

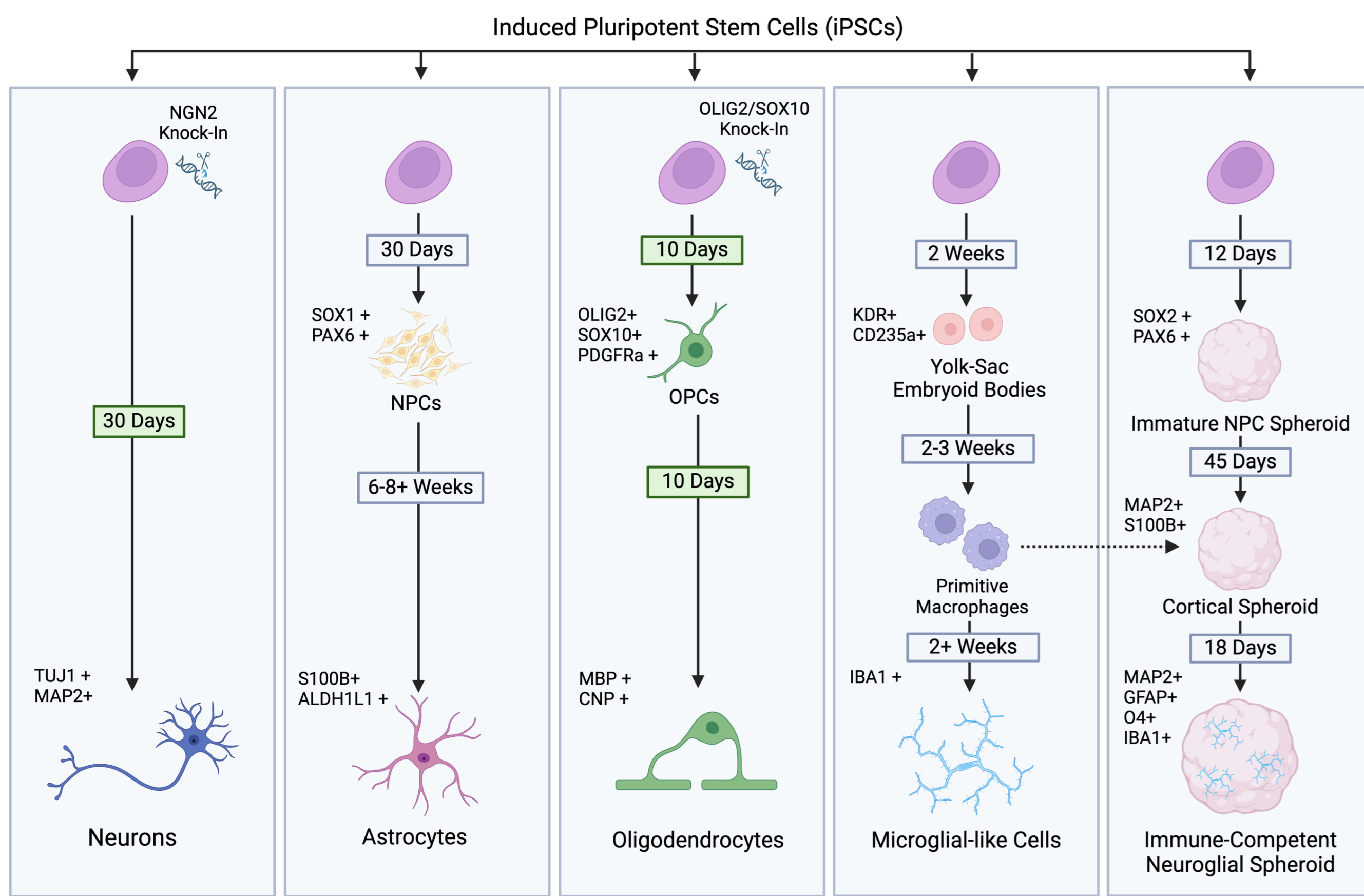
## Introduction

Human Herpesvirus 6 (HHV6) is an endemic pathogen, infecting 90% of the world's population at any given time. Acute infection manifests as rash and fever, but can progress to seizures, encephalitis, and even death. Following infection, HHV-6 establishes life-long latency in humans through genomic integration. The consequences of HHV-6 latency are currently unknown but are hypothesized to contribute to the pathogenesis of late-onset adult neurodegenerative disorders including multiple sclerosis and Alzheimer's disease.

