

MYT1L haploinsufficiency in human neurons and mice causes reversible autism-associated phenotypes

Bettina Weigel^{1,2,3}, Jana Tegethoff^{1,2,3}, Bryce Lim^{1,2,3}, Sarah Grieder^{1,2,3}, Bhuvaneswari Nagarajan^{1,2,3}, Christian Arnold⁴, Judith B. Zaugg⁴, Moritz Mall^{1,2,3}

¹⁾ Cell Fate Engineering and Disease Modeling Group, German Cancer Research Center (DKFZ) and DKFZ-ZMBH Alliance, 69120 Heidelberg, Germany

²⁾ HITBR Hector Institute for Translational Brain Research gGmbH, 69120 Heidelberg, Germany

³⁾ Central Institute of Mental Health, Medical Faculty Mannheim, Heidelberg University, 68159 Mannheim, Germany

⁴⁾ European Molecular Biology Laboratory, Structural and Computational Biology Unit, 69115 Heidelberg, Germany.

Research for a Life without Cancer

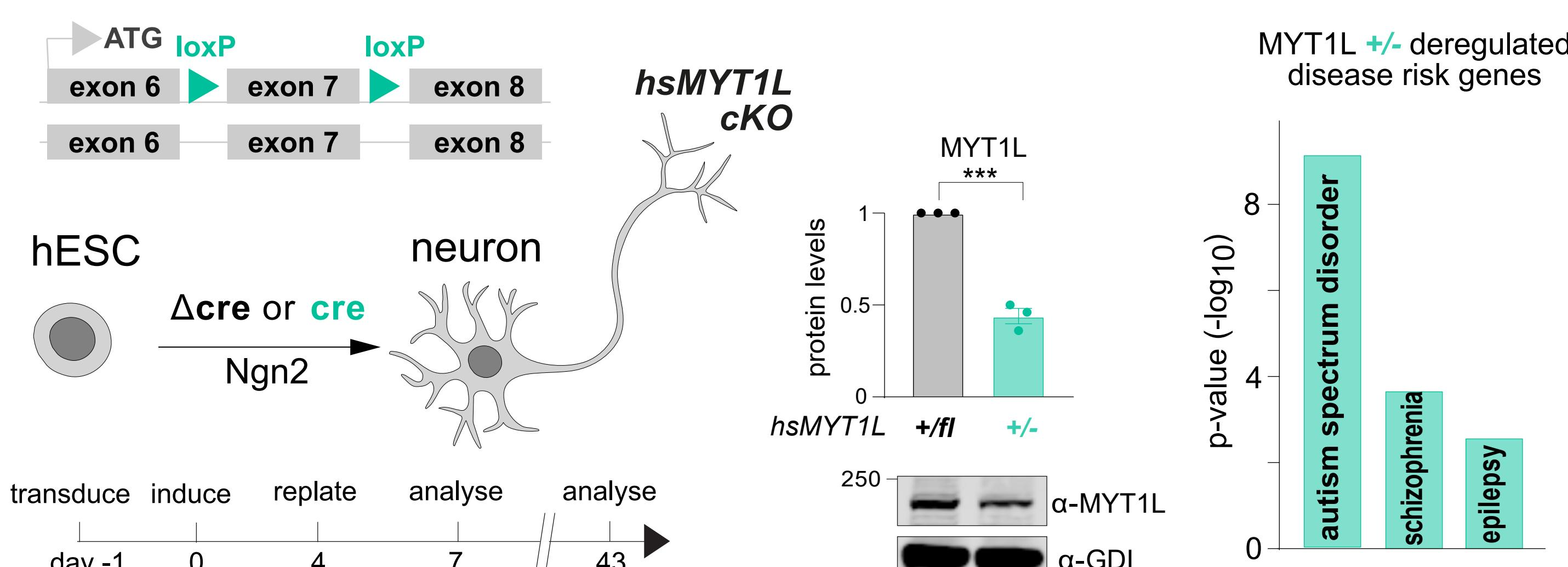
"The brain's utter centrality in the cockpit of our being means its malfunctions are unusually dangerous and far-reaching, attacking not just our health but our whole sense of self in particularly damaging and debilitating ways"

