

INTRODUCTION

- Frontotemporal lobar degeneration (FTLD) can be caused by autosomal dominant mutations in the *MAPT* gene (FTLD-Tau), such as P301L.
- Pathological tau is hyperphosphorylated and mislocalizes to the post-synaptic compartment where it can cause synaptic disruption.

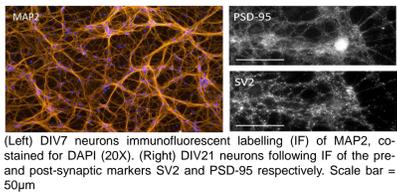
AIMS

1. Develop and characterise a neuronal model of tauopathy
2. Determine if there are any changes to network activity
3. Identify tau binding partners that might alter synaptic function

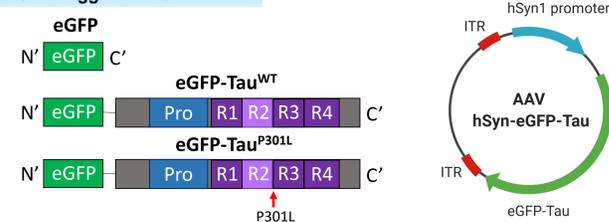
PRIMARY MOUSE CORTICAL NEURONS TRANSDUCED WITH AAVS CONTAINING eGFP-TAU CONSTRUCTS

Primary Cortical Neurons

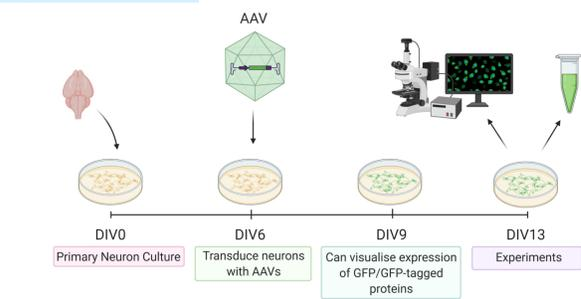
Primary cortical neurons from embryonic CD1 mice form extensive networks and active synapses in culture.



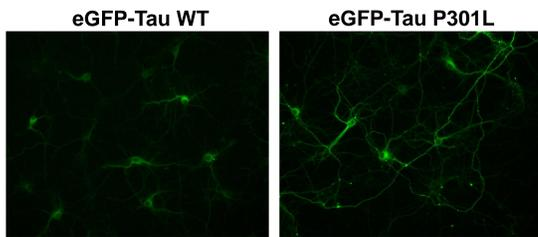
eGFP-Tagged Tau Constructs



Experimental Timeline

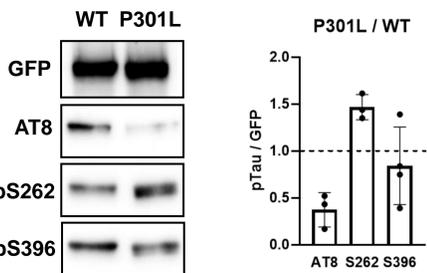


Transduced Neurons

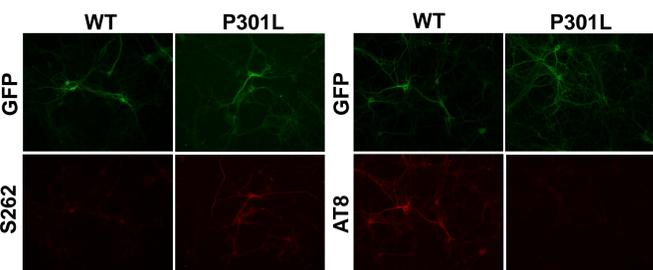


DIV14 primary cortical neurons overexpressing eGFP-tagged tau. Note the increased levels of tau in neurons expressing eGFP-Tau P301L

P301L TAU IS DIFFERENTIALLY PHOSPHORYLATED TO WT TAU



P301L tau is more phosphorylated at Ser262 but less phosphorylated at AT8. Immunoblotting for different phospho-epitopes of tau, AT8 represents phosphorylation at Ser202/Thr205. Results were normalised to GFP (total tau) and then compared against WT tau. Error bars are +/- standard deviation, n = 3.



ACKNOWLEDGEMENTS

This work was supported in part by grant MR/N0137941/1 for the GW4 BIOMED MRC DTP, awarded to the Universities of Bath, Bristol, Cardiff and Exeter from the Medical Research Council (MRC)/UKRI.

