

Functional biomarker of rare and familial diseases: NPC1 knockout neuronal networks phenotyped with HTS micro electrode arrays and artificial intelligence machine learning methods

Introduction

Background: There is an urgent need to develop fast and reliable disease models in rare diseases which can be used for screening in a standardized manner. One strategy is the development of more predictive pre-clinical in vitro models for gene-associated diseases such as Niemann-Pick Disease Typ C1. Niemann-Pick Type C disease (NPC) is an autosomal recessive neurodegenerative disease caused by a mutation in either the NPC1 gene (in 95% of cases) or the NPC2 gene (in 5%). The pathological changes in NPC1 are characterized by the excessive storage of unesterified cholesterol in lysosomes. NPC Patients show increasing loss of motor control, seizures and other neuropathological symptoms.

Objectives: Using cell cultures from NPC1 knockout mice, we aimed to identify a correlation between the disease-associated genotype and its functional *in vitro* phenotype.

Results

Structural/morphological effects of NPC1 knockout

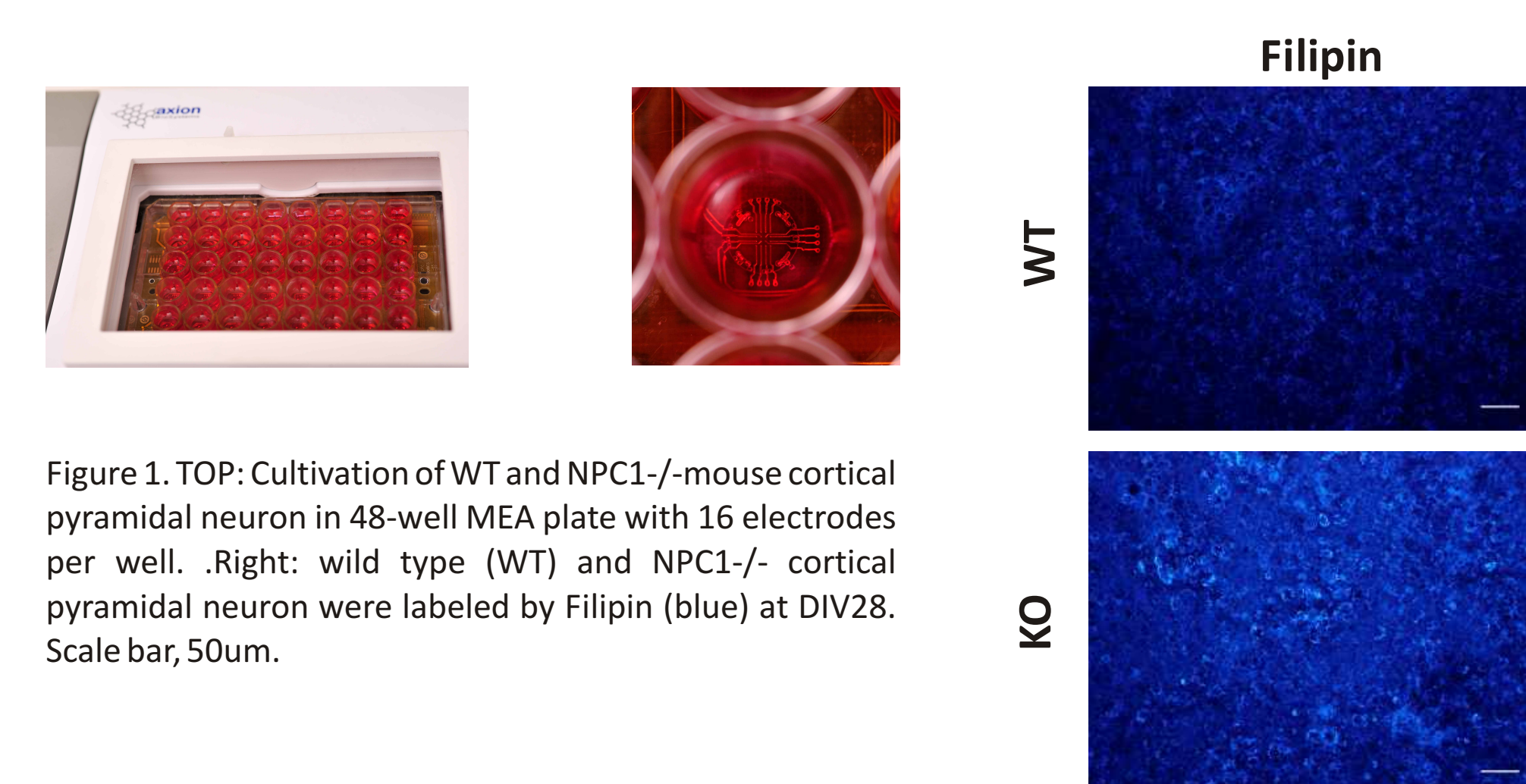


Figure 1. TOP: Cultivation of WT and NPC1-/- mouse cortical pyramidal neuron in 48-well MEA plate with 16 electrodes per well. Right: wild type (WT) and NPC1-/- cortical pyramidal neuron were labeled by Filipin (blue) at DIV28. Scale bar, 50um.

