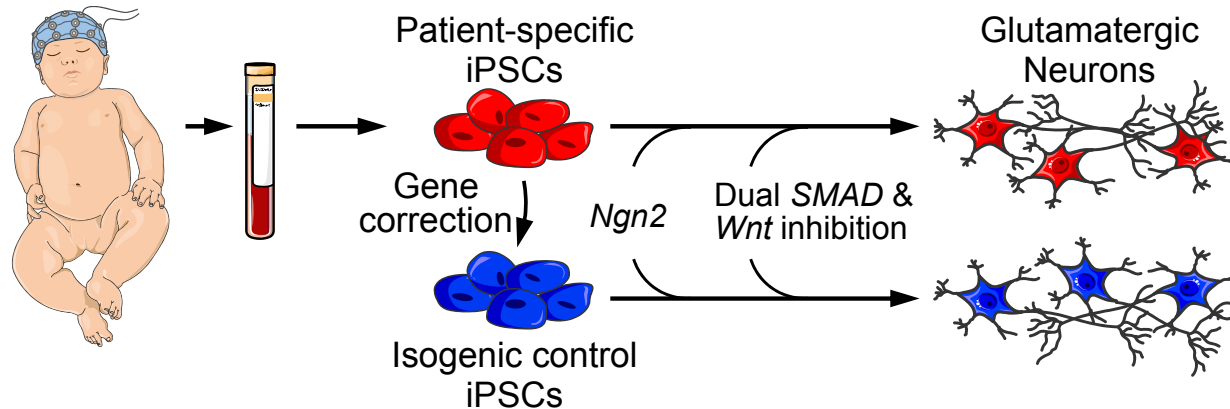


Mechanisms of maladaptive compensation in KCNQ2-related epilepsy in patient-specific iPSC-derived neurons

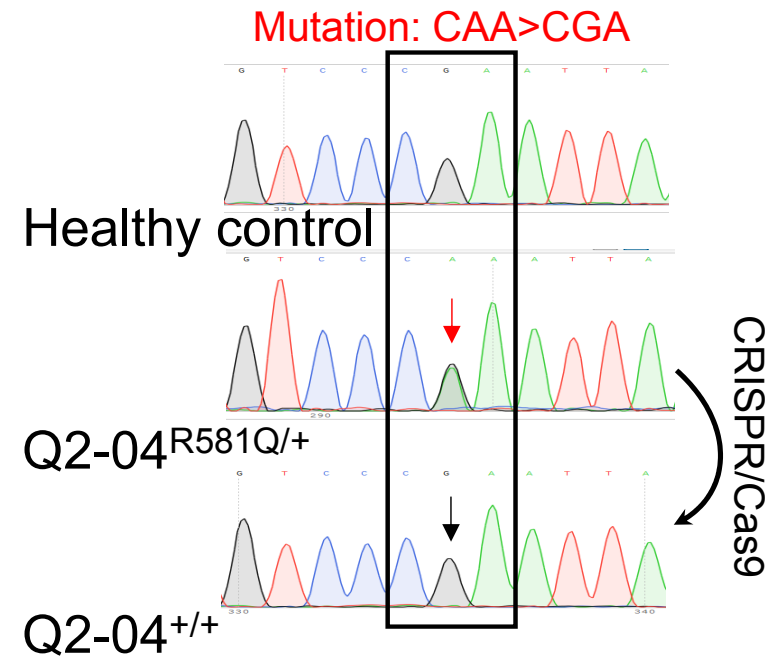
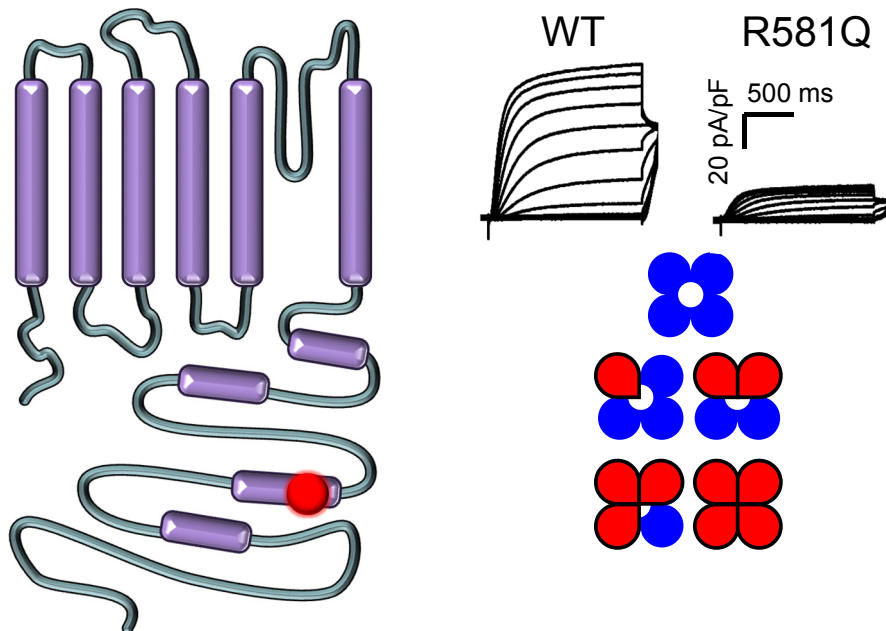
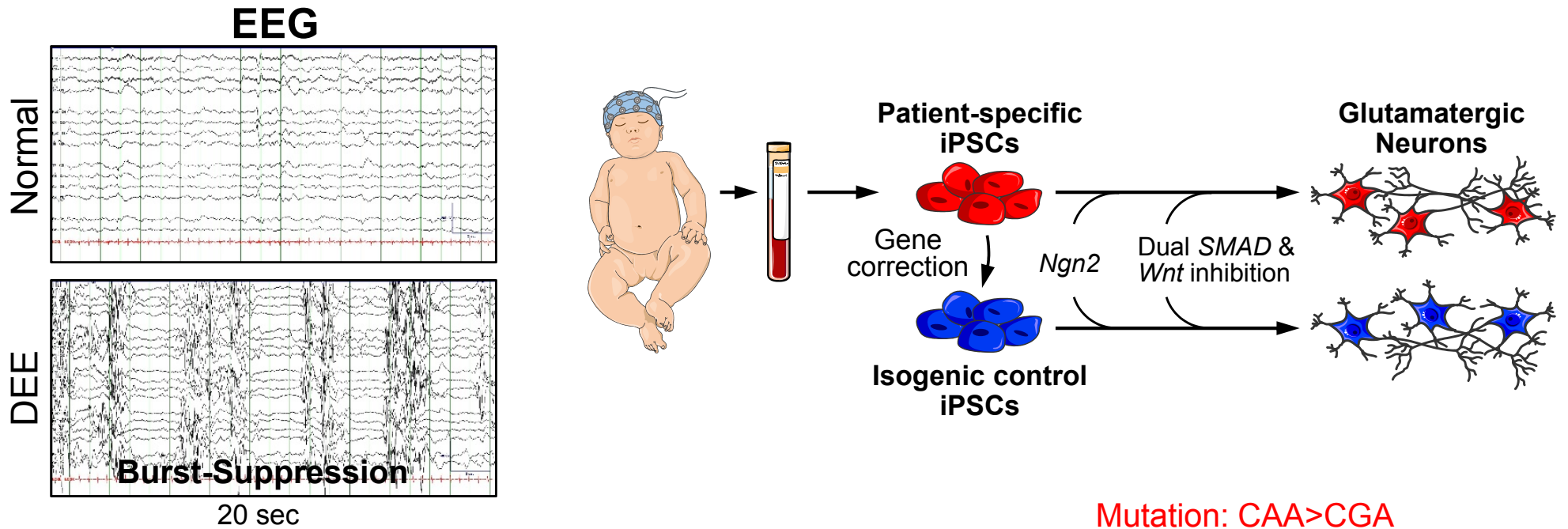
Dina Simkin¹

