

# Targeting NaV1.7: How Mechanism-Based Multidisciplinary Research Drives Precision Therapy in Inherited Pain Syndromes – A Translational Approach

Marie Hoeks<sup>1,3</sup>, Angelika Lampert<sup>1,3</sup> & Ralf Hausmann<sup>2,3</sup> - on behalf of all SCN<sup>AACHEN</sup> members

<sup>1</sup>Institute of Neurophysiology, RWTH Aachen University, Aachen, Germany; <sup>2</sup>Institute of Clinical Pharmacology, RWTH Aachen University, Aachen, Germany; <sup>3</sup>Scientific Center for Neuropathic Pain Aachen, SCN<sup>AACHEN</sup>, Uniklinik RWTH Aachen University, Aachen, Germany

## Rationale and medical need

Neuropathic pain affects the quality of life of around 8% of the population, but treatment has made only limited progress over the last two decades. Thus, an innovative approach is needed to meet this medical need. At the Scientific Center for Neuropathic Pain Aachen (SCN<sup>AACHEN</sup>) we have established an interdisciplinary analysis platform/pipeline to investigate individual sodium channel variants, with a focus on NaV1.7, identified in patients of our SCN<sup>AACHEN</sup>-patient registry suffering from small fiber neuropathy. We aim to address the treatment gap and to advance neuropathic pain research by interdisciplinary and translational research on neuropathic pain mechanisms.

### Pharmacology

- High throughput patch clamp
- Multi-electrode arrays
- Repurposing

### 3D structural modeling

- Molecular dynamics
- Drug docking
- *In silico* mutagenesis

### Nociceptor profiling

- Functional assesment with microneurography on patients
- PatchSeq of pig, human and iPSC derived sensory neurons
- Neurite RNAseq
- Nav subtype select. iPSC-nocic. classification by optogenetics

