

## Application Report

# hCa<sub>v</sub>1.2 recordings on Qube 384 using step, train and CiPA protocols

A reproducible high-throughput assay with high success rates for biophysical and pharmacological studies of the Ca<sub>v</sub>1.2 channel

### Summary

The development of screening assays for the Ca<sub>v</sub>1.2 channel has been challenging due to the tendency of Ca<sub>v</sub>1.2 expressing cell lines to exhibit declining current levels (rundown) during the experiment<sup>1,2</sup>. Here we report a robust Ca<sub>v</sub>1.2 assay yielding high success rates, low rundown and reliable pharmacology.

- Pharmacology and current-voltage relationship in accordance with literature values
- Success rates up to **94%**
- Stable currents with rundown as low as **(1.2 ± 0.9)%** per minute
- The CiPA protocol yields stable currents with rundown as low as **(1 ± 1)%** per minute

### Introduction

The L-type voltage-gated calcium channel Ca<sub>v</sub>1.2 is expressed in various mammalian tissues, including heart and smooth muscle<sup>3-5</sup>, and its dysfunction has been implicated in a range of cardiovascular and neurological diseases<sup>6,7</sup>. Opening of the Ca<sub>v</sub>1.2 channel results in an increase in the intracellular concentration of calcium ions (Ca<sup>2+</sup>), affecting a variety of cellular processes including muscle contraction, hormone secretion and neuronal transmission<sup>7-9</sup>, thus rendering the channel an important target in e.g. safety pharmacology screening.

Ca<sub>v</sub>1.2 channels require a strong depolarization to activate, display relatively long-lasting activity and can be blocked by low micromolar concentrations of e.g. dihydropyridines, phenylalkylamines and benzothiazepines<sup>8,9</sup>. Following activation, the channel displays both Ca<sup>2+</sup>- and voltage-dependent inactivation<sup>10</sup>. In this study, current traces from HEK-hCa<sub>v</sub>1.2 cells were recorded on the high-throughput platform Qube 384 using both single-hole and multi-hole (10 patch holes) QChips. The current-voltage relationship and pharmacology of three different compounds were determined with high success rates and low rundown.

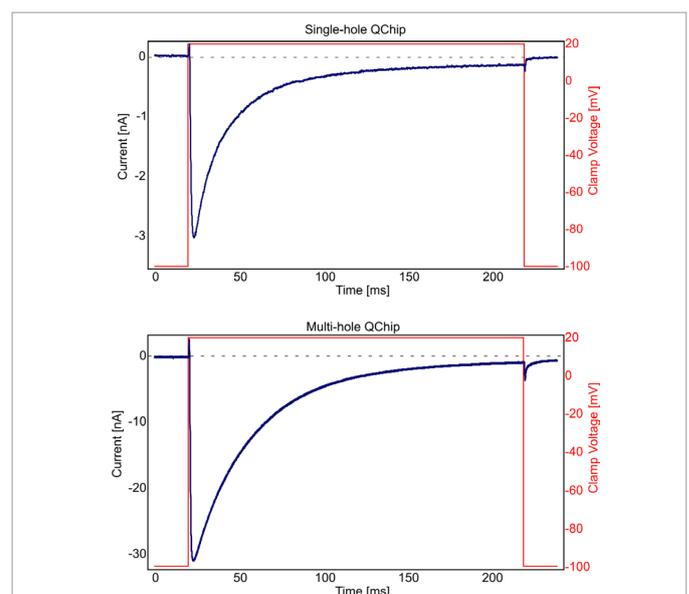
### Results and discussion

The following experiments were carried out:

1. Recordings of Ca<sup>2+</sup> currents
2. Current-voltage relationship of Ca<sub>v</sub>1.2 channels
3. Pharmacology of Ca<sub>v</sub>1.2 channels
4. Rundown analysis using multi-hole QChips
5. Recordings of Ca<sup>2+</sup> currents and rundown analysis using a step-ramp (CiPA) protocol

#### 1. Recordings of Ca<sup>2+</sup> currents

Cells were clamped to -100 mV and Ca<sup>2+</sup> currents were evoked by application of a depolarization step to +20 mV for 200 ms (Fig. 1).



