# Simultaneous Measurement of Cardiac Nav1.5 Peak and Late Currents in an Automated QPatch Platform

### ABSTRACT

High throughput in vitro profile of the cardiac Nav1.5 peak sodium current  $(I_{Na})$  is widely used in cardiac safety screening. However, there is no standardized high throughput method to measure late I<sub>Na</sub>. A variety of differing protocols used across industry and academia may contribute to variation in the data.

The objectives of this study were to (1) assess pharmacology and biophysical properties of veratridine- and ATX-II-induced late INa together with that induced by a mutation in the channel ( $\Delta$ KPQ-Nav1.5), (2) develop a protocol to allow simultaneous measurement of both peak and late INa under a single protocol using automated QPtach system.

The planar patch clamp technique (QPatch) was applied to record the peak and late  $I_{Na}$  from the Nav1.5 channel (human SCN5A gene) or  $\Delta$ KPQ-Nav1.5 mutant channel expressed in mammalian cells.

When measured at the maximal response during the ramp of the voltage protocol, the  $\Delta KPQ-Nav1.5$ mutant produced a small late  $I_{Na}$  (41.9 ± 5.4 pA). Veratridine and ATX-II induced concentration-dependent increases in the late  $I_{Na}$ . The amplitude of late  $I_{Na}$  were 1162.2 ± 258.5 pA and 392.4 ± 71.3 pA in the presences of 100 µM veratridine and 100 nM ATX-II, respectively. Veratridine inhibited the peak I<sub>Na</sub>  $(IC_{50} = 84.1 \pm 10.5 \mu M)$  and altered the biophysical properties of the I<sub>Na</sub>. ATX-II showed minimal effects on the peak I<sub>Na</sub> and preserved the biophysical properties of the I<sub>Na</sub>. In the presence of 100 nM ATX-II, potencies of 25 clinical I<sub>Na</sub> blockers on peak and late I<sub>Na</sub> were characterized. In addition, the IC<sub>50</sub> values of these clinical  $I_{Na}$  blockers on peak  $I_{Na}$  correlated well with and without ATX-II. The results also demonstrated that the potency of a compound blocking late I<sub>Na</sub> could be either overestimated or underestimated if the late  $I_{Na}$  was measured at the end of the depolarizing pulse versus during the ramp.

In conclusion, in the presence of ATX-II, both peak and late I<sub>Na</sub> could be assessed simultaneously under a single protocol. Our results suggest that late I<sub>Na</sub> may be best assessed using the maximum response obtained during the ramp after 200 ms depolarizing pulse at 40 mV.

### **METHODS**

### Automated Electrophysiology

WT Nav1.5 and  $\Delta$ KPQ-Nav1.5 mutant currents were recorded on the QPatch HTX automated patch clamp platform (Sophion). For WT Nav1.5 channels, ATX-II (Alomone labs) or veratridine (Sigma) was used as a chemical modifier to induce the late I<sub>Na</sub>. Test compounds were evaluated in the continued presence of these chemical modifiers. Test compounds were purchased from Sigma-Aldrich (St Louis, MO).

#### **Recombinant Cell Lines**

Wild type Nav1.5 channels were expressed in CHO cells (Charles Rivers, Cleveland, OH) and ∆KPQ-Nav1.5 mutant channels were expressed in HEK 293 cells (Icagen, Durham, NC).

#### Solutions

Extracellular solution (in mM):137 NaCl, 1.8 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>, 4 KCl, 10 HEPES, 10 Glucose, pH=7.4 (NaOH), ~295 mOsm. Intracellular solution (in mM): 92 CsF, 55 CsCl, 10 HEPES, 10 EGTA, 2 MgCl<sub>2</sub>, 2 MgATP pH=7.2 (KOH), ~300 mOsm.

#### Data analysis

Analyses of current (I) / Voltage (V)-relationships and concentration-dependent drug effects (Hill fit and IC<sub>50</sub>) were performed using QPatch Assay Software (Sophion, Denmark) and Graphpad Prism 7.0.4 (GraphPad Software, La Jolla, CA).

## WORLDWIDE RESEARCH & DEVELOPMENT

### RESULTS





Donglin Guo, Sunny Sun and Stephen Jenkinson Drug Safety Research & Development, La Jolla CA, USA