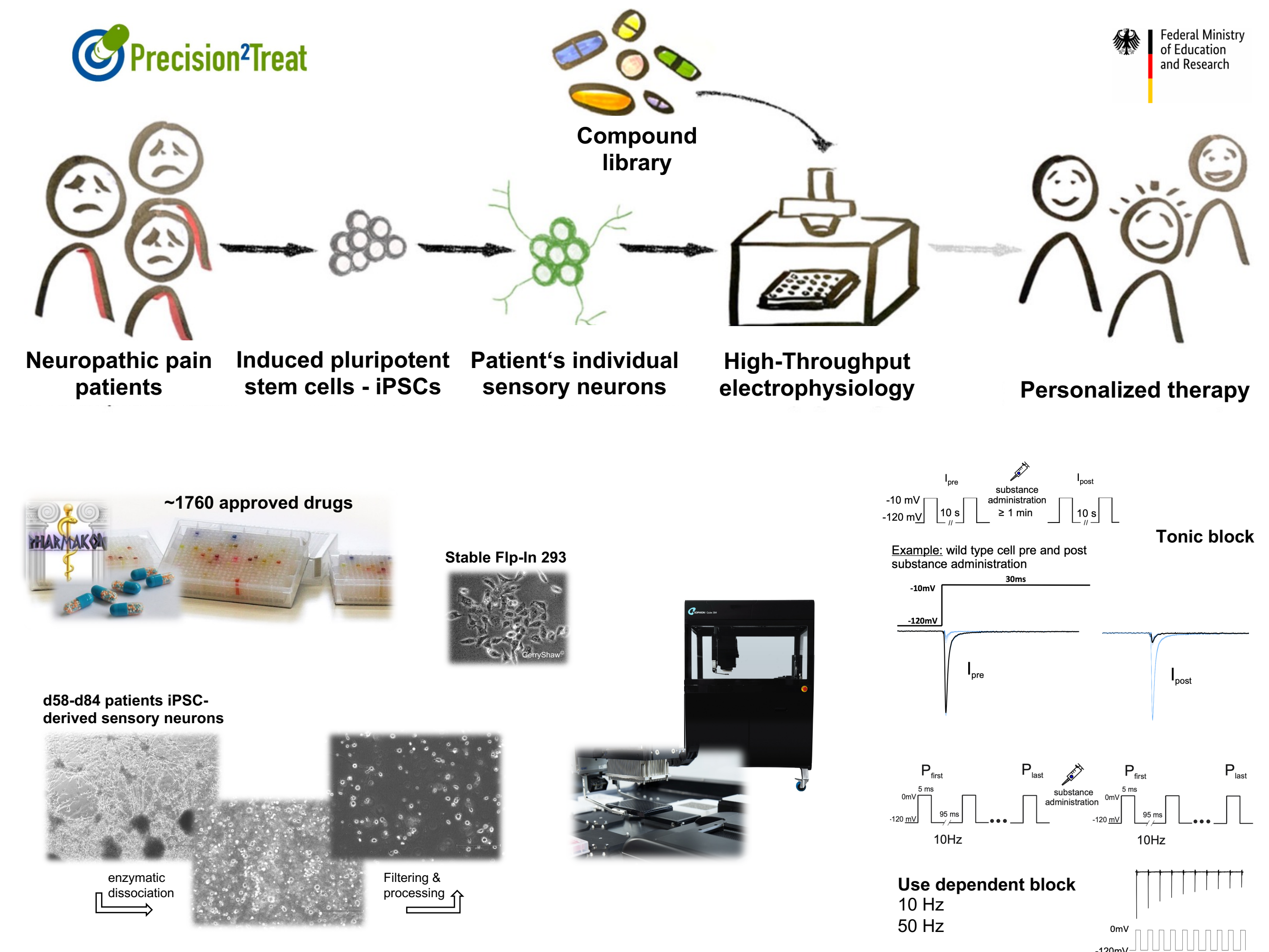
### Genetic and clinical phenotyping

- Cohort of 250+ small fiber neuropathy patients
- Genetic sequencing (WES and WGS)
- In-depth Clinical and biometric data (PainWatch)
- Brain Connectivity (fMRI)

### Digital Pain Patient

- Digital human nociceptor
- Markov Models of Nav
- Predictive models for disease progression and therapy

## Nav variant druggability testing



## Easy-to-Clone SCN9A cDNA

SCN9A plasmids are technically challenging as they frequently undergo rearrangements during propagation and result in very low yield after purification. Thus the efficiency of mutagenesis is usually low. To overcome this bottleneck a codon optimization strategy (Bertelli et al., 2018) was used to create an *easy-to-clone* SCN9A cDNA. SCN9A and SCN5A amino acid (aa) sequences were aligned and for each position in which the same aa was present, the corresponding codon of the SCN9A cDNA sequence was substituted by the SCN5A codon. For all other aa, the codon present in SCN9A was replaced by the synonymous codon being most similar to the corresponding SCN5A codon, if possible. The modified cDNA sequence was synthesized and subcloned into a modified pcDNA5.1/FRT/TO vector enabling Gateway®-cloning and FRT-mediated Flip-In® stable celline generation.