**Fig. 1:** Representative recordings of Ca<sup>2+</sup> current (blue) on a single-hole (left) and multi-hole (right) QChip following a 200 ms depolarization from -100 mV to 20 mV (red) in Ca<sub>v</sub>1.2 expressing cells.

The criteria for a successful recording were (average of first 10 sweeps):

**Single-hole QChip:**

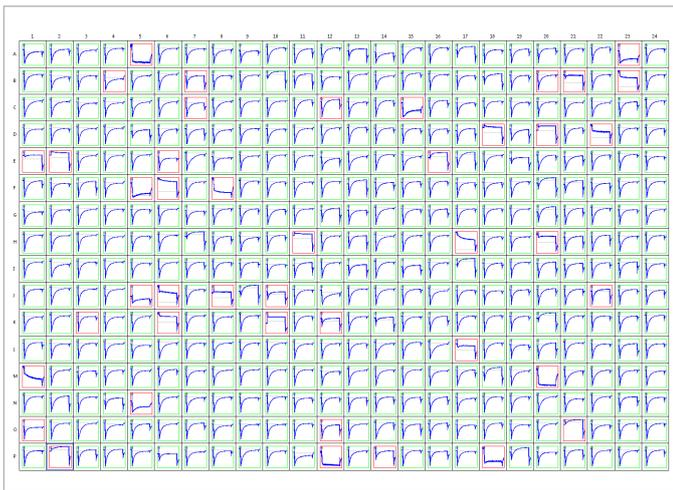
- $I_{\text{peak}} < -100 \text{ pA}$
- $R_{\text{membrane}} > 100 \text{ M}\Omega \text{ per cell}$
- $C_{\text{total}} > 4 \text{ pF per cell}$

**Multi-hole QChip:**

- $I_{\text{peak}} < -500 \text{ pA}$
- $R_{\text{membrane}} > 100 \text{ M}\Omega \text{ per cell}$
- $C_{\text{total}} > 4 \text{ pF per cell}$

Where  $I_{\text{peak}}$  is the peak current,  $R_{\text{membrane}}$  is the membrane resistance and  $C_{\text{total}}$  is the total capacitance.

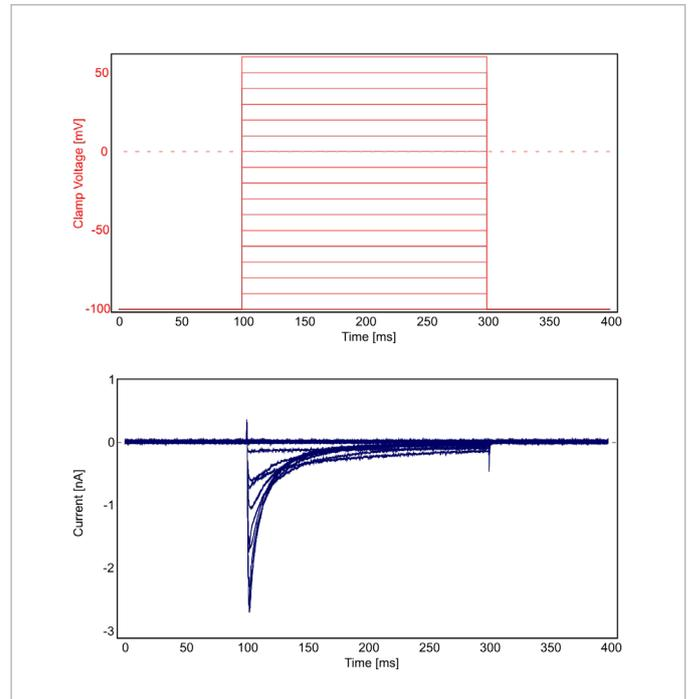
A plate-view of  $\text{Ca}_v1.2 \text{ Ca}^{2+}$  currents recorded on a multi-hole QChip is shown in Figure 2. The success rates were up to 89% and 94%, for single-hole and multi-hole QChips, respectively. The mean success rates and peak currents were  $(85 \pm 3)\%$  and  $(-2.9 \pm 0.4) \text{ nA}$  for single-hole QChips ( $N = 3$ ), and  $(82 \pm 8)\%$  and  $(-14 \pm 2) \text{ nA}$  for multi-hole QChips ( $N = 8$ ). The success rates and peak currents started declining at cell passage numbers above 20.



