

# Functional assessment of spontaneous and evoked activity in iPSC-derived Fragile X neurons

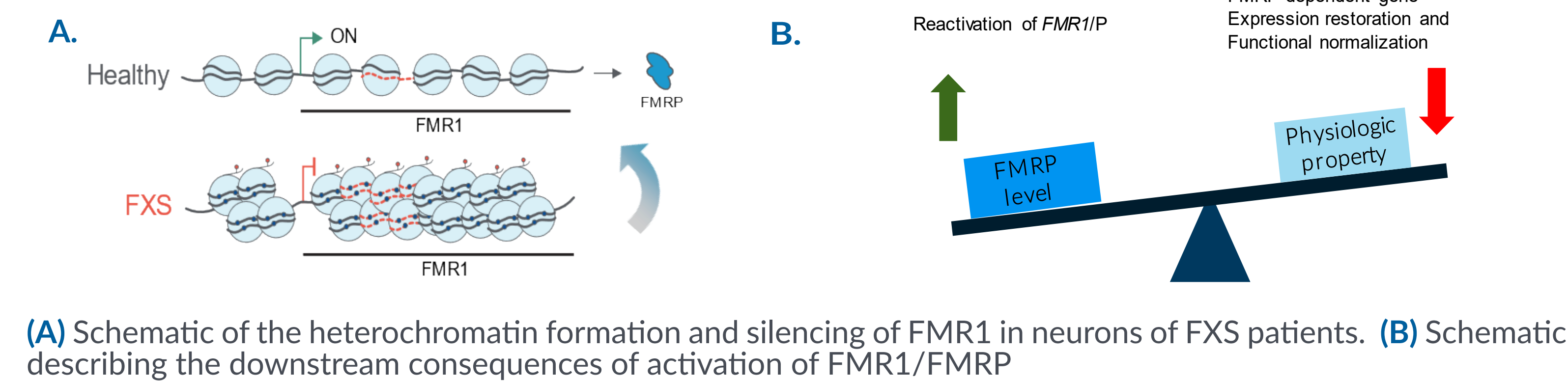
John D Graef, Wu H, Sun C, Lin L, Webb S, Villegas V, Ng C, Warren S\*, Cacace AM.  
Fulcrum Therapeutics. One Kendall Square, Suite B7102, Cambridge, Massachusetts, USA; \*Emory University. jgraef@fulcrumtx.com