Adil Wafa¹, Mennat Gharib¹, Kelly Marshall¹, Alfred L. George² & Evangelos Kiskinis¹

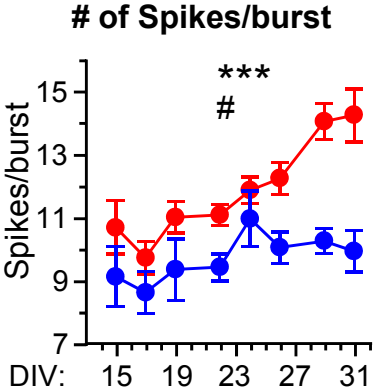
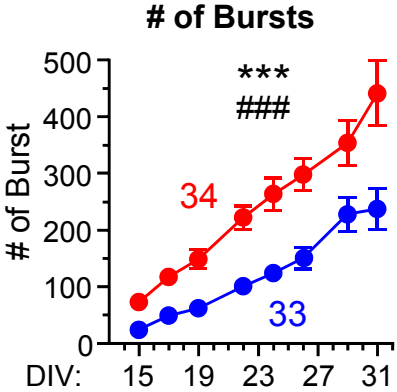
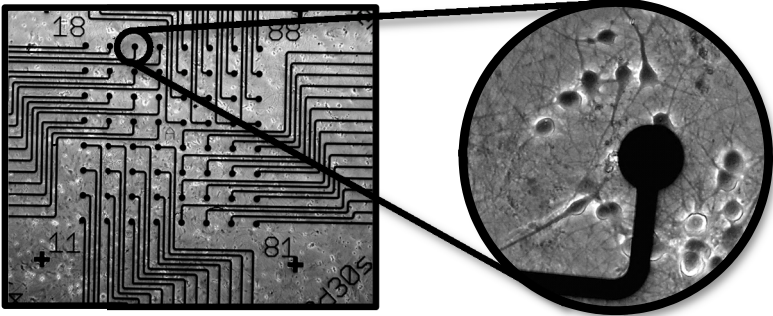
¹Department of Neurology; ²Department of Pharmacology,
Northwestern University, Feinberg School of Medicine, Chicago, IL



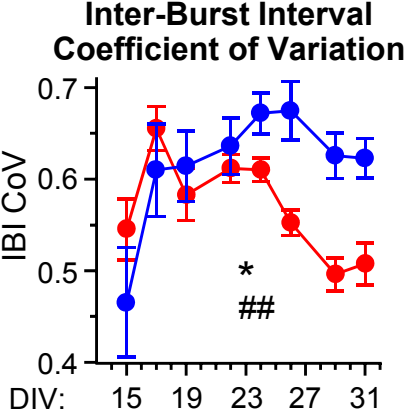
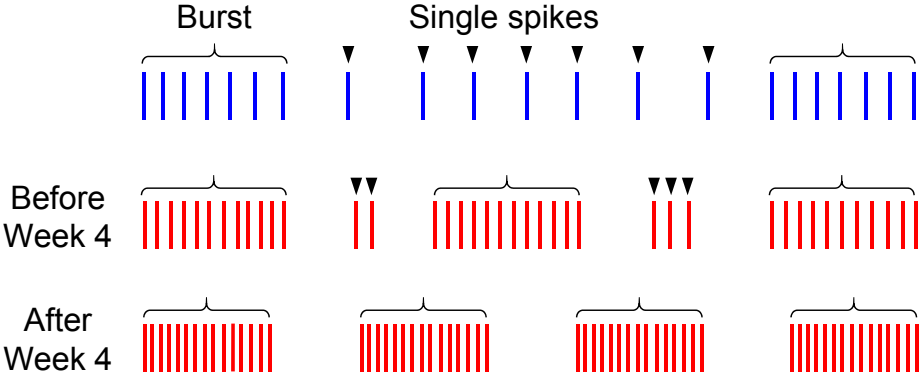
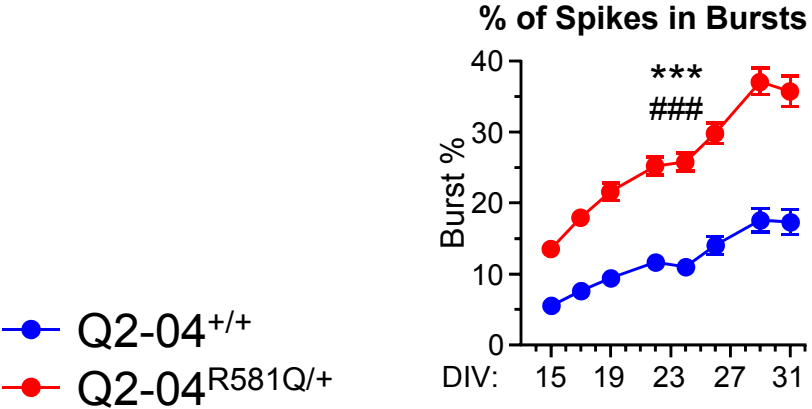
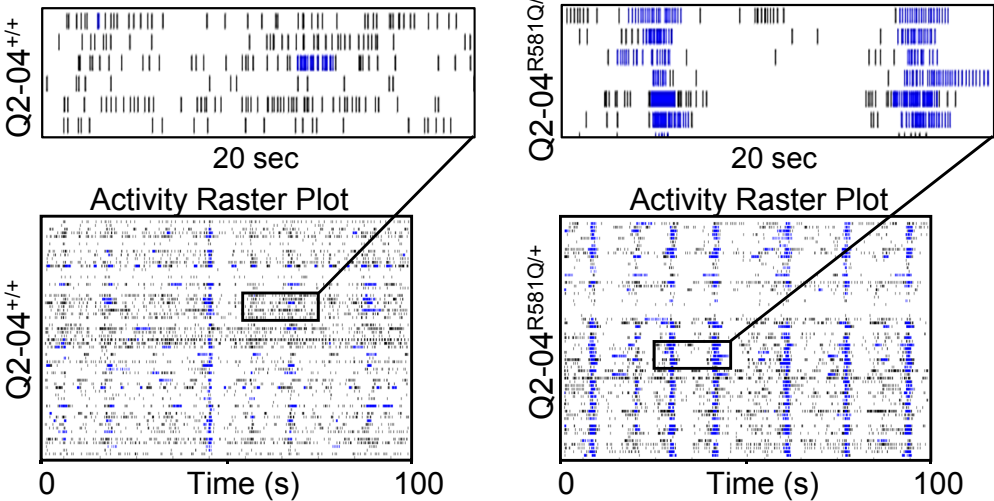
Developmental and epileptic encephalopathy (DEE) associated with KCNQ2-R581Q loss-of-function mutation



