

# STEMdiff™ Cerebral Organoid Kit: A New Tool for the Culture of 3-D Brain Organoids Derived from Human Pluripotent Stem Cells

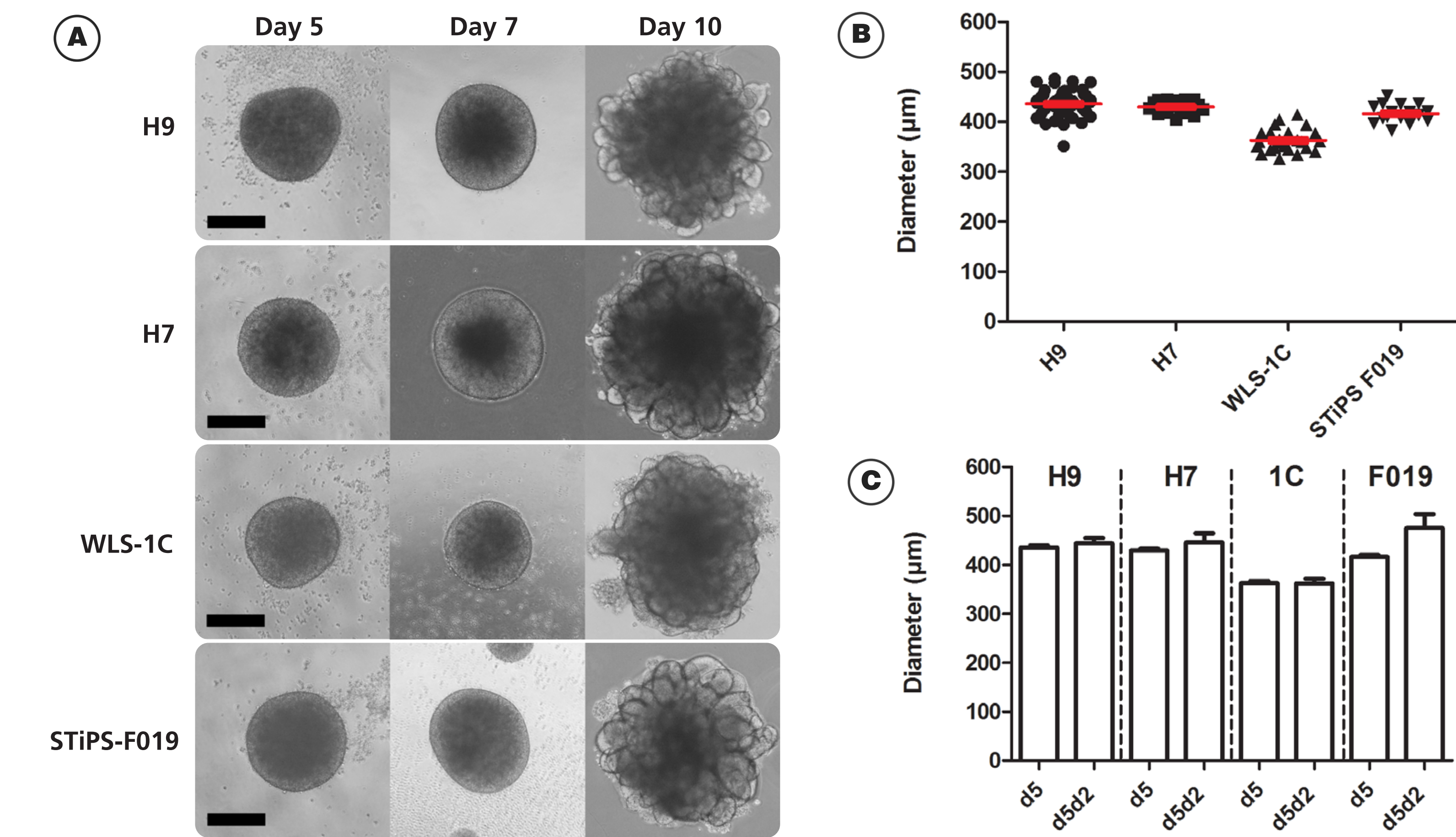
Leon H. Chew<sup>1</sup>, Vivian M. Lee<sup>1</sup>, Adam Añonuevo<sup>1</sup>, Mark Hills<sup>1</sup>, Sam Lloyd-Burton<sup>1</sup>, Terry E. Thomas<sup>1</sup>, Allen C. Eaves<sup>1,2</sup>, and Sharon A. Louis<sup>1</sup>  
<sup>1</sup>STEMCELL Technologies Inc., Vancouver, Canada    <sup>2</sup>Terry Fox Laboratory, BC Cancer Agency, Vancouver, Canada

