

# Use- and state-dependent Na<sub>v</sub>1.5 blockers on QPatch X and *in vivo* and *in vitro* assays.

1. SOPHION BIOSCIENCE A/S  
Baltorpevej 154  
DK - 2750 Ballerup  
DENMARK

2. SOPHION BIOSCIENCE, Japan  
1716-6, Shinmachi  
Takasaki-shi, Gumma 370-1301  
JAPAN

3. DAIICHI SANKYO CO.LTD  
1-16-13 Kitakasai Edogawa-ku  
Tokyo 134-8630  
JAPAN

info@sophion.dk  
www.sophion.com



1 M. KNIRKE JENSEN

2 TOMOKO SAKAKURA

3 YASUYUKI ABE

4 HIDEO TAKAMORI

5 KIIYOSHI TAKASUNA

6 YUJI TSURUBUCHI

7 MORTEN SUNESEN

## Introduction

The cardiac voltage dependent sodium channel (Na<sub>v</sub>1.5) is responsible for the upstroke and directed propagation of action potentials in the heart, and is therefore a central ion channel in safety assessment and drug discovery. It is often important to determine the mode of action of a drug candidate, and this requires high-quality recordings and careful experiment planning. Such experiments usually do not lend themselves well to testing of large drug libraries. We show that the QPatch X in multi-hole mode can be successfully used in such a screening scenario. Furthermore, since Na<sub>v</sub>1.5 inhibition *per se* is not necessarily arrhythmogenic or cardiotoxic as is hERG inhibition, integrated *in vitro* and *in vivo* studies are necessary to assess a potential risk of cardiotoxicity including proarrhythmia. QPatch data combines relevantly with the data from Langendorff-perfused rabbit hearts and the ECG assay and effectively discriminate the cardiovascular profiles of compounds. For drug screening, we established a protocol for QPatch X that both tests the decay of the sodium current (30 Hz pulsetrain), and the recovery of the current from this pulse-train induced decay. The easy assay set-up in the QPatch Assay Software allows the combination of these elements into a single protocol, thereby shortening experimental time and costs.

## Materials & Methods

### QPatch experiments

**Cells:** CHO Na<sub>v</sub>1.5 QCells, grown according to Sophion SOP. The cells were supplied by B'SYS (Switzerland). **Ringer's solutions:** Extracellular (in mM): 2 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>, 10 HEPES, 4 KCl, 145 NaCl, 10 glucose, 10 TEA-Cl. pH 7.4, 310 mOsm. Intracellular (in mM): 135 CsF, 1 EGTA (solubilized in CsOH), to a total Cs concentration of 140 mM, 10 HEPES, 10 NaCl. pH 7.3. Adjusted to approx. 310 mOsm with sucrose. **Compounds:** Lidocaine, Flecaïnide (Sigma – dissolved in ethanol to a stock solution x1000 of the highest final concentration. Quinidine (Sigma) – dissolved directly in extracellular Ringer's solution. TTX (Alomone labs, Israel), dissolved in H<sub>2</sub>O to a stock solution x1000 of the highest final concentration. **Experimental setup:** All compounds were tested in six concentrations with increasing concentration of compound applied to the same measurement site. Experiments were done on QPatch X in multi-hole mode. QPatch operation in single-hole mode is a classic patch clamp experiment where one cell is in whole-cell configuration, whereas multi-hole mode comprises 10 cells in whole-cell configuration. The multi-hole mode therefore measures the summed current of ten cells. The voltage protocol (Figure 1) consisted of a 30 Hz pulsetrain (50 depolarizations to -20 mV) followed by a interpulse for recovery at V<sub>hold</sub>, and a final depolarization at -20 mV. V<sub>hold</sub> was either -80 mV or -115 mV. The interpulse duration for recovery lasted 200 ms, increasing by 60% per sweep to max. 3355 ms (7 sweeps) in the screening experiments. Another, similar, voltage protocol with finer increments in time (25%, starting at 50 ms), thus giving more datapoints in this segment, was used to determine time constants of recovery after the pulsetrain more precisely ("extended voltage protocol").

Figure 1.