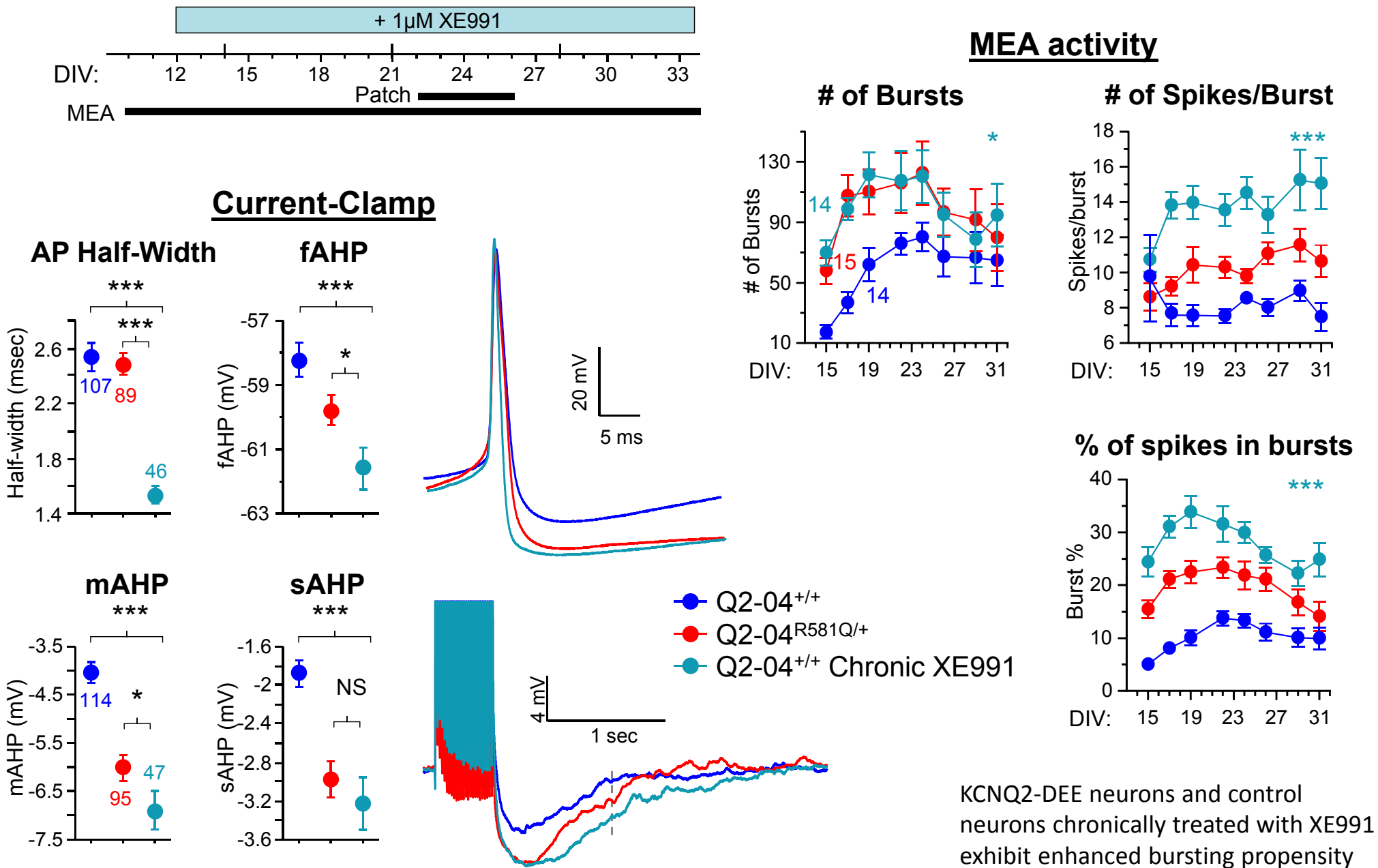
KCNQ2-DEE neurons exhibit enhanced spontaneous bursting on MEAs