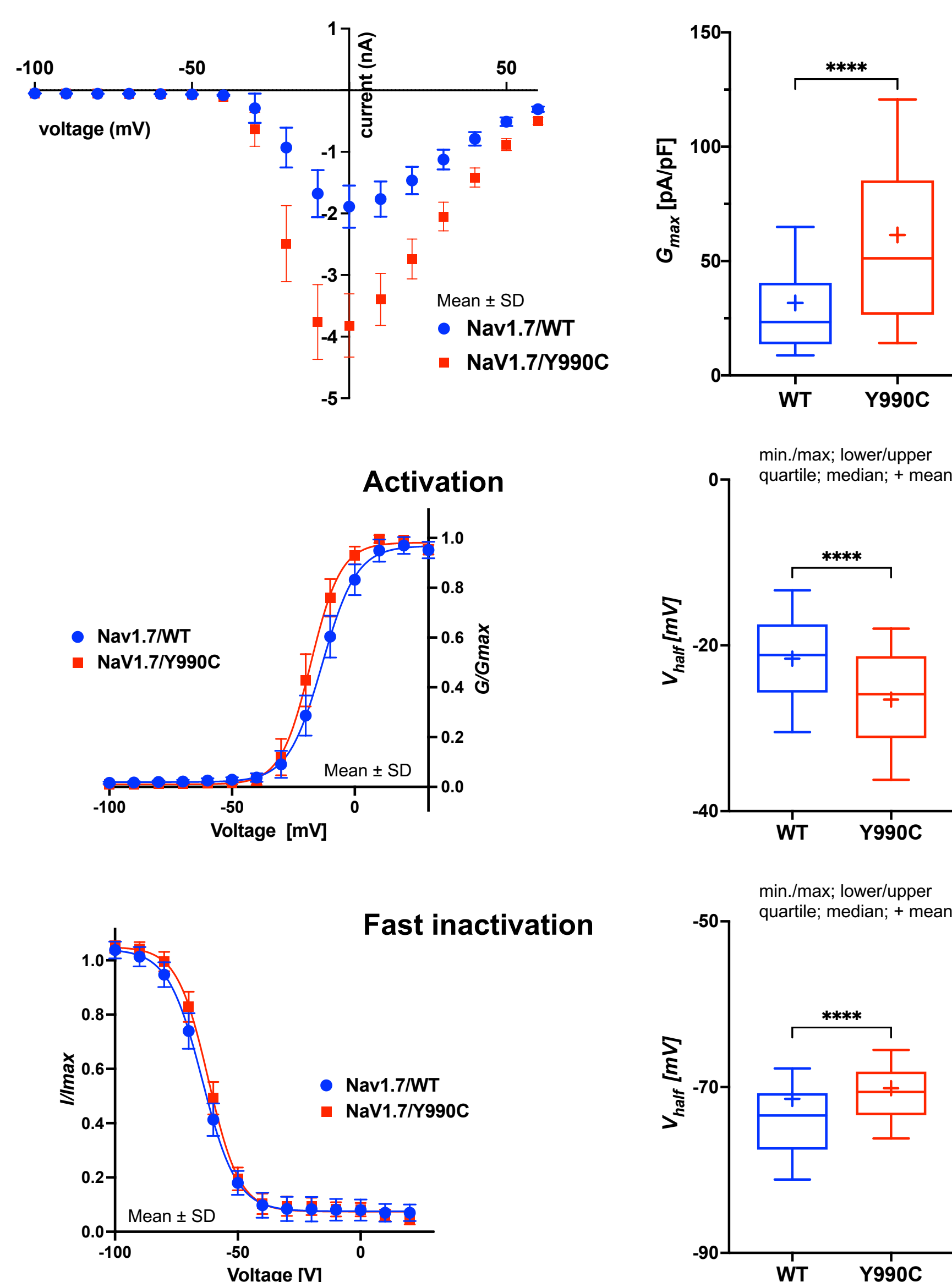
SCN5A	7P	R	G	T	S	S	F	R	R	F	T
	19cct	cgg	ggc	acc	agc	agc	ttc	cgc	agg	ttc	aca
SCN9A	6P	P	G	P	Q	S	F	V	H	F	T
	16ccc	cca	gga	cct	cag	agc	ttt	gtc	cat	ttc	aca

↓ codon optimization ↓

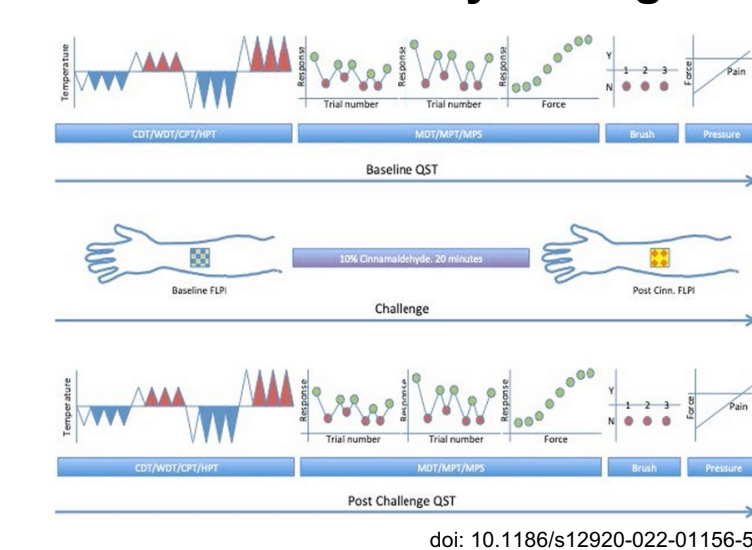
SCN9A cod.opt.	16cct	ccg	ggc	ccc	cag	agc	ttc	gtc	cat	ttc	aca
	6P	P	G	P	Q	S	F	V	H	F	T

