critical for their use as a therapeutic agent.

Introduction

The increasing number of clinical trials with human mesenchymal stem cells (MSCs) and progress in regenerative medicine^{1,2} has created a need for higher quantities of these cells at the desired quality. For autologous therapies, a minimum of 5×10^8 cells are required for one single dose². It has already been demonstrated that MSCs can successfully be grown on microcarriers in single-use stirred bioreactors at benchtop-scale instead of the commonly used planar, one- or multiple-layer flasks, such as CellSTACKs or Cell Factories³⁻⁶.



Medium, Methods and Equipment

Overview of Set-up Procedure

Day -1

Preparation of SoloHill Pronectin F (0.75% solid fraction) as well according to the manufacturer's recommendation.

Day 0

Installation of the Univessel® SU bioreactor. Transfer of the microcarriers into the bioreactor vessel and filling with 0.7 L of custom-made MSC medium. Equilibration of the optical sensor patches for pH and DO measurement for 4 hours.

Thawing and inoculation with a seeding density of 1.5×10^7 MSCs (pooled cells from cryo-preserved vials). 4-hour cell attachment phase without agitation. Filling up to 2 L working volume with medium. Starting agitation at 100 rpm.

Day 0-7

Sampling, analytics and recalibration of online pH and DO sensors. Instead of the sampling device depicted in Fig. 1, single-use manifold bags were used (see also⁵).

Day 4

Exchange of 50% of medium with fresh medium.

Day 7

Cell harvest by separation of the microcarrier-cell aggregates from the medium, enzymatic cell detachment, cell resuspension, vialing and freezing.

Cells, Medium and Microcarrier

Cryopreserved adipose-derived MSCs (Lonza Cologne GmbH, Germany) from a single consenting and informed donor (second passage, PDL = 10) were used for inoculation, with a recommended seeding number of 1.5×10^7 cells. The cells were cultivated in a serum-reduced (5% FBS) medium. The required amount of SoloHill ProNectin F microcarriers (0.75% solid fraction) was prepared according to the manufacturer's specifications.

Cultivation Setup

The cultivation process of the MSCs in the Univessel® SU includes 3 steps: (1) initial cell attachment, (2) cell expansion, and (3) cell harvest. The microcarriers along with 0.7 L of medium were transferred to the Univessel® SU and equilibrated for 4 hours at 37 °C, 5% CO₂, 85 rpm. Afterwards, the thawed and pooled MSCs were inoculated at an initial density of 2.8×10^3 cells cm⁻². No agitation was performed for 4 h to allow the cells to attach to the microcarriers. The Univessel® SU was then filled with medium to the maximal working volume of 2 L. The impeller speed was initially set to 100 rpm and increased to 135 rpm as cultivation progressed. The MSCs were grown at 37 °C, pH 7.2 and 0.1 vvm headspace aeration for 7 days. On day 4 of cultivation (cell density between 2 and 3×10^4 cells cm⁻²), a 50% medium exchange was performed to prevent nutrient limitation. Once the maximum cell density was reached (day 7 of cultivation), the cells were separated from the microcarriers by using a sieving procedure combined with enzymatic cell detachment and washing steps⁵. Finally, the microcarrier-free suspension was centrifuged at 2000 g for 8 min, followed by supernatant removal and cell resuspension in fresh culture medium before vialing and freezing occurred.

