

## Identification of novel activators of two-pore domain potassium (K2P) channels

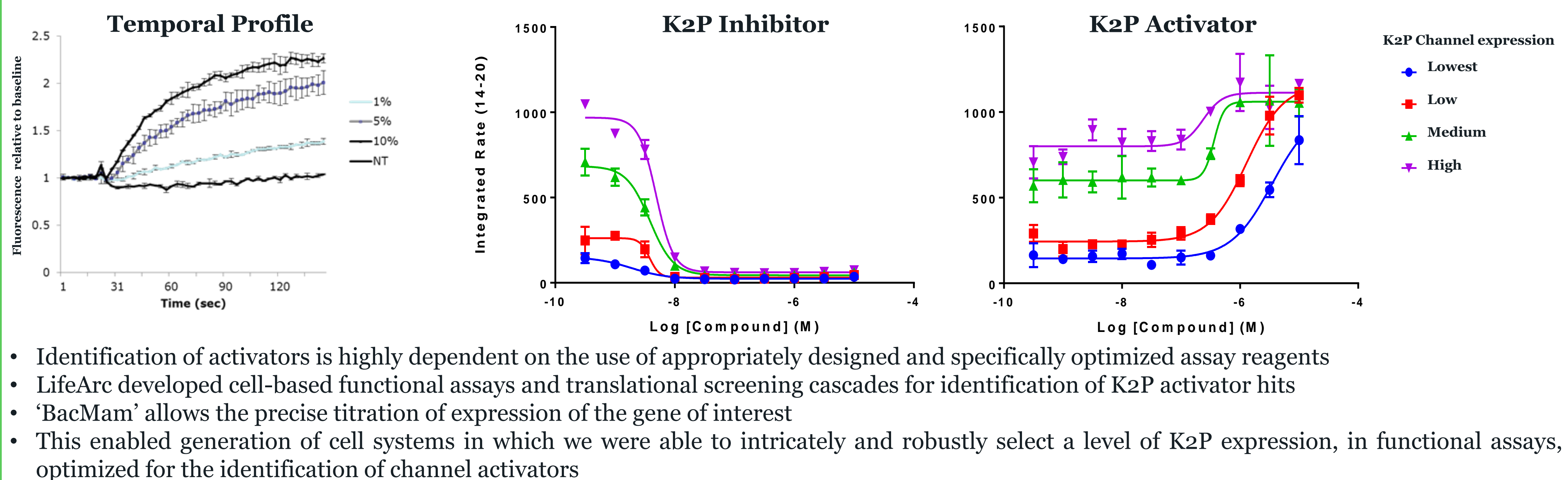
David McCoull, Lewis Byrom, Jonathan M Large, Jeff Jerman, Paul D Wright

LifeArc, Accelerator Building, Open Innovation campus, Stevenage, SG1 2FX, United Kingdom

