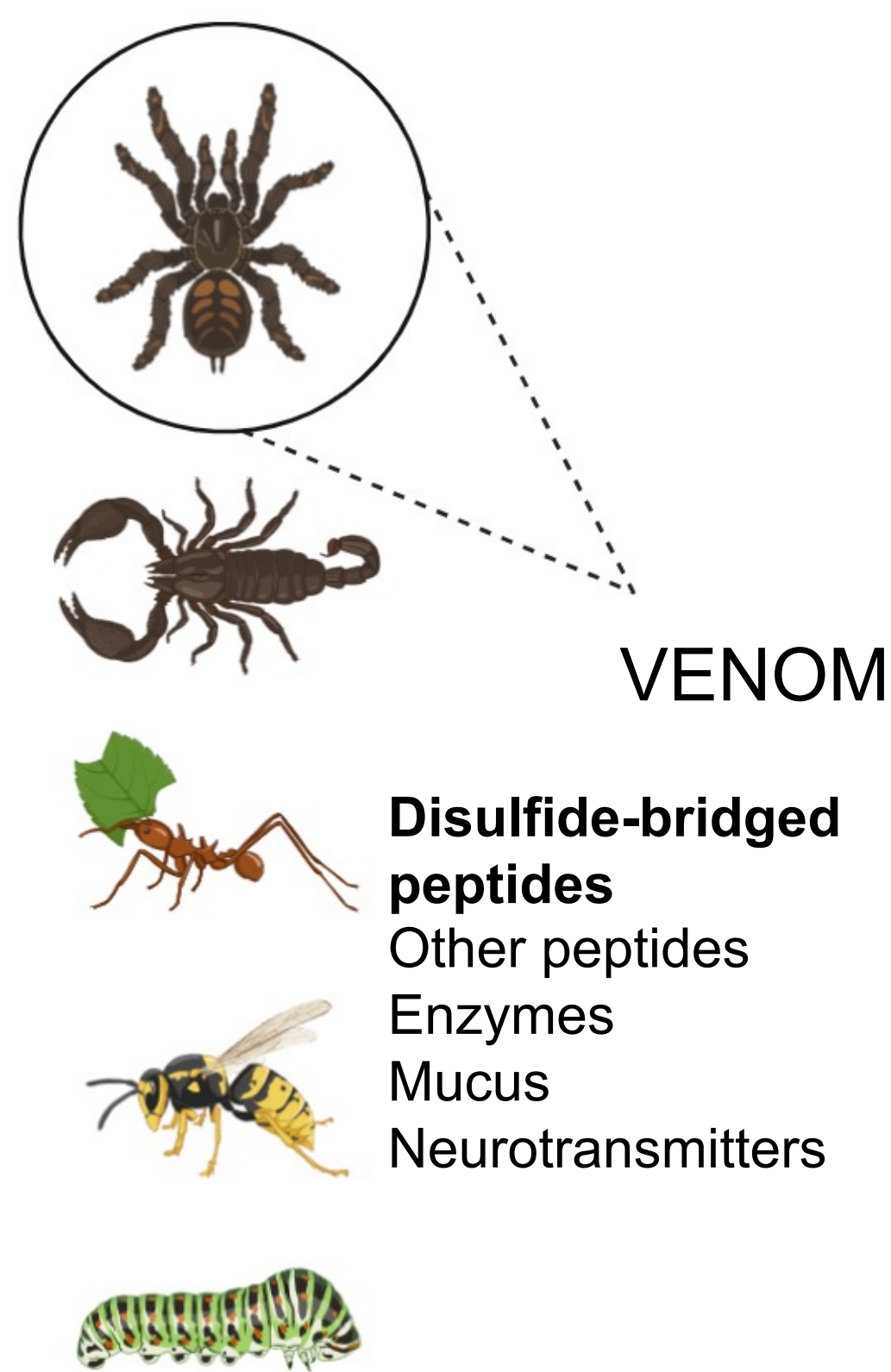


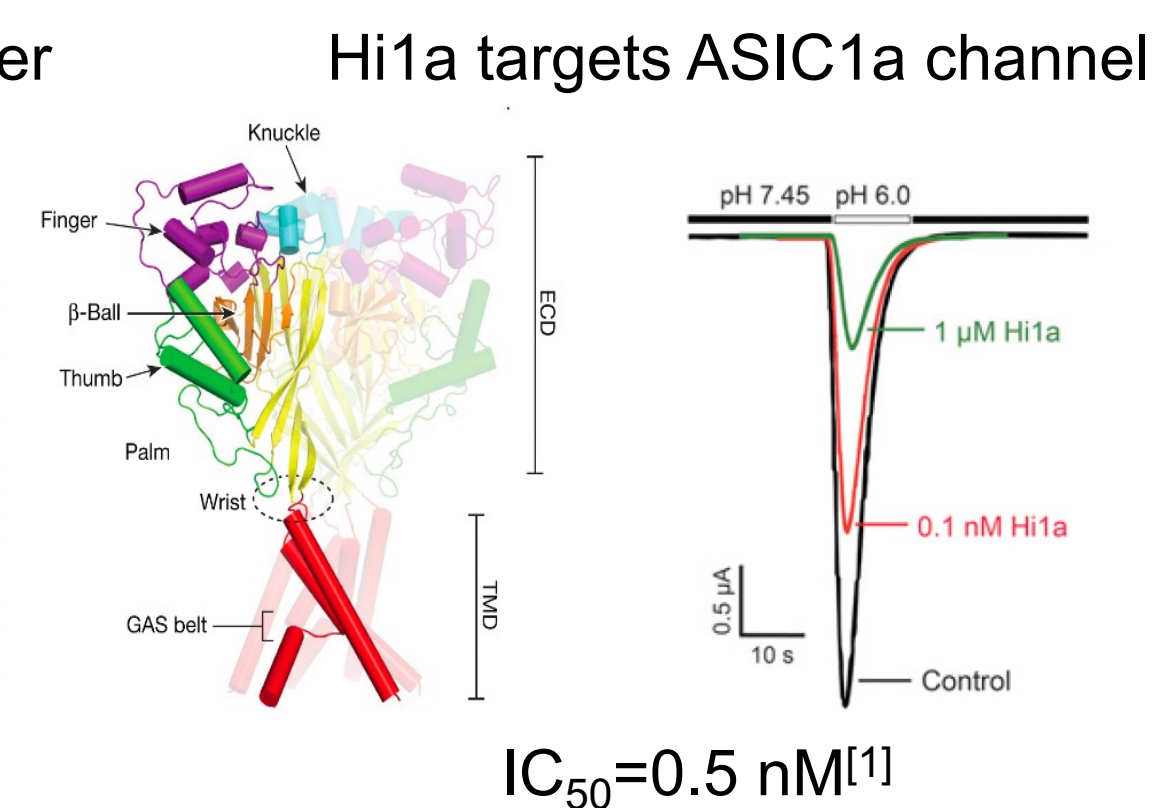
# From bites to insights: Utilization of arachnid venoms in ion channel studies and their potential in therapeutic development

**Daniela Rojas-Azofeifa**, Luca