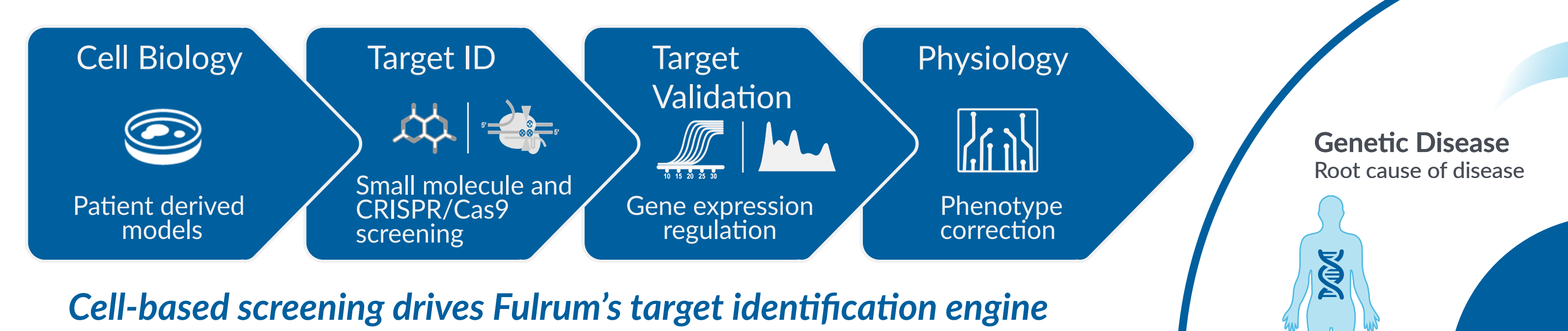
## ABSTRACT

- Fragile X Syndrome (FXS) is a neurodevelopmental disorder caused by >200 repeat CGG expansion in the 5'-UTR of the Fragile X mental retardation gene *FMR1*, leading to hypermethylation and silencing of the *FMR1* transcript, thereby reducing FMRP protein.
- To date, there has been no quantitative assessment of the relationship between varying levels of *FMR1* / FMRP expression and effects on spontaneous neuronal activity.
- We have developed patient-specific iPSC-derived neurons that retain the pathological repeat expansion, have decreased FMRP levels, and demonstrate increased spontaneous activity in *in vitro* neuronal networks as measured by multielectrode arrays (MEAs).
- Using CRISPR/Cas9, we have also generated a *FMR1* knockout line, as well as isogenic Fragile X cell lines with truncated numbers of CGG repeats that demonstrate increased spontaneous activity in the absence of FMRP and attenuated activity when FMRP is restored to near normal levels.
- We further show through *FMR1* mRNA transient transfections and titration of isogenic control neurons into a Fragile X neuronal network, that low levels of FMRP (<10%) partially attenuate FMRP-induced hyperexcitability, with higher levels of FMRP (>10%) leading to full normalization of increased levels of spontaneous activity seen in Fragile X neuronal networks.

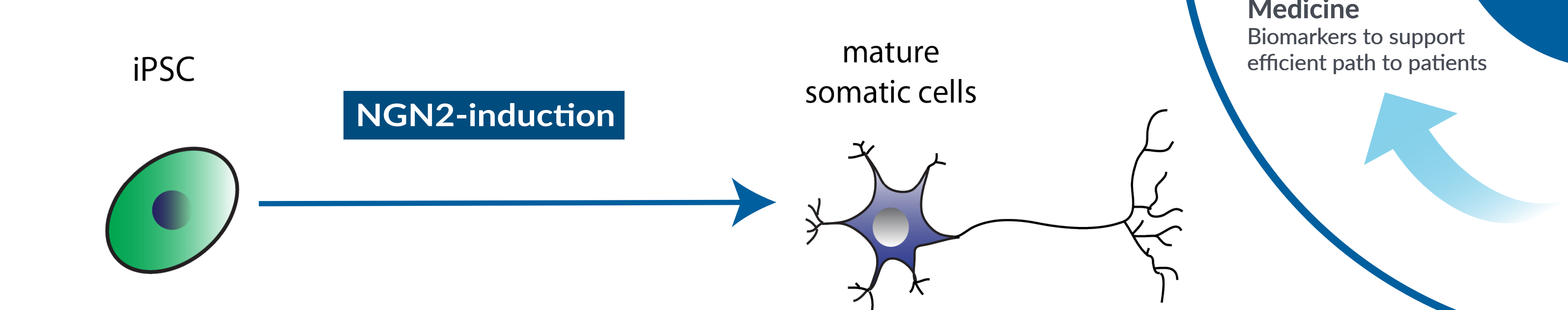
## 1. FRAGILE X SYNDROME (FXS)



