

A role for the primary cilium in network properties of neurons

Emma Dyke¹, Chantal Bijl nagte-Schoenmaker¹, Rachel Mijdam³, Lisa Rahm⁴, Marie Le Bihan^{1,2}, Ummi Ciptasari^{1,2}, Jindřiška Leischner Fialová⁵, Brooke Latour¹, Dirk Schubert⁶, Ronald Roepman¹, Nael Nadif Kasri^{1,2}

1) Department of Human Genetics, Radboudumc, Nijmegen, The Netherlands; 2) Donders Institute for Brain, Cognition and Behaviour, Nijmegen, The Netherlands; 3) Radboudumc, Neurology, Nijmegen, The Netherlands; 4) Radboudumc, Donders Institute for Brain, Cognition, and Behaviour, Human Genetics, Nijmegen, The Netherlands; 5) Department of Biology, University of Copenhagen, Copenhagen, Denmark; 6) Donders Institute for Brain, Cognition and Behavior, Department of Cognitive Neuroscience, Nijmegen, The Netherlands.



Radboudumc



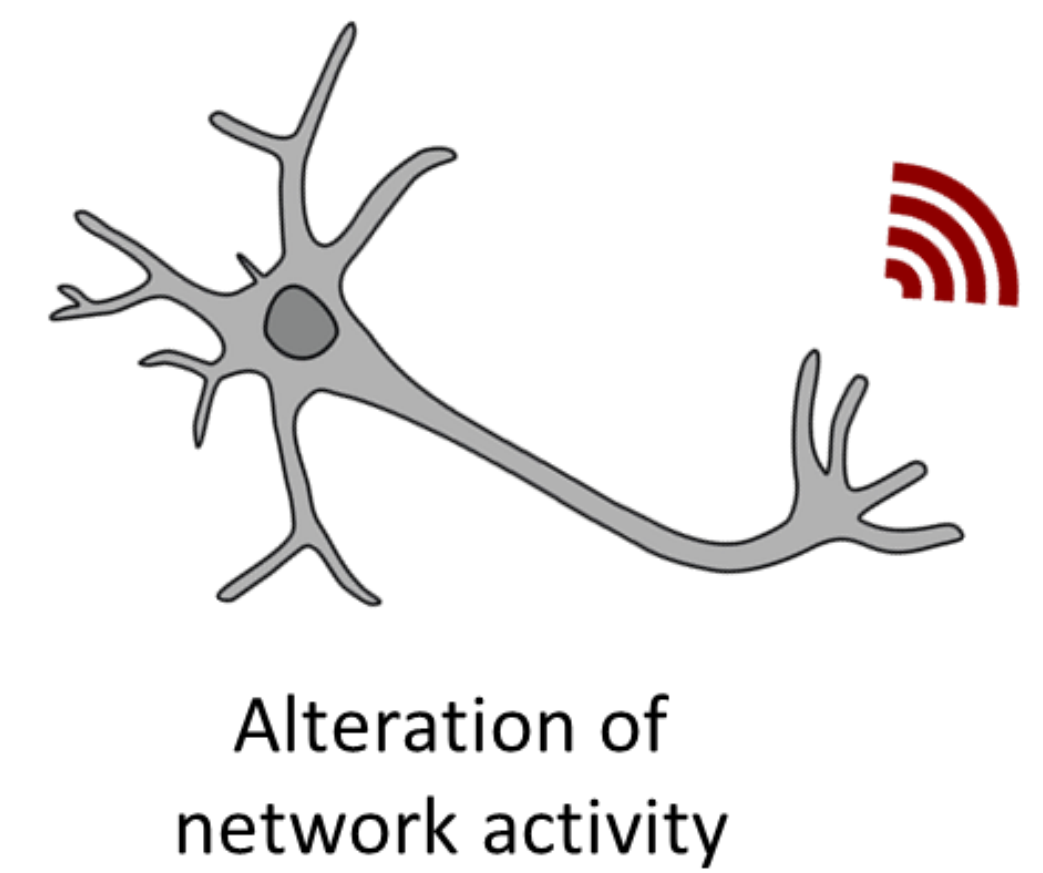
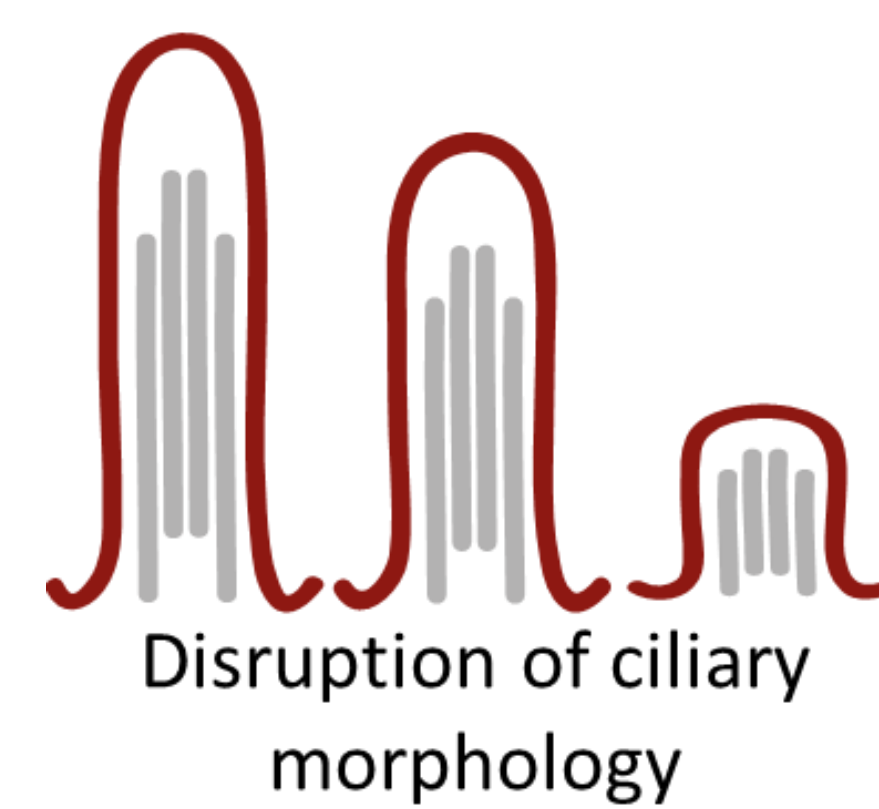
Introduction

