

Concentration-Response Evaluation of ToxCast Compounds for Multivariate Activity Patterns of Neural Network Function

Marissa B. Kosnik^{1*}, Jenna D. Strickland^{2†}, Skylar W. Marvel¹, Dylan J. Wallis¹, Kathleen Wallace³, Ann M. Richard⁴, David M. Reif¹, and Timothy J. Shafer³

¹Department of Biological Sciences, North Carolina State University, Raleigh, NC, USA, ²Axion Biosystems, Atlanta, GA, USA, ³Integrated Systems Toxicology Division, NHEERL, U.S. Environmental Protection Agency, Research Triangle Park, NC, USA, ⁴National Center for Computational Toxicology, U.S. Environmental Protection Agency, Research Triangle Park, NC, USA

*Current affiliation: Science for Life Laboratory, Department of Environmental Science and Analytical Chemistry, Stockholm University, Stockholm, Sweden

†Current affiliation: Department of Pharmacology and Toxicology, Michigan State University, E. Lansing, MI, USA

Summary

- Microelectrode arrays (MEAs) have been used to assess neuroactivity of ToxCast compounds.
- Screened 384 compounds in concentration response across 43 network activity parameters using MEAs.
- Identified 15 parameters crucial in characterizing neuroactivity of 237 compounds.
- For known neurotoxic compounds, these chemical-parameter potencies were more sensitive than for most ToxCast assays.
- Identified three clusters of chemical-parameter activity with varied bioactivity patterns and chemical structural features.

Introduction

- ToxCast evaluates activity of chemicals using high-throughput screening (HTS).¹
 - Over 1,000 assays.
 - Cannot assess breadth of neurotoxicity targets.
- Microelectrode Arrays (MEAs) can record activity in plated neural networks.
- Previous analysis: 1056 ToxCast chemicals screened at single concentration on MEAs.²
 - 326 altered mean firing rate (MFR).

