

## Application Report

# hCa<sub>v</sub>1.2 recordings using QPatch

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 the channel 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 of up to **98%**
- Stable currents with rundown as low as **2% per min**

### Introduction

The L-type voltage-gated calcium channel Ca<sub>v</sub>1.2 is expressed in various mammalian tissues, including heart, smooth muscle and brain<sup>3,4,5</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>7,8,9</sup>, thus rendering the channel an important pharmacological target.

Ca<sub>v</sub>1.2 channels require a strong depolarization to be activated, they 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 using QPatch employing both single-hole and multi-hole (10 patch holes per well) QPlates. The current-voltage relationship and pharmacology of three different compounds were determined with high success rates and low rundown.

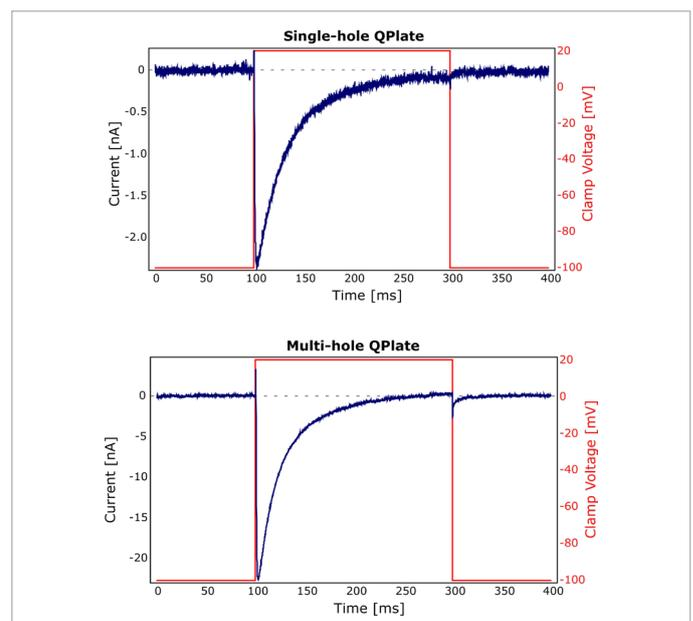
### Results and discussion

The following experiments were carried out:

- Recordings of Ca<sup>2+</sup> currents
- Current-voltage relationship of Ca<sub>v</sub>1.2 channels
- Pharmacology of Ca<sub>v</sub>1.2 channels
- Rundown analysis

#### 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 (top) and multi-hole (bottom) QPatch 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 QPlate:**

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

**Multi-hole QPlate:**

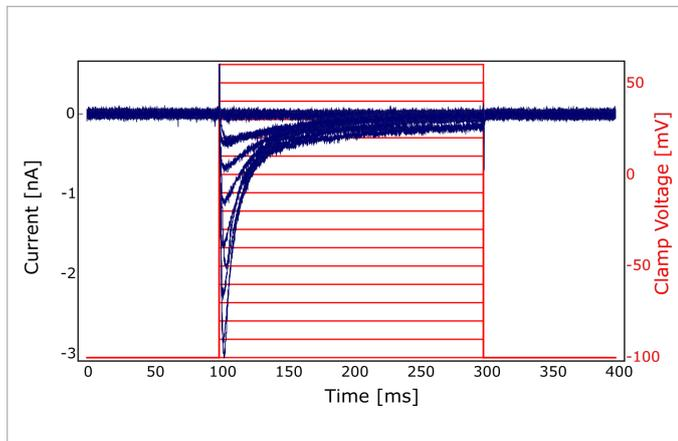
$I_{\text{peak}} < -500 \text{ pA}$   
 $R_{\text{membrane}} > 100 \text{ M}\Omega$  per cell  
 $C_{\text{total}} > 4 \text{ pF}$  per cell

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

The success rates were up to **71%** and **98%**, for single-hole and multi-hole QPlates, respectively. The mean success rates and peak currents were  $(58 \pm 8) \%$  and  $(1.2 \pm 0.3) \text{ nA}$  for single-hole QPlates ( $N = 6$  QPlates), and  $(88 \pm 5) \%$  and  $(14 \pm 3) \text{ nA}$  for multi-hole QPlates ( $N = 6$  QPlates). The success rates and peak currents started declining at cell passage numbers above 20.