The transcription repressor MYT1L

- neuron specific [1]
- life-long expressed [1]
- important for neuronal cell identity [2]
- mutations are strongly linked to mental disorders such as autism spectrum disorder (ASD), schizophrenia and intellectual disability [3,4]

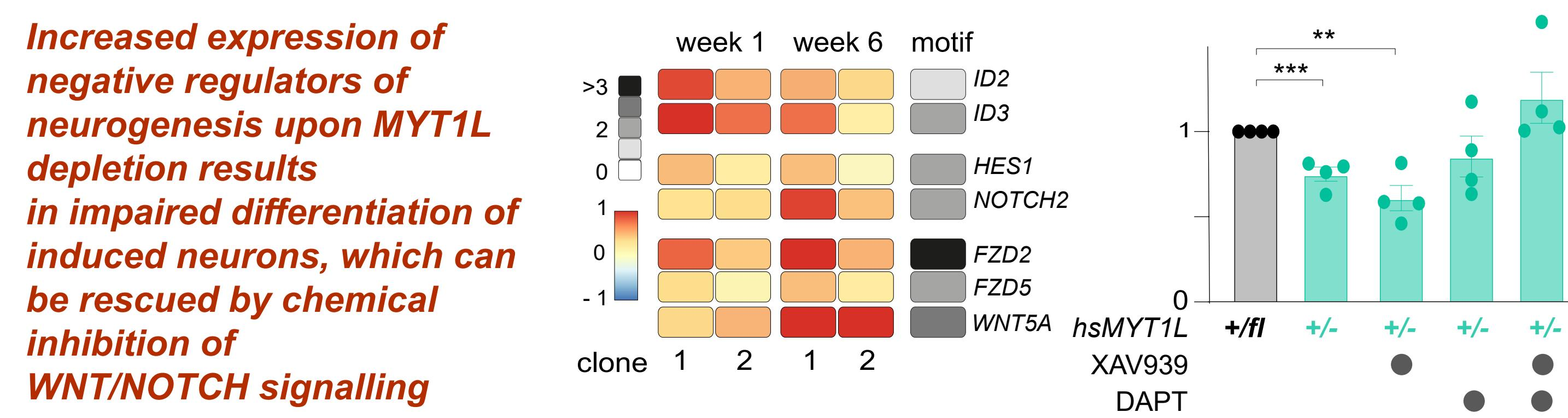
Can Myt1l mutations be causative for neuropsychiatric disorders?

Disease modelling with human induced neurons

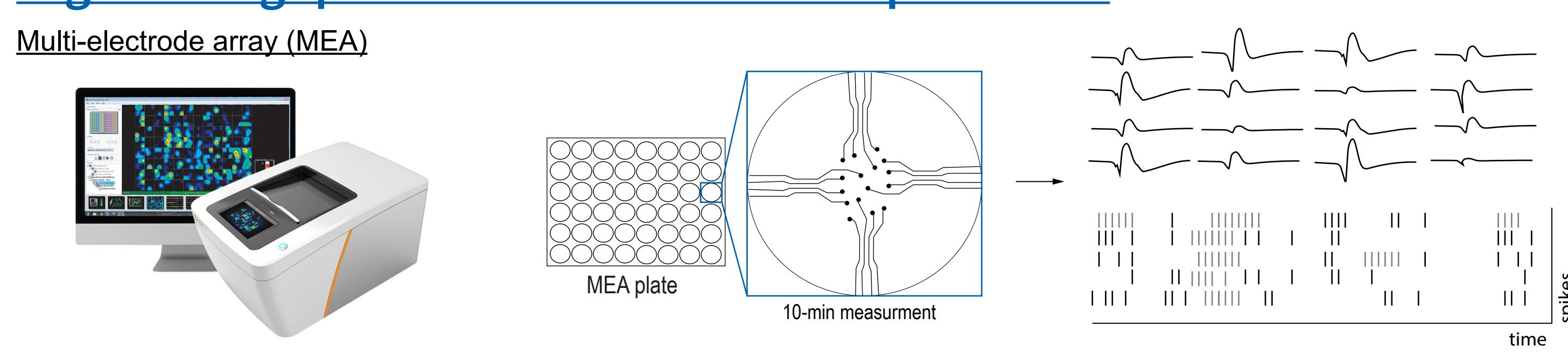


Human embryonic stem (hES) cells can be engineered to carry inducible loss-of-function mutations and can be differentiated into induced neurons (hIN) by overexpression of the proneural transcription factor Ngn2 [5].

