

# Development and Evaluation of Novel Solution Pairs to Enhance Seal Resistance in Automated Patch Clamp Experiments

Catherine Hodgson<sup>1</sup>, Edward Stevens<sup>1</sup>, Andrew James<sup>1</sup>, Anissa Bara<sup>2</sup>, Beatrice Badone<sup>2</sup>, Damian Bell<sup>2</sup>, Daniel Sauter<sup>2</sup>, Gary Clark<sup>1</sup>

<sup>1</sup>Metrion Biosciences Ltd, Granta Centre, Granta Park, Cambridge, CB21 6AL, UK

<sup>2</sup>Sophion Bioscience A/S, Industriparken 39, 2750 Ballerup, Denmark

## Introduction

Gigaohm seals, or 'gigaseals', are crucial for patch clamp electrophysiology, ensuring excellent electrical access to cells to enable high-quality recordings. These seals form through chemical bonds and electrostatic forces between the cell membrane and the glass pipette in manual patch clamp, or between the cell membrane and chip substrate in planar patch clamp. Planar patch clamp often employs 'seal enhancers' to increase seal resistances, with CaF<sub>2</sub> being the most commonly used. It is hypothesized that high extracellular Ca<sup>2+</sup> and intracellular F<sup>-</sup> concentrations lead to CaF<sub>2</sub> precipitate formation at the solution interface, promoting seal formation. However, CaF<sub>2</sub> as a seal enhancer has limitations. F<sup>-</sup> interacts with various internal components such as protein kinase A, adenylate cyclase, and lipid phosphatases, which can affect the biophysical properties of some ion channels. Additionally, F<sup>-</sup> is not ideal when recording from Ca<sup>2+</sup>-activated ion channels due to the resulting unknown concentrations of free intracellular Ca<sup>2+</sup>.

In an effort to overcome these limitations, Sophion developed new solution pairs in 2017 that foster seal formation (Patent: WO2018100206A1). Building on this technology, Metrion and Sophion collaborated to further determine whether other insoluble salts can act as seal enhancers and how these solution pairs affect the biophysical properties and pharmacology of the investigated ion channels.

## Conclusion

- BaSO<sub>4</sub> was identified as an equivalent seal enhancer to CaF<sub>2</sub>.
- The two solution pairs were characterized across two ion channels: hNa<sub>v</sub>1.5 and hCa<sub>v</sub>1.2.
- Intracellular F<sup>-</sup> caused depolarizing shifts in the voltage dependence of inactivation of hNa<sub>v</sub>1.5 where no such effects were observed with SO<sub>4</sub><sup>2-</sup> in the intracellular solution.
- No difference in pharmacological effects of inhibitory compounds against hNa<sub>v</sub>1.5 or hCa<sub>v</sub>1.2 was observed between the two solution pairs, CaF<sub>2</sub> and BaSO<sub>4</sub>.
- BaSO<sub>4</sub> is well-suited as a seal enhancer for recording from non-K<sup>+</sup>-conducting Ca<sup>2+</sup>-activated channels, such as TME-M16A. In particular as BaSO<sub>4</sub> allows more accurate estimation of free intracellular Ca<sup>2+</sup> concentration.

## Material and methods

Experiments were conducted at 22°C using a Sophion Bioscience Qube 384 with QChip 384 (single hole) and QChip 384X (multihole) consumables.

Analysis was conducted using Sophion Analyzer v9.0.42 and GraphPad Prism v10.2.2.

CHO-hNa<sub>v</sub>1.5 and HEK293-hCa<sub>v</sub>1.2 cell lines were provided by Metrion Biosciences.

