

Application Report

Validation of ligand-gated ion channels in multi-hole mode on QPatch

Comparison between single-hole and multi-hole mode to determine if the signal onset was comparable and if the effects of known compounds (agonists, antagonist and modulators) to see if any changes could be detected

Summary

Investigations of fast ligand-gated applications on QPatch in multi-hole mode. We evaluated rise times of rP2X3, nAChR $\alpha 1$ and $\alpha 7$ and found that they were similar when compared to single-hole recordings. We used known reference compounds to validate the pharmacology and the results for GABA_A, P2X3, nAChR $\alpha 1$ and GluR5 agreed with data obtained in single-hole recordings. We noted that running assays in multi-hole mode increased the success rates and we can conclude that QPatch in multi-hole mode is well suited for running fast ligand-gated ion channels without sacrificing data quality.

Introduction

This report summarizes the experiments that were done running fast ligand-gated assays on QPatch.

QPatch in multi-hole mode patches 10 cells per well instead of just one which has been shown to increase the overall success rate because cells with less or no expression of the ion channel of interest, will be patched in parallel with other cells that has better expression. The currents from all 10 holes are thus added and therefore gives rise to increased current amplitudes from each well.

The 10 holes take up more space in the flow channel in the QPlate chip and we were curious to know if these changes had any effect on the measured signal from the ten cells.

We set out to test different ligand-gated ion channels that are well known and already validated on QPatch to determine if the signal onset was comparable to QPatch in multi-hole mode. Furthermore, we investigated the effect of known compounds: agonists, antagonist and modulators to see if any changes in data quality could be detected.

Results

RBL2H3

This cell line has been widely used to answer different flow and pipetting related questions. The protocol used is shown in figure 1a and a typical response is shown in figure 1b.

RBL2H3 expresses an inward rectifier in large amounts and changing driving force for K will induce a "ligand"-gated response that can be analyzed.

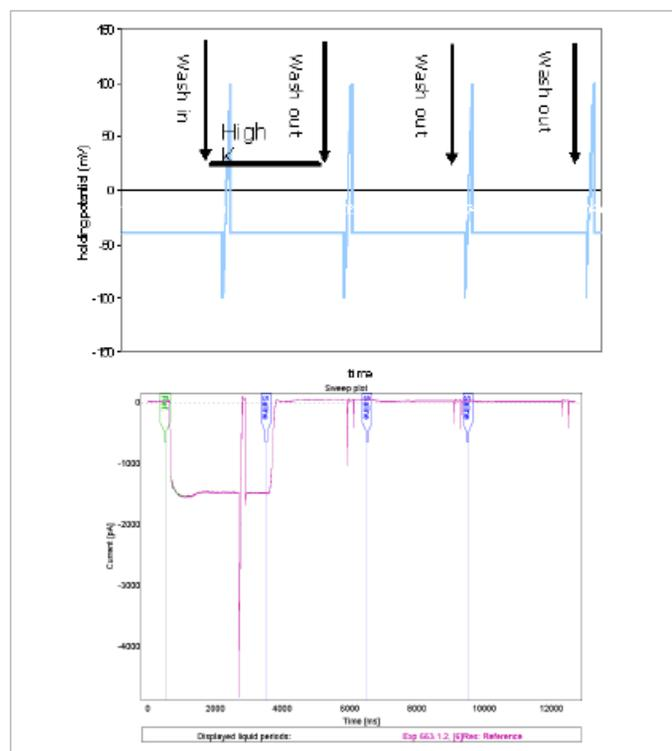


Fig. 1: Top: Voltage protocol including pipetting time stamps. High K is added for 3 sec. and then washed off again. The protocol includes 4 ramps where K concentration in the flow channel can be estimated. Bottom: Typical response from a single cell experiment. The bottom of the ramp is highlighted with black dots. Note that three washes are needed for complete washout.

In this set of experiments the exchange rate was evaluated in two ways. By adding a 10-90% rise time cursor and by fitting the "response" to a single exponential function (Figure 2).