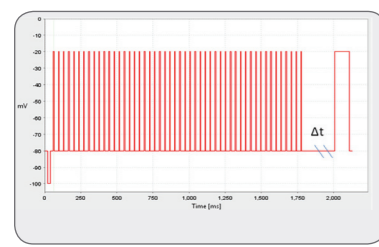


Figure 1. Voltage protocol for QPatch screening experiments. Here V<sub>hold</sub> of -80 mV is shown.

**Data analysis:** All data analysis for QPatch experiments was done using QPatch Assay Software 3.4 in combination with Origin 7.5. Time constants of recovery were calculated with a monoexponential fit. Data is represented as mean ± standard error.

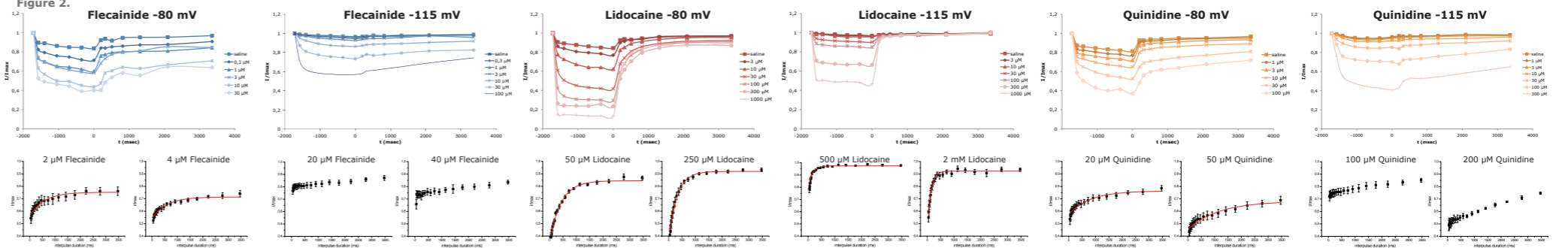
### Langendorff heart model

**Animals:** New Zealand White female rabbits (Oriental Yeast) at 8-13 weeks. Commercial diet (RC 4, Oriental Yeast) and chlorinated tap water was given ad libitum.

## Conclusion

The results obtained convincingly show that: 1) Flecaïnide is a use-dependent blocker of open-channels, as demonstrated by a slow decay of I<sub>Na</sub> (30 Hz pulsetrain), and a delayed recovery from the use-dependent block in the QPatch experiments. Flecaïnide produces ECG changes (marked PR and QRS prolongation, and slight QTc prolongation), and lethal arrhythmia (VT ~Vf) in both Langendorff-hearts and anesthetized rabbits at lower doses. 2) Lidocaine is a state-dependent blocker of inactivated channels, in that lidocaine produces a fast decay of I<sub>Na</sub> (30 Hz pulsetrain) and a fast recovery. Lidocaine produces slight ECG changes (PR and QRS prolongation), but no arrhythmia in Langendorff-hearts or anesthetized rabbits. 3) Quinidine demonstrates a slow decay of I<sub>Na</sub> (30 Hz pulsetrain), and delayed recovery from use-dependent block. Quinidine shows moderate PR and QRS prolongation, and severe morphological changes in ECGs. The present results suggest that: 1) The QPatch X can offer a time- and cost-effective unique screening scenario to select safe compounds with Na<sub>v</sub>1.5 blocking activity and no proarrhythmic activity, and 2) the pro-arrhythmic activity of the Na<sub>v</sub>1.5 blockers flecaïnide and quinidine might be attributable to their marked delay of recovery from use-dependent block.