## Biophysical Profiling of NaV1.7/Y990C

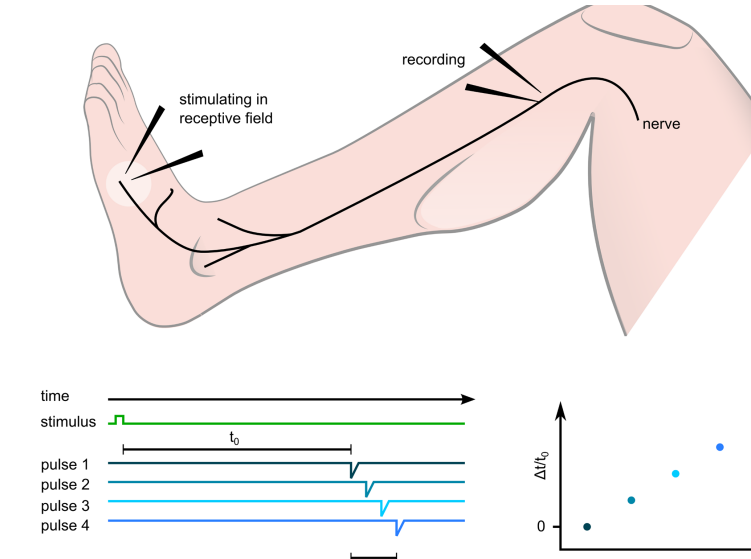
Biophysical profiling by APC of the pathogenic NaV1.7/Y990C variant associated with SFN showed a hyperpolarising shift of activation and a depolarizing shift in fast inactivation confirming a gain-of-function gating phenotype consistent with SFN. However, an initial screening of 90 approved „sodium channel“ drugs showed no different effect compared to the WT channel.



