

## Application Report

# Internal solution exchange on Qube 384

Internal addition of compounds targeting the chloride channel CIC-1. A robust assay with good pharmacology and high Z-score

### Summary

- A robust assay of CIC-1 with Z' constantly >0.5
- Biophysical and pharmacological characteristics as expected for CIC-1
- Qube offers the possibility to apply compounds both intra- and extracellularly
- Advanced analysis of compound effects on channel kinetics are available at a few clicks

### Introduction

The family of CLC proteins comprises both chloride channels and transporters that are expressed in a great variety of tissues. Subtype 1 of the CLC family (CIC-1) constitutes a voltage-gated chloride channel that is predominantly expressed in skeletal muscle fibers. In these cells, CIC-1 channels are key players in repolarizing the membrane potential and propagating action potentials that eventually regulate muscle contraction. Consequently, it is not surprising that malfunction of the channel is related to disease states. For example, the hereditary disease myotonia that is characterized by muscle membrane hyperexcitability was linked to mutations in the CLCN1 gene that codes for CIC-1 (Miryounesi, Ghafouri-Fard, & Fardaei, 2016). Aside from its relevance in muscle cells, a recent study revealed a role of CIC-1 in epilepsy (Chen et al., 2013).

CIC-1 has a complex gating, channels are time- and voltage-dependently activated at positive potentials and the channel is strongly inwardly rectifying. CIC-1 has two distinct gating mechanisms that are governed by a fast (protopore) and a slow (common) gate. The fast gate is about 10-fold faster than the common gate, a fact that has often been exploited by scientists to study the two gates individually. The channel is further sensitive to pH, anions and intracellular ATP (Jentsch, 2015).

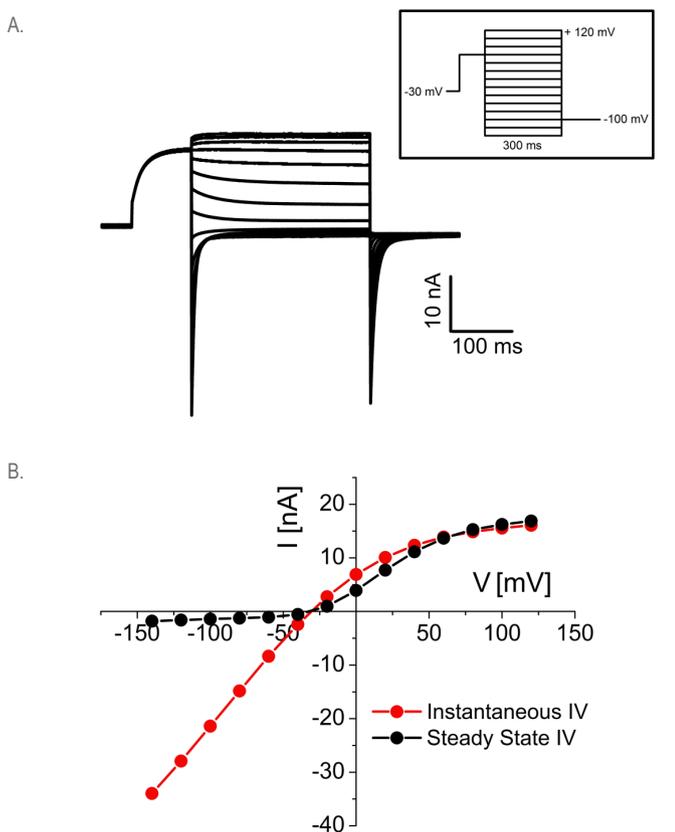
Albeit CIC-1 has been subject to many investigations over the past three decades, its pharmacology is only poorly developed with few moderately selective inhibitors and activators available at the moment. Many pharmacological tool agents targeting CIC-1 bind to the channel from the cytosolic side. This is not a problem as long as the compound is sufficiently membrane permeable. The classical CIC-1 inhibitor 9-anthracen carboxylic acid (9-AC) is highly lipophilic and hence passes the plasma membrane with ease. A fact that allows extracellular application of the compound in a patch clamp experiment. However, possible new compounds that also bind to CIC-1 from the cytosolic side may show different characteristics. Therefore, an assay that is limited to compound application to the external side may result in false negatives. Qube offers the possibility to exchange both extracellular and intracellular solution during an experiment making it well suited to study CIC-1 pharmacology.