The **primary cilium** – often referred to as the cells' antenna - is a small organelle involved in a diverse array of signalling pathways. Dysfunction of the cilium results in a spectrum of developmental disorders including the neurodevelopmental disorder, **Joubert syndrome**. Up to 90% of neurons in the mammalian cerebral cortex are ciliated, yet the function of this organelle in excitatory neurotransmission has not been fully established.

We seek to:

- Further characterise the role of the primary cilium in establishing a neuronal network,
- Generate a model in which to study the role *Joubert syndrome*-related proteins play in network formation and maturation.

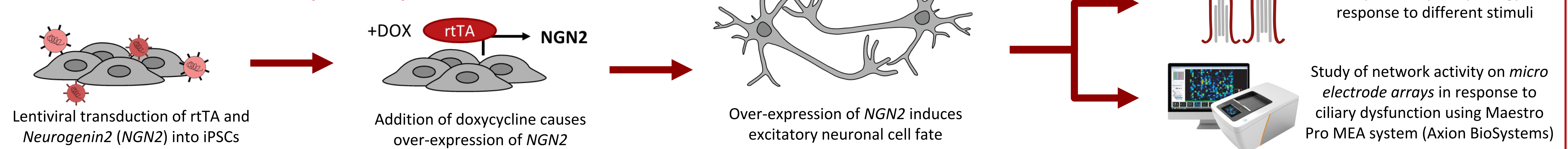
Discussion and concluding remarks



The primary cilium is linked to the excitatory properties of *iNeurons* with morphological changes preceding and following abnormal network activity.