**Fig. 2:** Plate overview of  $\text{Ca}_v1.2 \text{ Ca}^{2+}$  currents recorded on a multi-hole QChip.

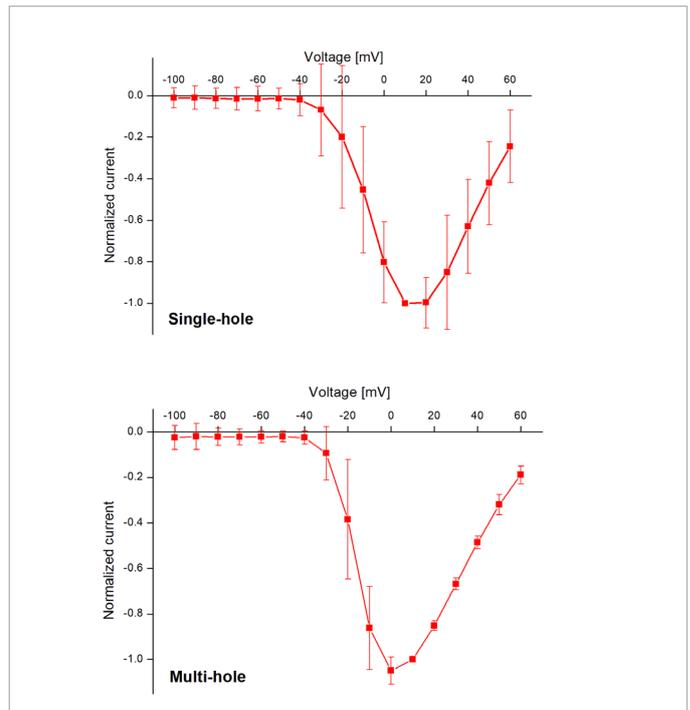
**2. Current-voltage relationship of  $\text{Ca}_v1.2$  channels**

The current-voltage relationship was quantified on single-hole and multi-hole QChips by applying a depolarization step protocol from  $-100 \text{ mV}$  to  $+60 \text{ mV}$  for  $200 \text{ ms}$  in increments of  $10 \text{ mV}$ . The time between sweeps was  $15 \text{ s}$  (Fig. 3).



**Fig. 3:** Representative measurement of current-voltage relationship using a single-hole QChip. Top: Depolarization step protocol in  $10 \text{ mV}$  increments from  $-100 \text{ mV}$  to  $+60 \text{ mV}$ . Bottom:  $\text{Ca}^{2+}$  current response of a cell expressing  $\text{Ca}_v1.2$  channels.

The peak current was extracted at each voltage step, normalized to the peak current at  $10 \text{ mV}$  and plotted as a function of voltage (Fig. 4).

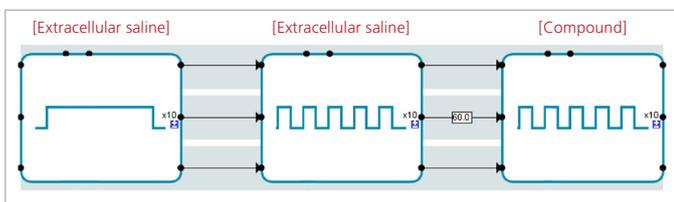


**Fig. 4:** Current-voltage plots of normalized current vs. voltage (mV) for a single-hole QChip (top) and a multi-hole QChip (bottom). The error bars are the SD.

The current-voltage relationship is in agreement with previous findings<sup>2</sup>. Note that the current-voltage plot recorded on multi-hole QChips is left-shifted due to a lack of  $R_s$  compensation.

