

BIBLIOGRAPHY

To order a copy of any document listed below please highlight the required article and email to us at:
biomedicals.support@upm.com

SCIENTIFIC PAPERS

- **Vallin, J. and J. Grantham, (2021).** Functional assessment of the V390F mutation in the CCT δ subunit of chaperonin containing tailless complex polypeptide 1. *Cell Stress and Chaperones*. Available from: <https://doi.org/10.1007/s12192-021-01237-x>.
- **Leppiniemi, J., et al., (2021).** Avidin-Conjugated Nanofibrillar Cellulose Hydrogel Functionalized with Biotinylated Fibronectin and Vitronectin Promotes 3D Culture of Fibroblasts. *Biomacromolecules*, 22(10): p. 4122-4137. Available from: <https://doi.org/10.1021/acs.biomac.1c00579>.
- **Zini, J., et al., (2021).** Drug diffusivities in nanofibrillar cellulose hydrogel by combined time-resolved Raman and fluorescence spectroscopy. *Journal of Controlled Release*, 334: p. 367-375. Available from: <https://www.sciencedirect.com/science/article/pii/S016836592100198X>.
- **Zhang, X., et al., (2021).** Effect of laminin, polylysine and cell medium components on the attachment of human hepatocellular carcinoma cells to cellulose nanofibrils analyzed by surface plasmon resonance. *Journal of Colloid and Interface Science*, 584: p. 310-319. Available from: <https://www.sciencedirect.com/science/article/pii/S0021979720312698>.
- **Posch, W., et al., (2021).** C5aR inhibition of non-immune cells suppresses inflammation and maintains epithelial integrity in SARS-CoV-2-infected primary human airway epithelia. *J Allergy Clin Immunol*. Available from: <https://pubmed.ncbi.nlm.nih.gov/33852936/>.
- **Koivunotko, E., et al., (2021).** Molecular Insights on Successful Reconstitution of Freeze-Dried Nanofibrillated Cellulose Hydrogel. *ACS Applied Bio Materials*. Available from: <https://doi.org/10.1021/acsabm.1c00739>.
- **Phiwchai, I., et al., (2020).** Liver Cancer Cells Uptake Labile Iron via L-type Calcium Channel to Facilitate the Cancer Cell Proliferation. *Cell Biochemistry and Biophysics*. Available from: <https://doi.org/10.1007/s12013-020-00951-0>.

- **Manninen, H., et al., (2020).** Taste compound – Nanocellulose interaction assessment by fluorescence indicator displacement assay. *Food Chemistry*, 318: p. 126511. Available from:
<https://www.sciencedirect.com/science/article/pii/S0308814620303733>.
- **Heuer, R.A., et al., (2020).** Three-Dimensional Otic Neuronal Progenitor Spheroids Derived from Human Embryonic Stem Cells. *Tissue Engineering Part A*. Available from:
<https://doi.org/10.1089/ten.TEA.2020.0078>.
- **Chang, H.-T., et al., (2020).** An engineered three-dimensional stem cell niche in the inner ear by applying a nanofibrillar cellulose hydrogel with a sustained-release neurotrophic factor delivery system. *Acta Biomaterialia*, 108: p. 111-127. Available from: <http://www.sciencedirect.com/science/article/pii/S1742706120301392>.
- **Kögler, M., et al., (2020).** Assessment of recombinant protein production in *E. coli* with Time-Gated Surface Enhanced Raman Spectroscopy (TG-SERS). *Scientific Reports*, 10(1): p. 2472. Available from: <https://doi.org/10.1038/s41598-020-59091-3>.
- **Mäkelä, R., et al., (2020).** Ex vivo modelling of drug efficacy in a rare metastatic urachal carcinoma. *BMC Cancer*, 20(1): p. 590. Available from:
<https://doi.org/10.1186/s12885-020-07092-w>.
- **Auvinen, V.V., et al., (2019).** Effects of nanofibrillated cellulose hydrogels on adipose tissue extract and hepatocellular carcinoma cell spheroids in freeze-drying. *Cryobiology*, 91: p. 137-145. Available from:
<https://www.ncbi.nlm.nih.gov/pubmed/31533026>.
- **Barnawi, R., et al., (2019).** β 1 Integrin is essential for fascin-mediated breast cancer stem cell function and disease progression. *International Journal of Cancer*, 145(3): p. 830-841. Available from: <https://onlinelibrary.wiley.com/doi/abs/10.1002/ijc.32183>.
- **Charbonneau, A.M., et al., (2019).** 3D Culture Histology Cryosectioned Well Insert Technology Preserves the Structural Relationship between Cells and Biomaterials for Time-Lapse Analysis of 3D Cultures. *Biotechnology Journal*, 0(0): p. 1900105. Available from: <https://onlinelibrary.wiley.com/doi/abs/10.1002/biot.201900105>.
- **Harjumäki, R., et al., (2019).** Quantified forces between HepG2 hepatocarcinoma and WA07 pluripotent stem cells with natural biomaterials correlate with in vitro cell behavior. *Scientific Reports*, 9(1): p. 7354. Available from:
<https://doi.org/10.1038/s41598-019-43669-7>.
- **Nugroho, R.W.N., et al., (2019).** Quantifying the interactions between biomimetic biomaterials – collagen I, collagen IV, laminin 521 and cellulose nanofibrils – by