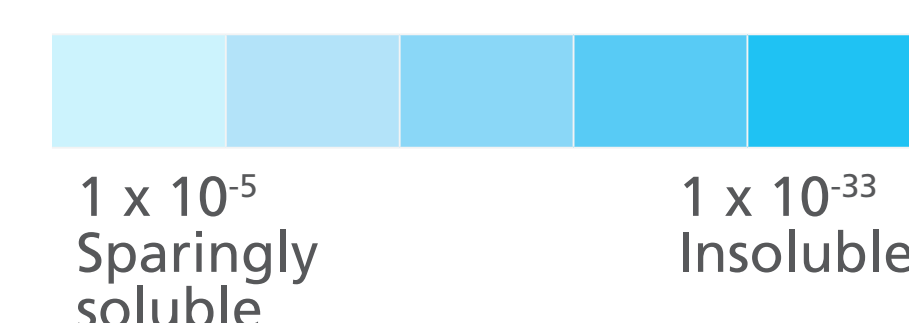
All compounds were tested at: 0.001, 0.01, 0.1, 1, 10 and 100 µM.

## Results

**Table 1: Correlation between salt pair solubility product constants (K<sub>sp</sub>) and gigaseal formation on Qube 384.** No correlation was found between the solubility product constants (K<sub>sp</sub>) of Ca<sup>2+</sup>, Ba<sup>2+</sup> and Sr<sup>2+</sup> salts (A) and their ability to foster gigaseal formation (B). Despite the PO<sub>4</sub><sup>3-</sup> salts having very low K<sub>sp</sub> values and SrCO<sub>3</sub> having a similar K<sub>sp</sub> value to CaF<sub>2</sub> and BaSO<sub>4</sub>, these salts failed to produce gigaohm seals. Moderate seal resistances with PO<sub>4</sub><sup>3-</sup> salts were transient and unstable. Median resistances calculated from 24 cells per salt pair.

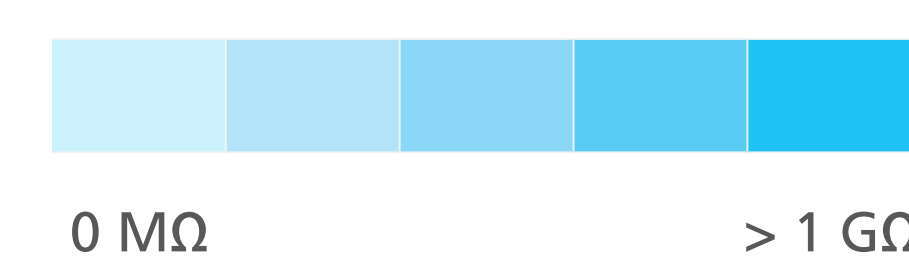
**A.**

K <sub>sp</sub>	Solubility product constant (K <sub>sp</sub> ) at 25°C; pH7.0				
	F <sup>-</sup>	SO <sub>4</sub> <sup>2-</sup>	CO <sub>3</sub> <sup>2-</sup>	PO <sub>4</sub> <sup>3-</sup>	C <sub>2</sub> O <sub>4</sub> <sup>2-</sup>
Ca <sup>2+</sup>	1.5 × 10 <sup>-5</sup>	1.1 × 10 <sup>-10</sup>	3.4 × 10 <sup>-9</sup>	1.0 × 10 <sup>-13</sup>	2.3 × 10 <sup>-11</sup>
Ba <sup>2+</sup>	1.5 × 10 <sup>-5</sup>	1.1 × 10 <sup>-10</sup>	3.4 × 10 <sup>-9</sup>	1.0 × 10 <sup>-13</sup>	2.3 × 10 <sup>-11</sup>
Sr <sup>2+</sup>	1.5 × 10 <sup>-5</sup>	1.1 × 10 <sup>-10</sup>	3.4 × 10 <sup>-9</sup>	1.0 × 10 <sup>-13</sup>	2.3 × 10 <sup>-11</sup>