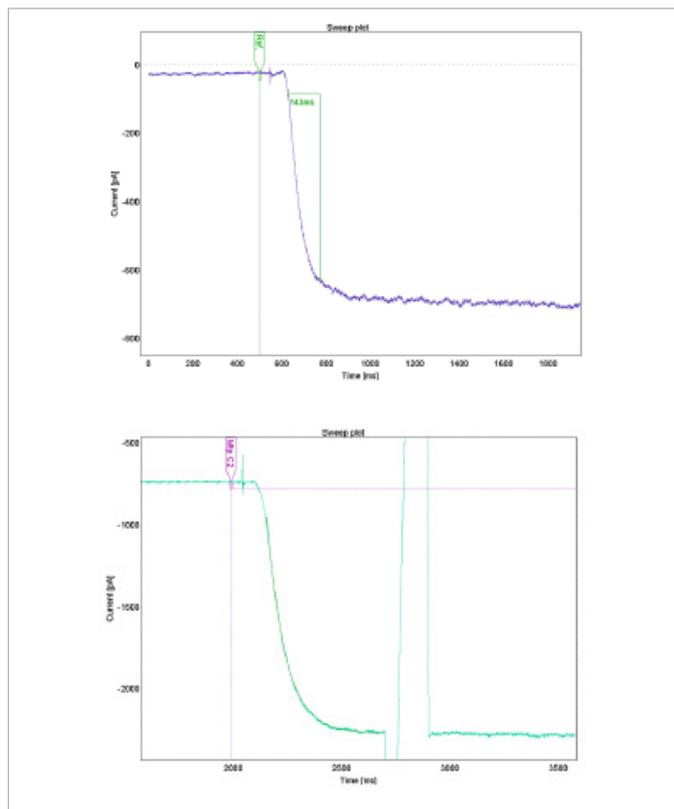


Fig. 2: Top: 10 – 90% rise time cursor on a single-hole experiment. Bottom: RBL response recorded in multi-hole mode and fitted to $A + B\exp(-t/\tau)$.

The data from two inserts collected from single-hole and multi-hole experiments are summarized in Table 1. It was found that the current amplitude was increased from 0.7 to 5.5 nA i.e. approx. 8-fold when comparing single-hole and multi-hole. This indicates that we have close to, but not full whole cell efficiency when running multi-hole.

When comparing the rise time parameters between the two systems it was found that there was no difference in this regard. The rise time was 170 ms for single-hole and 183 ms for multi-hole and tau was estimated to be 89 and 69 respectively.

Table 1: Current response rise time and activation time constant measured on RBL2H3 with single-hole and multi-hole measurement plates.

	Single-hole	Multi-hole
Current response (nA)	0.7 ± 0.3	5.05 ± 2.0
Rise time (10-90%) (ms)	170 ± 49	183 ± 27
"Activation" $1/C$ (ms)	89 ± 18	65 ± 14

TE671 nicotinic acetylcholine receptor $\alpha 1$

We tested rise times on the acetylcholine sensitive nicotinic $\alpha 1$ channel in TE671 cells. This cell line expresses the channel endogenously and is a more relevant assay since $\alpha 1$ is a "real" ligand-target especially used for counter screening.

This could potentially be an assay where a slower rise time could also be detected. This time we looked at the rise time as a function of increasing concentrations of acetylcholine. We ran the same assay on QPatch in multi-hole and single-hole mode and compared the 10-90% rise time in the same manner as in the RBL 2H3. The results are summarized in Figure 3.

The rise time decreases as a function on acetylcholine as expected - from approx. 150 ms to 25 ms on the highest concentrations. It was found that there was no difference between single-hole and multi-hole when comparing rise times in the two modes.

The success rate was found to be good in both systems: 90% and 93% sites with $\alpha 1$ currents were found for single-hole and multi-hole mode respectively. This success rate is counted from a total insert i.e. 16 sites or more also counting sites that do not prime, gets a cell positioned or fails in another way. Hence 90% means that 90% of all sites obtained whole-cell configuration with measurable $\alpha 1$ current in single-hole mode.

