

# Modulation of GABA<sub>A</sub> activity: Investigations in hiPSC-derived neural co-cultures and human ion channel assays

K. Rockley<sup>1</sup>, B. Kelly<sup>1</sup>, K. Jones<sup>1</sup>, R. Roberts<sup>1</sup>, M. Morton<sup>1</sup>

<sup>1</sup> Apconix, Alderley Edge, Cheshire, United Kingdom

## INTRODUCTION

A balance between inhibitory neurotransmission and neuronal excitation is critical for normal brain function.  $\gamma$ -aminobutyric acid (GABA) is the most abundant inhibitory neurotransmitter, which acts on GABA<sub>A</sub> receptors. Perturbation of GABA<sub>A</sub> signalling by drug-induced inhibition and potentiation are common mechanisms producing seizure and sedation, respectively. The introduction of commercially available human induced pluripotent stem cell (hiPSC-) derived neurons facilitates the *in vitro* study of neuronal function and, in our work, the detection of seizure liability during drug discovery. It is known that GABA<sub>A</sub> antagonists such as picrotoxin increase neuronal firing and induce a seizure-like phenotype in hiPSC-derived neurons, however further characterisation of GABA<sub>A</sub> modulation within these cell models is lacking. This study aimed to address this by assessing the effects of a selection of GABA modulators on the electrical activity of hiPSC-derived neuronal co-cultures, and the ion flux of  $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub>.

