



Modeling epilepsy-related SCN2A mutation L1342P with CRISPR/Cas9-edited human-induced pluripotent stem cell-derived cortical spheroids

Maria I. Olivero-Acosta¹, Zhefu Que¹, Conrad Otterbacher¹, Muhan Wang¹, Trang Nguyen¹, Ningren Cui¹, Xiaoling Chen¹, Jingliang Zhang¹, Kyle Wettschurack¹, Jiaxiang Wu¹, William C. Skarnes², Yang Yang¹.

¹Department of Medicinal Chemistry and Molecular Pharmacology (MCMP), Purdue Institute for Integrative Neuroscience (PIIN) Purdue University, West Lafayette, 47906 IN. ² The Jackson Lab.



Advancing pharmacogenomics to cure diseases of the nervous system and cancer

Introduction

The SCN2A gene encodes for sodium channel Nav1.2, a protein that mediates action potentials in neurons. SCN2A pathogenic mutations have been associated with **epilepsy**. An example is the **L1342P** mutation, identified in several patients with untreatable seizure episodes (Que, Olivero-Acosta et al., 2021).

3D Structure of Sodium Channel Nav1.2

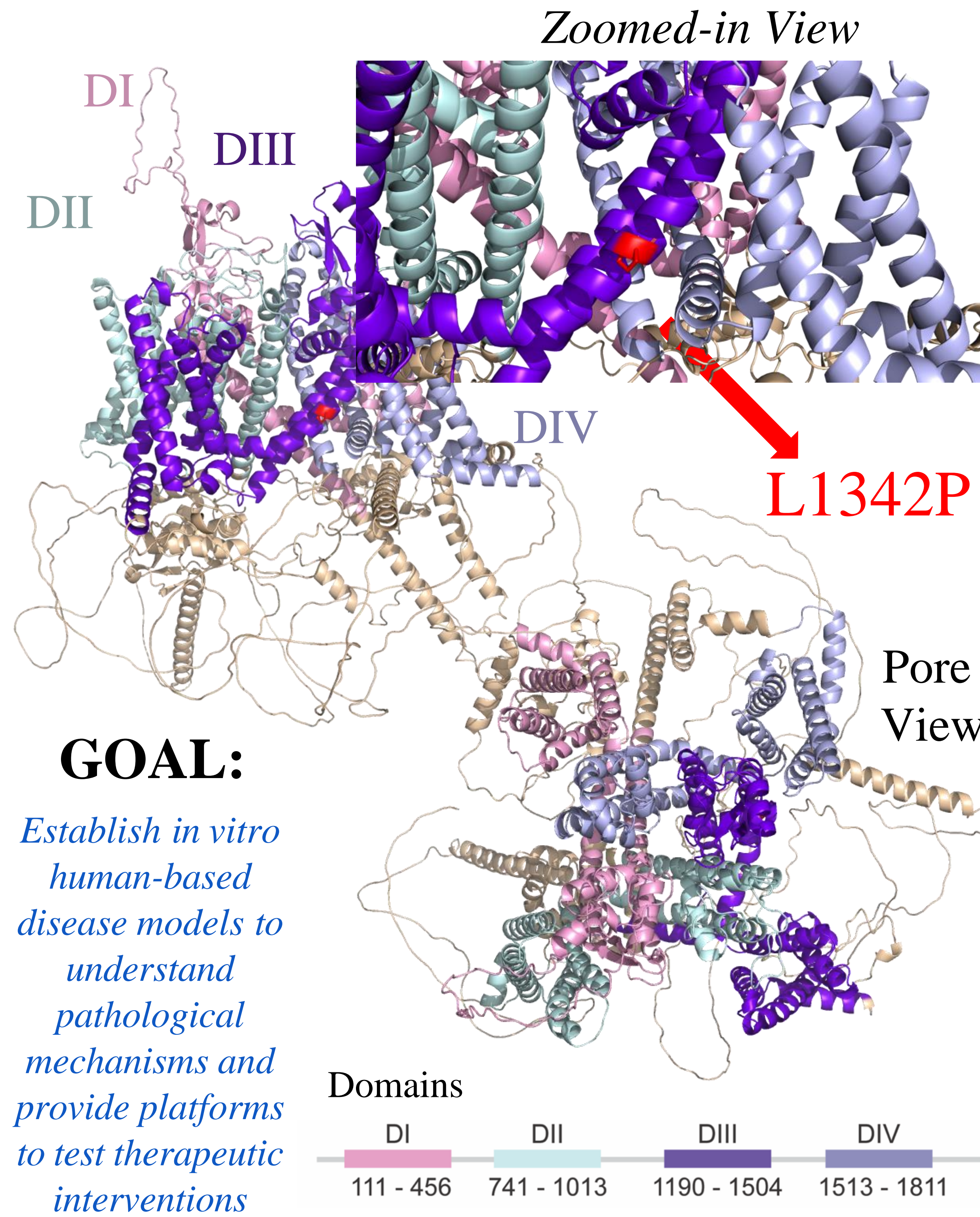
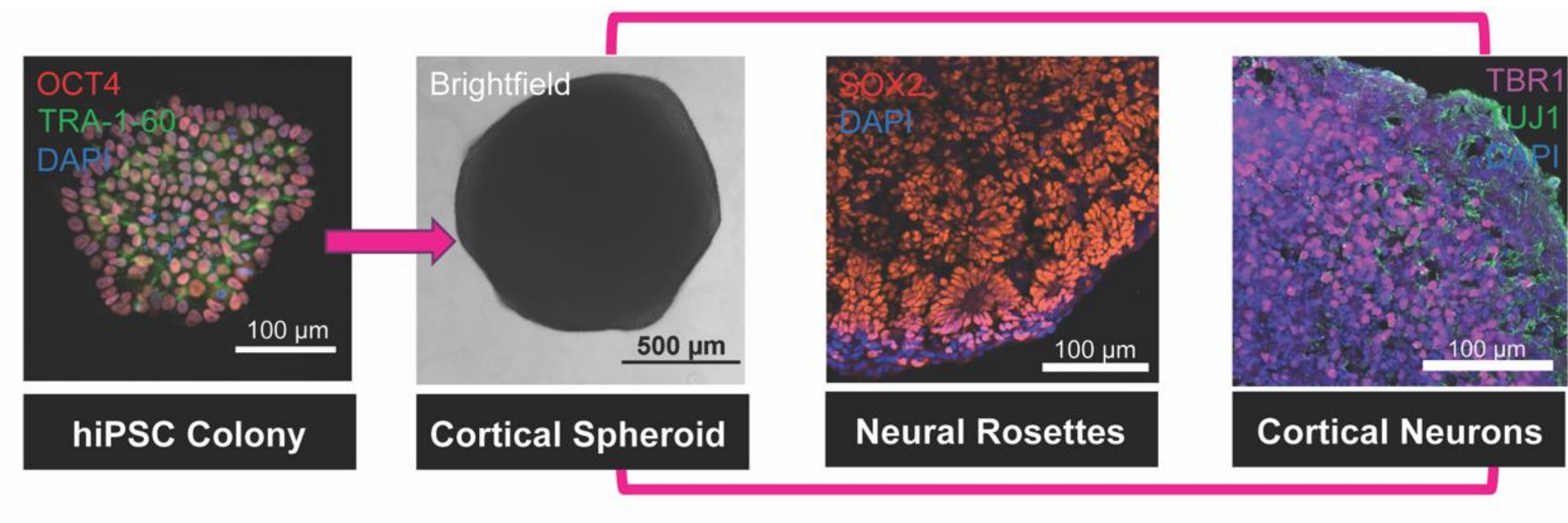