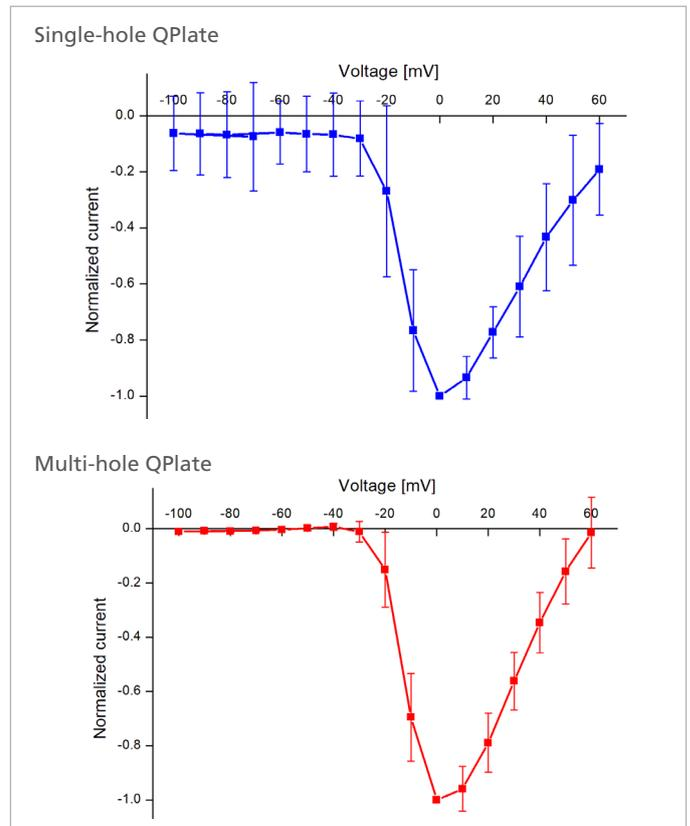
**2. Current-voltage relationship of  $\text{Ca}_v1.2$  channels**

The current-voltage relationship was quantified on single-hole and multi-hole QPlates 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. 2).



**Fig. 2:** Representative measurement of current-voltage relationship using a single-hole QPlate.  $\text{Ca}^{2+}$  current response of a cell expressing  $\text{Ca}_v1.2$  channels (blue) following a depolarization step protocol in  $10 \text{ mV}$  increments from  $-100 \text{ mV}$  to  $+60 \text{ mV}$  (red).

The peak current was extracted at each voltage step, normalized to the peak current at  $0 \text{ mV}$  and plotted as a function of voltage (Fig. 3).



**Fig. 3:** Current-voltage plots of normalized current versus voltage [mV] for a single-hole QPlate (top) and a multi-hole QPlate (bottom). Displayed is the averaged normalized current  $\pm$  SD.

**3. Pharmacology of  $\text{Ca}_v1.2$  channels**

The  $\text{Ca}_v1.2$  pharmacology was quantified on single-hole and multi-hole QPlates. Cells were clamped to  $-100 \text{ mV}$  and  $\text{Ca}^{2+}$  currents were evoked by application of a depolarization step to  $+20 \text{ mV}$  for  $200 \text{ ms}$  (see Fig. 1).

After 10 initial depolarization steps ( $15 \text{ s}$  between sweeps), a voltage protocol of 10 depolarization trains (5 pulses at  $2.5 \text{ Hz}$  and  $30 \text{ s}$  between sweeps) was applied three times. First, the baseline was established ( $0.1\% \text{ DMSO}$ ) (Fig. 4A). Subsequently, one of the three compounds, diltiazem, nifedipine or verapamil, was applied in 5 concentrations (non-cumulative). The concentration range was a 4-fold dilution with the highest concentration being:

