

March 2, 2022

## MSCgo™ Differentiation Media Scientific References

- **MSCgo™ Chondrogenic Differentiation Medium**
- **MSCgo™ Adipogenic Differentiation Medium**
- **MSCgo™ Osteogenic XF Differentiation Medium**
- **MSCgo™ Rapid Osteogenic XF Differentiation Medium**

# MSCgo™ Chondrogenic Differentiation Medium

1. J. Wang, et al. **Trehalose glycopolymers for cryopreservation of tissue-engineered constructs.** *Cryobiology*, 2021, <https://doi.org/10.1016/j.cryobiol.2021.11.004>.
2. V.A. Nikitina, et al. **Cytogenetic Characteristics of Diploid Lines of Mesenchymal Multipotent Stromal Cells.** *Cell Tiss. Biol.* 15, 604–615 (2021). <https://doi.org/10.1134/S1990519X21060146>
3. S. Horikoshi, et al. **Clumps of Mesenchymal Stem Cells/Extracellular Matrix Complexes Generated with Xeno-Free Chondro-Inductive Medium induce Bone Regeneration via Endochondral Ossification.** *Biomedicines* 2021, <https://doi.org/10.3390/biomedicines9101408>
4. C. Siyu, et al. **Single Cell Transcriptome Sequencing Reveals the Potential Mechanism of Heterogeneity in Immunoregulatory Function Between Mesenchymal Stromal Cells.** *research square* 2021, DOI: <https://doi.org/10.21203/rs.3.rs-823639/v1>  
Y. Galat, et al. **Crispr editing of the gli1 first intron abrogates gli1 expression and differentially alters lineage commitment.** *Stem Cells*, January 2021, <https://doi.org/10.1002/stem.3341>
5. 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>
6. S. Kikuchi, et al. **Development of a nasal mucosa-removal model for evaluating cell therapy.** *Regenerative Therapy*, Volume 16, 2021, P. 32–41, ISSN 2352-3204, <https://doi.org/10.1016/j.reth.2020.12.004>.
7. J. Min, et al. **Phenotype and biological characteristics of endometrial mesenchymal stem / stromal cells: A comparison between intrauterine adhesion patients and healthy women.** *AJRI*, 18 November 2020, <https://doi.org/10.1111/aji.13379>
8. E. Çerçi & H. Erdost, **Rapid, practical and safe isolation of adipose derived stem cells.** *Biotechnic & Histochemistry*, 23 Jun 2020. DOI: 10.1080/10520295.2020.1776895
9. E. Çerci & H.Erdost, **Stem Cell Identification and Clinical Practice.** *International Journal of Agricultural and Natural Sciences* E-ISSN:2651-3617 12(1): 17-19, 2019
10. O. Ben Menachem - Zidon, et al. **Systemically transplanted mesenchymal stem cells induce vascular-like structure formation in a rat model of vaginal injury.** *plos one*, June 13, 2019 <https://doi.org/10.1371/journal.pone.0218081>
11. M.Y. Meng, et. al. **Assessment of tumor promoting effects of amniotic and umbilical cord mesenchymal stem cells in vitro and in vivo.** *J Cancer Res Clin Oncol* (2019) 145: 1133. <https://doi.org/10.1007/s00432-019-02859-6>
12. R. Moloudi, et al. **Inertial-Based Filtration Method for Removal of Microcarriers from Mesenchymal Stem Cell Suspensions.** *Scientific Reports* volume 8, Article number: 12481 (2018)
13. J. Leber, et al. **Microcarrier choice and bead-to-bead transfer for human mesenchymal stem cells in serum-containing and chemically defined media.** *Process Biochemistry*, volume 59, Part B, August 2017, Pages 255-265
14. L. Pu, et al. **Compared to the amniotic membrane, Wharton’s jelly may be a more suitable source of mesenchymal stem cells for cardiovascular tissue engineering and clinical regeneration.** *Stem Cell Research & Therapy* 2017 8:72