Figure 1. Schematic representation of an AlphaFold predicted folding topology of sodium channel Nav1.2. Structures Rendered using PYMOL. Mutated residue L1342P is indicated in red.

In our recent work, we have demonstrated that hiPSC-derived 2D-neuronal monolayers carrying the CRISPR-Cas9-edited L1342P-mutant channel display a marked hyperexcitability phenotype (Que, Olivero-Acosta et al., 2021). However, the hyperexcitable L1342P mutation's impact on neurodevelopment remains unknown. Cortical spheroids (organoids) are *in-vitro* generated 3D cellular aggregates that resemble the features of the human cortex.

In this poster, we describe the generation of the first SCN2A Cortical Spheroid model, aiming to understand the impact of the L1342P mutation on neuron development and further probe at its characteristic hyperexcitability phenotype.



Starting material, can be a patient-derived or a CRISPR-Cas9 edited cell line.

Neural Cell aggregate can grow up to 4 mm.

Made up of Neural Progenitors, which eventually mature to become neurons.

Arrange themselves in patterns resembling aspects of the prenatal brain.

Figure 2. Representative immunofluorescence and brightfield images of hiPSC-derived cortical spheroids through different maturation stages.

Methods

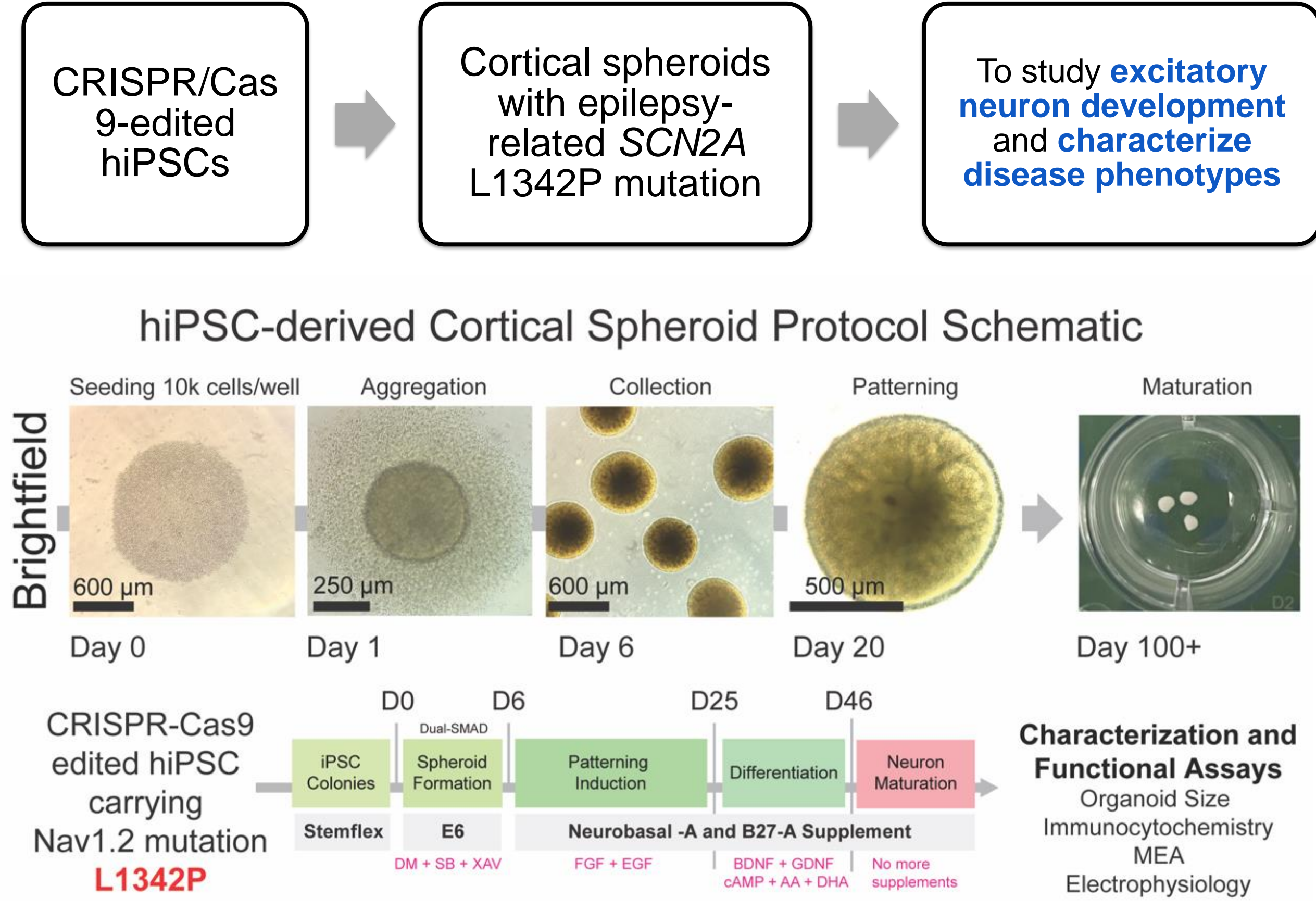


Figure 3. The procedure used to generate human induced pluripotent stem cell-derived cortical spheroids in the Yang Lab. Based on (Sloan et al., 2018).