- Understanding the effects of HHV-6 on neurons and glia requires a human model that is robust and scalable. Animal models do not recapitulate HHV-6 replication dynamics observed in humans, and primary tissue is difficult to obtain on the scale required to investigate viral infections of the brain.
- Human induced pluripotent stem cells (iPSCs) are an unlimited and accessible resource from which to generate large numbers of isogenic neurons and glia, offering a unique opportunity to study HHV-6 infection.
- Using novel optimized differentiation protocols, we can generate each of the main cell types in the CNS: neurons, astrocytes, oligodendrocytes, and most importantly the primary immune responder in the brain, microglia.
- Generating each of these cell types in individually, as well as in co-cultures, will enable capture of direct viral effects on each cell type, and allow investigation into the global inflammatory response and cell-cell crosstalk that occurs *in vivo* following infection.

## Differentiation of neurons and glia from iPSCs

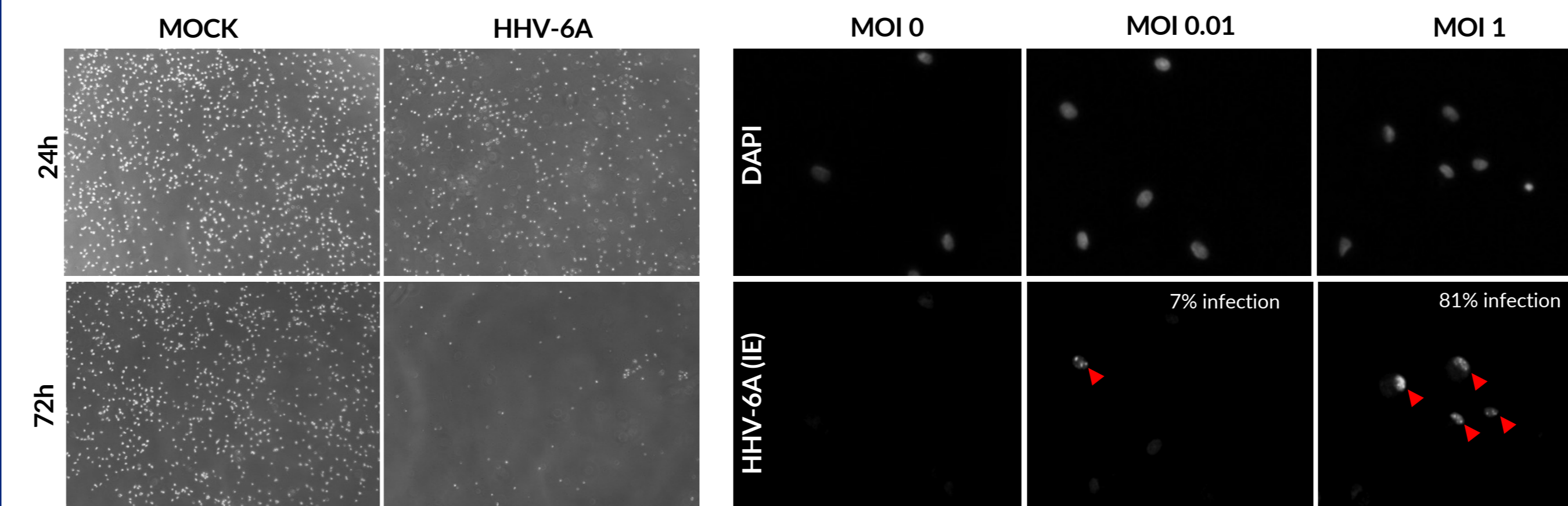
- Neurons and glia can be derived from stem cells through guided differentiation via small molecule supplementation following developmental timelines, or forced differentiation through iPSCs engineered to express a given protein when induced with an antibiotic that commits them to a particular fate.



