

# Automated Patch Clamp assay for proton-activated hTOP1 channel

Sarah Jane Lilley<sup>4</sup>, Joshua P. Kaplan<sup>1</sup>, Zachary Krieger<sup>2</sup>, Emily R. Liman<sup>1</sup>, Daniel Sauter<sup>3</sup>

## Introduction

Manual patch clamp is the gold standard for ion channel studies but is low-throughput and labor-intensive. Automated patch clamp (APC) systems, like the 8-well QPatch Compact (QPC) with rapid solution exchange capabilities, overcome these limitations, enabling efficient study of channels. Central to the QPC's functionality is a microfluidic flow channel that ensures precise and controlled delivery of solutions. This microfluidic design facilitates complete and rapid solu-

tion exchange around the cell, a critical feature for accurately studying channels activated by transient stimuli such as protons. OTOP1 is a proton-selective ion channel crucial for acid sensing in tissues like the inner ear and taste buds, acting as a sour taste receptor and ammonium sensor. Characterizing OTOP1's biophysical properties is vital to understand its physiological roles and therapeutic potential.

## Conclusion

This study demonstrates the successful implementation of an automated patch clamp assay using the QPatch Compact system to characterize proton-activated hTOP1 channels. The system enabled rapid and reproducible solution exchange, critical for studying transient proton-evoked responses. Our results confirm that hTOP1 mediates inward currents in response to acidic pH and ammonium, and that these currents are modulated by extracellular zinc in a

concentration-dependent manner. The fitted  $IC_{50}$  for  $Zn^{2+}$  inhibition aligns with previous findings, validating the assay's reliability. This automated approach offers a scalable and efficient platform for future pharmacological screening and mechanistic studies of proton-sensitive ion channels.

## Materials and methods

**Automated Patch Clamp:** All data were recorded using the QPatch Compact System (Sophion Bioscience A/S). All experiments were conducted using QPlate8X consumables.

**Cell preparation:** HEK293 cells stably expressing human OTOP1 channels were harvested either using detachin or trypsin. Cells were transferred to serum-free medium (EX-CELL<sup>®</sup> ACF CHO Medium, Sigma-Aldrich) supplemented with 25 mM HEPES. The cells were washed and resuspended in extracellular buffer immediately before the experiment.

**Solutions:** Extracellular solution used for priming (in mM): 145 NaCl, 4 KCl, 1  $MgCl_2$ , 2  $CaCl_2$ , 10 HEPES, 10 Glucose, pH 7.4.

During the experiment, NMDG-based extracellular solutions consisted of (in mM): 160 NMDG, 2  $CaCl_2$ , 10 CHES

(N-cyclohexyl-2-aminoethanesulfonic acid) for pH 10 / 10 HEPES (2-(4-(2-hydroxyethyl)-1-piperazinyl)ethanesulfonic acid) for pH 7.4 / 10 MES (2-Morpholinoethanesulfonic acid) for pH 6.0 / 10 or 100 Homo-PIPES (homopiperazine-1,4-bis(2-ethanesulfonic acid)) for pH 5.0 and 4.0.

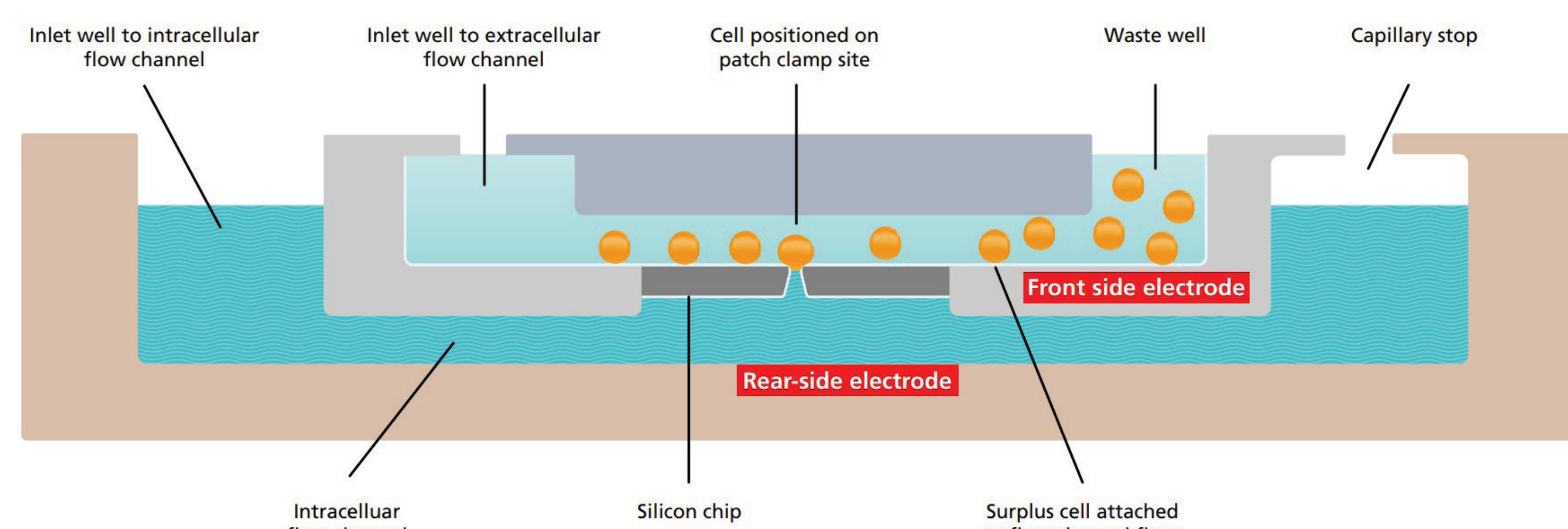
$NH_4Cl$  solution consisted of (in mM): 160  $NH_4Cl$ , 2  $CaCl_2$ , 10 HEPES.