## Introduction

The metazoan brain is a highly complex yet organized structure. 2-D neural cultures derived from human pluripotent stem cells (hPSCs) are useful models to study the nervous system, but they are limited in their capacity to recapitulate the complex organization of brain tissues. Lancaster et al.<sup>1,2</sup> have recently established an hPSC-based cerebral organoid culture system that models the major processes and structural features of human brain development. Cerebral organoids have important applications in studying human brain development and diseases such as autism, schizophrenia and Zika virus infection<sup>3</sup>. Based on the published formulations, we have developed the STEMdiff™ Cerebral Organoid Kit to enable generation of organoids in a highly reproducible and user-friendly manner. The kit contains 2 basal media and 5 supplements, which are combined separately to prepare complete media for each of the four distinct stages of organoid formation. Cerebral organoids generated with this kit develop progenitor populations that organize into distinct layers giving rise to mature cortical neurons, matching what is observed in the developing human cortex.

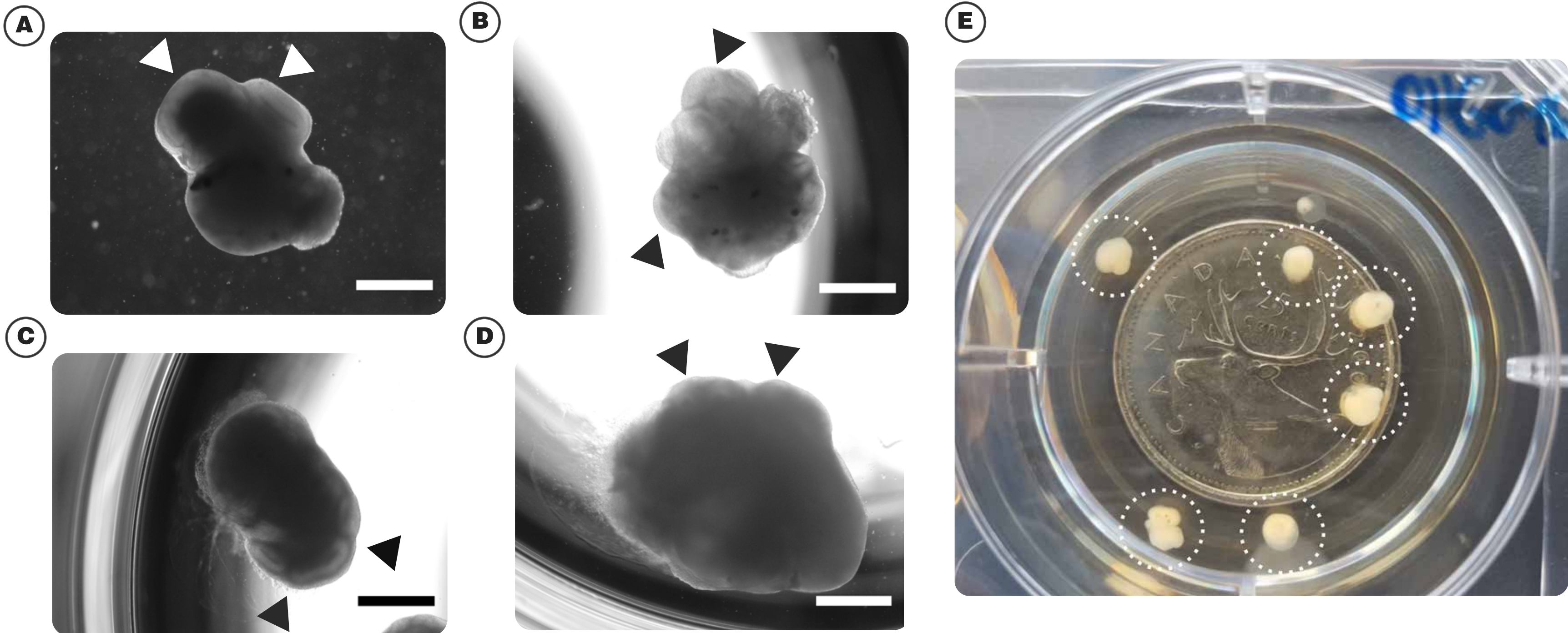
## Results

Figure 1. The STEMdiff™ Cerebral Organoid Kit Supports EB Formation and Neuroepithelia Expansion for Multiple hPSC Lines