In this report, we show a CIC-1 assay on Qube that shows biophysical characteristics as expected for CIC-1. We further introduce Sophion's Analyzer software that allows rapid analysis of large data sets to answer advanced electrophysiological questions, in the present case: What is the mode of action of a novel, unknown compound?

### Results and discussion

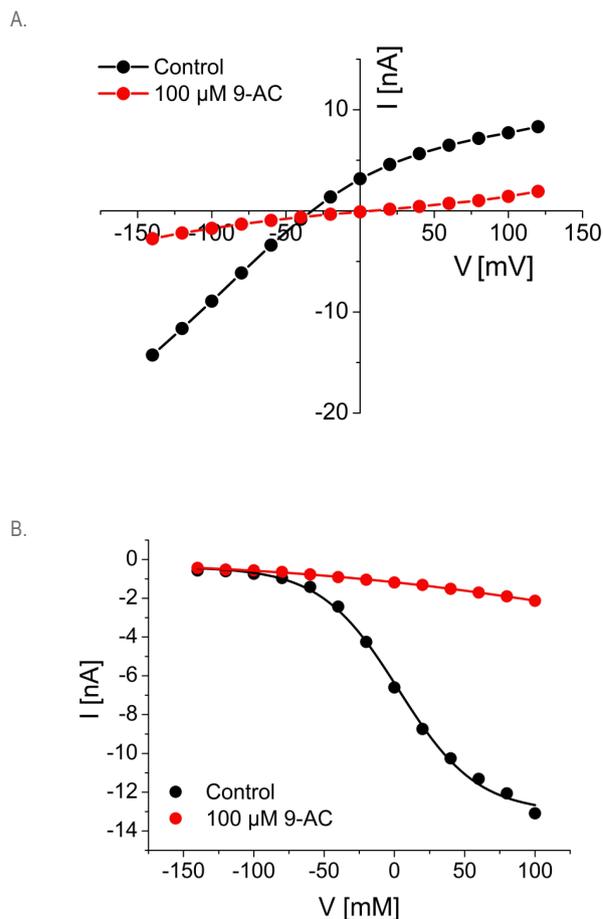
To characterize CIC-1 biophysically, we applied a series of voltage steps to the cells in the whole cell configuration. Typical current response traces are shown in Figure 1A. The respective current voltage relationships of the instantaneous (red) and steady state current (black) are shown in Figure 1B. A large inward current was recorded instantaneously, representing the current passing through all channels open at +60 mV. Due to the voltage sensitive nature of CIC-1, almost no inward current was detected in the steady state.

Different pharmacological tool agents are available that either act as gating modifier (i.e. shift the equilibrium between open and closed channels towards more depolarized potentials) or as pore blocker. CIC-1's intimate relationship between permeability and gating suggests that any pore blocking agent will also modify CIC-1 gating i.e. shift the half-activation potential ( $V_{1/2}$ ) towards more positive potentials. 9-AC acts as pore blocker and was shown to bind to the pore from the intracellular side (Estevez, Schroeder, Accardi, Jentsch, & Pusch, 2003).



**Fig. 1:** Typical CIC-1 whole cell voltage clamp recording using multi hole QChips. A: Current traces measured using the voltage protocol shown in the insert. B: Respective current voltage relationship. Instantaneous current was determined 1 ms after the voltage step. Steady state currents were measured at the end of the voltage step.

