

TRP'ING ON QPATCH IN MULTI-HOLE MODE



SOPHION BIOSCIENCE A/S
Baltorpvej 154
DK-2750 Ballerup
DENMARK
info@sophion.dk
www.sophion.com

SOPHION BIOSCIENCE, INC.
675 US Highway One
North Brunswick, NJ 08902
USA
Phone: +1 732 745 0221
www.sophion.com

SOPHION JAPAN
1716-6, Shimmachi
Takasaki-shi, Gumma 370-1301
JAPAN
Phone: +81 274 50 8388
www.sophion.com

RASMUS B. JACOBSEN

HERVØR L. OLSEN

SØREN FRIIS

RIKKE SCHRØDER

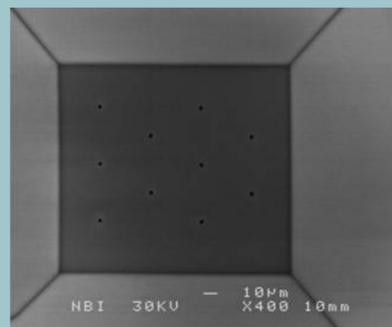
M. K. JENSEN

MORTEN SUNESEN

INTRODUCTION

True giga seal patch clamping can be performed in parallel with QPatch multi-hole technology. This new multi-hole functionality has been tested on three different transient receptor potential ion channels (TRPA1, TRPV1 & TRPM8). All three targets were tested in both single-hole mode and in multi-hole mode with different agonists and antagonists.

The advantages of testing compounds in multi-hole mode will be a minimization of biological variance and an increase in current amplitude, since the current response from several cells are summarized. The disadvantage is lack of serial resistance compensation and in some cases large leak currents.



MATERIALS AND METHODS

Cells: HEK - rTRPV1, HEK - hTRPA1, HEK - hTRPM8

Ringer's solutions:

TRPA1: IC (in mM): KF 120, KCl 20, HEPES 10, EGTA 10, pH 7.2, 300 mOsm EC (in mM): NaCl 155, KCl 5, MgCl₂ 1.6, HEPES 10, BaCl₂ 2, EGTA 5, glucose 10, pH 7.4, 314 Osm

TRPV1: IC (in mM): KF 120, KCl 20, HEPES 10, EGTA 10, pH 7.2, 300 mOsm EC (in mM): NaCl 145, KCl 5, MgCl₂ 1.75, HEPES 10, CaCl₂ 2, EGTA 10, glucose 10, pH 7.4, 310 Osm

TRPM8: IC (in mM): NaCl 4, KCl 130, MgCl₂ 1, HEPES 10, BAPTA (or EGTA) 10, pH 7.4 EC NaCl 150, KCl 4, CaCl₂ 0.5, MgCl₂ 1, HEPES 10, pH 7.4

Compounds:

A1: Antagonist: 4-(4-Chlorophenyl)-3-methylbut-3-en-2-oxime (AP-18), supercinnamaldehyde (CA), Ruthenium Red. All compounds from Sigma. Each compound was dissolved in DMSO. The final DMSO concentration did not exceed 0.1%.

V1: Agonist: Capsaicin (sigma M2028). Antagonist: Capsazepine (Sigma C191) and Tetrabutylammonium chloride (86870 Aldrich)

M8: Agonist: Menthol (Sigma-Aldrich product # M2780. Antagonist: Capsazepine, Cayman Chemical (Cat.no. 10007518).

Electrophysiology: Whole-cell patch-clamp experiments were performed on QPatch. Standard and experimental voltage protocols are presented in the figures.

Data analysis: Recorded ion channel whole-cell currents were stored in an integrated database (Oracle). Drug effects were analyzed as function of time (I-t plot) and concentration (dose-response relationship). Data analysis was accomplished with the QPatch Assay Software.

Acknowledgement: AstraZeneca Södertälje Sweden.

REFERENCE LIST

Andersson D A, Chase H W N, Bevan S, TRPM8 activation by menthol, icilin and cold is differentially modulated by intracellular pH, *J. Neurosci.* 2004; 24(23):5364-5369

Benetikt J, Teisinger J, Vyklicky L, Vlachova V, Ethanol inhibits cold-menthol receptor TRPM8 by modulating its interaction with membrane phosphatidylinositol 4,5-bisphosphate, *J. Neurochem.* 2007; 100:211-224

Chuang H, Neuhauser W M, Julius D, The super-cooling agent icilin reveals a mechanism of coincidence detection by a temperature-sensitive TRP channel, *Neuron* 2004, (43):859-869

Hui K, Gui Y, Feng Z, Biophysical properties of menthol-activated cold receptor TRPM8 channels, *Biochem. Biophys. Res. Comm.* 2005, (333):374-382

McKemy D, Neuhauser W, Julius D, Identification of a cold receptor reveals a general role for TRP channels in thermosensation, *Nature* 2002 (414):52-58

Voets T, Owsianik G, Nilius B, TRPM8, *Handb. Exp. Pharmacol.* 2007; (179):329-44.