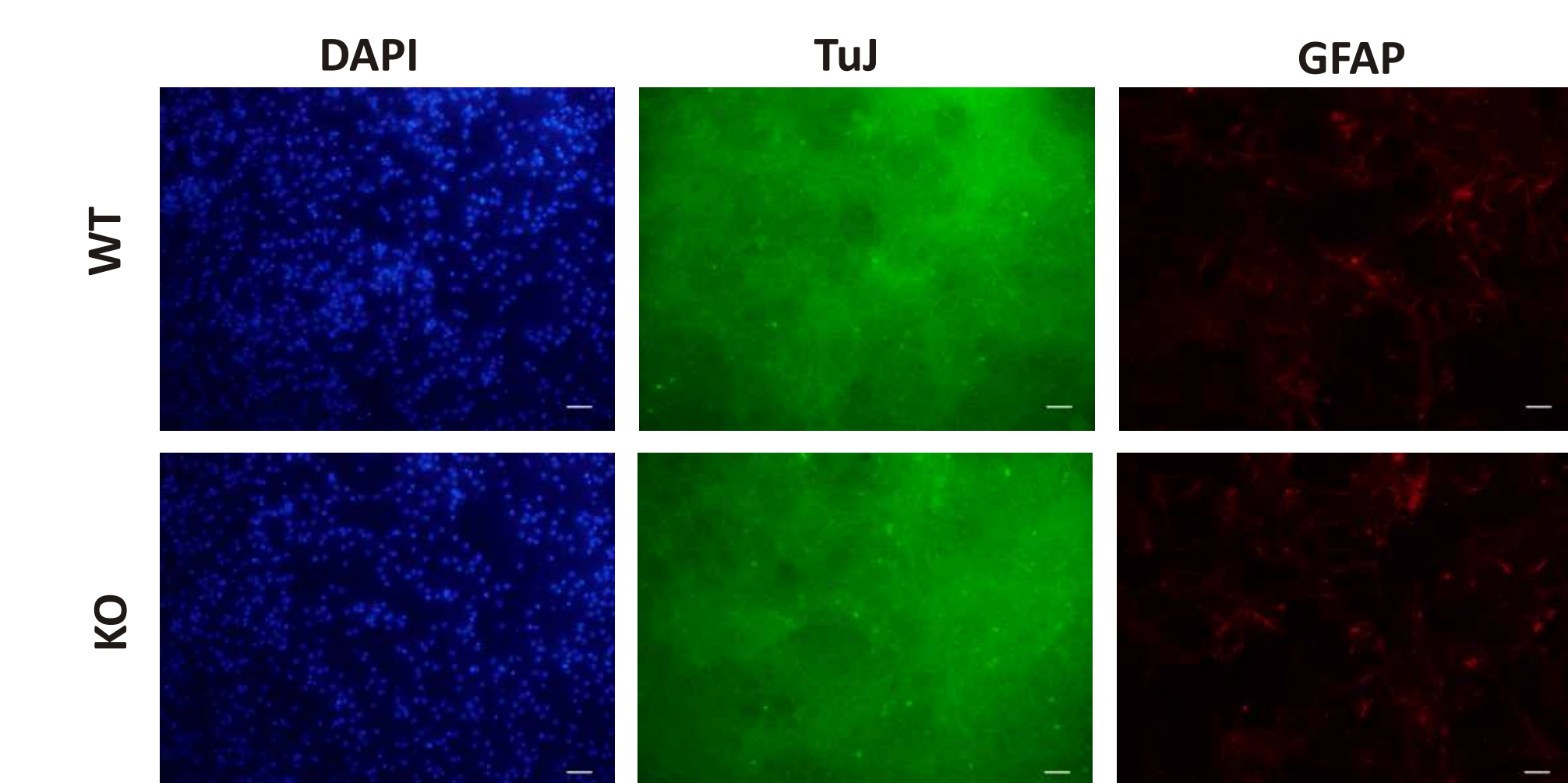


Figure 2: Immunostaining of WT and NPC1-/- cortical pyramidal neurons at 28 days in vitro (DIV). Tuj1 neurons (green), GFAP Glial cells (red), DAPI for nuclei (blue). Scale bar, 50um. Except for the differences in filipin staining (Figure 1), no general morphological differences are observed.