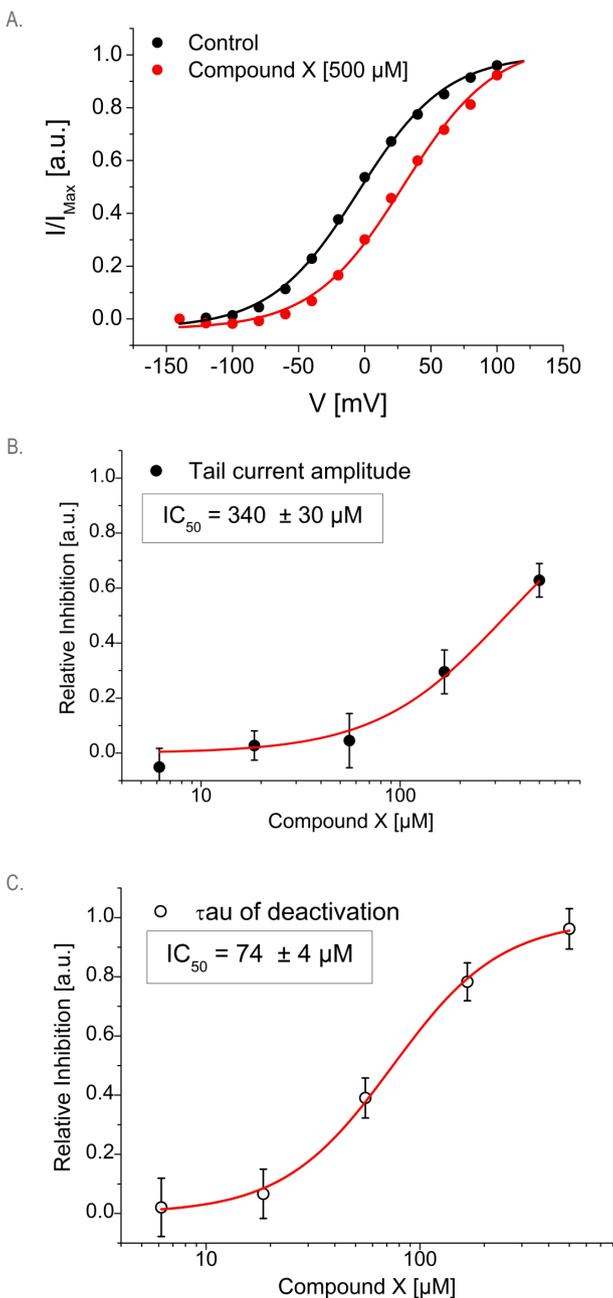
Figure 2 shows the effect of intracellularly applied 9-AC (100  $\mu$ M) on CIC-1 conductance. Almost no current was mediated through CIC-1 channels over the entire voltage spectrum following 9-AC exposure (Figure 2A). Analysis of the inactivation curve (Figure 2B) revealed that the compound in fact acts predominantly as pore blocker as even at +100 mV, almost no tail currents were detected.



**Fig. 2:** 9-AC is a CIC-1 pore blocker. A: Typical instantaneous IV curve recorded in control condition (black circles) and after internal application of 9-AC (100  $\mu$ M) (red circles). B: Tail current amplitude recorded at  $V = -100$  mV plotted against voltage of applied pre-pulse. Tail currents were almost completely inhibited in the presence of 9-AC.

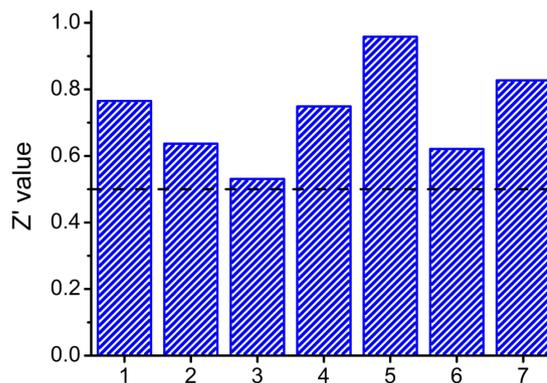
We used a compound (Compound X) that is known to act as gating modifier to demonstrate the assay's capability to differentiate between pore block and gating modification. To assess the compound's effect on CIC-1 gating, a set of voltage steps from -140 mV to +120 mV in 20 mV increments was applied to the cells and tail currents at -100 mV were used to generate deactivation curves (Figure 3A). The half activation potential ( $V_{1/2}$ ) in the control condition was calculated to  $V_{1/2}(\text{Control}) = -4$  mV. This value is more positive than values reported in the literature as the used internal solution contained a relatively low concentration of Cl<sup>-</sup> and  $V_{1/2}$  strongly depends on ionic strength (Accardi & Pusch, 2000). Exposure to Compound X (500  $\mu$ M) caused a shift in  $V_{1/2}$  by 27 mV to  $V_{1/2}(\text{Compound X}) = +23$  mV. To further investigate the mode of action of compound X, we recorded a dose response curve both on tail current amplitude and on deactivation kinetics represented