## METHODS

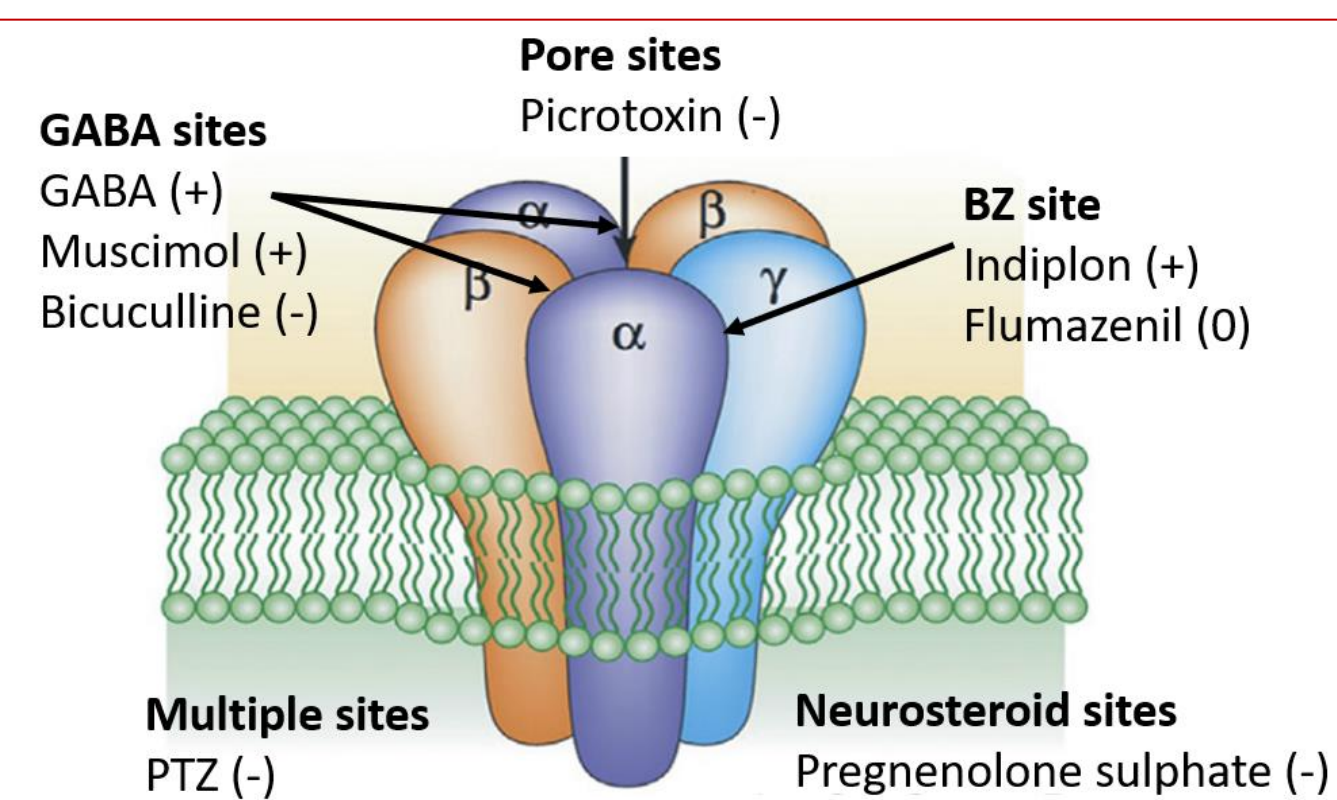
### hiPSC-DERIVED NEURONAL CO-CULTURES

- iCell Glutaneurons (80% glutamatergic/20% GABAergic neurons) were plated with astrocytes (85%:15%) and monitored using a microelectrode array (MEA) system (Maestro Edge, Axion).
- On DIV22 and DIV23, spontaneous electrical activity was recorded at baseline and 1 hour after exposure to GABA<sub>A</sub> modulators and solvent controls.
- Cells exposed to agonists were subsequently challenged with antagonists and spontaneous electrical activity was measured 15 minutes after application.

### HUMAN $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub> ION CHANNEL ASSAYS

- The activity of GABA modulators was assessed by automated patch-clamp (QPatch II, Sophion) using a CHO  $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub> cell line.
- All modulators except for ligands were co-applied with 30 $\mu$ M GABA.
- 6-point dose-response curves were generated for all modulators.
- For agonists, a 5-point dose-response curve plus subsequent antagonist challenge was generated.