## QPatch experiments



IC <sub>50</sub> (μM)	V <sub>hold</sub> -115 mV		V <sub>hold</sub> -80 mV	
	Resting block (1 <sup>st</sup> peak)	Use-dependent block (50 <sup>th</sup> peak)	Resting block (1 <sup>st</sup> peak)	Use-dependent block (50 <sup>th</sup> peak)
Flecaïnide	40.3 ± 10.4	22.6 ± 6.6	4.1 ± 0.5	2.2 ± 0.3
Lidocaine	1915.3 ± 434.9	586.1 ± 112.7	247.4 ± 61.4	44.1 ± 11.3
Quinidine	166.1 ± 18.6	97.0 ± 10.4	48.3 ± 7.7	18.7 ± 3.4
TTX	14.6 ± 1.4	7.5 ± 0.3	2.4 ± 0.3	0.81 ± 0.17

IC <sub>50</sub> (μM)	V <sub>hold</sub> -115 mV				V <sub>hold</sub> -80 mV			
	Conc. (μM)	tau (ms)	Conc. (μM)	tau (ms)	Conc. (μM)	tau (ms)	Conc. (μM)	tau (ms)
Flecaïnide	40	n/a	20	n/a	4	575	2	560
Lidocaine	2000	128	500	147	250	317	50	425
Quinidine	200	n/a	100	n/a	50	1151	20	795

Table 1. IC<sub>50</sub>s for all compounds tested were determined at two different points in the pulsetrain – part of the voltage protocol, namely the first and last (50<sup>th</sup>) peaks, corresponding to resting-state block and use-dependent block, respectively. All tests were conducted at a holding potential of either -80 mV, to be close to the true physiological conditions of the cell; or -115 mV to test at a potential where no ion channels are inactivated (V<sub>0</sub> for Na<sub>v</sub>1.5 is approximately -60 mV under the same experimental conditions as is used here – data not shown).

Table 2. Time constants of recovery from use-dependent block. Recovery was determined after the 30 Hz pulsetrain-induced decay. The extended voltage protocol with small time increments per sweep was used for this determination. The rank order of time constants obtained with the extended voltage protocol correlates well with the rank order of recovery from use-dependent block obtained with the screening protocol (see the upper graphs in Figure 2 (fits/time constants not shown) and compare with the small graphs, where the time constants are calculated from), where lidocaine has the quickest recovery, flecaïnide intermediate and quinidine the slowest.

## Langendorff heart model

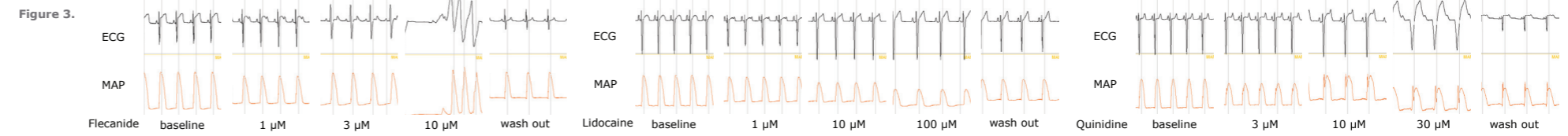


Table 3.	Conc. (μM)	HR (%)	PR (%)	QRS (%)	MAP <sub>90c</sub> (%)
Flecaïnide	1	-7.5	15.5	9.0	-2.1
	3	-15.8	54.0	80.2	-9.6
	10	Arrhythmia appeared			
Lidocaine	1	-4.0	4.8	0.7	-7.6
	10	-9.4	5.8	3.3	-12.2
	100	-24.4	29.7	21.2	-6.4
Quinidine	3	-11.8	14.5	4.0	10.8
	10	-29.8	31.9	20.2	9.5
	30	Severe morphology change of ECG			

Table 3. Summary of rabbit Langendorff-perfused heart preparations (n=3).

### Figure 3.

Flecaïnide at 1 μM showed only 15% prolongation of PR. At 3 μM, flecaïnide reduced the heart rate 16% and showed marked prolongation of PR (54%) and QRS (80%) but had almost no effect on MAP<sub>90c</sub> (reduction of 10%) (see also Table 3). At 10 μM, flecaïnide induced ventricular tachycardia, ventricular fibrillation or ventricular pause in all the preparations. Lidocaine showed no effect at concentrations of 1 and 10 μM. At 100 μM, lidocaine reduced the heart rate by 25% and showed a 30% prolongation of PR and a 21% prolongation of QRS. No effect on MAP<sub>90c</sub> was observed (see also Table 3). Quinidine at 3 μM showed 12% reduction of the heart rate, 15% prolongation of PR and 11% prolongation of QRS. At 10 μM, the changes in heart rate and PR were enhanced to 30% and 32%, respectively, and QRS was also prolonged 20%. MAP<sub>90c</sub> was not more enhanced at 10 μM than at 3 μM (about 10%). At 30 μM, ECG could not be analyzed because quinidine induced severe morphological changes of the ECG.