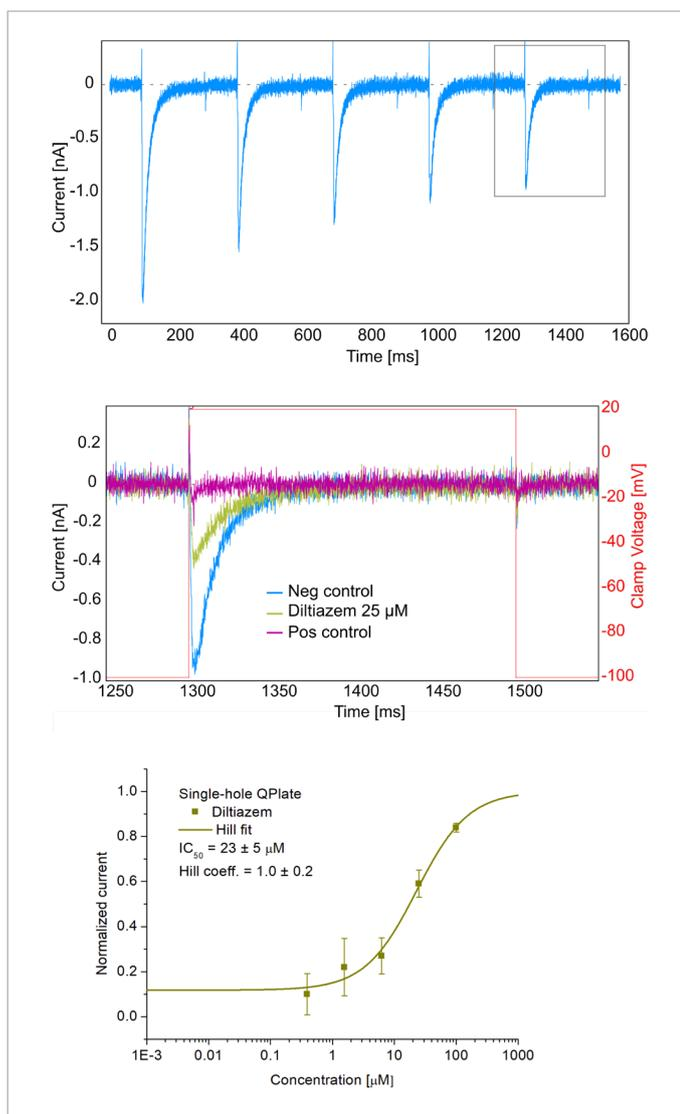
- Diltiazem  $100 \mu\text{M}$
- Nifedipine  $3 \mu\text{M}$
- Verapamil  $50 \mu\text{M}$

Finally,  $10 \mu\text{M}$  nifedipine was applied as a positive control.

The peak current of the last pulse of the depolarization train (ROI in Fig. 4 top) was extracted for the compound ( $I_{comp}$ ) and normalized to the corresponding peak current of the negative ( $I_-$ ) and positive ( $I_+$ ) controls (Fig. 4 middle), according to:

$$\text{Normalized current} = \frac{I_{comp} - I_-}{I_+ - I_-}$$

The normalized current was plotted as a function of compound concentration (Fig. 4 bottom). The  $IC_{50}$  values were extracted by fitting the Hill equation to the data and listed in Table 1 together with literature values.



**Fig. 4:** Figure 4:  $Ca_v1.2$  pharmacology measurement using a **single-hole QPlate**. Top: Representative  $Ca_v1.2$  current trace upon the application of a depolarization train (5 pulses from  $-100$  mV to  $+20$  mV at 2.5 Hz). Middle: Peak current of the last depolarization of the train (ROI in top graph) displayed for the negative control (0.1 % DMSO, blue), upon the addition of  $25 \mu M$  diltiazem (green), and for the positive control ( $10 \mu M$  nifedipine) (pink). Bottom: Dose-response relationship of diltiazem. Each data point is the average  $\pm$  SD of at least 5 cells and the  $IC_{50}$  value and Hill coefficient were estimated by fitting the Hill equation.