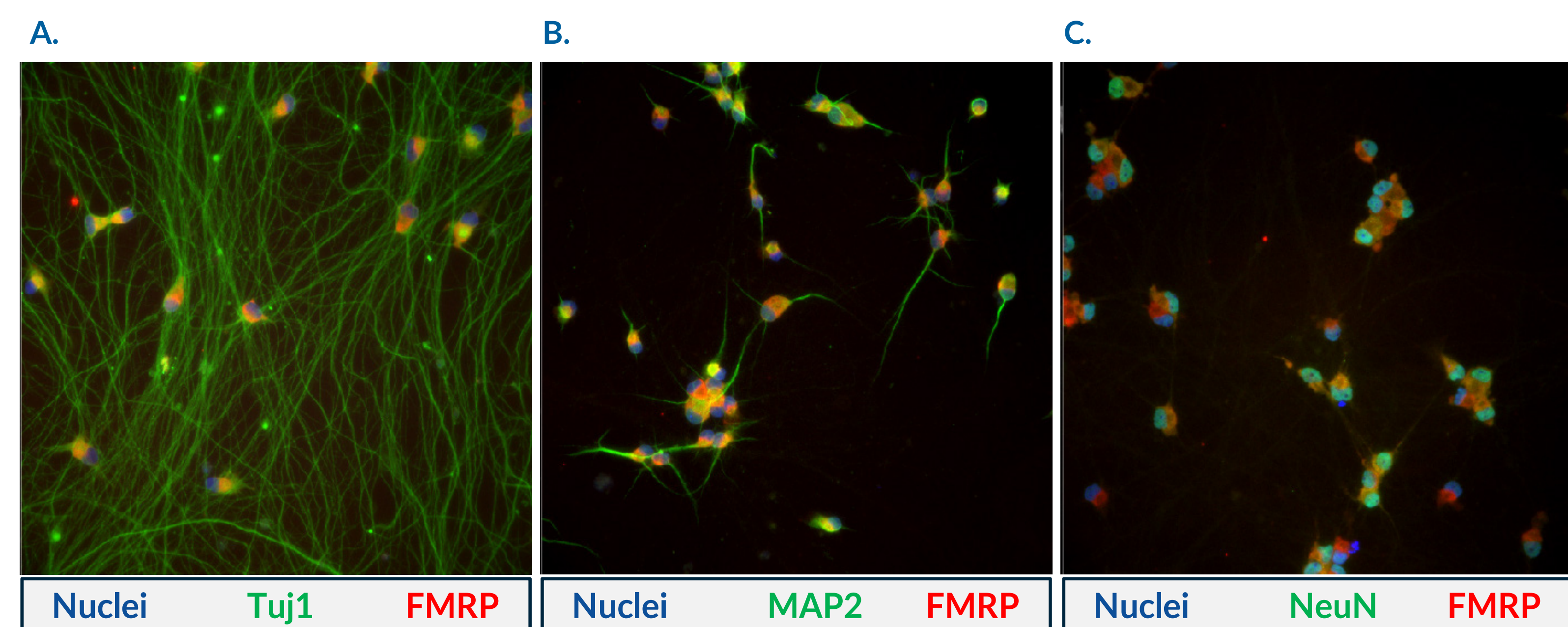
## 2. FULCRUM'S APPROACH TO TARGET ID AND VALIDATION