### 3. Pharmacology of $Ca_v1.2$ channels

The  $Ca_v1.2$  pharmacology was quantified on single-hole and multi-hole QChips. Cells were clamped to -100 mV and  $Ca^{2+}$  currents were evoked by application of a depolarization step to +20 mV for 200 ms (see Fig. 1). The experiment setup in the Sophion Viewpoint software is shown in Figure 5. After 10 initial depolarization steps (15 s between sweeps), 2 blocks of 10 depolarization trains (30 s between sweeps) were executed. Each train contained 5 pulses at 2.5 Hz (Fig. 5).



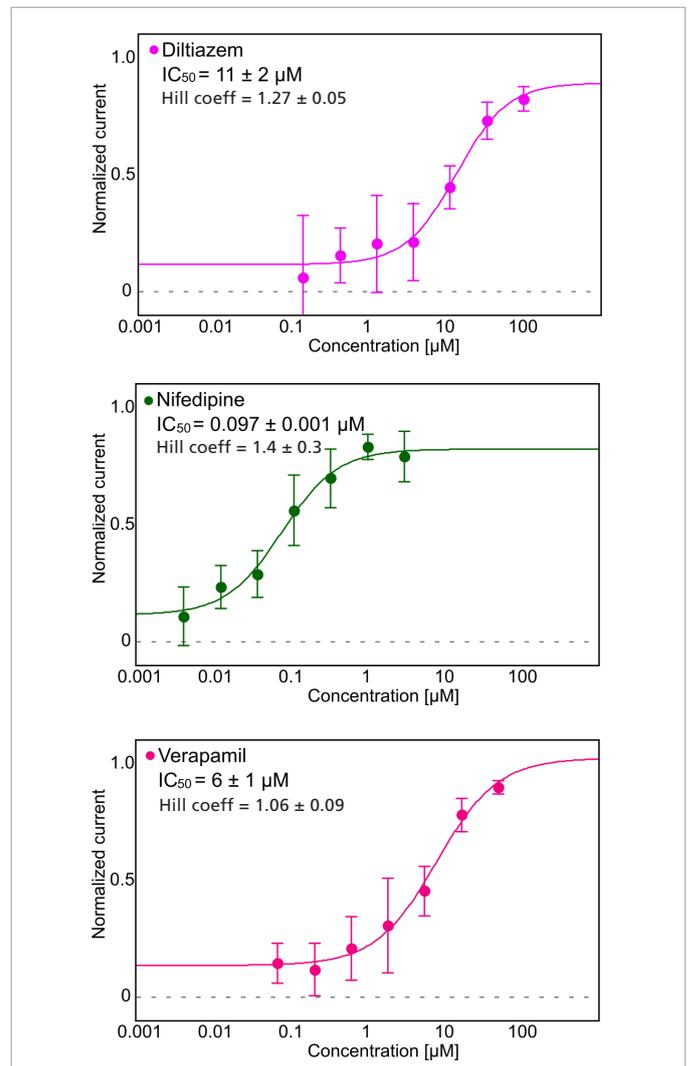
**Fig. 5:** Experiment setup in the Sophion Viewpoint software for  $Ca_v1.2$  pharmacology experiments. Voltage protocols: block 1: 10 single depolarization steps (left), block 2 and 3: 10 train depolarizations (middle and right). The compound was applied in block 3 (right).

The saline baseline was established in the second block of the experiment and in the third block of the experiment the three compounds, diltiazem, nifedipine and verapamil, were applied in 7 concentrations (non-cumulative). The concentration range was a 3-fold dilution with the highest concentration being:

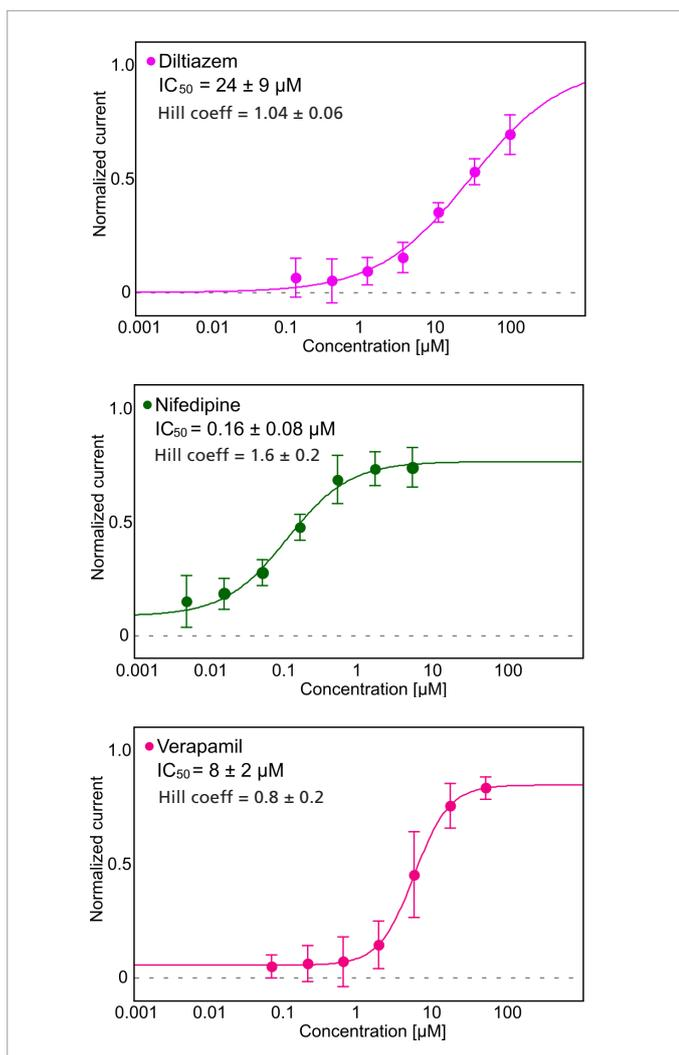
- Diltiazem 100  $\mu\text{M}$
- Nifedipine 3  $\mu\text{M}$
- Verapamil 50  $\mu\text{M}$

The peak current of the last train depolarization step in block 3 was normalized to the corresponding peak current in block 2 according to:  $1 - \left( \frac{I_{\text{peak}}(\text{compound})}{I_{\text{peak}}(\text{saline})} \right)$  and plotted as a function of concentration for the three compounds (Fig. 6 and Fig. 7).

The  $IC_{50}$  values were extracted by fitting the Hill equation to the data and listed in Table 1 together with literature values.



**Fig. 6:** Dose-response relationships measured using single-hole QChips for diltiazem (left), nifedipine (middle) and verapamil (right). The  $IC_{50}$  values and Hill coefficients were estimated by fitting the Hill equation to the data and the reported values are the average  $\pm$  SD of three QChips with a total of 48 experiments per compound.



**Fig. 7:** Dose-response relationships measured using multi-hole QChips for diltiazem (left), nifedipine (middle) and verapamil (right). The  $IC_{50}$  values and Hill coefficients were estimated by fitting the Hill equation to the data and the reported values are the average  $\pm$  SD of three QChips with a total of 48 experiments per compound.

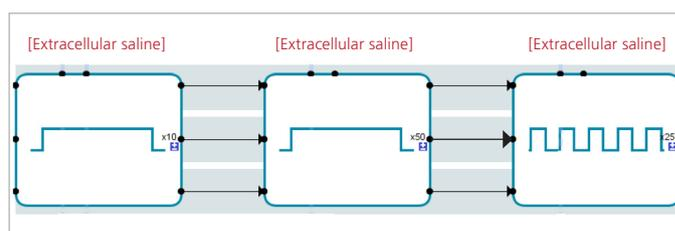
**Table 1:** Potency of the three tested compounds.  $IC_{50}$  values (average  $\pm$  SD) for the listed compounds, quantified on single-hole and multi-hole QChips, together with literature values.