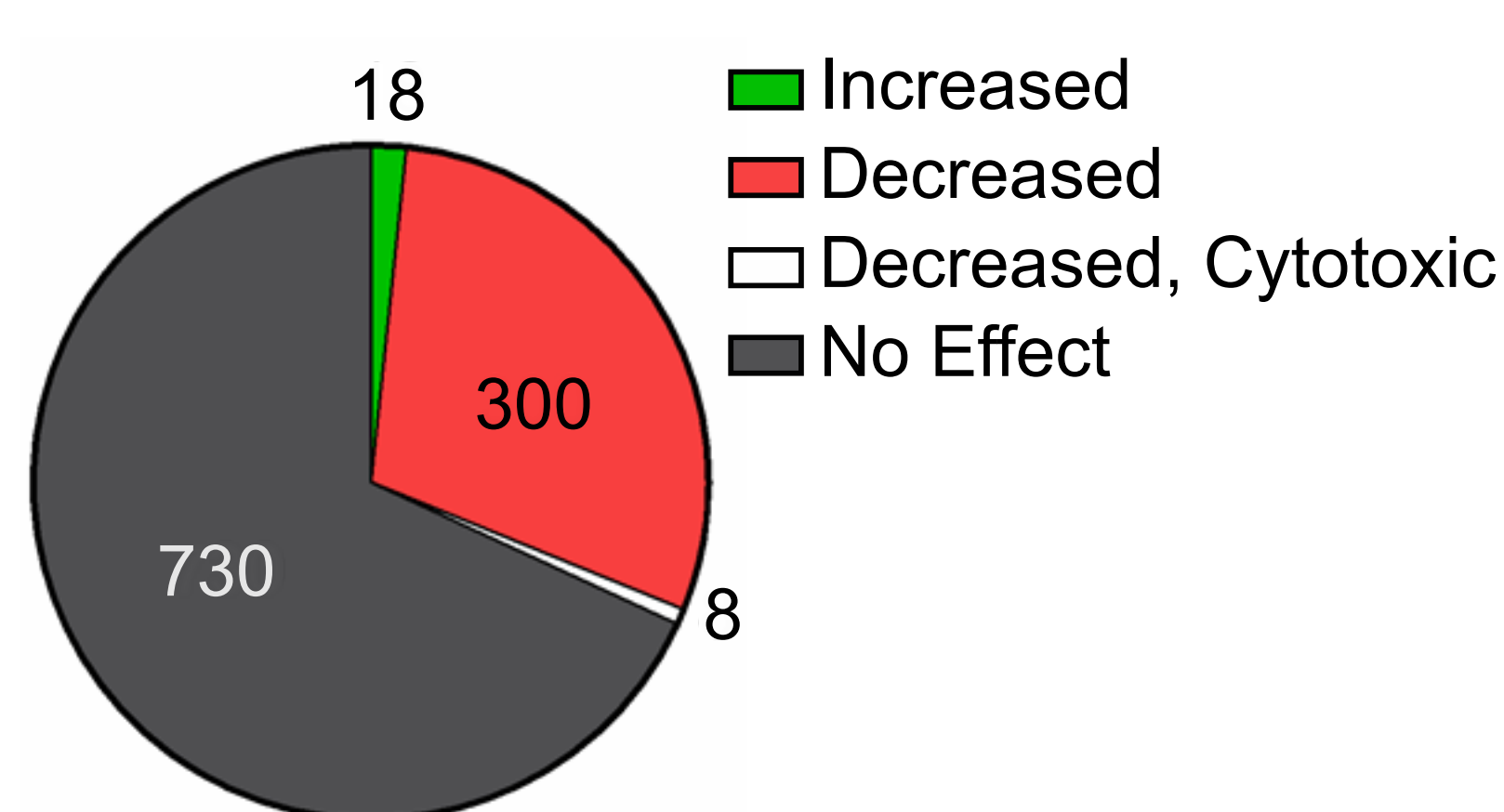


Figure 1. MFR activity in ToxCast compounds.²

Methods

Compound-parameter activity determined with Axion Maestro 1st generation MEA system and classic MEA M768-KAP-48 plates.