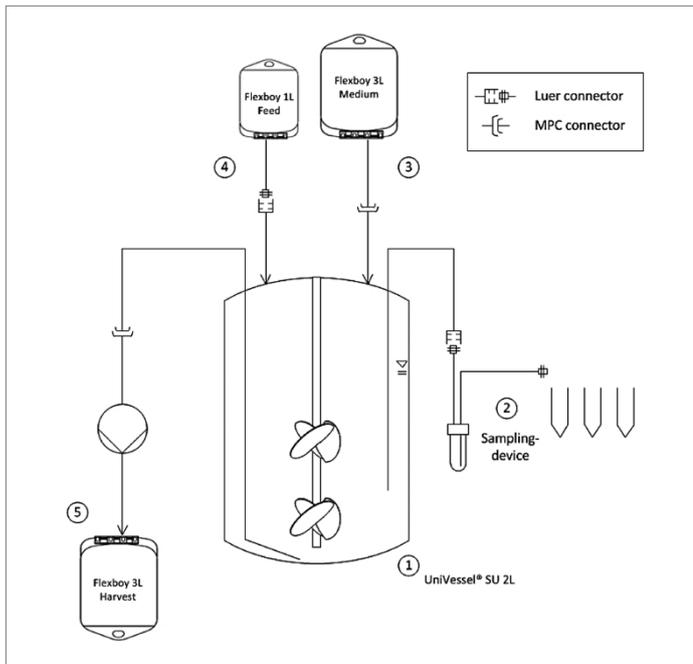


Figure 1: Experimental Setup of the Cultivation Process in the Univessel® SU 2 L.
Schematic overview of the experimental setup. The Univessel® SU 2 L (1) is additionally equipped with a sampling device (2), medium bag for medium feed after cell attachment (3), feeding bag for partial media exchange on day 4 (4), and harvest bag for cell harvest on day 7 (5).

Culture Conditions

Working volume	2 L
Microcarrier concentration	0.75% microcarrier solid fraction
Inoculated cell number	1.5×10^7 MSCs
Agitation speed	100 rpm (below NS1u) and 135 rpm (NS1)
Temperature	37 °C
Head aeration	0.1 vvm
Dissolved oxygen (DO)	> 20%
Cultivation time	7 days

Sampling and Quality Control

A daily sample of approximately 25 mL was taken with a sampling device to measure metabolites with the CedexBio (Roche Diagnostic) and/or the BioProfile (Nova Biomedical) as well as cell densities with the NucleoCounter® NC-100™ (Chemometec). Furthermore, a staining of the microcarrier-cell suspension with 4,6-diamidino-2-phenylindole (DAPI) was performed to evaluate carrier colonization and aggregate formation. In order to check the quality of the MSCs after cell harvest, flow cytometric investigations with fluorochrome-conjugated anti-human CD34, CD45, CD73, CD90 and CD105 (eBiosciences) were carried out. The results were compared to the surface marker profiles of the inoculated cells.

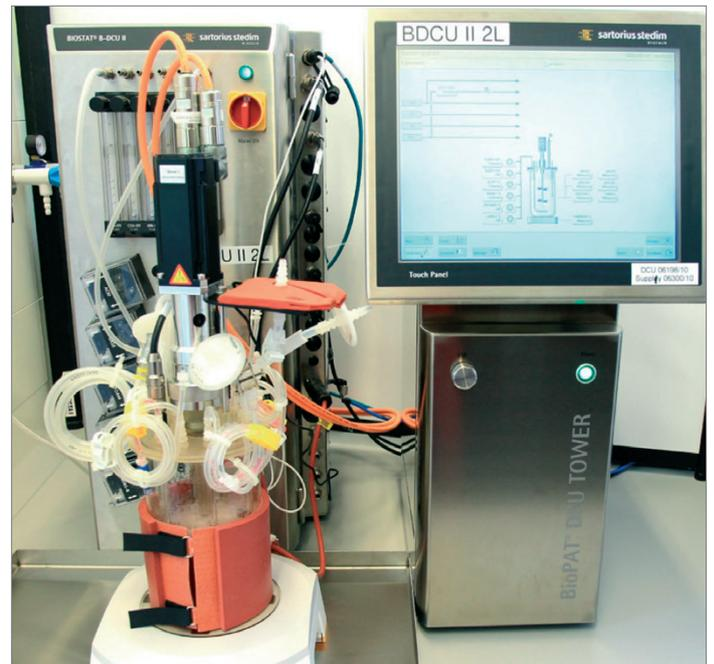


Figure 2: Experimental Setup at the Zurich University of Applied Sciences (ZHAW).

Results

Figure 4 exemplarily shows a typical time-dependent profile of the cell growth in the Univessel® SU bioreactor and control spinner flask. The comparability of the growth patterns in the two cultivation systems is evident. Within 7 days of expansion, a maximum cell number of $(5.3 \pm 0.5) \times 10^8$ MSCs was achieved in the Univessel® SU. This corresponds to an expansion factor of 35.4 ± 0.4 and a population doubling time of (25.6 ± 0.2) hr. On day 3, the cells were in a lag phase followed by an exponential growth phase until day 7. Glucose and lactate concentrations (data not shown) in the medium correlated well with cell growth. On day 4, a 50% media exchange was performed. DAPI staining of microcarrier-cell aggregates was performed on day 7, immediately before cell harvest. Notably, the microcarriers were completely covered with cells and formed aggregates of 2 mm in diameter. TrypLE Select-based detachment of the cells led to a harvest recovery of > 98% with cell viabilities exceeding 98%.