pH of solutions was adjusted with NMDG-OH and HCl.

Intracellular solution (in mM): 140 CsF, 1/5 EGTA/CsOH, 10 HEPES, 10 NaCl, pH 7.3

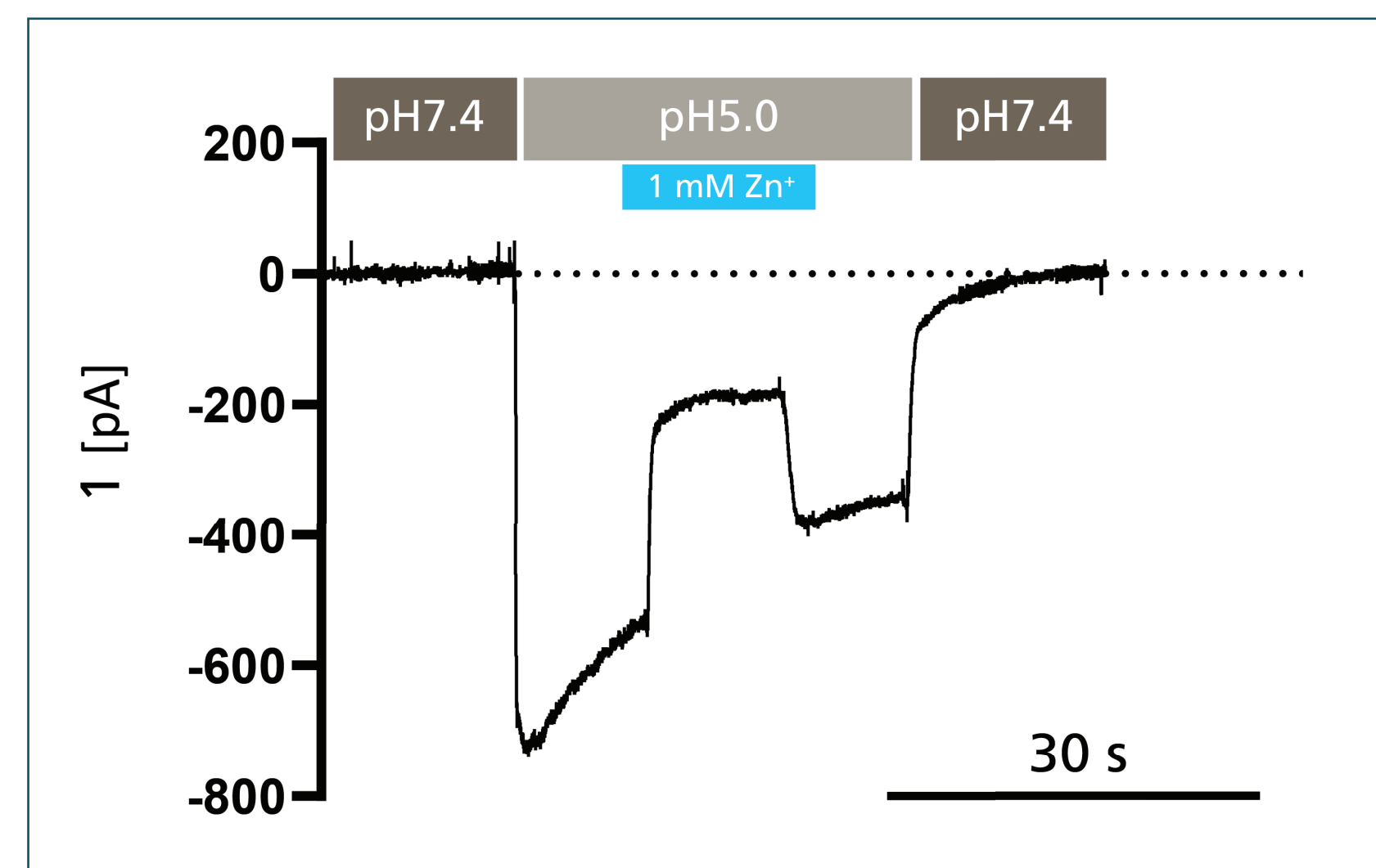
**Whole-cell protocol:** The whole-cell configuration was established using a standard whole-cell protocol for CHO-cells.