**Table 1:** Potency of the three tested compounds.  $IC_{50}$  values for the listed compounds, quantified on single-hole and multi-hole QPlates, together with literature values.

| Compound   | $IC_{50}$ ( $\mu M$ )<br>Single-hole | $N_{QPlate}$ | $IC_{50}$ ( $\mu M$ )<br>Multi-hole | $N_{QPlate}$ | $IC_{50}$ ( $\mu M$ )<br>Literature values |
|------------|--------------------------------------|--------------|-------------------------------------|--------------|--|
| Diltiazem  | $23 \pm 5$                           | 4            | $21 \pm 5$                          | 4            | 6 - $250^{11,12}$                          |
| Nifedipine | $0.32 \pm 0.03$                      | 4            | $0.11 \pm 0.08$                     | 4            | 0.06 - $0.10^{1,11}$                       |
| Verapamil  | $12 \pm 2$                           | 4            | $5.7 \pm 0.6$                       | 4            | 2.7 - $5.3^{1,11}$                         |

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

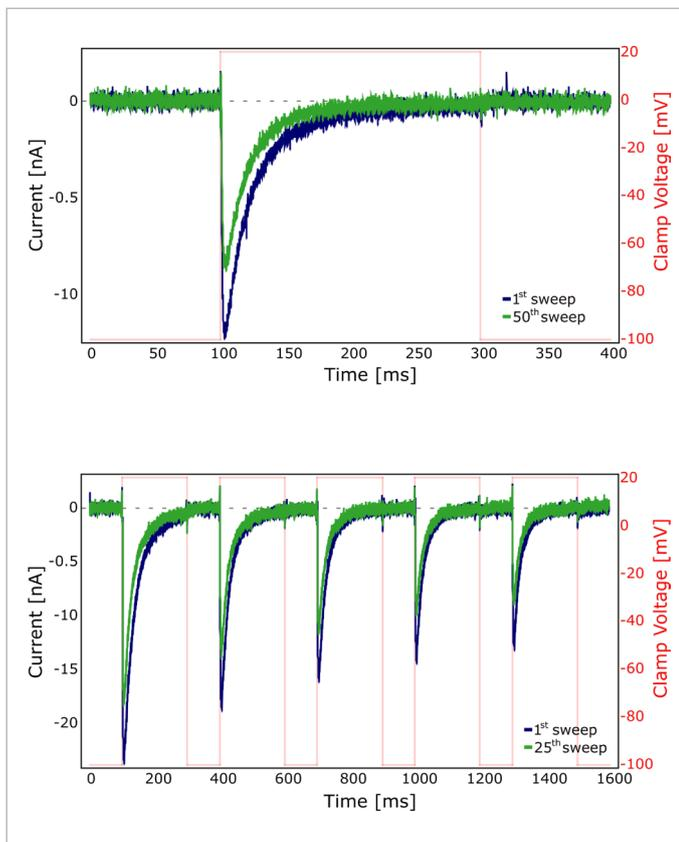
A concern with  $Ca_v1.2$  assays is the rundown of the current under continuous stimulation. Here, the rundown was quantified using single-hole and multi-hole QPlates. Cells were clamped to  $-100$  mV and  $Ca^{2+}$  currents were evoked by application of a depolarization step to  $+20$  mV for 200 ms (see Fig. 1). After 10 initial depolarization steps (15 s between sweeps), an additional 50 sweeps of single depolarization steps were applied (15 s between sweeps), followed by 25 depolarization trains that contained 5 pulses at 2.5 Hz (30 s between sweeps).

The rundown was quantified for the second voltage protocol by comparing the peak currents evoked by the 1<sup>st</sup> and 50<sup>th</sup> depolarization (Fig. 5, top):

$$\text{Rundown (\% per min): } \frac{I_{peak_1} - I_{peak_{50}}}{I_{peak_1} \cdot \Delta t}$$

where  $\Delta t$  is the time between sweep 1 and 50 in minutes. The mean rundown was  $(3.0 \pm 0.1)$  % per min and  $(1.7 \pm 0.2)$  % per min for multi-hole and single-hole QPlates, respectively.

The rundown was also quantified for the third voltage protocol, by comparing the 1<sup>st</sup> and 25<sup>th</sup> train depolarization (see Fig. 5, bottom). For the first depolarization within the train the mean rundown was  $(4.1 \pm 0.1)$  % per min and  $(3.5 \pm 0.2)$  % per min for multi-hole and single-hole QPlates, respectively.