Compound	$IC_{50}$ ( $\mu$ M) Single-hole	$N_{QChip}$	$IC_{50}$ ( $\mu$ M) Multi-hole	$N_{QChip}$	$IC_{50}$ ( $\mu$ M) Literature values <sup>11</sup>
Diltiazem	11 $\pm$ 2	3	24 $\pm$ 9	3	5.6 $\pm$ 0.2
Nifedipine	0.097 $\pm$ 0.01	3	0.16 $\pm$ 0.08	3	0.056 $\pm$ 0.002
Verapamil	6 $\pm$ 1	3	8 $\pm$ 2	3	2.69 $\pm$ 0.09

The  $IC_{50}$  values were in good agreement with  $IC_{50}$  values measured on other  $Ca_v1.2$  cell lines in-house (11  $\mu$ M  $\pm$  3  $\mu$ M, 0.052  $\mu$ M  $\pm$  n/a  $\mu$ M, 5  $\mu$ M  $\pm$  2  $\mu$ M for diltiazem, nifedipine and verapamil, respectively).

#### 4. Rundown analysis using multi-hole QChips

A concern in  $Ca_v1.2$  assays is the rundown of the current under continuous stimulation. Here, the rundown was quantified using multi-hole QChips. Cells were clamped to -100 mV and  $Ca^{2+}$  currents were evoked by application of a depolarization step to +20 mV for 200 ms (see Fig.1). The experiment setup in the Sophion Viewpoint software is shown in Figure 8. After 10 initial depolarization steps (block 1, 15 s between sweeps), an additional 50 sweeps of single depolarization steps were applied (block 2, 15 s between sweeps), followed by 25 depolarization trains, each containing 5 pulses at 2.5 Hz (block 3, 30 s between sweeps).



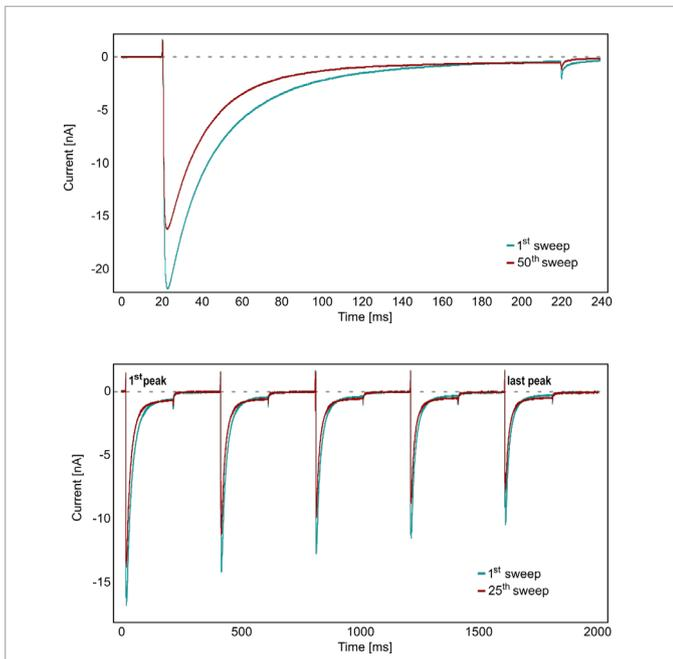
**Fig. 8:** Experiment setup for a  $Ca_v1.2$  rundown experiment in the Sophion Viewpoint software. Voltage protocols: block 1: 10 single depolarization steps (left), block 2: 50 depolarization steps (middle), and block 3: 25 train depolarizations.

The rundown was quantified for the second experiment block (see Fig. 8) by comparing the peak currents evoked by the 1<sup>st</sup> and 50<sup>th</sup> depolarization (Fig. 8):

$$\text{Rundown (\% per min): } \frac{I_{\text{peak}_1} - I_{\text{peak}_{50}}}{I_{\text{peak}_1} \cdot \Delta t} \cdot 100\%$$

where  $\Delta t$  is the time between sweep 1 and 50 in minutes. The mean rundown was (1.2  $\pm$  0.9)% per min.