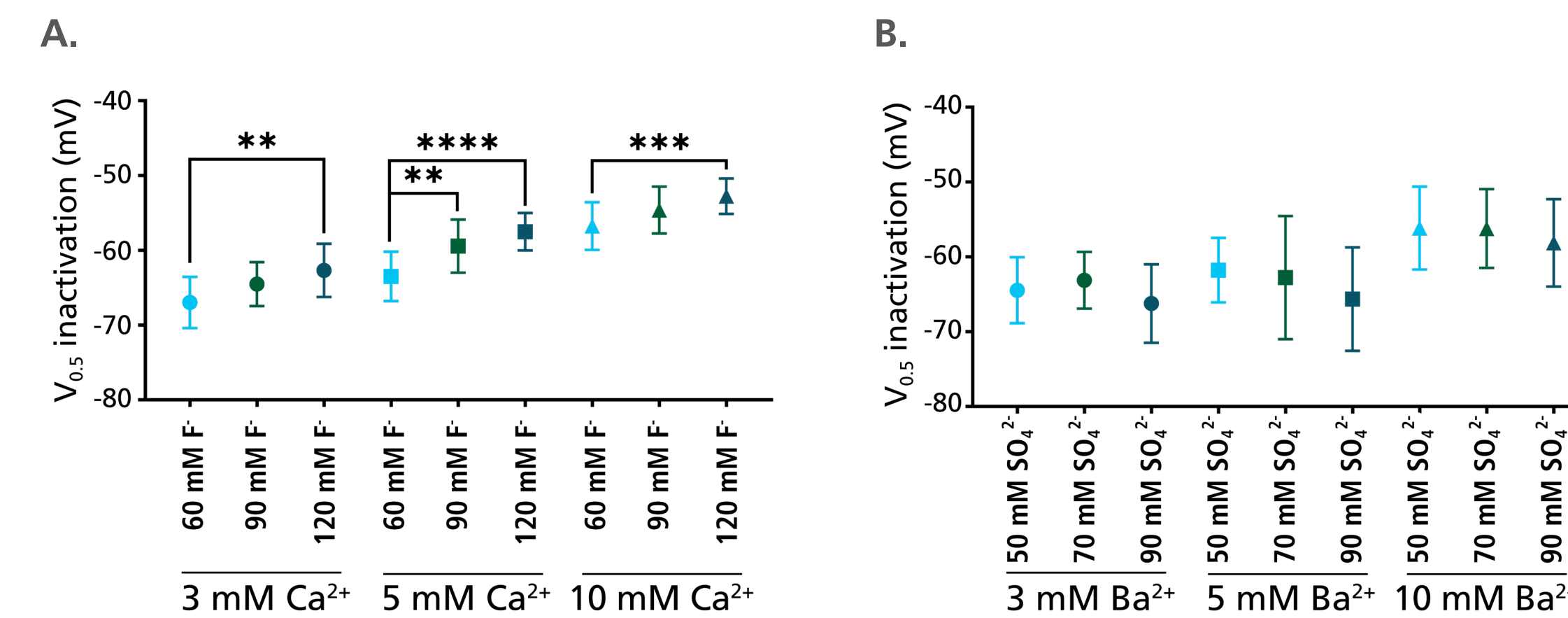
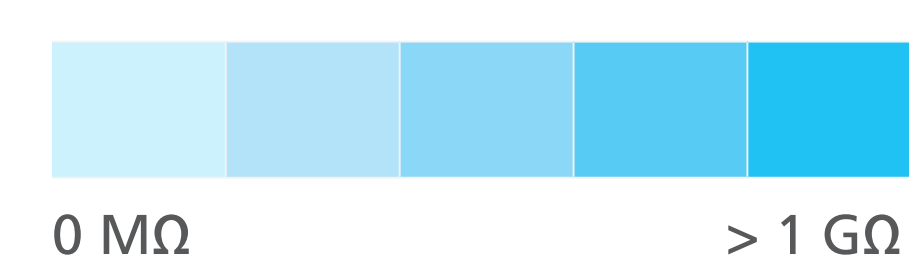
**B.**

GΩ	Median R <sub>membrane</sub> (GΩ)				
	120 mM F <sup>-</sup>	90 mM SO <sub>4</sub> <sup>2-</sup>	90 mM CO <sub>3</sub> <sup>2-</sup>	70 mM PO <sub>4</sub> <sup>3-</sup>	90 mM C <sub>2</sub> O <sub>4</sub> <sup>2-</sup>
10 mM Ca <sup>2+</sup>	0.5	0.5	0.5	0.5	0.5
10 mM Ba <sup>2+</sup>	0.5	0.5	0.5	0.5	0.5
10 mM Sr <sup>2+</sup>	0.5	0.5	0.5	0.5	0.5