## 3. HUMAN IPSC-DERIVED NEURONS

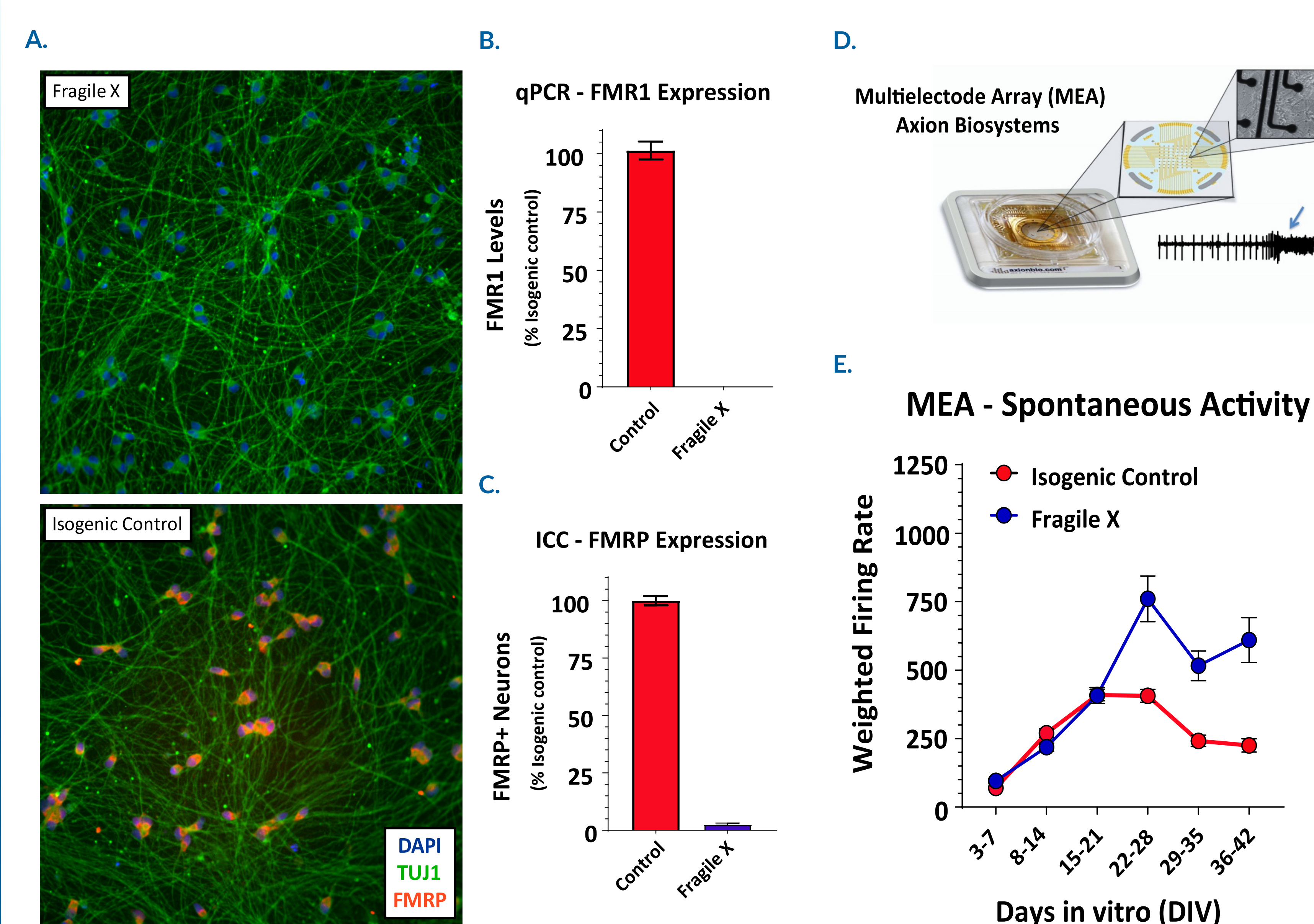


Patient-specific iPSCs are differentiated into excitatory glutamatergic neurons using a modified protocol adapted from Zhang et al. (2013), "Rapid Single-Step Induction of Functional Neurons from Human Pluripotent Stem Cells", *Neuron*.