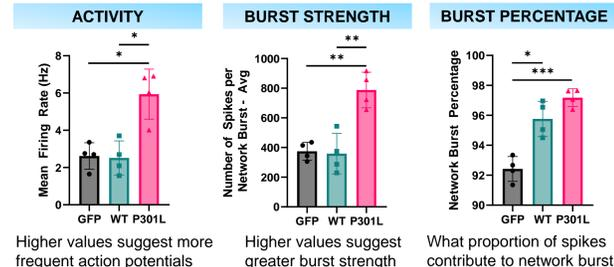
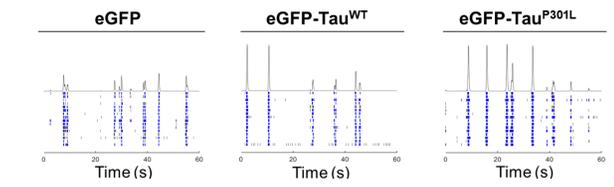
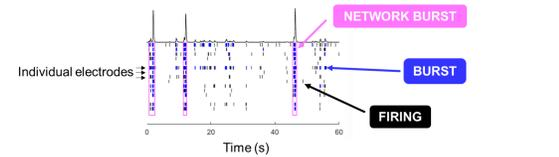
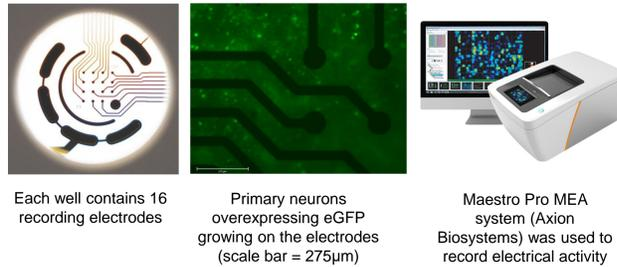
Figures were created with BioRender.com.

IS NETWORK FUNCTION ALTERED IN PRIMARY CORTICAL NEURONS OVEREXPRESSING eGFP-TAU P301L?

- Pathological tau is known to mislocalize to dendritic spines and disrupt synaptic function.
- Therefore we wanted to investigate whether network excitability was altered in this model.
- To do this, we used a multi-electrode array.

METHOD

- Primary neurons were plated onto a CytoView MEA plate at a density of 120,000 cells/well.
- Recordings were undertaken at DIV12.
- Baseline recordings were undertaken for 10 minutes.



Overexpression of P301L tau increases network activity. Brown-Forsythe and Welch ANOVA tests with multiple comparisons. p* < 0.05, p** < 0.01, p*** < 0.001, p**** < 0.0001. Error bars are +/- standard deviation, n = 4.

KEY FINDINGS

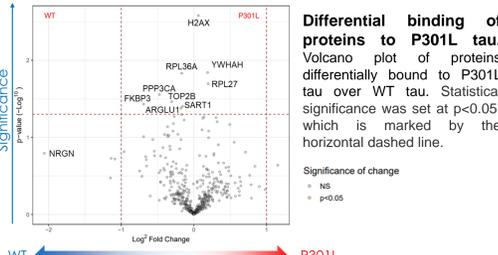
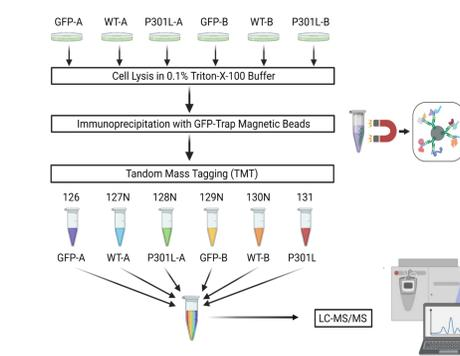
- Overexpression of P301L tau increases baseline excitability in primary cortical neurons
- This could result in plasticity changes which make synapses vulnerable.

CAN THE TAU INTERACTOME EXPLAIN THE INCREASE IN BASAL EXCITABILITY?

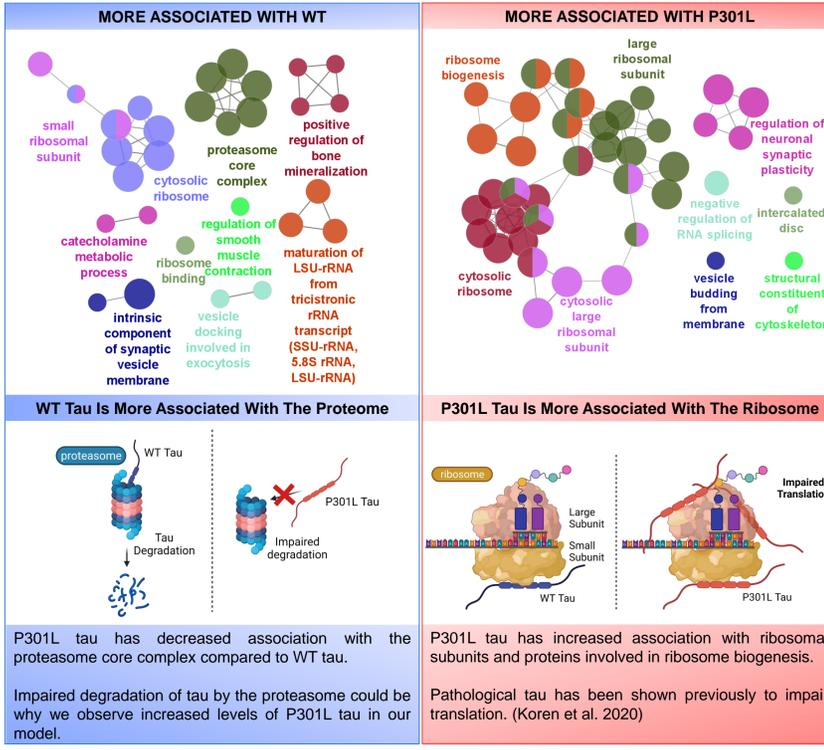
In order to investigate how tau interactions might be affecting neuronal activity, a proteomics study was undertaken to identify the tau interactome.