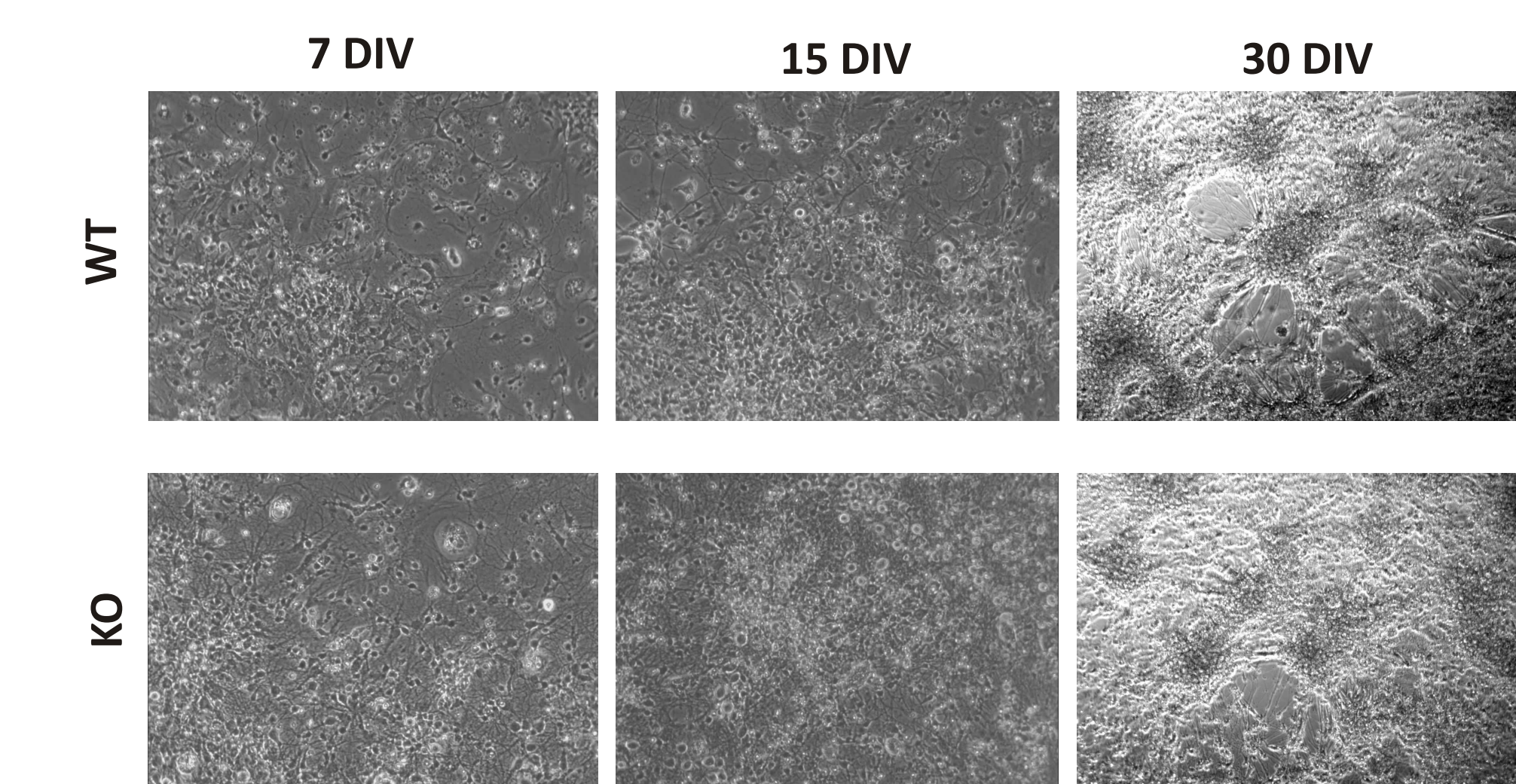


Figure 3: Phase contrast images of neuronal networks between 7 and 30 DIV. Cells growing under same conditions as in 48 well MEAs (coating, density). No general morphological differences are observed.

Summary and Conclusion

We demonstrate that NPC1 knockout cell culture exhibit a significantly different functional phenotype shown by different activity levels and neuronal communication. Moreover, the functional development into mature neuronal networks is affected by the knockout of the NPC1 gene. This specific phenotype can partially be rescued by compound addition.

In conclusion, we present a means to functionally phenotype gene-associated disease in vitro models by multi-variate and machine learning methods and thus provide functional biomarkers for screening and rank small molecules or natural compounds based on their rescue efficacy of disease genotype-associated functional phenotypes towards isogenic controls or wild type phenotypes.