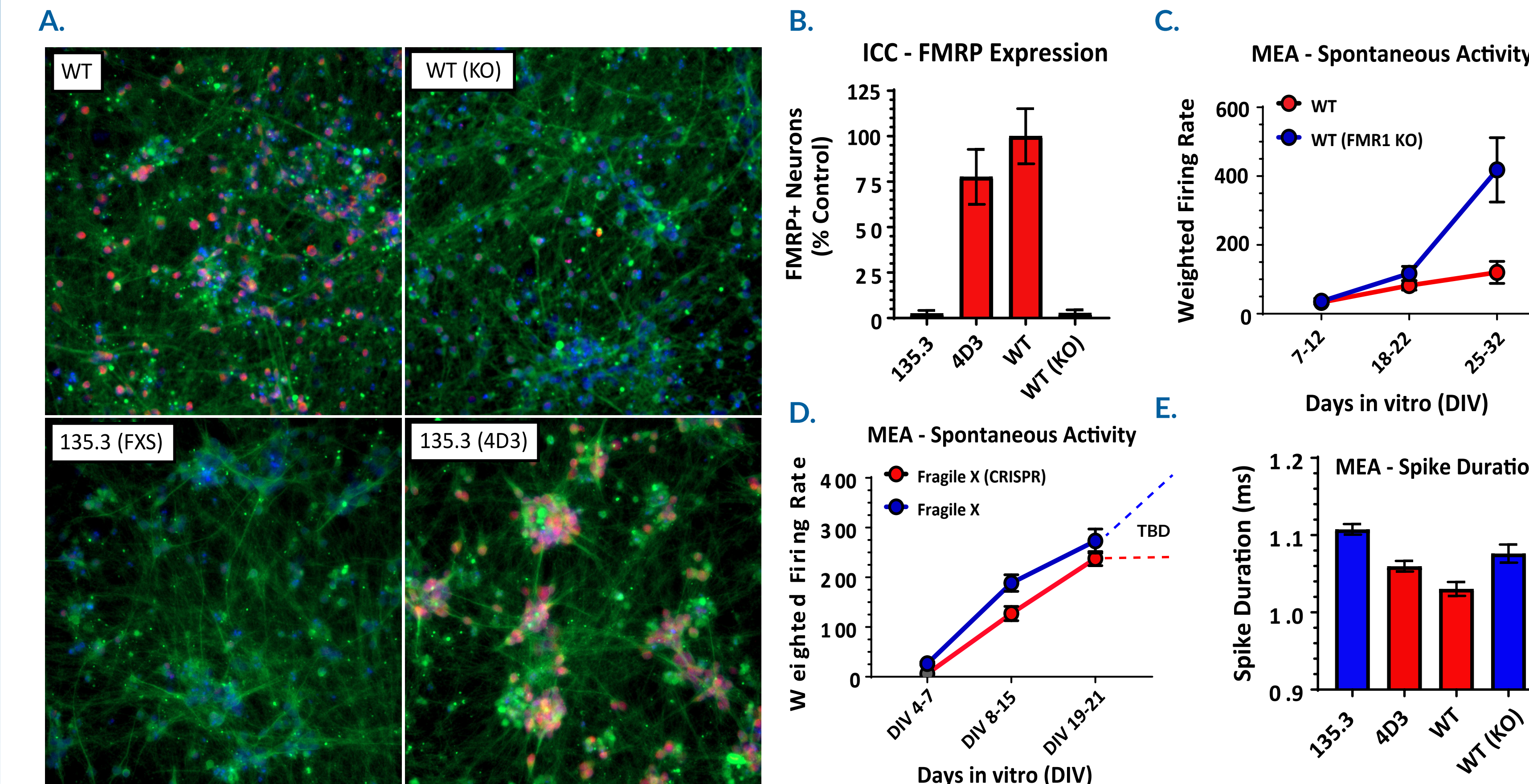
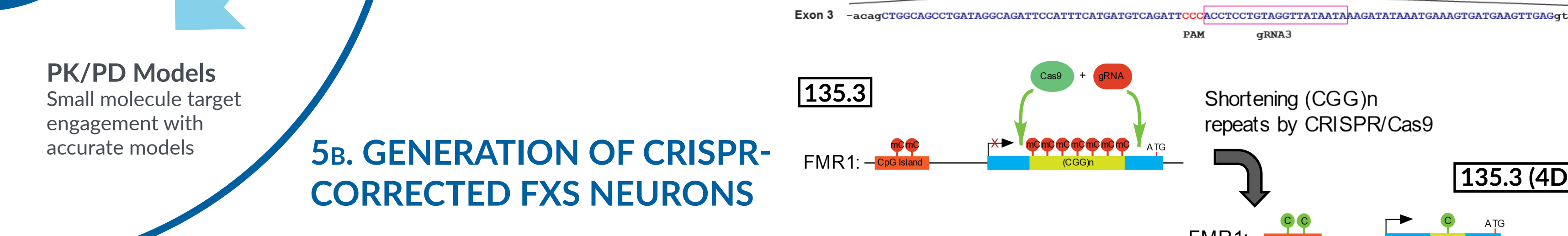
Immunofluorescent images from WT human iPSC-derived neurons. Cells were fixed and stained after 4 weeks in culture. Nuclei (blue) are stained with DAPI, neuron-specific markers (green) are stained using the following antibodies: (A) Tuj1, Aves, TUJ89947984; (B) MAP2, Aves, MAP717984; and (C) NeuN, Abcam, ab177487. FMRP (red) is stained using the 5C2 antibody from Biologend. As can be seen from the images, FMRP can only be detected in the soma and not in the neurites.