T. Rieder, Linda Blomster, Glenn F. King.  
Institute for Molecular Bioscience, University of Queensland, Australia  
uqdrojas@uq.edu.au

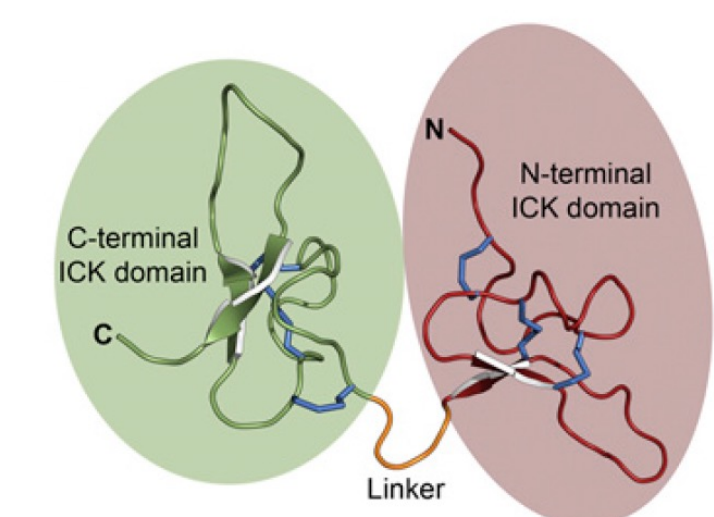
## Inspired by nature to create heart therapeutics