# MSCgo™ Adipogenic Differentiation Medium

1. Y. Xiong, et al. **Effects of different methods of demineralized dentin matrix preservation on the proliferation and differentiation of human periodontal ligament stem cells.** *Journal of Dental Sciences*, 2022, <https://doi.org/10.1016/j.jds.2022.01.007>.
2. E. Atanasova, et al. **Normal ex vivo mesenchymal stem cell function combined with abnormal immune profiles sets the stage for informative cell therapy trials in idiopathic pulmonary fibrosis patients.** *Stem Cell Res Ther*, 2022. <https://doi.org/10.1186/s13287-021-02692-0>
3. J. Wang, et al. **Trehalose glycopolymers for cryopreservation of tissue-engineered constructs.** *Cryobiology*, 2021, <https://doi.org/10.1016/j.cryobiol.2021.11.004>.
4. V.A. Nikitina, et al. **Cytogenetic Characteristics of Diploid Lines of Mesenchymal Multipotent Stromal Cells.** *Cell Tiss. Biol.* 15, 604–615 (2021). <https://doi.org/10.1134/S1990519X21060146>
5. C. Siyu, et al. **Single Cell Transcriptome Sequencing Reveals the Potential Mechanism of Heterogeneity in Immunoregulatory Function Between Mesenchymal Stromal Cells.** *research square* 2021, DOI: <https://doi.org/10.21203/rs.3.rs-823639/v1>
6. E. Atanasova, et al. **Autologous Mesenchymal Stem Cell Therapy for Idiopathic Pulmonary Fibrosis and Comprehensive Assessment of Circulating Immune Populations Cells.** *Mayo Clinic College of Medicine*. 2021 <https://doi.org/10.21203/rs.3.rs-570630/v1>
7. Y. Galat, et al. **Crispr editing of the gli1 first intron abrogates gli1 expression and differentially alters lineage commitment.** *Stem Cells*, January 2021, <https://doi.org/10.1002/stem.3341>
8. 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>
9. S. Kikuchi, et al. **Development of a nasal mucosa-removal model for evaluating cell therapy.** *Regenerative Therapy*, Volume 16, 2021, P. 32-41, ISSN 2352-3204, <https://doi.org/10.1016/j.reth.2020.12.004>.
10. J. Min, et al. **Phenotype and biological characteristics of endometrial mesenchymal stem / stromal cells: A comparison between intrauterine adhesion patients and healthy women.** *AJRI*, 18 November 2020, <https://doi.org/10.1111/aji.13379>
11. L. Xie, et al. **Alteration of circRNA and lncRNA expression profile in exosomes derived from periodontal ligament stem cells undergoing osteogenic differentiation.** *Archives of Oral Biology*, 2020, 104984, <https://doi.org/10.1016/j.archoralbio.2020.104984>.
12. G. Çiçek, et al. **Examination of Adipose Tissue-derived Mesenchymal Stem Cell Surface Markers in a Hypoxic Environment.** *Cell Tiss. Biol.* 14, 325–331 (2020). <https://doi.org/10.1134/S1990519X20050028>
13. E. Çerçi & H. Erdost, **Rapid, practical and safe isolation of adipose derived stem cells.** *Biotechnic & Histochemistry*, 23 Jun 2020. DOI:10.1080/10520295.2020.1776895
14. R. Moloudi, et al. **Inertial-Based Filtration Method for Removal of Microcarriers from Mesenchymal Stem Cell Suspensions.** *Scientific Reports* volume 8, Article number: 12481 (2018)
15. A. Mamchur, et al. **Adipose-Derived Stem Cells of Blind Mole Rat Spalax Exhibit Reduced Homing Ability: Molecular Mechanisms and Potential Role in Cancer Suppression.** *Stem Cells*, Volume 36, Issue 10, October 2018, Pages 1630-1642
16. L. Sun, et al. **Human gastric cancer mesenchymal stem cell-derived IL15 contributes to tumor cell EMT via up-regulation of Tregs ratio and PD-1 expression in CD4+T cell.** *Stem Cells and Development*, 2018
17. J. Leber, et al. **Microcarrier choice and bead-to-bead transfer for human mesenchymal stem cells in serum-containing and chemically defined media.** *Process Biochemistry*, volume 59, Part B, August 2017, Pages 255-265
18. L. Pu, et al. **Compared to the amniotic membrane, Wharton's jelly may be a more suitable source of mesenchymal stem cells for cardiovascular tissue engineering and clinical regeneration.** *Stem Cell Research & Therapy* 2017 8:72
19. M. Meng, et al. **Umbilical cord mesenchymal stem cell transplantation in the treatment of multiple sclerosis.** *American Journal of Translational Research*, 2018;10(1):212-223