## 4. FXS NEURONS HAVE REDUCED FMR1/FMRP AND INCREASED ACTIVITY



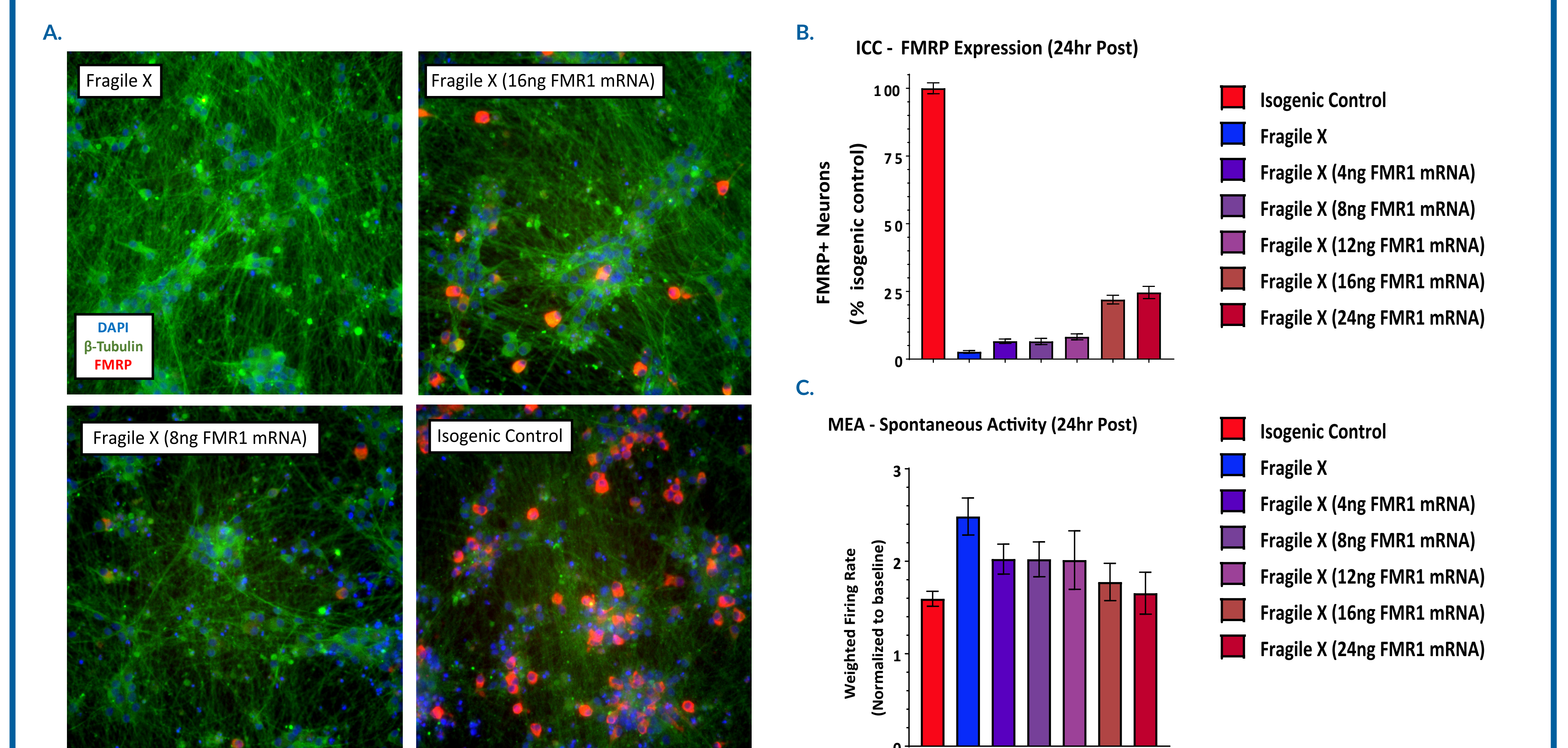
(A) Immunofluorescent images of iPSC-derived neurons from a Fragile X cell line (SW\_FXS) and an isogenic control line (SW\_C1\_2). Neurons were fixed 4 weeks post-induction. Nuclei (blue) are stained using DAPI, neurites (green) are stained with a Tuj1 antibody (Aves), and FMRP (red) is stained using a FMRP antibody (Biologend, 5C2). Quantification of *FMR1* mRNA levels are shown in panel B. Ct values are normalized to housekeeping gene *POP4*, then normalized to the mean of the control values. Quantification of the number of FMRP+ neurons are shown in panel C and normalized to the mean of the control values. Example image of a multielectrode array with an image of neurons plated directly onto the electrode array along with an example recording of spontaneous activity. (E) Quantification of the weighted firing rate (average spike rate within a well multiplied by the number of active electrodes in the well) for both Fragile X and isogenic control neurons over a period of 6 weeks. As can be seen from the graph, Fragile X neurons begin exhibiting hyperactivity starting 3 weeks post differentiation. Data represented as mean +/- SEM.