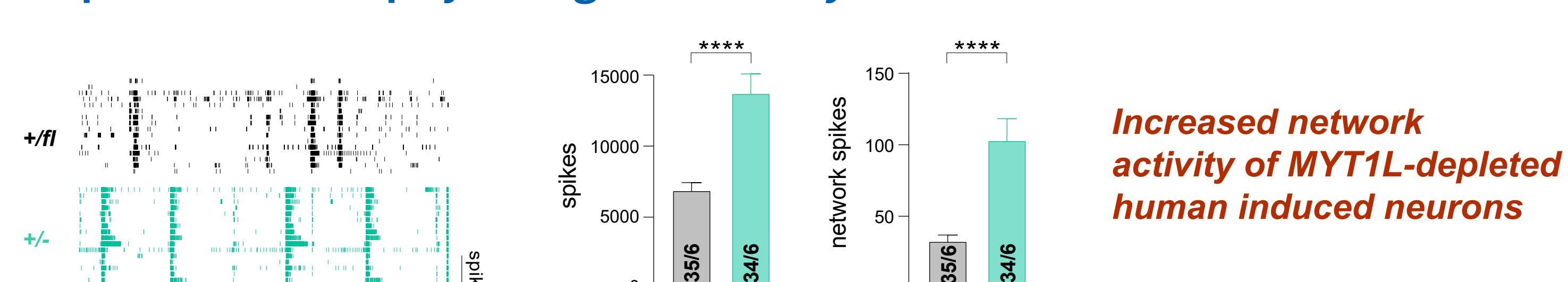
Impaired repression of developmental pathways



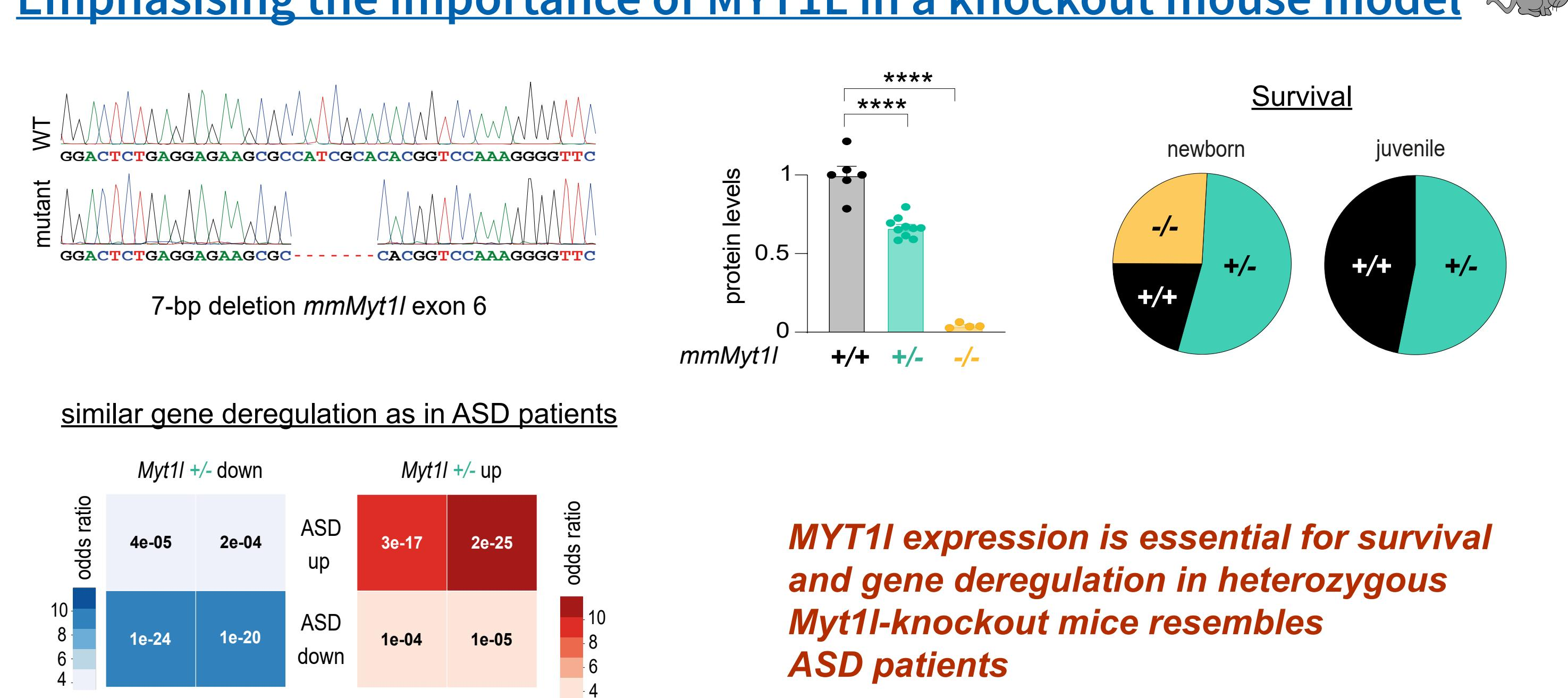
High-throughput measurement of field potentials