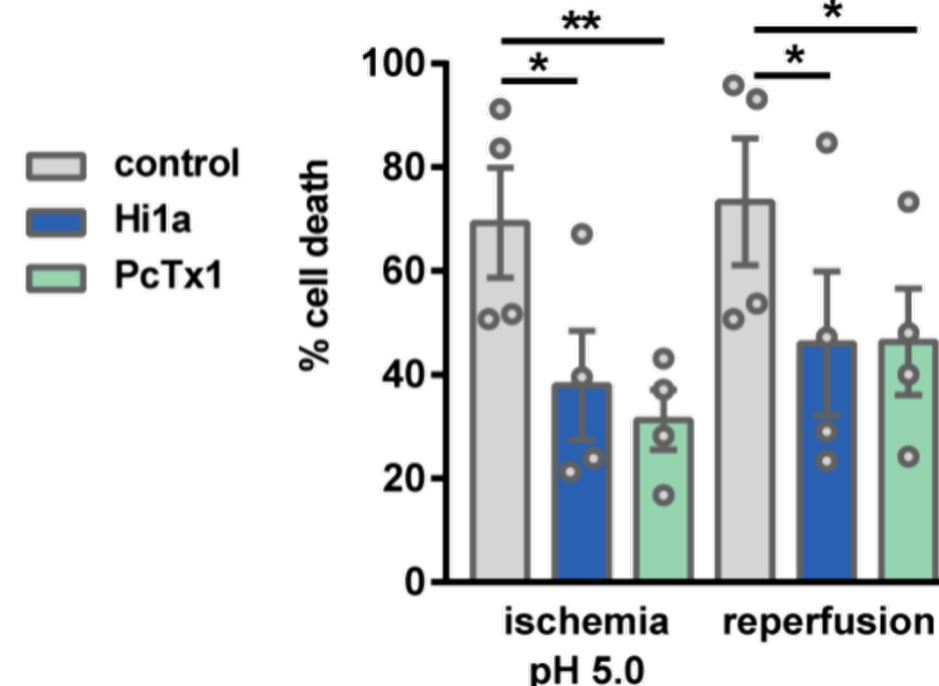
Australian funnel-web spider  
*H. infensa*.