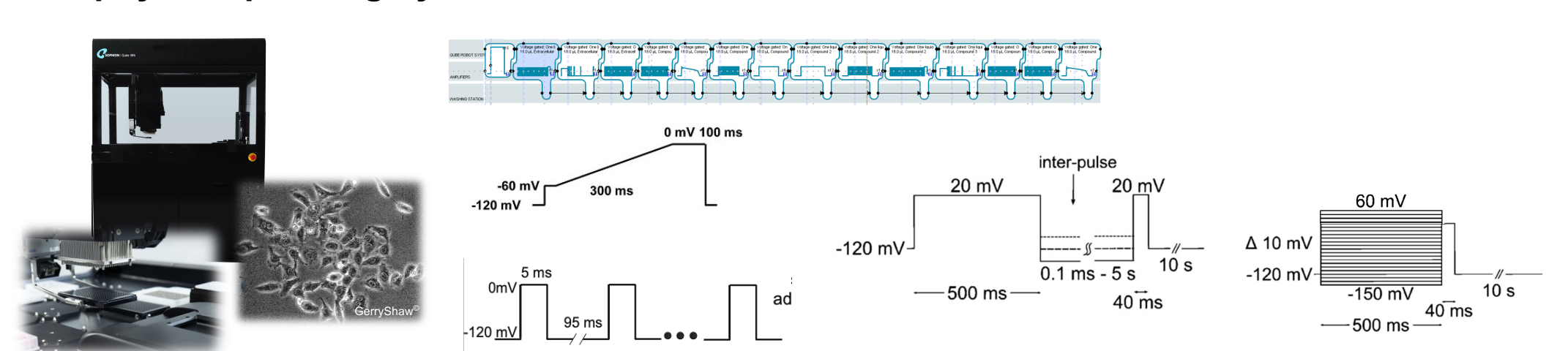
### Quantative sensory testing



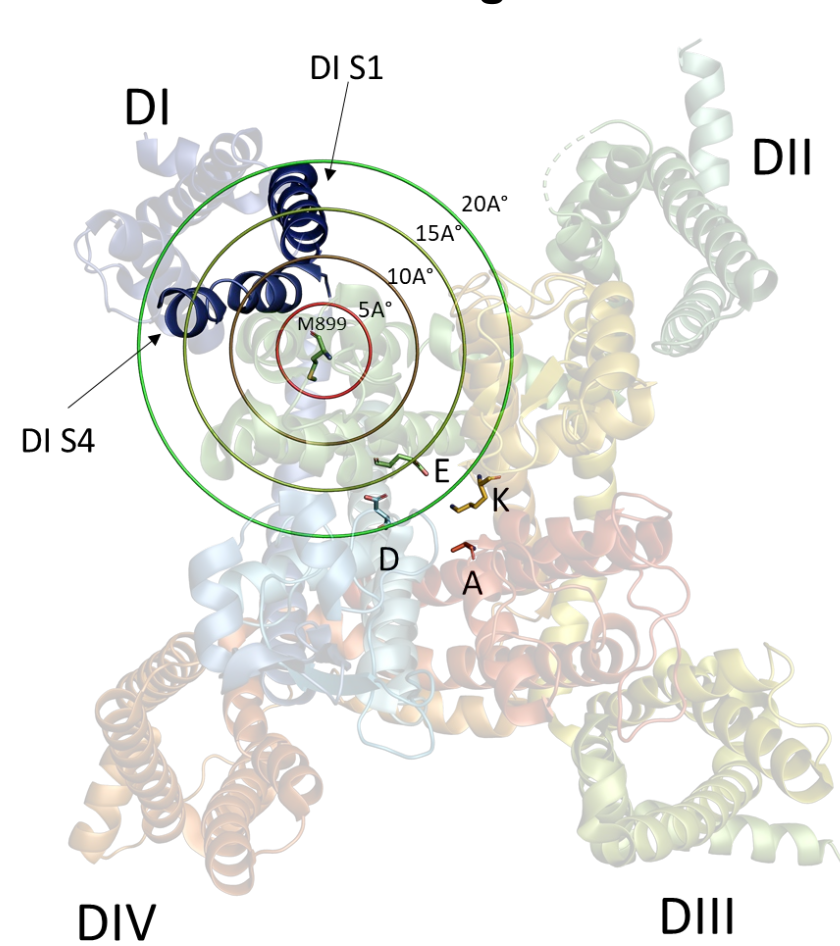
### Microneurography



