

# High-throughput assay of seizurogenic activity using multiwell microelectrode array technology

\*D. C. MILLARD, A. M. NICOLINI, S. A. CHVATAL, M. W. BROCK, K. K. COOK, J. D. ROSS;  
Axion BioSystems, Inc., 1819 Peachtree Road, Suite 350, Atlanta, GA, 30309

220.03

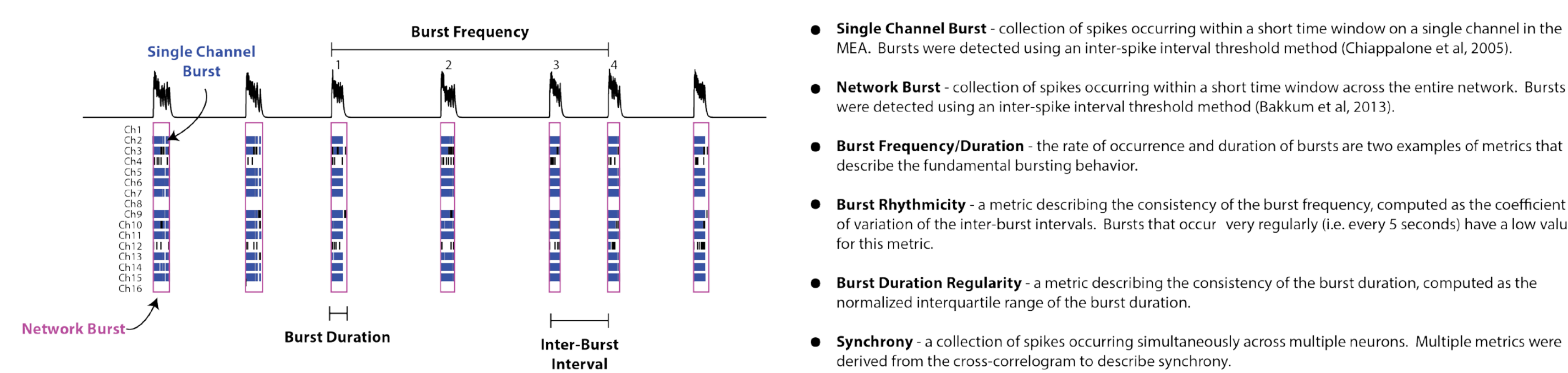


## I. Introduction

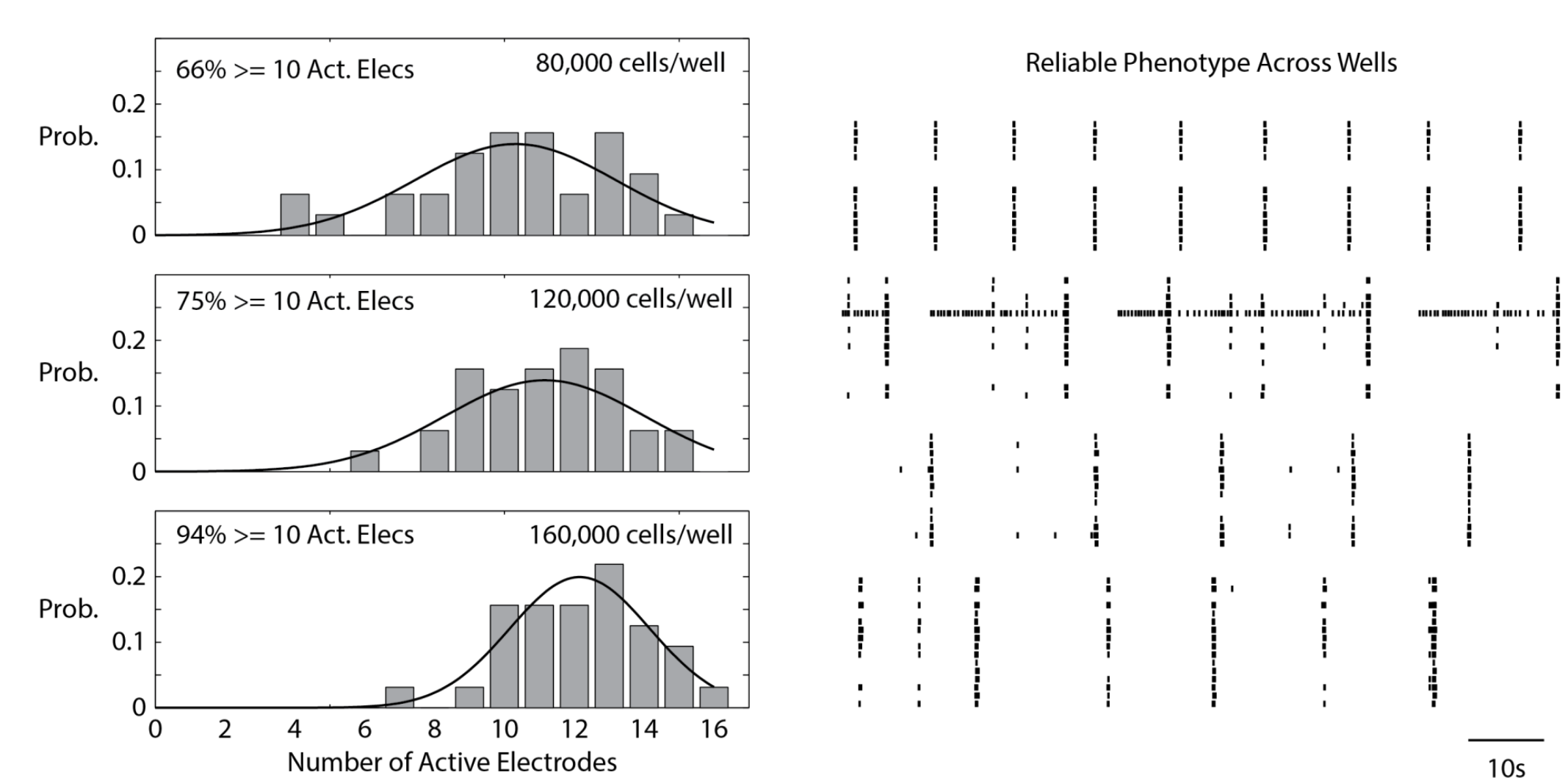
- The lack of advancement in anti-epileptic drugs (AEDs) over the last 30 years, along with the continued need for improved proconvulsant screening in drug safety, motivates the need for new assays of seizurogenic neural activity.
- Here, we present the development of an *in vitro* assay of seizurogenic activity based upon the Axion BioSystems Maestro multiwell MEA system, using previously published metrics for quantifying bursting and synchrony within networks of cryopreserved cortical neurons.
- Our results support the use of multiwell MEA technology for the high-throughput evaluation of complex neuronal networks *in vitro* to inform the development of AEDs, while also quantifying the proconvulsant risk of candidate pharmaceuticals in a pre-clinical setting.

