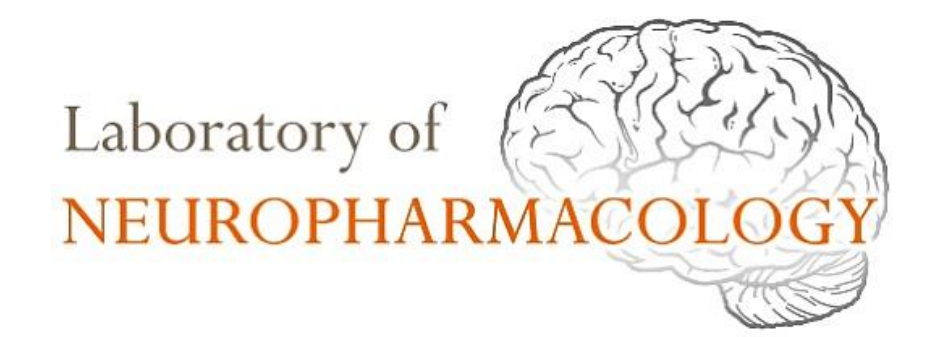




# Properties of Human Stem Cell-derived Neurons in Long-term Cell Culture as a Model for Antiseizure Drug Discovery

Hamed Salmanzadeh and Robert F. Halliwell

Thomas J. Long School of Pharmacy, University of the Pacific, Stockton, CA



## Introduction

Studies in our lab have addressed the functional properties of receptors and ion channels expressed in neurons derived from human stem cells (e.g., Halliwell *et al.*, 2021). Here we have investigated a pharmacologically diverse selection of antiepileptic agents that act *via* distinct receptors and ion channels on spontaneous spiking and induced epileptiform activity to determine their sensitivity, reliability and validity for drug investigations.

## Methods

**Cell culture.** The human stem cell line TERA2.cl.SP12 was cultured under differentiating conditions towards a neuronal phenotype using retinoic acid (10 $\mu$ M).

**Immunocytochemistry.** Stem cell-derived neurons were immunolabeled with the neural markers  $\beta$ III-tubulin, and MAP2.

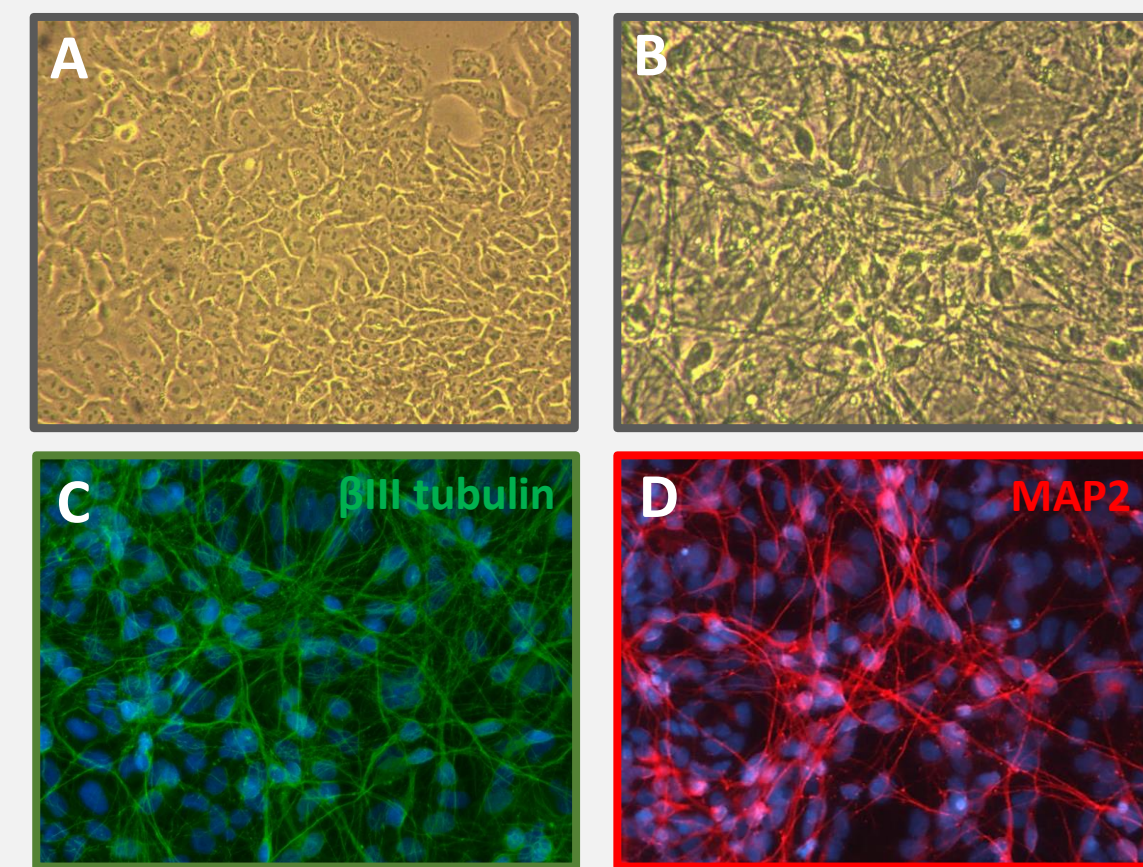
**Multi-Electrode Array Electrophysiology.** TERA2.cl.SP12 stem cells were cultured in 6-well MEA plates and allowed to differentiate. Activity of differentiated cells was recorded weekly for 1 year using the Maestro Edge MEA system (Axion Biosystems).