## Project Aims

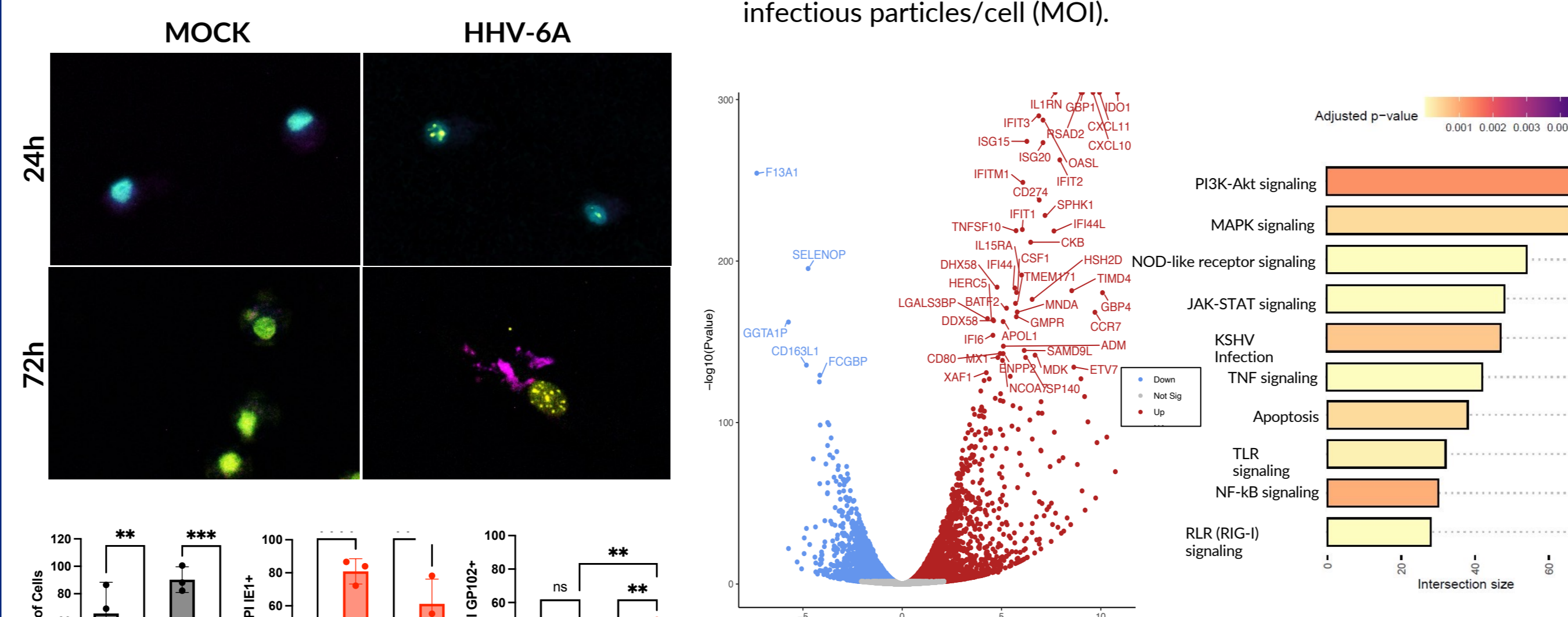
- 2D Monoculture infections:**
  - Differentiate neurons and glia from iPSCs
  - Infect 2D monocultures of each cell type with HHV-6A
  - Evaluate HHV-6A tropism within the CNS and the subsequent cellular response to viral presence and/or infection with immunofluorescence and RNASeq
- 3D Co-culture infections:**
  - Generate and infect 3D immune competent cortical spheroids comprised of neurons and glia
  - Analyze the subsequent inflammatory response on the level of single cells