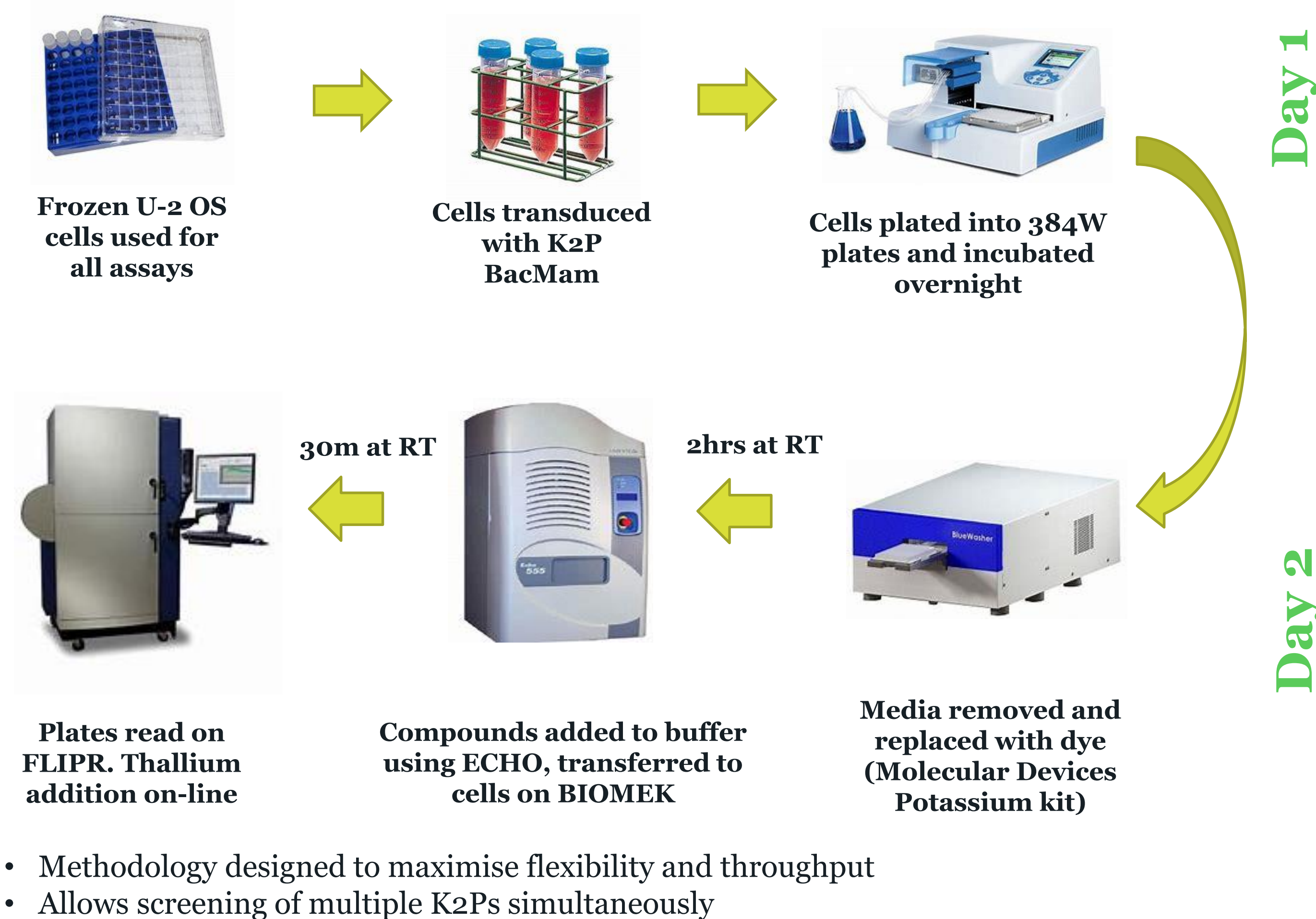
### OVERVIEW

- Two-pore domain potassium channels (K2Ps) carry background (or leak) potassium current
- Primarily act to maintain resting membrane potential.
- K2P channels are characterised by their four transmembrane domain, two-pore topology
- Genetic and functional evidence points to a role in multiple pathophysiology, including pain and migraine
- K2Ps have proved difficult to modulate with small molecules and there is a lack of useful specific pharmacological tools
- This has limited the interrogation of their precise physiological function and efforts to generate K2P based therapeutics
- LifeArc developed a novel system to identify K2P activators, with the aim of providing tools for research and ultimately novel therapeutics

