

March 22, 2021

MSC NutriStem® XF Medium Scientific References

- **Clinical Trials**
- **Extracellular Vesicles and Exosomes**
- **Publications**

Clinical Trials

1. Y. Matsuo, et al. **Isolation of adipose tissue-derived stem cells by direct membrane migration and expansion for clinical application.** *Human Cell* (2021). <https://doi.org/10.1007/s13577-021-00505-3>
2. T. Iwanaka, et al. **A model study for the manufacture and validation of clinical-grade deciduous dental pulp stem cells for chronic liver fibrosis treatment.** *Stem Cell Res Ther* 11, 134 (2020). <https://doi.org/10.1186/s13287-020-01630-w>
3. V.T. Hoang, et al. **Expansion of Human Mesenchymal Stromal/Stem Cells Using Standardized Xeno-Free, Serum-Free Culture Condition.** *blood*, Volume 134, Issue Supplement_1, November 13 2019, <https://doi.org/10.1182/blood-2019-132140>
4. C. Bozkurt, et al. **The Use of Allogeneic Mesenchymal Stem Cells in Childhood Steroid-Resistant Acute Graft-Versus-Host Disease: A Retrospective Study of a Single-Center Experience.** *tjh*, 2019, 2019.0090 ;36:186-192
5. D. Ben-David, et al. **Autologous cell-coated particles for the treatment of segmental bone defects-a new cell therapy approach.** *Journal of Orthopaedic Surgery and Research*, volume 14, Article number: 198 (2019)
6. L.A. McIntyre et. al. **Cellular Immunotherapy for Septic Shock. A Phase I Clinical Trial.** *American Journal of Respiratory and Critical Care Medicine*, Vol. 197, No. 3, 2018
7. K. Schlosser et al. **Effects of Mesenchymal Stem Cell Treatment on Systemic Cytokine Levels in a Phase 1 Dose Escalation Safety Trial of Septic Shock Patients.** *Critical Care Medicine*, doi: 10.1097/CCM.00000000000003657

Extracellular Vesicles and Exosomes

1. M. A. Skylar-Scott, et al. **Biomanufacturing of organ-specific tissues with high cellular density and embedded vascular channels.** *Sci. Adv.* 2019.
2. Y. Li, et al. **Mesenchymal Stem Cells and Acellular Products Attenuate Murine Induced Colitis.** *Stem Cell Research & Therapy*. 03 Sep, 2020 DOI:10.21203/rs.3.rs-65529/v1
3. K. Narbute, et al. **Time-Dependent Memory and Gait Improvement by Intranasally-Administered Extracellular Vesicles in Parkinson's Disease Model Rats.** *Cellular and Molecular Neurobiology*, 5 May 2020 <https://doi.org/10.1007/s10571-020-00865-8>
4. X. Wang, et al. **Exosomes influence the behavior of human mesenchymal stem cells on titanium surfaces.** *Biomaterials*, 24 October 2019, <https://doi.org/10.1016/j.biomaterials.2019.119571>
5. J.W Li, et al. **Mesenchymal stromal cells-derived exosomes alleviate ischemia/reperfusion injury in mouse lung by transporting anti-apoptotic miR-21-5p.** *European Journal of Pharmacology*, 2019
6. Z. Zhu, et al. **Exosomes derived from human umbilical cord mesenchymal stem cells accelerate growth of VK2 vaginal epithelial cells through MicroRNAs in vitro.** *Human Reproduction*, dey344, <https://doi.org/10.1093/humrep/dey344>, 2018
7. T.C. Peak, et al. **Exosomes secreted by placental stem cells selectively inhibit growth of aggressive prostate cancer cells.** *Biochemical and Biophysical Research Communications*, Volume 499, Issue 4, 23 May 2018, Pages 1004-1010.
8. A. Jarmalavičiūtė, et al. **Exosomes from dental pulp stem cells rescue human dopaminergic neurons from 6-hydroxy-dopamine-induced apoptosis.** *Cytotherapy*, 2015
9. U Pivoraitė, et. al. **Exosomes from Human Dental Pulp Stem Cells Suppress Carrageenan-Induced Acute Inflammation in Mice.** *Inflammation*, April 2015
10. K. Narbute, et al. **Intranasal Administration of Extracellular Vesicles Derived from Human Teeth Stem Cells Improve Motor Symptoms and Normalize Tyrosine Hydroxylase Expression in the Substantia Nigra and Striatum of the 6-Hydroxydopamine-Treated Rats.** *Stem Cells Translational Medicine*, pp. 1-10, 2019
11. U. Jonavičė, et al. **Extracellular vesicles can act as a potent immunomodulators of human microglial cells.** *Journal of Tissue Engineering and Regenerative Medicine*, 2019
12. S. Bobis-Wozowicz, et al. **Diverse impact of xeno-free conditions on biological and regenerative properties of hUC-MSCs and their extracellular vesicles.** *Journal of Molecular Medicine*, 2016
13. M. Pokrywczynska, et al. **Transdifferentiation of Bone Marrow Mesenchymal Stem Cells into the Islet-Like Cells: the Role of Extracellular Matrix Proteins.** *Archivum Immunologiae et Therapiae Experimentalis*, May 2015