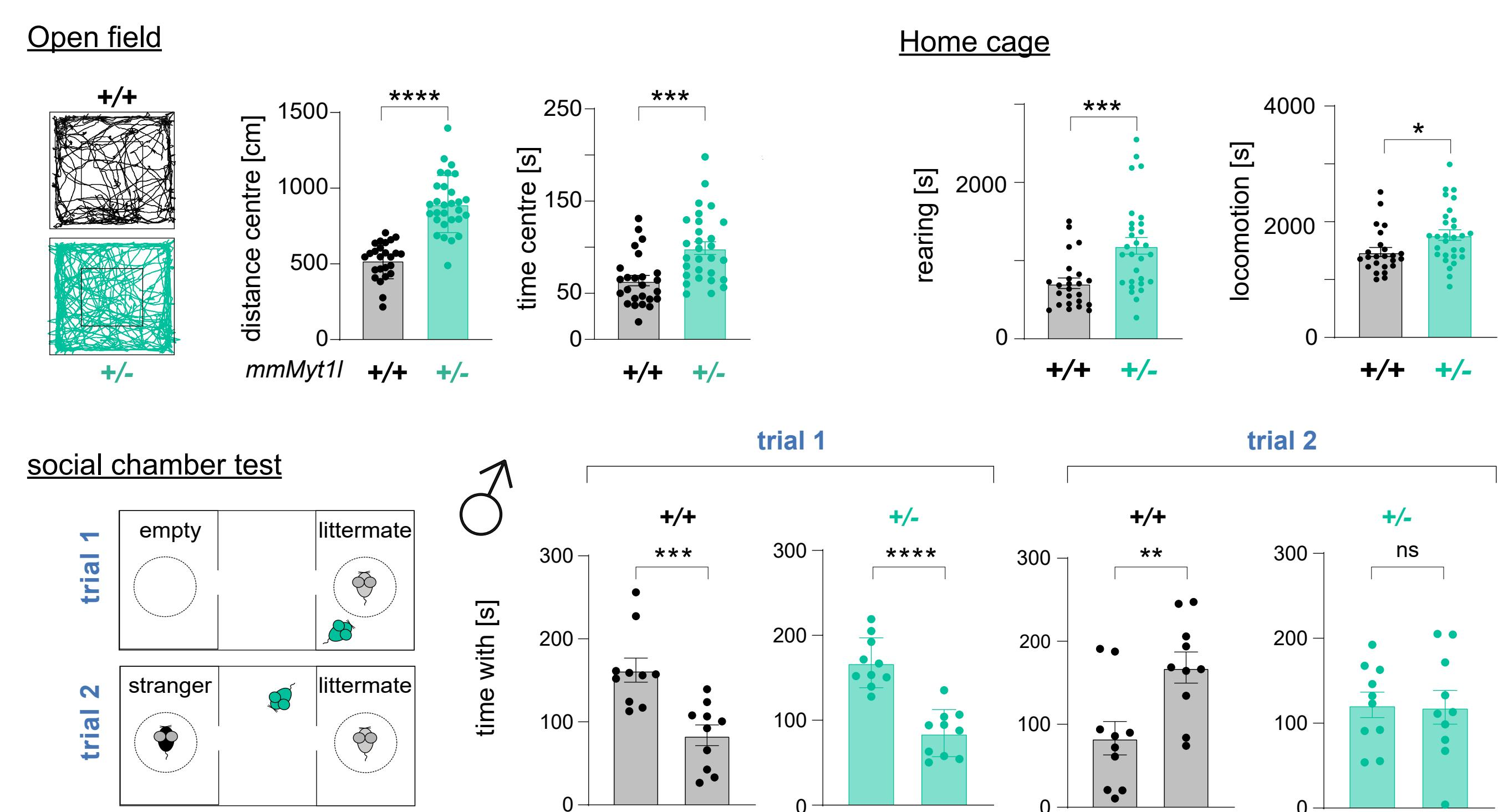
Impaired electrophysiological activity



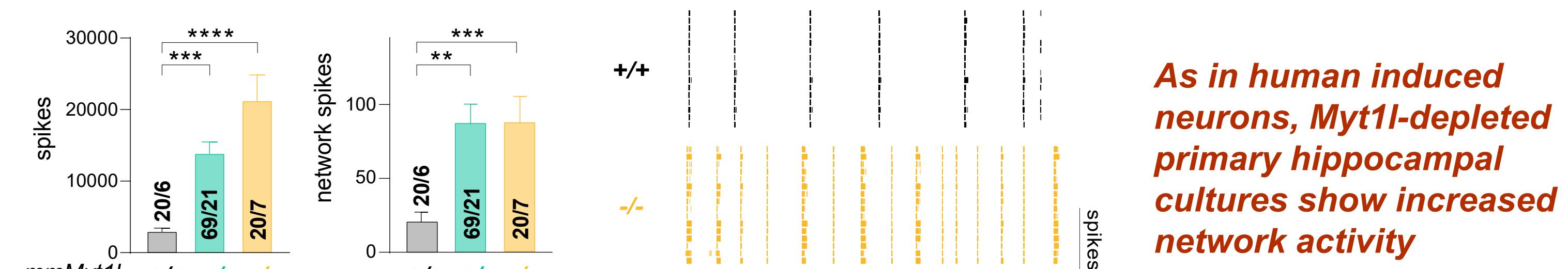