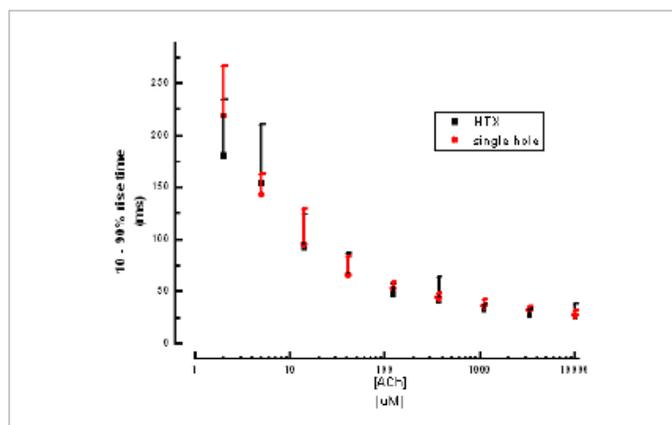


Fig. 3: Average 10-90% rise times \pm SD from at least 16 sites measured as a function of acetylcholine concentration in multi-hole and single-hole mode.

RLE rP2X3

This ATP-gated channel was the third cell line where the rise time was evaluated. This rat cell line has been used to validate ligand-gated assays on QPatch before where it was found to have a success rate of around 25%, not because of patch ability issues, but mainly due to low expression rates in the cell culture. It was therefore interesting not only because it is one of the fastest channels of the P2X family but also because we wanted to test if we could improve the overall success rate with QPatch in multi-hole mode. Figure 4 showing a raw data recording of the rise time as a response to 10 mM ATP and a summarizing figure of the rise time as a function of ATP concentration.

Again, the rise time decreases as a function agonist - from approx. 250 ms to 15 ms on the highest concentrations. It was found that there was no difference between single-hole and multi-hole mode when comparing rise times.

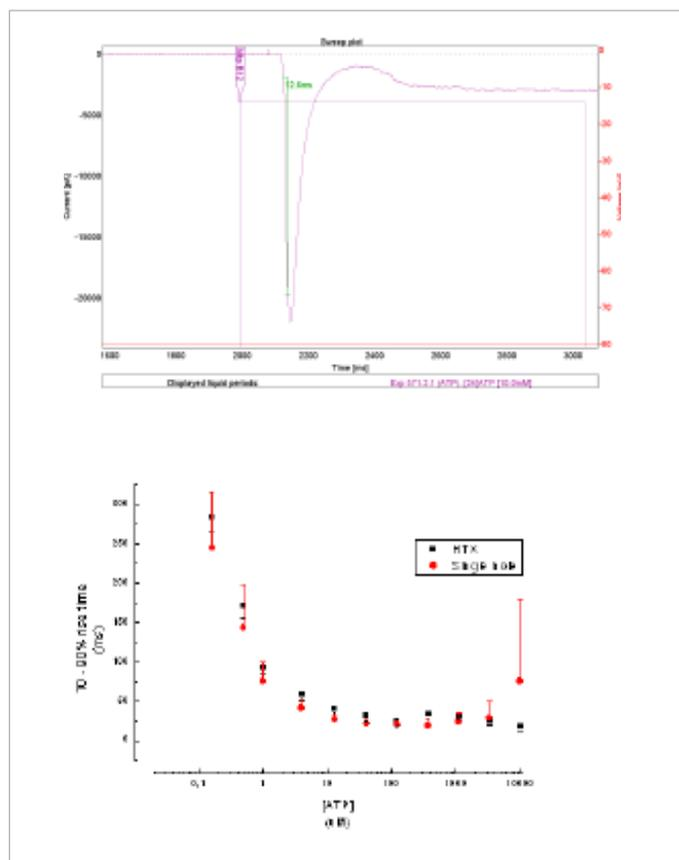


Fig. 4: Top: Leak subtracted 10 mM ATP induced P2X3 response with 10 – 90% rise time cursor. Rise time is 17 ms. Bottom: Average rise time +/- SD vs. ATP concentration.

Surprisingly, we found that the success rate was 81% in single-hole vs. 100% in multi-hole mode. As stated above we had expected the success rate to be close to 25% sites with current. This increase indicates that we have become better at culturing these cells or that this assay has improved over time. In any case we could again see that the success rate went up to 100% when running in multi-hole mode.

GH4C1 nicotinic acetylcholine receptor $\alpha 7$

The nicotinic acetylcholine $\alpha 7$ channel has already been shown to be measurable on QPatch (Friis, Mathes et al. 2009). This target is an interesting test case for QPatch in multi-hole mode for several reasons: First, because this channel also has a very fast onset and desensitization, but mainly because these channels are very hard to express in high enough quantities to get a usable signal (more than 50 pA) in single-hole mode.