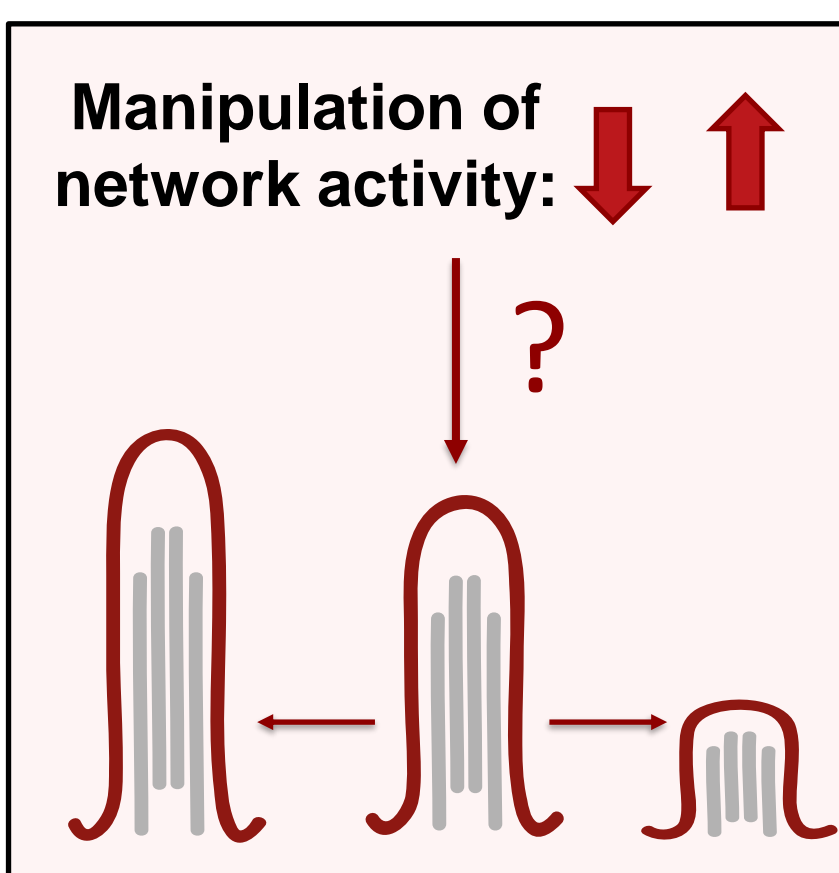
Model system

Generation of inducible neurons (*iNeurons*):



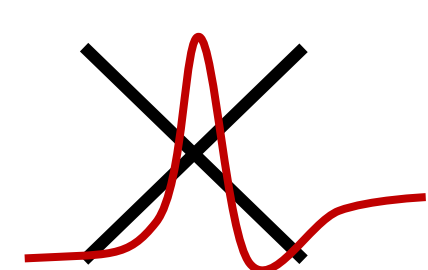
Cilia respond to changes in network properties

The primary cilia of *iNeurons* show bidirectional changes in volume in response to decrease or increase in activity:

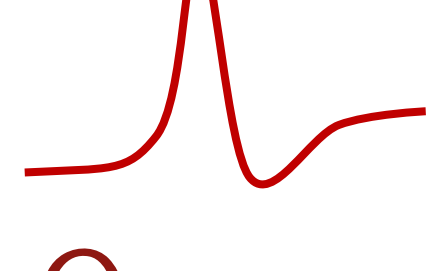


Recent work has shown that knockdown of ciliary protein expression alters the intrinsic excitability of cultured neurons. We sought to investigate if manipulation of network state could affect ciliary morphology.

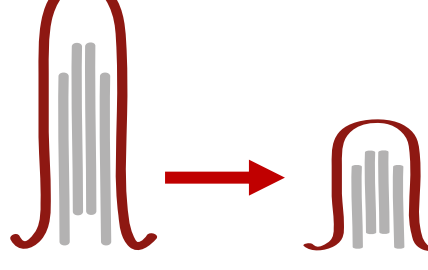
Tetrodotoxin (TTX):