Publications

1. M.G. Svahn, **ALLOGENEIC COMPOSITION**. US Patent App. 16/969,558, 2021
2. S. Bhat, et al. **Expansion and characterization of bone marrow derived human mesenchymal stromal cells in serum-free conditions**. *Sci Rep* 11, 3403 (2021). <https://doi.org/10.1038/s41598-021-83088-1>
3. N. Aydoğdu, et al. **Isolation, Culture, Cryopreservation, and Preparation of Umbilical Cord-Derived Mesenchymal Stem Cells as a Final Cellular Product Under Good Manufacturing Practices-Compliant Conditions**. *Methods in Molecular Biology*. (2020) Springer, New York, NY. https://doi.org/10.1007/7651_2020_332
4. N. Aydoğdu, et al. **Isolation, Culture, Cryopreservation, and Preparation of Skin-Derived Fibroblasts as a Final Cellular Product Under Good Manufacturing Practice-Compliant Conditions**. *Methods in Molecular Biology*. (2020) Springer, New York, NY. https://doi.org/10.1007/7651_2020_333
5. J. Jeriha, et al. **mRNA-Based Reprogramming Under Xeno-Free and Feeder-Free Conditions**. *Methods in Molecular Biology*, 22 June 2020. DOI https://doi.org/10.1007/7651_2020_302
6. S. Suresh, et al. **A nanocomposite hydrogel delivery system for mesenchymal stromal cell secretome**. *Stem Cell Res Ther* 11, 205 (2020). <https://doi.org/10.1186/s13287-020-01712-9>
7. V. Alonso-Camino & B. Mirsch, **Development of standard protocol for the cGMP production of human mesenchymal stem/stromal cells from different source tissues**. *Cytotherapy* Volume 22, Issue 5, Supplement, May 2020, Page S565-S566, <https://doi.org/10.1016/j.jcyt.2020.03.099>
8. A. López Fernández, et al. **Successful scale up expansion of Wharton's jelly mesenchymal stromal cells in different commercial xeno-free and serum-free media**. *Cytotherapy* Volume 22, Issue 5, Supplement, May 2020, Page S94, <https://doi.org/10.1016/j.jcyt.2020.03.162>
9. B. Brinkhof, et al. **ALCAM (CD166) as a gene expression marker for human mesenchymal stromal cell characterisation**. *Gene*: X 14 March 2020, 100031, <https://doi.org/10.1016/j.gene.2020.100031>
10. S. Guan, et al. **Indium-111-labeled CD166-targeted peptide as a potential nuclear imaging agent for detecting colorectal cancer stem-like cells in a xenograft mouse model**. *EJNMMI Res* 10, 13 (2020). <https://doi.org/10.1186/s13550-020-0597-3>
11. S. Mount, et al. **Physiologic expansion of human heart-derived cells enhances therapeutic repair of injured myocardium**. *Stem Cell Research & Therapy* (2019) 10:316
12. S. R. Cohen, et al. **Cellular Optimization of Nanofat: Comparison of Two Nanofat Processing Devices in Terms of Cell Count and Viability**. *Aesthetic Surgery Journal Open Forum*, <https://doi.org/10.1093/asjof/ojz028> Published: 30 September 2019
13. D. Boruckowski, et al. **Wharton's Jelly Mesenchymal Stem Cell Administration Improves Quality of Life and Self-Sufficiency in Children with Cerebral Palsy: Results from a Retrospective Study**. *Hindawi Stem Cells International* Volume 2019, Article ID 7402151, 13 pages, <https://doi.org/10.1155/2019/7402151>
14. G. Sagaradze, et al. **Conditioned Medium from Human Mesenchymal Stromal Cells: Towards the Clinical Translation**. *Int.J.Mol.Sci.* 2019, 20, 1656; doi:10.3390/ijms20071656
15. M. Valitsky, et al. **Cerebrospinal Fluid (CSF) Exchange with Artificial CSF Enriched with Mesenchymal Stem Cell Secretions Ameliorates Experimental Autoimmune Encephalomyelitis**. *Int.J.Mol. Sci.*, 2019, 20(7),1793; <https://doi.org/10.3390/ijms20071793>
16. Y. Ben, et al. **STABILIZED AMORPHOUS CALCIUM CARBONATE FOR TREATMENT OF NEUROLOGICAL, MUSCULAR AND INFERTILITY DISEASES OR CONDITIONS**. US Patent App. 16/069,762, 2019
17. J. Favaloro, et al. **Evaluation of human platelet lysate (HPL) as a substitute for foetal bovine serum (FBS) and recombinant proteins for the growth and maintenance of bone marrow derived mesenchymal stromal cells (BM-MSc)**. *Cytotherapy*, Volume 21, Issue 5, Supplement, May 2019, Pages S84-S85
18. G. Chew, et al. **Tide motion bioreactors for large-scale cultivation and expansion of human mesenchymal stem cells**. *Cytotherapy*, Volume 21, Issue 5, Supplement, May 2019, Page S85
19. V. Alonso-Camino and B. Mirsch, **Rapid expansion of Mesenchymal Stem/Stromal Cells using optimized media supplemented with human platelet lysate PLTMax® or PLTGold®, suitable for cGMP expansion at large scale**. *Cytotherapy*, Volume 21, Issue 5, Supplement, May 2019, Page S85
20. S. Pang, et al. **Dissecting the molecular pathways of apoptosis in mesenchymal stromal cell therapy**. *Cytotherapy*, Volume 21, Issue 5, Supplement, May 2019, Page S85
21. M. Valitsky, et al. **Cerebrospinal Fluid (CSF) Exchange with Artificial CSF Enriched with Mesenchymal Stem Cell Secretions Ameliorates Experimental Autoimmune Encephalomyelitis**. *Int. J. Mol. Sci.* 2019,

