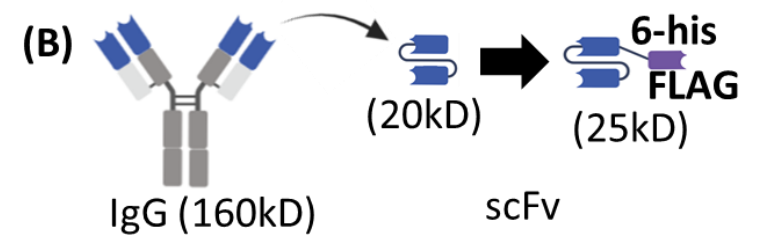
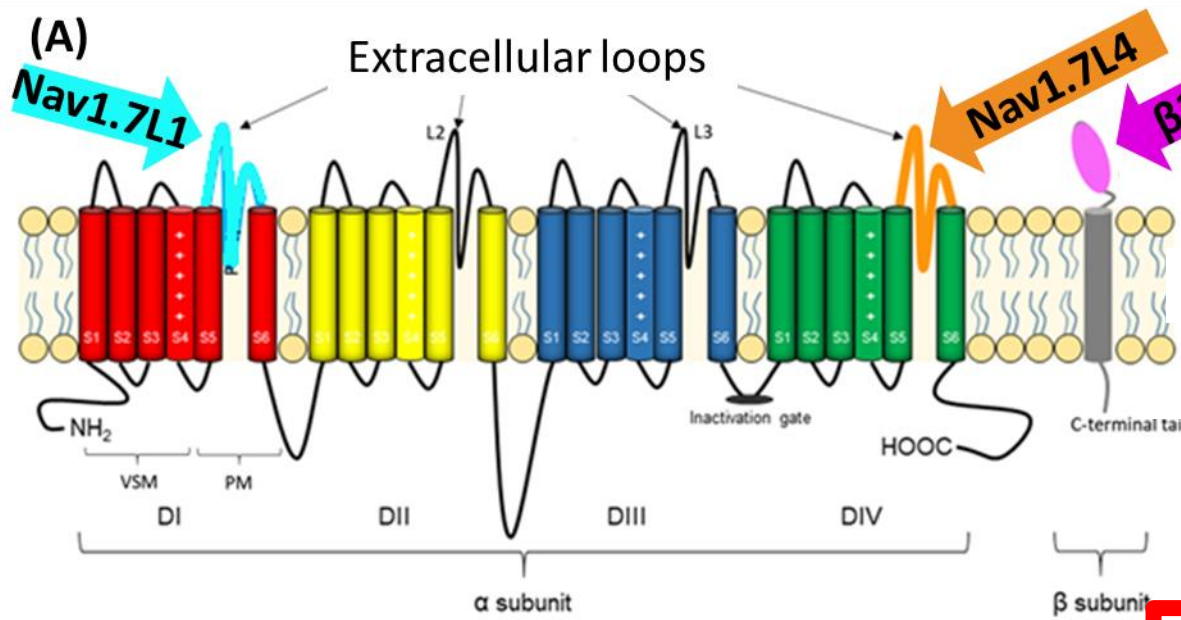
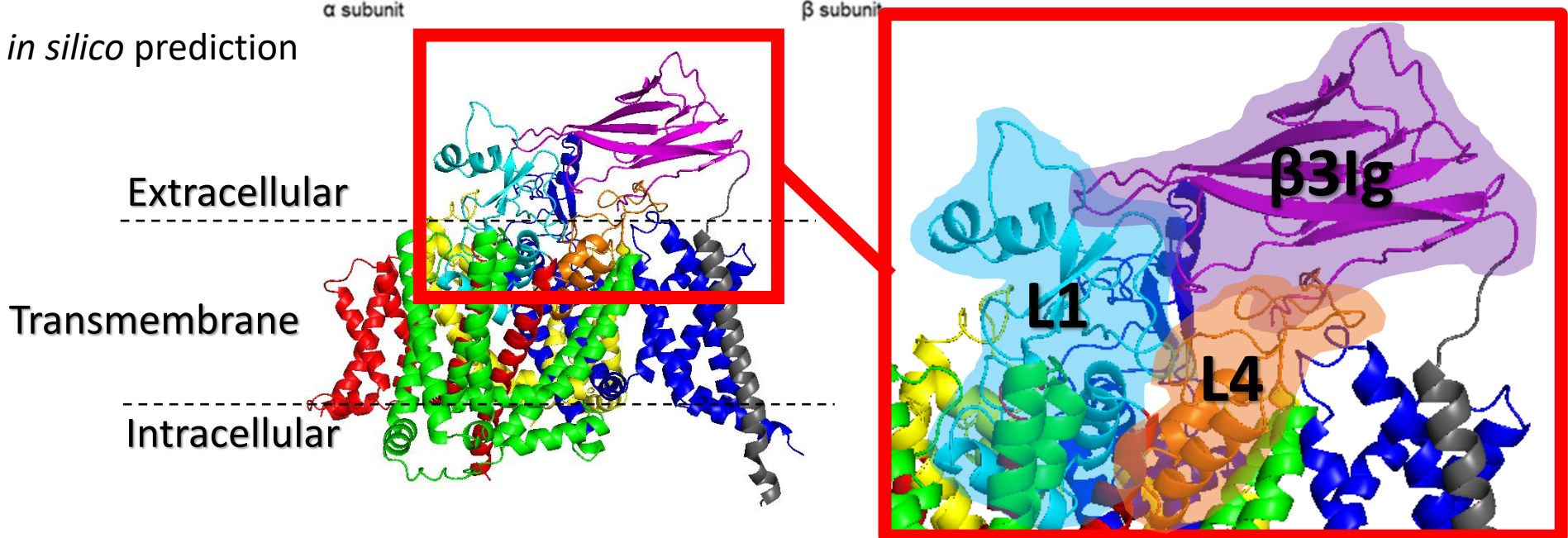