MEA Activity Raster Plot



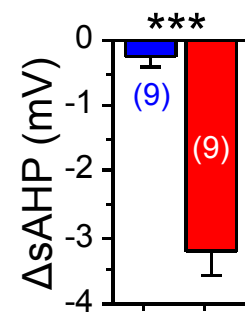
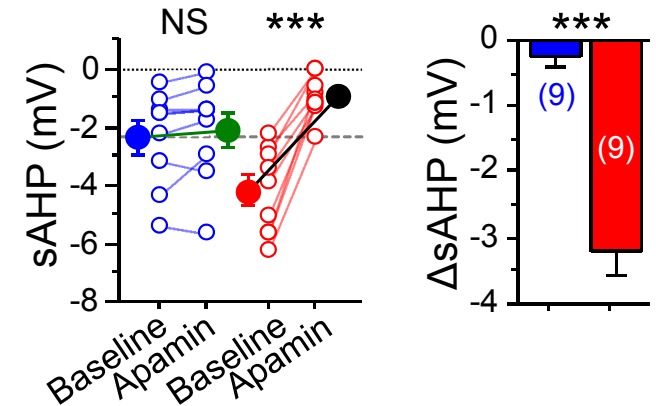
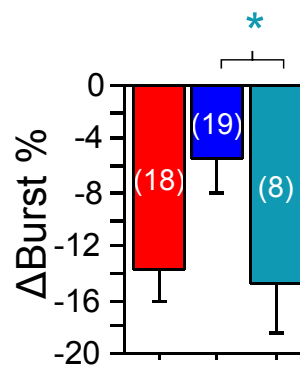
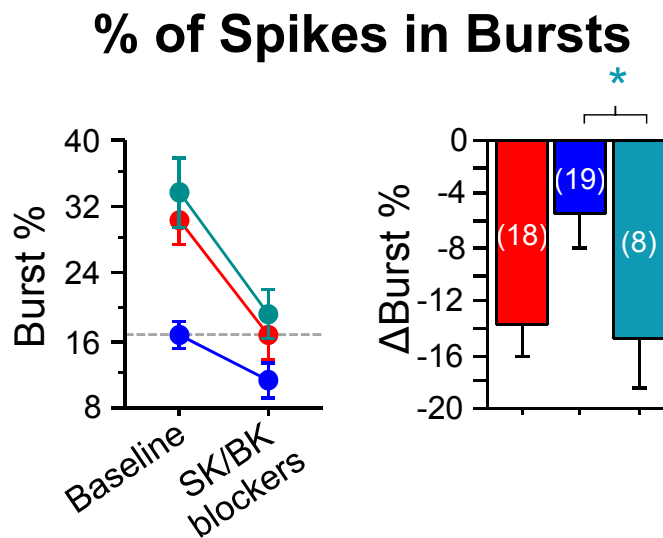
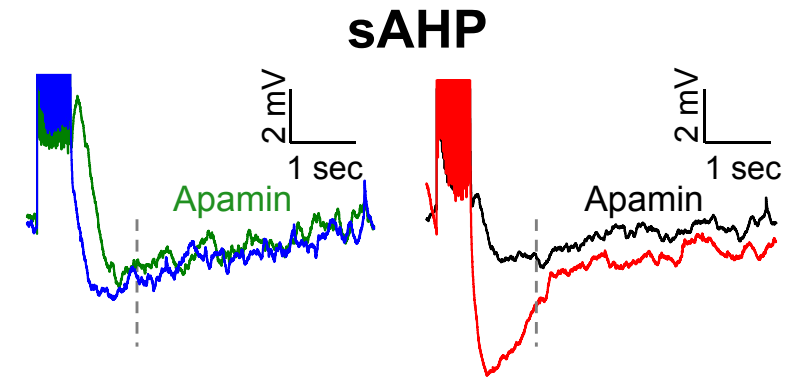
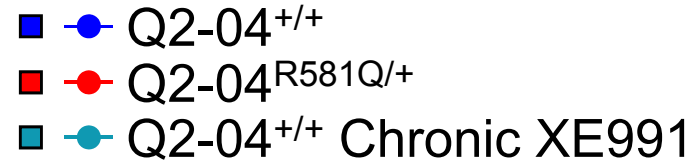
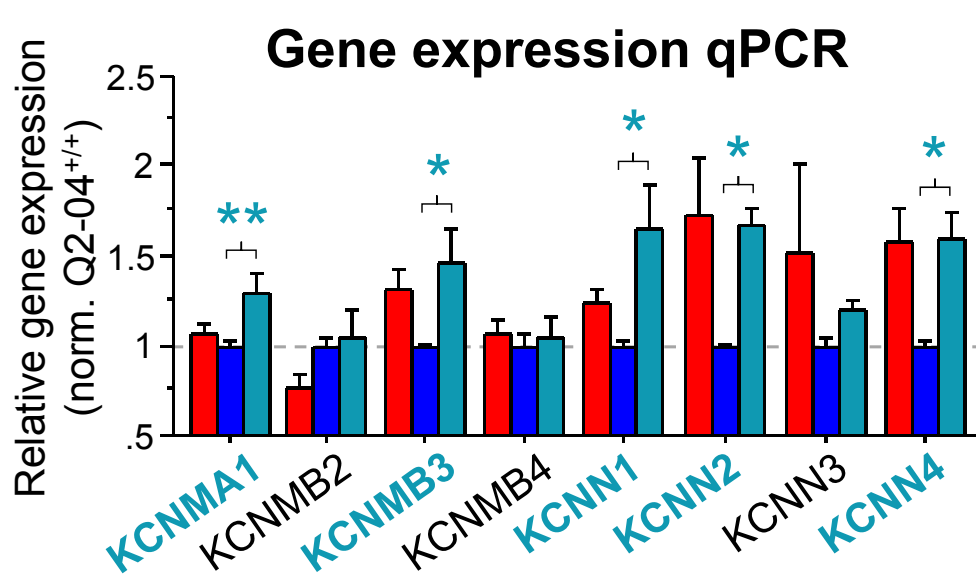
Chronic inhibition of M-current in control neurons phenocopies KCNQ2-DEE



KCNQ2-DEE neurons and control neurons chronically treated with XE991 exhibit faster AP repolarization and larger post-burst AHPs