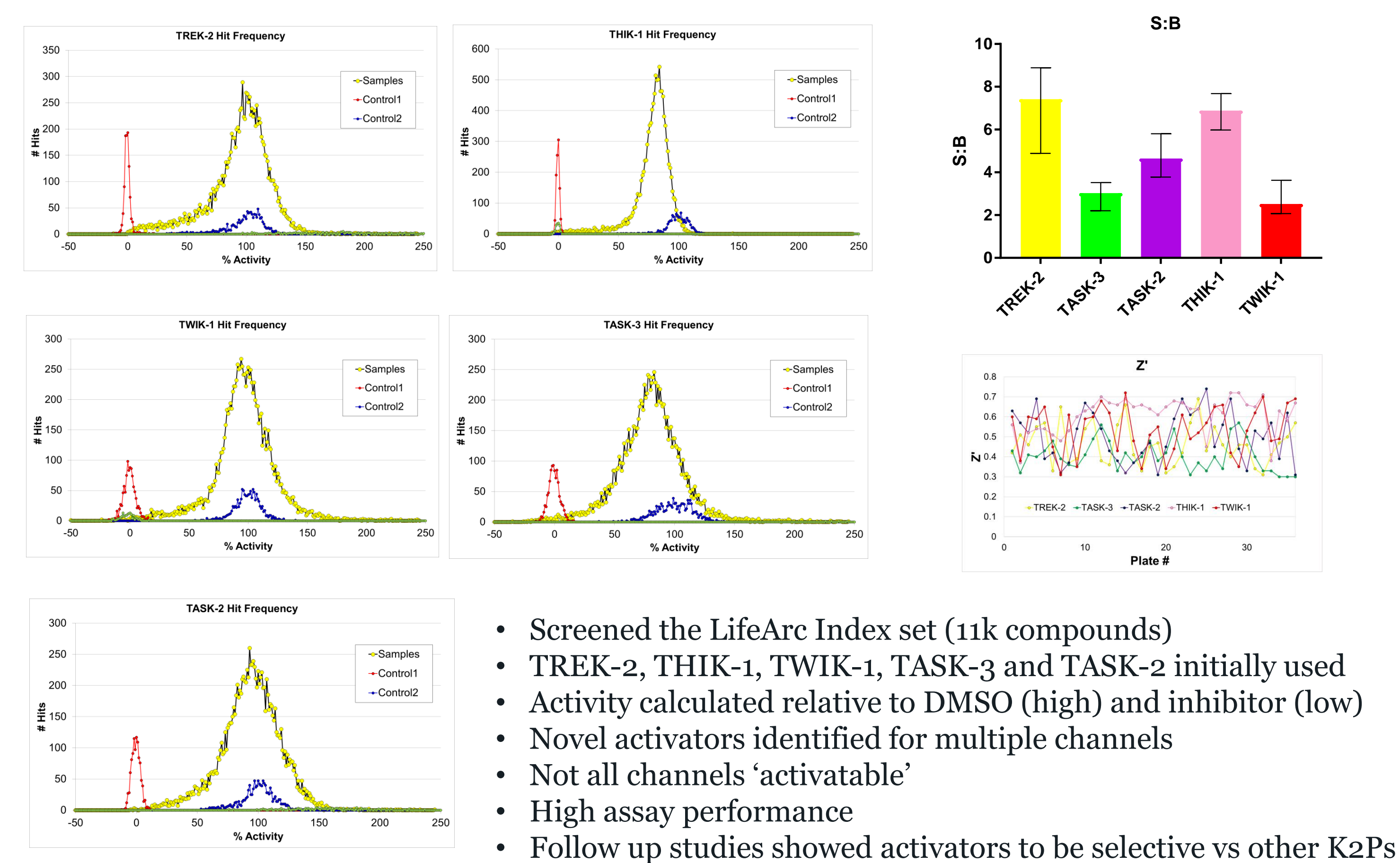
### IDENTIFYING ACTIVATORS REQUIRES BESPOKE ASSAYS/REAGENTS



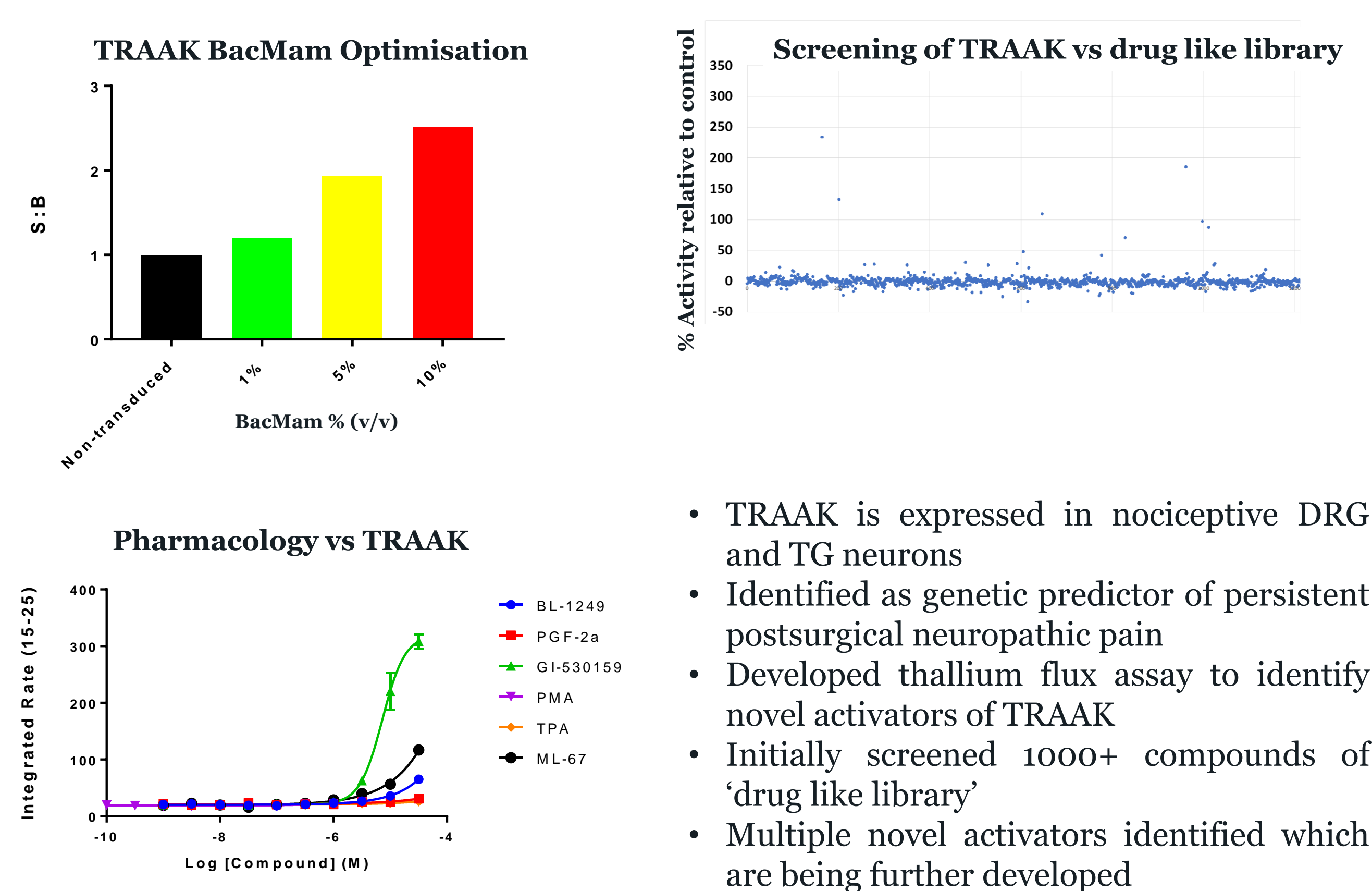
### ASSAY PROCESS AND METHODS



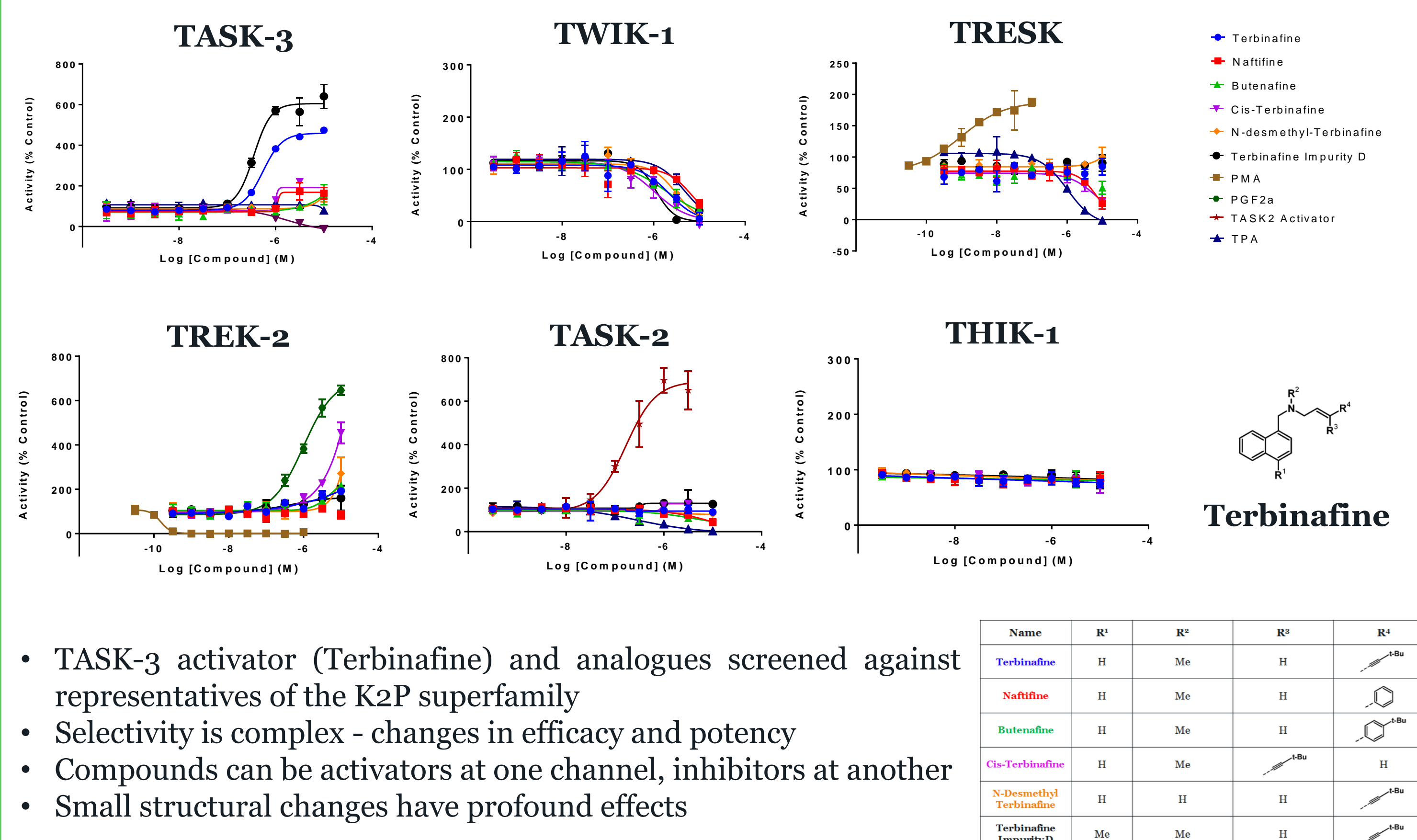
### SCREENING FOR K2P ACTIVATORS



### DEVELOPMENT OF A NOVEL TRAAK ASSAY



### SELECTIVITY OF K2P ACTIVATORS



### CONCLUSIONS

- Assays developed to assess 'ligandability' and facilitate the identification of novel activators of K2P channels
- LifeArc Index set screened and novel activators of multiple K2P channels observed
- Activators show selectivity across K2P channels but selectivity and SAR are complex
- Not all K2Ps are 'druggable' using assay system described