## Microglia are acutely vulnerable to infection with HHV-6A in iPSC-derived monocultures



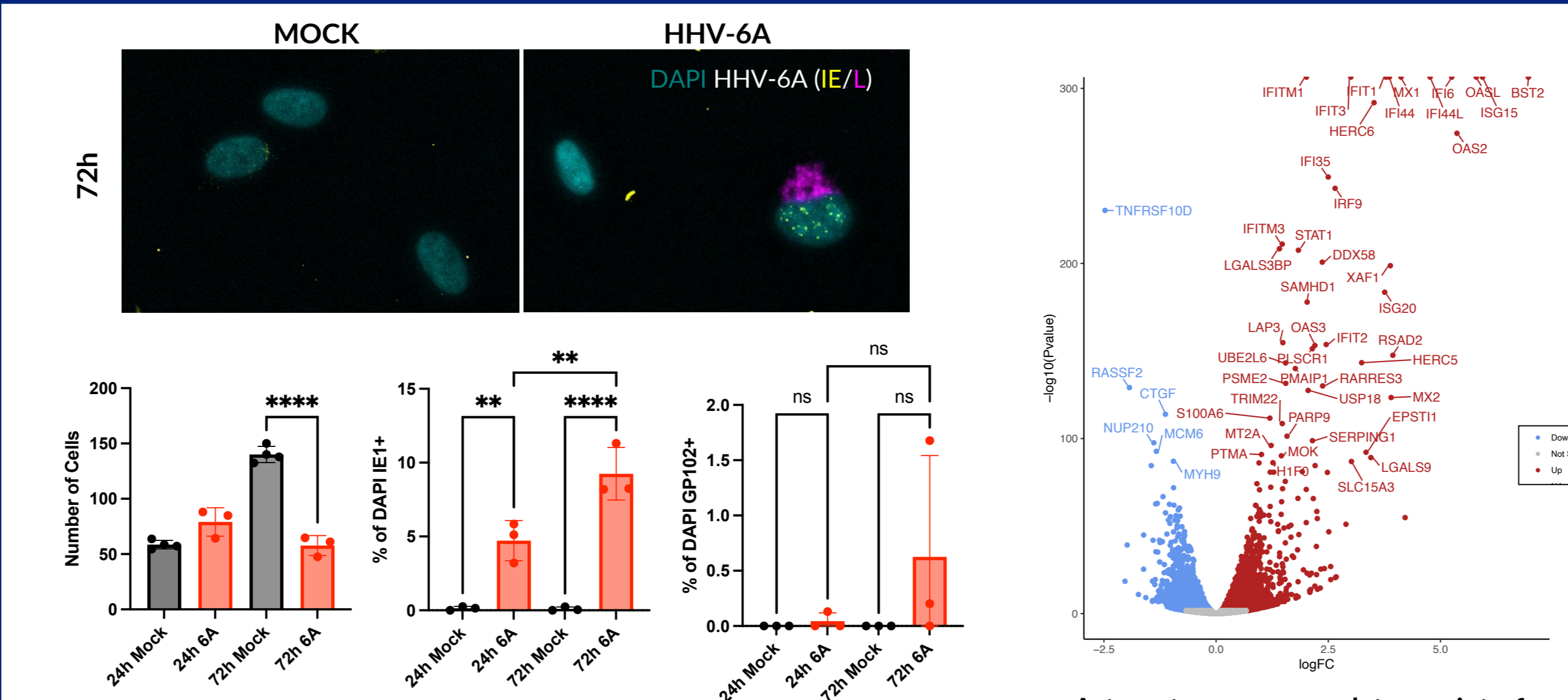
Live imaging of iPSC-derived microglia following infection with HHV-6A. Acute death observed at MOI 3 over a 72h time course relative to an uninfected control.

Immunofluorescent staining for immediate early (IE) HHV-6A viral proteins in microglia 24h post-infection across a range of viral doses. Expression of IE indicating initiation of viral replication observed following infection with the lowest dose attempted, 0.01 infectious particles/cell (MOI).