(A) Robust early organoid formation was observed in four hPSC lines (2 hESC; H9 and H7, 2 hiPSC; WLS-1C and STiPS-F019) (scale bar = 300 µm). **Day 5:** Embryoid Body Formation; Single-cell suspensions of hPSCs formed spherical aggregates with smooth edges. **Day 7:** Induction stage; EB edges smoothed and developed a translucent quality. **Day 10:** Expansion stage; Matrigel® embedded EBs displayed expanded neuroepithelia evident by bubbling of their surface. Over 80% of all embedded organoids exhibit this type of morphology indicative of neuroepithelia expansion (n = 2 per hPSC line, 8 to 24 organoids measured per hPSC line). (B) Diameter of EBs measured at day 5 shows that all EBs grew to a diameter over 300 µm (n = 2 per cell line, 16 – 48 EBs measured per hPSC line). (C) Measurement of EBs taken at Day 7 (d5d2) show that EBs do not shrink in size when switched to Induction Medium (n = 2 per cell line, 16 – 48 organoids measured per hPSC line).

Figure 2. Day 40 Cerebral Organoids Grow to Over 1 mm in Diameter and Exhibit Tissue Structures with Dense Cores



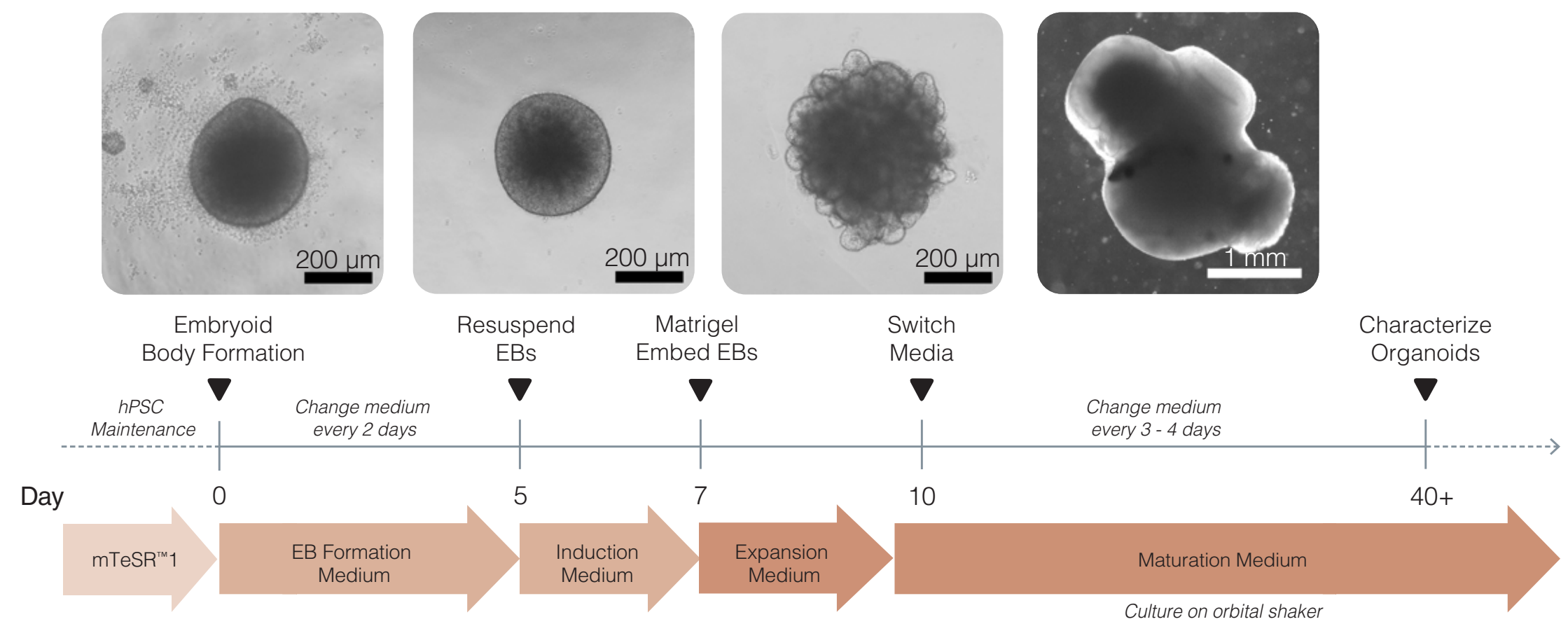
Cerebral organoids cultured in Maturation Medium at day 40 developed thick tissue-like structures and areas that exhibit layering (arrowheads). More than 50% of organoids from each experiment exhibited this type of morphology (n = 2 per cell line, 8 – 24 Organoids measured per hPSC line) (A) H9 (B) H7 (C) WLS-1C (D) STiPS-F019. (E) Size of day 60 organoids (H9). A Canadian quarter is used for scale. Cerebral organoids are outlined (white dotted circles).