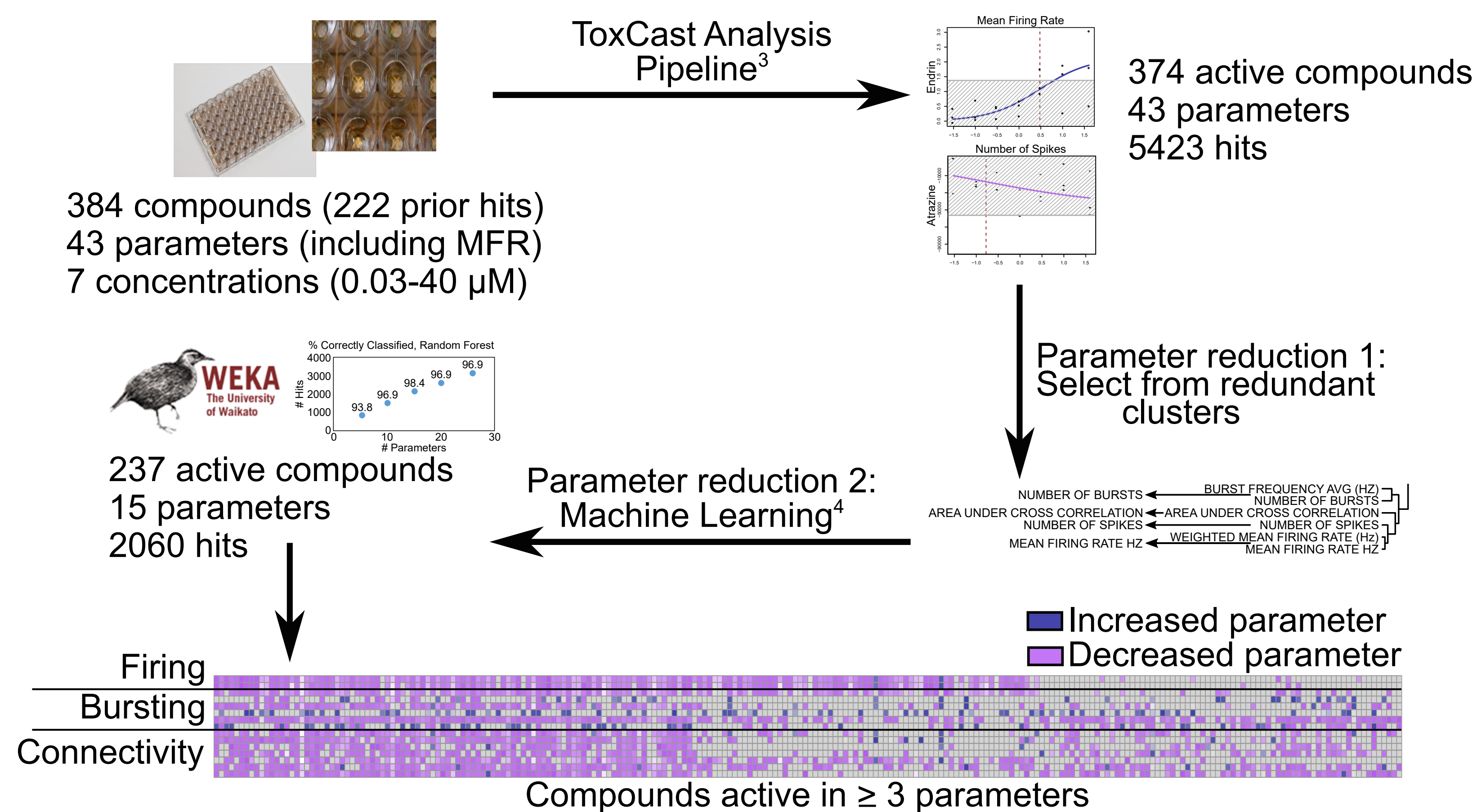


Figure 2. Methods to characterize active compound-parameter associations (heatmap). Darker colors = more potent.

Results

For known neuroactive compounds, MEA parameter potencies are more sensitive than ToxCast assays for most compounds

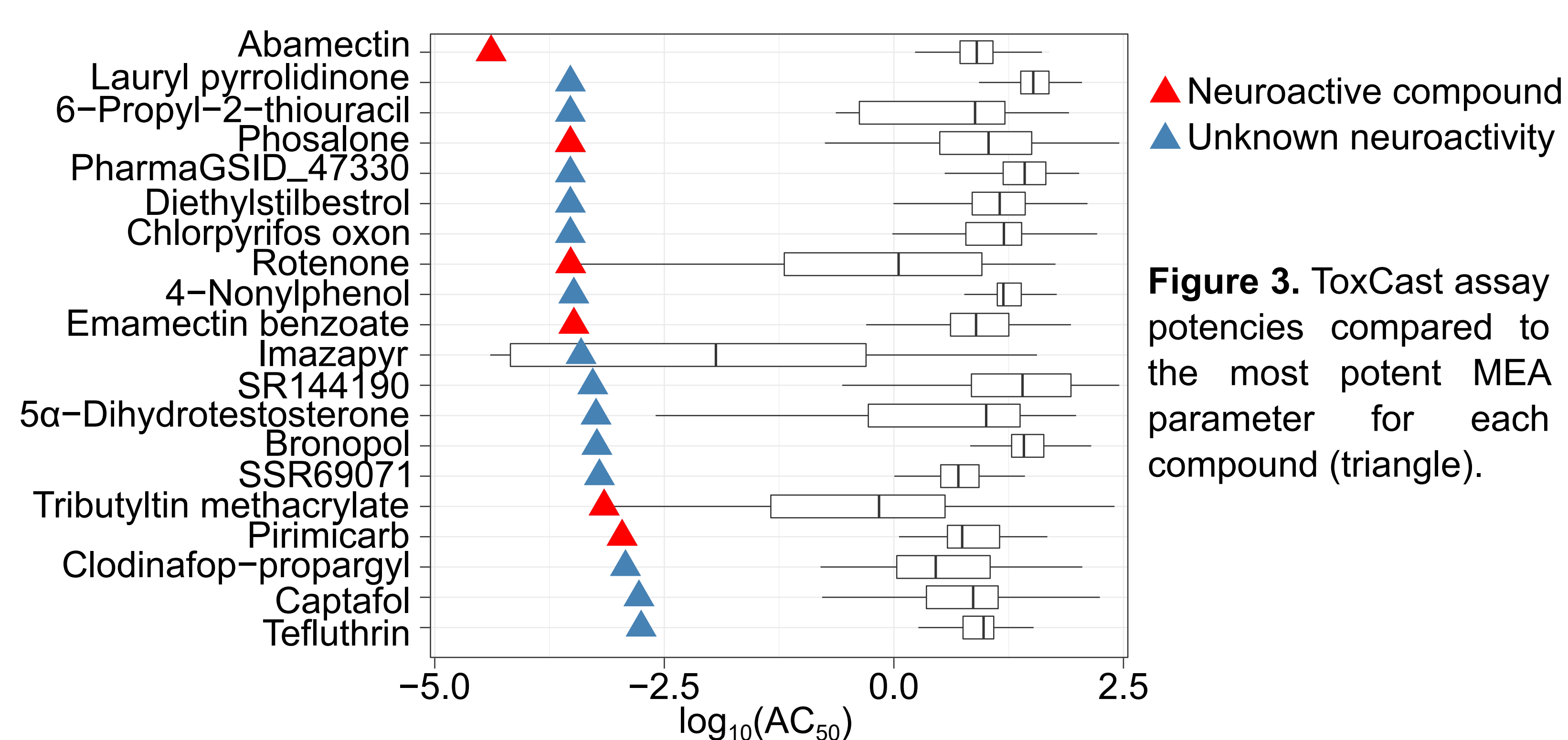


Figure 3. ToxCast assay potencies compared to the most potent MEA parameter for each compound (triangle).