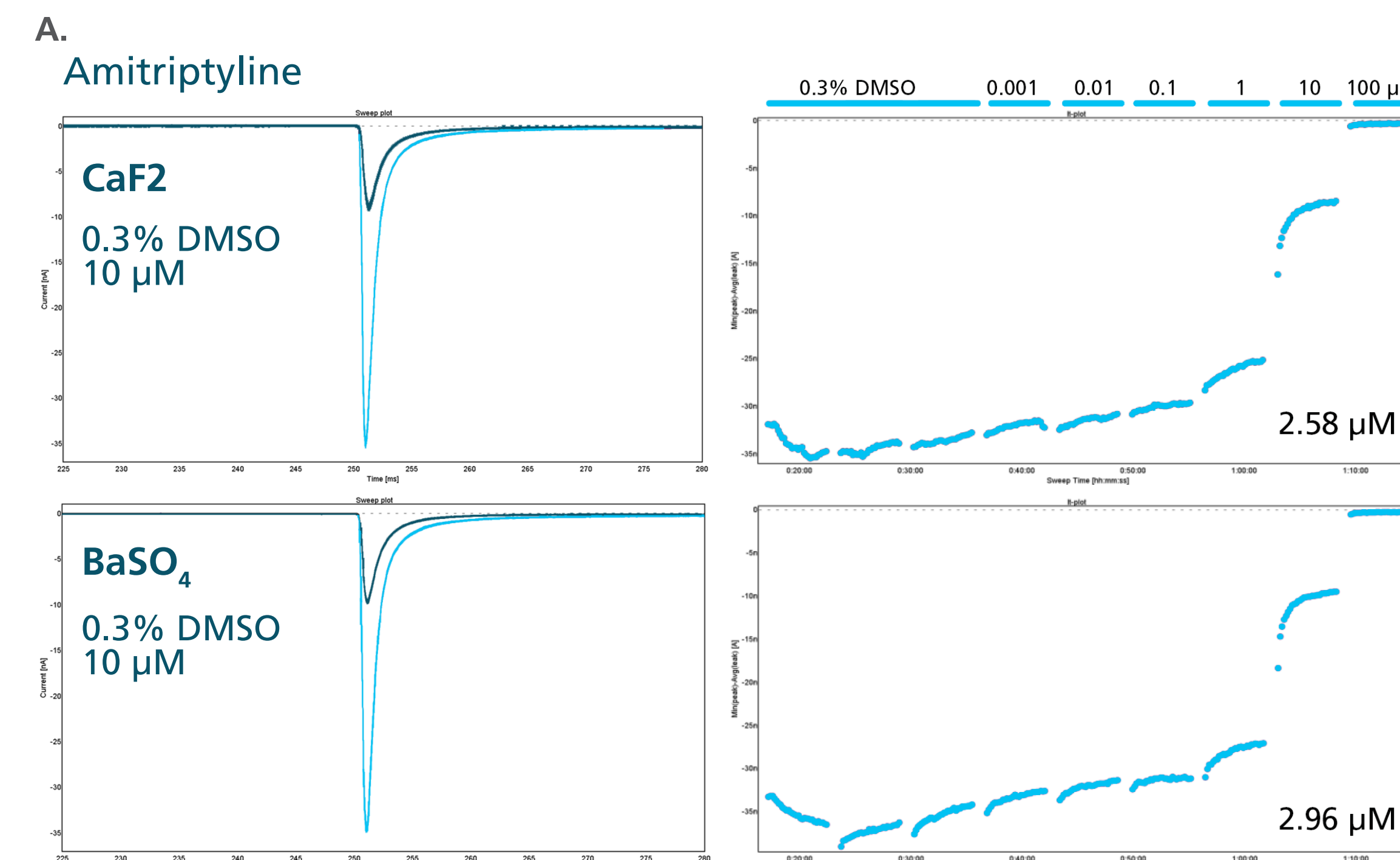


**Table 2: Resistances of seals formed using various concentrations of extracellular Ba<sup>2+</sup> and intracellular SO<sub>4</sub><sup>2-</sup>.** Gigaseals only formed with ≥ 3 mM Ba<sup>2+</sup> and in the presence of SO<sub>4</sub><sup>2-</sup>. Median resistances calculated from 24 or 48 cells per condition.