## II. Methods

- Process**
- Cell Source – Rat Cortical Neurons (QBM Cell Science)
  - Cell Density –  $8 \times 10^4$  to  $1.6 \times 10^5$  per well
  - Surface Coating – Polyethylenimine, Laminin
  - MEAs – 12-, 48- and 96-well (Axion BioSystems)
- Acquisition**
- Inclusion Criteria – At least 5 spikes recorded/minute for a given electrode (McConnell et al, 2012).
  - Settings – Signals acquired from 200-3000Hz, Spike detected at 6 x Std. Dev. of noise
  - Analysis – MatLAB, Axion BioSystems Neural Metric Tool
- Application**
- Compound Sensitivity – Compounds were prepared in DMSO such that the final [vehicle]  $\leq 0.1\%$
  - Compound Dosing – Compounds were dosed in both single dose and sequential dose formats
  - Electrical Stimulation – a  $\pm 800$ mV, 1ms duration, biphasic voltage stimulus was applied through 1 electrode per well across the MEA plate at 0.1-0.2 Hz.
  - Optogenetics – Chr2 transfected neurons were electrically stimulated as above, then optically stimulated using blue light (475nm) (all at 0.1 Hz). Neurons were then dosed with 100 $\mu$ M Picrotoxin and the stimulations were repeated.