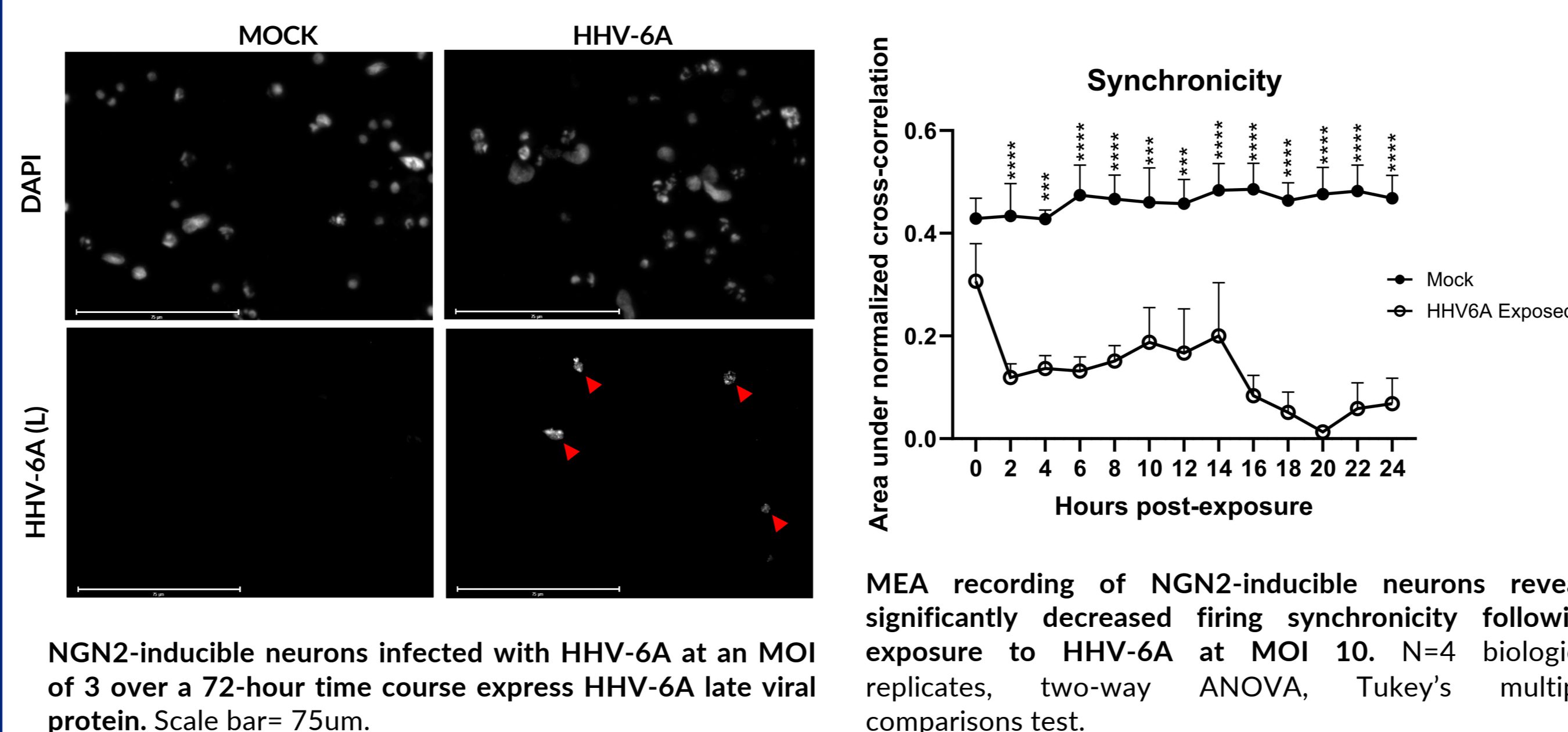
Microglia upregulate genes corresponding to immune activation and viral defense following exposure to HHV-6A. RNASeq performed on samples taken after infection at MOI 1 for 24h. Volcano plot generated in Galaxy showing significantly (p<0.05) upregulated (red) and downregulated (blue) genes, with many interferon-stimulated genes showing significant upregulation (left). GO terms associated with upregulated genes with significant (p<0.05) intersection following analysis with gProfiler (right). Terms include known pathways important to immune response, such as TLR and NF-κB signaling, which are known drivers of inflammation.