Methods

Primary culture: P1 mice were genotyped and *npc1* knockout and homozygous litters selected for cortical pyramidal neurons isolation at P2 according to the protocol by Brewer & Torricelli 2007⁽¹⁾. **Genotyping:** PCR with primer sets for knockout: 5'-GGTGTGGACAGCAAGTA-3' and 5'-TGAGCCCAAGCATAACTT-3' and for wildtype: 5'-TCTCACAGCCACAAGCTTC-3' and 5'-CTGTAGCTCATCTGCCATCG-3'. **MEA culture:** 48-well MEAs were coated with PEI/Laminin. Recording was performed with Maestro MEA System (Axion BioSystems). **Immunocytochemistry:** cells were PFA-fixed, permeabilized and incubated overnight with mouse anti-βIII tubulin, 1:1000 (Santa Cruz) and GFAP, 1:500 (Agilent) and goat anti-mouse Alexa 488, goat anti-Rabbit Cy3, 1:500. Filipin 0.5 mg/ml (Polysciences). **Data analysis:** multi-parametric data analysis of 204 spike train parameters was performed using NPWaveX software (NeuroProof). "Effect Score": Z' factor-based projection of multiple parameters into a single layer parameter.

NPC1 knockout affects functional phenotype

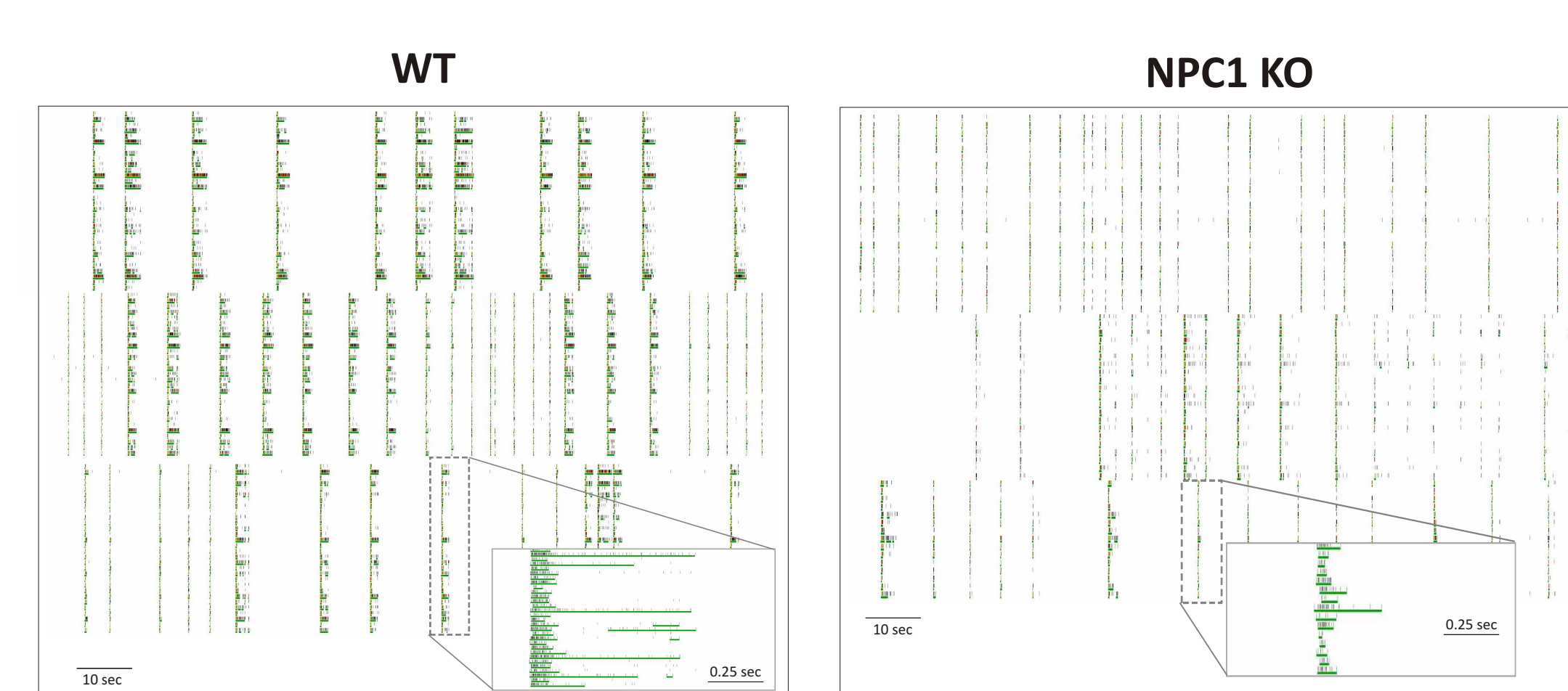
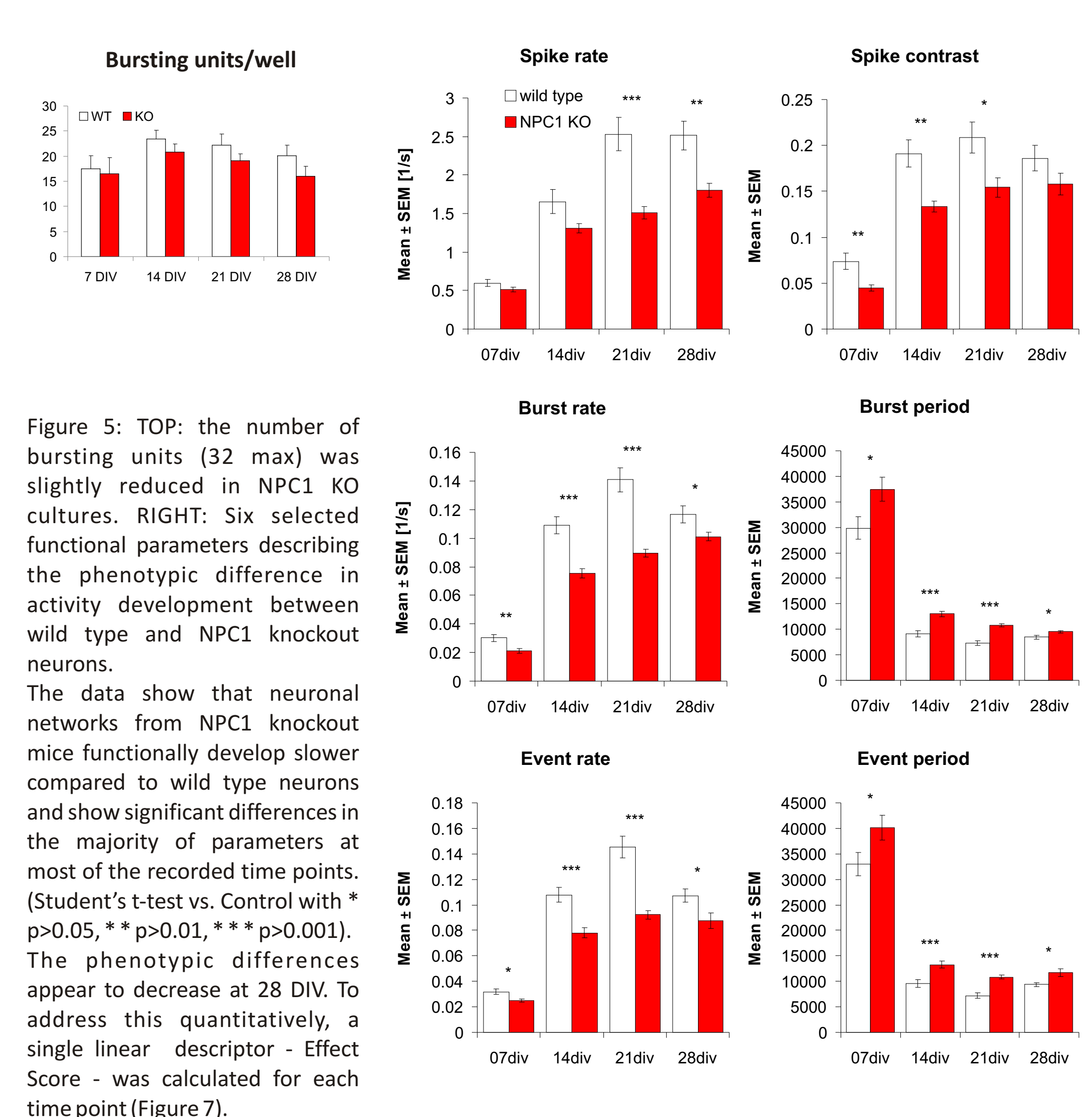