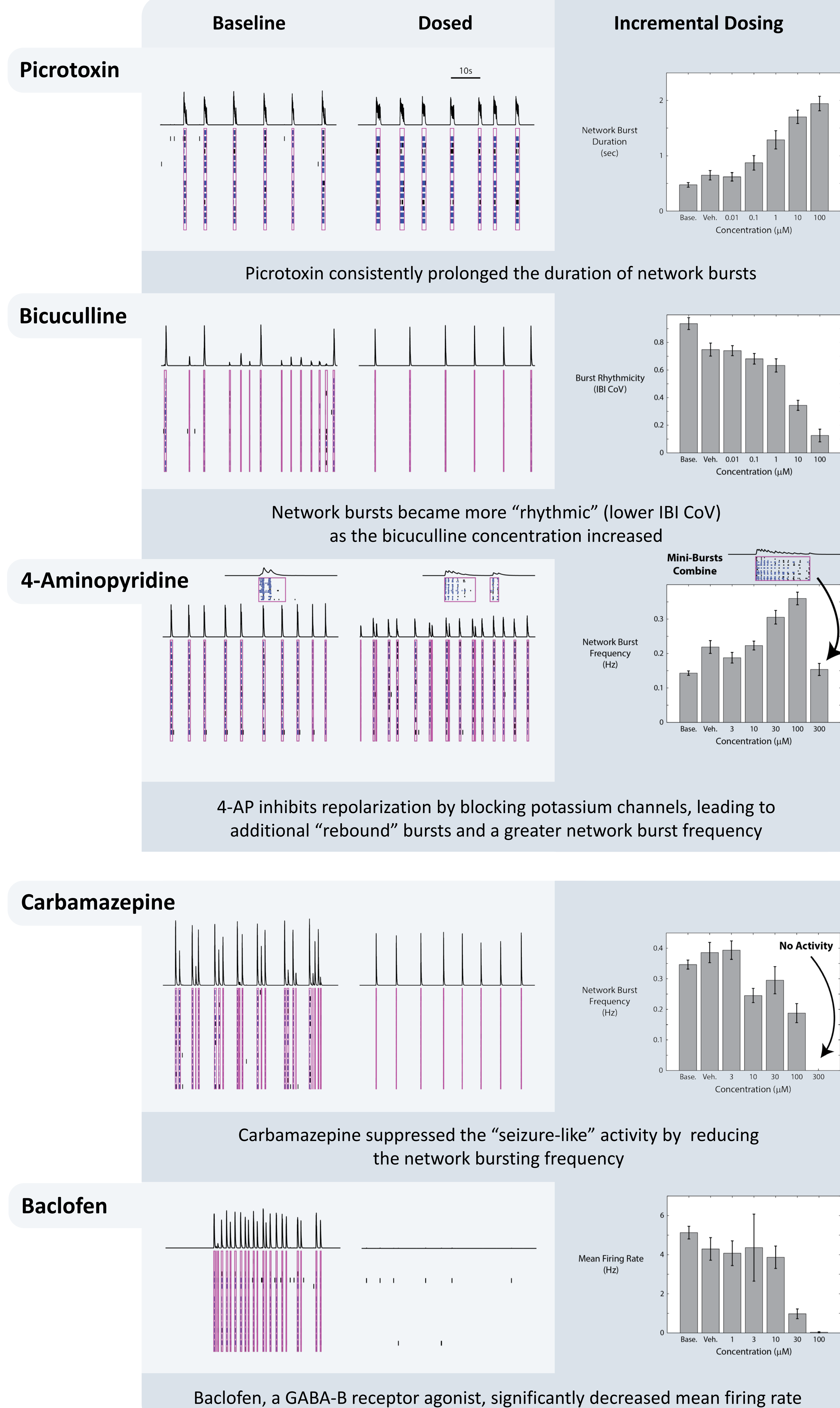


## III. Cell Density Variation

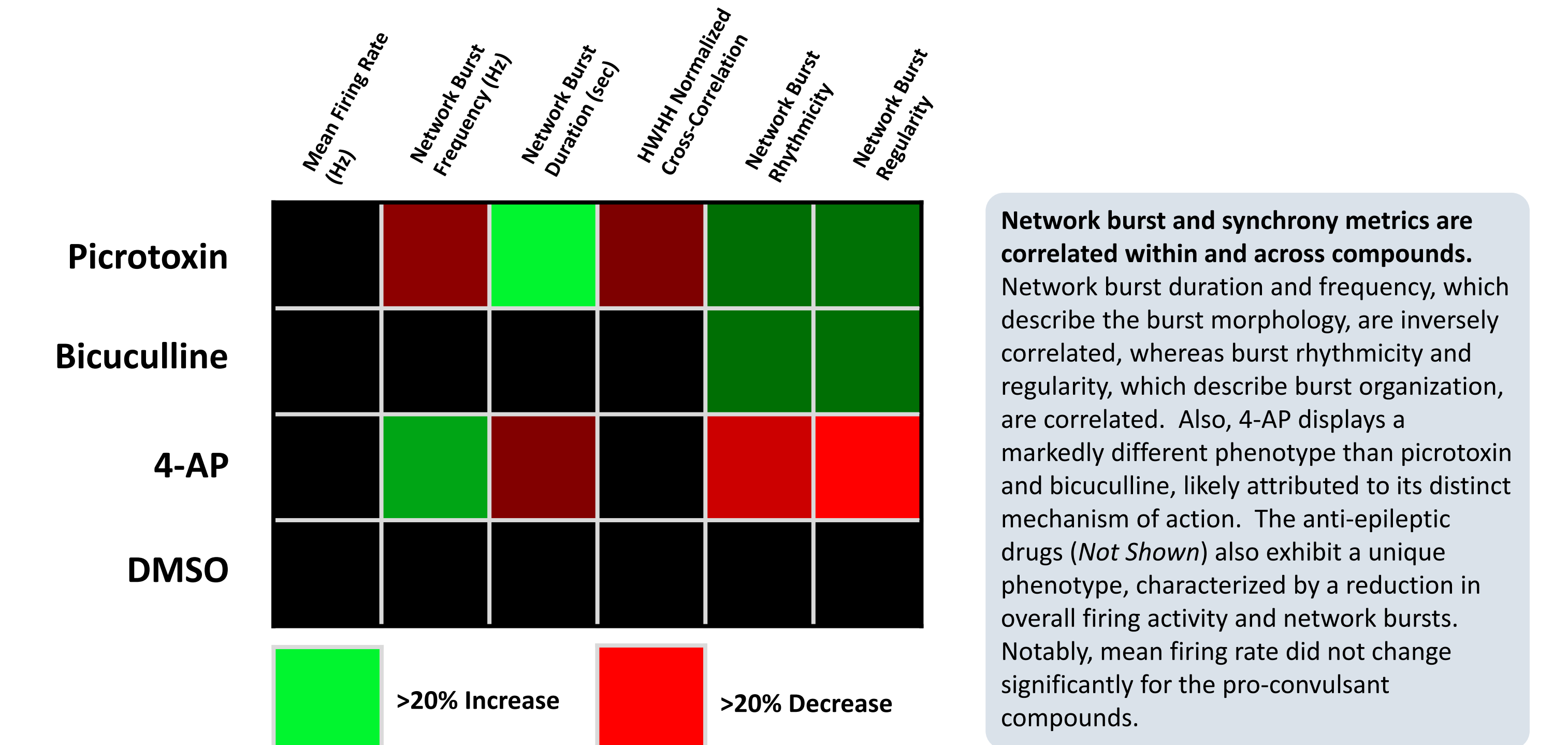


High density cultures ensure electrode coverage and reliable bursting phenotypes. Increased plating densities produce more active electrodes per well. At the highest plating density, 94% of wells had at least 10 active electrodes. Also, at this density the network bursting phenotype was consistent and regular, providing a uniform baseline across wells.

