

Application Report:

# CHO-K<sub>v</sub>1.5



## Voltage gated potassium channels



The voltage gated potassium channel K<sub>v</sub>1.5 is a homotetrameric protein present in the heart. It is a delayed rectifier, participating in the early phase of the heart action potential. This report shows data from CHO cells stably expressing K<sub>v</sub>1.5 tested on the QPatch platform. The cells are obtained through a collaboration with STZ (Germany).

AR\_PUBLIC15419-5

---

## Introduction

The voltage gated potassium channel  $K_v1.5$  is a homotetrameric protein present in the heart. It is a delayed rectifier, participating in the early phase of the heart action potential. This report shows data from CHO cells stably expressing  $K_v1.5$  tested on the QPatch platform.

## Materials & Methods

Intracellular saline (in mM): 5.374  $CaCl_2$ , 1.75  $MgCl_2$ , 3.125/10 KOH/EGTA, 10 HEPES, 120 KCl, 4  $Na_2$ -ATP, pH 7.2 with KOH, 285-296 mOsm. Extracellular saline (in mM): 2  $CaCl_2$ , 1  $MgCl_2$ , 10 HEPES, 4 KCl, 145 NaCl, 10 Glucose, pH 7.4 with NaOH, ~305 mOsm.

Cells: CHO cells stably expressing  $K_v1.5$  were obtained from STZ (Mannheim, Germany). Cells were cultured and harvested for QPatch experiments as described in the Sophion SOP. Data shown here is from STZ CHO- $K_v1.5$  clone 16.

## Results

Experiments were conducted to evaluate the IV-relationship of  $K_v1.5$  as well as dose-response for inhibitors.

Figure 1 shows the currents elicited at potentials ranging from -90 mV to +50 mV in a representative experiment with CHO-KV1.5. The corresponding IV plot for both maximum and steady-state current is shown in Figure 2.

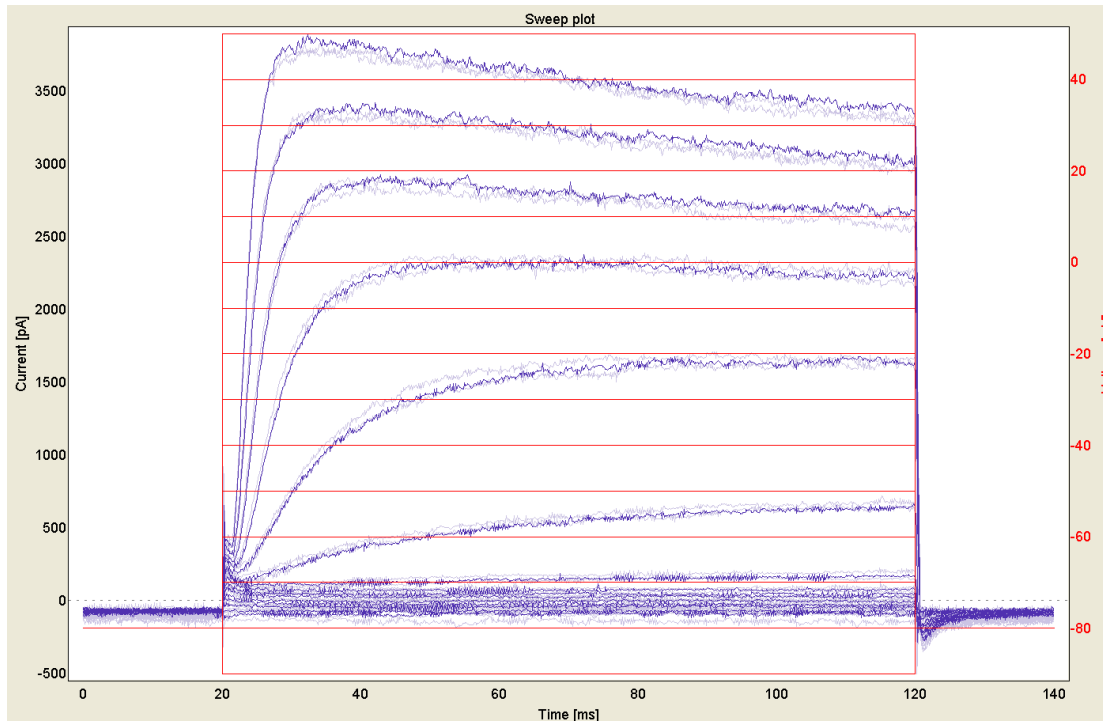


Figure 1. Kv1.5 raw data sweeps elicited in an IV-protocol with steps ranging from -90 to +50 mV.