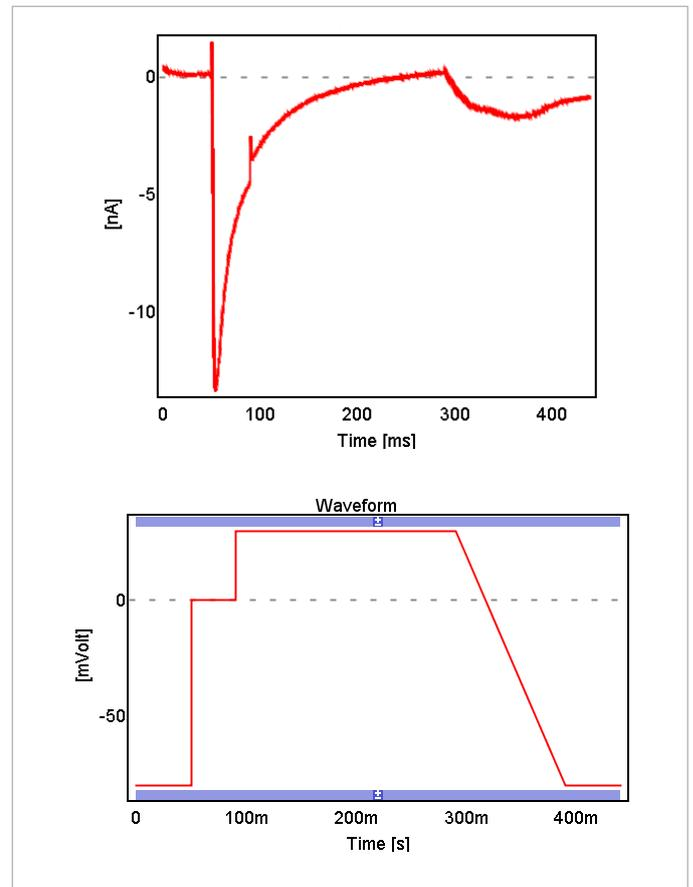
The rundown was also quantified for the third experiment block (see Fig. 8), by comparing the 1<sup>st</sup> and 25<sup>th</sup> train depolarization (see Fig. 9). For the first depolarization within the train the mean rundown was (2.0  $\pm$  0.5)% per min. For the last depolarization within the train the mean rundown was (2.5  $\pm$  0.5)% per min.



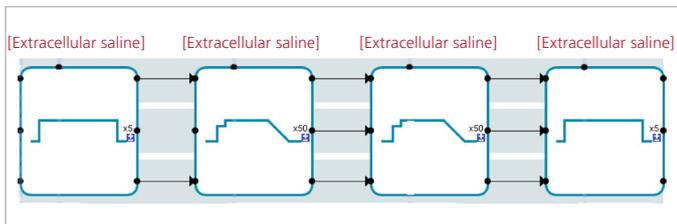
**Fig. 9:** Example current traces used for quantifying rundown. Top 1<sup>st</sup> (green) and 50<sup>th</sup> (red) depolarization in experiment block 2. Bottom: 1<sup>st</sup> (green) and 25<sup>th</sup> (red) depolarization train in experiment block 3. We quantified rundown for the first and last peak of the train.

## 5. Recordings of Ca<sup>2+</sup> currents and rundown analysis using a step-ramp (CiPA) protocol

Here we report the response of Ca<sub>v</sub>1.2 cells to a step-ramp voltage protocol, which is similar to the voltage protocol proposed in the Comprehensive *in vitro* Proarrhythmia Assay (CiPA) initiative. The objective of the CiPA initiative is to initiate a new paradigm for safety assessment, taking several ion channels – including Ca<sub>v</sub>1.2 – into account. For the Ca<sub>v</sub>1.2 CiPA voltage protocol, cells were clamped to -80 mV and Ca<sup>2+</sup> currents were evoked by application of a depolarization step to 0 mV for 40 ms, followed by an additional depolarization step to +20 mV for 200 ms and a 100 ms ramp to -80 mV (Fig. 10). In this study, a total of 100 sweeps (15 or 30 s between sweeps) of the CiPA protocol stimulations (block 2 and 3) were flanked by five 200 ms standard depolarization pulses from -100 mV to 20 mV (block 1 and 4), as indicated in the Sophion Viewpoint software (Fig. 11).



**Fig. 10:** Original recording on a multi-hole QChip of a Ca<sub>v</sub>1.2 Ca<sup>2+</sup> current (top) following a depolarization according to the CiPA protocol (bottom).