## Astrocytes decrease proliferation following infection with HHV-6A



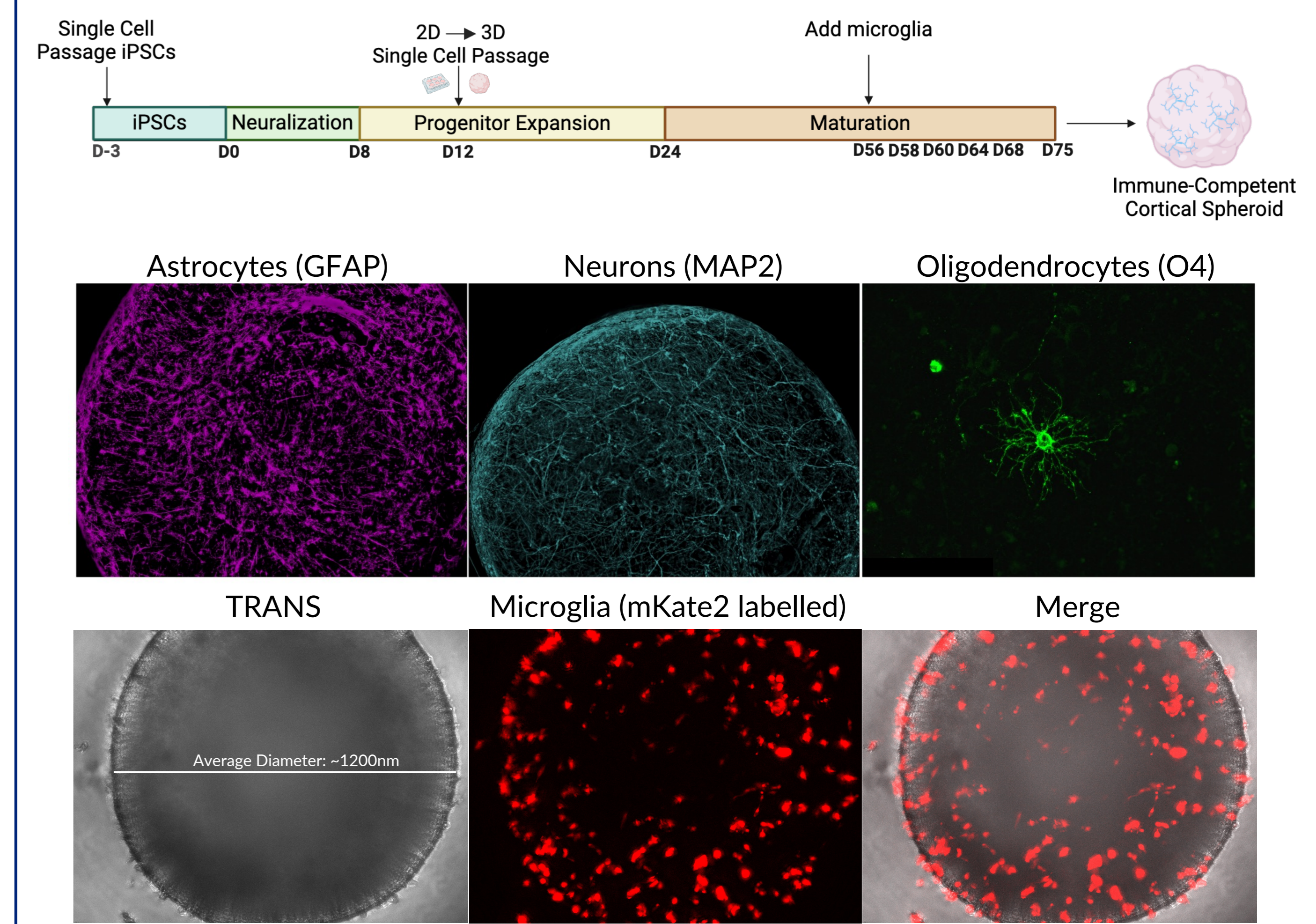
Astrocytes significantly express early (IE1)HHV-6A viral proteins at 24 and 72h, and do not proliferate at the same rate as an uninfected control by 72h. N=3 biological replicates, 9 FOV quantified/well. One-way ANOVA with Tukey's multiple comparisons test. Asterisks denote significance (P<0.05).