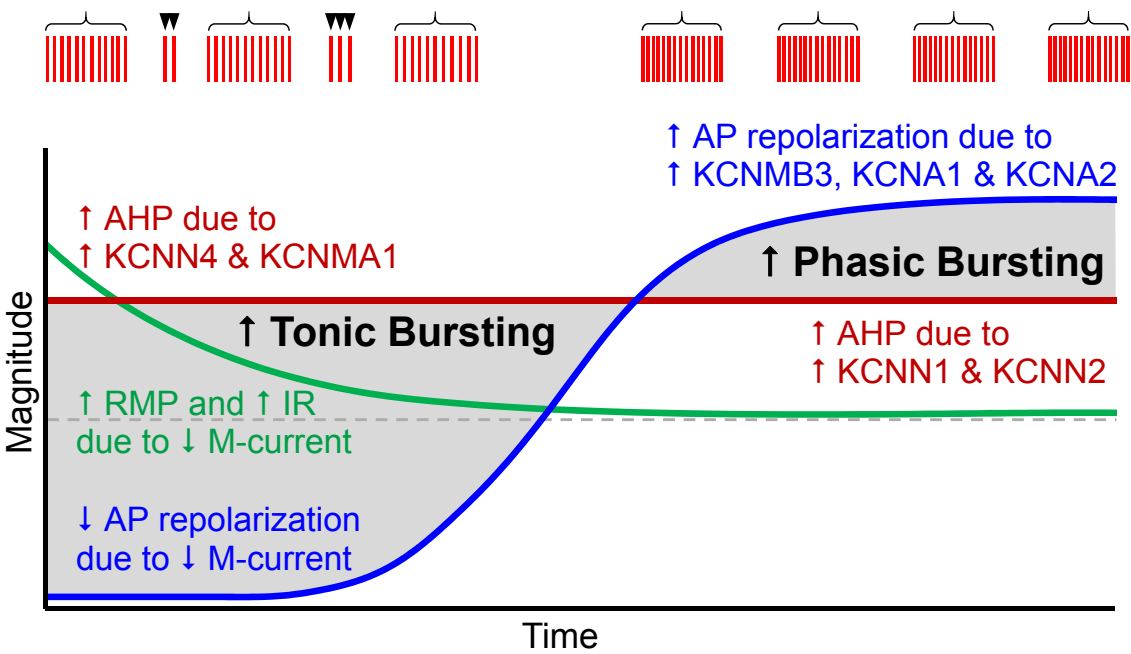
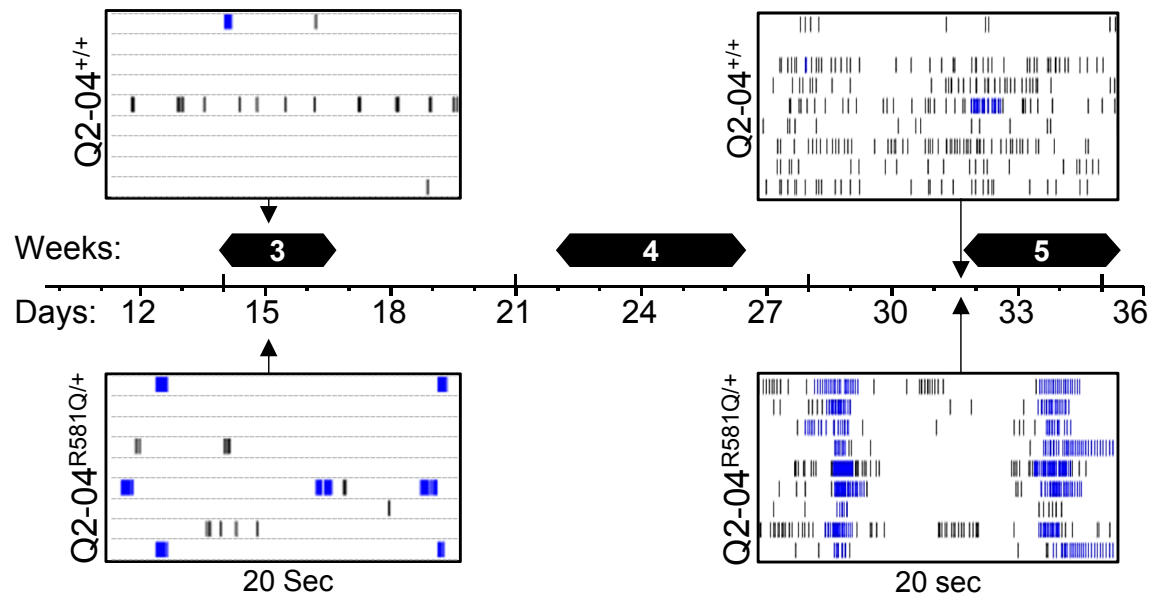
KCNQ2-DEE neurons and control neurons chronically treated with XE991 exhibit enhanced bursting propensity

Maladaptive enhancement of SK/BK channels leads to enhanced AHPs and bursting in KCNQ2-DEE neurons and XE991 treated control neurons



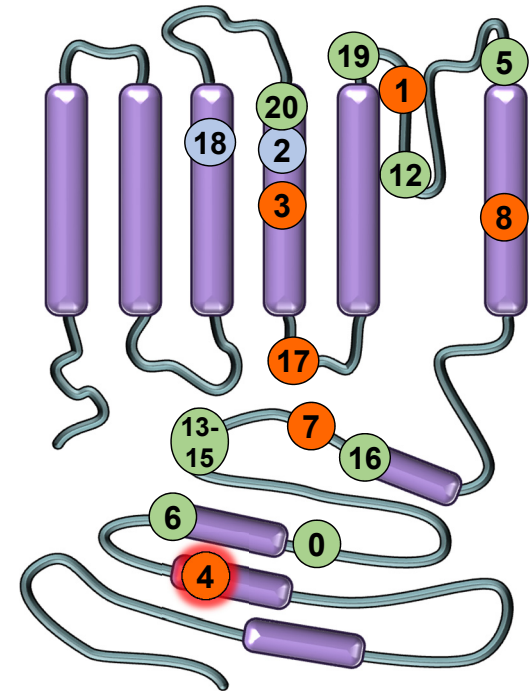
Apamin (500nM): SK channel blocker
Paxilline (20μM): BK channel blocker