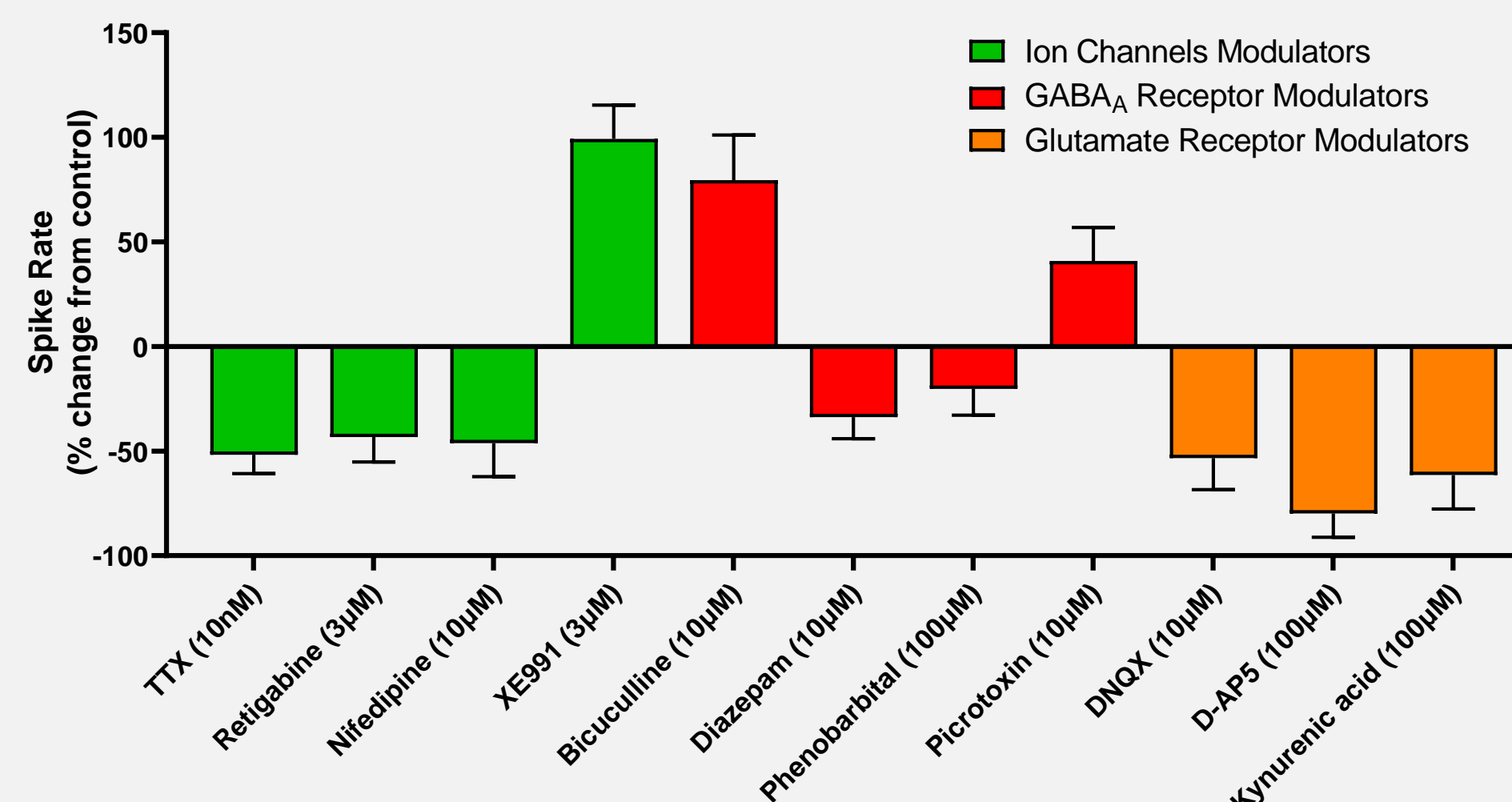
Human stem cell derived neurons differentiated for 120 days *in vitro* on a 64 electrode MEA well.

