

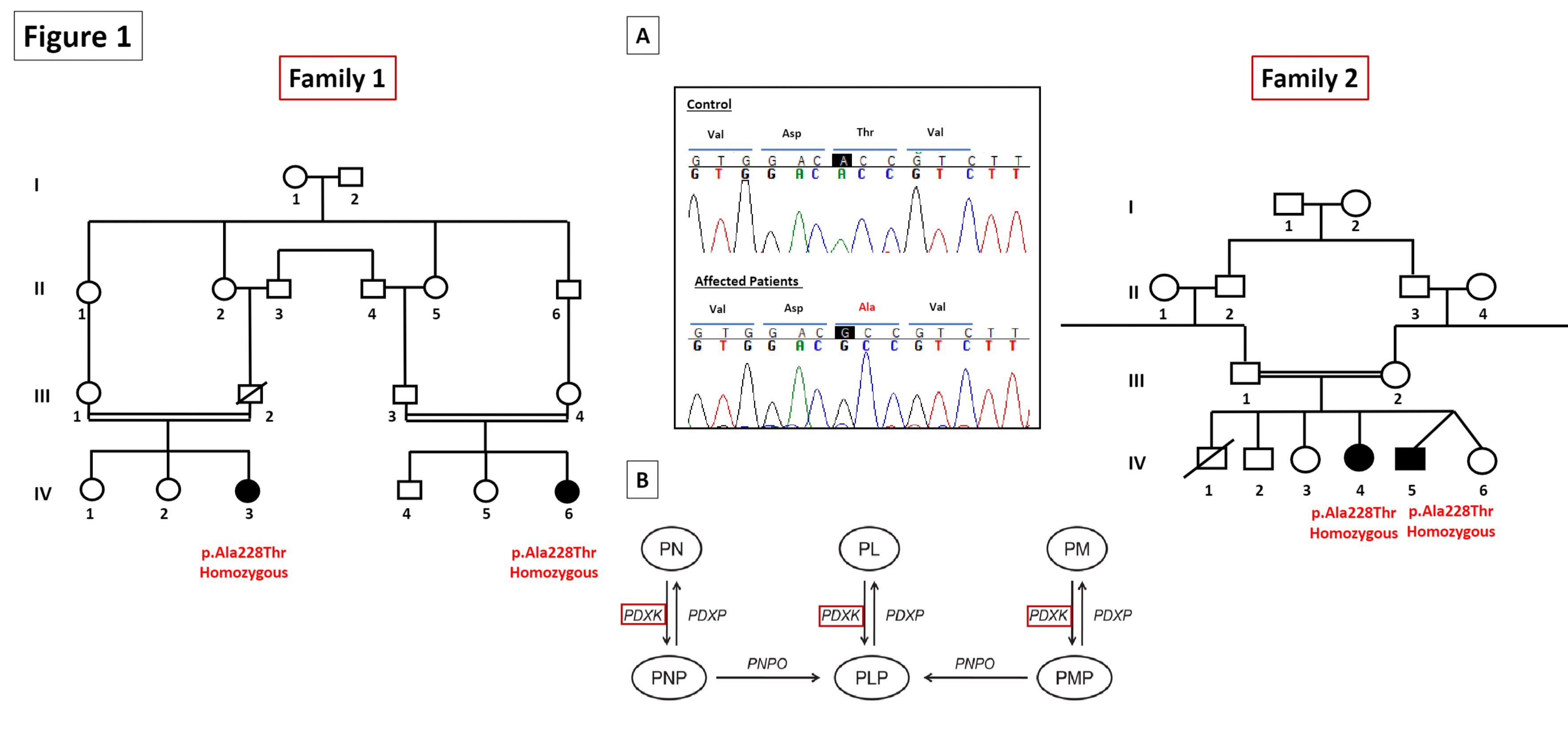
Pathophysiological mechanisms in a new form of Charcot-Marie-Tooth due to a mutation in *PDXK*