Figure 4: Three example Spike trains for NPC1 knockout and wild type at 21 DIV show a strong phenotypic difference in activity patterns. Spiking and bursting is reduced accompanied by weaker bursts (insert image). Green bars represent detected bursts. Two unit separation was performed.

NPC1 knockout delays functional development



Compound-mediated rescue of NPC1 knockout phenotype

Multi-parametric data complicate the ranking for identification of optimal assay time points. Our NeuroProof Effect Score allows ranking of phenotypic differences which are multi-parametrically described based on a single readout: a generic analysis of multiple parameters out of the 204 parameters selected by an optimization algorithm, which enables an objective, clear and easy-to-compare documentation of the phenotypic relationship of the complete phenotypic data.

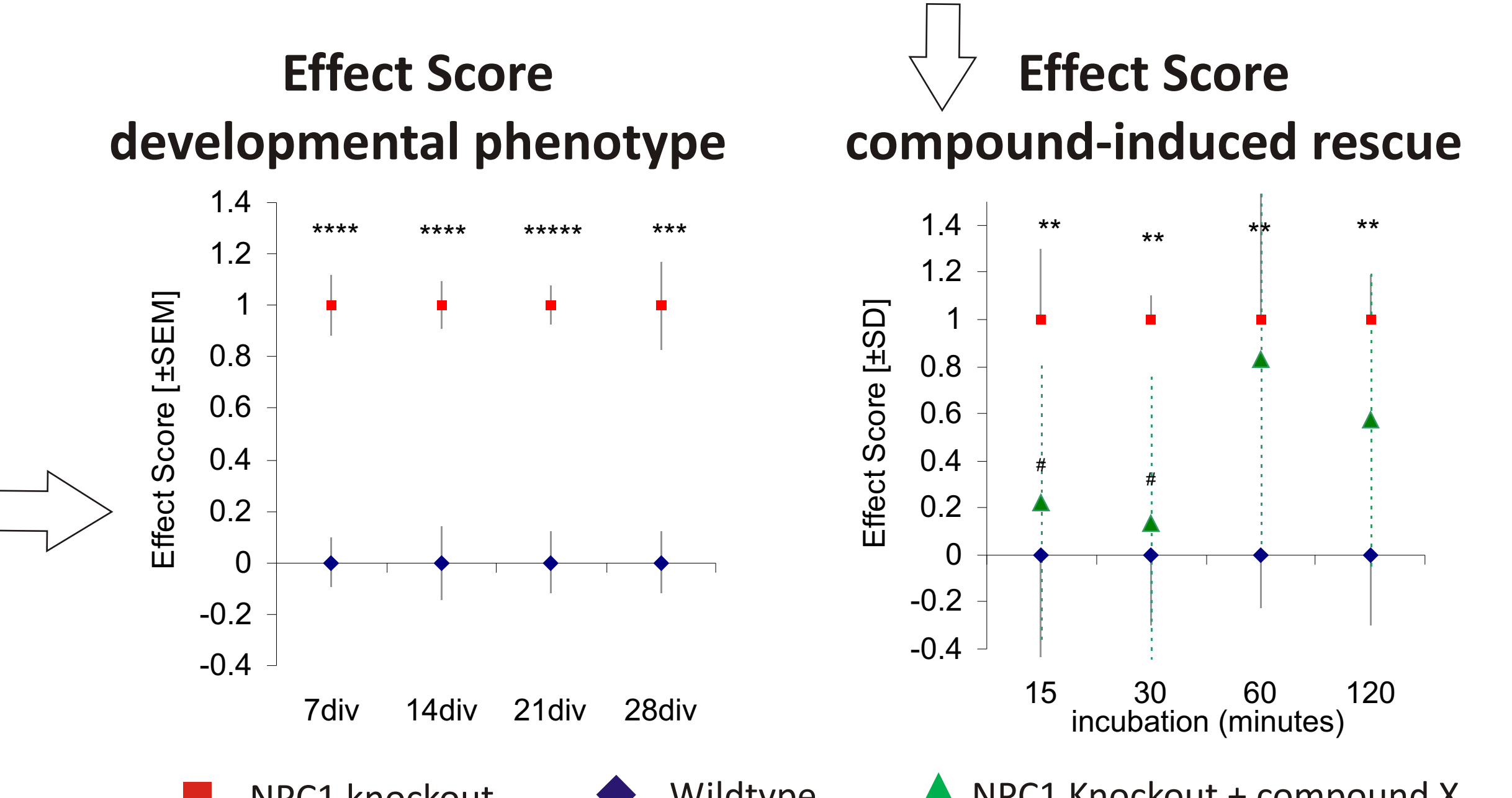
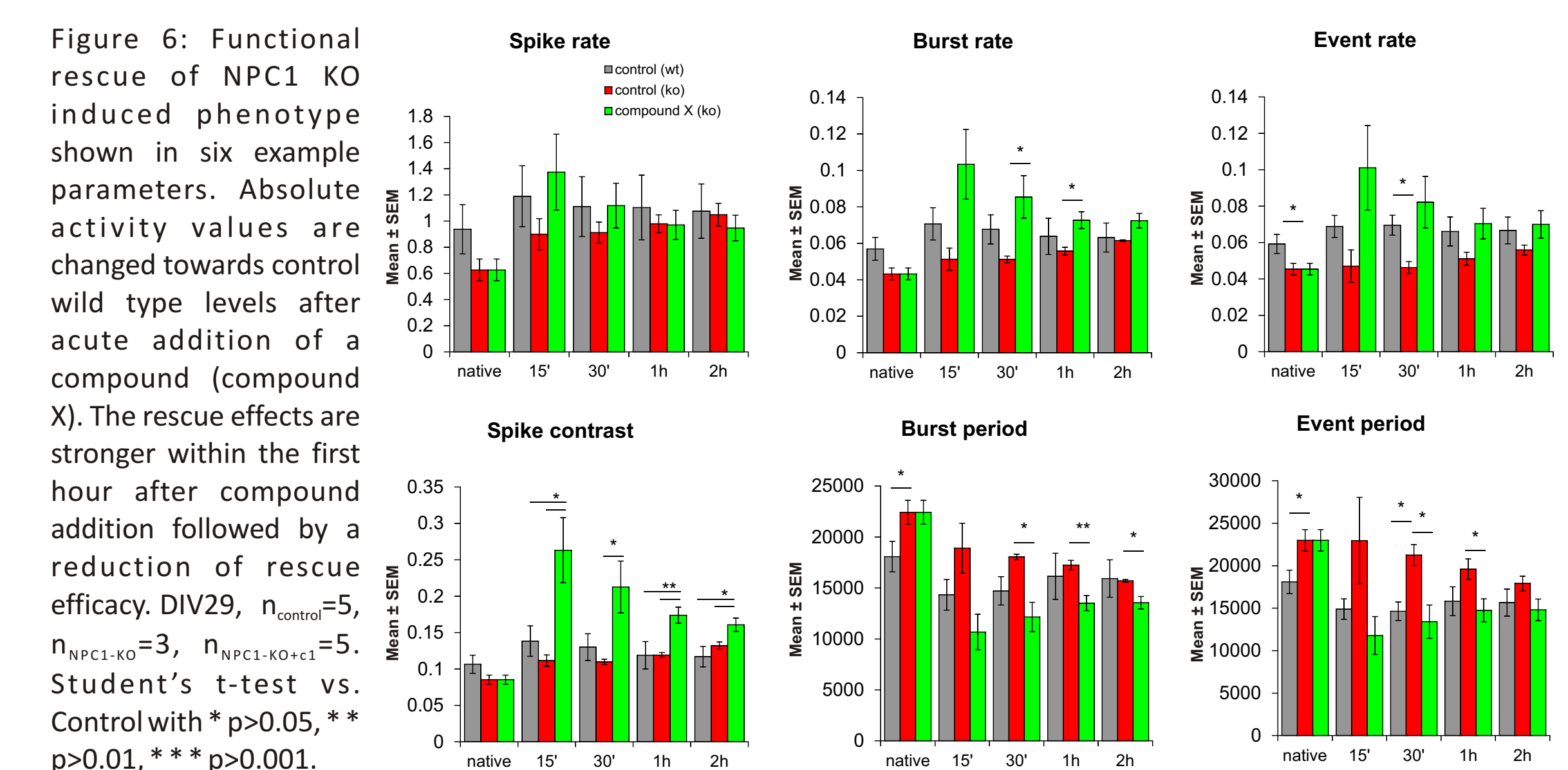


