

TRPM8

SyncroPatch 384

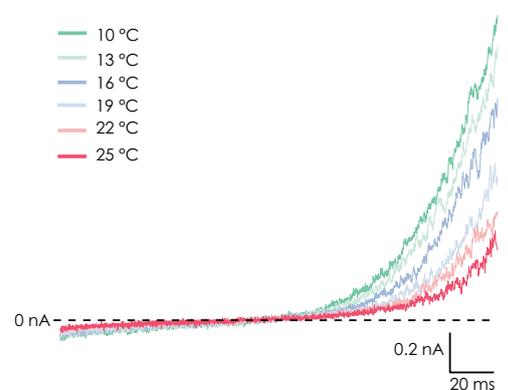
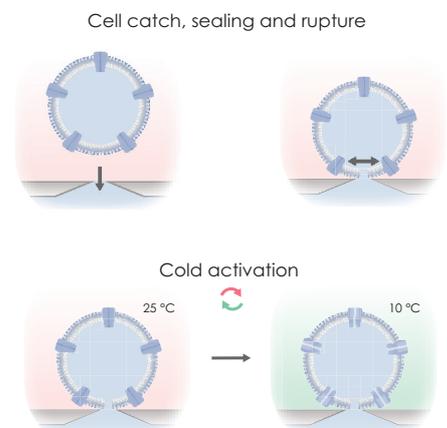
Cool application: activation of TRPM8 on the SyncroPatch 384

TRPM8 is located on the plasma membrane of a subset of unmyelinated (C-type) afferents in the dorsal root and trigeminal ganglion. Activated by the ligands menthol, icilin and eucalyptol, and also by low temperatures, TRPM8 plays a critical role in the detection of environmental cold temperatures and has been implicated in chronic and neuropathic pain. Therefore, it is a potential target in drug discovery, and there is a need for reliable methods to record TRPM8 activated using different stimuli to find TRPM8 modulators.

Nanonion's SyncroPatch 384 is equipped with a sophisticated temperature control that can record at temperatures as low as 10 °C, reliably activating TRPM8. The temperature can be adjusted at any time during the experiment for full flexibility of the experimental design.



Contact us today

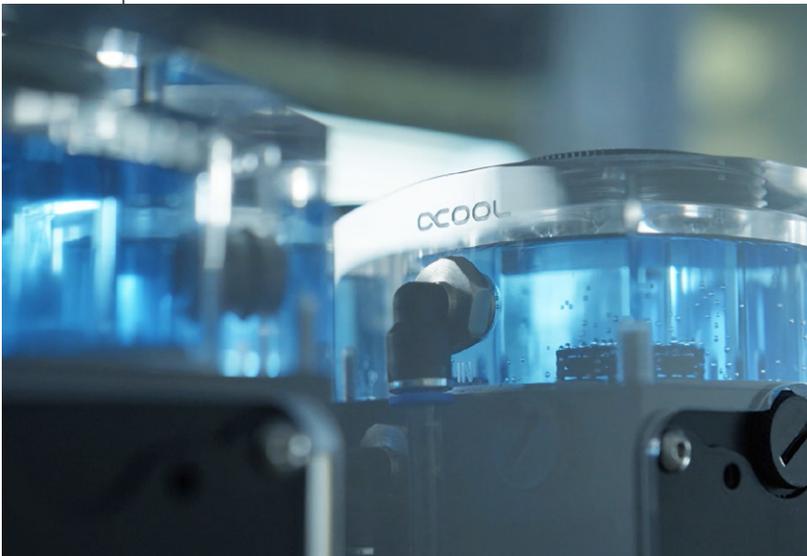


SyncroPatch 384

For all your electrophysiology needs

Sophisticated temperature control

- Full control of temperature at the measurement site – heating and cooling
- Temperature changes during an experiment
- Minimal temperature fluctuations: pre-cooling or pre-heating of solutions and compounds individually at each deck position



Features

- 384 amplifier channels
- Flexible throughput -> 32 to 384 wells in parallel
- Voltage & current clamp
- External & internal solution exchange
- Temperature control
- Single- & multi-hole chips, produced in-house
- Rs compensation
- Automated IV analysis
- Optical stimulation (optional)
- Easy & customizable analysis tools
- User & advanced modes of operation

Applications

- Voltage-, ligand-, mechanically & temperature-activated ion channels
- Action potential pharmacology
- Whole cell and perforated patch
- Cell lines, primary cells and stem cells
- CiPA methods