## MSCgo™ Osteogenic XF Differentiation Medium

## MSCgo™ Rapid Osteogenic XF Differentiation Medium

1. A.V. Kotova, et al. **Comparative Analysis of Dental Pulp and Periodontal Stem Cells: Differences in Morphology, Functionality, Osteogenic Differentiation and Proteome.** *Biomedicines* 2021. <https://doi.org/10.3390/biomedicines9111606>
2. J. Wang, et al. **Trehalose glycopolymers for cryopreservation of tissue-engineered constructs.** *Cryobiology*, 2021, <https://doi.org/10.1016/j.cryobiol.2021.11.004>.
3. V.A. Nikitina, et al. **Cytogenetic Characteristics of Diploid Lines of Mesenchymal Multipotent Stromal Cells.** *Cell Tiss. Biol.* 15, 604–615 (2021). <https://doi.org/10.1134/S1990519X21060146>
4. C. Siyu, et al. **Single Cell Transcriptome Sequencing Reveals the Potential Mechanism of Heterogeneity in Immunoregulatory Function Between Mesenchymal Stromal Cells.** *research square* 2021, DOI: <https://doi.org/10.21203/rs.3.rs-823639/v1>
5. N.I. Erukashvily, et al. **Fibrin Glue Implants Seeded with Dental Pulp and Periodontal Ligament Stem Cells for the Repair of Periodontal Bone Defects: A Preclinical Study.** *Bioengineering* 2021, <https://doi.org/10.3390/bioengineering8060075>
6. L. Zhang, et al. **LGALS3BP/Gal-3 promotes osteogenic differentiation of human periodontal ligament stem cells.** *Archives of Oral Biology*, 2021, <https://doi.org/10.1016/j.archoralbio.2021.105149>.
7. Y. Galat, et al. **Crispr editing of the gli1 first intron abrogates gli1 expression and differentially alters lineage commitment.** *Stem Cells*, January 2021, <https://doi.org/10.1002/stem.3341>
8. 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>
9. S. Kikuchi, et al. **Development of a nasal mucosa-removal model for evaluating cell therapy.** *Regenerative Therapy*, Volume 16, 2021, P. 32–41, ISSN 2352-3204, <https://doi.org/10.1016/j.reth.2020.12.004>.
10. J. Min, et al. **Phenotype and biological characteristics of endometrial mesenchymal stem / stromal cells: A comparison between intrauterine adhesion patients and healthy women.** *AJRI*, 18 November 2020, <https://doi.org/10.1111/aji.13379>
11. R. Moloudi, et al. **Inertial-Based Filtration Method for Removal of Microcarriers from Mesenchymal Stem Cell Suspensions.** *Scientific Reports* volume 8, Article number: 12481 (2018)
12. L. Sun, et al. **Human gastric cancer mesenchymal stem cell-derived IL15 contributes to tumor cell EMT via up-regulation of Tregs ratio and PD-1 expression in CD4+T cell.** *Stem Cells and Development*, 2018
14. J. Leber, et al. **Microcarrier choice and bead-to-bead transfer for human mesenchymal stem cells in serum-containing and chemically defined media.** *Process Biochemistry*, volume 59, Part B, August 2017, Pages 255–265
15. L. Pu, et al. **Compared to the amniotic membrane, Wharton’s jelly may be a more suitable source of mesenchymal stem cells for cardiovascular tissue engineering and clinical regeneration.** *Stem Cell Research & Therapy* 2017 8:72
16. M. Meng, et al. **Umbilical cord mesenchymal stem cell transplantation in the treatment of multiple sclerosis.** *American Journal of Translational Research*, 2018;10(1):212–223
17. S. Uchida, et al. **mRNA as a Tool for Gene Transfection in 3D Cell Culture for Future Regenerative Therapy.** *Micromachines* 2020, 11, 426; doi:10.3390/mi11040426 [www.mdpi.com/journal/micromachines](http://www.mdpi.com/journal/micromachines)
18. A. Picken, et al. **A Monte Carlo framework for managing biological variability in manufacture of autologous cell therapy from mesenchymal stromal cells therapies.** *Cytotherapy* 27 February 2020 <https://doi.org/10.1016/j.jcyt.2020.01.006>
19. Y.T. Hsiao, et al. **Neuromedin U (NMU) regulates osteoblast differentiation and activity.** *Biochemical and Biophysical Research Communications* Available online 11 February 2020, <https://doi.org/10.1016/j.bbrc.2020.02.003>
20. 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>
21. 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

**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