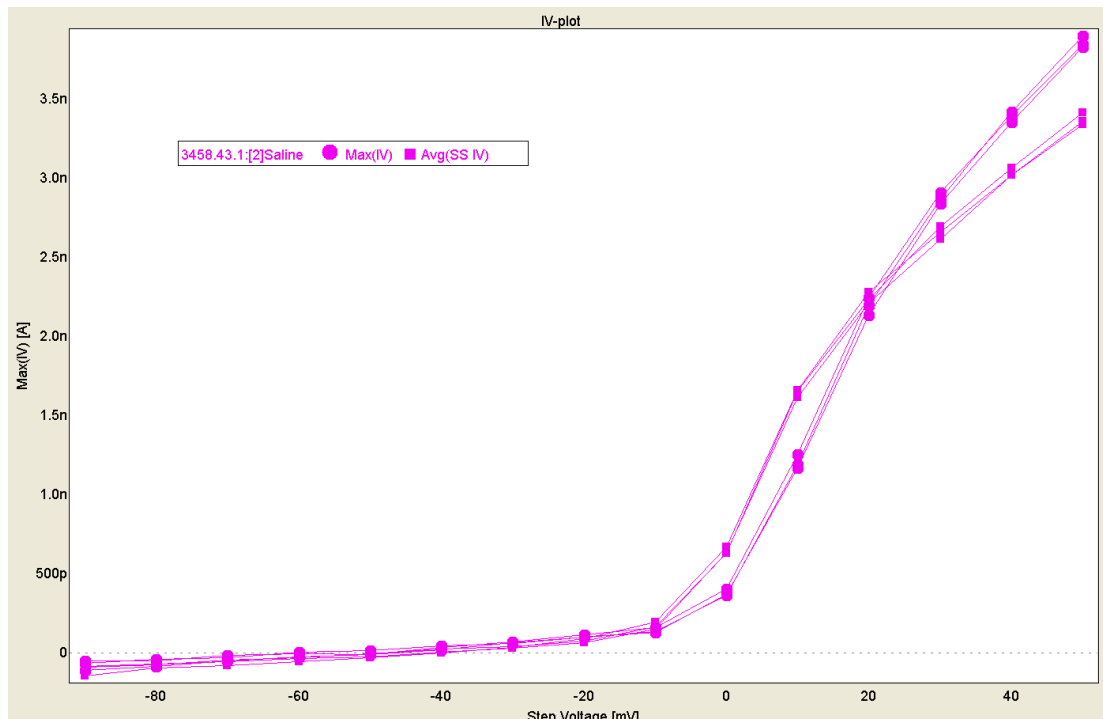


Figure 2. Current-voltage relationship (IV plot) of the data shown in Figure 1. Circles show the maximum elicited current, squares show the steady-state current.

The response of  $K_v1.5$  to a known blocker was also tested. Figure 3 shows the raw data traces of the steady-state response to six different concentrations of 4-aminopyridine. Figure 4 and Figure 5 show the corresponding current versus time (IT) plot and Hill fit, respectively. The resulting  $IC_{50}$  for 4-aminopyridine is  $63.8 \mu M$ .