Zeinab Hamze¹, Patrice Quintana¹, André Mégarbané², Valérie Delague¹

¹Aix Marseille Univ, INSERM, MMG, U 1251, Marseille, France

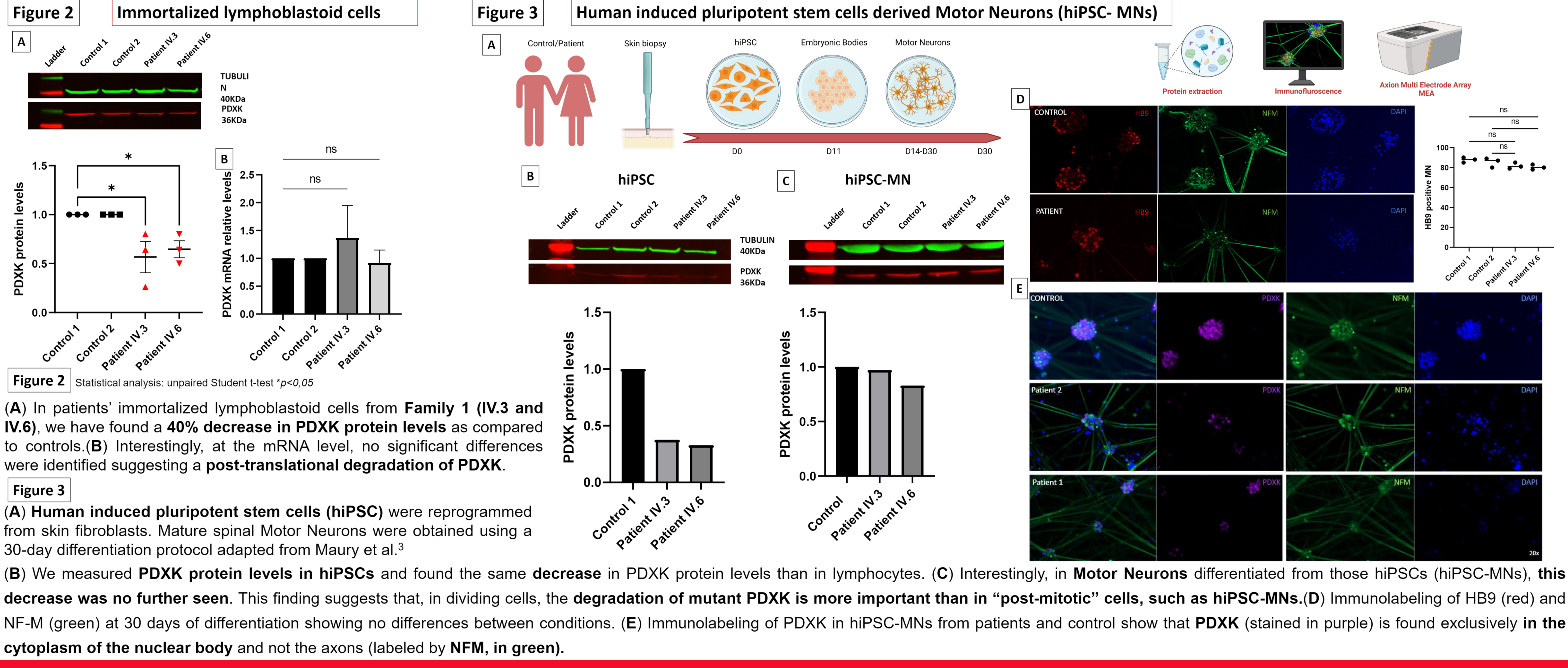
²Department of Human Genetics, Gilbert and Rose-Marie Chagoury School of Medicine, Lebanese American University, Lebanon