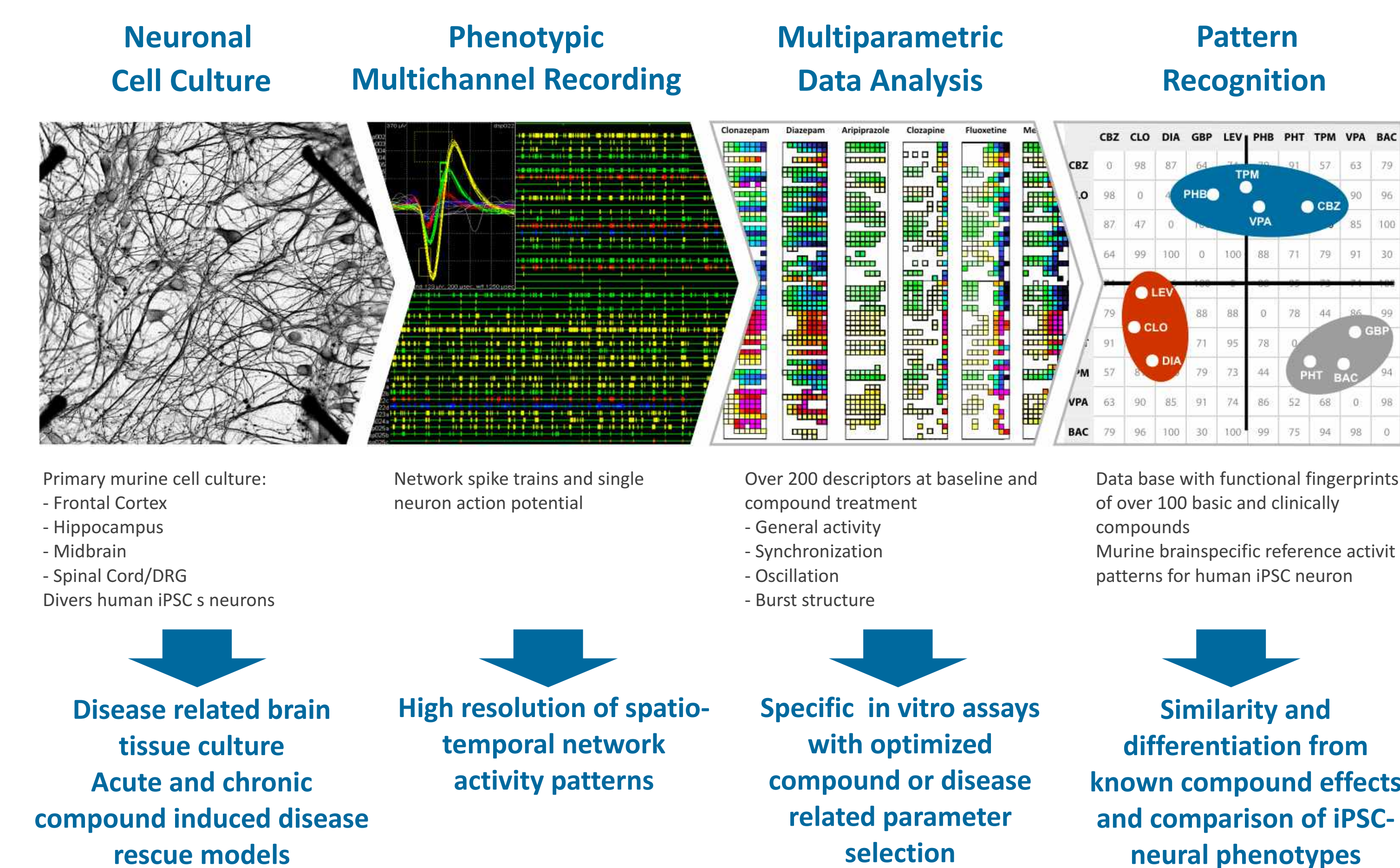
Figure 7: "Effect Score" calculation: Projection of up to 204 parameters into a single parameter allows projecting the complete functional fingerprint into one dimension and thus, LEFT: identifying the optimal time point showing the highest phenotypic difference (here 21 DIV). Wild type is set to "0", NPC1 knockout is set to "1", $n_{control}=14|14|15|12$, $n_{NPC1-KO}=12|17|15|12$. RIGHT: the evaluation of test compounds which rescue the disease-related phenotype towards wild-type conditions. Wild type is set to "0". NPC1 knockout is set to "1". Acute addition of compound X results in a partial rescue of KO-induced functional phenotype towards wild type conditions. Number of experiment: $n_{control}=5$, $n_{NPC1-KO}=3$, $n_{NPC1-KO+X}=5$. Student's t-test with vs. Control, # vs. NPC1 KO with $p>0.05$, $p>0.001$, $p>0.0001$, $p>0.00001$ (number of * (#)). The calculated Z-factor (describing the effect size) is optimized to find the best discrimination between WT and NPC1 KO. The "effect score" describes the rescue. For more information on z'-factor, see Kuemmel et al. 2010⁽²⁾ and Kozak et al. 2010⁽³⁾.