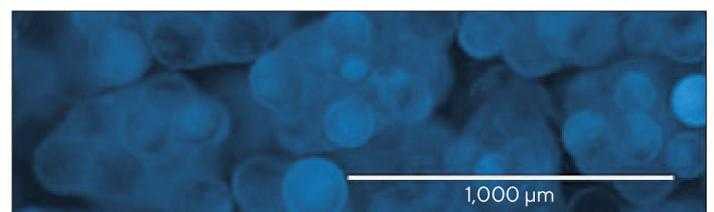


Figure 3: Polystyrene-Based Microcarriers Completely Covered With MSCs (White Bar = 1000 µm).

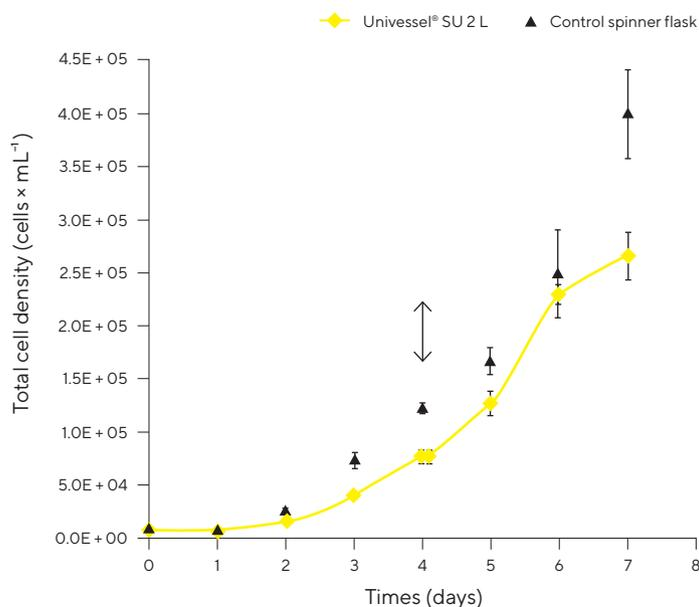


Figure 4: Growth Courses of MSCs Expanded in the UnivesseL® SU Bioreactor and Control Spinner Flask and DAPI-Staining on Day 7 of the MSC-Cultivation. The arrow indicates the 50% medium exchange.

Cell Quality Control

Flow cytometric analysis of the MSCs (presence of CD73⁺, CD90⁺, CD105⁺ and absence of CD34⁻, CD45⁻ surface markers) demonstrated that the phenotypic properties were maintained (data not shown).

Conclusion

The results demonstrate that the UnivesseL® SU bioreactor is suitable for the expansion of adipose-derived MSCs growing on microcarriers whilst maintaining the expected surface marker expression profile. Moreover, the UnivesseL® SU bioreactor can be used to generate clinically relevant doses for autologous cell therapy, typically in the order of 5×10^8 for a single dose.

References

- Wang J., Liao L., Tan J. Mesenchymal-stem-cell-based experimental and clinical trials: Current status and open questions. *Expert Opin Biol Ther.* 2011;11(7):893-909.
- Ren G., Chen X., Dong F., Li W., Ren X., Zhang Y., Shi Y. Mesenchymal stem cells and translational medicine: Emerging issues. *Stem Cells Transl Med.* 2012;1(1):51-58.
- Cierpka K., Elseberg CL., Niss K., Kassem M., Salzig D., Czermak P. MSC production in disposable bioreactors with regards to GMP and PAT. *Chem Ing Tech.* 2013;85(1-2):67-75.
- Jing D., Sunil N., Punreddy S., Aysola M., Kehoe D., Murrel J., Rook M., Niss K. Growth kinetics of human mesenchymal stem cells in a 3-l single-use, stirred-tank bioreactor. *Biopharm Int.* 2013;26(4):28-38
- Schirmaier C., Jossen V., Kaiser SC., Jüngerkes F., Brill S., Safavi-Nab A., van den Bos C., Eibl D., Eibl R.. Scale-up of adipose tissue-derived mesenchymal stem cell production in stirred single-use bioreactors under low-serum conditions. *Eng Life Sci.* 2014;14(3):292-303.
- Jossen V., Kaiser SC., Schirmaier C., Herrmann J., Tappe A., Eibl D., Siehoff A., van den Bos C., Eibl R.. Modification and qualification of a stirred single-use bioreactor for the improved expansion of human mesenchymal stem cells at benchtop scale. *Pharm Bioprocess.* 2014;2(4):311-322.

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For more information, visit
www.sartorius.com