**Voltage protocol:** A constant holding potential  $V=-80$  mV was applied unless noted otherwise.

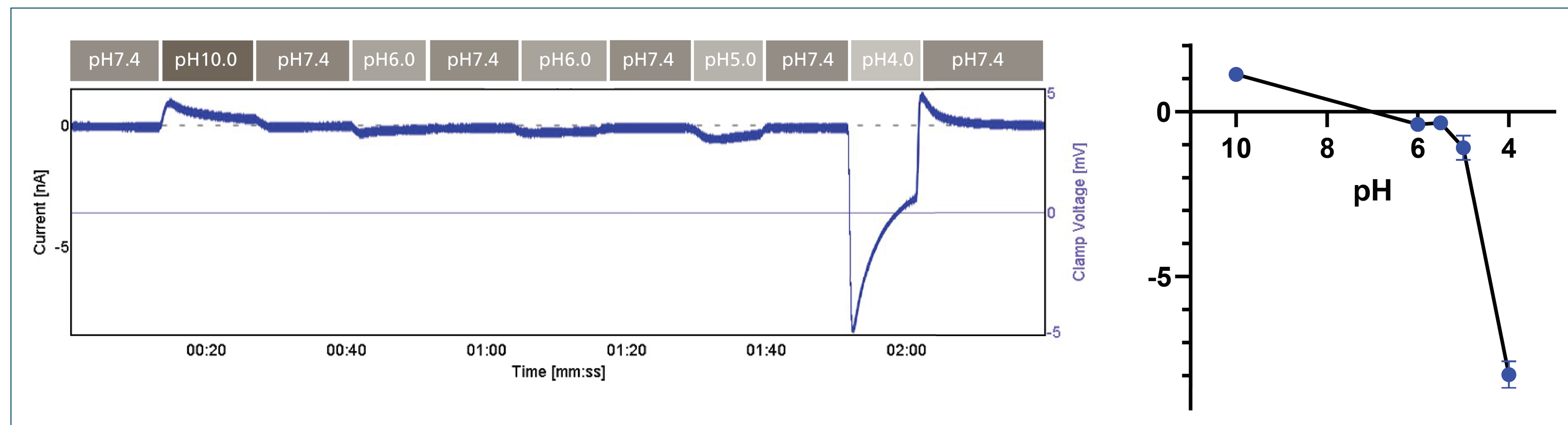


**Fig. 1:** Schematic representation of a microfluidic flow channel. A cell suspension is injected into the external microfluidic channel, while negative pressure is applied to capture a single cell at the patch hole. A controlled pressure sequence then ruptures the membrane patch at this site, establishing stable electrical access to the cell. Each consumable is equipped with built-in Ag/AgCl electrodes, ensuring precise and reproducible current and voltage recordings.