## 5A. GENERATION OF FMRP KO NEURONS

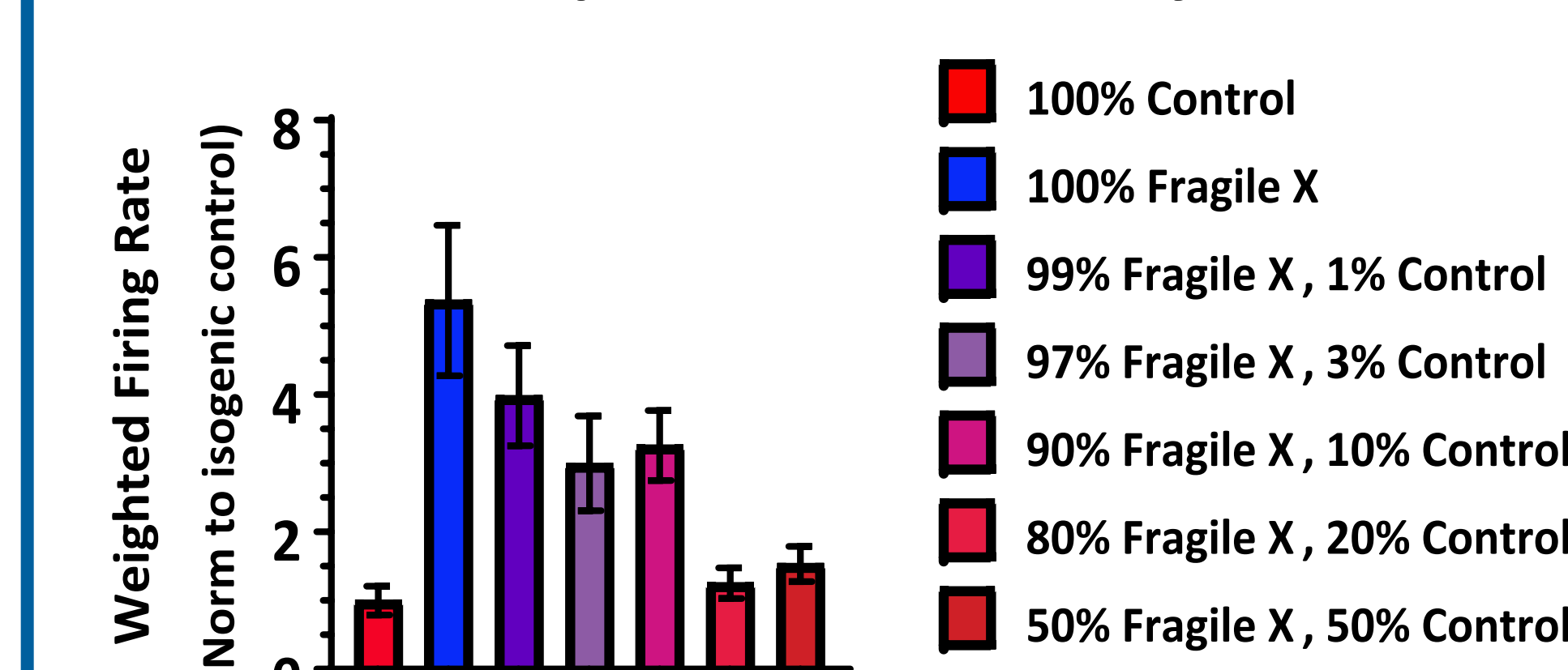


(A) Immunofluorescent images of iPSC-derived neurons from a healthy-normal line (WT) and a Fragile X line (135.3) where FMRP has been removed from the WT line through a premature stop codon in Exon 3 of *FMR1* (WT KO) and FMRP has been restored in the Fragile X line by removal of CGG repeats using CRISPR/Cas9 (135.3, clone 4D3). Quantification of the number of FMRP+ neurons are shown in panel B. (C) Quantification of spontaneous activity for WT and KO neurons over 4 weeks in culture showing increased activity in neurons lacking FMRP. (D) Quantification of spontaneous activity for both Fragile X and CRISPR-corrected neurons over 3 weeks in culture showing decreased activity in neurons re-expressing FMRP. (E) Average spike duration of spontaneous spikes for all four lines. Data represented as mean +/- SEM.