Developmental Time Course of KCNQ2-DEE Neurons



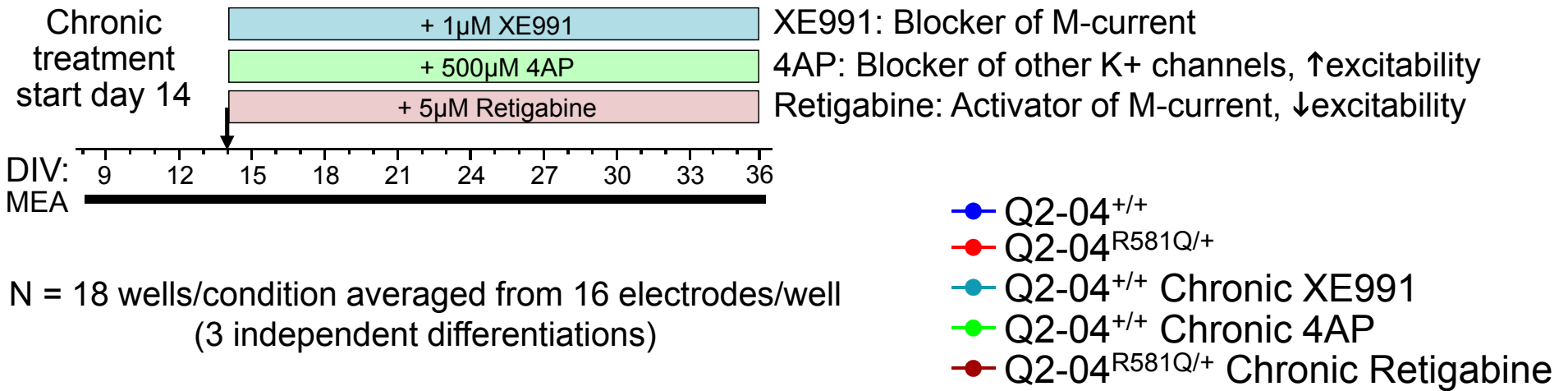
Unanswered questions:

1) Is maladaptive Ca²⁺-dependent K⁺ channel upregulation and bursting a result of reduced M-current specifically or the resulting hyperexcitability?

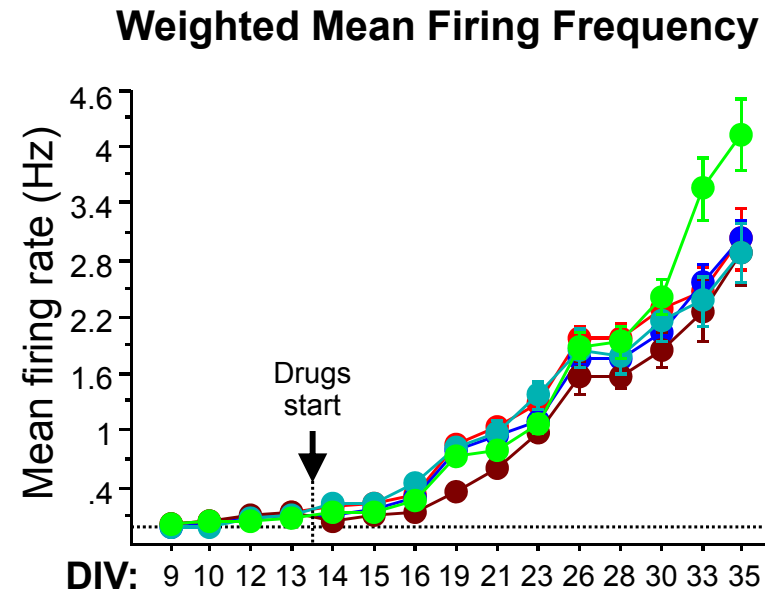
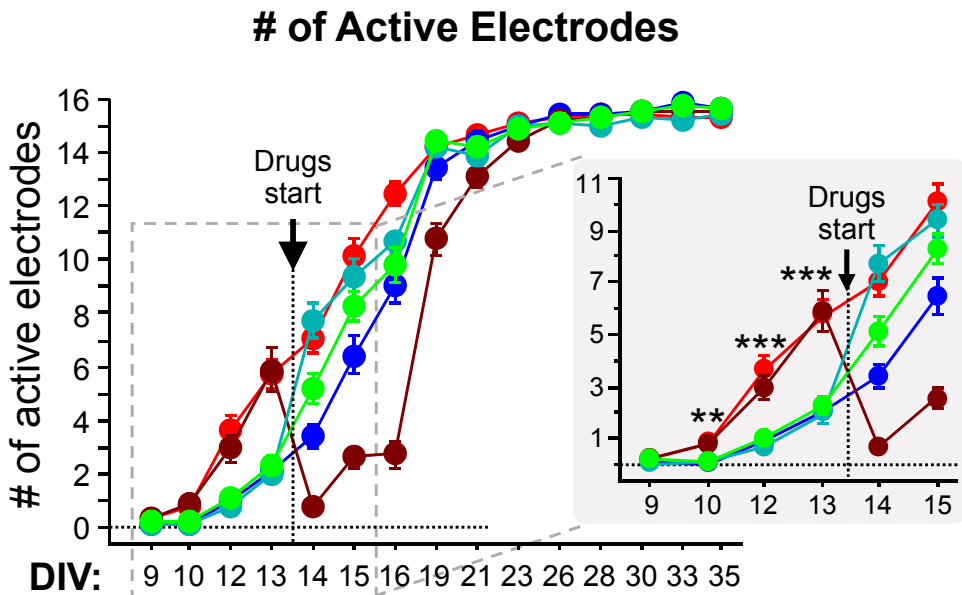


2) Is this bursting a common phenotype amongst KCNQ2-DEE patient neurons with other KCNQ2 mutations?

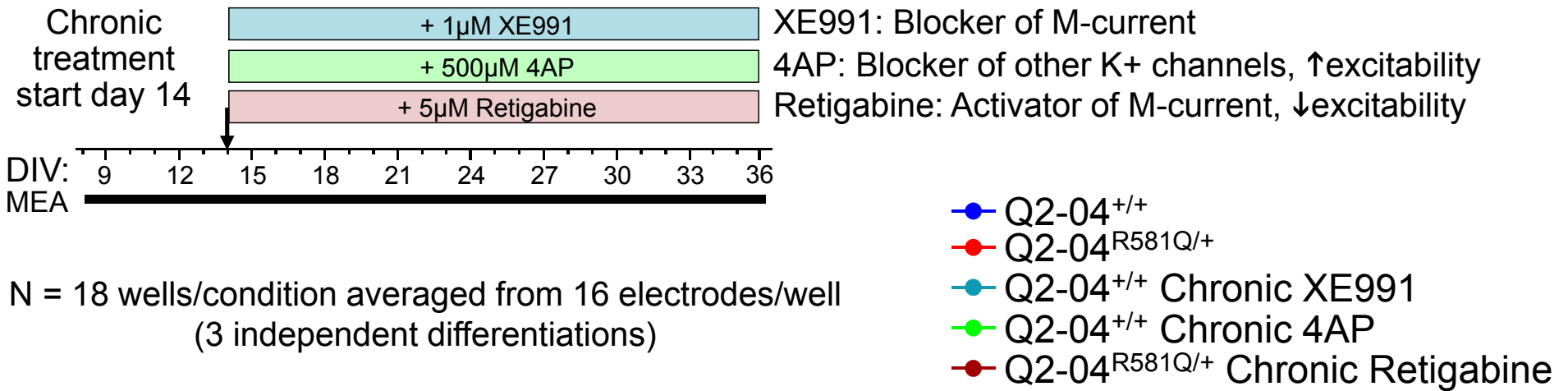
Chronic inhibition of K⁺ channels and increased neuronal activity with 4AP does not phenocopy KCNQ2-DEE maladaptive phenotype