Xiao Feng¹, Monica Segura-Castell², Luise Schultz², Konstantin Jügel², Olaf H.-U. Schröder² Jiankai Luo¹, Benjamin M. Bader²

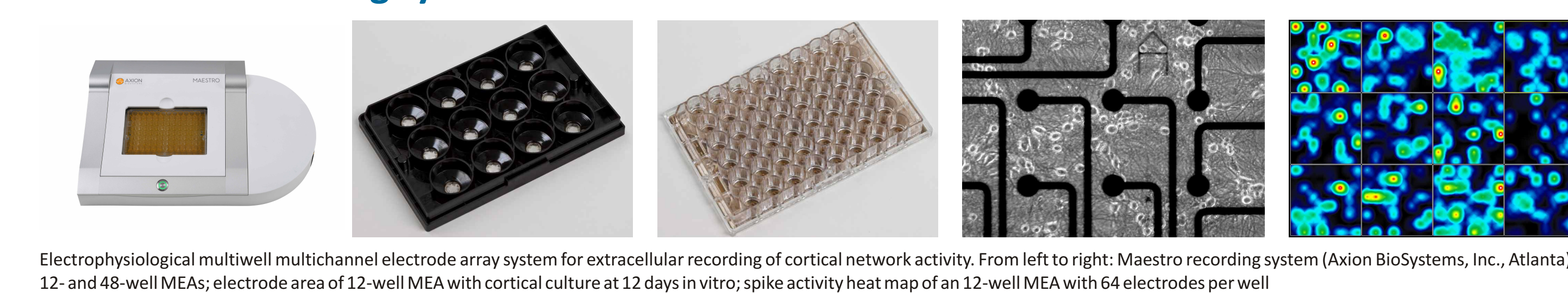
1 University Medicine Rostock, Albrecht-Kossel-Institut for Neuroregeneration, Rostock, Germany; 2 NeuroProof GmbH, Rostock, Germany.

NeuroProof Technology

Phenotypic Screening with MEA-Neurochips



MAESTRO Recording System



Multiparametric Characterization of Neuronal Network Activity

Read out:

- Extracellular action potentials on a single neuron and network activity level
- Spatio-temporal activity changes as well as synchronicity and oscillation in time scales of spikes and bursts

Each specific spike train is described by 200 parameters in 4 categories:

