

GrowDex® Supports Human iPSC-derived Neural Progenitor Differentiation *in vitro*

Andrew Schwab, PhD

ORISE Contractor, Research Triangle Park, NC, USA

INTRODUCTION

Three-dimensional (3D) *in vitro* culture of neuronal cells is necessary to recapitulate the complex cytoarchitecture of the central nervous system. As complex, multicellular models develop, there is a need for defined substrates that can allow for the differentiation of iPSC-derived neurons and astrocytes.

Here we establish methodologies for 3D cell culture of human iPSC-derived neural progenitors within nanofibrillar cellulose (NFC) hydrogel, GrowDex®.

MATERIALS

- DMEM/F12 media (Cat No. 11330032, ThermoFisher)
- B27 Supplement (Cat No. 17504044, ThermoFisher)
- GrowDex®, 1.5% (Cat No. 100 103 005, UPM)
- 96-well plate (Cat No. 89626, Ibidi)
- Anti β III Tubulin mAB (Cat No. G7121, Promega)
- Anti GFAP mAB (Cat No. 12389, Cell Signalling Technology)
- Matrigel (Cat No. 354277, Corning)

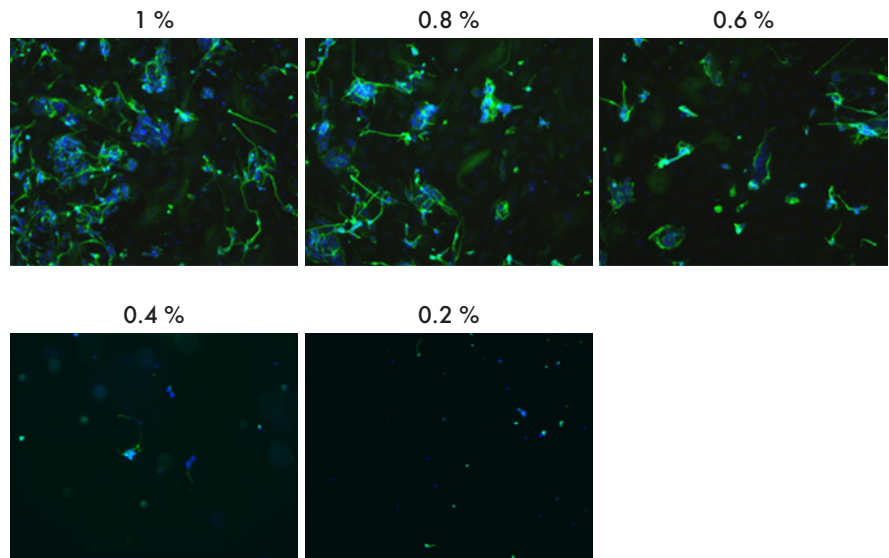
METHODS

- Human induced pluripotent stem cell (iPSC) derived neural progenitors (Ebert et al. 2013) were cultured for one week embedded in GrowDex and on Matrigel.
- Studied GrowDex concentrations were 1.0, 0.8, 0.6, 0.4 and 0.2 wt%.
- GrowDex hydrogel volume was 100 μ l in a Ibidi 96-well plate.
- Matrigel volume was 100 μ l and used at a final concentration of 0.08 mg/ml. The media was aspirated and the cells were plated on top of the Matrigel layer.
- Cell concentration for all conditions was 50,000 cells per well .
- Cells were fed every other day with DMEM/F12 containing 2% B27.
- Half of the media was replaced as to not disturb the hydrogel layer.
- The formation of neural cells inside the hydrogels was evaluated by immunocytochemical staining against β III Tubulin and GFAP.

RESULTS

After one week of differentiation, neural cells were observed by immuno-cytochemistry. Representative 10x images of neurons and astrocytes are shown in Fig. 1. Qualitatively, neural growth increased as GrowDex concentration increased from 0.2% to 1.0%. Cultures in GrowDex resulted in 3D neural structures and outgrowth, whereas cells on Matrigel resulted in neuronal monolayer.

GrowDex



Matrigel

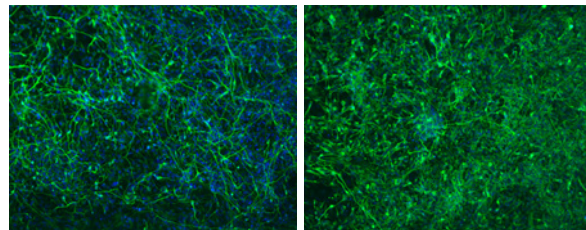


Figure 1. Human iPSC-derived neural progenitor cell aggregates (EZ spheres) differentiate into neurons and astrocytes within GrowDex. Representative images of cells cultured in different GrowDex concentrations (0.2% to 1.0%) showing cell nuclei (blue), differentiation of neural progenitors into β III Tubulin positive neurons (green), and GFAP positive astrocytes (red). In GrowDex cultures, cells were embedded in GrowDex whereas in Matrigel the cells were seeded on top for culture (2 representative images).

CONCLUSIONS

Compared to 2D, differentiation of neural progenitor cells in a 3D environment represents a more physiological approach to model organ-level responses. Neural progenitors grown on Matrigel differentiated into a monolayer of neurons whereas neural progenitor cells grown in GrowDex resulted in neuronal 3D structures. Based on the results, it can be concluded that GrowDex is a good substrate for 3D culture of human iPSC-derived neural progenitor cells and GrowDex concentrations of 0.8% and above are recommended. The biocompatible nature of GrowDex and ease of use in a 96-well plate format, provides a promising alternative scaffold for 3D cell culture. Notably, these results demonstrate that GrowDex could easily be adapted for use in 3D microfluidic neural cell culture models where a defined, animal-free, and tunable matrix is required. As 3D culturing becomes more common, GrowDex will be an important tool that may improve the success of new alternative models.

REFERENCES

Ebert, A.D., et al., EZ spheres: a stable and expandable culture system for the generation of pre-rosette multipotent stem cells from human ESCs and iPSCs. *Stem Cell Res*, 2013. 10(3): p. 417-427.



UPM

UPM Biomedicals

Alvar Aallon katu 1
P.O. Box 380

00101 Helsinki, Finland
biomedicals@upm.com

www.upmbiomedicals.com

www.upm.com

To order products online: www.revivity.com

To request a quote or place an order: www.revivity.com/contact-us

For technical support: LS.ReagentsTechSupport@revivity.com

revivity

An authorized distributor

Revvity, Inc.
940 Winter Street
Waltham, MA 02451
USA
www.revivity.com