Results

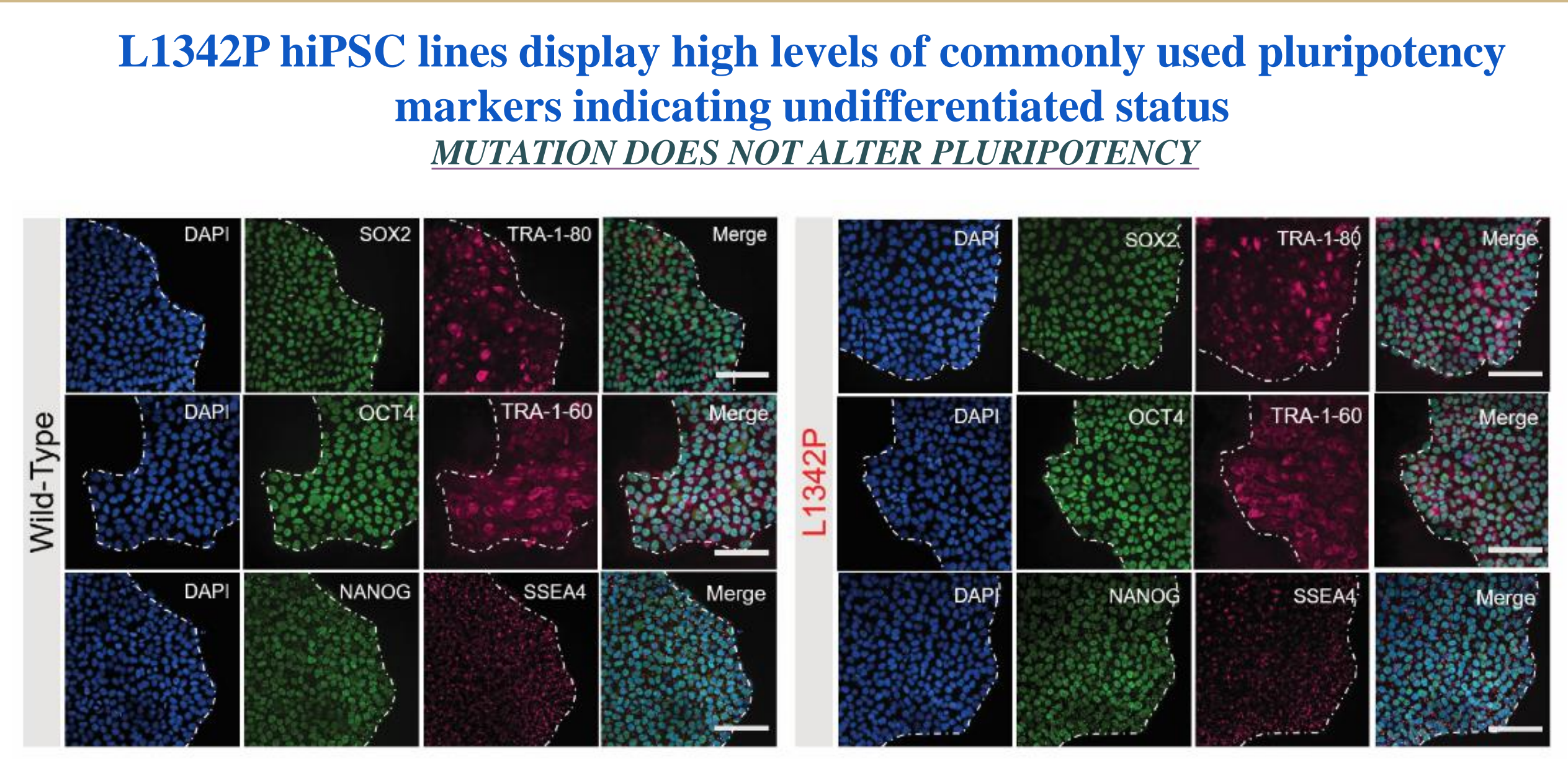


Figure 4. Representative images of pluripotent colonies. Markers include DAPI (nuclei), SRY-Box Transcription Factor 2 (SOX2), Tra-1-80, OCT4, Tra-1-60, NANOG and SSEA4. Scale bar set to 100 µm.

