

Validation of astrocytic cAMP signalling to study therapeutic targets for Alzheimer's disease

Lucy Granat¹, Sarah Jolly¹, Jamie Bilsland², Fiona. E. Ducotterd¹

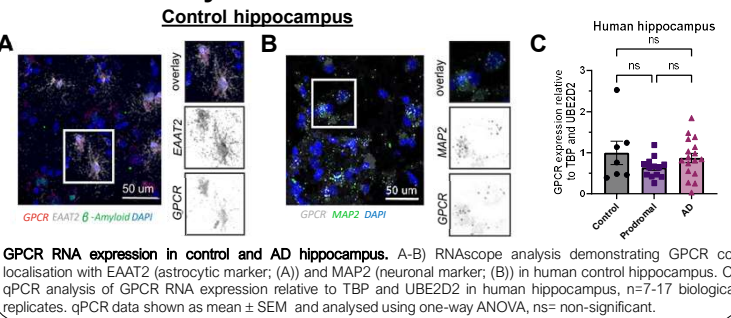
¹ Alzheimer's Research UK UCL Drug Discovery Institute, London, United Kingdom, ² AstronuTx, London, United Kingdom



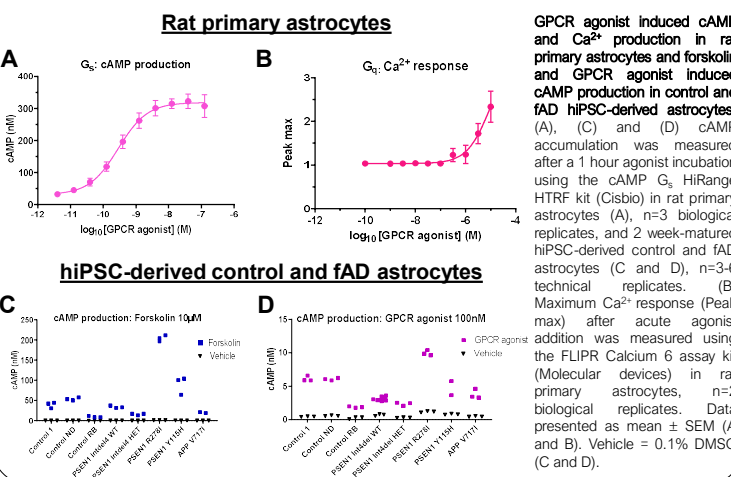
1. Introduction

Astrocytes play a fundamental role in pathological processes associated with neurodegenerative diseases, including neuroinflammation, impaired glutamate uptake, reduced neurotrophic support and defective metabolism. Activation of cAMP signalling in astrocytes elevates glycolytic rate, increases glutamate transporter and neurotrophic factor expression, and suppresses the immune response. Molecules that modulate astrocytic cAMP signalling are therefore potential therapeutic targets for neurodegenerative diseases such as Alzheimer's disease (AD). To support the identification and validation of astrocyte-centric targets, we have optimised a suite of *in vitro* assays including cAMP, Ca²⁺, and metabolic assays, multi-electrode array (MEA) and RNA-seq in rodent, human foetal and hiPSC-derived control and familial AD (fAD) astrocyte models. We have tested the adenylyl cyclase activator forskolin and a GPCR agonist in these assays to validate and increase our confidence in astrocytic cAMP signalling as an AD drug target.

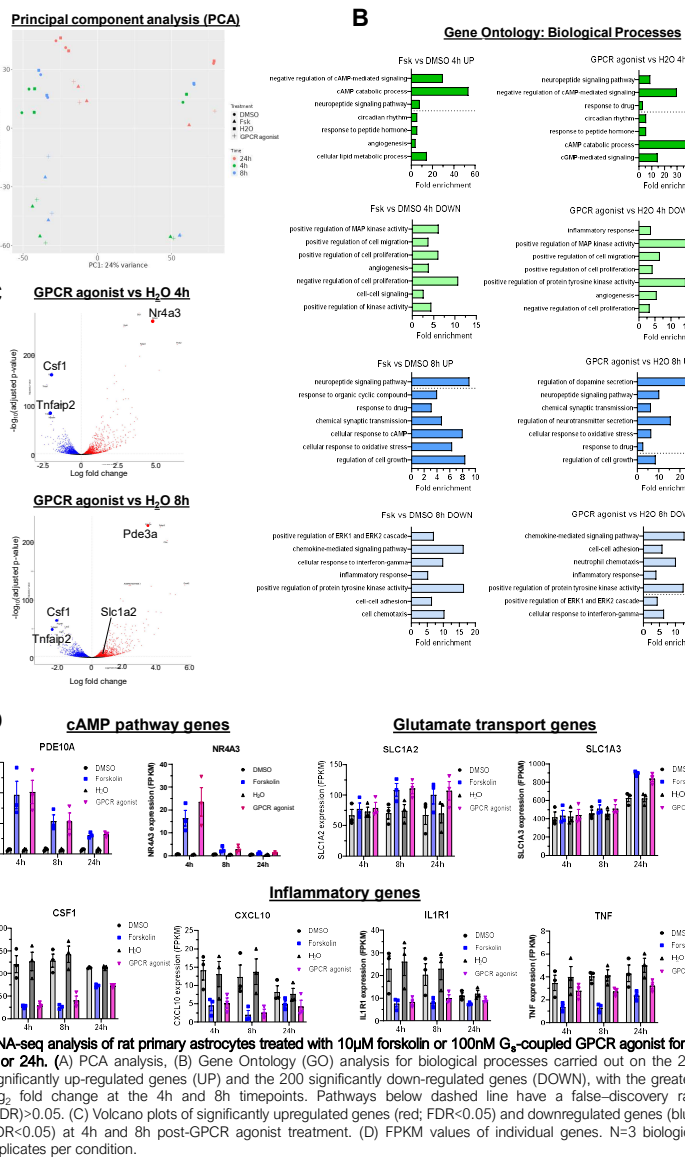
2. Levels of a G_s-coupled GPCR expressed by human astrocytes and neurons are unaltered in AD