Early-onset of functional maturation

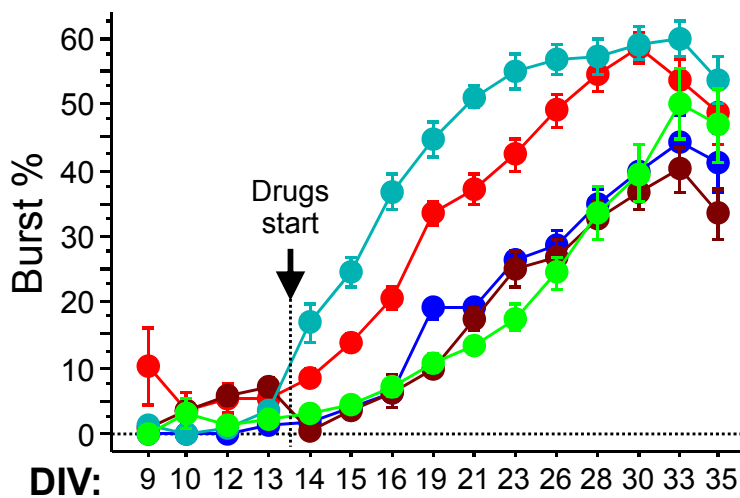


Chronic inhibition of K⁺ channels and increased neuronal activity with 4AP does not phenocopy KCNQ2-DEE maladaptive bursting phenotype