The channel was stimulated with 1 and 10 mM acetylcholine and the rise time was compared to single-hole data with the same protocol. Figure 5 a & b below shows raw data sweeps obtained with GH4 C1 stimulated with acetylcholine.

The dataset was acquired twice – first time with cells that had been frozen for about 1½ years. These vials contained cells from the previous optimization project on this cell line. A second data set was done with cells that were not optimized.

The results from the two data sets are summarized in Table 2.

Table 2: Current amplitude and success rate evaluated with two different batches of GH4 C1 cells in single-hole and multi-hole mode.

	Single-hole	Multi-hole	Cell type
Peak current @ 1 mM ACh (pA)	-114 ± 32	-460 ± 328	Optimized
% sites with $\alpha 7$ current	25%	81%	Optimized
Peak current @ 1 mM ACh (pA)	-33 ± 23	-136 ± 69	Non-optimized
% sites with $\alpha 7$ current	6	28	Non-optimized

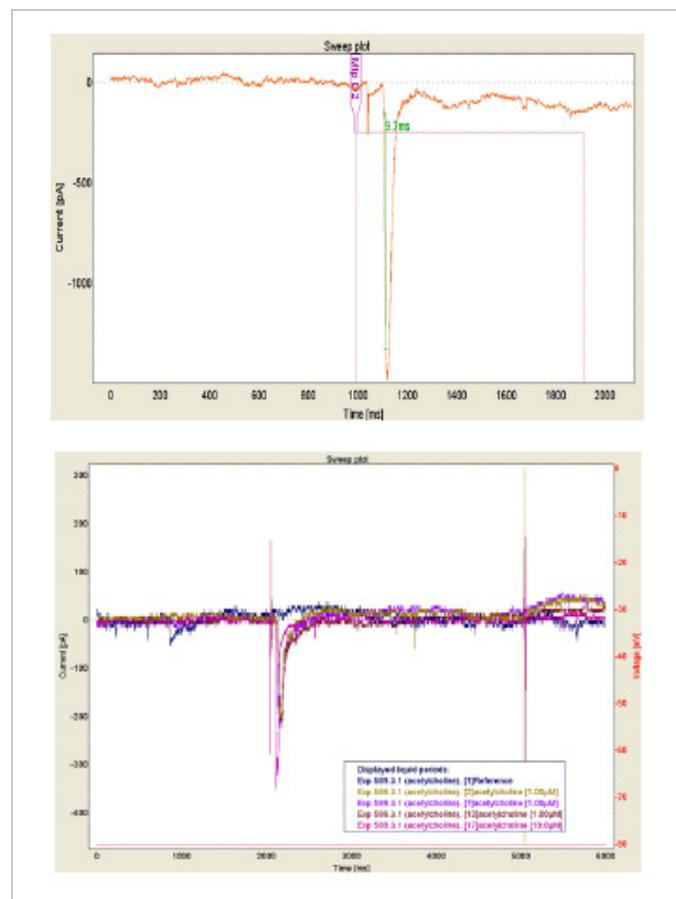


Fig. 5: Top: Leak subtracted non-optimized cells stimulated with 3x1 mM and 1x 10 mM ACh. Bottom: Leak subtracted optimized cells stimulated with 10 mM ACh.

Running the experiments with QPatch technology increases the success rate with both cell types. It was found that the optimized cell line performed better than the non-optimized cell line as expected. The success rate with the optimized cell is increased from 25% to 81%, which for this particular cell line is considered very good. When running non-optimized cells the success rate is also increased but this time from 6% to 28%. The rise time was found to be the same when comparing the two systems (data not shown).