### COMPOUND SELECTION

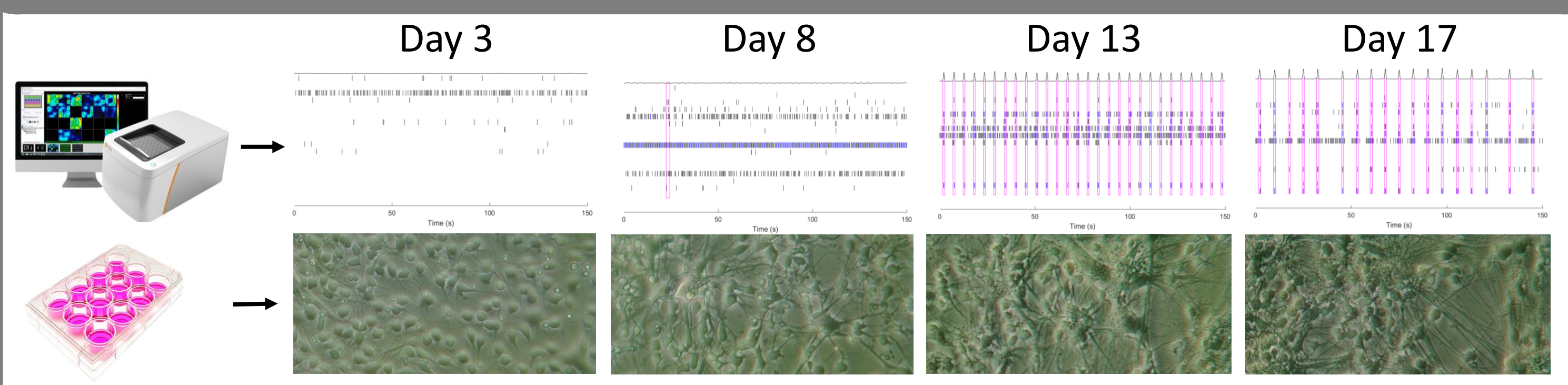


## RESULTS

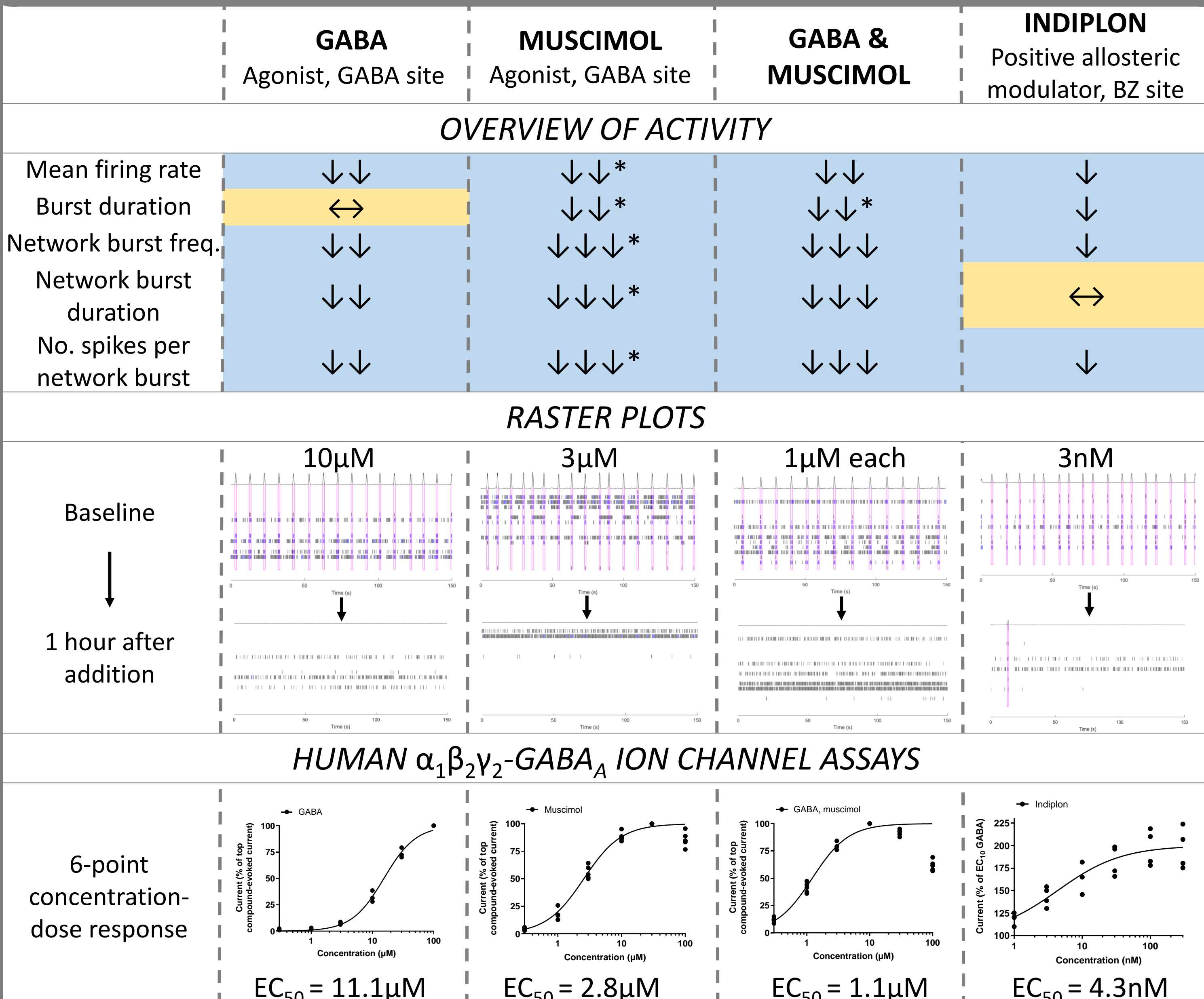
### MEA PARAMETERS

<b>Firing rate</b> - Weighted mean firing rate based on electrodes with activity greater than minimum spike rate, set by the neural statistics calculator.	↑↑↑ ≥100% ↑↑ 50 to 99% ↑ 20 to 50%
<b>Burst duration</b> - Average time between the first and last spike in a burst.	↔ within +/-20%
<b>Network burst freq.</b> - Total number of electrode bursts divided by recording time.	↓↓ -20 to -50% ↓↓↓ -50 to -99%
<b>Network burst duration</b> - Average time between the first and last spike in a network burst.	↓↓↓ ≥100% ***p<0.001 **p<0.01 *p<0.05
<b>No. spikes per network burst</b> - Average number of spikes in a network burst.	