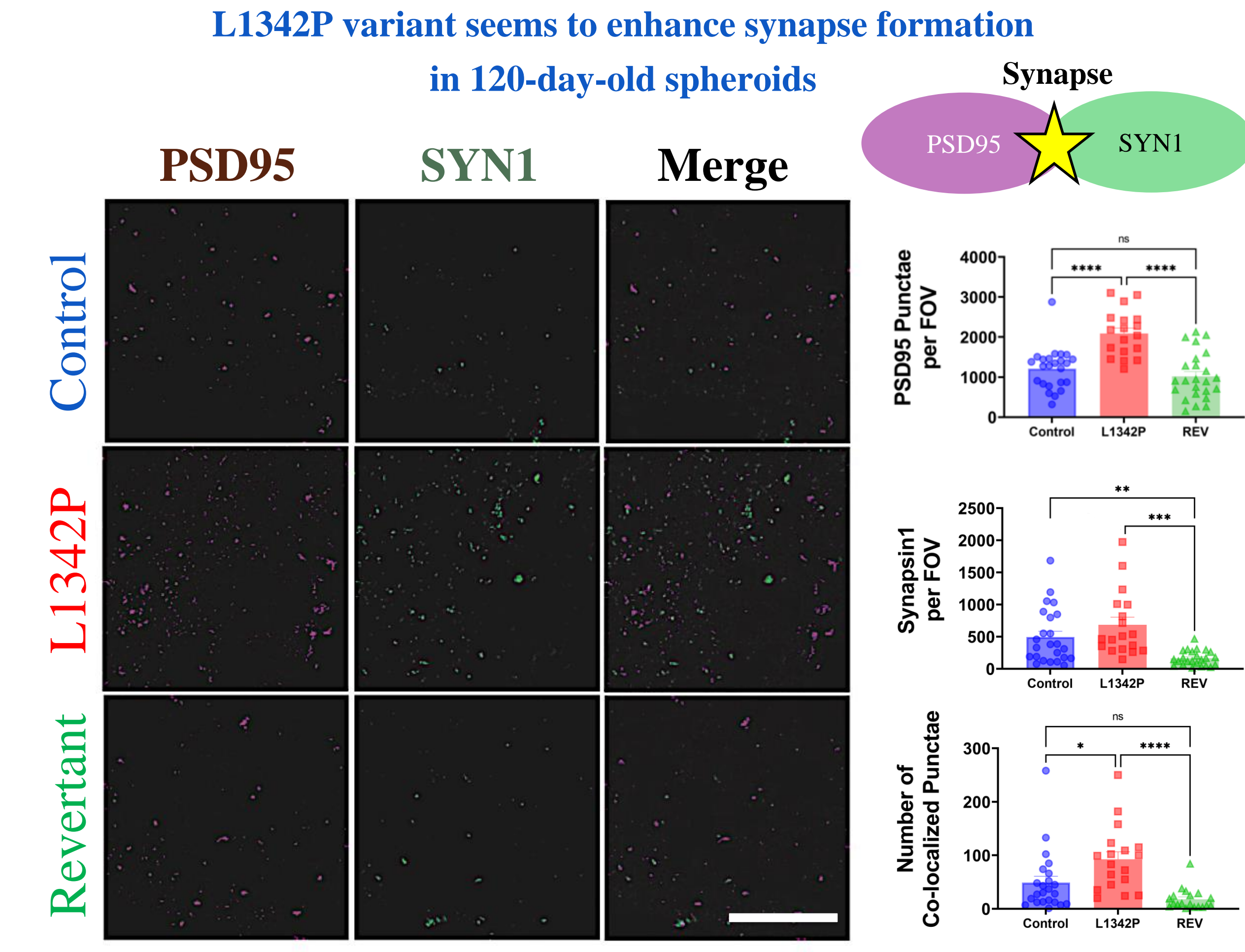


Figure 5. Preliminary data shows the cortical spheroids carrying the L1342P-SCN2A Mutation display increased synapse formation at Day 120. Markers include Postsynaptic density protein-95 (PSD95) in magenta, Synapsin1 (green). Analysis performed using Zen Blue. Magnification 63X. Each dot represents one field of view. At least 2 organoids per genotype. One-Way ANOVA. Scale bar set to 50 µm.