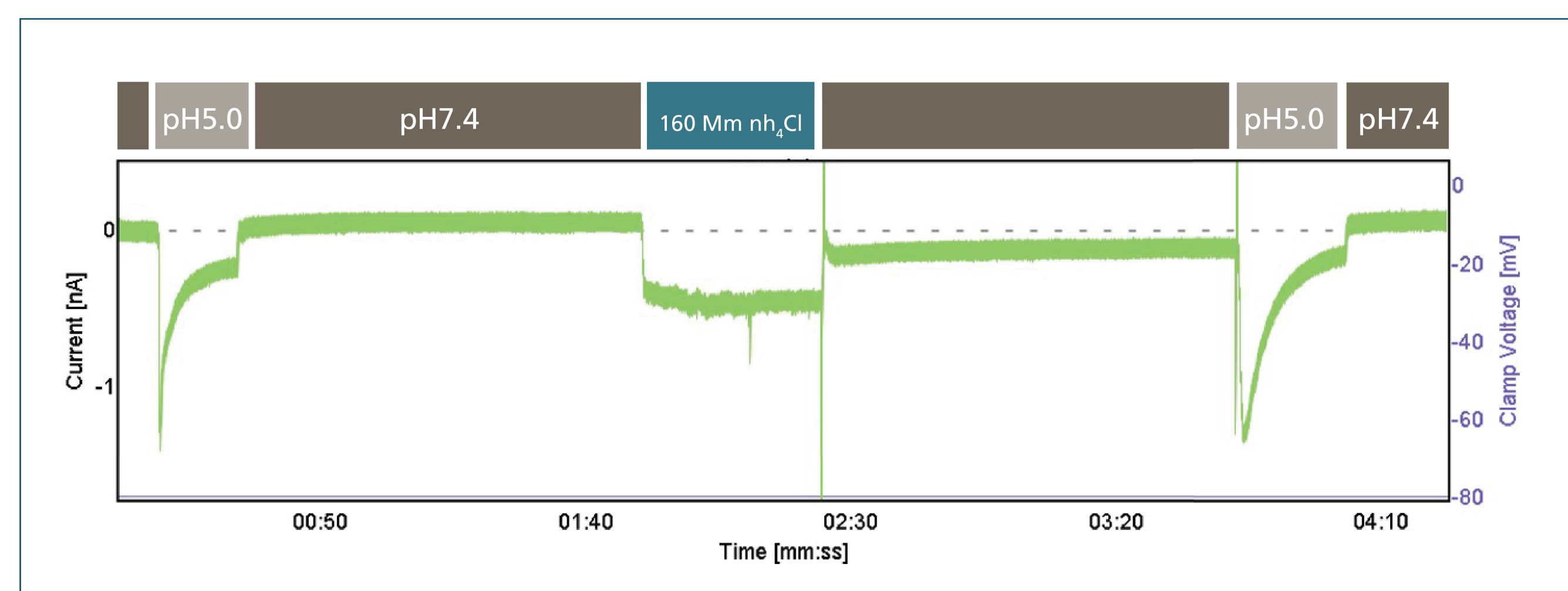
## Results



**Fig. 2:** Acidic pH evokes an inward current inhibited by zinc in HEK293-hTOP1 cells. Representative current trace recorded from HEK293-hTOP1 cells. External solution was rapidly exchanged from pH7.4 to pH5.0 using QPC's microfluidic flow channel. Subsequent application of 1 mM  $Zn^{2+}$  resulted in a significant reduction of the acid-evoked current. Wash out of  $Zn^{2+}$  recovered most of the acid-activated current.