Results

MEA parameters distinguish three clusters of compounds with distinct bioactivity patterns

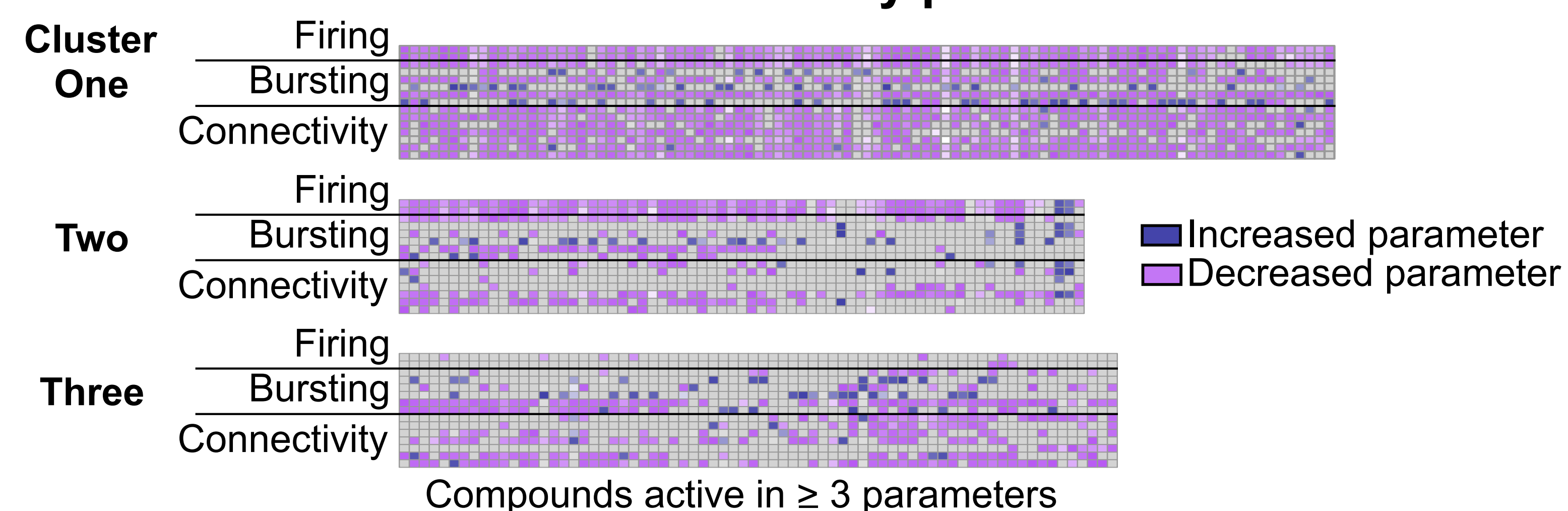


Figure 4. Bioactivity patterns of three MEA clusters determined through k-means clustering of compound-parameter potencies. Darker colors = more potent.

MEA clusters of compounds are enriched for different chemical structural features