Oseguera A J, et al, On the mechanism of TBA block of the TRPV1 channel, *Biophysical Journal*, 2007; 92:3901-3914

Pedersen S F, Owsianik G, Nilius B, TRP channels: An overview, *Cell Calcium*, 2005; 38; 233-252

Clapham D E, TRP channels as cellular sensors, *Nature*; 426: 517-524

SUMMARY

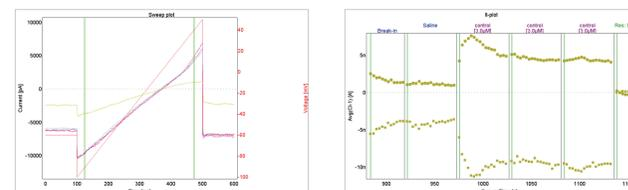
Transient receptor potential channels are activated and regulated by a wide range of stimuli and several different physiological pathways are involved. Automated patch clamping on TRP channels is therefore not trivial. To obtain high quality data assays need to be specially designed to test these targets. For most voltage gated targets an online leak protocol will be sufficient for running experiments in multi-hole mode. However, this is not possible for the TRP channels and other methods need to be applied. Recording of TRP channels in multi-hole mode shows a relative large leak component. By adding a high concentration of reference compound at the end of the experiment this large leak

current can automatically be subtracted. This method requires a specific blocker which blocks the TRP current completely.

By testing the three TRP targets in multi-hole mode the total current level increases by a factor of 10. It is therefore an advantage for targets with low expression on this platform. The agonists and antagonists tested in this study have comparable effects on the three targets using multi-hole and single-hole technology. Setting up the optimal assay the QPatch patch clamp technologies are suitable for getting insight to the behaviour of TRP channels.

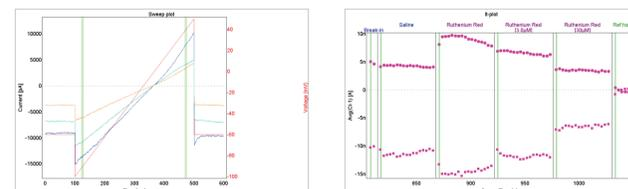
TRPA1 MULTI-HOLE

Supercinnamaldehyde



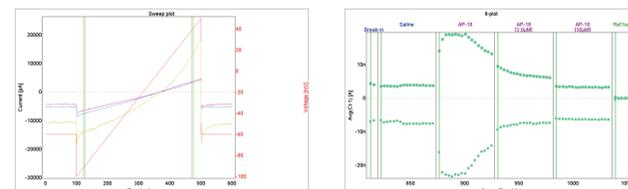
A voltage ramp from -100 mV to +50 mV (duration 400 ms) was executed every three seconds from a holding potential of -60 mV. Supercinnamaldehyde (30 μM) were added through out and the outward current was measured at +40 mV and inward current at -90 or -100 mV. Data were sweep subtracted (the fourth last recorded sweep was subtracted from the others) before further data handling. The data were sampled with a frequency at 2000 Hz and filtered at 800 Hz.

Ruthenium Red



The current was activated by 30 μM of supercinnamaldehyde through out the experiment and a voltage ramp from -100 mV to +50 mV (duration 400 ms) was executed every three seconds from a holding potential of -60 mV and increasing concentrations of Ruthenium Red (3 & 30 μM) were added. Data were sweep subtracted (the fourth last recorded sweep was subtracted from the others) before further data handling. The data were sampled with a frequency at 2000 Hz and filtered at 800 Hz.

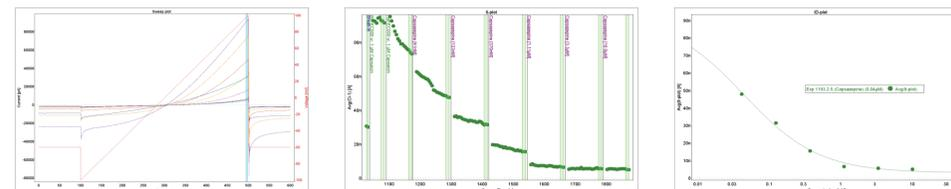
AP-18



The current was activated by 30 μM of supercinnamaldehyde through out the experiment and a voltage ramp from -100 mV to +50 mV (duration 400 ms) was executed every three second from a holding potential of -60 mV and increasing concentrations of AP-18 (3 & 30 μM) were added. Data were sweep subtracted (the fourth last recorded sweep was subtracted from the others) before further data handling. The data were sampled with a frequency at 2000 Hz and filtered at 800 Hz.