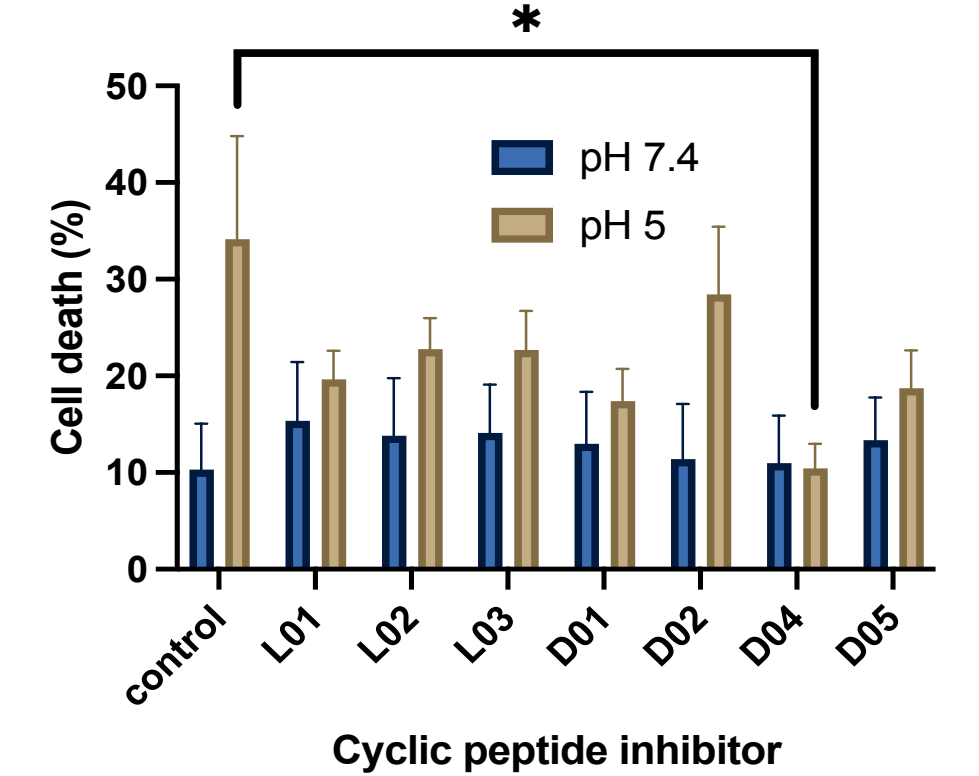
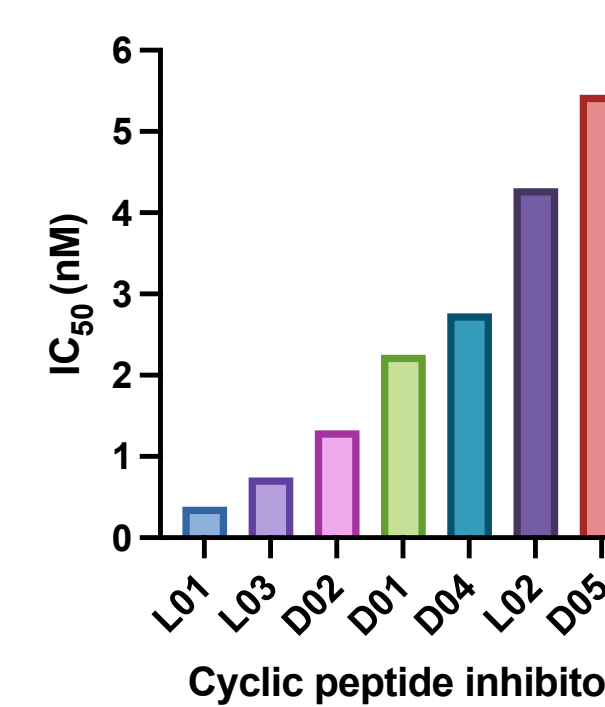
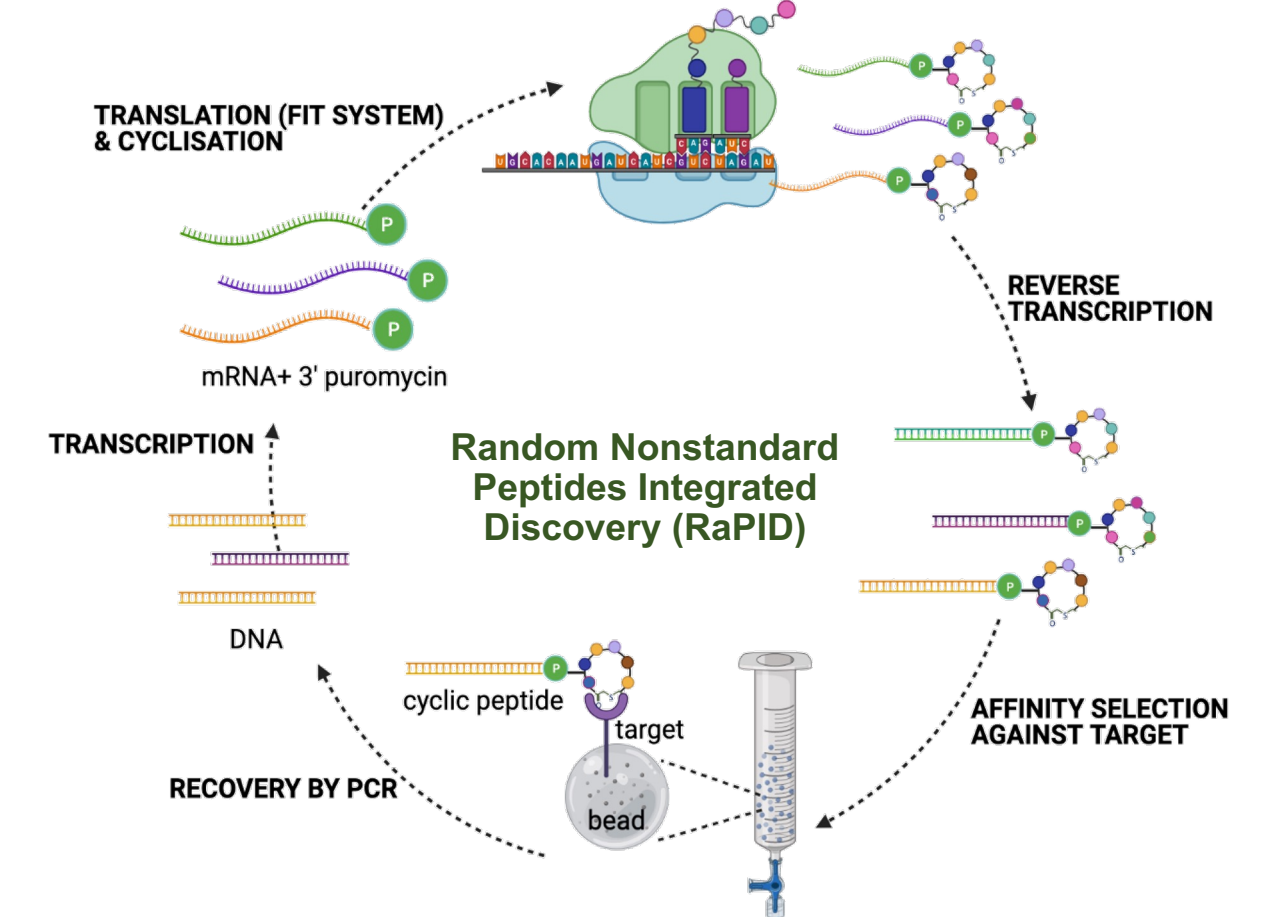
Hi1a peptide from the venom [1]



