

## Cold activation of TRPM8 using the External Perfusion System and the Port-a-Patch

The electrophysiology team at Nanion Technologies GmbH, Munich.

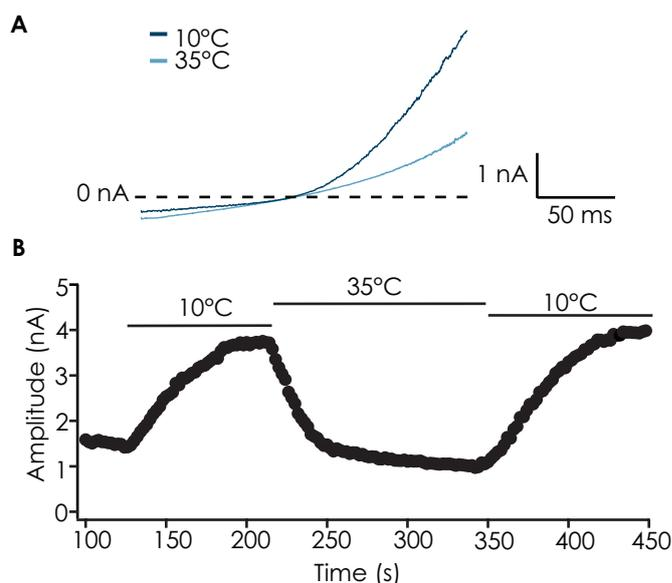
### Summary

TRPM8 is a member of the transient receptor potential channel (TRP) family. TRPM8 is known to be a thermo-sensitive channel, activated by cold temperatures (below  $\sim 25^{\circ}\text{C}$ ) and ligands such as menthol, Eucalyptol and icilin<sup>1-4</sup>. It belongs to the melastatin subfamily of TRP channels<sup>5</sup> and shows an outward rectification with a relatively high permeability for calcium ions and little selectivity between monovalent cations. Menthol, a secondary alcohol produced by the peppermint herb, *Mentha piperita*, is widely used in the food and pharmaceutical industries as a cooling/soothing compound and odorant. It induces  $\text{Ca}^{2+}$  influx in a subset of sensory neurons from dorsal root and trigeminal ganglia, where the TRPM8 channel is specifically expressed<sup>2-4</sup>. It is known that TRPM8 plays a role cold allodynia after inflammation<sup>5</sup> and it has been implicated as a possible pharmaceutical target to treat chronic pain and migraine<sup>6</sup>. Additionally, TRPM8 channels have been implicated to play a role in cancer<sup>7,8</sup>, in particular prostate cancer<sup>7</sup> and pancreatic cancer<sup>8</sup>, and may provide novel clinical biomarkers and therapeutic targets for these types of cancer.

Here we present data of hTRPM8 collected on the Port-a-Patch using cooled solution via the External Perfusion System. Cold activated hTRPM8 was blocked by capsaizepine with an  $\text{IC}_{50}$  in good agreement with the literature<sup>1,9</sup>.

### Results

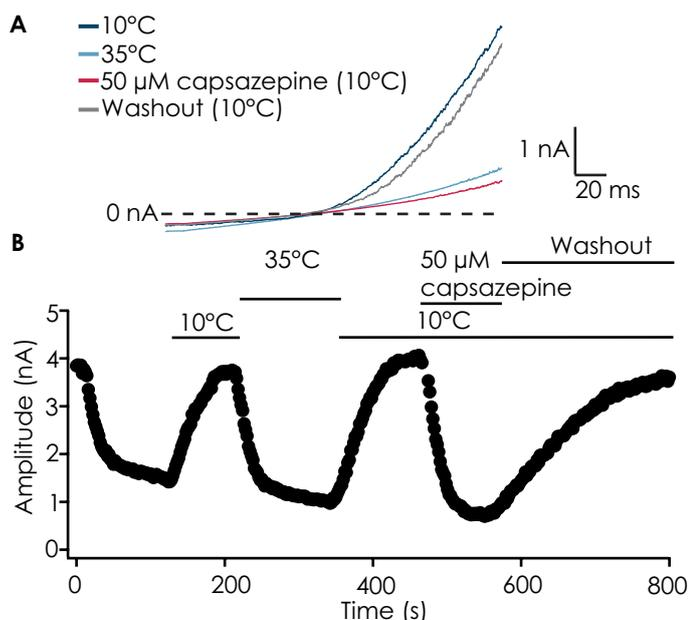
TRPM8 can be activated by a number of ligands including the cooling compounds menthol and icilin<sup>1-4</sup>, as well as cold temperatures<sup>1-4</sup>. TRPM8 is a cation channel with voltage-dependence of activation exhibiting a strong outward rectification at depolarized membrane potentials and fast, voltage-dependent closure at negative potentials<sup>9,10</sup>. TRPM8-mediated currents are suppressed by high temperatures<sup>2,10</sup>. Figure 1 shows activation of TRPM8 expressed in CHO cells by continuous perfusion of external solution at  $10^{\circ}\text{C}$  and subsequent suppression by elevated temperature ( $35^{\circ}\text{C}$ ).