## Conclusion

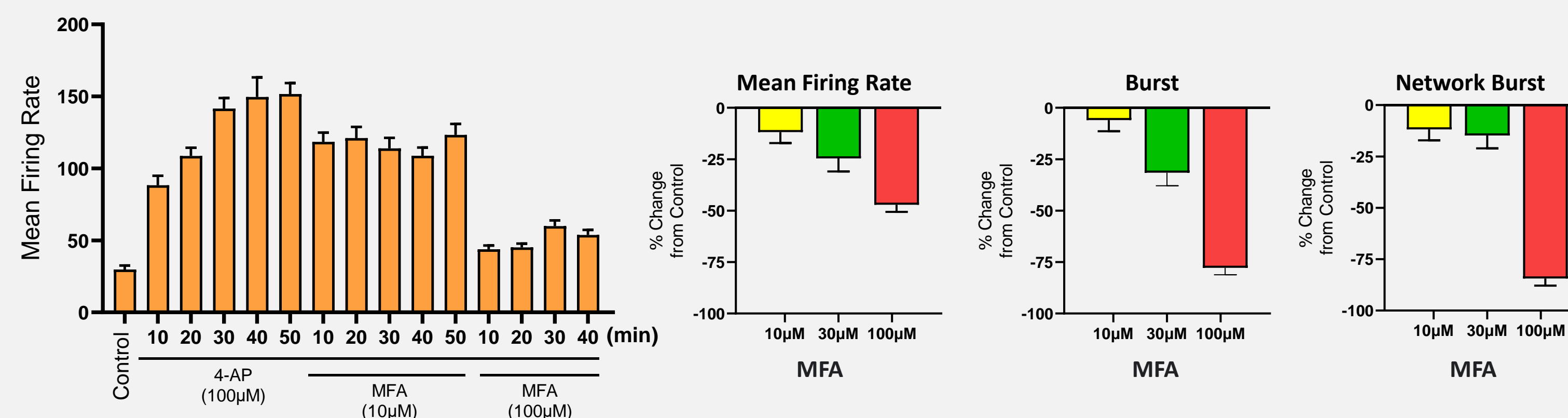
2D cultures of human stem cell-derived neurons are now a valuable tool in drug discovery. Here, we addressed the development and maturation of the endogenous electrochemical and pharmacological properties of neurons differentiated over 1 year *in vitro*. Our data show that neurons develop an array of voltage and ligand- ion channels and synaptic neurotransmission is mediated *via* the major inhibitory and excitatory receptors. We also demonstrate that these neural networks are sensitive to established and potentially novel (e.g. MFA) anticonvulsants, supporting their real value in neuropharmacological investigations.