GENETIC ANALYSIS



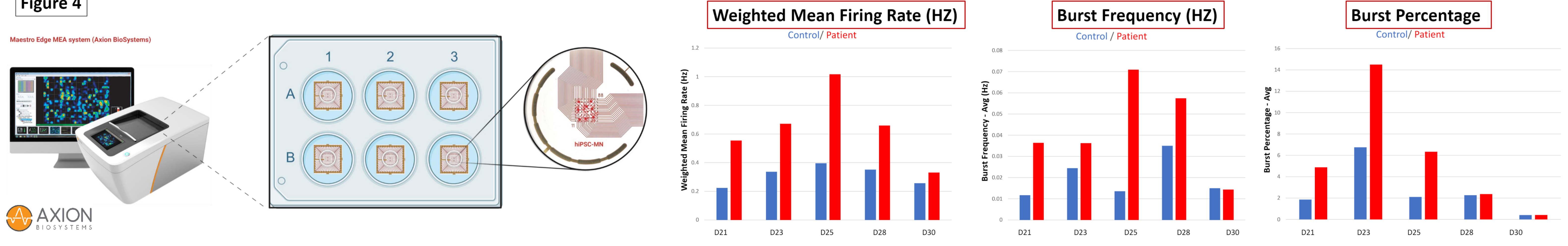
Charcot-Marie-Tooth (CMT) disease is the **commonest inherited group of neuromuscular diseases**. This group of diseases, affecting the **peripheral nervous system**, is characterized by wide clinical and **genetic heterogeneity**, with around 100 genes identified to date¹. Here, we present **4 new patients**, from two **consanguineous** Middle Eastern families, presenting with a rare subtype of CMT, for whom we identified, by Whole Exome Sequencing (WES), a **homozygous missense p.Ala228Thr mutation in the *PDXK* gene (A)**. The same mutation was described previously by Chelban et al 2019, who first identified *PDXK* as responsible for CMT². However, while the patients described by Chelban et al. are affected with axonal CMT associated with visual loss, our patients have **mixed axonal/demyelinating CMT** and no visual impairment. *PDXK* encodes for **Pyridoxal Kinase**, a cytoplasmic protein that converts the inactive B6 vitamers pyridoxine (PN), pyridoxal (PL), and pyridoxamine (PM) into catalytically active PLP (pyridoxal 5-phosphate) (**B**). PLP is an essential cofactor for more than 70 human enzymes implicated in diverse, essential, biological pathways, including amino acid and neurotransmitter metabolism.

FUNCTIONAL ANALYSIS



ELECTRICAL ACTIVITY OF PATIENT'S hiPSC-MNs

DATA ANALYSIS



We have used Maestro Edge MEA system (Axion BioSystems) to measure the global electrical activity of our hiPSC-MNs at different differentiation stages (D21, D23, D25, D28 and D30). This technology allows the recording of the action potentials of hiPSC-MNs plated in 6-well plates for 10 min inside a 37°C chamber at 5% CO₂. We measured three different parameters, i.e. **weighted mean firing rate, burst frequency and burst percentages**, which characterize the general activity of the motor neurons. In this preliminary experiment (n=1), we found an **increase in all three parameters**, in both patients' hiPSC-MNs as compared to controls, at all differentiation time points. The increase seem to peak at differentiation day 25, but **the values come back to normal while MNs gets more mature (D25 to D30)**. These results are preliminary and need to be completed, but they are in line with the differences observed by WB.

CONCLUSION

Here, we describe 4 new patients from 2 unrelated consanguineous families affected with a rare subtype of CMT, harboring a **homozygous p.Ala228Thr mutation in the *PDXK* gene**. Interestingly, our patients present a different phenotype than the ones described earlier with this mutation by Chelban et al, as they have mixed axonal/demyelinating CMT and no visual loss. Measurement of the global general activity using **MEA technology** suggest increased global electrical activity in patients' hiPSC-MNs before they reach maturity. These preliminary results will be completed by the assessment of the role for *PDXK* in myelination, by knock-down of the gene in an *in vitro* myelination model based on the coculture of mouse Dorsal Root Ganglion neurons and Schwann cells. This study is very encouraging toward using **hiPSC-derived motor neurons** to study the pathogenicity of mutations in this specific subtype of CMT, and in CMT in general.

REFERENCES

(1) Pipis, M., Rossor, A. M., Laura, M., & Reilly, M. M. (2019). Next-generation sequencing in Charcot-Marie-Tooth disease: Opportunities and challenges. *Na*
 (2) Chelban, V., Wilson, M. P., Warman Chardon, J., Vandrovcova, J., Zanetti, M. N., Zamba-Papanicolaou, E., Efthymiou, S., Pope, S., Conte, M. R., Abis, G., Liu, Y., Tribollet, E., Haridy, N. A., Botfa, J. A., Rytén, M., Nicolaou, P., Minaidou, A., Christodoulou, K., Kernohan, K. D., ... Tucci, A. (2019). *PDXK* mutations cause polyneuropathy responsive to pyridoxal 5'-phosphate supplementation. *Annals of Neurology*, 86(2), 225-240.
 (3) Maury, Y., Côme, J., Piskrowski, R. A., Salah-Mohellibi, N., Chevalere, V., Peschanski, M., Martinat, C., & Nedelec, S. (2015). Combinatorial analysis of developmental cues efficiently converts human pluripotent stem cells into multiple neuronal subtypes. *Nature Biotechnology*, 33(1), Article 1. <https://doi.org/10.1038/nbt.3049>