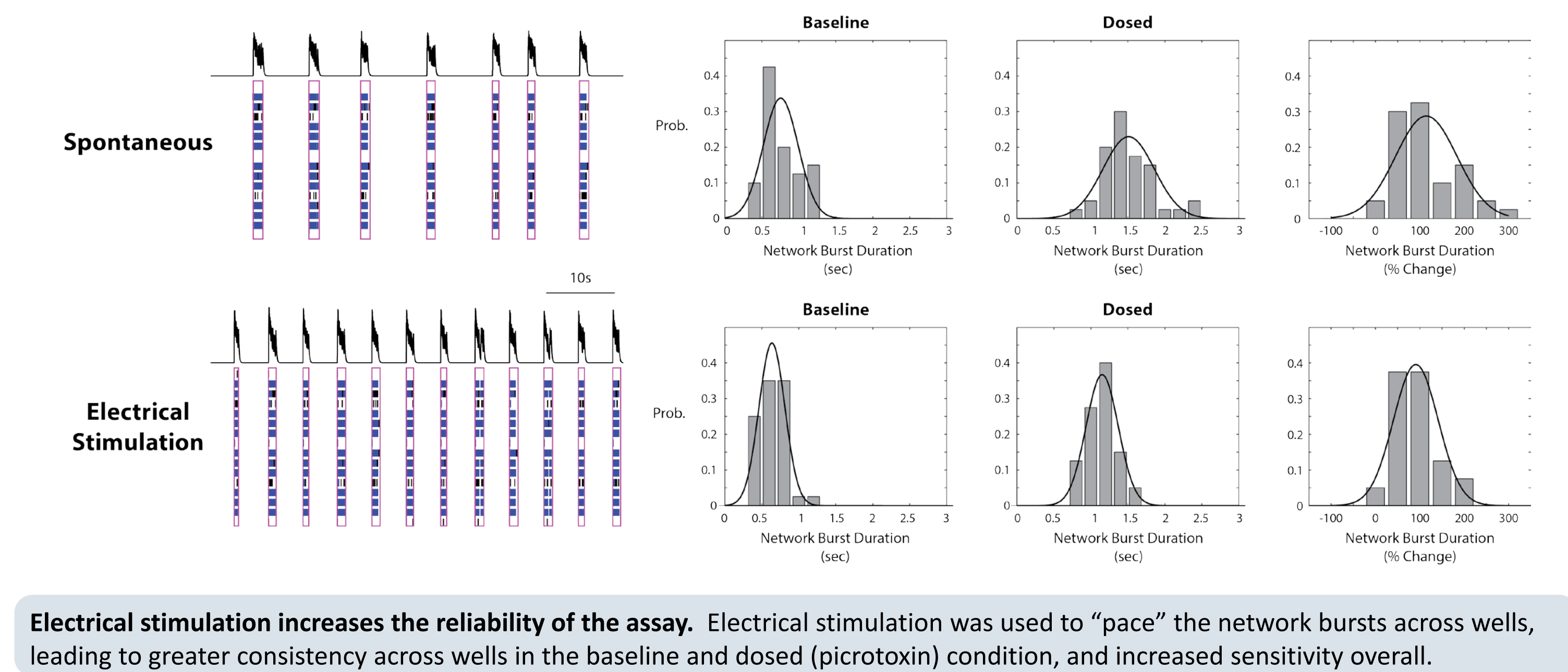
## IV. Neural Metrics



## V. Neural Metrics Continued



## VI. Electrical Stimulation



## VII. Conclusion

The network activity of dissociated cortical cultures, quantified through burst and synchrony metrics, was extremely sensitive to known pro-convulsant compounds, and electrical stimulation further increased the reliability across wells. These results support the use of multiwell MEA technology for the high-throughput evaluation of complex neuronal networks *in vitro* to evaluate the pro-convulsant risk of candidate compounds.

### References

- Bakum, D. J., Radivojevic, M., Frey, U., Franke, F., Heilmann, A., & Takahashi, H. (2013). Parameters for burst detection. *Frontiers in Computational Neuroscience*, 7, 193. doi:10.3389/fncom.2013.00193
- Chiappalone, M., Novellino, A., Vajda, L., Vato, A., Marinoni, S., & van Pelt, J. (2005). Burst detection algorithms for the analysis of spatio-temporal patterns in cortical networks of neurons. *Neurocomputing*, 65, 466-468. doi:10.1016/j.neurocom.2004.10.094
- Delfanti, E., Novellino, A., Whelan, M., Vogel, S., Ramirez, T., van Raveesteyn, B., & Landisoddi, R. (2013). Feasibility Assessment of Micro-Electrode Chip Assay as a Method of Detecting Neurotoxicity *in vitro*. *Frontiers in NeuroEngineering*, 4(April), 6. doi:10.3389/fneng.2011.00006
- Helliday, D. M., Rosenberg, J. R., Breese, P., & Conway, B. A. (2006). Neural spike train synchronization indices: definitions, interpretations, and applications. *IEEE Transactions on Bio-Medical Engineering*, 53(6), 1056-66. doi:10.1109/TBME.2006.873392
- Kreuz, T., Chicharro, D., Houghton, C., Andrzejak, R. G., & Mormann, F. (2013). Monitoring spike train synchrony. *Journal of Neurophysiology*, 109(5), 1457-72. doi:10.1152/jn.00873.2012
- Leghney, C. R., & Saltman, M. (1983). Bursts and recurrences of bursts in the spike trains of spontaneously active striate cortex neurons. *Journal of Neurophysiology*, 53(4), 926-39. Retrieved from http://www.ncbi.nlm.nih.gov/pubmed/6998798
- McConnell, E.R., et al. Evaluation of multi-well microelectrode arrays for neurotoxicity screening using a chemical training set. *Neurotoxicology*, 2012, 33(5): 1048-57
- Palva, A. R. C., Park, J., & Principe, J. C. (2010). A comparison of binned spike train measures. *Neural Computing and Applications*, 19(3), 405-415. doi:10.1007/s00221-009-0307-6
- Wagenaar, D. A., Pine, J., & Potter, S. M. (2006). An extremely rich repertoire of bursting patterns during the development of cortical cultures. *BMC Neuroscience*, 7, 11. doi:10.1186/1471-2202-7-11

### Acknowledgements

Dr. Tony Krantis – QBM Cell Science / Dr. Chris Strock – Cyprotex / Dr. Khuram Chaudry and Matt Cato – GSK