Pharmacological Data

To further evaluate these ligand-gated assays on QPatch in multi-hole mode, we performed several tests with known agonists, antagonists and modulators of these channels. The protocols used for these experiments are all standard protocols for ligand-gated assays. The cycling interval is 150-200 seconds between each round, and this is also the length of preincubation used when compounds have been tested together with an agonist. In some assays, enzymatic wash is used between each cycle – yielding 5 test pt. per well. Assays without enzymatic wash yielded 6 pt. per well.

rP2X3

ATP, $\alpha\beta$ - met-ATP and TNP-ATP were tested on rP2X3. The results can be seen in Figure 6. The number of repeats of each experiment type is in no case less than 4. The success rates in these experiments were found to be equivalent to the initial rise time experiments. Figure 6 shows raw data sweeps obtained with GH4 C1 stimulated with acetylcholine

Table 3: EC₅₀ values for the agonists ATP and ab-met-ATP and IC₅₀ for the antagonist TNP-ATP obtained on P2X3 compared between multi-hole, single-hole and literature values.

XC50 of known compounds on rP2X3 (uM)			
	Single-hole	Multi-hole	Lit. values
ATP	0.7	1.5	1.8 (Wildman, Brown <i>et al.</i> 1999; Wildman, King <i>et al.</i> 1999)
$\alpha\beta$ -met-ATP	1.5	0.3	2.7 (Pratt, Brink <i>et al.</i> 2005)
TNP_ATP @ 1 μ M $\alpha\beta$ - met-ATP	0.021	0.029	0.009 (Virginio, Robertson <i>et al.</i> 1998)

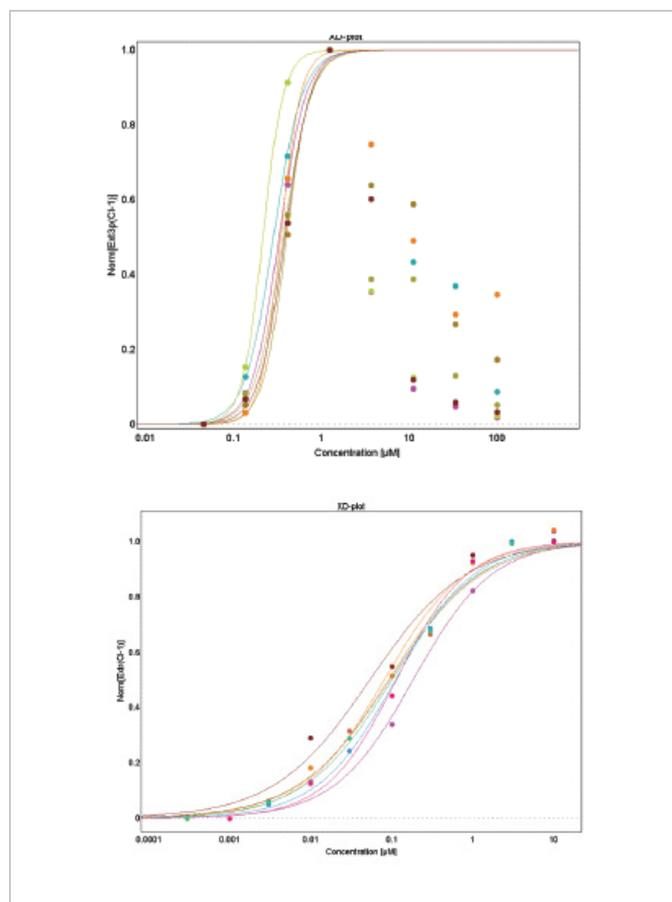


Fig. 6: Top: Normalized $\alpha\beta$ - met-ATP rP2X3 peak responses vs. concentration from 4 different cells. The responses from concentrations eliciting increasing responses are fitted to the Hill equation to estimate EC₅₀. Bottom: Normalized TNP-ATP induced block. Cells stimulated with 1 μ M $\alpha\beta$ - met-ATP. Data combined from 4 different cells.

The data that was collected with multi-hole on QPatch was found to be comparable to reference data from both single-hole on QPatch and literature references.

GABA_A

GABA $\alpha 1\beta 2\gamma 2$ expressed in HEK293 cells was also one of the cell lines tested in this validation study. The reason for this is that this particular cell line has previously been shown to have an effective success rate of around 40% mainly due to the low expression of the ion channel.

GABA is also a widely used ion channel in transient expression assays due to the many possible combinations of subunits. The success rate for transfections is typically also around 40% and hence this would be an assay that would be well-suited for running on QPatch in multi-hole mode.

The results are summarized in Table 4 and Figure 7.

Table 4: EC₅₀ values for GABA, the positive modulator diazepam and IC₅₀ for bicuculline compared between multi-hole, single-hole and literature values.