*In vitro* effect of Hi1a in the heart [2]

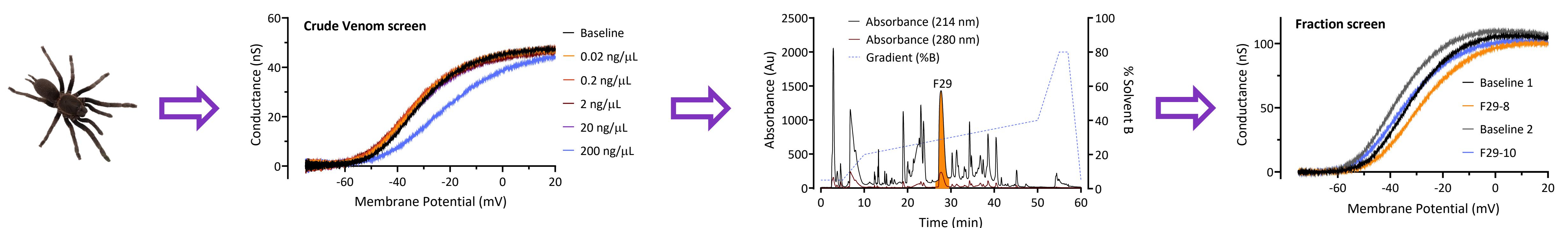


We produced smaller and less complex peptides, with a high potency for ASIC1a and easier to produce