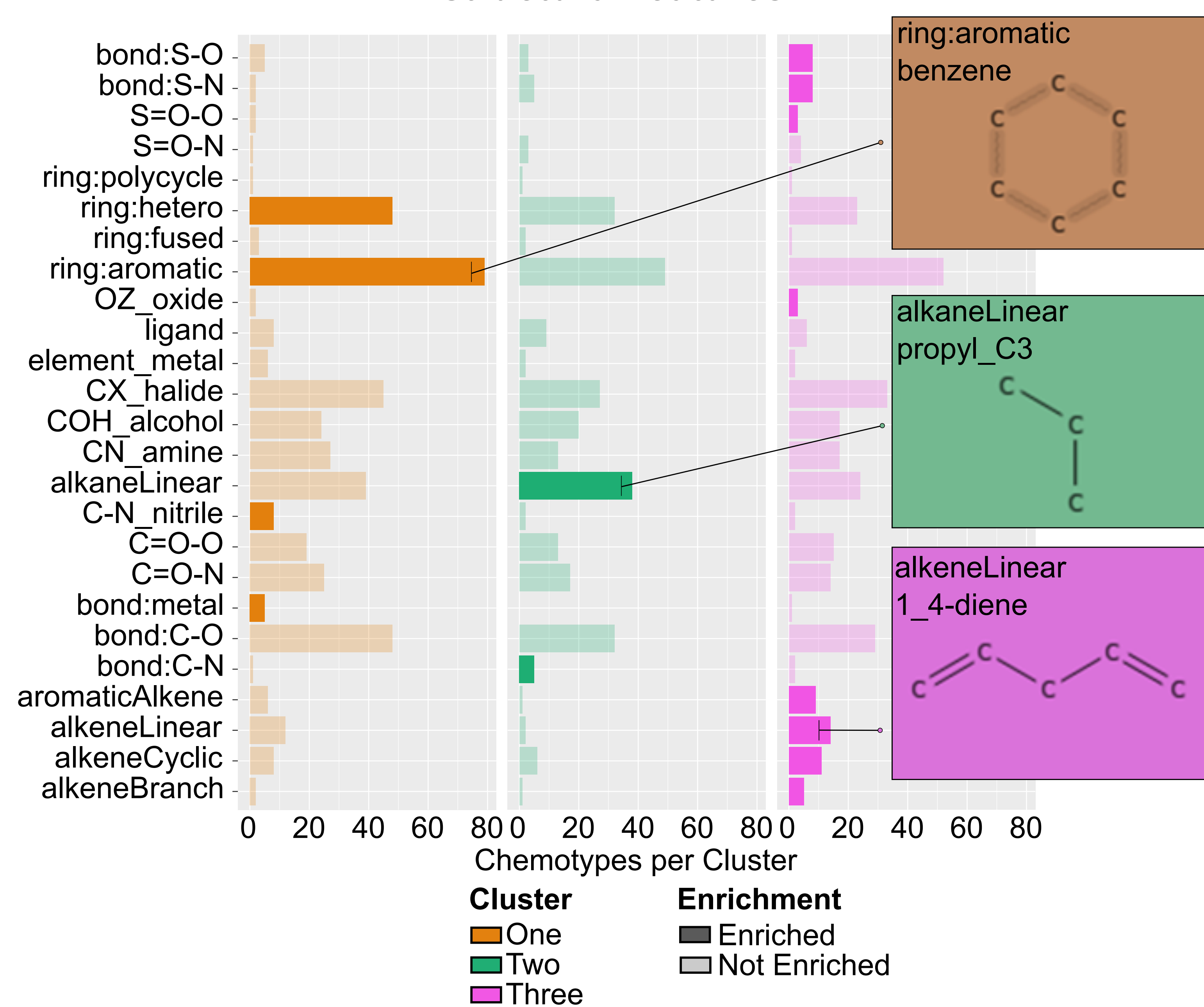


Figure 5. Chemical features (identified as chemotypes⁵) enriched within each cluster using hypergeometric test. Inset images depict example structures within respective chemotype.

Conclusions

- We developed a robust assessment for neuroactivity across 15 network parameters (Figure 2).
- For known neuroactive compounds, MEA parameters were more sensitive overall than other ToxCast assays (Figure 3).
- MEA chemical-parameter potencies can be used to group compounds based on bioactivity (Figure 4).
- Biological activity groups were consistent with the underlying structure of the compounds (Figure 5).
- Multivariate MEA activity patterns can efficiently screen for diverse chemical neuroactivities.
- MEA activity patterns may have predictive value related to chemical structural features.

References & Funding

This work was supported by EPA CRADA 644-11, National Institutes of Health [ES025128, ES030007] and the US Environmental Protection Agency [STAR R835802] This poster does not reflect EPA policy.

- Dix, D.J. et al. (2007) The toxcast program for prioritizing toxicity testing of environmental chemicals. *Toxicol. Sci.*
- Strickland, J.D. et al. (2018) Screening the ToxCast phase II libraries for alterations in network function using cortical neurons grown on multi-well microelectrode array (mwMEA) plates. *Arch. Toxicol.*
- Filer, D.L. et al. (2017) Tcpl: The ToxCast pipeline for high-throughput screening data. *Bioinformatics*
- Frank, E. et al. (2016) The WEKA Workbench. Online Appendix for "Data Mining: Practical Machine Learning Tools and Techniques"
- Chemotyper application (<https://chemotyper.org/>), ToxPrint feature set (<https://toxprint.org/>)