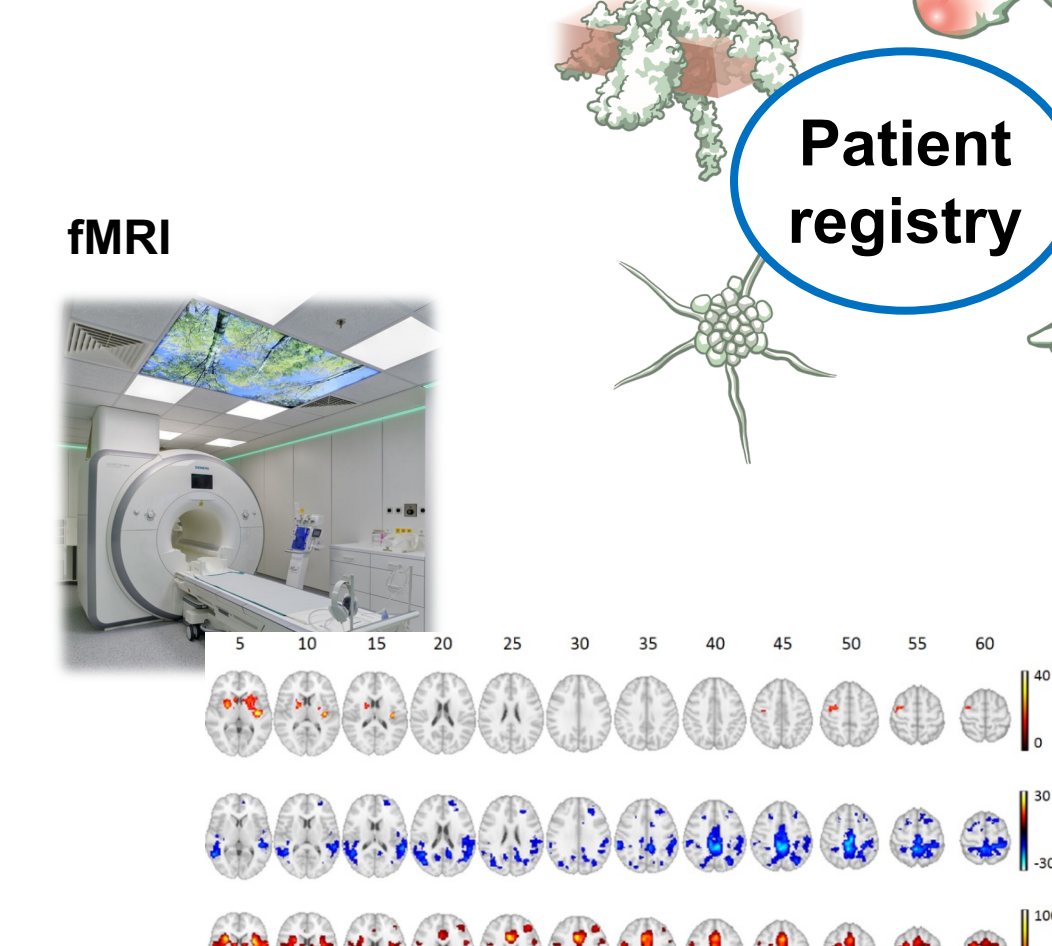
### Biophysical profiling by APC



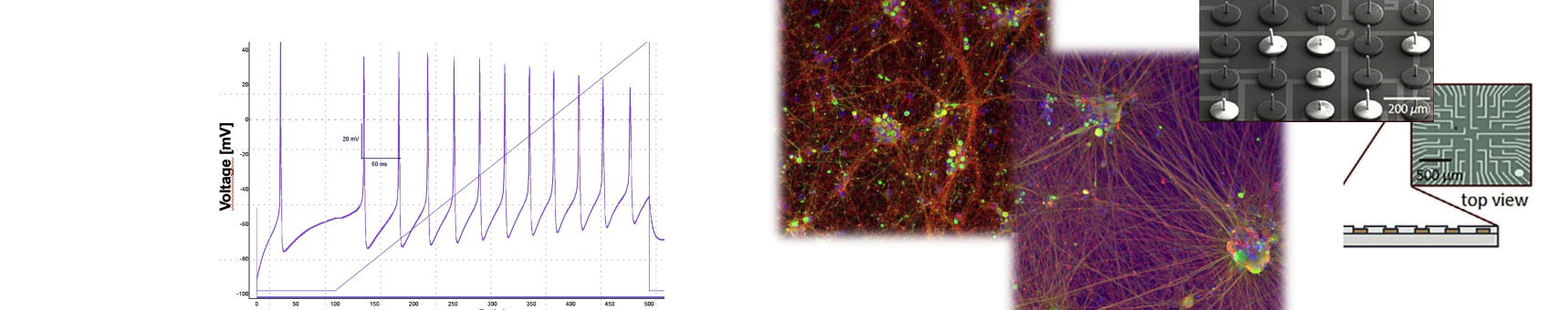
### Structural modelling & MDS