**Fig. 5:** Example current traces used for quantifying rundown. Top: 1<sup>st</sup> (blue) and 50<sup>th</sup> (green) depolarization in voltage protocol 2. Bottom: 1<sup>st</sup> (blue) and 25<sup>th</sup> (green) depolarization train in voltage protocol 3. Rundown was quantified for the first depolarization of the train.

## Conclusion

Ca<sub>v</sub>1.2 Ca<sup>2+</sup> currents were recorded using the QPatch on both single-hole and multi-hole QPlates. 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 rundown as low as **~2% per min** and high success rates were achieved (up to **71%** and **98%** for single-hole and multi-hole QPlates, respectively) with cells at passage numbers below 20.

## Methods

Experiments in this study were performed on HEK-hCa<sub>v</sub>1.2 cells, kindly provided by SB Drug Discovery. The cell line is stably transfected with the Ca<sub>v</sub>1.2 channel and cultured in 30 μM verapamil according to SB Drug Discovery guidelines. HEK-hCa<sub>v</sub>1.2 cells were harvested according to Sophion standard procedures.

## References:

1. Kuryshev, Y. A., Brown, A. M., Duzic, E. & Kirsch, G. E. Evaluating State Dependence and Subtype Selectivity of Calcium Channel Modulators in Automated Electrophysiology Assays. *ASSAY Drug Dev. Technol.* **12**, 110–119 (2014).
2. Balasubramanian, B. et al. Optimization of Ca<sub>v</sub>1.2 screening with an automated planar patch clamp platform. *J. Pharmacol. Toxicol. Methods* **59**, 62–72 (2009).
3. Sinneger-Brauns, M. J. et al. Isoform-specific regulation of mood behavior and pancreatic β cell and cardiovascular function by L-type Ca<sup>2+</sup> channels. *J. Clin. Invest.* **113**, 1430–1439 (2004).
4. Moosmang, S. et al. Dominant role of smooth muscle L-type calcium channel Ca<sub>v</sub>1.2 for blood pressure regulation. *EMBO J.* **22**, 6027–6034 (2003).
5. Zhang, J. et al. Role of Ca<sub>v</sub>1.2 L-type Ca<sup>2+</sup> channels in vascular tone: effects of nifedipine and Mg<sup>2+</sup>. *AJP Heart Circ. Physiol.* **292**, H415–H425 (2006).
6. Zamponi, G. W. Targeting voltage-gated calcium channels in neurological and psychiatric diseases. *Nat. Rev. Drug Discov.* **15**, 19–34 (2015).
7. Liao, P. & Soong, T. W. Ca<sub>v</sub>1.2 channelopathies: from arrhythmias to autism, bipolar disorder, and immunodeficiency. *Pflüg. Arch. - Eur. J. Physiol.* **460**, 353–359 (2010).
8. Hofmann, F., Flockerzi, V., Kahl, S. & Wegener, J. W. L-Type Ca<sub>v</sub>1.2 Calcium Channels: From In Vitro Findings to In Vivo Function. *Physiol. Rev.* **94**, 303–326 (2014).
9. Reuter, H. Calcium channel modulation by neurotransmitters, enzymes and drugs. *Nature* **301**, 569–574 (1983).
10. Hennessey, J. A. et al. A CACNA1C Variant Associated with Reduced Voltage-Dependent Inactivation, Increased Ca<sub>v</sub>1.2 Channel Window Current, and Arrhythmogenesis. *PLoS ONE* **9**, e106982 (2014).
11. Elkins, R. C. et al. Variability in high-throughput ion-channel screening data and consequences for cardiac safety assessment. *J. Pharmacol. Toxicol. Methods* **68**, 112–122 (2013).
12. Zhang, H. Y., Liao, P., Wang, J. J., Yu, D. J. & Soong, T. W. Alternative splicing modulates diltiazem sensitivity of cardiac and vascular smooth muscle Ca<sub>v</sub>1.2 calcium channels: Alternative splicing affects diltiazem block. *Br. J. Pharmacol.* **160**, 1631–1640 (2010).

## Author:

Kadla R Rosholm, Application scientist