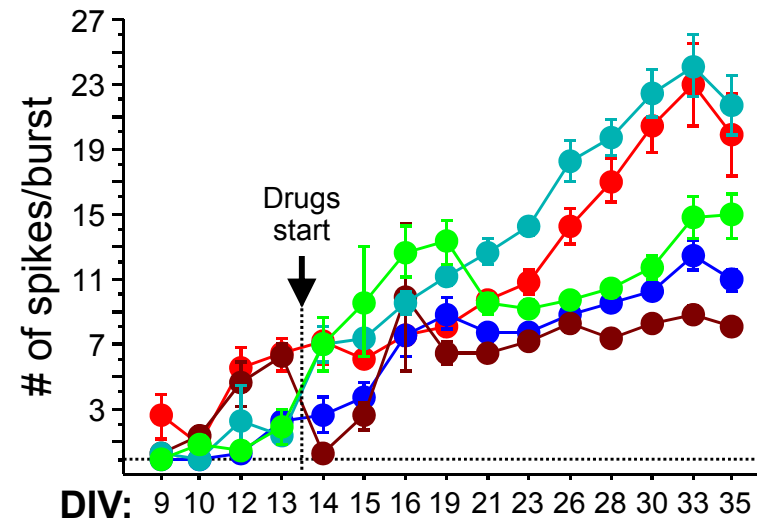


Bursting Propensity

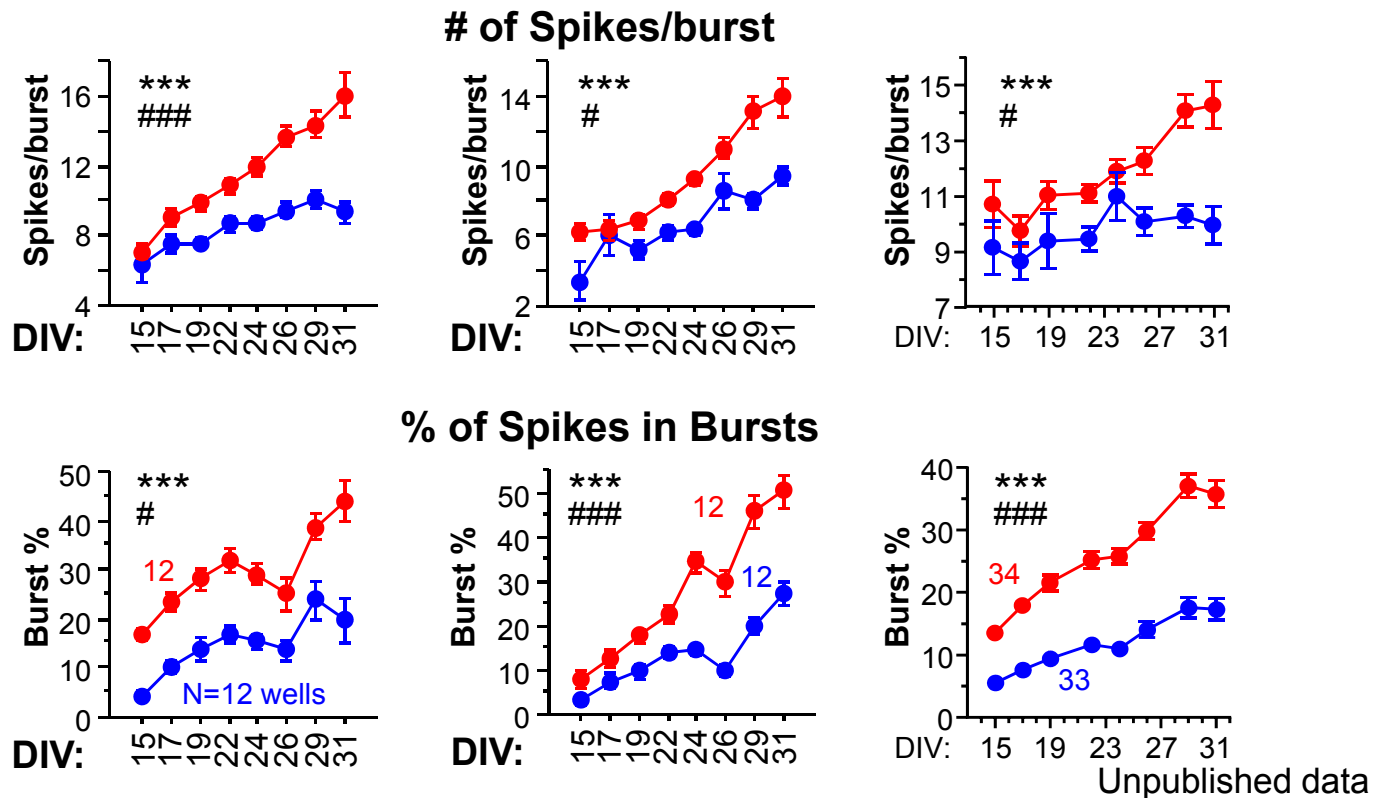
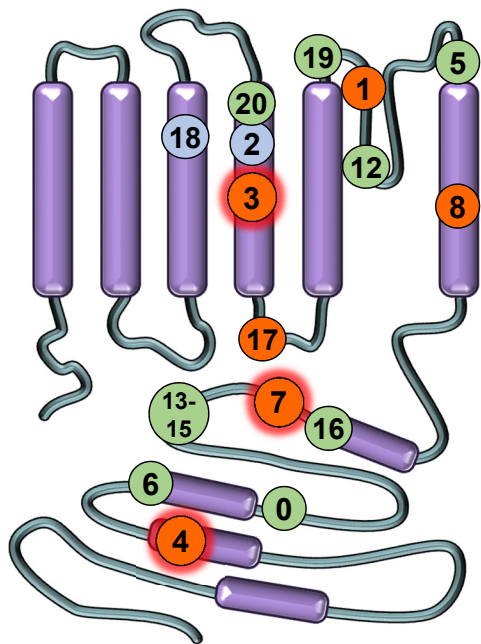
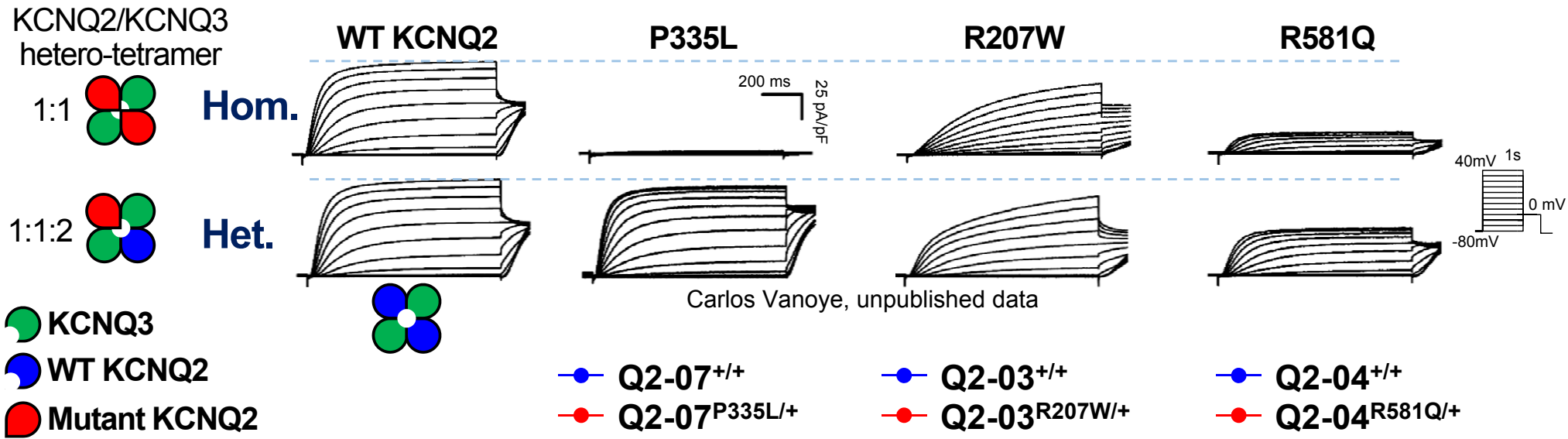
% of Spike in Bursts



of Spikes per Burst



Common bursting phenotype amongst KCNQ2-DEE patient-specific iPSC-derived excitatory neurons with different KCNQ2 mutations



Acknowledgments

Northwestern Medicine
Feinberg School of Medicine

Evangelos Kiskinis & Lab

Kelly Marshall

Adil Wafa

Mennat Gharib

Al George & Lab

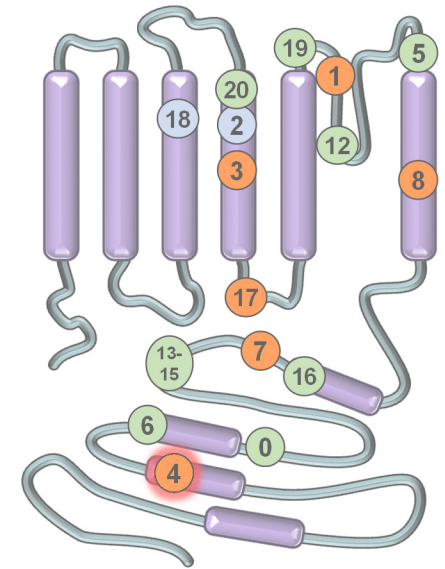
 Ann & Robert H. Lurie
Children's Hospital of Chicago

John Millichap

Dyshomeostatic modulation of Ca²⁺-activated K⁺ channels in a human neuronal model of KCNQ2 encephalopathy



Dina Simkin^{1,2}, Kelly A Marshall¹, Carlos G Vanoye², Reshma R Desai², Bernabe I Bustos¹, Brandon N Piyevsky¹, Juan A Ortega¹, Marc Forrest^{3,4}, Gabriella L Robertson¹, Peter Penzes^{3,4}, Linda C Laux⁵, Steven J Lubbe¹, John J Millichap⁵, Alfred L George Jr^{2*}, Evangelos Kiskinis^{1,3*}



NYSCF
The New York
Stem Cell Foundation
NYSCF – Robertson Investigator



National Institute of
Neurological Disorders
and Stroke

U54 NS108874

Falk Medical
Research Trust