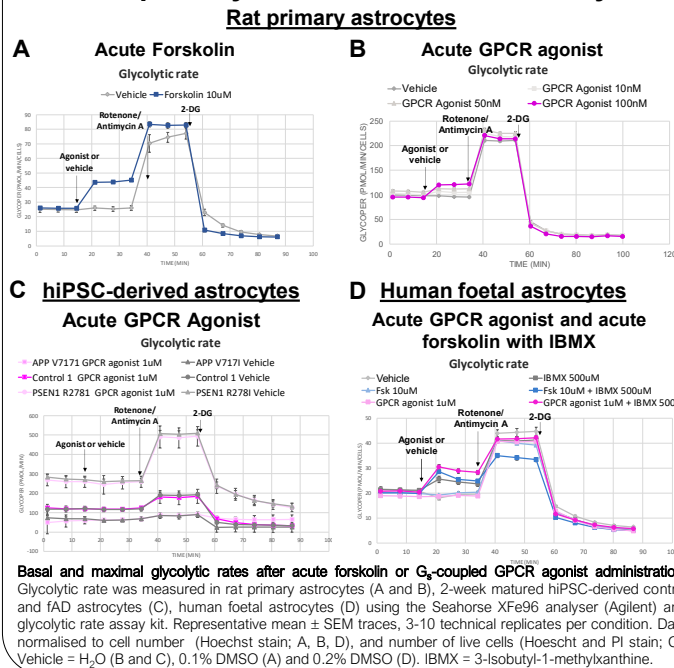
3. Forskolin and a GPCR agonist induce cAMP activation in astrocytes *in vitro*



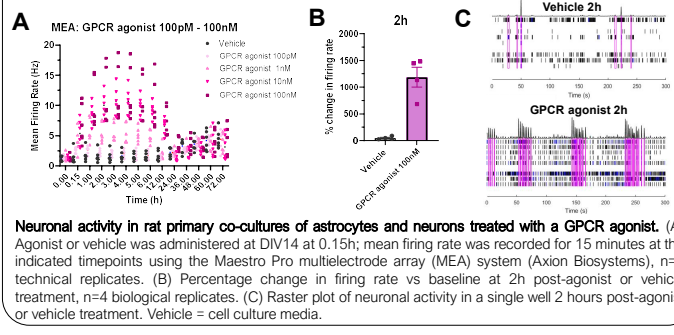
4. cAMP activation in rodent primary astrocytes reduces the expression of inflammatory genes



5. cAMP activation increases glycolytic rate in rodent primary and human foetal astrocytes



6. A GPCR agonist enhances neuronal activity in rodent co-cultures of astrocytes and neurons



7. Summary

We have demonstrated that RNA levels of a G_s-coupled GPCR are not altered in the AD brain. Both forskolin and the GPCR agonist elevate glycolytic rate in rat and human foetal astrocytes, suggesting that the mechanism is cAMP-mediated. Interestingly, the GPCR agonist enhances neuronal firing *in vitro* in neuron-astrocyte co-cultures. Finally, RNA-seq analysis identified cAMP pathway and inflammatory genes as major targets of both forskolin and GPCR agonist administration. Overall, our data suggest that activation of astrocytic cAMP signalling has exciting therapeutic potential as a treatment for AD.



Work supported by Alzheimer's Research UK Grant ARUK-2021 DDI-UCL.

Contact: f.ducotterd@ucl.ac.uk