by Tau of a monoexponential fit to the tail currents (Figure 3B and C) (see methods). Compound X exhibited a ~5-times higher potency on deactivation kinetics (Figure 3C) compared with open channel inhibition (Figure 3B). This suggests that the compound stabilizes the closed state of CIC-1 rather than inhibiting the conducting pore.



**Fig. 3:** Compound X inhibits CIC-1 currents by modulating channel gating. A: Representative inactivation curve recorded at the tail current in the presence ( $V_{1/2}(9\text{-AC}) = +23\text{ mV}$ ) and absence ( $V_{1/2}(\text{Control}) = -4\text{ mV}$ ) of cytosolic compound X (500  $\mu\text{M}$ ). B: Noncumulative dose response curve obtained from the tail current amplitude of a -20 mV pre-pulse and normalized to control period and full block period ( $n = 6\text{-}9$  per concentration). C: Noncumulative dose-dependent effect of compound X on CLC-1 deactivation kinetics represented by Tau (Monoexponential fit to the tail current of a +120 mV pre-pulse).

Finally, we validated the assay using  $Z'$  analysis.  $Z'$  values were determined from a column with 0.1% DMSO control versus a column with internally applied 100  $\mu\text{M}$  9-AC.  $Z'$  was  $> 0.5$  in 7 individual experiments (Figure 4).



**Fig. 4:**  $Z'$  values were calculated from normalized current inhibition values obtained from column 23 (0.1% DMSO control) and column 24 (9-AC 100  $\mu\text{M}$  applied intracellularly) of a 384-well plate.

## Conclusion

As illustrated above by means of the challenging CIC-1 assay, Qube offers a high degree of flexibility in the assay design. In addition to this, the related Qube software, Analyzer, makes it very easy to rapidly set up advanced analyzing templates that, once generated, can be applied to every newly generated data. Taken together this makes Qube an ideal platform for automated high-throughput patch clamp studies.