		IC <sub>ε0</sub> (μM)		
	Late I <sub>Na</sub> with ATX-II	Peak I <sub>Na</sub> with ATX-II	Peak I <sub>na</sub> withou	t ATX-II
5-OH Propafenone	1.3 ± 0.1 (5)	0.8 ± 0.1 (9)	0.4 ± 0.0	) (9)
Amitriptyline	2.3 ± 0.2 (11)	2.7 ± 0.4 (12)	1.1 ± 0.2	2 (7)
Bupivacaine	2.1 ± 0.3 (8)	3.8 ± 0.8 (9)	4.5 ± 0.5	5 (9)
Carbamazepine	25.4 ± 2.1 (9)	173.6 ± 34.8 (7)	76.3 ± 10	.0 (7)
Clomipramine	4.1 ± 0.6 (11)	6.7 ± 0.8 (12)	7.2 ± 0.7	' (9)
Desipramine	2.5 ± 0.3 (9)	2.8 ± 0.5 (9)	2.8 ± 0.4	I (12)
Disopyramide	89.1 ± 5.0 (8)	84.3 ± 10.3 (9)	102.6 ± 25	.5 (5)
Flecainide	1.9 ± 0.3 (6)	1.9 ± 0.3 (5)	1.8 ± 0.1	(8)
Imipramine	3.6 ± 0.5 (9)	3.8 ± 0.6 (12)	1.2 ± 0.1	(8)
Lamotrigine	34.5 ± 3.2 (8)	36.3 ± 6.8 (9)	31.3 ± 3.6	5 (7)
Maprotiline	1.8 ± 0.2 (8)	1.9 ± 0.3 (7)	1.6 ± 0.2	2 (9)
Mesoridazine	3.6 ± 0.3 (6)	2.8 ± 0.5 (6)	3.1 ± 0.3	3 (9)
Mexiletine	12.2 ± 1.1 (12)	23.9 ± 3.3 (10)	21.9 ± 2.6	ኝ (10)
Nortriptyline	3.3 ± 0.6 (7)	2.9 ± 0.5 (10)	1.7 ± 0.1	(11)
Procainamide	554.4 ± 85.2 (9)	515.9 ± 103.8 (5)	487.5 ± 54	.2 (12)
Procaine	82.5 ± 7.1 (6)	81.7 ± 6.8 (5)	91.7 ± 7.3	3 (12)
Propafenone	1.9 ± 0.2 (9)	2.4 ± 0.2 (12)	1.4 ± 0.1	(9)
Quinidine	16.8 ± 1.4 (8)	18.4 ± 0.7 (6)	14.9 ± 2.4	ł (9)
Quinine	11.4 ± 1.5 (8)	19.5 ± 2.5 (9)	19.0 ± 3.1	(12)
Ranolazine	16.7 ± 1.7 (5)	75.0 ± 14.6 (5)	79.5 ± 5.5	5 (12)
Risperidone	26.2 ± 5.9 (10)	31.8 ± 6.4 (10)	14.8 ± 3.0	) (9)
Ropivacaine	5.6 ± 0.6 (5)	10.2 ± 1.5 (5)	11.5 ± 1.4	¥ (11)
Tetracaine	1.5 ± 0.4 (6)	1.1 ± 0.2 (6)	1.1 ± 0.1	(11)
Thioridazine	2.8 ± 0.2 (10)	4.5 ± 0.5 (9)	6.9 ± 1.0	) (12)
Venlafaxine	461.4 ± 38.6 (6)	79.7 ± 6.5 (6)	36.5 ± 3.2	2 (8)

**Table 1.**  $IC_{50}$  values for sodium channel blockers on peak and late  $I_{Na}$ . Data represent mean ± SEM. Values in parenthesis represent the number of individual experiments performed

Figure 7. Correlations of compound IC<sub>50</sub> values for (A) peak and late  $I_{Na}$  in the presence of ATX-II, and on (B) peak  $I_{Na}$  in the absence or presence of ATX-II

## **CONCLUSIONS**

- Under control conditions no significant late I<sub>Na</sub> was detected in cells expressing the WT Nav1.5 channel when measured during the ramp. A small late  $I_{Na}$  (41.9 ± 5.4 pA) was observed in cells expressing the  $\Delta$ KPQ-Nav1.5 mutant channel that was insufficient for high throughput profiling of compounds
- Veratridine significantly enhanced the late  $I_{Na}$  (EC<sub>50</sub> = 27  $\mu$ M). However, it also inhibited the peak  $(IC_{50} = 84 \mu M)$  and altered the voltage-dependency of the  $I_{Na}$ , demonstrating that it is not an ideal enhancer for assessing either the peak or late  $I_{Na}$
- ATX-II produced a concentration-dependent increase in the late  $I_{Na}$  (392.4 ± 71.3 pA (ATX-II 100 nM)), but had minimal effect on the peak I<sub>Na</sub> and preserved the biophysical properties of I<sub>Na</sub>. ATX-II therefore had a profile that was suitable as an enhancer for measuring both peak and late  $I_{Na}$  simultaneously
- The presence of ATX-II did not significantly affect the pharmacology of the peak I<sub>Na</sub> for a series of sodium channel inhibitors when compared with results obtained in the absence of ATX-II
- We suggest that late  $I_{Na}$  should be measured at the maximum response generated during the ramp after a 200 ms depolarizing pulse at +40 mV (which allows for the completely inactivation of peak  $I_{Na}$ ) in the presence of ATX-II
- Our data demonstrate that in the presence of ATX-II (100 nM), peak and late  $I_{Na}$  can be measured simultaneously, allowing for a rapid and cost effective assessment of both currents