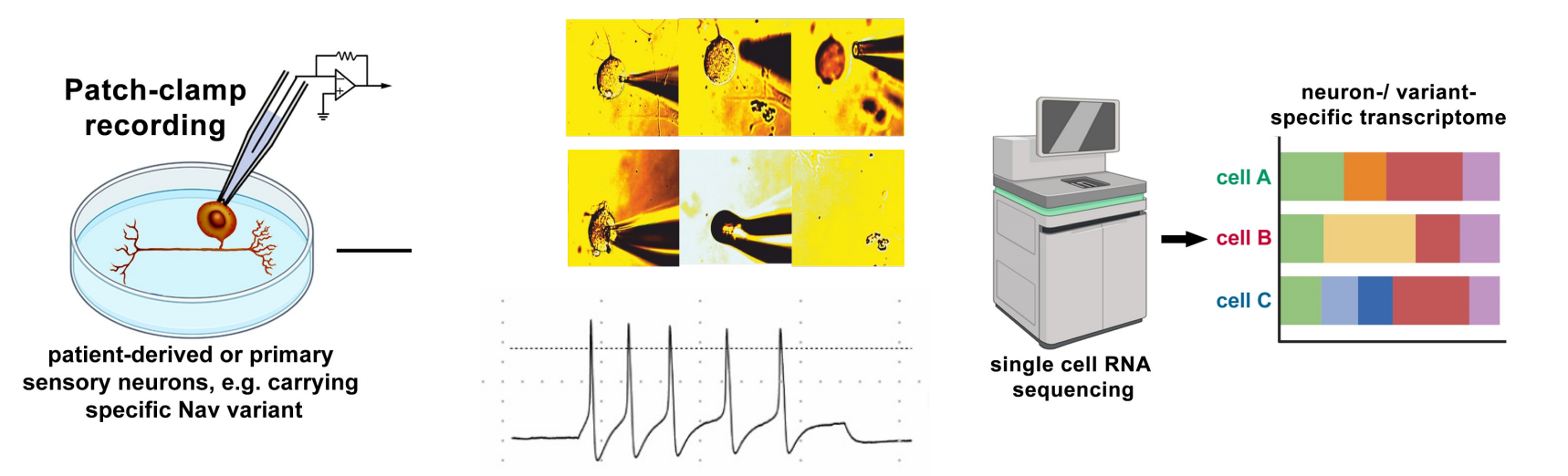
### fMRI



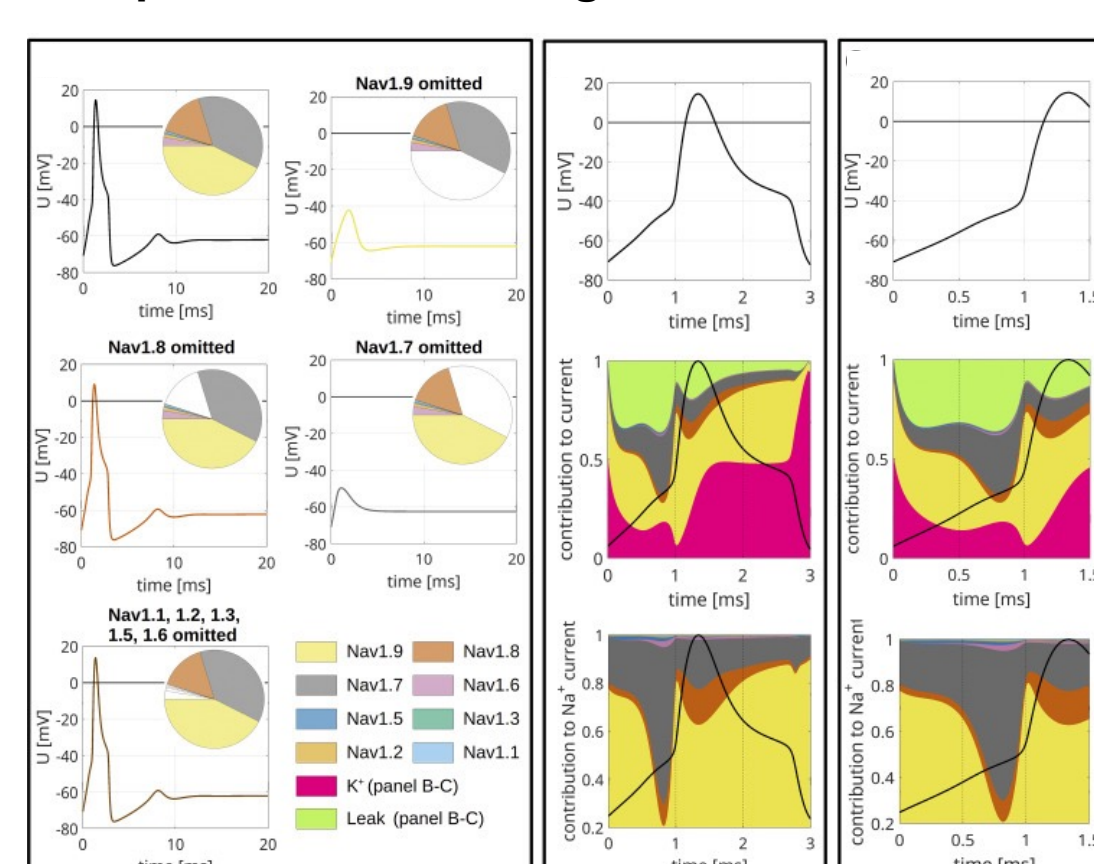
### iPSC derived sensory neurons – MEA and APC CC