## Summary

### STEMdiff™ Cerebral Organoid Kit

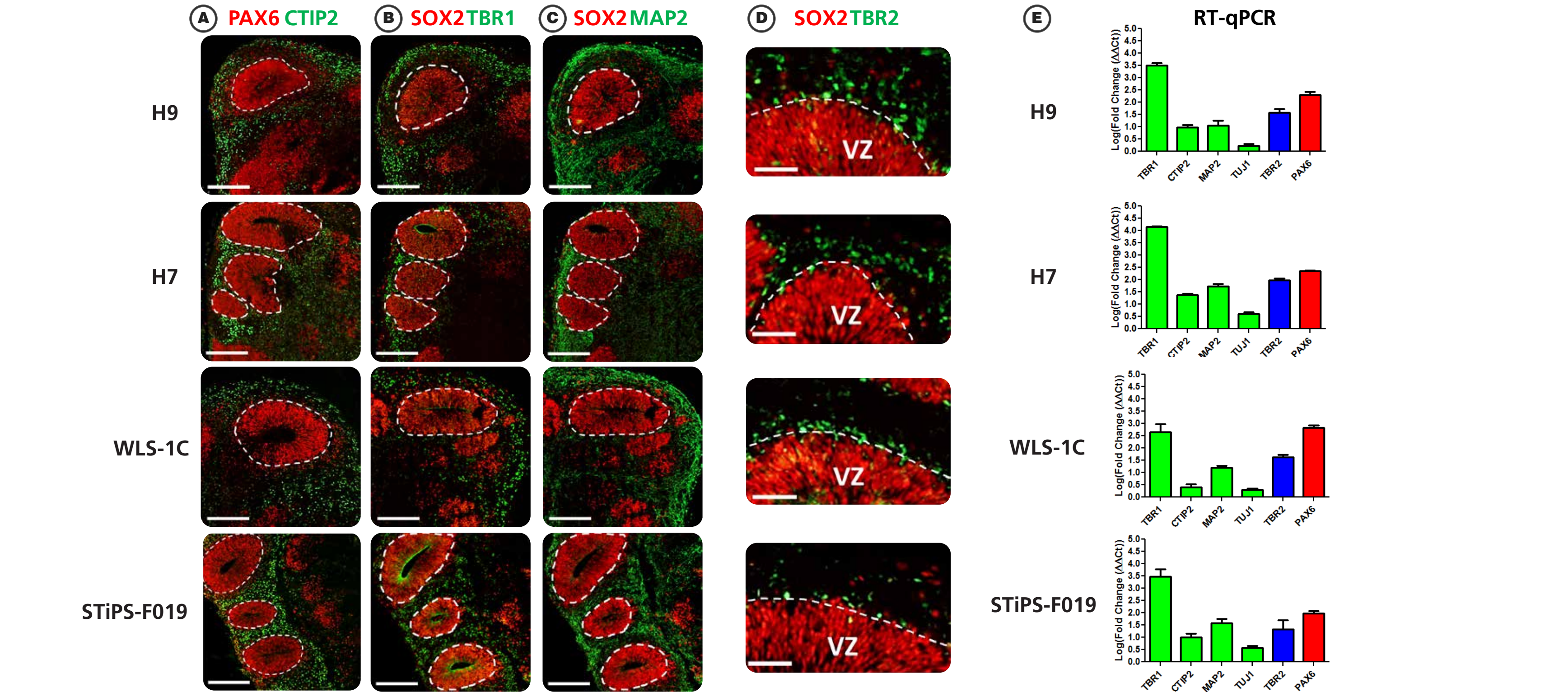
- Supports highly efficient generation of hPSC-derived cerebral organoids from multiple hPSC lines
- Generates organoids with cortical architecture consisting of both neural progenitor and mature neuron populations
- Reproduces cerebral organoids that are consistent with published results by Lancaster et al.<sup>1,2</sup>