CHO-hNa <sub>v</sub> 1.5	Median R <sub>membrane</sub> (GΩ)			
	0 mM SO <sub>4</sub> <sup>2-</sup>	50 mM SO <sub>4</sub> <sup>2-</sup>	70 mM SO <sub>4</sub> <sup>2-</sup>	90 mM SO <sub>4</sub> <sup>2-</sup>
Single-hole experiments	0	0	0	0
Ba <sup>2+</sup> (mM)	0	0	0	0
	0.01	0	0	0
	0.03	0	0	0
	0.1	0	0	0
	0.3	0	0	0
	1	0	0	0
	3	0	0	0
	5	0	0	0
	10	0	0	0



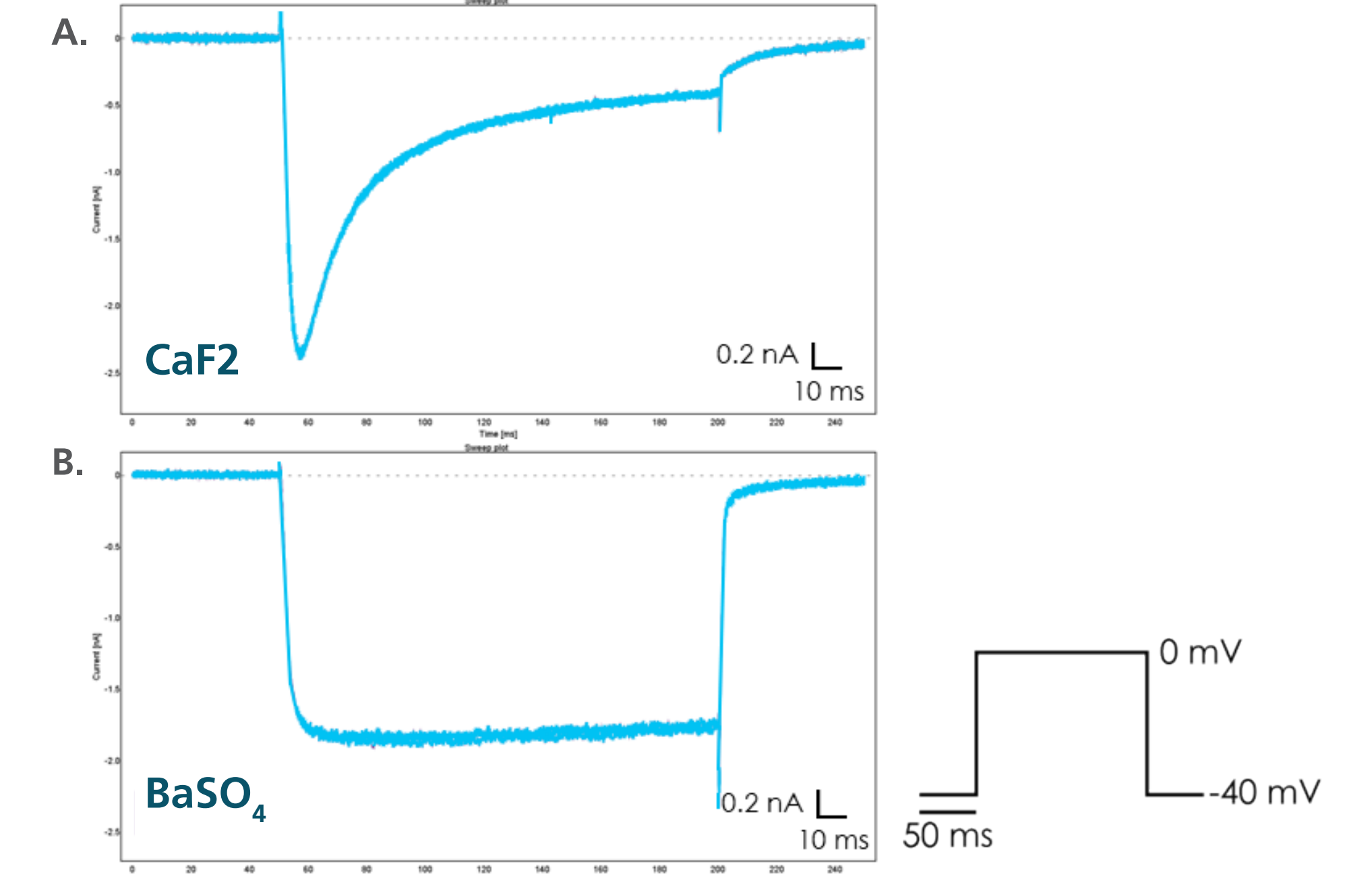
**Fig. 1: Effects of CaF<sub>2</sub> and BaSO<sub>4</sub> on hNa<sub>v</sub>1.5 channel biophysics.** hNa<sub>v</sub>1.5 V<sub>0.5</sub> inactivation with different cation and anion concentrations (mean ± S.D.; N ≥ 11). Increasing concentrations of intracellular F<sup>-</sup> caused a depolarizing shift in V<sub>0.5</sub> inactivation (A). In contrast, increasing concentrations of SO<sub>4</sub><sup>2-</sup> had no effect on hNa<sub>v</sub>1.5 V<sub>0.5</sub> inactivation (B). One-way ANOVAs conducted within each cation group followed by Tukey's Honestly Significant Difference post-hoc tests: \*\* = p < .01; \*\*\* = p < .001; \*\*\*\* = p < .0001.