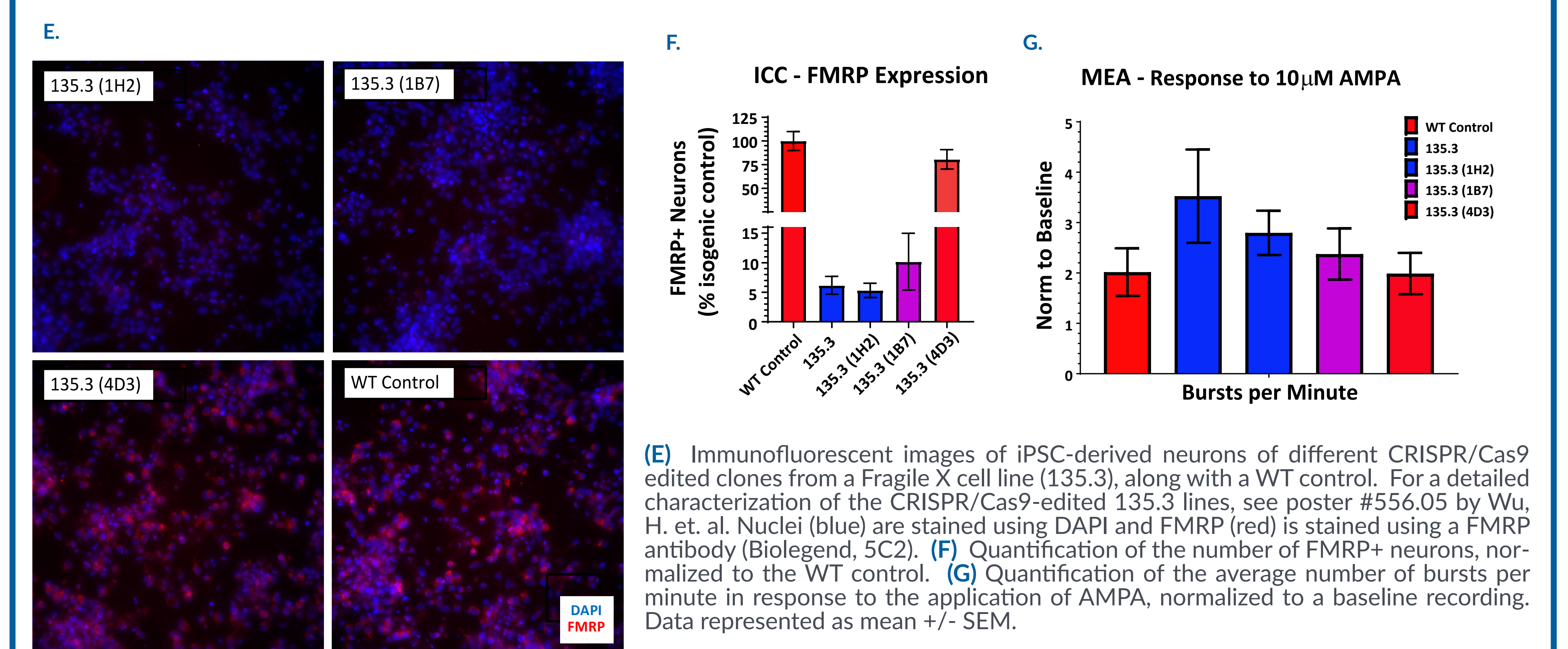
## 6. HOW MUCH FMRP IS NEEDED TO NORMALIZE ACTIVITY?



## MEAs - Spontaneous Activity



## DIFFERENT CRISPR/CAS9-EDITED CLONES RESULT IN DIFFERENT FMRP LEVELS



## CONCLUSIONS

- We have developed patient-specific iPSC-derived neurons that retain the pathological repeat expansion, have decreased FMRP levels, and demonstrate increased spontaneous activity in *in vitro* neural networks as measured by multielectrode arrays.
- Using CRISPR/Cas9, we have also generated a *FMR1* knockout line, as well as a Fragile X cell line with a truncated number of CGG repeats that demonstrates increased spontaneous activity in the absence of FMRP and attenuated activity when FMRP is restored to near normal levels.
- We further show through *FMR1* mRNA transient transfections and titration of isogenic control neurons into a Fragile X neuronal network, that low levels of FMRP (<10%) partially attenuate FMRP-induced hyperexcitability, with higher levels of FMRP (>10%) leading to full normalization of increased levels of spontaneous activity seen in Fragile X neuronal networks.