METHOD

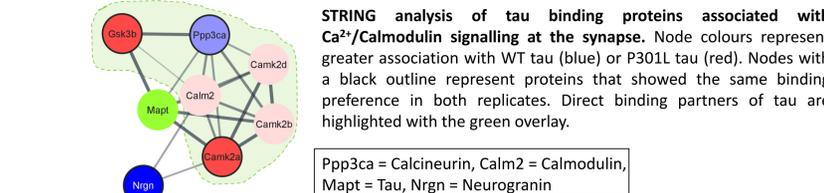
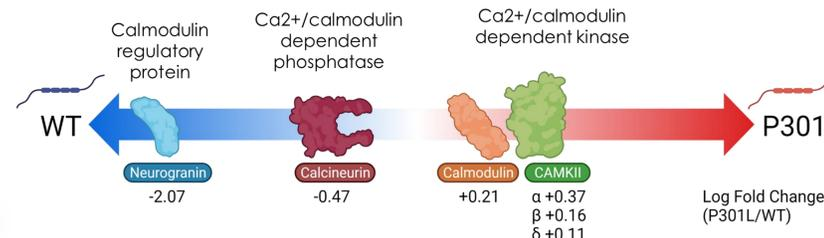
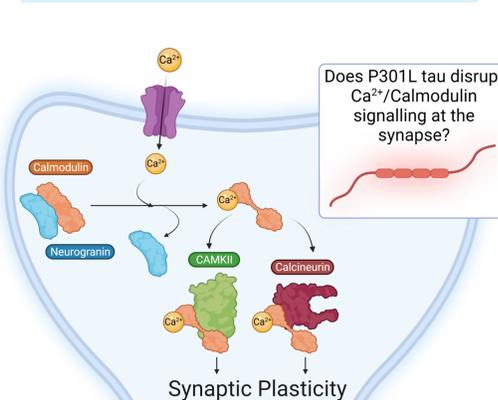
- A GFP-Trap was used to immunoprecipitate eGFP/eGFP-tagged tau and any proteins associated with it
- At the Bristol Proteomics Facility, samples were labelled by TMT to enable quantitative analysis and analysed by LC-MS/MS



Differential association of proteins with WT and P301L tau. ClueGO Biological Process, Cellular Compartment and Molecular function analysis of proteins that were more associated with (left) WT tau over P301L tau or (right) P301L tau over WT tau in both repeats. Node colours represent functionally grouped networks. Node size corresponds to p-value corrected with Bonferroni step-down.



Tau Associates with Proteins Involved in Ca²⁺/Calmodulin Signalling at the Synapse



CONCLUSION

In this model of tauopathy:

- P301L tau alters association with key Ca²⁺ dependent synaptic binding partners and enhances network excitability which could result in synaptic vulnerability
- Increased synaptic activity correlates with changes in tau phosphorylation, most notably an increase at pSer262

FUTURE WORK

- Link changes in synaptic proteome to enhanced excitability
- Investigate what effect pathological tau has on the Ca²⁺/Calmodulin signalling pathway

REFERENCES

Bindea G, Mlecnik B, Hackl H, Charoentong P, Tosolini M, et al. ClueGO: a Cytoscape plug-in to decipher functionally grouped gene ontology and pathway annotation networks. *Bioinformatics*. 2009, 25(8):1091-3.
Bindea G, Galon J, Mlecnik B, CluePedia Cytoscape plugin: pathway insights using integrated experimental and in silico data. *Bioinformatics*. 2013, 29(5):661-3.
Hole K, Staniaszek L, Balan Gayathri M, Mason J, Brown J & Williams R. (-)-Epicatechin inhibits progressive tau pathology in Tg4510 mice independent of direct actions at GSK3β. *Frontiers in Neuroscience*, 2021
Koren S, Galvis-Escobar S & Abisambra J. Tau-mediated dysregulation of RNA: Evidence for a common molecular mechanism of toxicity in frontotemporal dementia and other tauopathies. *Neurobiology of Disease*, 2020