

1

BIOREACTORS

Maintaining organoids in culture allows them to mature and develop, but their growth and lifespan are limited because the larger they get the less access they have to nutrient diffusion, especially for cells in the inner core. Bioreactors—sterile vessels that gently stir contents, maintain temperature, pH, and oxygen levels, and disperse nutrients and biomolecular signals—help researchers control the environmental conditions of brain organoid cultures to maximize their size and lifespan. Researchers are exploring 3D printed and miniature bioreactors as application-specific, low cost, high-throughput, and reproducible alternatives to conventional bioreactors.^{1,2}



2

MICROFLUIDICS

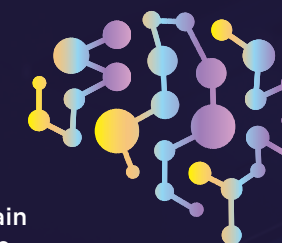
Microfluidic chips enable more precise fluid flow control to further optimize oxygen and nutrient exchange, and generate chemical gradients that guide cell differentiation and migration—addressing the limitation of large fluid volumes used to grow brain organoids in conventional bioreactors.³ These chips can also accommodate complex, patterned co-cultures of various cell types—including nerve, blood vessel, and glial cells—using selective membranes and multiple chambers and channels, which simulate the structure and function of human brains and reduce heterogeneity between samples.^{4,5} In doing so, researchers can create a cellular microenvironment that mimics different brain regions, the blood-brain barrier, and complex cell-cell interactions.



3

BIOMATERIALS

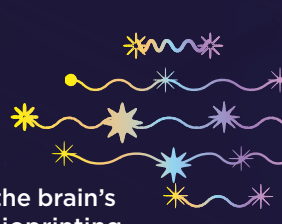
Researchers take advantage of a wide range of biological and synthetic biomaterials to mimic brain tissue's structure and molecular cues. For example, mini brains' growth and cellular patterning depends on the interaction of cells with extracellular matrix—a 3D biological network of molecules and minerals that supports tissue structure and provides biochemical cues to guide cell differentiation, migration, organization, and patterning.⁶ Similarly, scaffolds made of synthetic polymers augment the extracellular environment of cultured stem cells to support 3D growth and organization.^{7,8}



4

3D BIOPRINTING

3D bioprinting is a powerful method for mimicking the brain's tissue architecture with computer modeling and bioprinting technologies (e.g. inkjet, laser-assisted, extrusion-based, or stereolithography).⁹ The ink used for bioprinting organoids contains stem cells plus various biomaterials such as extracellular matrix components that provide cells with appropriate biomolecular cues, architectural scaffolding, and mechanical support. Bioprinting mini brains enables researchers to control the spatial positioning of cell aggregates and ensure that the assembled tissue retains appropriate cell-cell interactions that mimic brain development.¹⁰ Researchers also create structured scaffolds and patterned substrates to encourage stem cells to develop and interact within a topographical map.



5

MICROELECTRODES

As researchers develop increasingly complex brain organoids that exhibit advanced features such as rudimentary brainwaves and the ability to detect basic visual inputs, there is a growing need to characterize and monitor their neural activity. Researchers use specially designed microelectrode arrays—grids of tiny electrodes—to measure the electrical activity of 3D brain organoids.¹¹ Innovative 3D microelectrode arrays use flexible nanowires, novel materials, and unique customizable shapes to overcome the limitations of those traditionally used for 2D cultures by enabling simultaneous recordings across a 3D network, maximizing contact area, and minimizing physical disruption of the organoids.^{12,13,14}



MAJOR ADVANCES IN MINI BRAIN BIOENGINEERING

Sponsored By

SARTORIUS



Empower Your Research With Incucyte® Live-Cell Analysis

The Incucyte® Live-Cell Analysis System speeds scientific discovery by combining lab-tested protocols and reagents with powerful, automated image acquisition and analysis. Gain dynamic insights into the health, morphology, movement and function of cell models, all from the stable environment of a tissue culture incubator. Select an Incucyte® for your research at

www.sartorius.com/incucyte

Specifications subject to change without notice. © 2022. All rights reserved. Incucyte and all names of Sartorius products are registered trademarks and the property of Sartorius AG.

Simplifying Progress

SARTORIUS

SARTORIUS

The Sartorius Group is a leading international partner of biopharmaceutical research and the industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services Division concentrates on serving the needs of laboratories performing research and quality control at pharma and biopharma companies and those of academic research institutes. The Bioprocess Solutions Division with its broad product portfolio focusing on single-use solutions helps customers to manufacture biotech medications and vaccines safely and efficiently.

REFERENCES

1. M.A. Phelan et al., "Mini and customized low-cost bioreactors for optimized high-throughput generation of tissue organoids," *Stem Cell Investig*, 5:33, 2018.
2. X. Qian et al., "Brain-region-specific organoids using mini-bioreactors for modeling ZIKV exposure," *Cell*, 165(5):1238-54, 2016.
3. H. Castiglione et al., "Human brain organoids-on-chip: advances, challenges, and perspectives for preclinical applications," *Pharmaceutics*, 14(11):2301, 2022.
4. Y. Zhu et al., "Engineering human brain assembloids by microfluidics," *Adv Mater*, Jan 12:e2210083, 2023.
5. A.N. Cho et al., "Microfluidic device with brain extracellular matrix promotes structural and functional maturation of human brain organoids," *Nat Commun*, 12(1):4730, 2021.
6. J.M. Unagolla, A.C. Jayasuriya, "Recent advances in organoid engineering: a comprehensive review," *Appl Mater Today*, 29:101582, 2022.
7. G.D. Mahumane et al., "3D scaffolds for brain tissue regeneration: architectural challenges," *Biomater Sci*, 6(11):2812-37, 2018.
8. M.N. Collins et al., "Emerging scaffold- and cellular-based strategies for brain tissue regeneration and imaging," *In Vitro Models*, 1:129-50, 2022.
9. F. Salaris, A. Rosa, "Construction of 3D in vitro models by bioprinting human pluripotent stem cells: challenges and opportunities," *Brain Res*, 1723:146393, 2019.
10. R. Samanipour et al., "A review on 3D printing functional brain model," *Biomicrofluidics*, 16(1):011501, 2022.
11. A. Pelkonen et al., "Functional characterization of human pluripotent stem cell-derived models of the brain with microelectrode arrays," *Cells*, 11(1):106, 2022.
12. T. Kazi, J. Liu, "Emerging bioelectronics for brain organoid electrophysiology," *J Mol Biol*, 434(3):167165, 2022.
13. Q. Huang et al., "Shell microelectrode arrays (MEAs) for brain organoids," *Sci Adv*, 8(33):eabq5031, 2022.
14. D. Soscia et al., "A flexible 3-dimensional microelectrode array for in vitro brain models," *Lab Chip*, 20(5):901-11, 2020.

MAJOR ADVANCES IN MINI BRAIN BIOENGINEERING

Sponsored By

SARTORIUS



More Colors.
More Insights.
More Possibilities.

See more information in every sample with the new Incucyte® SX5 featuring patent-pending optics. Do more with up to 5 colors specifically designed for live-cell analysis.

www.sartorius.com/sx5

Specifications subject to change without notice. © 2022. All rights reserved. Incucyte and all names of Sartorius products are registered trademarks and the property of Sartorius AG.



Simplifying Progress

SARTORIUS