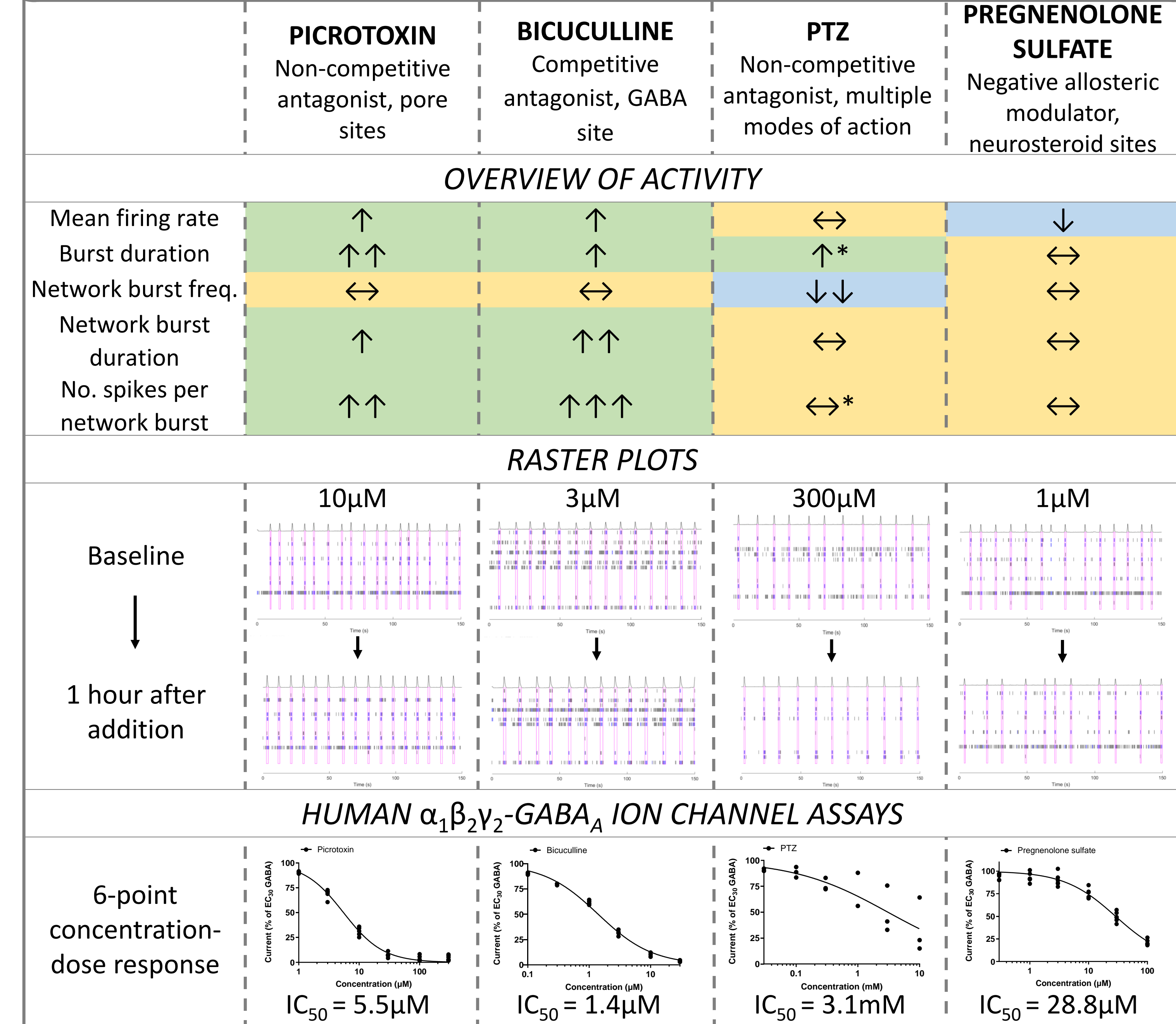
### DEVELOPMENT OF SPONTANEOUS ELECTRICAL ACTIVITY