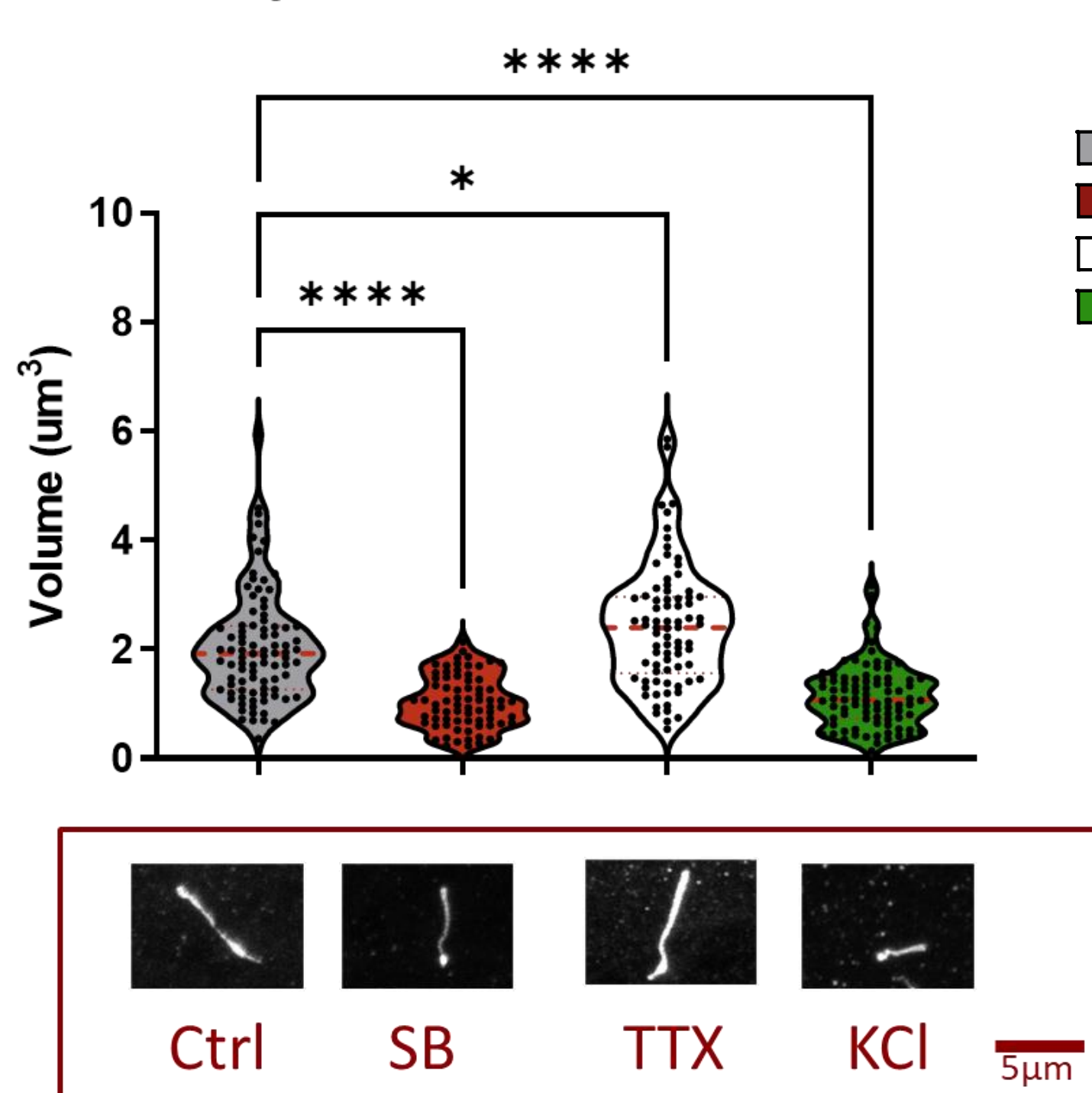
Potassium chloride (KCl):