Emphasising the importance of MYT1L in a knockout mouse model



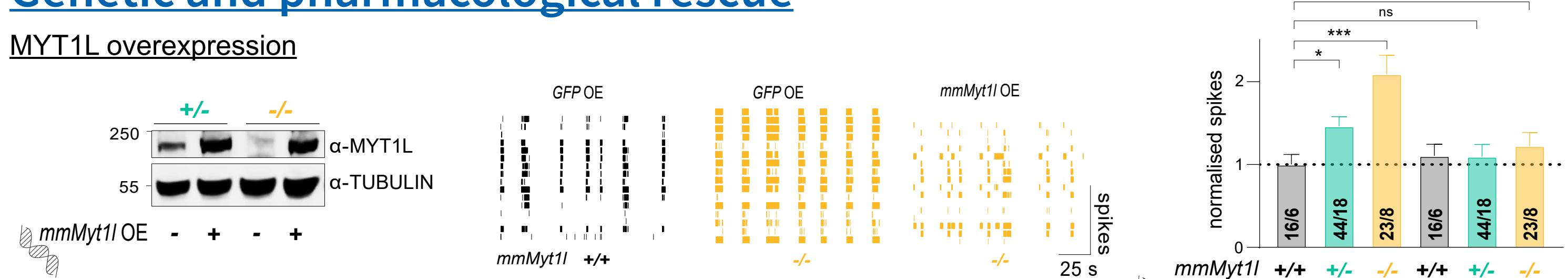
Behavioural abnormalities in Myt1l-mutant mice