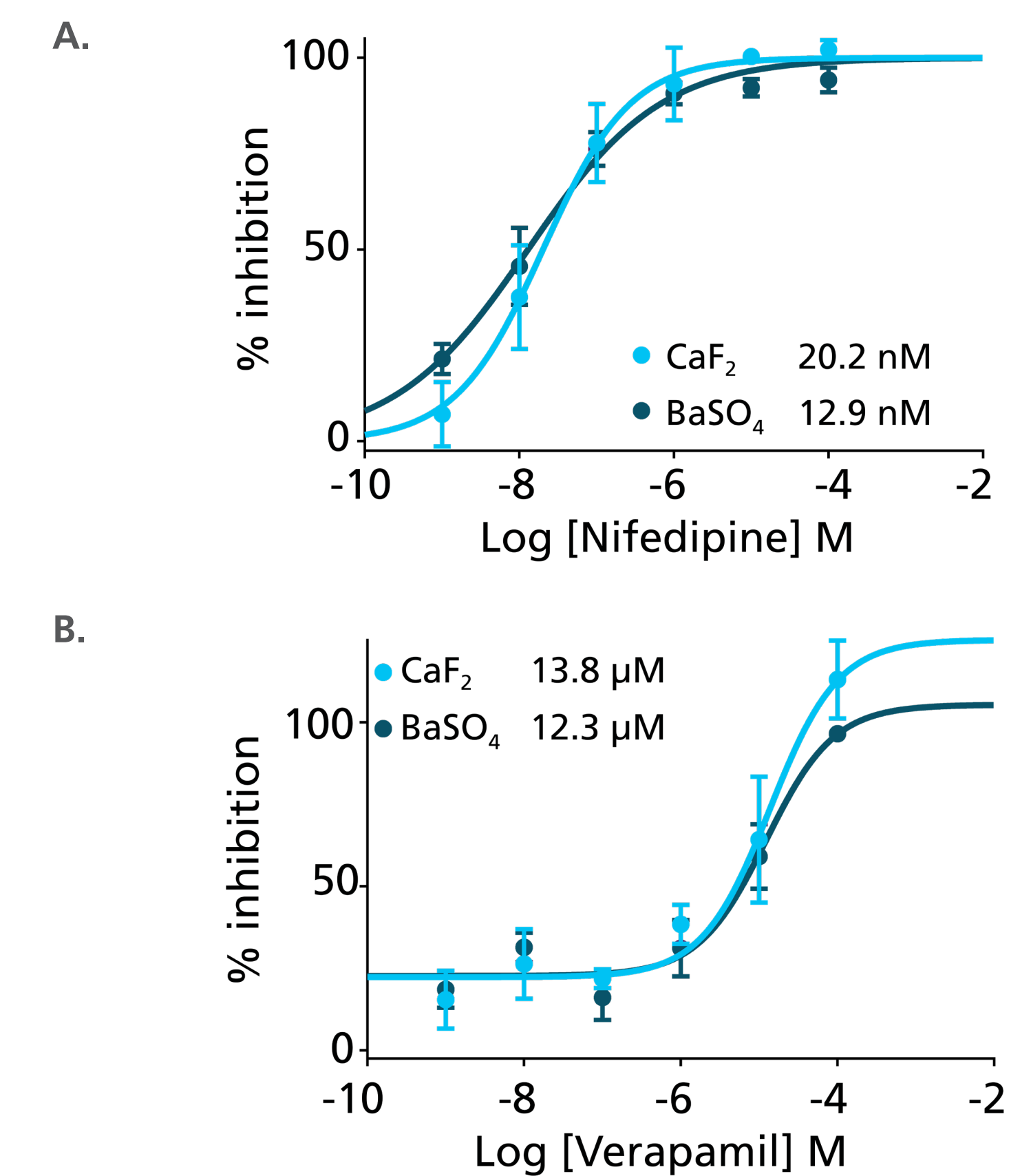
**B.**

CHO- hNa <sub>v</sub> 1.5	IC <sub>50</sub> (µM)	
	3 mM Ca <sup>2+</sup> + 120 mM F <sup>-</sup>	50 mM Ba <sup>2+</sup> + 70 mM SO <sub>4</sub> <sup>2-</sup>
Multihole experiments		
Amitriptyline	2.6	3.0
Tetracaine	3.7	5.5
Bepridil	12.0	14.0
Verapamil	16.4	18.0
Dapoxetine	9.1	11.6
Lidocaine	>100	>100
Phenytoin	>100	>100

**Fig. 2: CaF<sub>2</sub> versus BaSO<sub>4</sub> - hNa<sub>v</sub>1.5 pharmacology.** (A) Representative sweep plots (left) and current-time (I-t) plots (right) for hNa<sub>v</sub>1.5 inhibition by amitriptyline. There was no difference in cumulative inhibition of hNa<sub>v</sub>1.5 by increasing concentrations of amitriptyline between CaF<sub>2</sub> and BaSO<sub>4</sub>. (B) Screening of a range of inhibitory compounds showed no difference in hNa<sub>v</sub>1.5 pharmacology between CaF<sub>2</sub> and BaSO<sub>4</sub>. Concentration-response curves for amitriptyline against hNa<sub>v</sub>1.5 using CaF<sub>2</sub> or BaSO<sub>4</sub> as the seal enhancer (mean ± S.D.; N = 12 wells per concentration for CaF<sub>2</sub>, N = 8 wells per concentration for BaSO<sub>4</sub>).