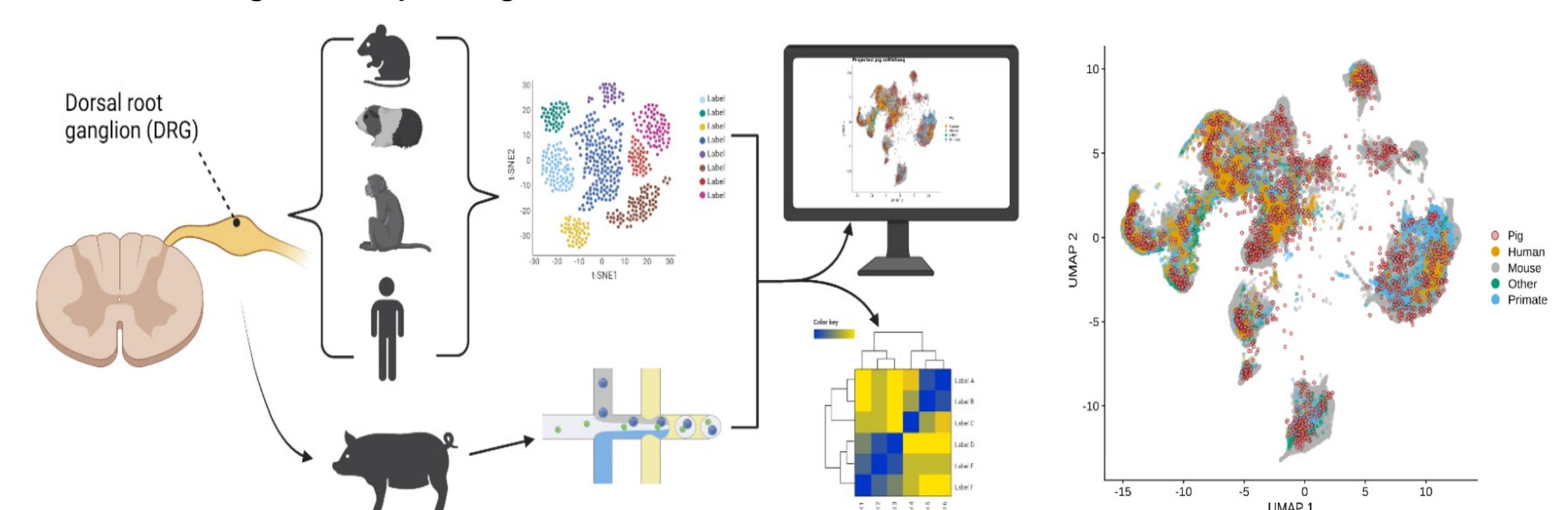
### Patch-Seq (scRNA-seq.)



### Computational Modelling



### Genome and single-cell sequencing



## Summary and Outlook

The SCN<sup>AACHEN</sup> patient registry, NGS, high-throughput electrophysiology, drug screening, and iterative modelling provides a scalable strategy for (patho)mechanistic annotation and precision pharmacology of NaV1.7 variants. It's translation potential extends beyond the example of NaV1.7/Y990C, accelerating rational therapy development for the spectrum of NaV1.7-mediated pain disorders.

### Selected References

Bertelli S et al., Gain of function of sporadic/familial hemiplegic migraine-causing SCN1A mutations: Use of an optimized cDNA. *Cephalalgia*. 2019 Apr;39(4):477-488. Köster PA et al., Nociceptor sodium channels shape subthreshold phase, upstroke, and shoulder of action potentials. *J Gen Physiol*. 2025;157:e202313526. Körner J et al., Molecular architecture of human dermal sleeping nociceptors. *bioRxiv PREPRINT* doi:10.1101/2024.12.20.629638.

### Funding