Results

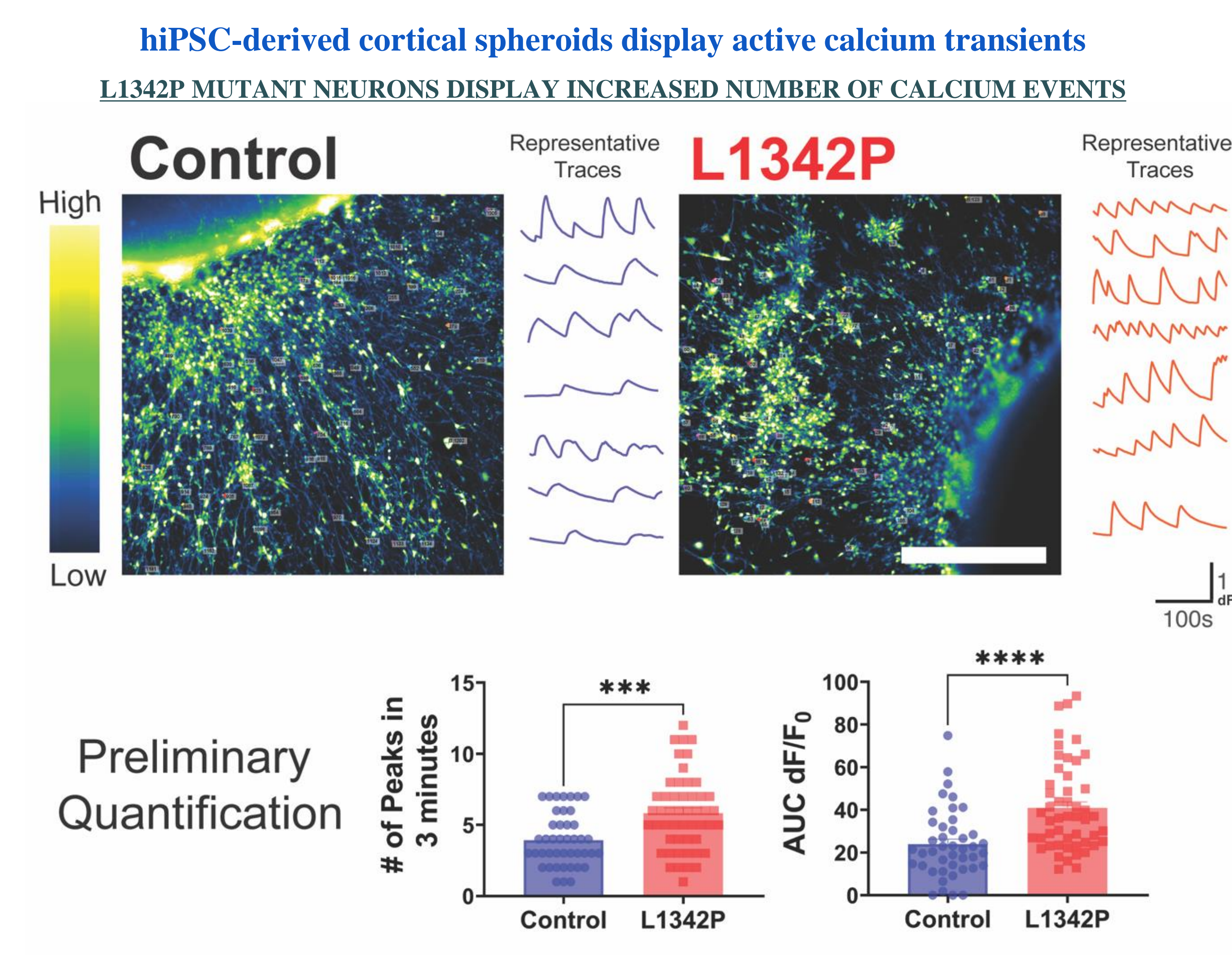


Figure 6. Mature cortical spheroid derived neurons carrying the L1342P-variant display active calcium transients, increased peak frequency and area under the curve. Pseudocolored fields of views containing neurons loaded with Fluo-4. Each dot represents an active neuron. Data are reported as mean ± error (SEM). Scale bar is set to 500 µm. Data analyzed by Student's t test; *p < 0.05.