## Methods



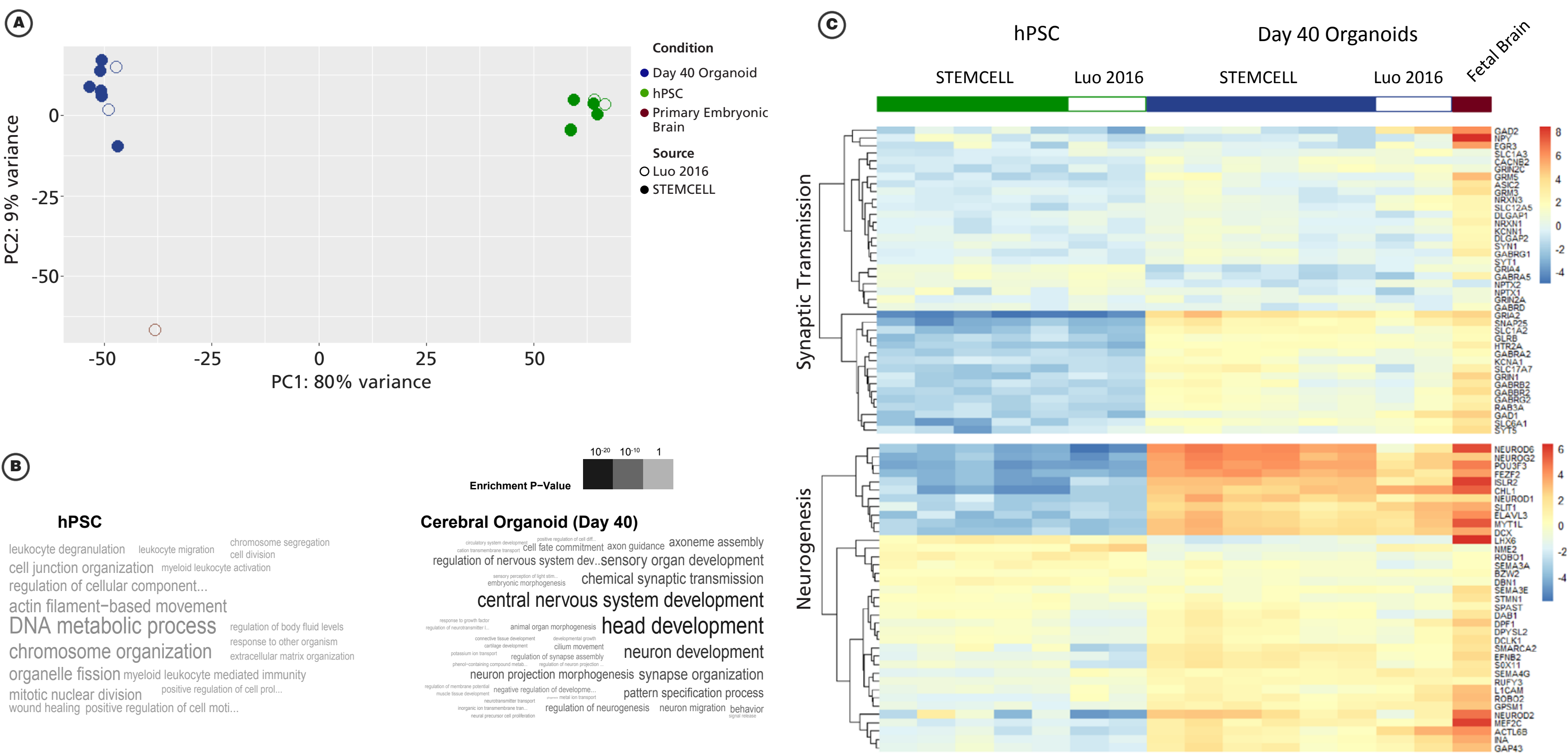
Human pluripotent stem cells (either embryonic or induced pluripotent stem cells) maintained in mTeSR1™ were dissociated into single-cell suspensions using Gentle Cell Dissociation Reagent (GCDR) and seeded at a density of 9,000 cells/well in a U-Bottom 96-well Ultra Low Attachment Plate (Corning) in EB Formation Medium + 10 µM Rho-Kinase Inhibitor (ROCKi). EBs were fed every 2 days with EB Formation Medium without ROCKi. After 5 days, EB diameters were measured from images taken using the Incucyte system (Essen BioScience) and they were transferred to Induction Medium in a 24-well Ultra low Attachment Plate (Corning®). EBs were cultured for an additional 2 days and were then embedded in liquid Matrigel® (Growth Factor Reduced, Corning®) followed by transfer to a non-tissue culture treated 6-well plate (12 -16 organoids/well). Embedded organoids were maintained in Expansion Medium for 3 days. On Day 10, organoids were switched to Maturation Medium and cultured on an orbital shaker set at 57-95 RPM (Infors HT). Organoids were fed every 3 – 4 days with Maturation Medium. On Day 40, organoids were analyzed for neuronal (TBR1, CTIP2, MAP2, beta III Tubulin), intermediate progenitor (ASCL1, TRB2) and neural progenitor (PAX6, SOX2) markers using either RT-qPCR or immunostaining following cryosectioning.