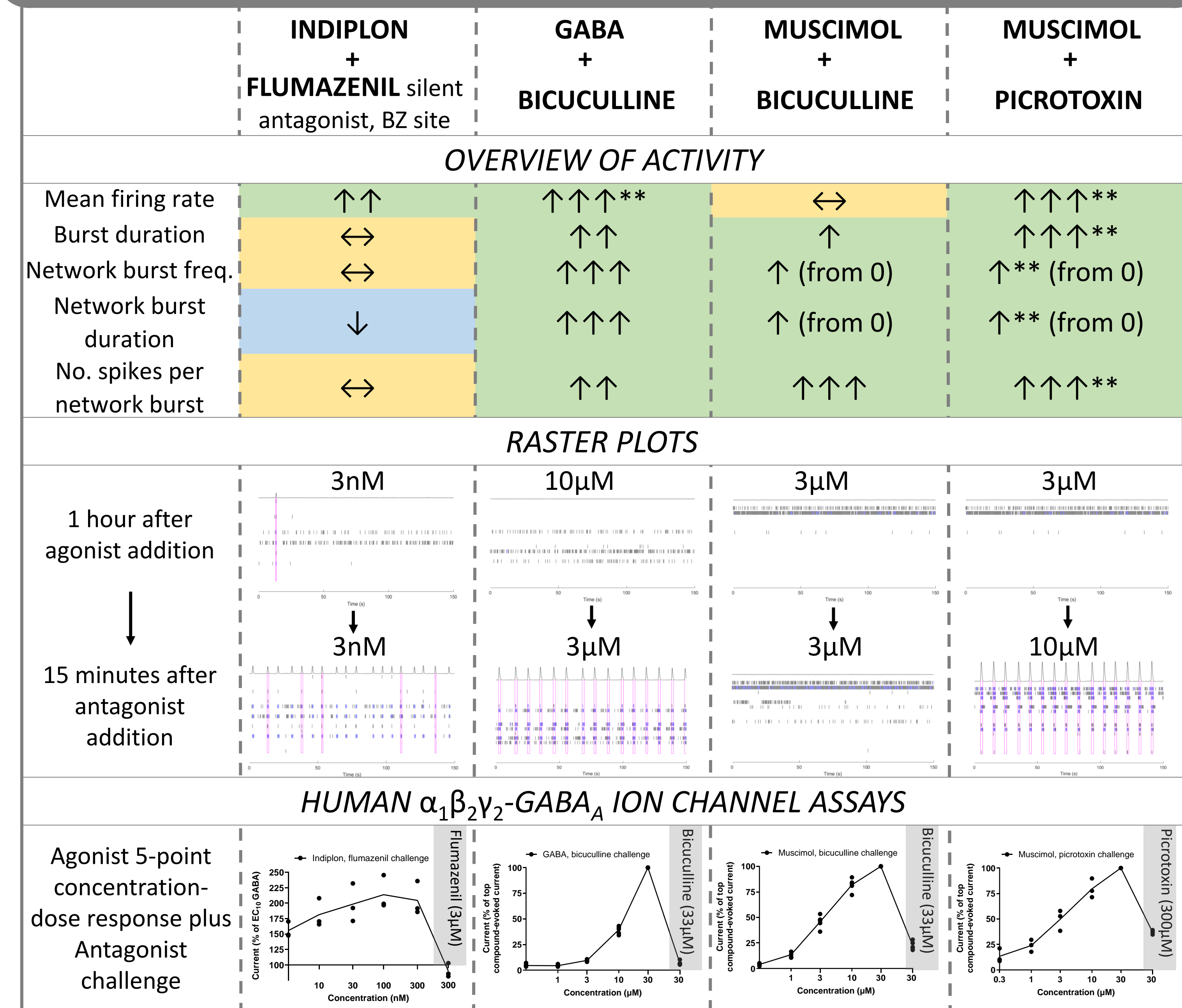
## 1 AGONISTS INCREASE GABA<sub>A</sub> RESPONSE, DECREASE POPULATION ACTIVITY



## 2 ANTAGONISTS DECREASE GABA<sub>A</sub> RESPONSE, MIXED POPULATION ACTIVITY



## 3 REVERSAL OF AGONIST-INDUCED SEDATION BY ANTAGONISTS



## DISCUSSION AND CONCLUSIONS

- Agonists GABA and muscimol induced a sedation like phenotype in hiPSC-derived neuronal co-cultures and increased  $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub> current (fig.1) while antagonists bicuculline and picrotoxin induced a seizure like phenotype in hiPSC-neuronal co-cultures and reduced  $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub> current (fig.2).
- PTZ is used *in vivo* to induce seizure. This often involves chronic repeat-dose application, suggesting PTZ may not translate well to single-dose *in vitro* studies (fig.2).
- Pregnenolone sulfate (PS) did not induce seizure in hiPSC-derived neuronal co-cultures, yet inhibited  $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub> current. This suggests the expression of other subtypes in neuronal cells, possibly GABA<sub>C</sub> which is considerably less sensitive to PS than GABA<sub>A</sub>.
- In ion channel assays, bicuculline blocked GABA- and muscimol-induced current (fig.3). In hiPSC-derived neuronal co-cultures however, muscimol-induced sedation was not reversed by bicuculline. It is known that bicuculline cannot compete with muscimol at GABA<sub>C</sub>, further suggesting its expression in hiPSC-derived neuronal co-cultures.
- Indiplon, a marketed sleeping aid, induced sedation in hiPSC-derived neuronal co-cultures and increased  $\alpha_1\beta_2\gamma_2$ -GABA<sub>A</sub> current (fig.3). It was competitively antagonized by flumazenil, a clinical antidote to indiplon overdose. As a silent antagonist, Flumazenil was inactive alone in both assays.
- These studies have further characterised modulation of GABA<sub>A</sub> activity within hiPSC-derived neuronal co-cultures by recapitulating expected clinical outcomes. This further validates the model as a translationally relevant screen for seizure detection which also shows promise for sedation.