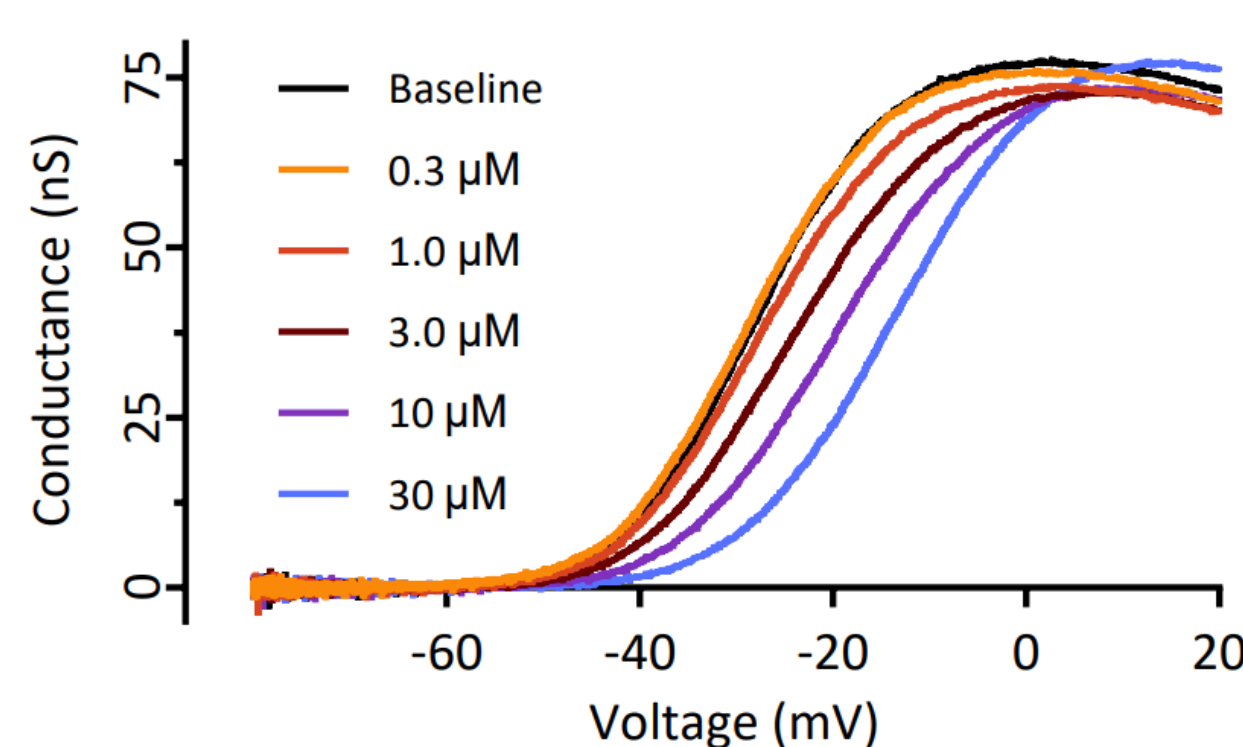


## Novel peptide modulators of K<sub>v</sub>7.2-7.3

Mutations in the voltage-gated potassium channel K<sub>v</sub>7.2-7.3 are associated with a wide spectrum of early-onset epileptic disorders. Patients with a **gain of function** mutation in the encoding KCNQ2 genes show severe developmental delay with prominent language impairment and autistic features, often accompanied by infantile- to childhood-onset epilepsy.[3] To treat this phenotype, selective negative modulators of the ion channel that shift the activation potential back towards normal activity are needed. Spider venoms provide a rich source of peptides that target ion channel activity.[1]



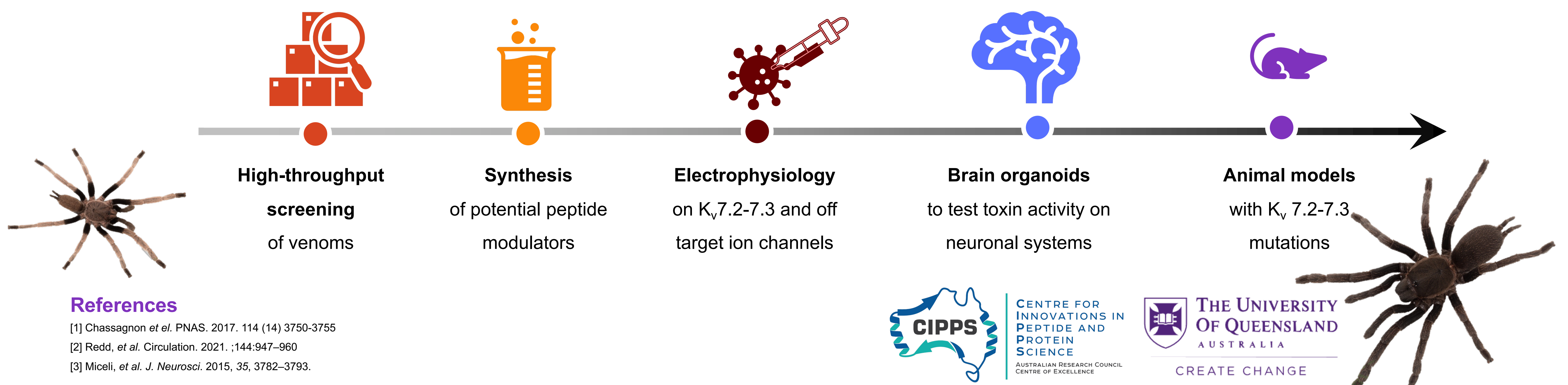
Effect of the venom peptide modulator obtained from the fraction



Several challenges:

- Screening venoms using manual patch clamp.
- Hundreds of venoms to screen
- Folding of peptides
- Off-target problems

## Outlook for finding modulators of K<sub>v</sub>7.2-7.3



### References

- [1] Chassagnon *et al.* PNAS. 2017. 114 (14) 3750-3755  
 [2] Redd, *et al.* Circulation. 2021. ;144:947-960  
 [3] Miceli, *et al.* J. Neurosci. 2015, 35, 3782-3793.



CREATE CHANGE