- 20(7), 1793; <https://doi.org/10.3390/ijms20071793>
22. A. Yamasaki, et al. **Osteochondral regeneration using constructs of mesenchymal stem cells made by bio three-dimensional printing in mini-pigs.** *Journal of Orthopaedic Research*, 2018
 23. N.B. Vu, et al. **Off-the-Shelf Mesenchymal Stem Cell Technology.** *Stem Cell Drugs - A New Generation of Biopharmaceuticals.* (2018) https://doi.org/10.1007/978-3-319-99328-7_7
 24. J.K. Ledwon, et al. **Osteogenic Differentiation Of Msc As A Model Study Of The Postnatal Progressive Crouzon Syndrome.** *Plastic and Reconstructive Surgery - Global Open.* 6(4S):116, APR 2018
 25. D.Lisini, et al. **Adipose tissue-derived mesenchymal stromal cells for clinical application: An efficient isolation approach.** *Current Research in Translational Medicine*, <https://doi.org/10.1016/j.retram.2018.06.002>
 26. S. Khan, et al. **cGMP-Compatible Large-Scale Production of Mesenchymal Stem Cells (MSCs) In Xeno- and Serum-Free Media for Allogeneic Cell Therapies.** *Cytotherapy*, Volume 20, Issue 5, Supplement, Page S42, 2018
 27. A.C. Volz and P.J. Kluger, **Completely serum-free and chemically defined adipocyte development and maintenance.** *Cytotherapy*, April 2018 Volume 20, Issue 4, Pages 576-588
 28. C. Elseberg, et al. **The Challenge of Human Mesenchymal Stromal Cell Expansion: Current and Prospective Answers.** *New Insights into Cell Culture Technology*, Dr. Sivakumar Joghi Thatha Gowder (Ed.), InTech.
 29. C. Ceccaliti, et al. **Optimization of Injectable Thermosensitive Scaffolds with Enhanced Mechanical Properties for Cell Therapy.** *Macromolecular Bioscience*, 2017
 30. D. Boruckowski, et al. **Third-party Wharton's jelly mesenchymal stem cells for treatment of steroid-resistant acute and chronic graft-versus-host disease: a report of 10 cases.** *Turkish Journal of Biology*, 40: 493-500, 2016
 31. H. Jianxia, et al. **Long term effect and safety of Wharton's jelly-derived mesenchymal stem cells on type 2 diabetes.** *Experimental and Therapeutic Medicine*, Volume 12 Issue 3, 2016
 32. L. Berger, et al. **Tumor Specific Recruitment and Reprogramming of Mesenchymal Stem Cells in Tumorigenesis.** *STEM CELLS* Volume 34, Issue 4, Version of Record online: 31 DEC 2015
 33. K.Y. Tan, et al. **Serum-free media formulations are cell line-specific and require optimization for microcarrier culture.** *Cytotherapy*, 2015
 34. C. Zhen, et al. **Chondrogenesis of Human Adipose-Derived Stem Cells by In Vivo Co-graft with Auricular Chondrocytes from Microtia.** *Aesthetic plastic surgery* 39.3 (2015): 431-439.
 35. S.H. Mei, et al. **Isolation and large-scale expansion of bone marrow-derived mesenchymal stem cells with serum-free media under GMP-compliance.** *Cytotherapy*, Volume 16, Issue 4, Supplement, April 2014
 36. M. Genser-Nir, et al. **Toward a serum-free, xeno-free culture system for optimal growth and expansion of hMSC suited to therapeutic applications.** *BMC Proc.* 2013; 7(Suppl 6): P6.
 37. M. J. McVey, et al. **Microparticles as biomarkers of lung disease-enumeration in biological fluids using lipid bilayer microspheres.** *American Journal of Physiology-Lung Cellular and Molecular Physiology* (2016): ajplung-00369.
 38. Y. Lopez, et al. **Identification of Optimal Conditions For Generating MSCs For Preclinical Testing: Comparison of Three Commercial Serum-Free Media And Low-Serum Growth Medium.** From 18th ISCT Annual Meeting, Seattle, USA, 2012.
 39. R. D. Foster et al. **GorodetskyFibrin microbeads loaded with mesenchymal cells support their long-term survival while sealed at room temperature.** *Tissue Eng Part C Methods* 2011, *TISSUE ENGINEERING: Part C*, Volume 17, Number 7, 2011

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