## Neuronal activity is significantly altered following HHV-6A exposure



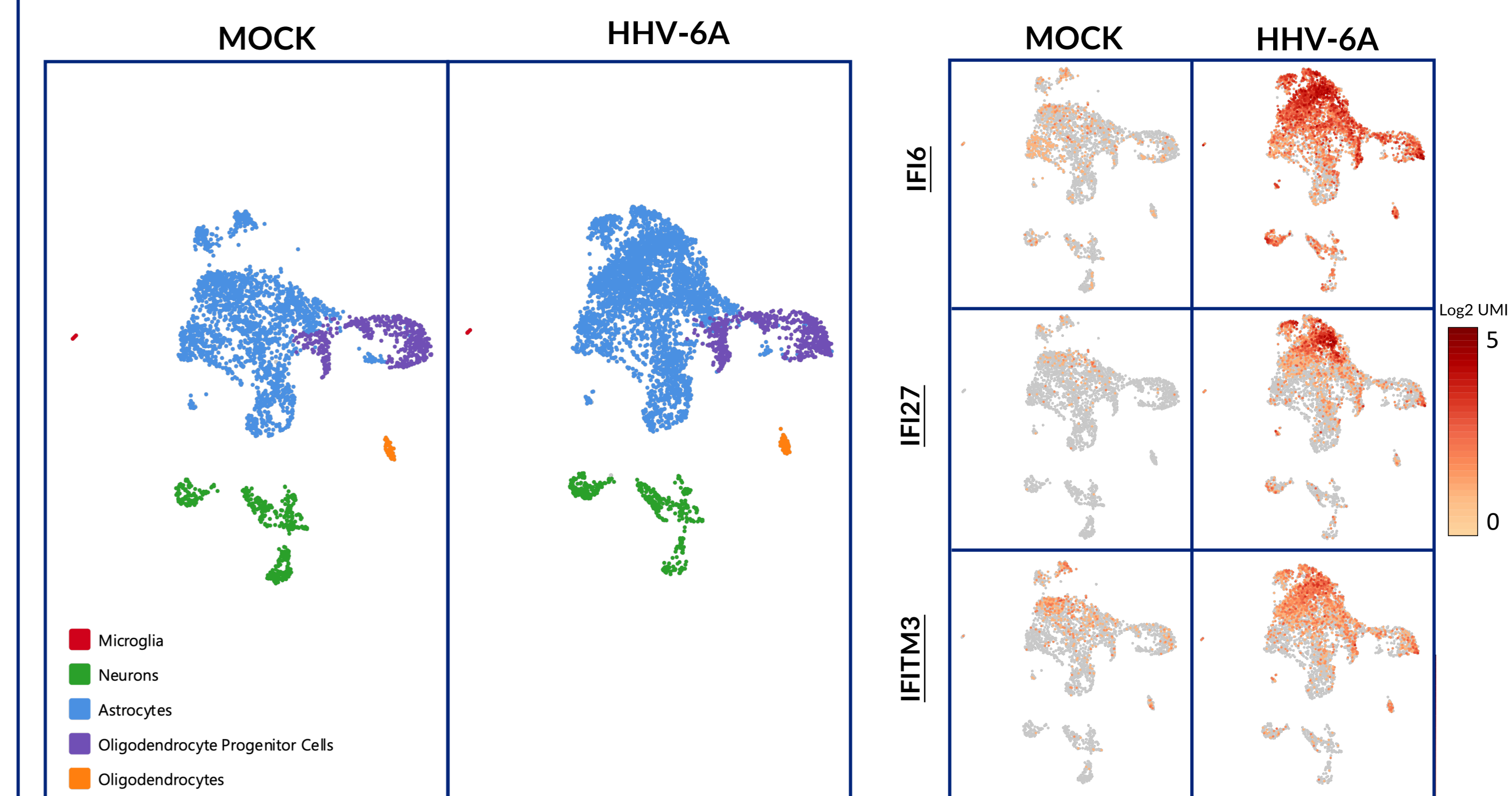
MEGA recording of NGN2-inducible neurons reveals significantly decreased firing synchronicity following exposure to HHV-6A at MOI 10. N=4 biological replicates, two-way ANOVA, Tukey's multiple comparisons test.

## Generation of immune-competent cortical spheroids



Immune-competent cortical spheroids express markers of neurons and glia after 75 days in culture. Following addition to culture medium, mKate2-labelled microglia efficiently integrate into the tissue.

## scRNASeq of immune-competent cortical spheroids infected with HHV-6A



Immune-competent cortical spheroids show induction of interferon responses following infection with HHV-6A. Sample UMAP plots of the expression of upregulated genes relative to an uninfected control (scale= log<sub>2</sub>-transformed UMI counts).

## Summary

- Stem cells can be manipulated to model the CNS, thereby enabling investigation of insults such as viral infection
- iPSC-derived neurons and glia are vulnerable to infection by HHV-6A, with infection resulting in acute cytotoxicity in microglia
- Organotypic cortical spheroid co-cultures can also be generated for infection studies, with the addition of microglia rendering cultures immune-competent
- Preliminary analyses of scRNASeq performed on infected co-cultures suggests strong induction of interferon-mediated response to infection