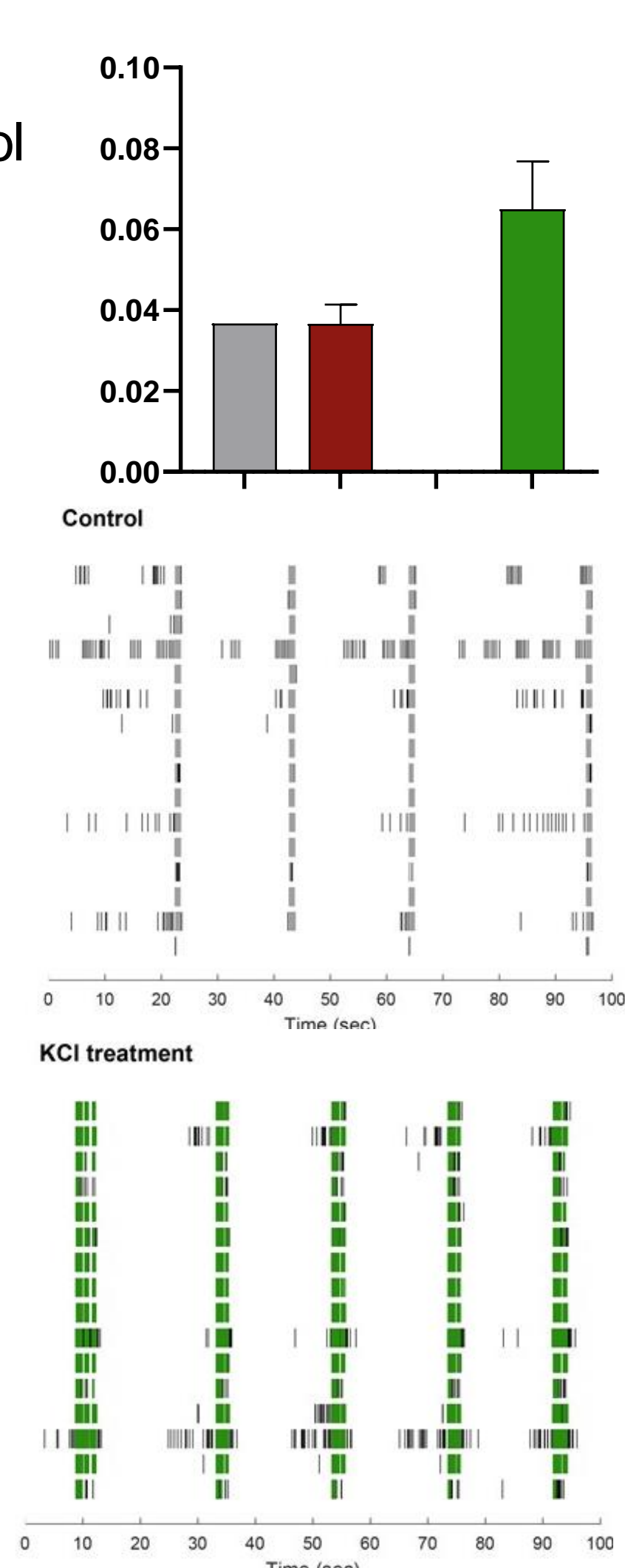
SB399885 (SB):