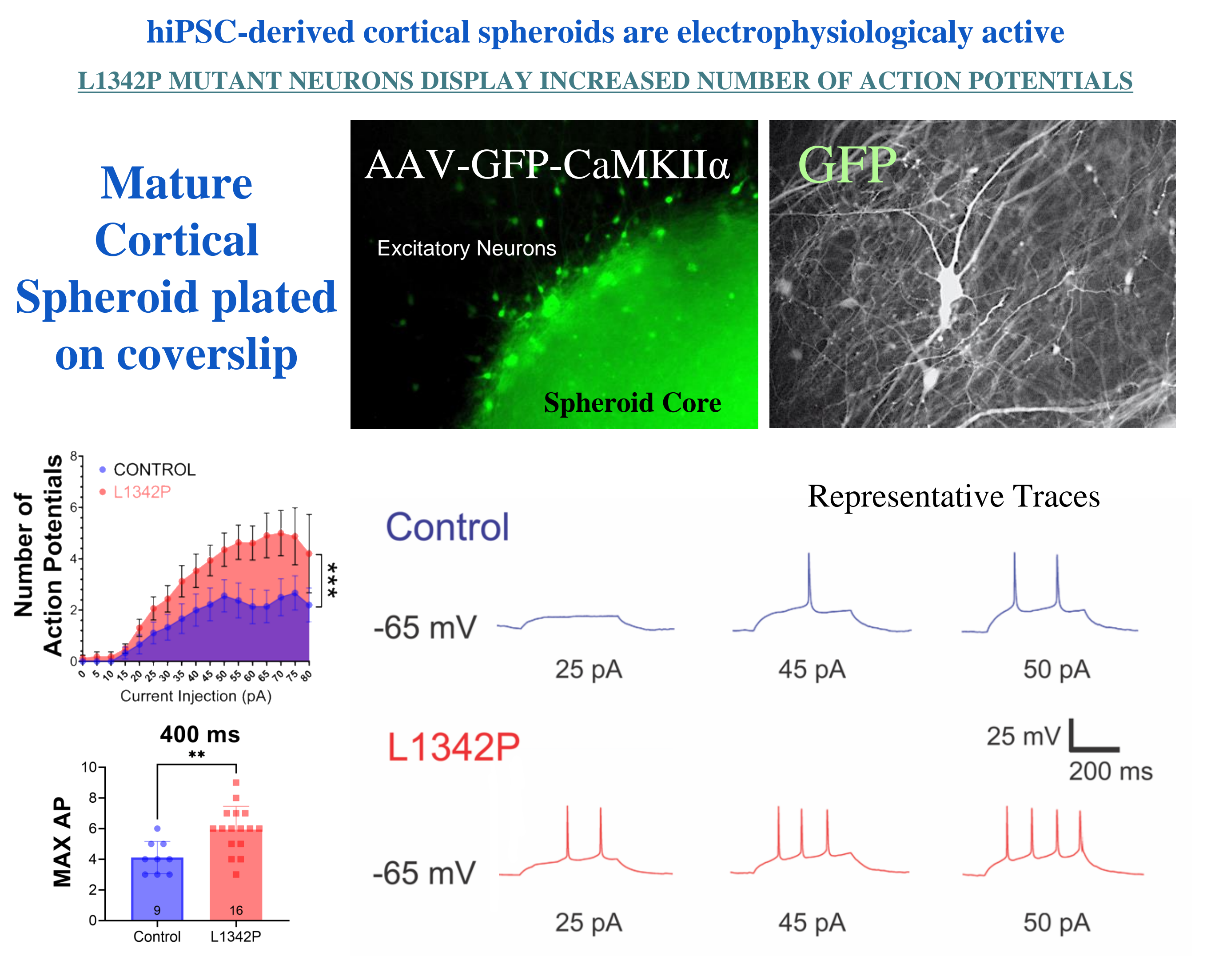


Figure 7. hiPSC-derived neurons with Nav1.2-L1342P variant display increased excitability. The L1342P variant enhances the repetitive firing of hiPSC-derived neurons. Plot showing AP number per epoch in response to graded inputs from 0- to 80-pA current injection (400-ms duration). Representative sustained AP firings from hiPSC-derived Nav1.2-L1342P (red) cortical neurons or isogenic control (blue). Data were collected from two differentiated batches, with two clones used for each genotype. Data analyzed by repeated-measures two-way ANOVA and Student's t test; *p < 0.05.

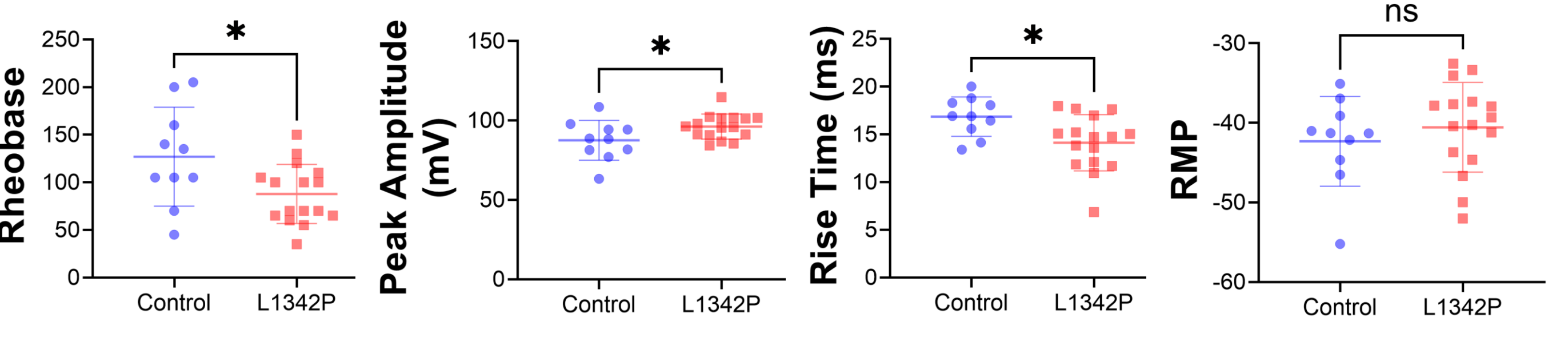


Figure 8. The L1342P variant increases the intrinsic excitability of hiPSC-derived neurons. Data analyzed by Student's t test; *p < 0.05.