**Figure 1:** **A** Whole cell current responses from induced CHO cells expressing TRPM8 to a ramp protocol ( $-100\text{ mV}$  to  $+100\text{ mV}$  over  $200\text{ ms}$ ) at  $10^{\circ}\text{C}$  (dark blue) and  $35^{\circ}\text{C}$  (light blue). TRPM8 was activated at  $10^{\circ}\text{C}$  but not at  $35^{\circ}\text{C}$ . **B** Timecourse of TRPM8 activation with cold solution ( $10^{\circ}\text{C}$ ) and block by warm solution ( $35^{\circ}\text{C}$ ).

# Application Note

TRPM8 was activated by cooled solution and this could be blocked by either increasing temperature, or by applying the blocker capsazepine via the external solution. Capsazepine was applied at 10°C. Figure 2 shows the time course of the experiment and corresponding example traces.



**Figure 2:** **A** Traces from an example cell showing activation by cooled solution (10°C, dark blue), block by either heated solution (35°C; light blue), or capsazepine (50 µM; red trace) and washout after capsazepine application (grey). TRPM8 current could be almost completely recovered upon washout of capsazepine. **B** Timecourse of the experiment.

## References

1. Behrendt, H.J., *et al.*, 2004. *Br. J. Pharmacol.* 141, 737–745
2. McKemy, D.D., *et al.*, 2002. *Nature* 416: 52–58
3. Peier, A.M., *et al.*, 2002. *Cell* 108: 705–715
4. Andersson, D.A., *et al.*, 2004. *J. Neurosci.* 24(23):5364–5369
5. Weyer, A.D. & Leyto, S.G. 2017. *Pharmaceuticals.* 10: 37
6. González-Muñiz, R., *et al.* 2019. *Int. J. Mol. Sci.* 20: 2618
7. Hantute-Ghesquier, A. *et al.* 2018. *Pharmaceuticals.* 11: 58
8. Yee, N.S. *et al.*, 2012. *Scientifica.* 2012: Article 415158
9. Malkia, A., *et al.*, 2009. *Mol. Pain.* 6: 62
10. Nilius, B., *et al.* 2005. *J. Physiol.* 567.1: 35–44

## Methods

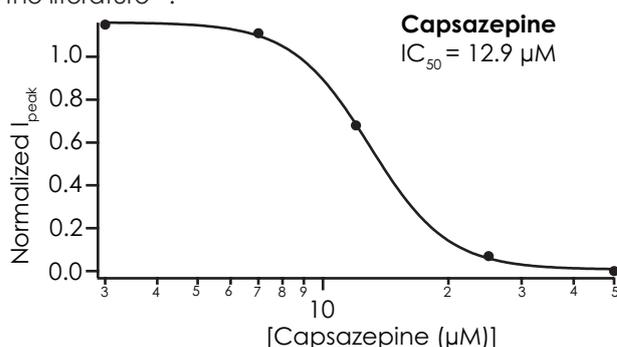
### Cells

CHO cells stably expressing hTRPM8

Nanon Technologies GmbH  
Ganghoferstr. 70A  
80339 Munich, Germany

phone +49 89 219 095-0  
fax +49 89 218997960  
www.nanon.de • info@nanon.de

A concentration response curve for capsazepine using cooled solution (10°C) as the activator could also be constructed. Increasing concentrations of capsazepine at 10°C were applied and the concentration response curve is shown in Figure 3. The  $IC_{50}$  for capsazepine block of TRPM8 was 12.9 µM (n = 1), in excellent agreement with the literature<sup>1,9</sup>.



**Figure 3:** Concentration response curve for capsazepine using cooled solution to activate TRPM8 reveals an  $IC_{50}$  of 12.9 µM, in good agreement with the literature<sup>1,9</sup>.

In conclusion, hTRPM8 expressed in CHO cells can be reliably activated using the External Perfusion System of the Port-a-Patch with Temperature Control which enables cooling of the external solution to 10°C. Cold-activated TRPM8 was also blocked by capsazepine as expected<sup>1,9</sup>. The Port-a-Patch is therefore a useful tool for identifying TRPM8 inhibitors as potential therapeutics for chronic pain and some cancers.

### Cell culture

CHO cells stably expressing hTRPM8 were cultured using standard culture conditions for the Port-a-Patch and induced using 1 mg/ml tetracycline for at least 24–36 hours prior to measurements.

### Electrophysiology

Whole cell patch clamp recordings were conducted according to Nanion's standard procedure for the Port-a-Patch. Currents were elicited by 200 ms voltage ramps from -100 mV to +100 mV,  $V_{hold} = -80$  mV. External solution was cooled to 10°C using the Temperature Control of the External Perfusion System to activate hTRPM8. Capsazepine concentrations were made up in external solution and applied to the cell at 10°C using the External Perfusion System of the Port-a-Patch.