Ciliary volume of treated *iNeurons*



Network burst frequency

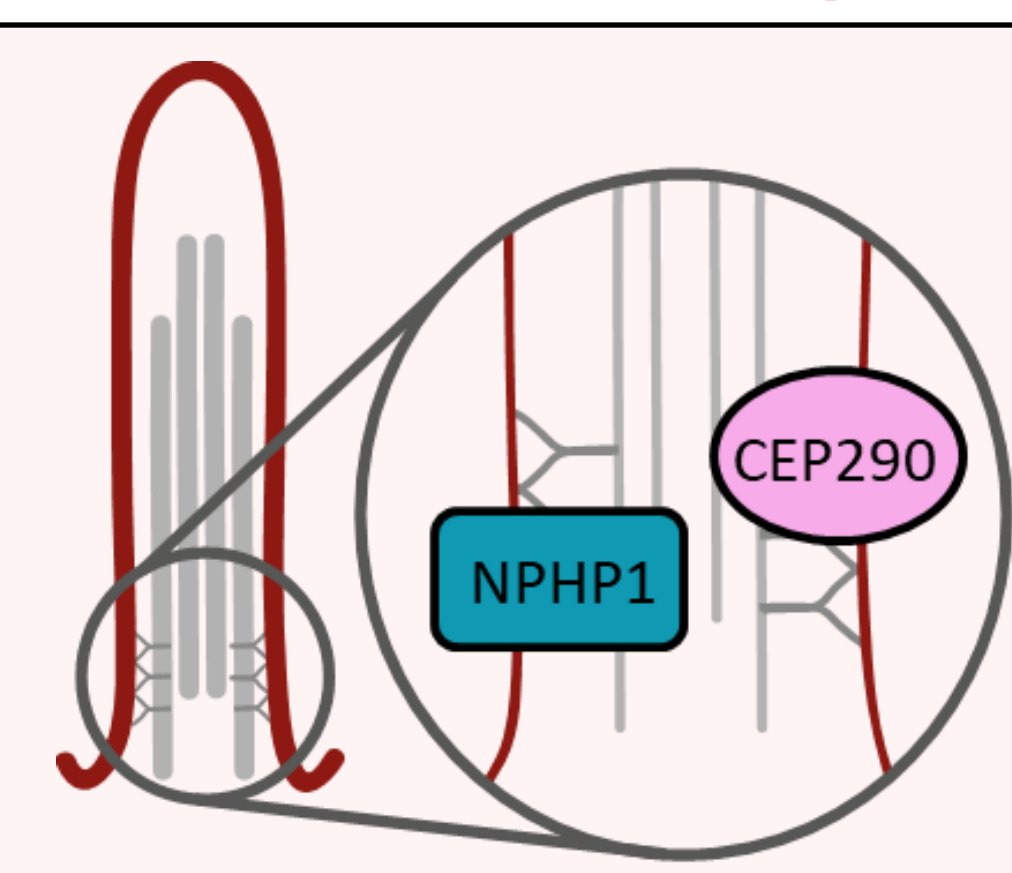


Mean inter burst interval

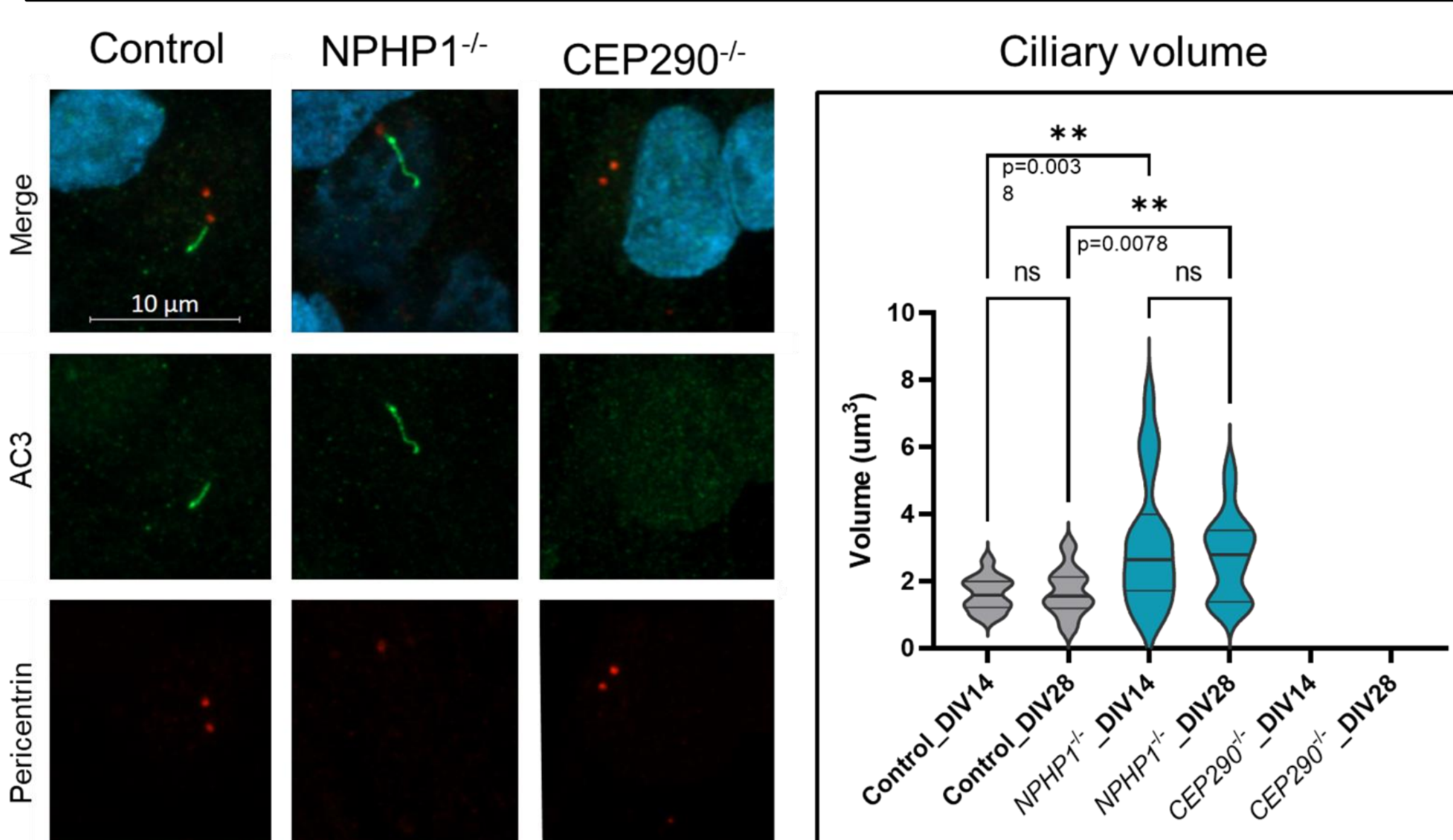


Alteration of ciliation changes activity

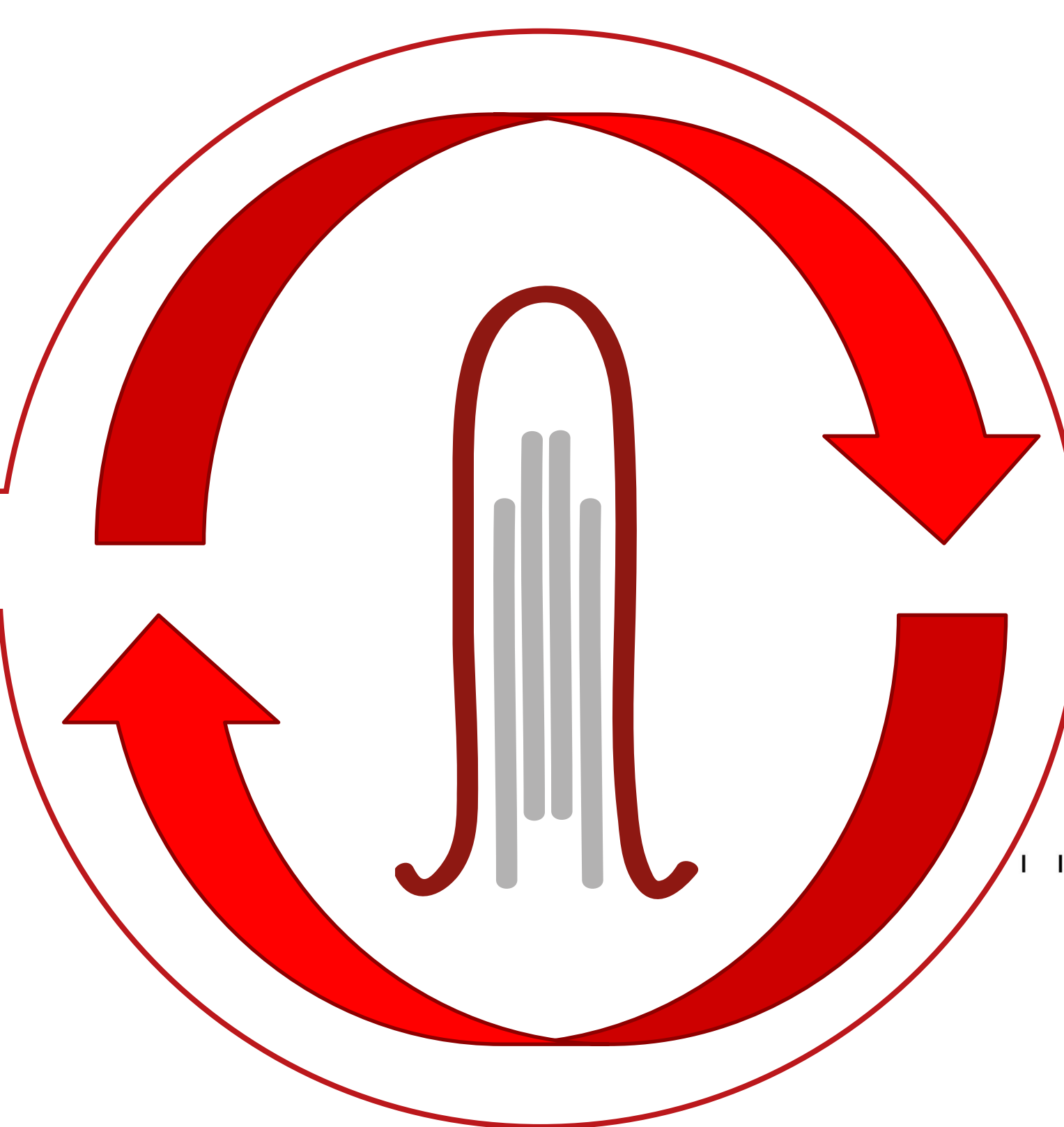
Loss of transition zone proteins alters ciliary morphology:



NPHP1 and CEP290 form components of important complexes localised to the transition zone. This part of the cilium is essential for regulating ciliary composition.