XC50 of known compounds on GABA (uM)			
	Single-hole	Multi-hole	Lit. values
GABA	5.6	12	9-18 (Curtis, Duggan <i>et al.</i> 1970)
Bicuculline	1.2	1.5	1-3 (Curtis, Duggan <i>et al.</i> 1970)
Diazepam @ 1 μM GABA	ND	0.14	0.15 (Kapur and McDonald 1996)
% sites with current	37	93	

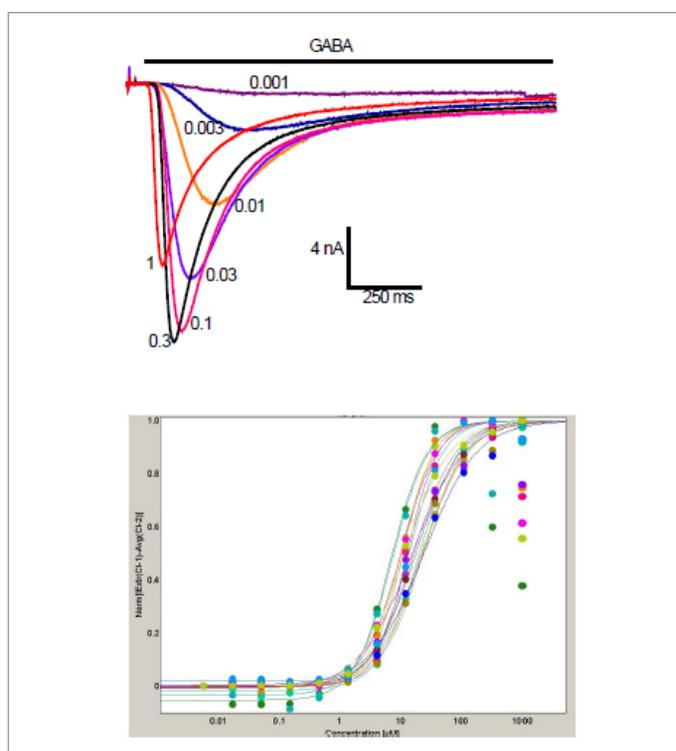


Fig. 7: Top: Leak subtracted GABA α1β2γ2 currents recorded as response to increasing concentrations of GABA (mM). Bottom: Normalized GABA responses from 16 individual sites fitted to the Hill equation. Each cell was exposed to 12 different concentrations of GABA.

The success rate was indeed increased from 37% (as expected) in single-hole mode to 93% sites with GABA current seen in multi-hole mode. The pharmacological data recorded on GABA was found to be very close to what has been seen in single-hole mode and what was found in the literature.

Nicotinic acetylcholine Receptor α1

This channel is affected by local anesthetics and a number of these were tested. The cells were again preincubated with the compound before a dose of acetylcholine (1 mM) was added together with the test concentration of compound. The results are shown in Table 5 and Figure 8.

Table 5: EC₅₀ for agonist acetylcholine and IC₅₀ for four known antagonists and success rates obtained with TE 671 on single hole, multi-hole (HTX) and literature values.

XC50 of known compounds on Nicotinic α1 (uM)			
	Single-hole	Multi-hole	Lit. values
Acetylcholine	5.6	12	8.5 (Shao, Mellor <i>et al.</i> 1998)
Tetracaine	ND	1.6	13 (Gentry and Lukas 2001)
Lidocaine	ND	<100	52 (Gentry and Lukas 2001)
Gallamine	5.3	2.2	0.9 (Paul Kindler <i>et al.</i> 2002; Purohit, Tate <i>et al.</i> 2007)
Rocuronium	ND	0.13	0.33 (Paul Kindler <i>et al.</i> 2002; Purohit, Tate <i>et al.</i> 2007)
% sites with current	90	93	NA