Figure 3. Day 40 Cerebral Organoids Contain Progenitor and Neuron Populations that Organize into Distinct Layers



Immunostaining was performed on 12 µm cryosections of 4% paraformaldehyde fixed organoids for the following markers: (A) PAX6/CTIP2, (B) SOX2/TBR1, (C) SOX2/MAP2, (D) SOX2/TBR2 (scale bar = 200 µm). Neural progenitors (SOX2<sup>+</sup>, PAX6<sup>+</sup>) localized in distinct apical regions surrounding a central ventricle. (E) RT-qPCR analysis showed upregulation of both neural progenitor and mature neuron transcripts as Log(Fold Change ΔΔCt) (Average ± SEM n = 2 per cell line, ≥ 3 organoids per analysis). Data is normalized to 18S/TBP and compared to undifferentiated hPSC control.

Figure 4. RNA-Seq Analysis of Day 40 Organoids Reveals Marker Expression of the Early Developing Human Brain



(A) Principal component analysis of hPSC and cerebral organoid transcriptomes. Cerebral organoids generated using the STEMdiff™ Cerebral Organoid Kit (filled blue circles) cluster together, and cluster with previously published cerebral organoids (open blue circles). The first principal component accounts for the majority of variance seen (PC1; 80 %) and distinguished the cerebral organoid samples from the hPSCs (green circles). The second principal component accounts for only 9 % of the variation, and highlights the modest expression differences between cultured organoids and primary embryonic fetal brain samples (19 post-conceptual weeks, brown circles). (B) Gene ontology terms enriched in Day 40 cerebral organoids with respect to hPSCs reveals transcripts related to brain development and function. (C) Heatmap of expression levels for genes associated with synaptic transmission function and neurogenesis in Day 40 organoids. These data show that gene expression of cerebral organoids generated from the STEMdiff™ Cerebral Organoid Kit are similar to published results.

## References

1. Lancaster, M.A., Renner, M., Martin, C.A., Wenzel, D., Bicknell, L.S., Hurler, M.E., Homfray, T., Penninger, J.M., Jackson, A.P., Knoblich, J.A. (2013). Cerebral organoids model human brain development and microcephaly. Nature, 501(7467), 373–9.
2. Lancaster, M.A. & Knoblich, J.A. (2014). Generation of cerebral organoids from human pluripotent stem cells. Nature Protocols, 9(10), 2329–2340.
3. Kelava, I., & Lancaster, M.A., (2016) Dishing out mini-brains: Current progress and future prospects in brain organoid research. Developmental Biology, 420(2), 199-209
4. Luo, C., Lancaster, M.A., Castanon, R., Nery, J.R., Knoblich, J.A., Ecker, J.R. (2016) Cerebral Organoids Recapitulate Epigenomic Signatures of the Human Fetal Brain. Cell Reports, 17(12), 3369-3384