Results

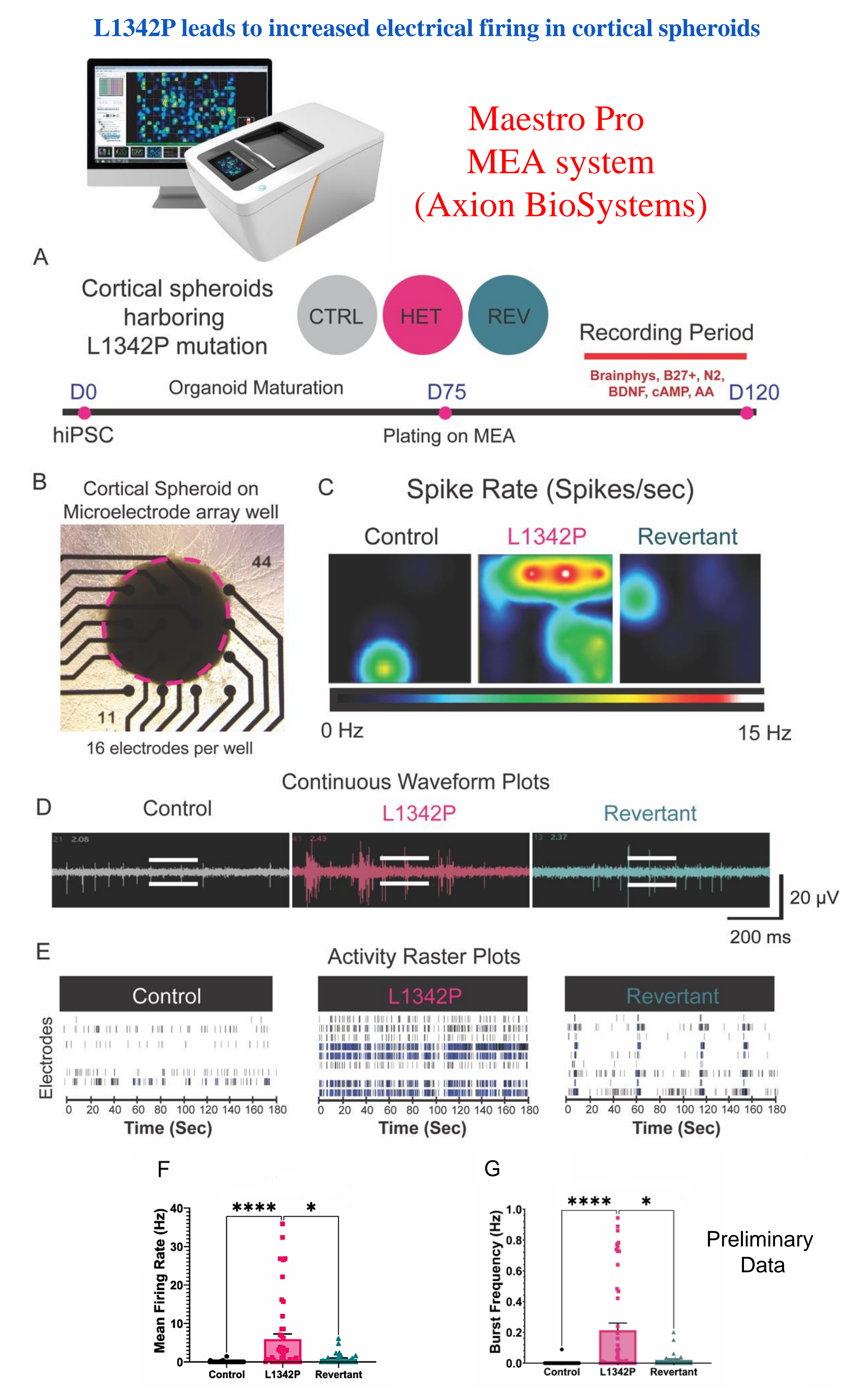


Figure 9. Cortical Spheroids carrying the L1342P Mutation display enhanced network excitability. (A) Experimental Design. (B) Description of organoids used in study and representative image of organoid plated on a 16-electrode MEA well. (C) Activity Heat-maps (D) Representative raw spikes (E) Representative spike raster plots Bursting events are depicted by a cluster of ticks in blue. (F) Mean Firing Rate (G) Burst Frequency. Each dot represents an active electrode. Data are reported as mean ± error (SEM). Data pooled from Control: n = 48, L1342P: n = 48, Revertant: n = 32. Kruskal-Wallis test was performed with *p < 0.05; **p < 0.01; ****p < 0.001

Conclusions

- L1342P Mutation displays characteristic hyperexcitability phenotype.

References

Que, Z., Olivero-Acosta, M. I., Zhang, J., Eaton, M., Tukker, A. M., Chen, X., Wu, J., Xie, J., Xiao, T., Wettschurack, K., Yunis, L., Shafer, J. M., Schaber, J. A., Rochet, J. C., Bowman, A. B., Yuan, C., Huang, Z., Hu, C. D., Trider, D. L., Yang, Y. (2021). Hyperexcitability and Pharmacological Responsiveness of Cortical Neurons Derived from Human iPSCs Carrying Epilepsy-Associated Sodium Channel Nav1.2-L1342P Genetic Variant. *The Journal of Neuroscience : The Official Journal of the Society for Neuroscience*, 41(49), https://doi.org/10.1523/JNEUROSCI.0564-21.2021

Sloan, S. A., Andersen, J., Pasca, A. M., Birey, F., & Pasca, S. P. (2018). Generation and assembly of human brain region-specific three-dimensional cultures. *Nature Protocols*, 13(9), https://doi.org/10.1038/s41596-018-0032-7

Acknowledgements

The authors acknowledge support from the Purdue University Institute for Drug Discovery (PIDD), Purdue Autism Research Center (PARC), and Institute for Integrative Neuroscience. This work was also supported by National Institutes of Health National Institute of Neurological Disorders and Stroke (NINDS) Grants R01 NS117585 and R01 NS123154 (to Y.Y.), NINDS Grant R03 NS108229 (to J.-C.R.), National Institute of Environmental Health Sciences Grant R01 ES031401 (to A.B.B), National Cancer Institute Grant R01 CA212403 (to C.-D.H.), and National Institute of Allergy and Infectious Diseases Grant R01 AI150847 (to D.J.T.). Y.Y. is supported by funds from Ralph W. and Grace M. Showalter Research Trust, and Purdue Big Idea Challenge 2.0 on Autism. M.I.O.A. is supported by Fulbright scholarship program. **The Yang lab thanks the FamiliesSCN2A foundation for the Action Potential Award and Axion for Travel Grant awards.**