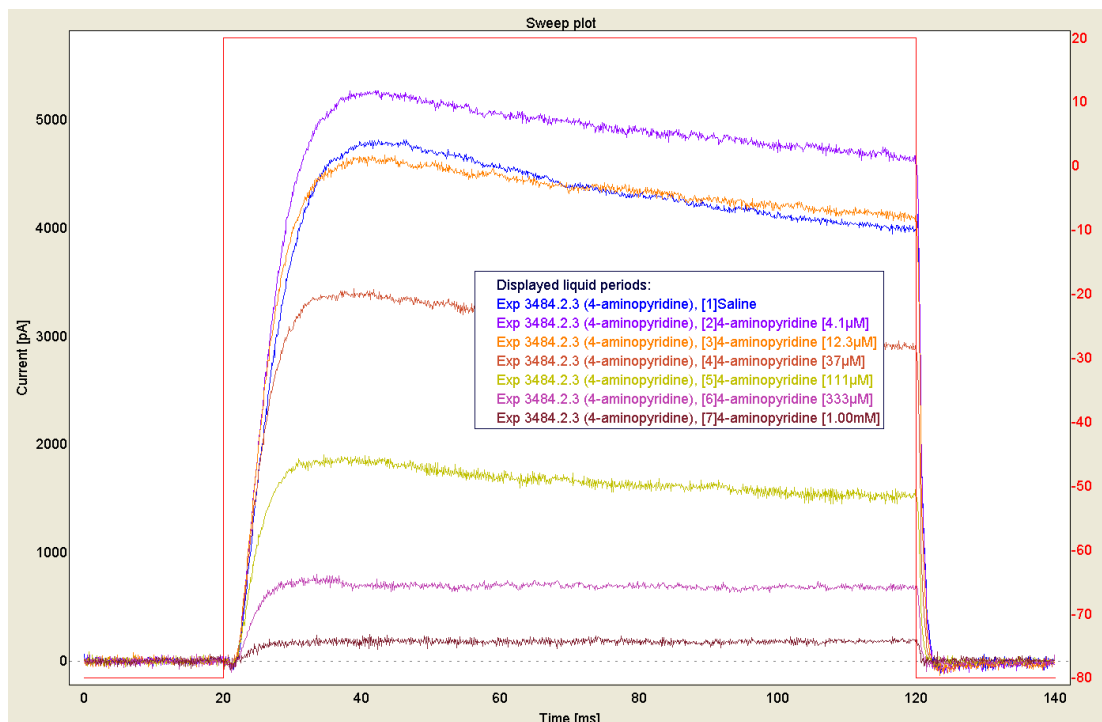


Figure 3. Six point cumulative dose-response experiment with 4-aminopyridine.