colloidal probe microscopy. *Colloids and Surfaces B: Biointerfaces*, 173: p. 571-580.

Available from:

<http://www.sciencedirect.com/science/article/pii/S0927776518306854>.

- **Sheard, J.J., et al., (2019).** Optically Transparent Anionic Nanofibrillar Cellulose Is Cytocompatible with Human Adipose Tissue-Derived Stem Cells and Allows Simple Imaging in 3D. *Stem Cells International*, 2019: p. 12. Available from: <https://doi.org/10.1155/2019/3106929>.
- **Zaderer, V., et al., (2019).** Turning the World Upside-Down in Cellulose for Improved Culturing and Imaging of Respiratory Challenges within a Human 3D Model. *Cells*, 8(10): p. 1292. Available from: <https://www.ncbi.nlm.nih.gov/pmc/articles/PMC6830077/>.
- **Broguiere, N., et al., (2018).** Growth of Epithelial Organoids in a Defined Hydrogel. *Advanced Materials*, 30(43): p. 1801621. Available from: <https://onlinelibrary.wiley.com/doi/abs/10.1002/adma.201801621>.
- **Laurén, P., et al., (2018).** Pectin and Mucin Enhance the Bioadhesion of Drug Loaded Nanofibrillated Cellulose Films. *Pharmaceutical Research*, 35(7): p. 145. Available from: <https://doi.org/10.1007/s11095-018-2428-z>.
- **Nuutila, K., et al., (2018).** Inhibition of Skin Wound Contraction by Nanofibrillar Cellulose Hydrogel. *Plastic and Reconstructive Surgery*, 141(3). Available from: https://journals.lww.com/plasreconsurg/Fulltext/2018/03000/Inhibition_of_Skin_Wound_Contraction_by.18.aspx.
- **O'Donnell, N., et al., (2018).** Cellulose-based scaffolds for fluorescence lifetime imaging-assisted tissue engineering. *Acta Biomaterialia*, 80: p. 85-96. Available from: <http://www.sciencedirect.com/science/article/pii/S1742706118305634>.
- **Azoidis, I., et al., (2017).** Three-dimensional cell culture of human mesenchymal stem cells in nanofibrillar cellulose hydrogels. *MRS Communications*, 7(3): p. 458-465. Available from: <https://www.cambridge.org/core/article/threedimensional-cell-culture-of-human-mesenchymal-stem-cells-in-nanofibrillar-cellulose-hydrogels/D907876D60057E3A8E4739900CDAD403>.
- **Laurén, P., et al., (2017).** Nanofibrillar cellulose-alginate hydrogel coated surgical sutures as cell-carrier systems. *PLOS ONE*, 12(8): p. e0183487. Available from: <https://doi.org/10.1371/journal.pone.0183487>.
- **Paukkonen, H., et al., (2017).** Nanofibrillar cellulose hydrogels and reconstructed hydrogels as matrices for controlled drug release. *International Journal of*

Pharmaceutics, 532(1): p. 269-280. Available from:
<http://www.sciencedirect.com/science/article/pii/S0378517317308487>.