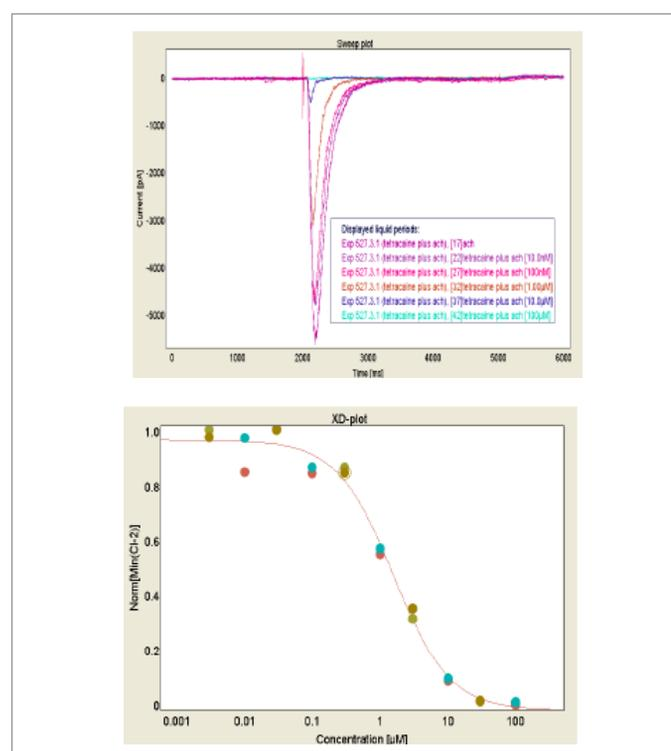


Fig. 8: Top: α1 current stimulated with 1 μM ACh in the presence of increasing concentrations of tetracaine (0.01-100 μM). Bottom: Resulting concentration response plot from 4 sites, IC₅₀ = 1.6 μM for tetracaine

Again, we found excellent agreement between data obtained on multi-hole and single-hole for all these compounds tested. The success rate was not increased tremendously but this is because this cell line performs very well in both single-hole and multi-hole mode – around 90% sites with current in each.

GluR5

This cell line was also tested in multi-hole mode and compared to single-hole data obtained previously. The channel is expressed in HEK 293 cells. The results of the tests with its natural agonist glutamate, the analogue kainate and an antagonist are shown in Table 6 and Figure 9.

Table 6: EC₅₀ values for agonists and IC₅₀ values for antagonists tested on GluR5 and compared between multi-hole, single-hole and literature values.

XC50 of known compounds on GluR5 (uM)			
	Single-hole	Multi-hole	Lit. values
Glutamate	226	344	630 (Lerma, Paternain <i>et al.</i> 2001)
Kainate	119	299	33-177 (Lerma, Paternain <i>et al.</i> 2001)
CNQX	1.9	0.8	0
% sites with current	65	91	

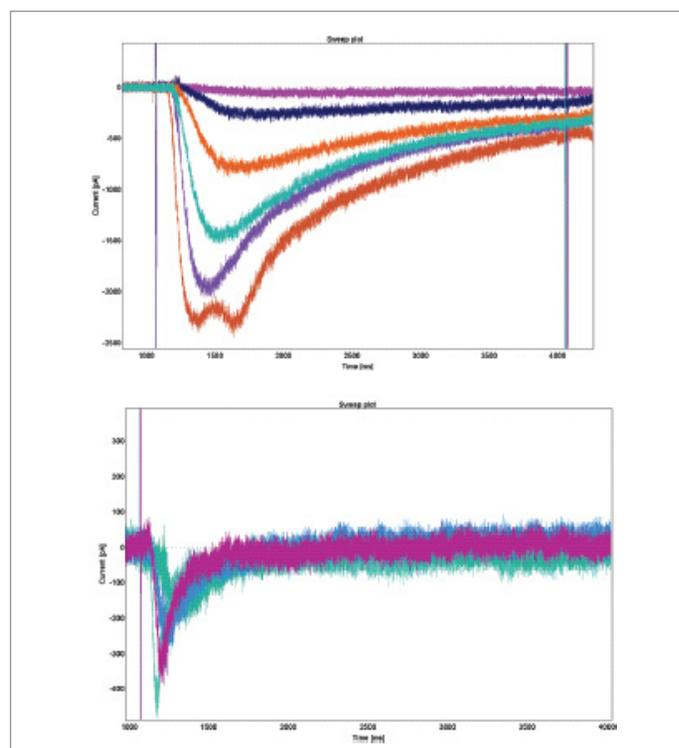


Fig. 9: Top: 1: Leak subtracted GluR5 currents recorded as response to increasing concentrations of kainate. Bottom: Leak subtracted GluR5 currents recorded as response to increasing concentrations of glutamate.