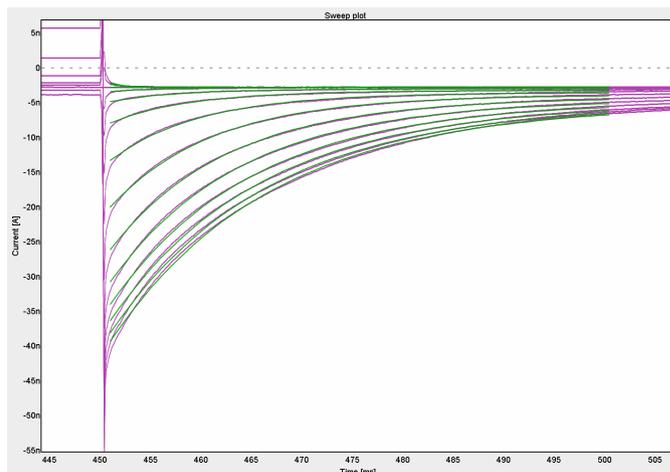
## Methods

### Cells and Cell Culture

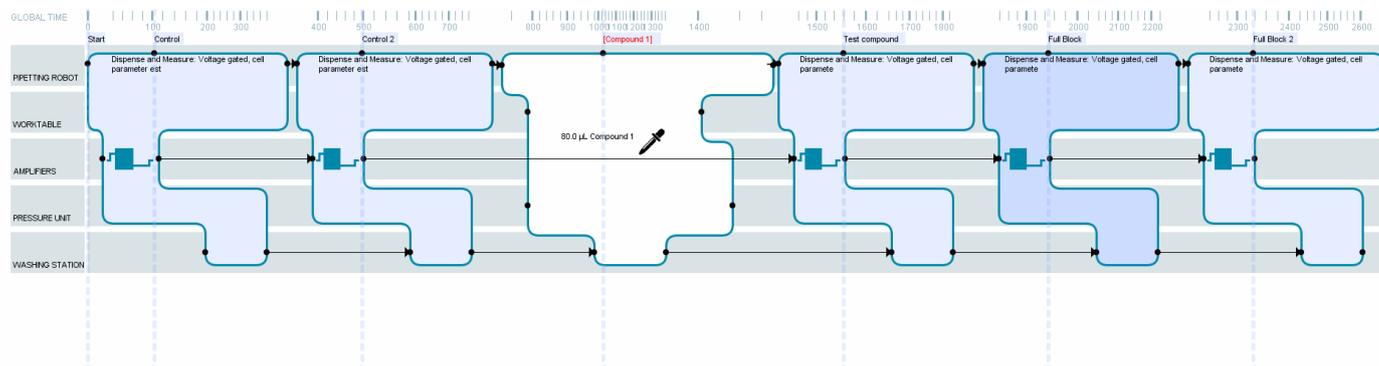
CHO cells stably expressing hCIC-1 were provided by ChanTest. Cells were cultured following the manufacturer's guidelines and harvested according to Sophion's standard procedure for CHO cells including Trypsin.

### Electrophysiology

Whole cell patch clamp recordings were performed on Qube following Sophion's standard procedure using multi hole QChips (10 holes per well). CIC-1 currents were elicited using a 100 ms long voltage pre step to +60 mV followed by a series of voltage steps (300 ms) ranging from -140 mV to +120 mV in 20 mV increments. Tail currents were analyzed at -100 mV.



**Fig. 5:** Monoexponential fit to the tail current. The function was fitted to an interval between 1 and 51 ms following the voltage step. Compound effects on deactivation kinetics were assessed by analyzing tau. Compound effects on tail current amplitude were analyzed at 1 ms after a +120 mV voltage pre-pulse.



**Fig. 6:** Experimental setup of the CIC-1 assay. Each block represents the addition of either extracellular (blue) or intracellular (white) solution. Each addition is followed by a voltage- (or current-clamp) recording. A small representation of the applied protocol is shown in each block. Two baseline recordings were followed by an internal solution exchange where the test compound was applied. Compounds were further applied extracellularly before the next measurement was taken. The experiment was finished with two subsequent applications of 100  $\mu$ M 9-AC that was used as reference compound.

Cells provided by:



### Author:

Daniel Sauter, Application scientist

Sophion Bioscience A/S, Baltorpvej 154, 2750 Ballerup, Denmark  
Phone: +45 4460 8800 Fax: +45 4460 8899, E-mail: info@sophion.com

sophion.com

### References:

- Accardi, A., & Pusch, M. (2000). Fast and slow gating relaxations in the muscle chloride channel CIC-1. *J. Gen. Physiol.*, 116(3), 433–444. <http://doi.org/10.1085/jgp.116.3.433>
- Chen, T. T., Klassen, T. L., Goldman, A. M., Marini, C., Guerrini, R., & Noebels, J. L. (2013). Novel brain expression of CIC-1 chloride channels and enrichment of CLCN1 variants in epilepsy. *Neurology*, 80(12), 1078–1085. <http://doi.org/10.1212/WNL.0b013e31828868e7>
- Estevez, R., Schroeder, B. C., Accardi, A., Jentsch, T. J., & Pusch, M. (2003). Conservation of chloride channel structure revealed by an inhibitor binding site in CIC-1. *Neuron*, 38(1), 47–59. <http://doi.org/S0896627303001685> [pii]
- Jentsch, T. J. (2015). Discovery of CLC transport proteins: cloning, structure, function and pathophysiology. *The Journal of Physiology*, 0(Mdc), n/a-n/a. <http://doi.org/10.1113/jphysiol.2014.270043>
- Miryounesi, M., Ghafouri-Fard, S., & Fardaei, M. (2016). A novel missense mutation in CLCN1 gene in a family with autosomal recessive congenital myotonia. *Iranian Journal of Medical Sciences*, 41(5), 456–458.