- **Paukkonen, H., et al., (2017).** Hydrophobin-nanofibrillated cellulose stabilized emulsions for encapsulation and release of BCS class II drugs. *European Journal of Pharmaceutical Sciences*, 100: p. 238-248. Available from:
<http://www.sciencedirect.com/science/article/pii/S0928098717300532>.
- **Pilapong, C., S. Siriwongnanon, and Y. Keereeta, (2017).** Development of targeted multimodal imaging agent in ionizing radiation-free approach for visualizing hepatocellular carcinoma cells. *Sensors and Actuators B: Chemical*, 245: p. 683-694. Available from:
<http://www.sciencedirect.com/science/article/pii/S0925400517302320>.
- **Reijonen, V., et al., (2017).** Multicellular dosimetric chain for molecular radiotherapy exemplified with dose simulations on 3D cell spheroids. *Physica Medica*, 40: p. 72-78. Available from:
<http://www.sciencedirect.com/science/article/pii/S1120179717302296>.
- **Rinner, B., et al., (2017).** MUG-Mel2, a novel highly pigmented and well characterized NRAS mutated human melanoma cell line. *Sci Rep*, 7(1): p. 2098. Available from:
<http://www.ncbi.nlm.nih.gov/pubmed/28522871>.
- **Hakkarainen, T., et al., (2016).** Nanofibrillar cellulose wound dressing in skin graft donor site treatment. *Journal of Controlled Release*, 244: p. 292-301. Available from:
<http://www.sciencedirect.com/science/article/pii/S0168365916304990>.
- **Mertaniemi, H., et al., (2016).** Human stem cell decorated nanocellulose threads for biomedical applications. *Biomaterials*, 82: p. 208-220. Available from:
<http://www.sciencedirect.com/science/article/pii/S014296121501008X>.
- **Lou, Y.-R., et al., (2015).** Silica bioreplication preserves three-dimensional spheroid structures of human pluripotent stem cells and HepG2 cells. *Scientific Reports*, 5: p. 13635. Available from: <http://dx.doi.org/10.1038/srep13635>.
- **Laurén, P., et al., (2014).** Technetium-99m-labeled nanofibrillar cellulose hydrogel for in vivo drug release. *European Journal of Pharmaceutical Sciences*, 65: p. 79-88. Available from:
<http://www.sciencedirect.com/science/article/pii/S0928098714003704>.
- **Lou, Y.-R., et al., (2014).** The Use of Nanofibrillar Cellulose Hydrogel As a Flexible Three-Dimensional Model to Culture Human Pluripotent Stem Cells. *Stem Cells and*

Development, 23(4): p. 380-392. Available from:
<http://www.ncbi.nlm.nih.gov/pmc/articles/PMC3920753/>.

- **Malinen, M.M., et al., (2014).** Differentiation of liver progenitor cell line to functional organotypic cultures in 3D nanofibrillar cellulose and hyaluronan-gelatin hydrogels. *Biomaterials*, 35(19): p. 5110-5121. Available from:
<http://www.sciencedirect.com/science/article/pii/S0142961214002555>.
- **Bhattacharya, M., et al., (2012).** Nanofibrillar cellulose hydrogel promotes three-dimensional liver cell culture. *Journal of Controlled Release*, 164(3): p. 291-298. Available from:
<http://www.sciencedirect.com/science/article/pii/S0168365912005391>.
- **Kolakovic, R., et al., (2012).** Spray-dried nanofibrillar cellulose microparticles for sustained drug release. *International Journal of Pharmaceutics*, 430(1): p. 47-55. Available from:
<http://www.sciencedirect.com/science/article/pii/S0378517312002724>.