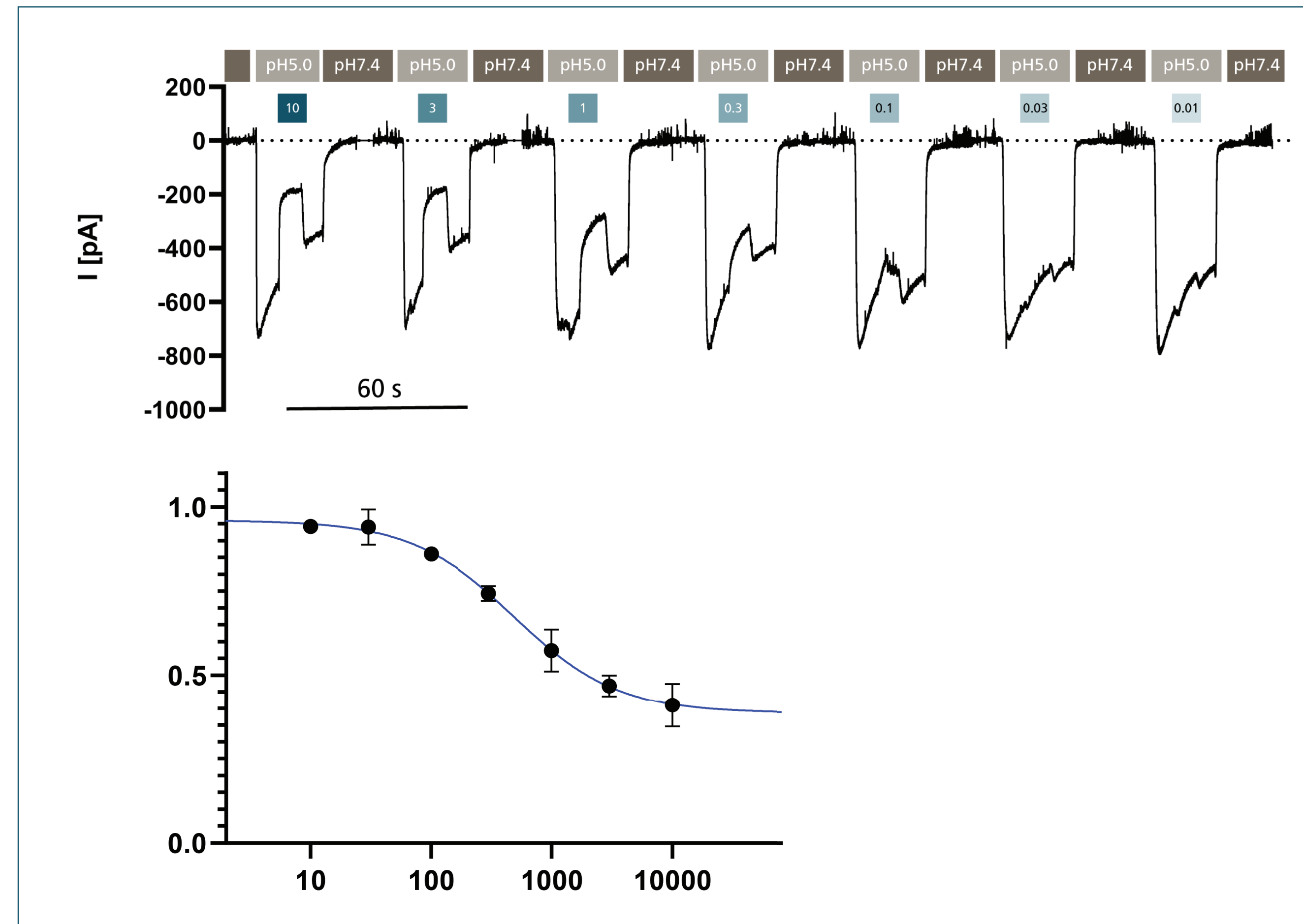
**Fig. 3:** pH-dependence of hTOP1 current in HEK293 Cells. Representative current trace (left) and average pH-response plot (right) illustrating hTOP1 activity. Cells were voltage-clamped at  $V_m=0$  mV to reduce proton driving force. External solutions with pH values ranging from 10.0 to 4.0 were applied to stimulate the cells. The bottom panel shows the average peak current responses ( $N=5$  cells, mean  $\pm$  SEM) across various pH stimulations.



**Fig. 4:** hTOP1 mediates ammonium-evoked inward currents. Representative current trace demonstrating the electrophysiological response of hTOP1-expressing cells to 160 mM  $NH_4Cl$ . Prior to and following  $NH_4Cl$  stimulation, cells were challenged with a pH 5.0 buffer solution to confirm hTOP1 functional expression. This data corroborates the initial findings from the Liman lab, which first identified hTOP1 as the molecular entity responsible for mediating ammonium-evoked inward currents (Liang et al., 2023, *Nature Communications*).

## References

1.Liang, Z., Wilson, C. E., Teng, B., Kinnamon, S. C., & Liman, E. R. (2023). *The proton channel OTOP1 is a sensor for the taste of ammonium chloride*. **Nature Communications**, 14, Article 6163. <https://doi.org/10.1038/s41467-023-41637-4>



**Fig. 5:** Concentration-dependent inhibition of hTOP1 by zinc. (Top) Representative whole-cell current trace demonstrating the inhibitory effect of varying  $Zn^{2+}$  concentrations on acid-evoked hTOP1 currents. Currents were activated by an external solution at pH5.0. Effect of  $Zn^{2+}$  was assessed by washing in varying concentrations of  $Zn^{2+}$  during the pH5.0 stimulation pulse followed by a wash out step. This protocol was repeated for seven different  $Zn^{2+}$  concentrations (100 nM to 10  $\mu$ M). (Bottom) Average concentration-response curve ( $N=6$  cells; mean  $\pm$  SEM) for  $Zn^{2+}$  inhibition of hTOP1. Residual current was calculated as the ratio of current immediately prior to  $Zn^{2+}$  washout to the current immediately after washout. The fitted  $IC_{50}$  for  $Zn^{2+}$  was 489  $\mu$ M (95% CI: 364 to 695  $\mu$ M).