**Fig. 3: CaF<sub>2</sub> versus BaSO<sub>4</sub> - hCa<sub>v</sub>1.2 kinetics.** hCa<sub>v</sub>1.2 exhibits Ca<sup>2+</sup>-dependent inactivation when CaF<sub>2</sub> is used as the seal enhancer (A). BaSO<sub>4</sub> as the seal enhancer (using Ba<sup>2+</sup> as a surrogate carrier ion) (B) confers loss of the Ca<sup>2+</sup>-dependent inactivation of hCa<sub>v</sub>1.2 observed with CaF<sub>2</sub>. Example sweep plots derived from Sophion Analyzer v9.0.42.



**Fig. 4: CaF<sub>2</sub> versus BaSO<sub>4</sub> - hCa<sub>v</sub>1.2 pharmacology.** Mean ± S.D. concentration-response curves for two common inhibitors against hCa<sub>v</sub>1.2, nifedipine (A) and verapamil (B) (CaF<sub>2</sub>: N = 2-5 wells per concentration; BaSO<sub>4</sub>: N = 6-12 wells per concentration). Compound potencies (IC<sub>50</sub> values) did not differ between CaF<sub>2</sub> and BaSO<sub>4</sub>.

## Sophion Qube 384



Sophion's automated patch clamp system **Qube 384** for high performance and high-throughput ion channel characterization and screening.