**Figure 1.** Retinoic acid (RA) promoted neuronal differentiation of TERA2.cl.SP-12. (A) and (B) are phase contrast images showing cells at day 0 (A) and day 80 (B) of differentiation. Immunofluorescence images show fixed stem cell-derived neurons labeled with (C)  $\beta$ -III tubulin (green) and (D) MAP2 (red) at 80 days differentiation.

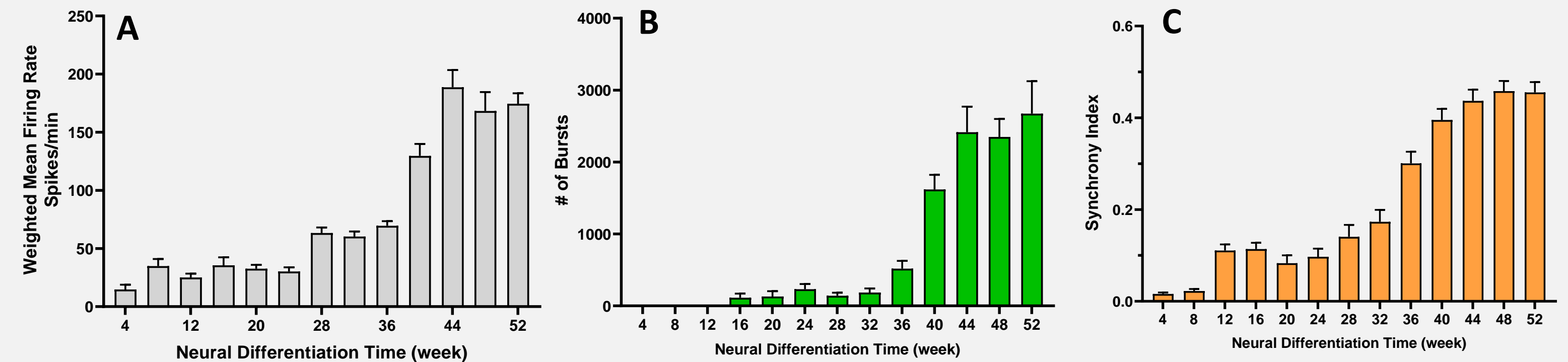


**Figure 3.** Pharmacological profile of spontaneous firing recorded from stem-cell derived neurons. Addition of ion channel and receptor modulators to the media showed that neurons expressed voltage- and ligand- gated ion channels and that synaptic transmission was mediated *via* glutamate and GABA<sub>A</sub> receptors in these neural networks.