Introduction



$\beta 3$ -subunit is co-expressed with Nav1.7 in pain-sensing neurons, but its binding site remains unknown. We used *in silico* modelling to predict the $\beta 3$ binding site. We used scFvs against Nav1.7 domain I extracellular loop (L1) and domain IV extracellular loop (L4) to test this model.

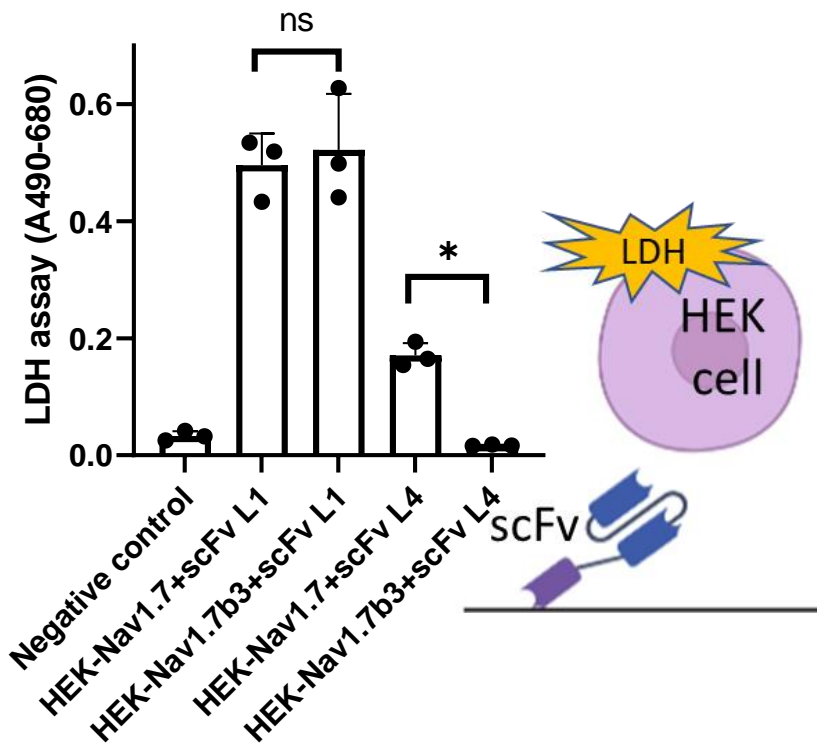
(C) *in silico* prediction



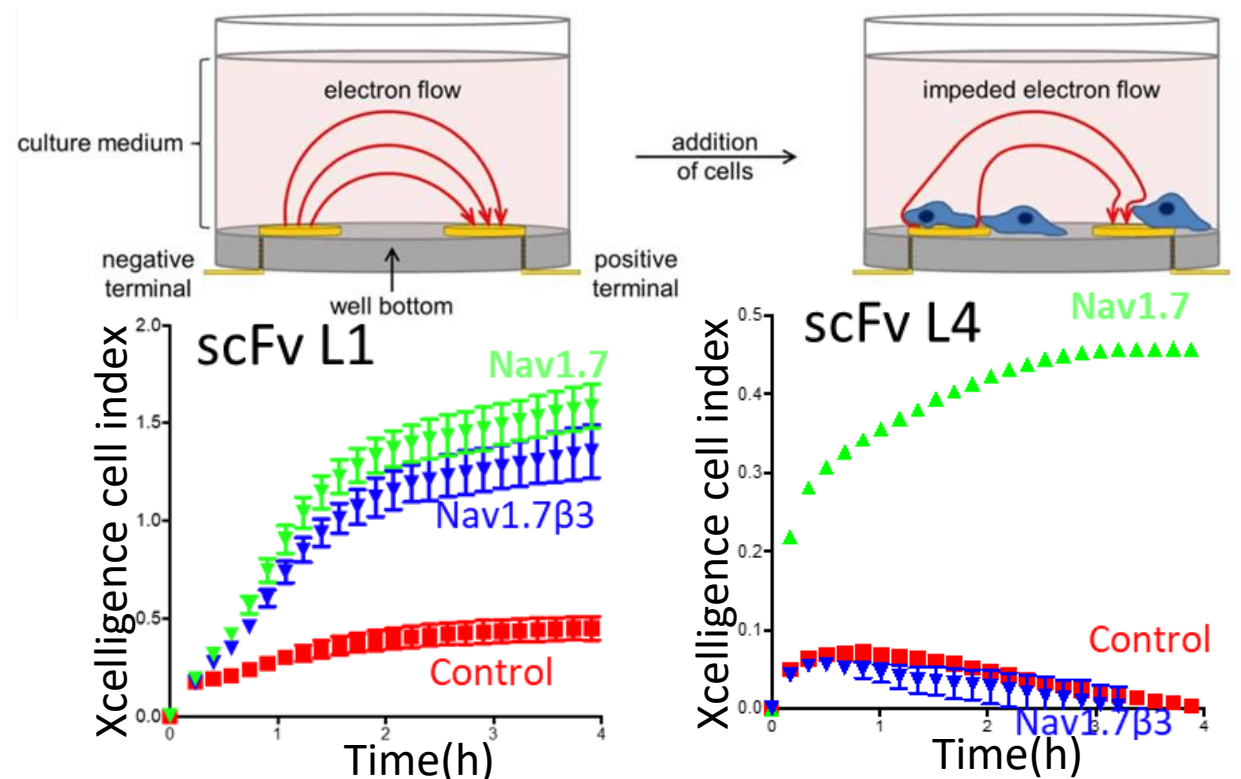
Hypothesis: L1 scFv should bind in the presence of $\beta 3$ Ig domain, but L4 scFv should not.