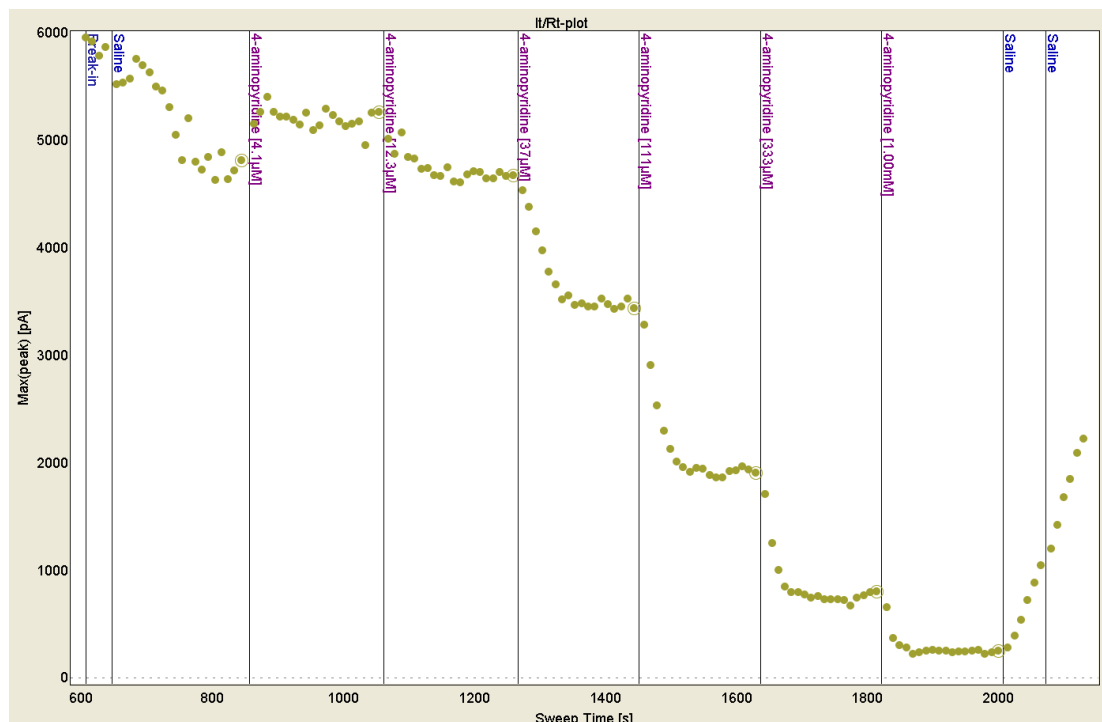


Figure 4. IT plot of  $KV1.5$  channel response to increasing concentrations of 4-aminopyridine.

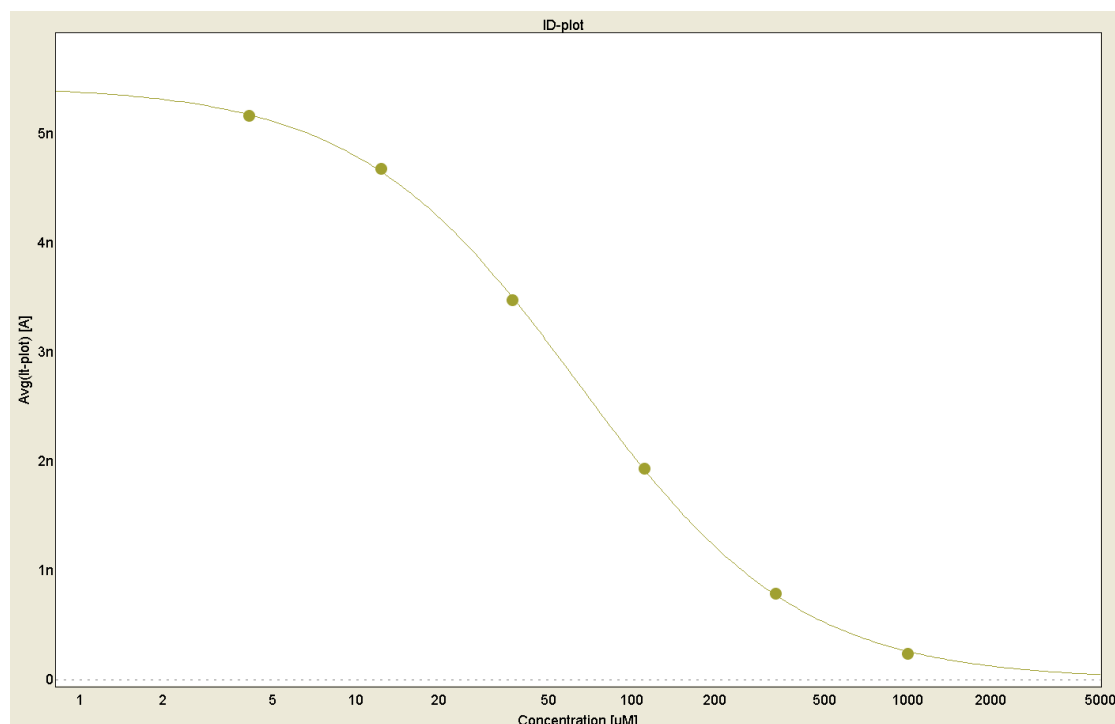


Figure 5. Dose-response plot, with Hill fit, of steady-state current level at six concentrations of 4-aminopyridine.

## Conclusion

IV characteristics and dose-response experiments with a  $K_v1.5$  channel blocker was successfully obtained using QPatch.  $K_v1.5$  shows its characteristic outward rectification and an  $IC_{50}$  for 4-aminopyridine within range of reported literature values (e.g. Gutman et al., *Pharmacological Reviews* 57:473-508, 2005, 270  $\mu$ M).