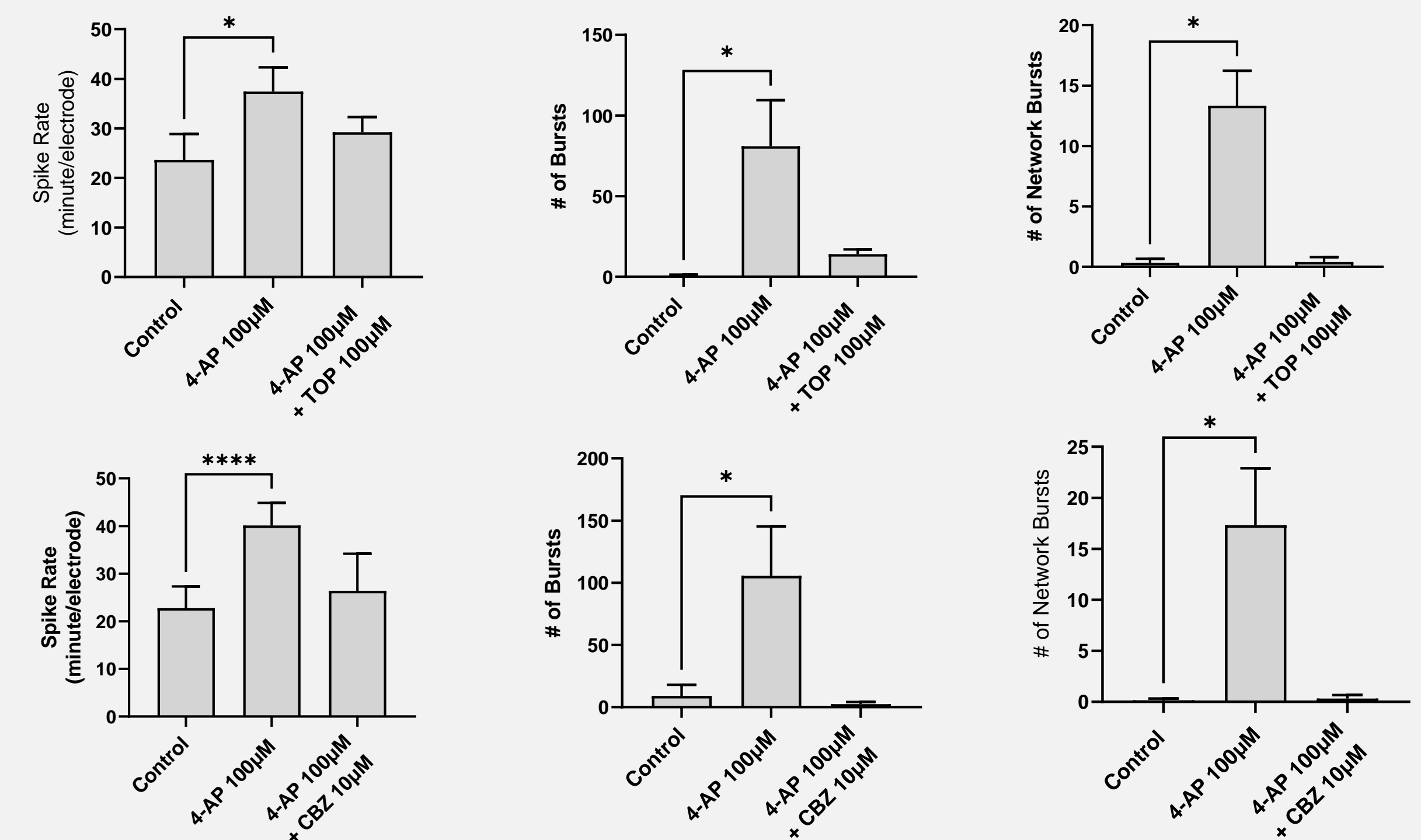


**Figure 5.** The NSAID, mefenamic acid (MFA) is also a GABA<sub>A</sub> receptor modulator. Here, we tested its potential value as a novel anticonvulsant against 4-AP induced epileptiform activity. The histograms (left) show that MFA reduced unit firing, bursting and network burst firing evoked by 4-AP in a highly concentration-dependent manner. The cells were 16 weeks differentiation.

## Results



**Figure 2.** Development and maturation of spontaneous firing of stem cell-derived neurons over one year of differentiation. (A) Time course of the average firing rate from 4 to 52 weeks differentiation and (B) development of spontaneous burst and (C) synchrony firing during long-term culture.



**Figure 4.** Impact of the anticonvulsants, topiramate (TOP) and carbamazepine (CBZ) on epileptiform activity induced by 4-aminopyridine (4-AP). The histograms show that 4-AP significantly increased spiking, burst firing, network bursts and synchronized epileptiform activity. Topiramate and carbamazepine reduced neural activity to control levels. The cells were 16 weeks differentiation.