With this cell line we again saw a marked increase in success rate when evaluating sites with kainate induced currents. The current response was found to be markedly increased when using kainate as the agonist compared to using glutamate which has also been seen on single-hole mode on QPatch. The response was found to be 10-fold increased when running multi-hole.

With this assay QPatch multi-hole mode could facilitate the use of glutamate which is the natural ligand for this channel since the currents increased to a level where it is usable. However, it is still recommended to use kainate since the signal is larger and more stable with this ligand.

The results XC50's found on this channel with multi-hole are comparable to the data found with single-hole and when compared to literature values.

Conclusion

This study investigates the use of fast ligand-gated applications on QPatch in multi-hole mode in order to evaluate these types of demanding channels with this technology.

We evaluated the rise times of rP2X3, nAChR α 1 and α 7 and found that these characteristics were the same when compared to single-hole recordings. In addition, we also tested several known reference compounds to validate the pharmacology on these channels. The results found on GABA_A, P2X3, nAChR α 1 and GluR5 were all found to be in very good agreement with data obtained in single-hole recordings.

It was seen that running these assays in multi-hole mode increased the success rates in all cases. In most assays the success rate was at least 90%. For non-optimized nAChR α 7 it was found to be 28%, but still increased from 6%.

We therefore conclude that QPatch in multi-hole mode is indeed very well suited for running fast ligand-gated ion channels – it increases the success rate without sacrificing any of the data quality found with single-hole recordings.

Methods

Cells

The cell lines used were RBL 2H3 (endogenous Kir); TE671 (endogenous nicotinic α 1); HEK 293 GABA_A; RLE P2X₃; HEK GluR5; GH4 C1 nicotinic α 7.

References:

- Curtis DR, Duggan AW et al. 1970. Bicuculline and central GABA receptors. *Nature* 228(5272):676-7
- Friis S, Mathes C et al. 2009. Characterization of compounds on nicotinic acetylcholine receptor α 7 channels using higher throughput electrophysiology. *J Neurosci Methods* 177(1):142-8
- Gentry CL and Lukas RJ. 2001. Local anesthetics noncompetitively inhibit function of four distinct nicotinic acetylcholine receptor subtypes. *J Pharmacol Exp Ther* 299(3):1038-48
- Kapur J and Macdonald RL. 1996. Pharmacological properties of gamma-aminobutyric acidA receptors from acutely dissociated rat dentate granule cells. *Mol Pharmacol* 50(3):458-66
- Jerina J, Paternain AV et al. 2001. Molecular physiology of kainate receptors. *Physiol Rev* 81(3):971-98
- Paul M, Kindler CH et al. 2002. The potency of new muscle relaxants on recombinant muscle-type acetylcholine receptors. *Anesth Analg* 94(3): 597-603; table of contents
- Pratt EB, Brink TS et al. 2005. Use-dependent inhibition of P2X₃ receptors by nanomolar agonist. *J Neurosci* 25(32):7359-65
- Purohit PG, Tate RJ et al. 2007. The role of the amino acid residue at α 1:189 in the binding of neuromuscular blocking agents to mouse and human muscle nicotinic acetylcholine receptors. *Br J Pharmacol* 150(7):920-31
- Shao Z, Mellor IR et al. 1998. Potentiation and inhibition of nicotinic acetylcholine receptors by spermine in the TE671 human muscle cell line. *J Pharmacol Exp Ther* 286(3):1269-76
- Virginio C, Robertson G et al. 1998. Trinitrophenyl-substituted nucleotides are potent antagonists selective for P2X₁, P2X₃, and heteromeric P2X_{2/3} receptors. *Mol Pharmacol* 53(6):969-73
- Wildman SS, Brown SG et al. 1999. Selectivity of diadenosine polyphosphates for rat P2X receptor subunits. *Eur J Pharmacol* 367(1):119-23
- Wildman SS, King BF et al. 1999. Modulatory activity of extracellular H⁺ and Zn²⁺ on ATP-responses at rP2X₁ and rP2X₃ receptors. *Br J Pharmacol* 128(2):486-92