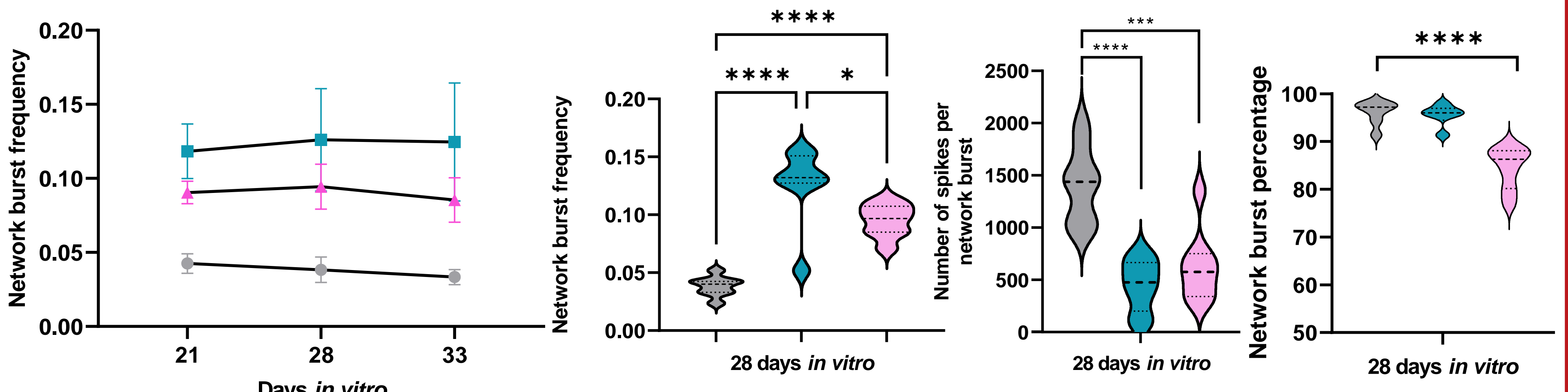
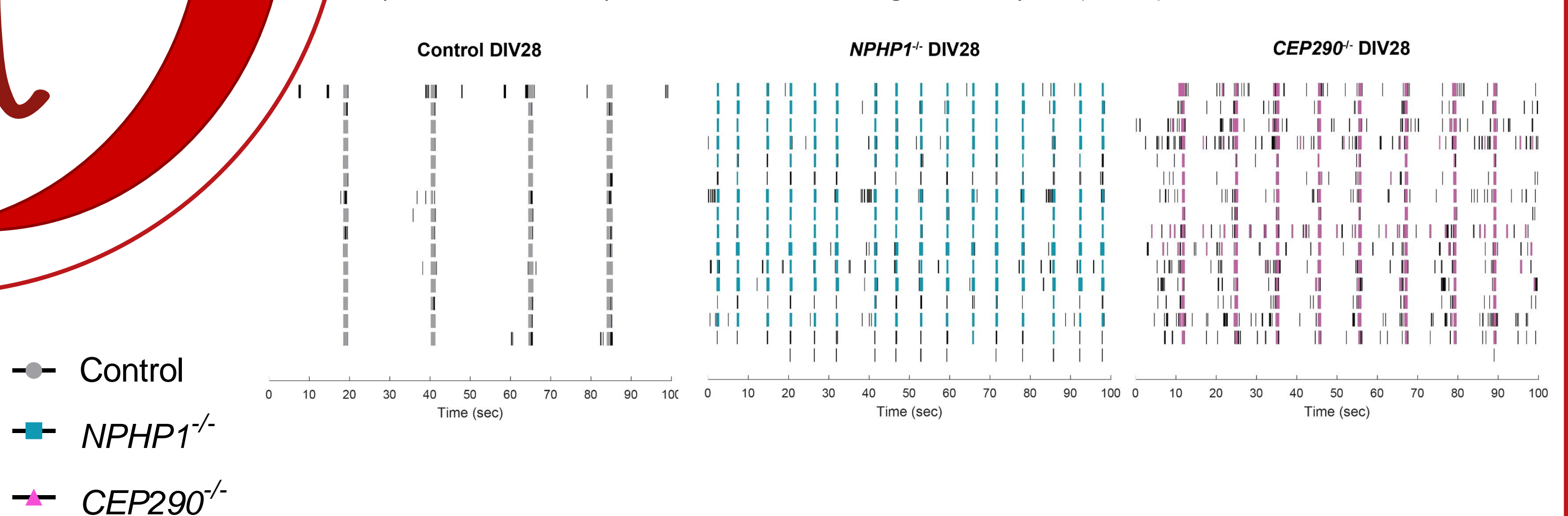


Immunostaining of *iNeurons* derived from *NPHP1*^{-/-} and *CEP290*^{-/-} iPSCs



NPHP1^{-/-} and *CEP290*^{-/-} derived *iNeurons* show increased network activity:

Representative rasterplots from MEA recordings of 28 day old (DIV28) *iNeurons*:



Loss of transition zone proteins CEP290 and NPHP1 leads to loss of cilia, and increase in ciliary volume respectively. Despite the opposite morphologies the cell lines display similar network dynamics indicating that loss of transition zone proteins alters the basal network activity of *iNeurons*.