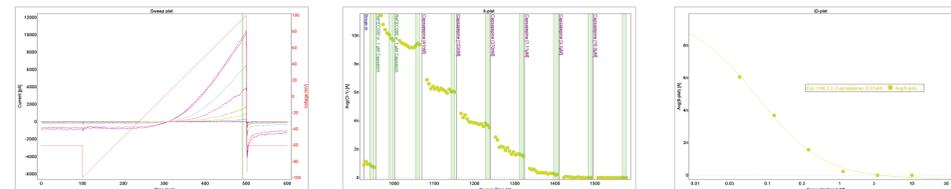
TRPV1 MULTI-HOLE

Capsazepine



TRPV1 SINGLE-HOLE

Capsazepine

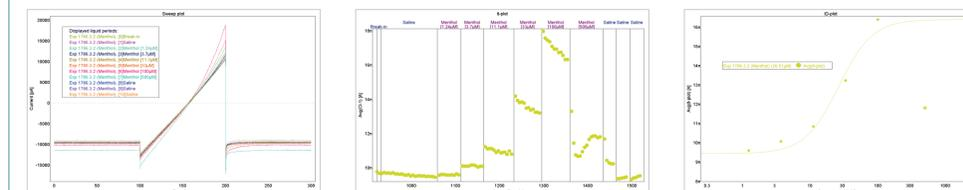


The current was activated with 1 μM capsaicin through out the experiment and a voltage ramp from -100 mV to +100 mV (duration 600 ms) was executed every fifth second from a holding potential of -60 mV. Increasing concentrations of

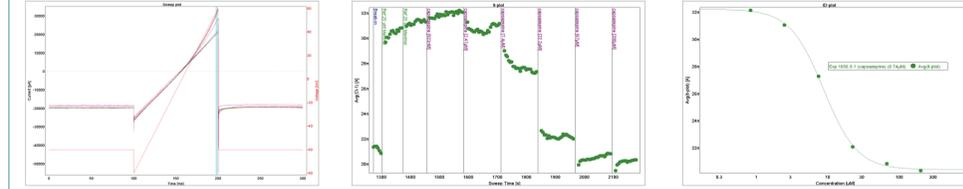
capsazepine were added and the outward current measured at +90. The data was sampled with a frequency at 5 kHz and filtered at 1 kHz.

TRPM8 MULTI-HOLE

Menthol

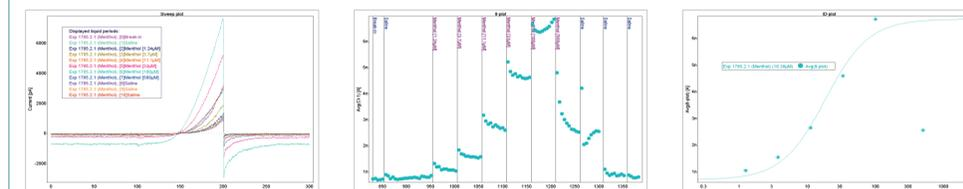


Capsazepine

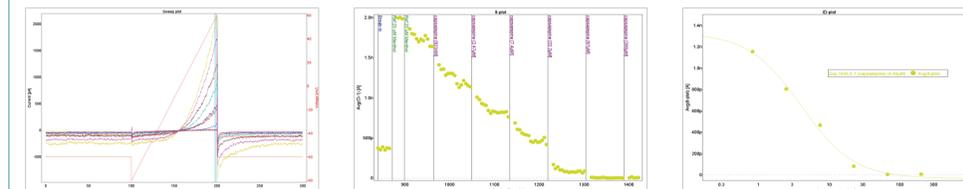


TRPM8 SINGLE-HOLE

Menthol



Capsazepine



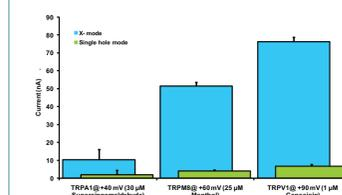
Menthol

A voltage ramp from -80 mV to +60 mV (duration 100 ms) was executed every fifth second from a holding potential of -60 mV. Increasing concentrations of menthol were added and the outward current measured at +60 mV. The data were sampled with a frequency at 10 kHz and filtered at 3 kHz.

Capsazepine

The current was activated with 25 μM Menthol through out the experiment and a voltage ramp from -80 mV to +60 mV (duration 100 ms) was executed every fifth second from a holding potential of -60 mV. Increasing concentrations of capsazepine were added and the outward current measured at +60 mV. The data were sampled with a frequency at 10 kHz and filtered at 3 kHz.

COMPARISON OF SINGLE-HOLE AND MULTI-HOLE RESULTS



Average outward TRP current amplitude in single-hole mode (green) and in multi-hole mode (blue). The current was measured at +40, +60 and +90 mV for TRPA1, TRPM8 and TRPV1, respectively.

		Single-hole mode	Multi-hole mode
TRPA1	Ruthenium Red	n.d.	20.19±5.37 μM (n=7)
	AP-18	n.d.	4.53±1.76 μM (n=10)
TRPV1	Capsaicin	n.d.	7.59±6.15 nM (n=6)
	Capsazepine	30.7±7.77 nM (n=4)	43.3±12.6 nM (n=8)
TRPM8	Menthol	99.35±50.43 nM (n=6)	21.47±1.12 nM (n=10)
	Capsazepine	6.03±1.31 μM (n=3)	15.56±3.04 μM (n=5)

EC₅₀ and IC₅₀ values for different TRP agonists and antagonists were obtained for both multi-hole, and in single-hole mode recordings.