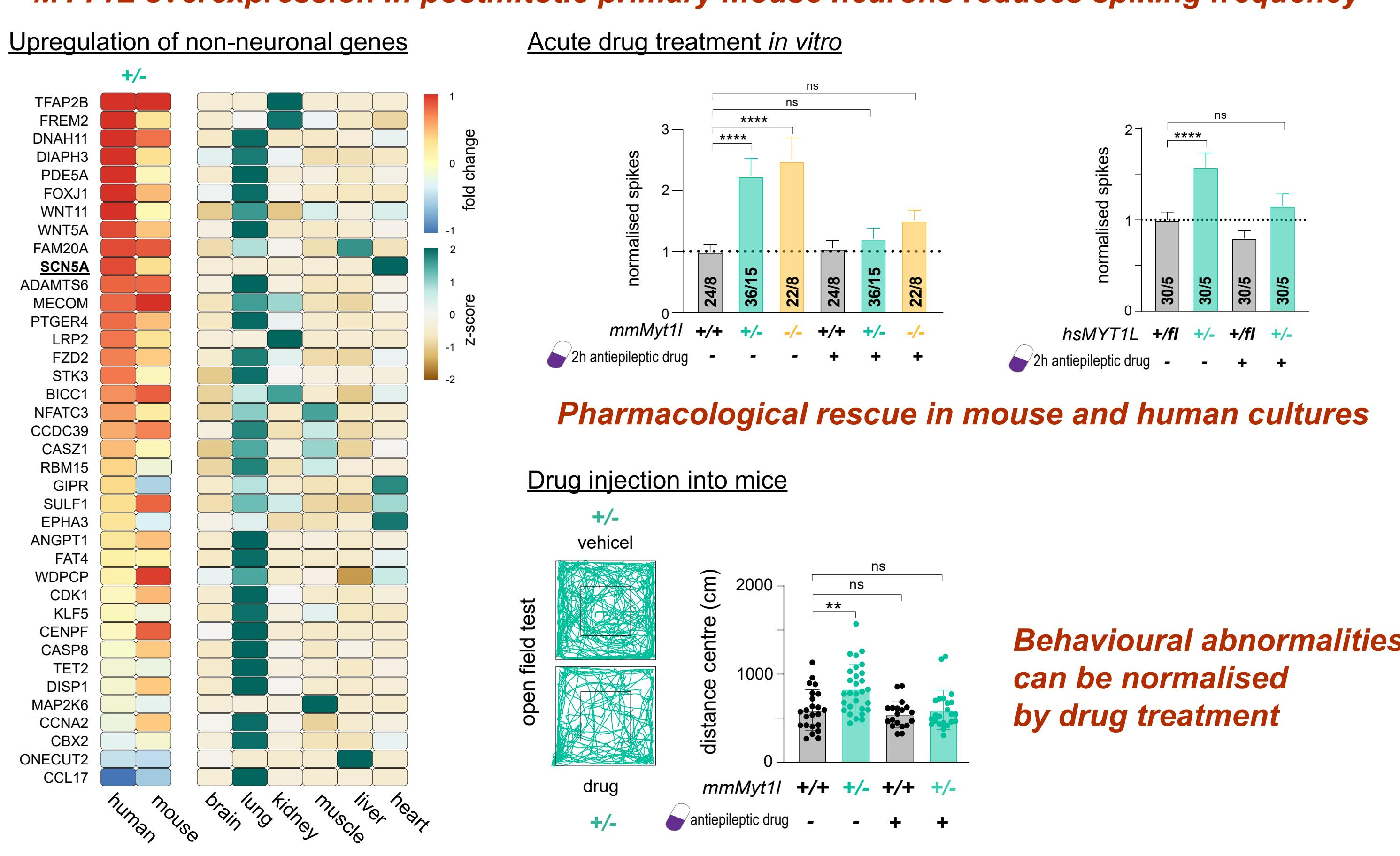
Conserved electrophysiological phenotypes in both models



Genetic and pharmacological rescue



MYT1L overexpression in postmitotic primary mouse neurons reduces spiking frequency



Key points

MYT1L mutations in human and mouse models result in gene deregulation that resembles ASD patients

Myt1l-mutant mice are hyperactive and present with male-specific social deficits

MYT1L-deficient primary mouse and human induced neurons show increased electrophysiological activity, which can be normalised by MYT1L overexpression in postmitotic primary mouse neurons

Application of an FDA-approved antiepileptic drug rescues electrophysiological and behavioural phenotypes

Bar graphs display mean values with number of MEA wells and cells from indicated biological replicates or data points from individual animals and iN batches, error bars = SEM, * p < 0.05, ** p < 0.01, *** p < 0.001, **** p < 0.0001, ns = not significant.