Methods & Results

Steady-state cell binding assay



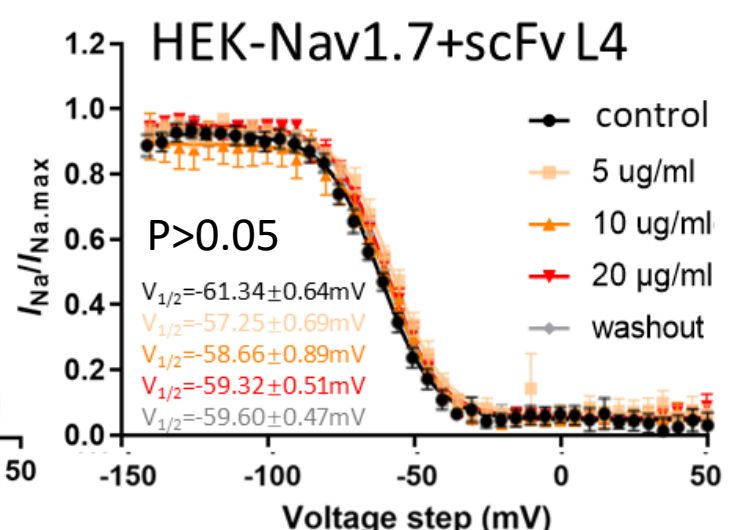
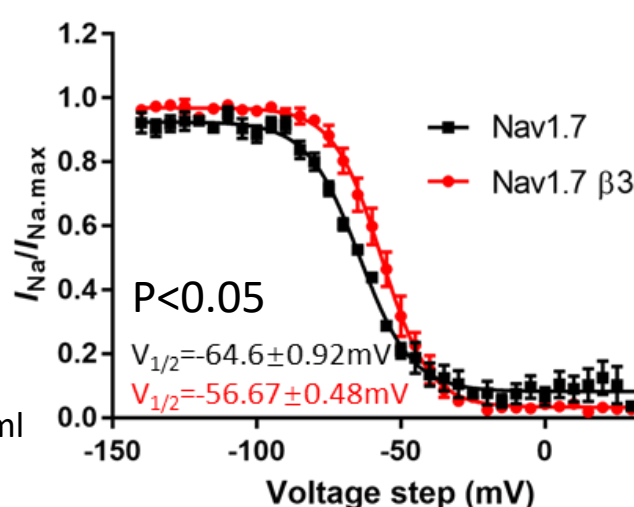
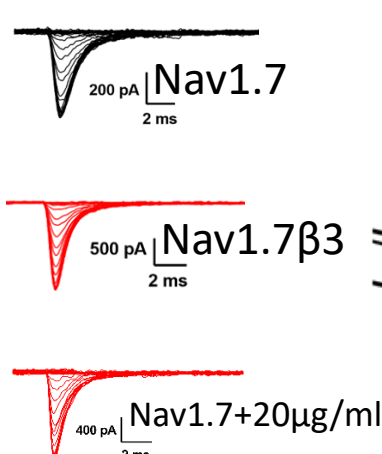
Real-time cell binding assay



Conclusion: These results supported the *in silico* modeling of $\beta 3$ binding to Nav1.7.

Effect of $\beta 3$ or scFv on Nav1.7 inactivation

Nanon Patchliner automated patch-clamp



Conclusion: Unlike $\beta 3$, the L4 scFv had no effect on Nav1.7 inactivation.