### Conclusions on Langendorff model

All the representative three Na<sub>v</sub>1.5 blockers, flecaïnide, lidocaine and quinidine, induced decrease in heart rate and prolongation of PR and QRS with minor changes on MAP<sub>90c</sub>. However, there were apparent differences of the degree of PR or QRS prolongation, and arrhythmogenic activity, namely flecaïnide showed the most potent inhibition on cardiac conduction velocity and arrhythmogenic activity, while lidocaine was the most weak and not arrhythmogenic.

## In vivo ECG anesthetized rabbit model

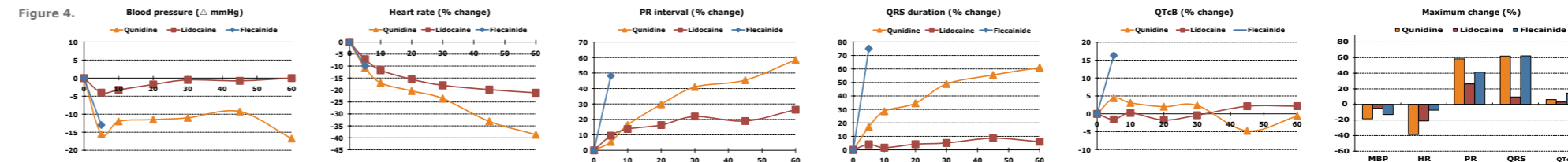


Table 4.	MBP	HR	PR	QRS	QTc	Arrhythmia incidence (%)	Onset of arrhythmia (min)
Flecaïnide	-13.0	-7.4	41.4	62.3	14.5	100%	6
Lidocaine	-5.0	-21.2	26.3	9.6	3.1	0%	>60
Quinidine	-18.5	-38.7	58.4	62.1	6.2	0%	>60

Table 4. Mean of individual maximum change (% change; BP: delta mmHg) in anesthetized rabbits (n=3-4).

### Figure 4.

Effect of Na<sub>v</sub>1.5 blockers on cardiovascular systems in ketamine and xylazine anesthetized rabbits (n=3-4). Flecaïnide at 1 mg/kg/min for 60 min i.v. infusion promptly decreased the blood pressure and heart rate, and prolonged the PR and QRS intervals and QTc with individual maximum changes of about -15 mmHg, -10%, 40%, 60%, and 15%. All the animals showed arrhythmia (VT followed by Vf), followed by death within 20 min after the start of administration (N=3). Lidocaine at 1 mg/kg/min for 60 min i.v. infusion gradually decreased the heart rate and prolonged the PR interval with individual maximum changes of about -20% and 25%, while it had little or no effect on blood pressure, QRS or QTc. No arrhythmia was observed (N=4). Quinidine at 1 mg/kg/min for 60 min i.v. infusion gradually decreased the heart rate and prolonged PR and QRS intervals with individual maximum changes of about -40%, 60% and 60%, while it introduced a slight decrease in blood pressure (about -20 mmHg at 60 min) with little or no effect on QTc. No arrhythmia was observed but severe morphological changes of ECG were observed in 2 out of 4 animals.

### Conclusions on in vivo rabbit model

The IC<sub>50</sub>s of Na<sub>v</sub>1.5 block of flecaïnide, quinidine and lidocaine are here reported to be 2 μM, 18 μM, and 44 μM (at the physiologically relevant V<sub>0</sub> of -80 mV, and determined at the last peak in the pulsetrain). In the present study, flecaïnide and quinidine showed stronger changes of ECG and arrhythmogenic activity than lidocaine. These differences are thus expected given that the same test concentration was used in this *in vivo* assay. The other Na<sub>v</sub>1.5 blocking profile such as an effect on recovery from use-dependent block might be related to their *in vivo* effects.