**Fig. 11:** CiPA experiment setup for a  $Ca_v1.2$  rundown measurements in the Sophion Viewpoint software. Voltage protocols: block 1 and 4: 5 single depolarization steps (15 s between sweeps), block 2: 50 CiPA depolarization steps (30 s between sweeps), block 3: 50 CiPA depolarization steps (15 s between sweeps).

According to the previously defined success criteria, the CiPA protocol execution had a 67% success rate and rundown of  $(1.0 \pm 0.7)\%$  per min (30 s in between sweeps) and  $(1 \pm 1)\%$  per min (15 s in between sweeps).

## Conclusion

$Ca_v1.2$   $Ca^{2+}$  currents were recorded on the high-throughput platform Qube 384 on both single-hole and multi-hole QChips. The reported current-voltage relationship and pharmacology experiments were in accordance with literature values and highly reproducible. The current levels were stable over time with an average rundown of  $(1.2 \pm 0.9)\%$  per min and high success rates were achieved (up to **89%** and **94%** for single-hole and multi-hole QChips, respectively) with cells at passage numbers below 20. In addition, the CiPA protocol exhibited stable currents albeit with slightly lower success rates. Conclusively, Qube is an ideal platform for running  $Ca_v1.2$  experiments for screening as well as characterization.

## Material and methods

Experiments in this study were performed on HEK-h $Ca_v1.2$  cells, kindly provided by SB Drug Discovery. The cell line is stably transfected with the  $Ca_v1.2$  channel and cultured in 30  $\mu$ M verapamil according to SB Drug Discovery guidelines. HEK-h $Ca_v1.2$  cells were harvested according to Sophion standard procedures.



Sophion Bioscience A/S, Baltorpevej 154, 2750 Ballerup, Denmark  
Phone: +45 4460 8800, E-mail: info@sophion.com

sophion.com

6 Sophion

## Author:

Melanie Schupp, application scientist, Sophion Bioscience A/S  
Kadla R Rosholm, application scientist, Sophion Bioscience A/S

## References:

- Balazubramanian, B et al. Optimization of  $Ca_v1.2$  screening with an automated planar patch clamp platform, *J Pharmacol Tox Met* 2009;**59**:62-72
- Kuryshv YA et al. Evaluating State Dependence and Subtype Selectivity of Calcium Channel Modulators in Automated Electrophysiology Assays. *Assay Drug Dev Technol* 2014;**12**(2): 110–119
- Sinnegger-Brauns MJ et al. Isoform-specific regulation of mood behavior and pancreatic  $\beta$  cell and cardiovascular function by L-type  $Ca^{2+}$  channels. *J Clin Invest* 2004;**113**:1430-9
- Moosmang S et al. Dominant role of smooth muscle L-type calcium channel  $Ca_v1.2$  for blood pressure regulation. *EMBO J* 2003;**22**:6027-34
- Zhang J et al. Role of  $Ca_v1.2$  L-type  $Ca^{2+}$  channels in vascular tone: effects of nifedipine and  $Mg^{2+}$ . *Am J Physiol Heart Circ Physiol* 2007;**292**:H415-25
- Zamponi, GW Targeting voltage-gated calcium channels in neurological and psychiatric diseases. *Nat Rev Drug Discovery* 2007;**15**:19-34
- Liao P & Soong TW  $Ca_v1.2$  channelopathies: from arrhythmias to autism, bipolar disorder, and immunodeficiency. *Eur J Physiol* 2016;**460**:353-359
- Hofman F et al. L-type  $Ca_v1.2$  calcium channels: from in vitro findings to in vivo function. *Physiol Rev* 2014;**94**:303-326
- Reuter H Calcium channel modulation by neurotransmitters, enzymes and drugs. *Nature* 1983;**301**:569–574
- Hennessey JA et al. A CACNA1C variant associated with reduced voltage-dependent inactivation, increased  $Ca_v1.2$  channel window current, and arrhythmogenesis. *PLOS One* 2014;**9**,e106982
- Elkins CR et al. Variability in high-throughput ion-channel screening data and consequences for cardiac safety assessment *J Pharmacol Tox Met* 2013;**68**:112-122